

## Contents by Research Area:

- Hypertension
- Thrombosis and Hemostasis
- Atherosclerosis
- Myocardial Infarction
- Ischemia/Reperfusion Injury
- Arrhythmias
- Heart Failure



Wild Garlic  
*Allium ursinum*  
A source of Allicin

# Cardiovascular Research

## Contents

	<i>Page</i>
Hypertension	3
Thrombosis and Hemostasis	6
Atherosclerosis	8
Myocardial Infarction	10
Ischemia/Reperfusion Injury	12
Arrhythmias	14
Heart Failure	16
List of Acronyms	18
Related Literature	19
Cardiovascular Research Products	20
Further Reading	31

## Introduction

Cardiovascular disease is the leading cause of death globally, accounting for about 30% of deaths worldwide. The term refers to a collection of diseases which affect the heart and blood vessels. These conditions can be genetic in nature or triggered by infection, but the major influences on the development of cardiovascular disease are environmental factors including a high saturated fat intake and a sedentary lifestyle.

One of the most common cardiovascular diseases is atherosclerosis – a condition which is characterized by the build-up of fatty deposits within arterial walls and which can trigger heart attack or stroke through the formation of a blood clot. Damage to the heart muscle caused by a heart attack can alter the contractility of the heart, leading to irregular heartbeats (arrhythmias). A reduced ability of the heart to pump blood around the body can eventually lead to the development of heart failure. Due to the close association of cardiovascular disorders, early treatment of these disorders is preferable.

Surgical interventions such as coronary artery bypass and artificial pacemakers are available to treat patients with cardiovascular disease, though the treatments themselves may worsen the overall function of the heart. For many patients, pharmacological intervention is therefore preferable to invasive surgery, particularly for the treatment of early atherosclerosis and high blood pressure. For example, a major breakthrough in the treatment of cardiovascular disease was the development of statins for the treatment of raised cholesterol. However, therapy for other diseases such as heart failure and advanced atherosclerosis is lacking and demands further research.

Our latest cardiovascular research guide aims to discuss current and future therapeutic targets in seven major cardiovascular diseases, all of which are the focus of significant preclinical research. Tocris provides a wide range of pharmacological tools for these targets, a selection of which are highlighted in our 'Key Products' boxes within each section. A full product listing can be found on page 20.

## Key Cardiovascular Research Products

Box Number	Title	Page
Box 1	Vasoconstrictor Key Products	5
Box 2	Vasodilator Key Products	5
Box 3	Thrombosis and Hemostasis Key Products	7
Box 4	Atherosclerosis Key Products	9

Box Number	Title	Page
Box 5	Myocardial Infarction Key Products	11
Box 6	Ischemia/Reperfusion Injury Key Products	13
Box 7	Arrhythmia Key Products	15
Box 8	Heart Failure Key Products	17

# Hypertension

Products by Category	Page
$\alpha_1$ and $\alpha_2$ Adrenergic Receptors	20
$\beta$ Adrenergic Receptors	20
Aldosterone Receptors	21
Angiotensin-converting Enzyme	21
Angiotensin II Receptors	21
Calcium Channels	22
Cyclic GMP	23
Endothelin Receptors	24
Muscarinic Receptors (mAChRs)	25
Natriuretic Peptide Receptors	25
Nitric Oxide	26
NKCC Cotransporter	26
Phosphodiesterases	26
Potassium Channels	27
Prostanoid Receptors	28
Rho-kinase	29
Soluble Guanylyl Cyclase	29

## Hypertension

Hypertension is defined as a chronic elevation in blood pressure with a systolic pressure over 140 mmHg and a diastolic pressure over 90 mmHg. The majority of hypertension is primary – that is, an increase in blood pressure with no underlying cause – yet pathologies that affect the kidney or endocrine system may also trigger hypertension. This is known as secondary hypertension. The exact mechanism of primary hypertension is yet to be elucidated, though dysfunctions in mechanisms that regulate vascular tone, both directly and indirectly, have been identified as having a major influence on hypertension.

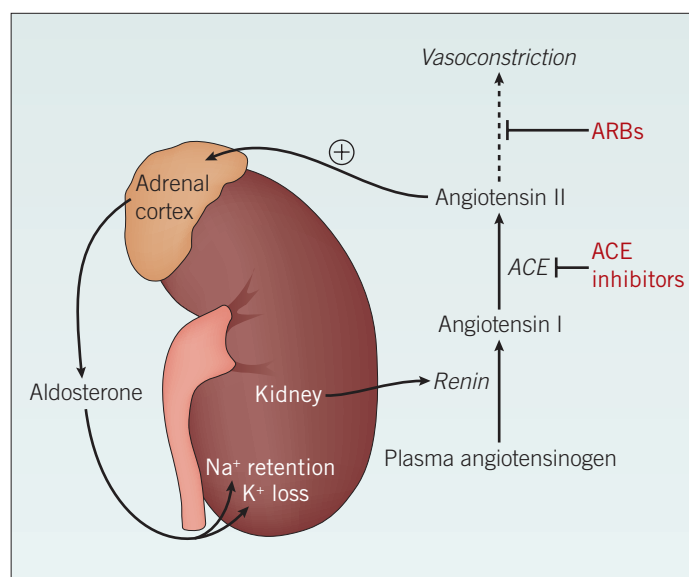
In hypertension, increased arterial pressure is detected by specialized mechanoreceptors called baroreceptors, present in the aortic arch and the carotid sinuses. Baroreceptors are innervated by nerves that synapse in the nucleus tractus solitarius (NTS), an area within the medulla oblongata that regulates blood pressure through the modulation of parasympathetic and sympathetic transmission. In the event of a rise in blood pressure, the baroreceptor firing rate increases; this stimulates the activation of sympathetic neurons that originate in the NTS and synapse in the outer arterial wall, or adventitia. Activation of these sympathetic neurons induces vasoconstriction through the release of noradrenaline and subsequent activation of  $G_q$  and the downstream  $IP_3$  signal transduction pathway. As a result, drugs that target  $\alpha$  adrenergic receptors modulate blood pressure. The precise effect on vascular tone is dependent on the  $\alpha$  adrenergic receptor subtype;  $\alpha_1$  adrenergic receptors stimulate the release of noradrenaline from sympathetic nerve terminals, whilst  $\alpha_2$  adrenergic receptors inhibit the release of noradrenaline, acting as a feedback mechanism to modulate its release from sympathetic nerve terminals.

In addition to sympathetic mechanisms, targeting the renin-angiotensin-aldosterone system (RAAS) is a proven and effective strategy in hypertension. The activation of the RAAS in response to a fall in blood pressure leads to the release of renin from the juxtaglomerular apparatus in the kidney (Figure 1). Renin cleaves angiotensinogen, which undergoes further cleavage to produce the highly potent vasoconstrictor, angiotensin II. Angiotensin II binding to the membrane-bound GPCR, angiotensin II receptor 1 ( $AT_1$ ), induces vasoconstriction directly through the potentiation of noradrenaline release from sympathetic nerve terminals within blood vessel walls.

The downstream effects of  $AT_1$  receptor activation are counterbalanced in part by the activation of  $AT_2$  receptors. Due to the potent vasoconstrictor properties of  $AT_1$  receptor activation, drugs which act as antagonists at this receptor, such as valsartan (Cat. No. 4216) and losartan (Cat. No. 3798), are effective antihypertensives through their indirect vasodilator activity. Inhibition of angiotensin-converting enzyme (ACE) blocks production of angiotensin II and therefore exhibits antihypertensive effects.

Angiotensin II also acts indirectly by stimulating the secretion of vasopressin (AVP) from the pituitary gland and increasing the release of aldosterone from the adrenal cortex. Both AVP and aldosterone augment water reabsorption in the kidney, thereby increasing blood volume and therefore blood pressure.

**Figure 1 | The renin-angiotensin-aldosterone system**



Activation of the RAAS in response to a fall in blood pressure stimulates the release of renin from the kidney. This leads to the production of the potent vasoconstrictor, angiotensin II. Angiotensin II also induces aldosterone release from the adrenal cortex, triggering increased water reabsorption. Together these mechanisms counter the decrease in blood pressure. Abbreviations: ACE – angiotensin-converting enzyme; ARB – angiotensin II receptor blocker

Aldosterone receptor antagonists including spironolactone (Cat. No. 2968) and eplerenone (Cat. No. 2397) also exert antihypertensive effects due to their inhibitory actions on water reabsorption. The resultant reduction in circulating blood volume in turn lowers blood pressure. Other diuretics such as furosemide (Cat. No. 3109) and bumetanide (Cat. No. 3108) also lower circulating blood volume and therefore blood pressure, though they act through the inhibition of the  $\text{Na}^+/\text{K}^+/\text{2Cl}^-$  cotransporter (NKCC). The NKCC is an integral membrane pump which drives calcium ion and magnesium ion reabsorption in the renal medulla, resulting in water reabsorption from the loop of Henlé.

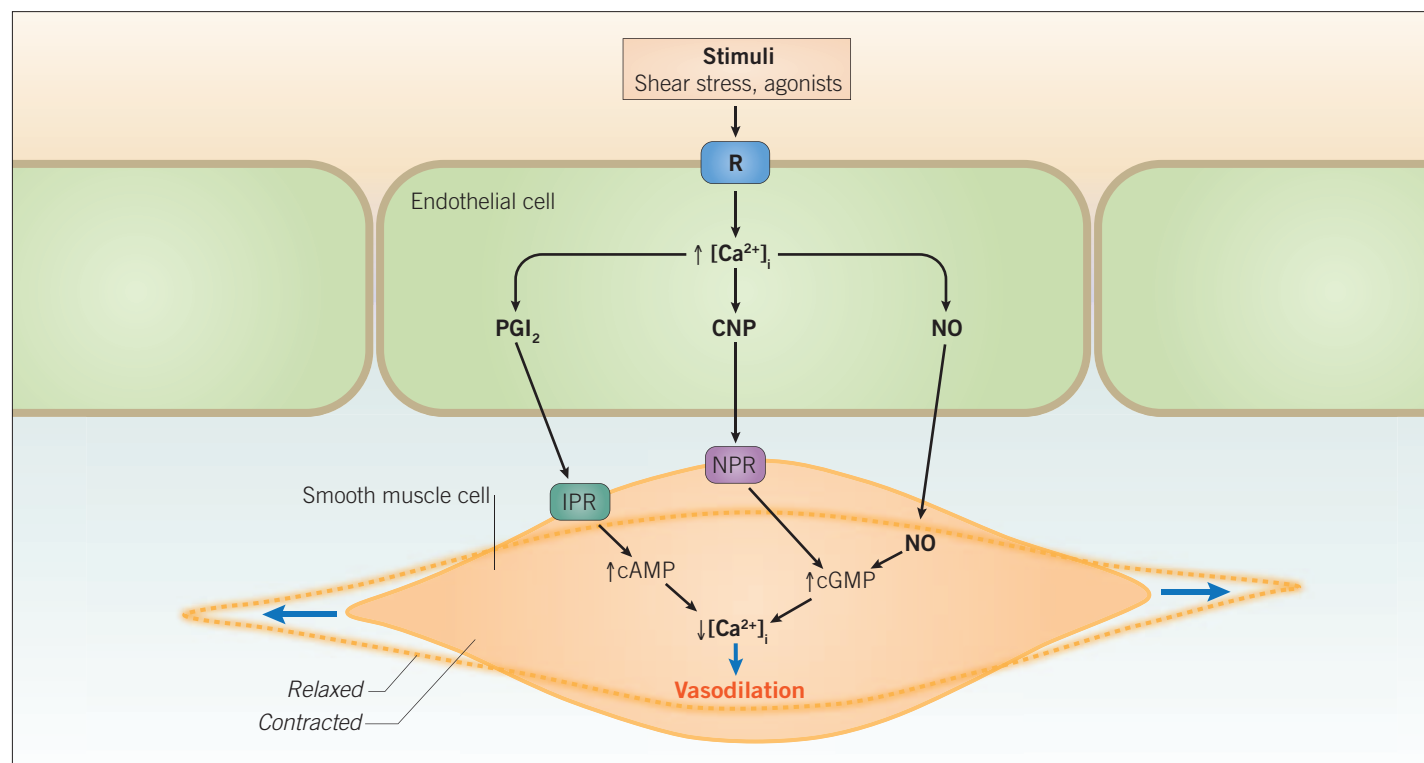
In addition to indirect control of vascular tone by the sympathetic nervous system and RAAS, direct control mechanisms within the blood vessel wall are also a valid therapeutic target in hypertension. Key regulators of blood pressure within the vasculature include nitric oxide (NO), endothelin 1 (ET-1) and prostacyclin ( $\text{PGI}_2$ ) (Figure 2). Other major vasodilators including acetylcholine and bradykinin also directly alter vascular tone by inducing the production of endothelial nitric oxide.

Endothelin receptors can be divided into two classes – endothelin A ( $\text{ET}_A$ ) receptors and endothelin B ( $\text{ET}_B$ ) receptors.  $\text{ET}_B$  receptors are highly expressed in the endothelium whereas  $\text{ET}_A$

receptors are absent, yet both receptor subtypes are present on the underlying vascular smooth muscle cells. Activation of  $\text{ET}_A$  receptors by ET-1 leads to vasoconstriction whilst the effects of  $\text{ET}_B$  receptor activation are cell type-specific; endothelial cell  $\text{ET}_B$  receptor ( $\text{ET}_{B1}$ ) activation leads to vasodilation through the production of NO and  $\text{PGI}_2$ , yet smooth muscle cell  $\text{ET}_B$  receptor ( $\text{ET}_{B2}$ ) activation causes vasoconstriction.

A further therapeutic target in hypertension is NO since its bioavailability is often impaired in hypertensive patients; this is a hallmark of endothelial dysfunction. NO is a key endogenous vasodilator that is secreted in response to endothelial membrane receptor stimulation by agonists such as acetylcholine, bradykinin and 5-HT, as well as shear stress. Activation of endothelial cell membrane receptors by agonist stimulation or shear stress results in an increase in intracellular calcium ion concentration. This increased calcium ion availability activates calmodulin (CaM), a calcium binding protein. The  $\text{Ca}^{2+}$ -calmodulin complex is vital in removing the caveolin-mediated inhibition of endothelial nitric oxide synthase (eNOS), enabling eNOS enzyme activity. The principal reaction of eNOS is to convert L-arginine to L-citrulline, generating nitric oxide as a by-product. NO production and release from endothelial cells triggers an increase in cyclic GMP

**Figure 2 | Nitric oxide-mediated vasodilation**



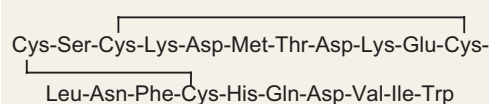
Endothelial cell surface receptor activation by vasodilatory stimuli triggers a rise in intracellular  $\text{Ca}^{2+}$  ( $[\text{Ca}^{2+}]_i$ ). Downstream intracellular signaling pathways lead to the generation of vasodilatory mediators including  $\text{PGI}_2$ , CNP and NO. By binding to their respective receptors on the underlying vascular smooth muscle cell membranes, these mediators prompt a decrease in intracellular calcium within vascular smooth muscle cells through the actions of second messengers including cAMP and cGMP. Decreased intracellular calcium triggers vascular smooth muscle cell relaxation, and subsequent vasodilation. Abbreviations: cAMP – cyclic adenosine monophosphate; cGMP – cyclic guanosine monophosphate; CNP – C natriuretic peptide; IPR – I prostanoid receptor; NO – nitric oxide; NPR – natriuretic peptide receptor;  $\text{PGI}_2$  – prostacyclin; R – receptor.



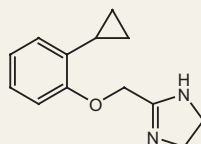
## Hypertension – continued

**Box 1: Vasoconstrictor Key Products**

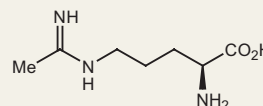
A full list of targets and related products are listed on pages 20-30



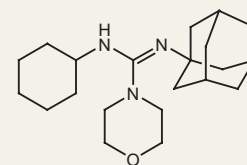
**Sarafotoxin S6a (1899)**  
Potent endothelin receptor agonist;  
vasoconstrictor



**Cirazoline (0888)**  
Selective  $\alpha_1$  agonist



**L-NIO (0546)**  
Potent eNOS inhibitor



**PNU 37883 (2095)**  
Vascular  $K_{ir6}$  ( $K_{ATP}$ ) channel blocker

concentration in the underlying smooth muscle cells through the activation of soluble guanylyl cyclase (sGC), which in turn lowers the intracellular calcium ion concentration, prompting smooth muscle cell relaxation and resulting in vasodilation.

Due to the integral involvement of calcium ions in the control of vascular tone, drugs which prevent the influx of calcium ions through calcium channels following cellular depolarization also directly induce vasodilation. Commonly used calcium channel blockers for the treatment of hypertension include nifedipine (Cat. No. 1075) and diltiazem (Cat. No. 0685). Inhibiting the sensitivity of downstream mediators involved in vasodilation, such as the Rho-associated protein kinase p160ROCK, to calcium ions also prompts vasodilation. An example of a compound that produces vasodilation through this alternate mechanism is Y-27632 (Cat. No. 1254).

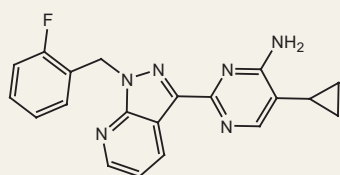
Other vasodilators secreted by endothelial cells that reduce intracellular calcium within underlying smooth muscle cells include prostacyclin, which acts on I prostanoïd (IP) receptors, and C-natriuretic peptide (CNP) through its actions on natriuretic peptide receptors (NPR). Compounds that target these receptors, directly modulate the vasodilatory properties of these endogenous mediators. The activity of these endogenous vasodilators can also be altered by targeting second messenger signaling molecules, namely cAMP and cGMP. Examples of compounds that increase cytoplasmic cAMP levels include the  $\beta$  adrenergic receptor agonist formoterol (Cat. No. 1448), whilst cGMP levels can be therapeutically increased by the activation of sGC, using compounds such as BAY 41-2272 (Cat. No. 4430) and A 350619 (Cat. No. 2753).

Vasodilation is also achieved by selectively increasing the membrane permeability to  $K^+$  ions, leading to cellular hyperpolarization and subsequent relaxation. Activators of ATP-sensitive potassium channels ( $K_{ir6}$ ;  $K_{ATP}$ ), such as cromakalim (Cat. No. 1377) and nicorandil (Cat. No. 2147), possess potent vasodilator activity as a result of this mechanism, though the NO donor activity of nicorandil in addition to its  $K_{ir6}$  channel activating properties further increases its potency as a vasodilator.

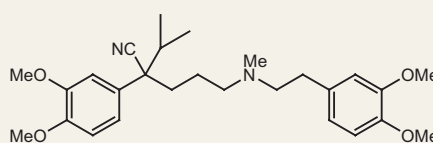
Despite the considerable advances in identifying relevant therapeutic targets involved in hypertension, the side effect profiles resulting from the ubiquitous expression of many of these targets limits their long term use. Rather than directly lowering circulating blood pressure, future perspectives for the treatment of hypertension tend to focus upon ameliorating endothelial dysfunction – through mechanisms such as increasing nitric oxide bioavailability – in order to reduce the target organ damage associated with hypertension and improve prognosis. Since endothelial dysfunction can be triggered by free radical-mediated damage, several clinical trials studied the effects of vitamin E administration in hypertensive patients. These studies reported little or no beneficial effects of the antioxidant, yet there is emerging data of an antihypertensive effect of statins, independent of their lipid-lowering activity. Further potential therapeutic targets in hypertension include increasing L-arginine levels, and preventing the neutral endopeptidase-mediated proteolysis of natriuretic peptides, though additional research focus is required in order to develop novel, effective treatments for hypertension.

**Box 2: Vasodilator Key Products**

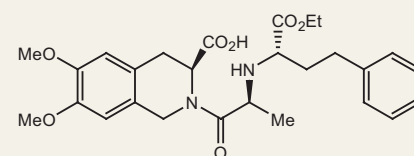
A full list of targets and related products are listed on pages 20-30



**BAY 41-2272 (4430)**  
Soluble guanylyl cyclase (sGC) activator



**Verapamil (0654)**  
 $Ca^{2+}$  channel blocker (L-type)



**Moexepiril (2691)**  
Angiotensin-converting enzyme (ACE) inhibitor

# Thrombosis and Hemostasis

Products by Category	Page
5-HT Receptors	20
Cell Adhesion Molecules	23
Cyclooxygenases	24
IGF-1 Receptor	24
Lipoxygenases	24
PDGF Receptors	26
Phosphodiesterases	26
Phospholipases	27
PI 3-Kinase	27
Prostanoid Receptors	28
Protease-activated Receptors	28
Purinergic P2 Receptors	28
TGF- $\beta$ Receptors	30

## Thrombosis and Hemostasis

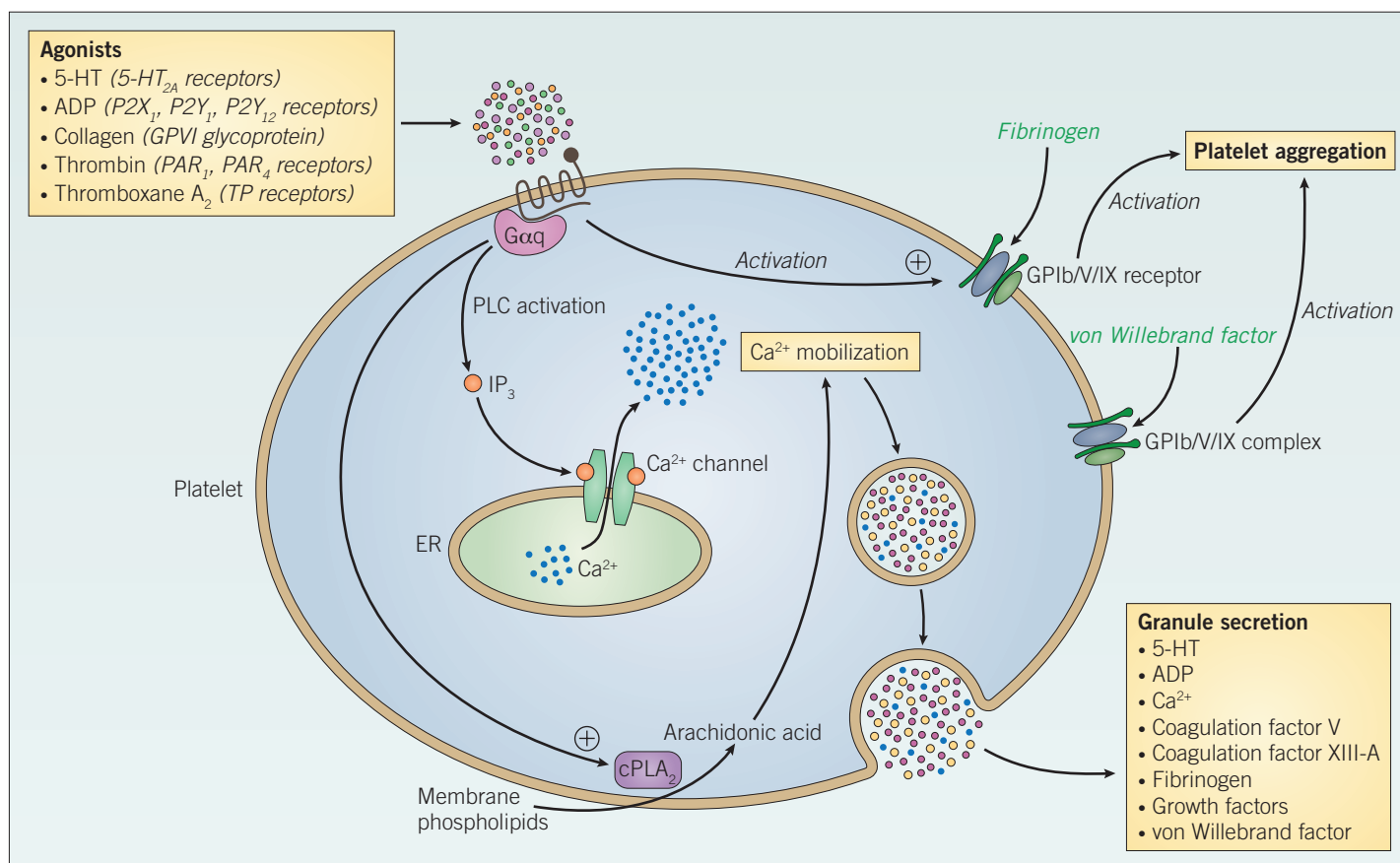
Thrombosis is a crucial hemostatic process for preventing excessive blood loss following injury, yet aberrant thrombosis can trigger pathological conditions including myocardial

infarction and stroke. Therefore the initiation of thrombosis is tightly controlled under physiological conditions.

Platelets are a central component of thrombosis and exhibit a rapid, exponential activation in the event of tissue damage. Produced in the bone marrow, platelets are anucleate cell fragments of megakaryocytes. Despite having no nucleus, platelets possess two different types of granules within the cytoplasm – alpha granules and dense granules – and also express a number of different receptors on their plasma membranes (Figure 3). Both alpha and dense granules contain a variety of bioactive mediators including ADP, calcium and 5-HT as well as growth factors such as platelet-derived growth factor (PDGF), insulin-like growth factor-1 (IGF-1) and transforming growth factor (TGF)  $\beta$ 1. Upon activation, platelets undergo degranulation; this releases granule contents into the surrounding environment and promotes the activation and aggregation of neighboring platelets.

Receptors present on the platelet plasma membrane include the purinergic receptors P2X<sub>1</sub>, P2Y<sub>1</sub> and P2Y<sub>12</sub>; the 5-HT receptor 5-HT<sub>2A</sub>; the thromboxane A<sub>2</sub> (TXA<sub>2</sub>) receptor TP; and the thrombin receptors (protease-activated receptors) PAR<sub>1</sub> and

**Figure 3 | Platelet signaling and activation**



Agonist stimulation of platelets triggers activation of glycoprotein Ib/V/IX; generation of arachidonic acid from membrane phospholipids by the actions of cytosolic phospholipase A<sub>2</sub> (cPLA<sub>2</sub>); and mobilization of intracellular calcium. This in turn triggers the release of granules containing platelet-activating mediators that subsequently act on neighboring platelets, amplifying platelet activation and aggregation.

## Thrombosis and Hemostasis – continued

PAR<sub>4</sub>. Platelet signaling may also be activated by exposure to collagen via the glycoprotein receptor, GPVI. Key downstream mediators of these receptors include the plasma membrane phospholipid phosphatidylinositol 4,5-bisphosphate (PIP<sub>2</sub>), which is cleaved by phospholipase C to form inositol triphosphate (IP<sub>3</sub>) and diacylglycerol (DAG). IP<sub>3</sub> receptor (InsP3R) activation triggers the efflux of calcium ions from intracellular calcium stores such as the endoplasmic reticulum, leading to a rise in intracellular calcium (Figure 3).

The mobilization of calcium causes the exocytosis of cytoplasmic granules and the activation of platelet aggregation-inducing glycoproteins such as GPIIb/IIIa. GPIIb/IIIa functions as a cell surface receptor for fibrinogen and promotes the activation of platelets. The pivotal involvement of GPIIb/IIIa in platelet aggregation renders it a useful therapeutic target for anti-coagulant therapy. GPIIb/IIIa inhibitors, such as abciximab and echistatin (Cat. No. 3202), are currently used as prophylactic therapy during angioplasty to prevent thrombus formation. Inhibitors of P2Y<sub>12</sub> – for example, clopidogrel (Cat. No. 2490) and ticlopidine (Cat. No. 3931) – also indirectly prevent platelet aggregation by blocking the activation of GPIIb/IIIa.

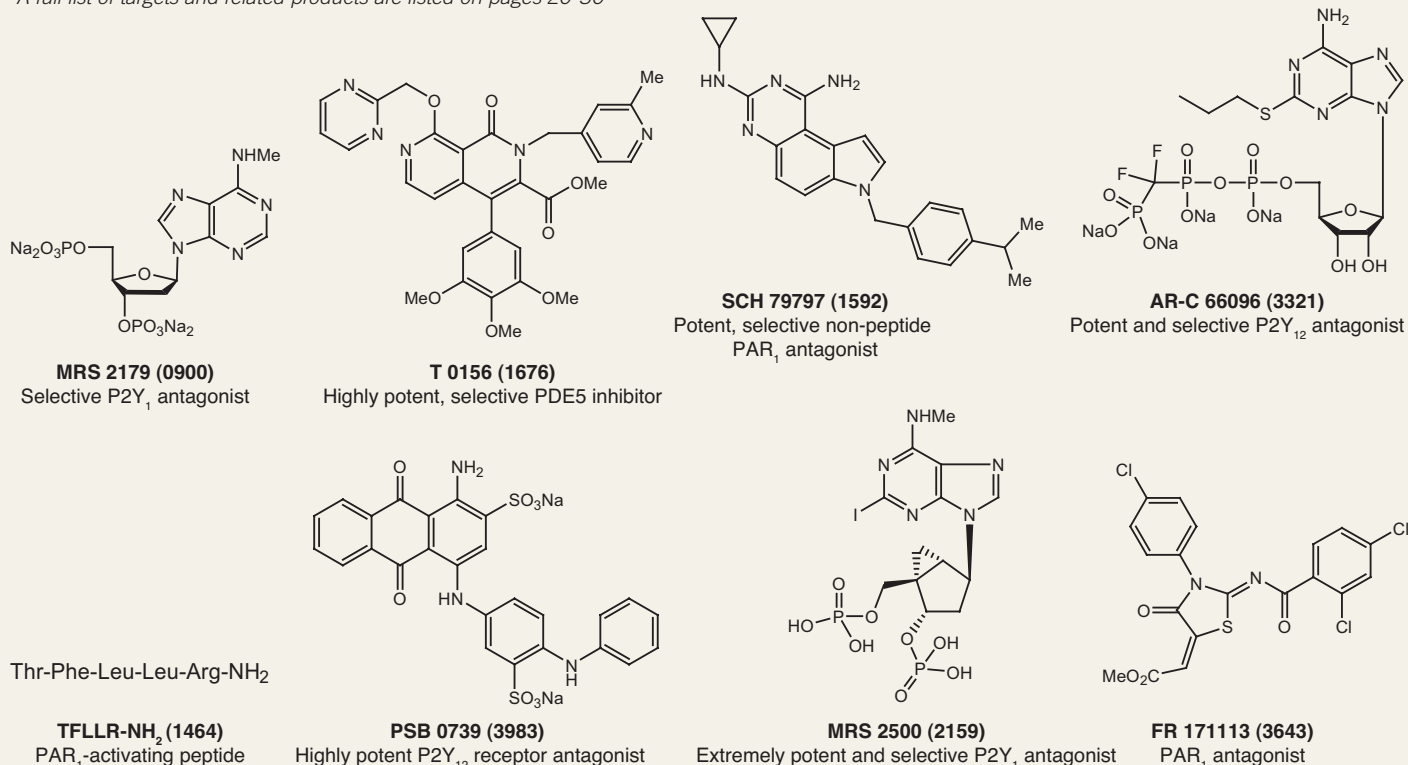
In addition to GPIIb/IIIa, the integrin receptor complex GPIb/V/IX is also integral in thrombosis. Interaction of the GPIb/V/IX complex with von Willebrand factor (vWF) on exposed collagen at the site of vascular damage promotes platelet

adhesion. This is achieved through GPIb/V/IX-mediated phosphorylation of intracellular signaling proteins, including pp60<sup>src</sup>, PLC<sub>γ</sub> and the Fc receptor γ-chain (FcRγ), which initiates and augments the adhesion of platelets with the extracellular matrix and also with neighboring platelets. Phosphorylation of FcRγ by GPIb/V/IX induces FcRγ-GPIb/V/IX functional coupling; this promotes platelet activation through immunoreceptor tyrosine-based activation motif (ITAM)-driven signaling.

Agonist stimulation of platelets also initiates the production of arachidonic acid (Cat. No. 2756) from membrane phospholipids, a reaction which is catalyzed by cytosolic phospholipase A<sub>2</sub> (cPLA<sub>2</sub>). Arachidonic acid can be utilized by both cyclooxygenases (COX) to form prostaglandin H<sub>2</sub> (PGH<sub>2</sub>), and also by lipoxygenases (LOX) to form the lipid mediator hydroperoxy-eicosatetraenoic acid (HPETE). PGH<sub>2</sub> is further metabolized to form prothrombotic eicosanoids including TXA<sub>2</sub>. Inhibition of the PGH<sub>2</sub> synthesis pathway using COX inhibitors such as aspirin (Cat. No. 4092), celecoxib (Cat. No. 3786) and diclofenac (Cat. No. 4454), and the TXA<sub>2</sub> synthase inhibitor, dipyridamole (Cat. No. 0691), is an effective therapeutic strategy for preventing thrombosis. PLA<sub>2</sub> inhibitors including AACOCF<sub>3</sub> (Cat. No. 1462) also exhibit antithrombotic effects by inhibiting the activity of the arachidonic acid pathway.

### Box 3: Thrombosis and Hemostasis Key Products

A full list of targets and related products are listed on pages 20-30



# Atherosclerosis

Products by Category	Page
<b>Cell Adhesion Molecules</b> .....	<b>23</b>
<b>Chemokine Receptors</b> .....	<b>23</b>
<b>Cholesterol Regulation</b> .....	<b>23</b>
<b>Cyclooxygenases</b> .....	<b>24</b>
<b>Cytokine Receptors</b> .....	<b>24</b>
<b>Elastases</b> .....	<b>24</b>
<b>Matrix Metalloproteinases</b> .....	<b>25</b>
<b>Phospholipases</b> .....	<b>27</b>
<b>PPAR</b> .....	<b>28</b>
<b>Protease-activated Receptors</b> .....	<b>28</b>
<b>Urokinase</b> .....	<b>30</b>
<b>Urotensin II</b> .....	<b>30</b>

## Atherosclerosis

Atherosclerosis is a disease of the arterial system that is characterized by the accumulation of fatty deposits within arterial walls (Figure 5). These fatty deposits, known as ‘atherosclerotic plaques’, also contain cellular debris and in advanced plaques, calcium deposits (in the form of hydroxyapatite) are often present. The plaque contents are highly thrombogenic and so are separated from the circulating blood by a fibrous cap, composed mainly of vascular smooth muscle cells and extracellular matrix proteins such as collagen, which are synthesized by these cells (Figure 4).

The exact cause of atherosclerosis remains unknown, but the initial process – termed ‘atherogenesis’ – involves the

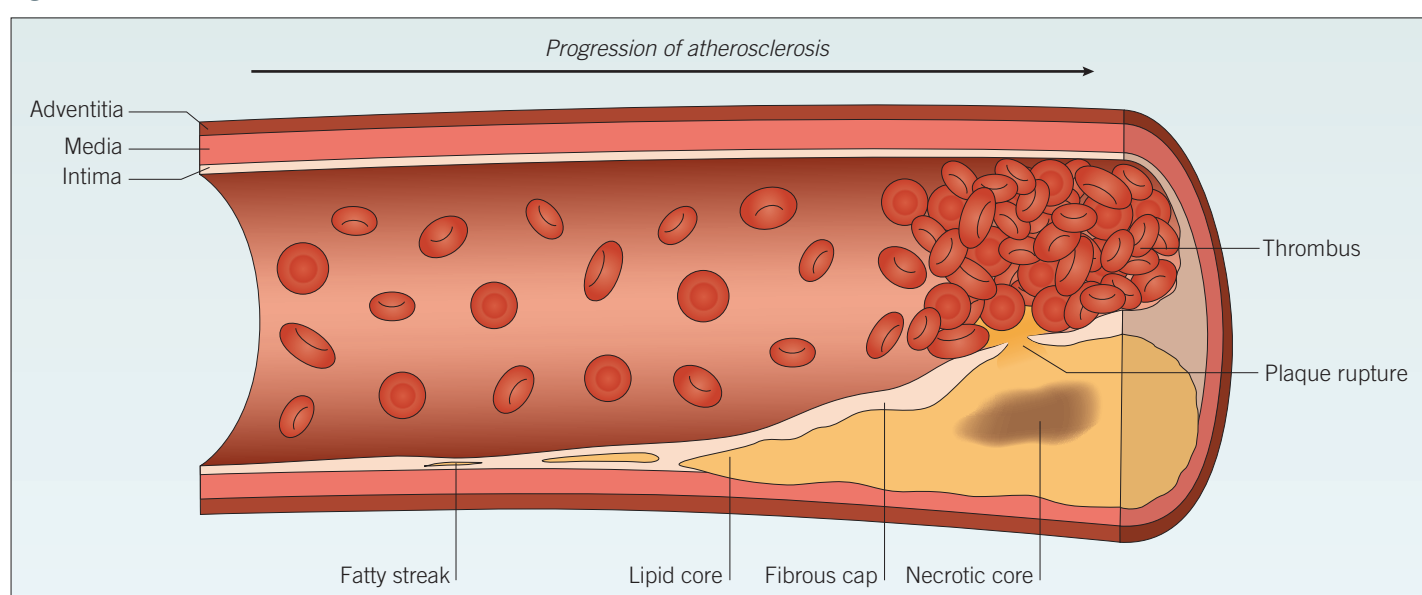
transmigration of monocytes/macrophages across the endothelium into the intima, where they scavenge oxidized lipid (ox-LDL), forming lipid-laden ‘foam cells’. Expression of adhesion molecules by endothelial cells promotes this transmigration. Foam cells undergo apoptosis, forming a lipid-rich core within the developing plaque. The formation of a plaque within the arterial wall triggers an inflammatory response, propagated by the secretion of growth factors and chemokines by resident plaque cells.

Circulating macrophages, recruited to the atherosclerotic plaque through chemotaxis, are a major source of matrix metalloproteinases (MMPs). These zinc-dependent enzymes degrade the fibrous cap, predisposing the plaque to rupture; they may also degrade the elastic laminae between layers of the arterial wall, facilitating the migration of smooth muscle cells from the tunica media to the fibrous cap, thus helping strengthen it.

A thick smooth muscle-cell rich fibrous cap with a small lipid core is characteristic of a ‘stable’ plaque, that is, one which is less likely to rupture. Plaques with a thin fibrous cap and a large, necrotic lipid core are termed ‘unstable’ or ‘vulnerable’ plaques and are more susceptible to rupture. Any defect in the fibrous cap exposes the thrombogenic plaque contents to the circulating blood, triggering thrombosis. Ninety percent of these plaque ruptures are clinically silent, yet the remaining proportion trigger myocardial infarction or stroke, and can be fatal.

In addition to the recruitment of macrophages within an atherosclerotic plaque, other immune mediators are also

**Figure 4 | Atherosclerosis**



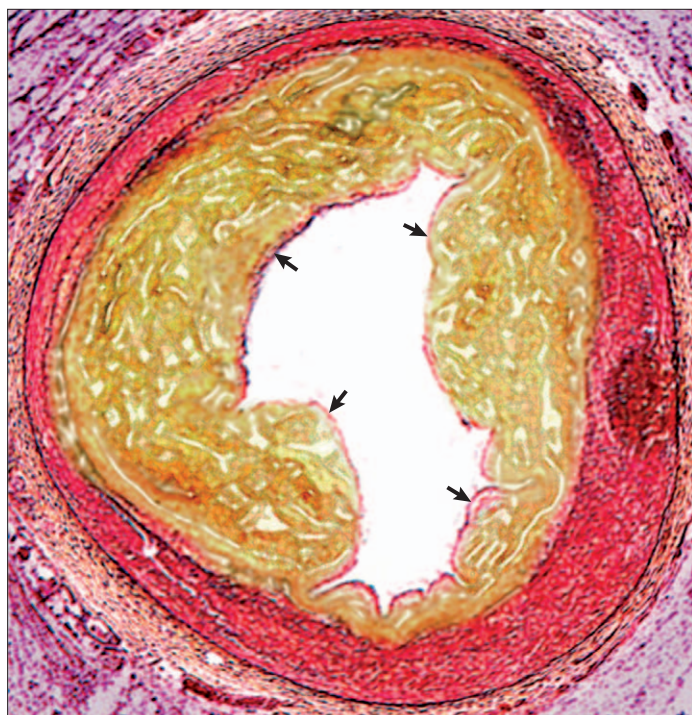
The development of atherosclerosis within a blood vessel wall starts as a fatty streak, composed of a small cluster of lipid-laden macrophages, which progresses over time to form a lipid pool within the vessel wall. Cell debris from apoptotic foam cells within the lipid pool forms a necrotic core. The fibrous cap separates the lipid pool from the blood, but rupture of this cap exposes the thrombogenic plaque contents to circulating blood, triggering thrombus formation.



## Atherosclerosis – continued

present during atherosclerosis. The expression of a number of pro-atherogenic cytokines, including interleukin (IL)-1 and IL-6, are known to be upregulated in atherosclerotic plaques,

**Figure 5 | Section through a coronary artery with an atherosclerotic plaque**



Lipid accumulation and inflammatory cell invasion (depicted in yellow) within an arterial wall (pink) causes vessel narrowing and stenosis. A thin fibrous cap (arrows) separates the plaque contents from the circulating blood. Rupture of the fibrous cap exposes these thrombogenic contents to the blood, triggering thrombosis.

and therefore targeting the receptors of these cytokines may hinder the progression of atherosclerosis. Whilst activation of the immune response within an atherosclerotic plaque may be detrimental, driving the progression and eventual rupture of a plaque, it may equally be beneficial in resolving the inflammation and clearing necrotic foam cells from the core of the plaque. Therefore, targeting the immune response in atherosclerosis may be a 'double-edged sword'.

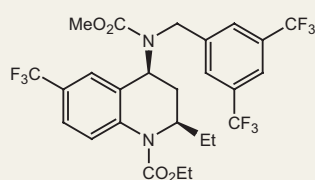
Current therapeutics in atherosclerosis predominantly aim to reduce blood cholesterol, thereby limiting the expansion of an atherosclerotic plaque. The gold standard lipid-lowering drugs are the statins, but other compounds including torcetrapib (Cat. No. 4184), probucol (Cat. No. 2775) and CI 976 (Cat. No. 2227) also exhibit antihypercholesterolemic activity, though their side effect profiles limit their use in the clinic.

Aside from lipid-lowering, a further potential avenue in the treatment of atherosclerosis is the coagulation cascade since it is the formation of an occlusive thrombus, not plaque rupture *per se*, which triggers the pathologies associated with atherosclerosis. Potential targets within the coagulation cascade include thrombin, urokinase plasminogen activator (uPA) and tissue plasminogen activator (tPA). Other mediators involved in platelet activation are also potential therapeutic targets in the prevention and treatment of atherothrombosis; this topic is discussed in greater detail in the 'Thrombosis and Hemostasis' section, page 6.

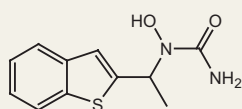
Additional targets with therapeutic potential in atherosclerosis include phospholipases and urotensin II, since the presence of both may trigger foam cell formation, furthering the progression of atherosclerosis.

#### Box 4: Atherosclerosis Key Products

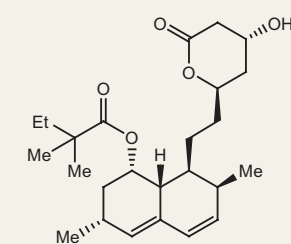
A full list of targets and related products are listed on pages 20-30



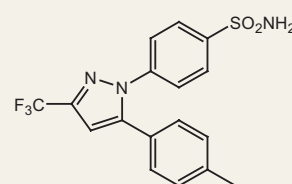
**Torcetrapib (4184)**  
Inhibitor of cholesteryl ester transfer protein (CETP)



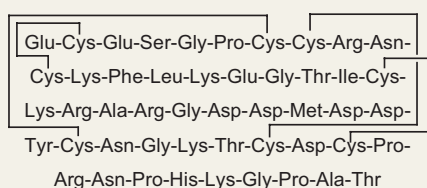
**Zileuton (3308)**  
Orally active 5-LOX inhibitor



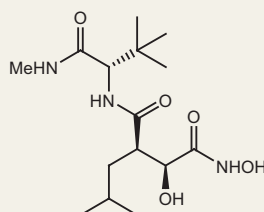
**Simvastatin (1965)**  
HMG-CoA reductase inhibitor



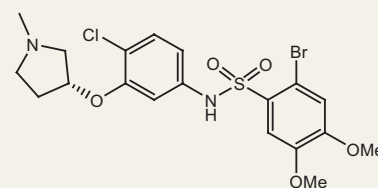
**Celecoxib (3786)**  
Selective COX-2 inhibitor



**Echistatin  $\alpha_1$  isoform (3202)**  
Potent, irreversible  $\alpha_v\beta_3$  integrin antagonist



**Marimastat (2631)**  
Broad spectrum MMP inhibitor



**SB 657510 (3571)**  
Selective urotensin-II receptor antagonist

# Myocardial Infarction

Products by Category	Page
<b><math>\beta</math> Adrenergic Receptors</b>	<b>20</b>
<b>Adenosine Receptors</b>	<b>21</b>
<b>Angiotensin-converting Enzyme</b>	<b>21</b>
<b>Calcium Channels</b>	<b>22</b>
<b>Fatty Acid Oxidation</b>	<b>24</b>
<b>Nitric Oxide</b>	<b>26</b>
<b>PI 3-Kinase</b>	<b>27</b>
<b>Prostanoid Receptors</b>	<b>28</b>
<b>Stem Cells</b>	<b>29</b>

## Myocardial Infarction

Myocardial infarction (MI) – more commonly referred to as a heart attack – is an acute event caused by the interruption of blood supply to regions of the heart, leading to myocardial necrosis. Infarction of a substantial area of the myocardium can disrupt normal conductance of the heart, leading to cardiac arrest.

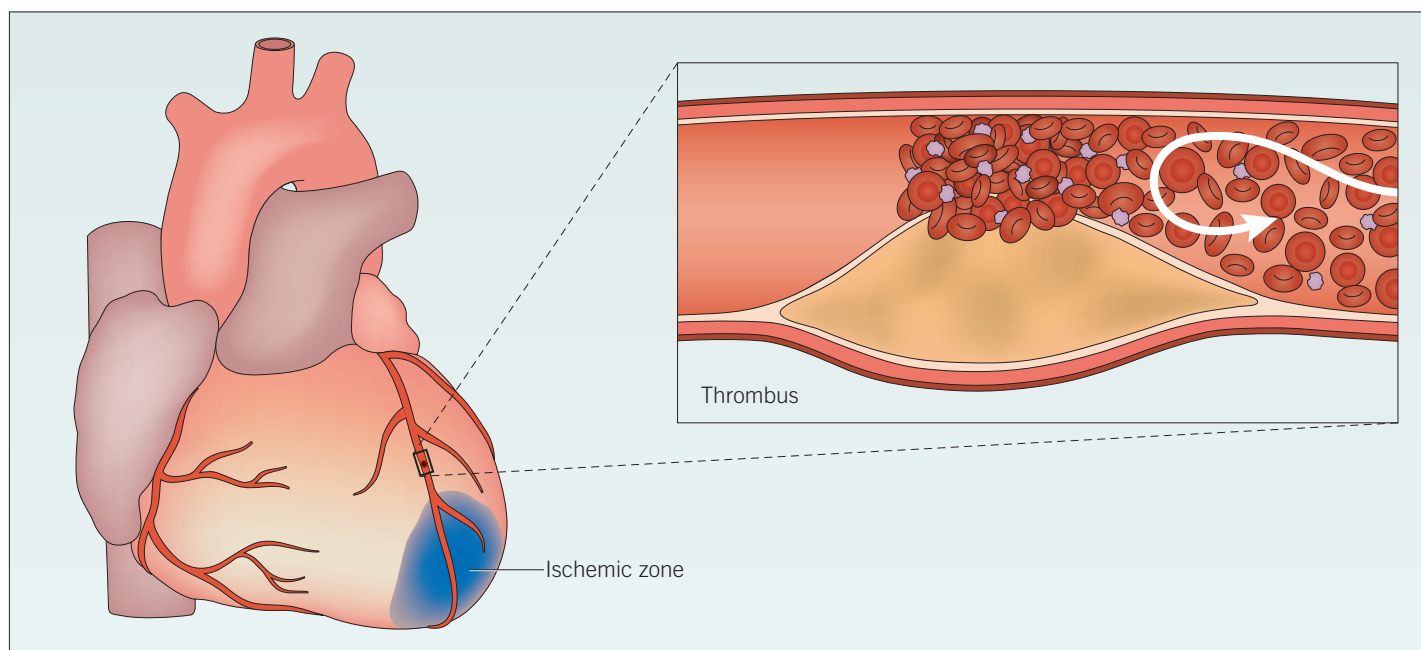
MI is immediately preceded by the presence of an occlusive thrombus within a coronary artery, blocking blood flow to the downstream tissue (Figure 6). The most common cause of an occlusive thrombus within a coronary artery is the rupture of an atherosclerotic plaque (see pages 8-9 for further details on Atherosclerosis). However, the occlusion of a coronary artery

may also result from coronary embolism. This can occur in patients following stent placement, angioplasty, and coronary artery bypass grafting.

One of the few warning symptoms for MI is the occurrence of angina pectoris – a severe, cardiac-induced chest pain which may also radiate down the left arm. Angina is caused by a lack of oxygen to the myocardium due to coronary artery obstruction or spasm, and can be classified as either ‘unstable’ or ‘stable’ angina. Patients with stable angina experience ‘predictable’ chest pain during exertion which resolves following rest or the administration of the NO donor, nitroglycerin. There is little damage to the myocardium during stable angina. However, an episode of unstable angina – that is, chest pain which occurs at rest or in patients with no history of stable angina – may induce myocardial necrosis, albeit at a reduced level to that observed during acute MI. The most common blood biomarkers used to diagnose acute MI or unstable angina are cardiac troponins T (cTnT) and I (cTnI), two components of cardiac muscle whose serum levels rise as a result of myocardial necrosis.

Immediate pharmacological treatment of angina is achieved by the administration of nitroglycerin, but longer term therapy involves either increasing blood supply to the heart using vasodilators such as calcium channel blockers, long-acting nitrates and nicorandil (Cat. No. 2147), or by reducing metabolic demand of the heart through decreasing heart rate by administering  $\beta$ -blockers or ivabradine. A further pharmacological

**Figure 6 | Myocardial infarction**



When an atherosclerotic plaque ruptures, blood flow is greatly impeded and thrombosis may occur. Consequently, the artery is occluded and the supply of oxygen to the heart is restricted. Ischemia results, and if it is sustained the heart muscle tissue may become damaged and die. The likelihood of plaque rupture is influenced by a number of factors, including elevated blood pressure and degree of inflammation. This damaged tissue does not function fully; acute myocardial ischemia results in ionic and metabolic disturbances that affect the membrane and action potentials of myocytes. Arrhythmias and heart failure may occur as a result of the slower conduction of electrical impulses.

## Myocardial Infarction Research – continued

mechanism for preventing angina is to increase ATP production whilst maintaining the same oxygen consumption. One of the first antianginal agents, perhexiline, exerts these effects by inhibiting the mitochondrial enzyme, carnitine palmitoyl-transferase-1 (CPT1). Newer antianginal agents such as ranolazine (Cat. No. 3118) and etomoxir (Cat. No. 4539) also alter fatty acid oxidation, increasing metabolic efficiency and preventing the occurrence of angina.

In addition to vasodilators, pharmacological treatment of acute MI also includes antiplatelet and fibrinolytic agents such as recombinant human tPA, uPA and streptokinase in combination with heparin (Cat. No. 2812) in order to break down the occluding thrombus and restore blood supply to the downstream myocardium. Angioplasty is the preferred method of restoring vessel patency, however, as it carries a lower risk of reocclusion and stroke. Reducing cardiac output by targeting the RAAS is also a useful therapeutic strategy following MI, particularly in patients at risk of developing heart failure (see pages 16–17 for further information on heart failure). Adenosine receptor activation is also an effective therapeutic strategy for reducing myocardial injury following MI. All four adenosine receptor subtypes –  $A_1$ ,  $A_{2A}$ ,  $A_{2B}$  and  $A_3$  – have been shown to be cardioprotective in the ischemic heart.

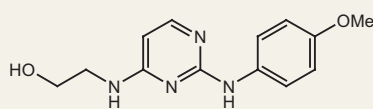
Research continues to identify novel targets for the treatment of MI. One such target is the cathepsin family, a group of enzymes which degrade myofibrillar proteins during MI and promote ventricular remodeling. The beneficial effects of cathepsin inhibition may be two-fold – in addition to reducing ventricular remodeling following MI, cathepsin inhibition may also

prevent atherosclerotic plaque rupture, the principal trigger of MI. A further target in reducing ventricular remodeling and preventing MI-induced heart failure is PI 3-kinase, though its ubiquitous expression limits its use as a therapeutic target.

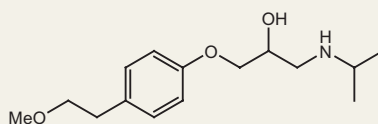
More recent therapy for the treatment of myocardial infarction centers around the use of stem cells to repair damaged myocardium. Following an infarction, damaged myocardium cannot regenerate and so is replaced by non-contractile scar tissue. This alters both the contractility and the conductance of the myocardium and may subsequently lead to the development of an arrhythmia or heart failure. The injection of multipotent cardiac stem cells to the infarcted area of the heart following a myocardial infarction has shown promise in facilitating regeneration of damaged myocardium, but their availability is limited. As a result, research efforts are currently focused on producing cardiomyocytes by differentiating more readily available stem cell populations, such as undifferentiated skeletal myoblasts or bone marrow-derived adult stem cells. Inducing cardiomyogenic function in these stem cell populations has been achieved using a range of methods. These include cardiac preconditioning, whereby stem cells are differentiated in media previously used to culture primary cardiomyocytes; and also by using small molecule inhibitors such as cardiogenol C (Cat. No. 3851) and XAV 939 (Cat. No. 3748) which modulate stem cell signaling pathways including the Wnt/ $\beta$ -catenin pathway. Facilitating the repair and regeneration of damaged myocardium following infarction, together with preventing aberrant remodeling, represent promising future therapeutic avenues within the field of myocardial infarction research.

**Box 5: Myocardial Infarction Key Products**

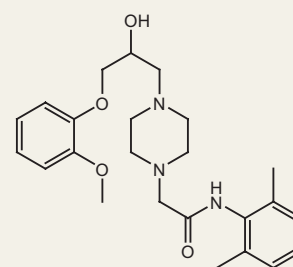
A full list of targets and related products are listed on pages 20-30



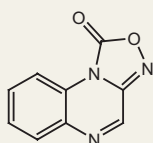
**Cardiogenol C (3851)**  
Induces cardiomyogenesis in ESCs



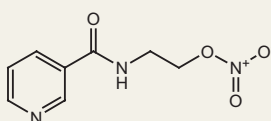
**Metoprolol (3256)**  
Selective  $\beta_1$  antagonist



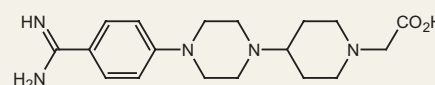
**Ranolazine (3118)**  
Antianginal, activates pyruvate dehydrogenase



**ODQ (0880)**  
Selective inhibitor of NO-sensitive guanylyl cyclase



**Nicorandil (2147)**  
 $K_{ir}6$  ( $K_{ATP}$ ) channel opener and NO donor



**GR 144053 (1263)**  
Glycoprotein IIb/IIIa (integrin  $\alpha_{IIb}\beta_3$ ) receptor antagonist. Antithrombotic



# Ischemia/Reperfusion Injury

Products by Category	Page
<b>Adenosine Receptors</b>	<b>21</b>
<b>Apoptosis</b>	<b>22</b>
<b>Calcium Signaling</b>	<b>22</b>
<b>Mitochondrial Calcium Uniporter</b>	<b>25</b>
<b>Mitochondrial Permeability Transition Pore</b>	<b>25</b>
<b>Na<sup>+</sup>/Ca<sup>2+</sup> Exchanger</b>	<b>25</b>
<b>Na<sup>+</sup>/H<sup>+</sup> Exchanger</b>	<b>25</b>
<b>Oxidative Phosphorylation</b>	<b>26</b>
<b>Potassium Channels</b>	<b>27</b>
<b>SERCA</b>	<b>29</b>
<b>Sodium Channels</b>	<b>29</b>

## Ischemia/Reperfusion Injury

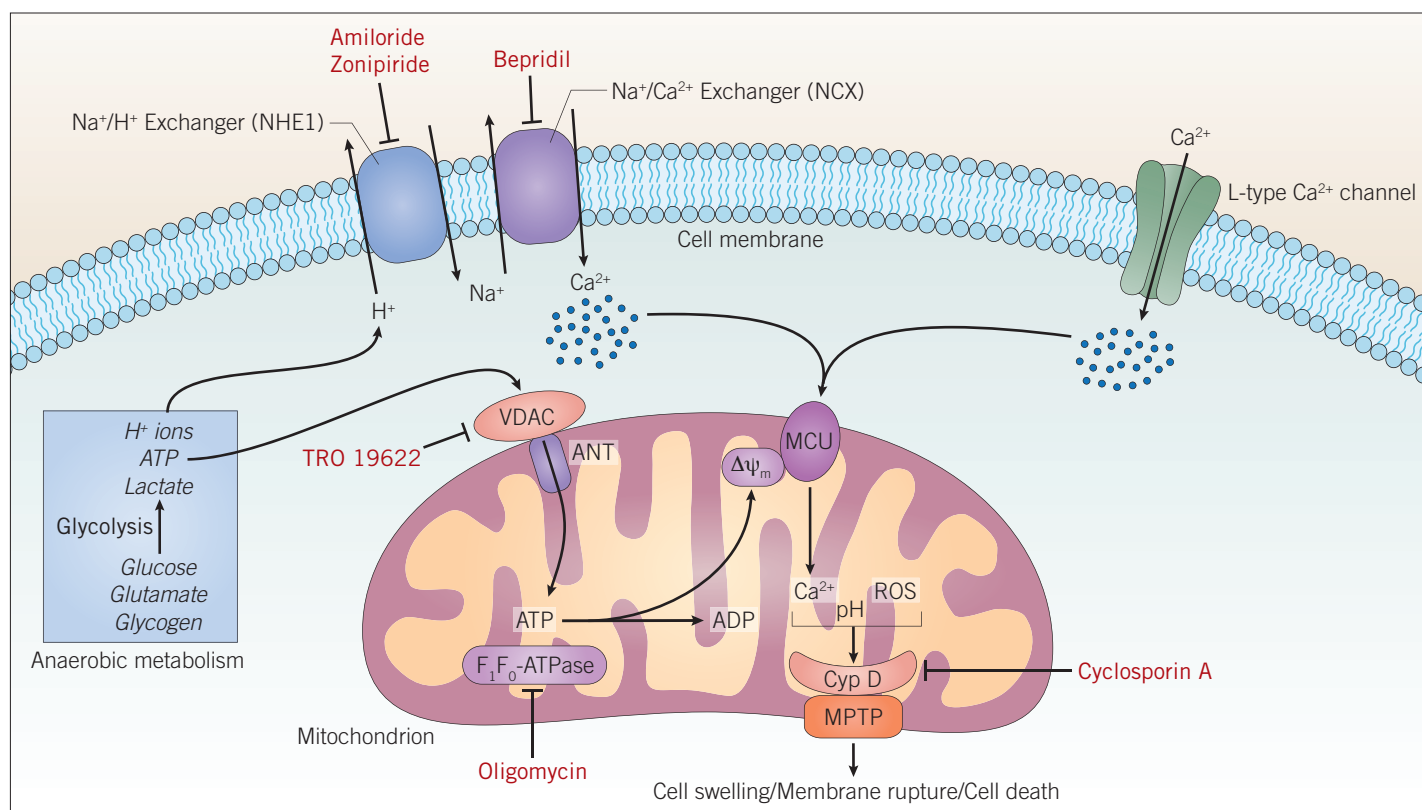
Myocardial ischemia, also known as cardiac ischemia, is defined as the deprivation of oxygen and nutrients to the heart. This phenomenon occurs during a myocardial infarction, when an occlusive thrombus within a coronary artery prevents blood

supply to the myocardium, but can occur during cardiac surgery as a result of pharmacological intervention to temporarily stop the heart. Reperfusion restores blood supply to ischemic tissue, but this is paradoxically associated with further tissue damage.

Under ischemic conditions, the lack of oxygenated blood supply to the myocardium means that the energy demands of the heart cannot be met. Numerous metabolic changes result from cardiac ischemia, and if it is prolonged, ischemia may result in irreversible injury. Hypoxic ischemic conditions and the rapid decline in oxygen availability blocks ATP production by oxidative phosphorylation, leading to reduced contractility of the heart. A prolonged reduction in contractility can itself induce ventricular dysfunction.

During anaerobic metabolism, glucose, glycogen and glutamate are broken down and lactate, alanine, succinate and hydrogen ions build up. A build-up of hydrogen ions lowers the pH of the intracellular and extracellular environment. The acidic environment inside and outside cardiomyocytes gradually affects ion homeostasis, leading to increases in intracellular Na<sup>+</sup> concentration. Increased Na<sup>+</sup> concentration prompts a resultant

**Figure 7 | Ionic disturbances in I/R injury within a cardiac muscle cell**



A rise in intracellular Ca<sup>2+</sup> concentration ([Ca<sup>2+</sup>]<sub>i</sub>) is evident in ischemia and early reperfusion. This increase has been shown to precede irreversible cardiac injury. Decrease in ATP production, as a result of ischemia, lowers the intracellular pH. This change results in the increased activity of Na<sup>+</sup>/H<sup>+</sup> and Na<sup>+</sup>/Ca<sup>2+</sup> exchangers, thus increasing [Ca<sup>2+</sup>]<sub>i</sub>. ATP, generated by glycolysis, is used by F<sub>1</sub>F<sub>0</sub>-ATPase to create a mitochondrial membrane potential (Δψ<sub>m</sub>). This is utilized by the mitochondrial calcium uniporter; an increase in mitochondrial Ca<sup>2+</sup> concentration, combined with ROS activity and a normalized pH, prompts the opening of the mitochondrial pore during reperfusion. Abbreviations: ANT – adenine nucleoside translocator; ATP – adenosine triphosphate; Cyp D – Cyclophilin D; MCU – mitochondrial calcium uniporter; MPTP – mitochondrial permeability transition pore; ROS – reactive oxygen species; VDAC – voltage-dependent anion channel.



## Ischemia/Reperfusion Injury – continued

increase in  $\text{Ca}^{2+}$  ions in a phenomenon known as ‘ $\text{Ca}^{2+}$  overloading’. Apoptosis and necrosis of cardiomyocytes induced by  $\text{Ca}^{2+}$  overloading can cause irreparable damage to the heart.

The reperfusion of an ischemic myocardium following a period of ischemia is also a focus of research. Restoration of blood supply to the ischemic zone triggers tissue damage via the release of intracellular enzymes, sarcolemmal rupture,  $\text{Ca}^{2+}$  influx and cardiomyocyte hypercontracture. Rupture of the sarcolemmal membranes leads to the movement of  $\text{Na}^+$  ions through gap junctions between adjacent cells and the induction of reverse  $\text{Na}^+/\text{Ca}^{2+}$  exchange, propagating the damage to neighbouring myocytes. This process of initial ischemia-related damage followed by further damage induced by reperfusion is known as ischemia/reperfusion (I/R) injury.

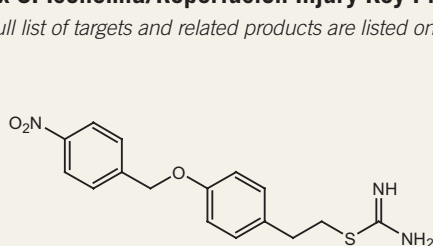
The cell death that underlies I/R injury is characterized by features typical of apoptosis, autophagy and necrosis. A key regulator of both apoptotic and necrotic cell death is the mitochondrial permeability transition pore (MPTP) (Figure 7). Low pH, induced during ischemia, inhibits MPTP opening; only upon reperfusion does pH return to normal. Elevated matrix  $\text{Ca}^{2+}$  concentrations and reactive oxygen species (ROS) are the primary activators of the MPTP during I/R. Inhibition of ion exchangers that influence cytosolic  $\text{Ca}^{2+}$  levels, such as the  $\text{Na}^+/\text{H}^+$  exchanger (NHE) and  $\text{Na}^+/\text{Ca}^{2+}$  exchanger (NCX), have been shown to reduce I/R injury. One such inhibitor – zoniporide (Cat. No. 2727) – selectively inhibits NHE1 and provides cardioprotection from ischemic injury *in vivo*. In addition to agents targeting the NHE and NCX, inhibitors of MPTP, such as cyclosporin A (Cat. No. 1101), help protect against reperfusion injury.

A further cardioprotective strategy currently under investigation for I/R injury is ischemic preconditioning. This method aims to reduce the damage associated with I/R injury by subjecting the vascular system to brief, sublethal periods of ischemia. An advantage of this method is that the same protective effects can be obtained even when inducing ischemia in a tissue distinct from the heart, such as the upper or lower limbs; this is known as remote preconditioning. The cardioprotective tissue response to ischemic preconditioning is thought to involve a number of biological targets, including adenosine receptors. Activation of these receptors prior to ischemia or during reperfusion has been shown to confer cardioprotection; for example, the subtype-selective  $\text{A}_3$  agonist, IB-MECA (Cat. No. 1066) exhibits cardioprotective properties in a rat model.

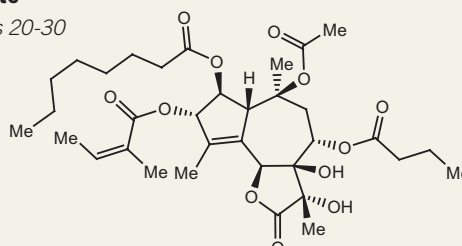
Newer cardioprotective targets include GSK-3 $\beta$ , a multifunctional kinase that has also been linked to protection against I/R damage; the mitochondrial calcium uniporter (MCU), which has been linked to the cardioprotective response to ischemic preconditioning; and the sarco/endoplasmic reticulum  $\text{Ca}^{2+}$ -ATPase (SERCA). Enhancement of SERCA activity has been shown to reduce infarct size and preserve cardiac function in a rodent model of transient myocardial ischemia. Administration of interventional drugs at the initiation of reperfusion is limited since they must be introduced within 10 minutes of reperfusion starting, yet therapeutic targeting of intracellular processes invoked during both ischemia and reperfusion remain a promising strategy for the prevention and/or limitation of both the occurrence and the extent of I/R injury.

### Box 6: Ischemia/Reperfusion Injury Key Products

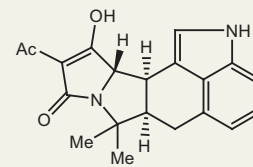
A full list of targets and related products are listed on pages 20-30



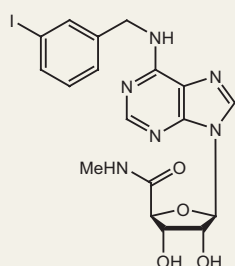
**KB-R7943 (1244)**  
Na<sup>+</sup>/Ca<sup>2+</sup> exchange inhibitor (reverse mode)



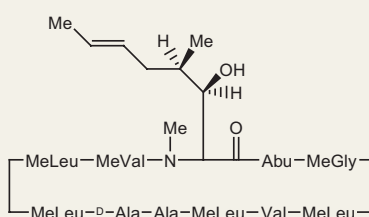
**Thapsigargin (1138)**  
Potent inhibitor of SERCA



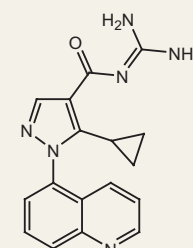
**Cyclopiazonic acid (1235)**  
Inhibitor of SERCA



**IB-MECA (1066)**  
 $\text{A}_3$  selective agonist



**Cyclosporin A (1101)**  
Inhibits formation and opening of the MPTP



**Zoniporide (2727)**  
Selective NHE1 inhibitor

# Arrhythmias

Products by Category	Page
<a href="#">β Adrenergic Receptors</a>	<a href="#">20</a>
<a href="#">Calcium Channels</a>	<a href="#">22</a>
<a href="#">Potassium Channels</a>	<a href="#">27</a>
<a href="#">Sodium Channels</a>	<a href="#">29</a>

## Arrhythmias

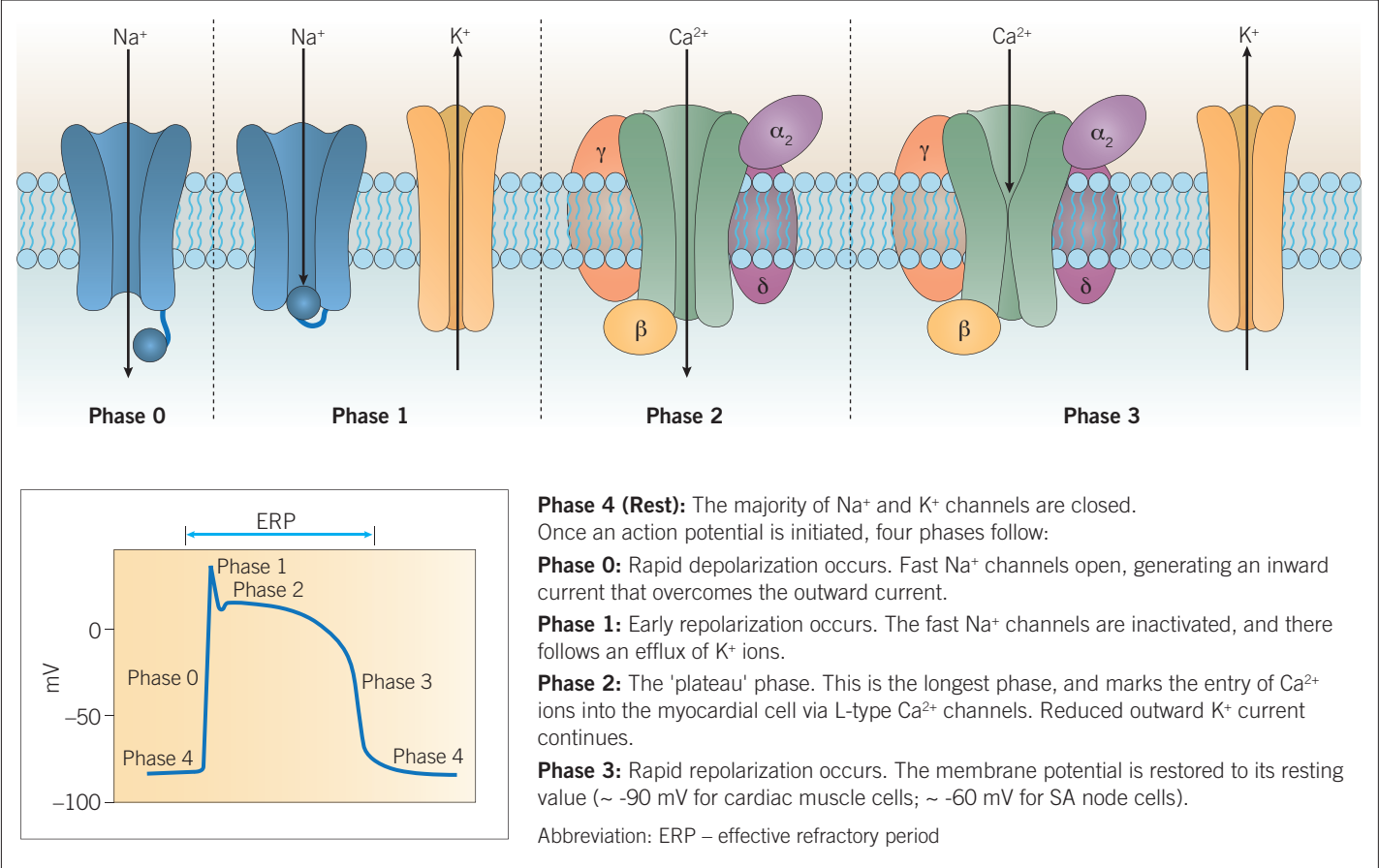
An arrhythmia (also known as cardiac dysrhythmia) is defined as an irregular heartbeat, and results from abnormal electrical activity in the heart. There are various different types of arrhythmia, often resulting in a heartbeat that is too fast (tachycardia) or too slow (bradycardia). Atrial and ventricular fibrillation, which are the most common cardiac arrhythmias, account for 10-20% of all deaths among adults in the Western world. The incidence of atrial fibrillation increases with age; it not only affects cardiac function, but also increases the risk of stroke and may worsen heart failure.

A normal heartbeat is driven by various stages of membrane depolarization and repolarization in single heart cells, propagating from the sinoatrial (SA) node to the atrium and the ventricle. There are two types of action potentials: the fast response action potential, which occurs in cardiac muscle and Purkinje

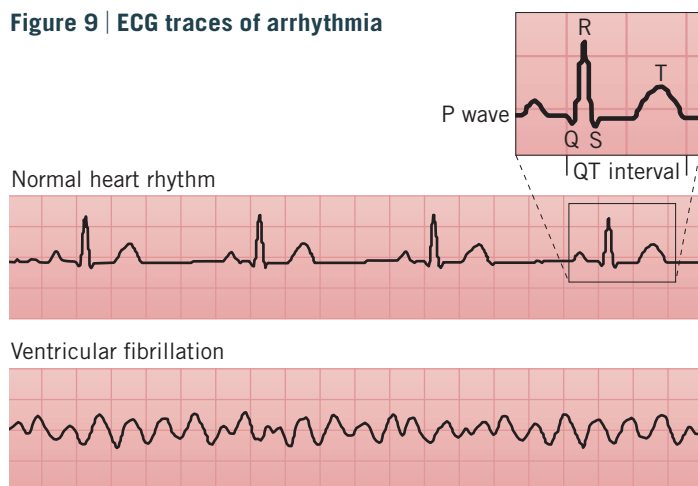
fibers; and the slow response, evident in the SA and atrioventricular (AV) nodes. Action potentials enable rapid changes in heart rate. Contraction of the cardiac muscle occurs in response to depolarization. By generating action potentials and setting off a wave of depolarization, the SA node thus acts as a pacemaker, setting the rate of contraction of the heart. Dysfunction of the SA node may therefore result in an irregular heartbeat. The mechanism of normal cardiac action potential generation in cardiac muscle cells is depicted in Figure 8. The effective refractory period (ERP) is a mechanism that helps protect the heart from arrhythmias, by preventing the generation of new action potentials during the propagation of an existing one. Antiarrhythmics such as quinidine (Cat. No. 4108) can be used to prolong the ERP, preventing premature activation. However, quinidine also prolongs the QT interval (Figure 9) and can induce Torsades de pointes (TdP; a type of ventricular tachycardia that can be transient or lead to lethal ventricular fibrillation).

In addition to SA node dysfunction, arrhythmias can also occur as a result of abnormalities in the electrophysiology of heart cells or in cell-to-cell (impulse) propagation, which takes place through gap junctions. These enable conduction of a wave of depolarization between cells. Ion channels are responsible for the conduction of coordinated electrical impulses, and

Figure 8 | Action potentials in cardiac muscle cells



## Arrhythmias – continued

**Figure 9 | ECG traces of arrhythmia**

Arrhythmias can be detected by electrocardiography (ECG or EKG), which measures the electrical activity of the heart. A normal ECG trace will have a consistent, regular form, representing the different intervals involved in cardiac rhythm. This includes the QT interval, during which the left and right ventricles depolarize and repolarize. During ventricular fibrillation, a type of arrhythmia, the heart does not contract in an ordered fashion; the absence of normal heart rhythm is apparent when comparing ECG traces.

consequently dysregulation of their activity has been linked to the development of arrhythmias. For example, mutations in genes encoding the  $K_{V11.1}$  (hERG) potassium channel, the sodium  $Na_v1.5$  channel, and the calcium  $Ca_v1.2$  channel have been linked to long QT syndrome (LQTS). In LQTS, the QT interval is extended and repolarization is delayed; this increases

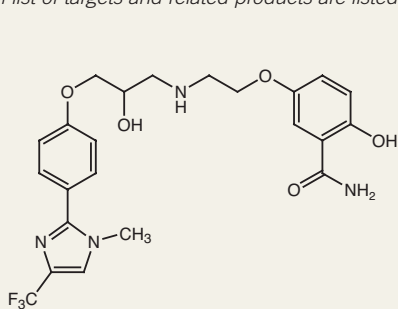
the risk of Torsades de pointes. Cardiac ion channel blockade represents the traditional action of antiarrhythmic drugs. Directly or indirectly altering ion channel conductance changes the characteristics of cardiac action potentials and decreases atrial fibrillation.

In addition to ionic imbalances, arrhythmogenic stimuli in the heart include: metabolic substances (e.g. phospholipids and eicosanoids), thrombosis, atheromas and coronary artery spasm (angina). Arrhythmias can also result from myocardial ischemia.

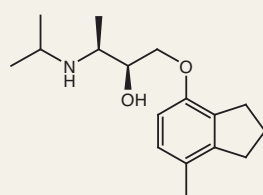
Antiarrhythmic drug therapy aims to restore normal cardiac rhythm and conduction, and to prevent more serious arrhythmias from occurring. Vaughan Williams created one of the most widely used classification schemes for antiarrhythmic drugs. The scheme divides antiarrhythmic drugs into five classes (I-V), each of which concerns a different target. Class I compounds are sodium channel blockers; class II are beta-blockers; class III contains agents that target potassium channels; and class IV agents block calcium channels. Class V includes drugs that act via an unknown mechanism. Within these classes exist further subclasses, which exhibit slightly different properties at different points of the cardiac action potential, and which may affect the overall duration of the action potential. For specific types of arrhythmia, different agents may also be used: for example, adenosine (Cat. No. 3624) and verapamil (Cat. No. 0654) may be used to treat supraventricular tachycardia. The main advances in arrhythmic therapy have been made in the use of electronic devices, such as artificial pacemakers, and direct current cardioversion.

**Box 7: Arrhythmia Key Products**

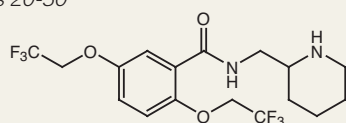
A full list of targets and related products are listed on pages 20-30



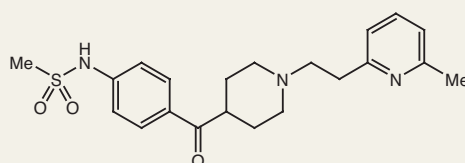
**CGP 20712 (1024)**  
Highly potent and selective  $\beta_1$  antagonist



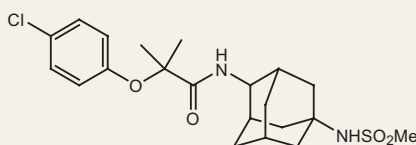
**ICI 118,551 (0821)**  
Very selective  $\beta_2$  antagonist



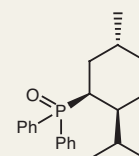
**Flecainide (1470)**  
Cardiac  $Na^+$  channel blocker. Antiarrhythmic



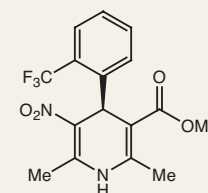
**E-4031 (1808)**  
 $K_{V11.1}$  (hERG) channel blocker; inhibits rapid delayed rectifier  $K^+$  current ( $I_{Kr}$ )



**JNJ 303 (3899)**  
Potent and selective  $I_{Ks}$  blocker



**DPO-1 (2533)**  
Blocker of  $K_{V1.5}$  channel and  $I_{Kur}$  current



**(S)-(-)-Bay K 8644 (1546)**  
 $Ca^{2+}$ -channel activator (L-type)

# Heart Failure

Products by Category	Page
<b><math>\beta</math> Adrenergic Receptors</b>	20
<b>Aldosterone Receptors</b>	21
<b>Angiotensin-converting Enzyme</b>	21
<b>Angiotensin II Receptors</b>	21
<b>GRK2</b>	24
<b>Na<sup>+</sup>/K<sup>+</sup> ATPase</b>	25
<b>Nitric Oxide</b>	26
<b>NKCC Cotransporter</b>	26
<b>PI 3-Kinase</b>	27

## Heart Failure

Heart failure, also known as congestive heart failure or CHF, is an inability of the heart to pump sufficient blood around the body. Heart failure typically occurs secondary to an existing pathology that alters cardiac function. Examples of syndromes that can precede heart failure include myocardial infarction, arrhythmia or infection. These can also cause dilated cardiomyopathy, a condition which accounts for around one third of all cases of heart failure.

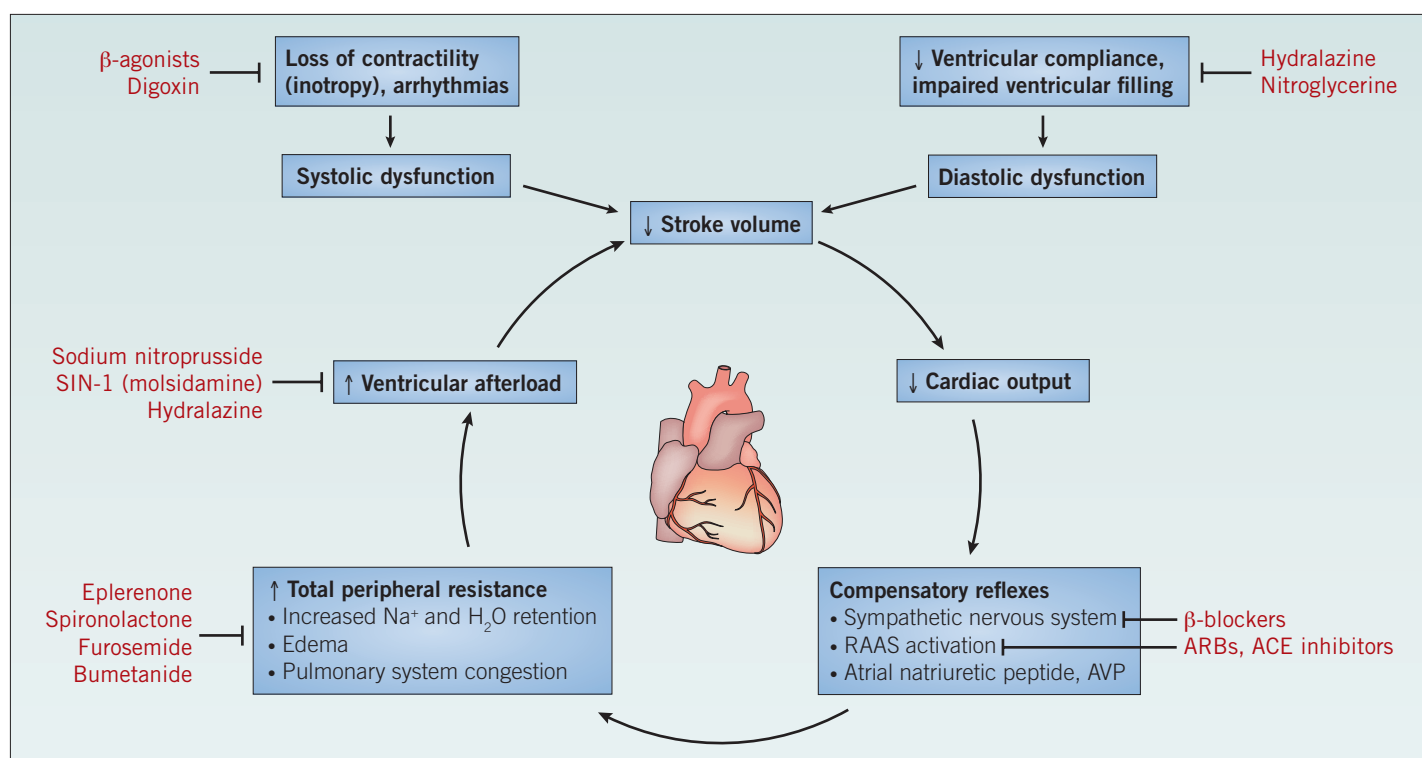
The pathogenesis of heart failure is cyclical and progressive; endogenous mechanisms, which are activated during heart failure in an attempt to counteract the symptoms, actually worsen

cardiac function. Cardiac dysfunction, either systolic or diastolic, triggers a decrease in stroke volume and a resultant decrease in cardiac output. In healthy individuals the body responds to decreases in cardiac output by initiating the renin-angiotensin-aldosterone system to promote fluid retention, and also by activating the sympathetic nervous system to cause peripheral vasoconstriction. Under normal circumstances this counteracts the imbalance in stroke volume, restoring cardiac output to normal levels.

In patients with heart failure the increase in blood volume, together with the heightened peripheral resistance and elevated levels of circulating catecholamines, causes an increased load on the already weakened ventricles with each contraction, and the stroke volume does not return to normal levels. Repeated cycles of this process further weaken the ventricle walls, prompting ventricular hypertrophy and a decreased force of contraction (Figure 10).

Surgical intervention is available for patients with heart failure – current options include the implantation of a left ventricular assist device – but these are invasive and are not suitable for all patients. Pharmacological intervention is common and there are a range of drugs available to target the different stages of the heart failure mechanism. However, none of these current therapeutic options are able to reverse the pathology of heart failure and act only to slow the progression of the disease.

**Figure 10 | The mechanism of heart failure**



Abbreviations: ACE – angiotensin-converting enzyme; ARBs – angiotensin II receptor blockers; AVP – arginine vasopressin; RAAS – renin-angiotensin-aldosterone system.



## Heart Failure – continued

The two most common pharmacological therapies for heart failure are to increase intracellular calcium concentration within myocytes by activating second messenger signaling pathways, and also to block or counteract the neurohormonal compensatory reflexes through the inhibition of the RAAS. Pharmacological agents which trigger a rise in intracellular calcium include the positive inotrope digoxin (Cat. No. 4583), an Na<sup>+</sup>/K<sup>+</sup> ATPase blocker;  $\beta$  adrenergic receptor agonists; and phosphodiesterase inhibitors.

However, indirectly targeting the signaling pathways involved in cardiac contractility also induces mechanism-related adverse effects. More recent therapy using small molecule drugs – such as the cardiac myosin activator, omecamtiv mecarbil – looks to be a promising and more effective strategy for improving contractility in patients with heart failure.

Blocking the neurohormonal reflexes in the failing heart is a more efficacious therapeutic strategy for heart failure compared to increasing contractility, and is achieved using ACE inhibitors such as perindopril (Cat. No. 4302), or ARBs including valsartan (Cat. No. 4216).

The administration of diuretics reduces symptoms associated with heart failure, such as peripheral edema, but it does not reverse or halt the disease pathology. Examples of diuretics used in the treatment of heart failure include furosemide (Cat. No. 3109), an NKCC co-transporter inhibitor, and spironolactone (Cat. No. 2968), an aldosterone receptor antagonist.

Vasodilators such as nicorandil (Cat. No. 2147), SIN-1 (Cat. No. 0756) and hydralazine are also used in the treatment of heart failure since they induce peripheral vasodilation, thereby reducing ventricular afterload.

In addition to the existing compounds, future therapeutic targets include the collagenase enzymes MMP-2 and MMP-9. The expression of these metalloproteinases is increased in heart failure whilst the expression of their endogenous inhibitors, tissue inhibitors of metalloproteinases (TIMPs), has been shown to be downregulated in the same tissue. In support of this, experimental inhibition of MMP-9 reduces ventricular dilatation in a model of heart failure.

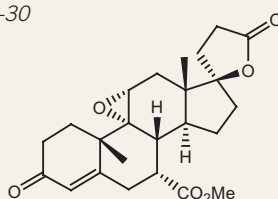
A further therapeutic target in heart failure, central to disease progression, is the activity of G protein-coupled receptor kinases (GRKs), in particular the cardiomyocyte-expressed GRK2. GRK2 activation during heart failure leads to the desensitization of  $\beta$  adrenergic receptors ( $\beta$ ARs) and therefore reduces contractility and depresses cardiac function. Circulating levels of GRK2 and GRK5 are upregulated in the early stages of heart failure, whilst both cardiac isoforms of  $\beta$ ARs –  $\beta_1$ ARs and  $\beta_2$ ARs – have been shown to be downregulated, or non-functional. Cardiomyocyte-specific overexpression of GRK2 triggered  $\beta$ AR uncoupling and led to a reduction in contractility, whereas the expression of an inactive form of GRK2 in either cardiac or adrenal tissue caused an increase in contractility in response to adrenergic stimulation. Therefore, cardiomyocyte- or adrenal-specific inhibition of GRK2 may represent a new therapeutic target for heart failure.

**Box 8: Heart Failure Key Products**

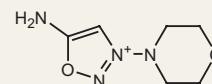
A full list of targets and related products are listed on pages 20-30

Trp-Lys-Lys-Glu-Leu-Arg-Asp-Ala-Tyr-Arg-  
Glu-Ala-Gln-Gln-Leu-Val-Gln-Arg-Val-Pro-  
Lys-Met-Lys-Asn-Lys-Pro-Arg-Ser

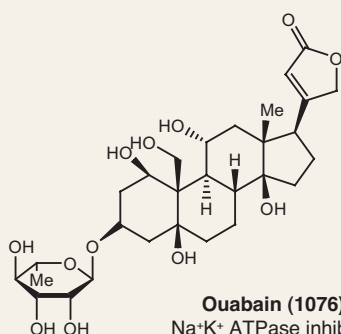
**GRK2i (3594)**  
GRK2 inhibitory polypeptide;  
G $\beta\gamma$  antagonist



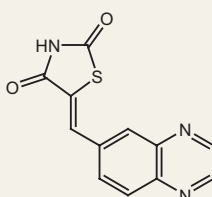
**Eplerenone (2397)**  
Selective mineralocorticoid  
receptor antagonist



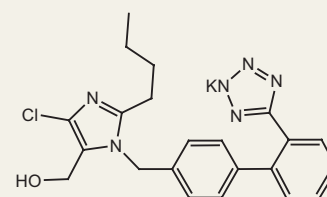
**SIN-1 (0756)**  
Water-soluble NO donor



**Ouabain (1076)**  
Na<sup>+</sup>/K<sup>+</sup> ATPase inhibitor



**AS 605240 (3578)**  
Potent and selective  
PI 3-kinase  $\gamma$  (PI3K $\gamma$ ) inhibitor



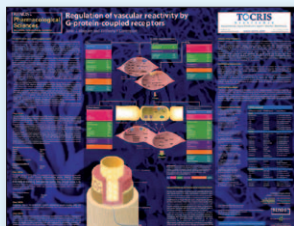
**Losartan (3798)**  
Selective, nonpeptide AT<sub>1</sub> antagonist

# List of Acronyms

Acronym	Definition
ACE	Angiotensin-converting enzyme
ACh	Acetylcholine
ANT	Adenine nucleotide translocase
ARB	Angiotensin II receptor blocker
AV	Atrioventricular
AVP	Arginine Vasopressin
$\beta$ AR	$\beta$ -androgen receptor
BP	Blood pressure
CaM	Calmodulin
cAMP	Cyclic adenine monophosphate
cGMP	Cyclic guanosine monophosphate
CHF	Congestive heart failure
CK	Creatine kinase
COX	Cyclooxygenase
cPLA <sub>2</sub>	Cytosolic phospholipase A <sub>2</sub>
CPT1	Carnitine palmitoyltransferase 1
cTnI	Cardiac troponin I
cTnT	Cardiac troponin T
DAG	Diacylglycerol
DNA-PK	DNA protein kinase
ECG/EKG	Electrocardiography/Elektrokardiogramm
eNOS	Endothelial nitric oxide synthase
ERP	Effective refractory period
ET	Endothelin
FcR $\gamma$	Fc receptor $\gamma$ -chain
GPVI	Glycoprotein VI
GRK	G protein-coupled receptor kinase
GSK-3 $\beta$	Glycogen synthase kinase-3 $\beta$
GTP	Guanosine triphosphate
hERG	Human ether-à-go-go-related gene
HPETE	Hydroperoxyeicosatetraenoic acid
IGF-1	Insulin-like growth factor-1
IL	Interleukin
InsP3R	IP <sub>3</sub> receptor
IP <sub>3</sub>	Inositol triphosphate
IPR	I prostanoid receptor
ITAM	Immunoreceptor tyrosine-based activation motif
K <sub>ATP</sub>	ATP-sensitive potassium channel

Acronym	Definition
K <sub>ir</sub>	Inward rectifying potassium channel
K <sub>v</sub>	Voltage-gated potassium channel
LOX	Lipoxygenase
LQTS	Long QT syndrome
MCU	Mitochondrial calcium uniporter
MLCK	Myosin light chain kinase
MMP	Matrix metalloproteinase
MPTP	Mitochondrial permeability transition pore
mTOR	Mammalian target of rapamycin
NA	Noradrenalin
NCX	Na <sup>+</sup> /Ca <sup>2+</sup> exchanger
NHE	Na <sup>+</sup> /H <sup>+</sup> exchanger
NKCC	Na <sup>+</sup> /K <sup>+</sup> /2Cl <sup>-</sup> cotransporter
NO	Nitric oxide
NTS	Nucleus tractus solitarius
ox-LDL	Oxidized low-density lipoprotein
PAR	Protease-activated receptor
PBR	Peripheral benzodiazepine receptor
PDGF	Platelet-derived growth factor
PGH <sub>2</sub>	Prostaglandin H <sub>2</sub>
PGI <sub>2</sub>	Prostacyclin
PI 3-K	Phosphoinositide 3-kinase
PIP <sub>2</sub>	Phosphatidylinositol 4,5-bisphosphate
PKG	Protein kinase G
PLC	Phospholipase C
PMN	Polymorphonuclear leukocyte
RAAS	Renin-angiotensin-aldosterone system
ROS	Reactive oxygen species
SA	Sinoatrial
SERCA	Sarco/endoplasmic reticulum Ca <sup>2+</sup> -ATPase
sGC	Soluble guanylyl cyclase
TGF-BRI	Transforming growth factor-B type I receptor
TIMP	Tissue inhibitor of metalloproteinase
tPA	Tissue plasminogen activator
TXA <sub>2</sub>	Thromboxane A <sub>2</sub>
uPA	Urokinase plasminogen activator
VDAC	Voltage-dependent anion channel
vWF	von Willebrand factor

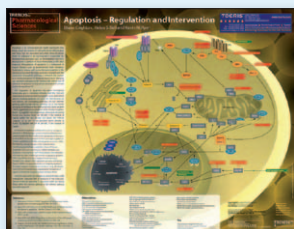
# Related literature from Tocris that you may be interested in:



## **Regulation of Vascular Reactivity by G-protein-coupled Receptors**

J. Maguire and A. Davenport, University of Cambridge

Cardiovascular disease remains one of the major causes of morbidity and mortality in the Western world and therefore this therapeutic area continues to be of great interest to researchers. This poster highlights the key GPCRs regulating vascular reactivity.



## **Apoptosis – Regulation and Intervention**

D. Crighton *et al.* Beatson Institute for Cancer Research

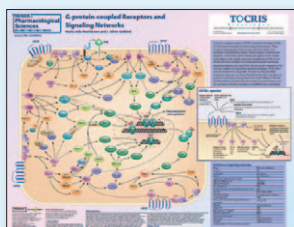
Apoptosis is an orchestrated cell death mechanism that brings about removal of cells without inflammation and stress. The key signaling pathways associated with the regulation of apoptosis are summarized in this poster.



## **7-TM Receptor Signaling**

T. Kenakin, GlaxoSmithKline Research and Development  
R. Lefkowitz and J. Violin, Duke University Medical Center  
M. Bouvier and G. Oligny-Longpré, Université de Montréal

Seven-transmembrane (7-TM) receptors are now recognized as complex processors of information that can bind to molecules and cytosolic interactants on the cell membrane. The 7-TM poster highlights the multiple behaviors of 7-TMs including G-protein-dependent and -independent signaling as well as the concept of collateral efficacy.



## **G-protein-coupled Receptors & Signaling Networks**

M. Marinissen, Universidad Autonoma de Madrid  
J. Gutkind, National Institutes of Health

This poster reviews G-protein-coupled receptors (GPCRs), their physiological roles and the cellular responses they mediate. The network of intracellular signaling pathways involved in GPCR function are also explored.



## **P2X and P2Y Receptors**

K. Jacobson, National Institutes of Health

P2X and P2Y receptors are widely distributed in the body. In particular, P2X<sub>1</sub>, P2Y<sub>1</sub> and P2Y<sub>12</sub> have been isolated in platelets. This review covers the subtypes and structures of P2 receptor families and the pharmacological probes used to study them.

To download or request copies, please visit [www.tocris.com/requestliterature](http://www.tocris.com/requestliterature)

# Cardiovascular Research Compounds from Tocris

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>5-HT Receptors</b>				
<i>Agonists</i>	3428	DMT	5-HT <sub>2A</sub> agonist; also endogenous $\alpha_1$ ligand	10 mg
	2643	DOI	Mixed 5-HT <sub>2A/2C</sub> agonist	10 mg 50 mg
	2201	PNU 22394	5-HT <sub>2C</sub> agonist and 5-HT <sub>2A/2B</sub> partial agonist	10 mg 50 mg
	2592	TCB-2	Potent, high affinity 5-HT <sub>2A</sub> agonist	10 mg 50 mg
<i>Antagonists</i>	0523	4F 4PP	Selective 5-HT <sub>2A</sub> antagonist	10 mg 50 mg
	0870	MDL 11,939	5-HT <sub>2A</sub> antagonist	10 mg 50 mg
	1742	R-96544	Potent, selective 5-HT <sub>2A</sub> antagonist	10 mg 50 mg
	2865	Risperidone	5-HT <sub>2A</sub> antagonist	10 mg 50 mg
	3739	Sarpogrelate	Selective 5-HT <sub>2A</sub> antagonist	10 mg 50 mg
<b><math>\alpha_1</math> adrenergic Receptors</b>				
<i>Agonists</i>	0888	Cirazoline	Selective $\alpha_1$ agonist	10 mg
	2838	(R)-(-)-Phenylephrine	$\alpha_1$ agonist	100 mg
<i>Antagonists</i>	2685	Carvedilol	$\alpha_1$ and $\beta$ adrenergic receptor antagonist	50 mg
	0545	Ifenprodil	$\alpha_1$ antagonist; also NMDA antagonist and $\sigma$ ligand	10 mg 50 mg
	0623	Prazosin	$\alpha_1$ and $\alpha_{2B}$ antagonist; MT <sub>3</sub> antagonist	100 mg
<b><math>\alpha_2</math> adrenergic Receptors</b>				
<i>Agonists</i>	0690	Clonidine	$\alpha_2$ agonist. Also I <sub>1</sub> ligand	100 mg
	2749	Dexmedetomidine	Potent, highly selective $\alpha_2$ agonist; active isomer of medetomidine (Cat. No. 2023)	10 mg 50 mg
	0885	Guanabenz	$\alpha_2$ agonist; also I <sub>2</sub> selective ligand	100 mg
	1030	Guanfacine	$\alpha_{2A}$ agonist	10 mg 50 mg
	2638	ST 91	$\alpha_2$ agonist, putative $\alpha_{2C}$ agonist	10 mg 50 mg
	2466	UK 14,304 tartrate	$\alpha_2$ agonist; water-soluble form of UK 14,304 (Cat. No. 0425)	10 mg 50 mg
<i>Antagonists</i>	2666	JP 1302	Potent and selective $\alpha_{2C}$ antagonist	10 mg 50 mg
<b><math>\beta</math> Adrenergic Receptors</b>				
<i>Agonists</i>	0435	Cimaterol	$\beta$ agonist	10 mg 50 mg
	0515	Dobutamine	$\alpha_1$ , $\beta_1$ and $\beta_2$ agonist	50 mg
	1448	Formoterol	Potent and selective $\beta_2$ agonist	10 mg 50 mg
	1747	Isoproterenol	Standard selective $\beta$ agonist	100 mg
<i>Antagonists</i>	2685	Carvedilol	$\alpha_1$ and $\beta$ adrenergic receptor antagonist	50 mg
	1024	CGP 20712	Highly potent and selective $\beta_1$ antagonist	10 mg 50 mg
	0821	ICI 118,551	Very selective $\beta_2$ antagonist	10 mg 50 mg
	0832	ICI 89406	$\beta$ antagonist	10 mg 50 mg
	3256	Metoprolol	Selective $\beta_1$ antagonist	50 mg
	0829	Pronethalol	$\beta$ antagonist	100 mg



## Cardiovascular Research Compounds – continued

Class	Cat. No.	Product Name	Primary Action	Unit Size
	0624	Propranolol	$\beta$ antagonist	100 mg
	0952	Sotalol	$\beta$ antagonist	10 mg 50 mg
<b>Adenosine Receptors</b>				
<i>Agonists</i>	1705	2-Chloro-N <sup>6</sup> -cyclopentyladenosine	Potent, selective A <sub>1</sub> agonist	10 mg 50 mg
	1104	2-Cl-IB-MECA	Highly selective A <sub>3</sub> agonist	10 mg 50 mg
	4472	BAY 60-6583	Potent A <sub>2B</sub> receptor agonist; cardioprotective	10 mg 50 mg
	1063	CGS 21680	A <sub>2A</sub> agonist	10 mg 50 mg
	1066	IB-MECA	A <sub>3</sub> selective agonist	5 mg 25 mg
<i>Antagonists</i>	0439	DPCPX	A <sub>1</sub> selective antagonist	100 mg
	1217	MRS 1220	Highly potent, selective hA <sub>3</sub> antagonist	5 mg 25 mg
	2752	MRS 1754	Selective A <sub>2B</sub> antagonist	10 mg 50 mg
	2009	PSB 1115	Selective human A <sub>2B</sub> receptor antagonist; water-soluble	10 mg 50 mg
	3198	PSB 603	Highly selective A <sub>2B</sub> antagonist	10 mg 50 mg
	2270	SCH 58261	Potent, highly selective A <sub>2A</sub> antagonist	10 mg 50 mg
	1036	ZM 241385	Potent, highly selective A <sub>2A</sub> antagonist	10 mg 50 mg
<b>Aldosterone Receptors</b>				
<i>Antagonists</i>	3281	Canrenone	Mineralocorticoid receptor antagonist	50 mg
	2397	Eplerenone	Selective mineralocorticoid receptor antagonist	10 mg 50 mg
	2970	RU 26752	Mineralocorticoid receptor antagonist	10 mg
	1672	RU 28318	Potent, selective mineralocorticoid receptor antagonist	10 mg 50 mg
	2968	Spironolactone	Mineralocorticoid receptor antagonist	50 mg
<b>Angiotensin-converting Enzyme</b>				
<i>Inhibitors</i>	2578	Benazepril	Angiotensin-converting enzyme (ACE) inhibitor	50 mg
	4455	Captopril	ACE inhibitor; also inhibits LTA <sub>4</sub> hydrolase	50 mg
	2691	Moexipril	Angiotensin-converting enzyme (ACE) inhibitor	10 mg 50 mg
	4302	Perindopril	Angiotensin-converting enzyme (ACE) inhibitor	50 mg
	2931	Spinorphin	Endogenous peptide inhibitor of ACE; also potent P2X <sub>3</sub> antagonist	1 mg
<i>Substrates</i>	1563	Angiotensin I (human, mouse, rat)	Endogenous precursor to angiotensin II (Cat. No. 1158)	1 mg
	1764	Hemopressin (rat)	Endogenous endopeptidase substrate; potent hypotensive <i>in vivo</i>	1 mg
<b>Angiotensin II Receptors</b>				
<i>Agonists</i>	2569	CGP 42112	Selective, high affinity AT <sub>2</sub> ligand	1 mg
	3615	Novokinin	Orally active AT <sub>2</sub> agonist	1 mg
<i>Antagonists</i>	3798	Losartan	Selective, non-peptide AT <sub>1</sub> antagonist	50 mg
	1361	PD 123319	Potent, selective non-peptide AT <sub>2</sub> antagonist	10 mg 50 mg
	4216	Valsartan	High affinity, selective AT <sub>1</sub> antagonist	10 mg 50 mg

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>Apoptosis</b>				
<i>Other</i>	2098	Apoptosis Activator 2	Promotes apoptosome formation and activates caspase-9/caspase-3 pathway; selectively induces tumor cell apoptosis	10 mg 50 mg
	2172	AZ 10417808	Selective non-peptide caspase-3 inhibitor	10 mg 50 mg
	2160	Bax channel blocker	Inhibits Bax-mediated mitochondrial cytochrome c release	10 mg 50 mg
	1786	Bax inhibitor peptide P5	Inhibitor of Bax-mediated apoptosis	1 mg
	1785	Bax inhibitor peptide V5	Inhibitor of Bax-mediated apoptosis	1 mg
	1787	Bax inhibitor peptide, negative control	Negative control peptide for Bax inhibitor peptides V5 and P5 (Cat. Nos. 1785 and 1786)	1 mg
	3590	Gambogic acid	Apoptosis inducer; activates caspases and inhibits Bcl-2 family proteins	10 mg 50 mg
	1541	HA14-1	Bcl-2 inhibitor; induces apoptosis	10 mg 50 mg
	3794	iMAC2	Suppressor of mitochondrial apoptosis	10 mg 50 mg
	2636	Ivachtin	Potent caspase-3 inhibitor	1 mg 10 mg
	2581	PAC 1	Activator of procaspase-3; proapoptotic	10 mg 50 mg
	1758	PETCM	Activator of caspase-3	50 mg
	2775	Probucol	Antioxidant, anti-inflammatory and hypocholesterolemic agent	100 mg
	4038	TW 37	Bcl-2 inhibitor; induces apoptosis	10 mg 50 mg
	2166	Z-DEVD-FMK	Cell-permeable, irreversible caspase-3 inhibitor	1 mg
	2163	Z-VAD-FMK	Cell-permeable, irreversible caspase inhibitor	1 mg
<b>Calcium Channels</b>				
<i>Activators</i>	1544	(±)-Bay K 8644	Ca <sup>2+</sup> -channel activator (L-type)	10 mg 50 mg
	1546	(S)-(-)-Bay K 8644	Ca <sup>2+</sup> -channel activator (L-type)	10 mg 50 mg
	1403	FPL 64176	Potent activator of Ca <sup>2+</sup> channels (L-type)	10 mg 50 mg
<i>Blockers</i>	0685	Diltiazem	Ca <sup>2+</sup> channel blocker (L-type)	1 g
	2004	Isradipine	Ca <sup>2+</sup> channel blocker (L-type)	10 mg 50 mg
	2198	Mibefradil	Ca <sup>2+</sup> channel blocker (T-type)	10 mg 50 mg
	1075	Nifedipine	Ca <sup>2+</sup> channel blocker (L-type)	100 mg
	0600	Nimodipine	Ca <sup>2+</sup> channel blocker (L-type)	100 mg
	2268	NNC 55-0396	Highly selective Ca <sup>2+</sup> channel blocker (T-type)	10 mg
	1439	Ruthenium Red	Non-selective Ca <sup>2+</sup> channel blocker (N- and P-type)	100 mg
	0654	Verapamil	Ca <sup>2+</sup> channel blocker (L-type)	1 g
<b>Calcium Signaling</b>				
<i>Inhibitors</i>	3954	<i>trans</i> -Ned 19	NAADP antagonist; inhibits Ca <sup>2+</sup> release	10 mg
	1329	Ryanodine	Ca <sup>2+</sup> release inhibitor	1 mg
	1147	SKF 96365	STIM1-mediated Ca <sup>2+</sup> influx inhibitor	10 mg 50 mg
	1280	(-)-Xestospongine C	Inhibits IP <sub>3</sub> -mediated Ca <sup>2+</sup> release	10 µg

## Cardiovascular Research Compounds – continued

Class	Cat. No.	Product Name	Primary Action	Unit Size
Other	1234	A23187	Calcium ionophore	10 mg
	2786	BAPTA	Selective calcium chelator	100 mg
	2787	BAPTA AM	Cell-permeable Ca <sup>2+</sup> chelator	25 mg
	2220	FURA-2AM	Fluorescent Ca <sup>2+</sup> indicator	1 mg
	1704	Ionomycin calcium salt	Calcium ionophore	1 mg
	2092	Ionomycin free acid	Calcium ionophore	1 mg
<b>Cell Adhesion Molecules</b>				
Antagonists	1263	GR 144053	Glycoprotein IIb/IIIa (integrin $\alpha_{IIb}\beta_3$ ) receptor antagonist; antithrombotic	10 mg 50 mg
	2524	A 205804	Selective inhibitor of E-selectin and ICAM-1 expression	10 mg 50 mg
Inhibitors	4228	A 286982	Potent inhibitor of the LFA-1/ICAM-1 interaction	10 mg 50 mg
	3202	Echistatin, $\alpha 1$ isoform	$\alpha_v\beta_3$ and glycoprotein IIb/IIIa (integrin $\alpha_{IIb}\beta_3$ ) inhibitor	100 $\mu$ g
	2748	KF 38789	Selective inhibitor of P-selectin-mediated cell adhesion	10 mg
	2877	MNS	Selective inhibitor of Src and Syk	50 mg
	2710	OGT 2115	Antiangiogenic; heparanase inhibitor	1 mg 10 mg
	4227	RWJ 50271	Inhibitor of LFA-1/ICAM mediated cell adhesion	10 mg 50 mg
	2812	Heparin	Anticoagulant	100 mg
Other	4034	PM 102	Antagonist of heparin (Cat. No. 2812)	1 mg
<b>Chemokine Receptors</b>				
Antagonists	3299	AMD 3100	Highly selective CXCR4 antagonist	10 mg 50 mg
	2423	DAPTA	Chemokine receptor 5 (CCR5) antagonist	1 mg
	2595	J 113863	Potent CCR1 chemokine receptor antagonist	1 mg 10 mg 50 mg
	3756	Maraviroc	Selective CCR5 antagonist	10 mg 50 mg
	2725	SB 225002	Potent and selective CXCR2 antagonist	10 mg 50 mg
	2724	SB 265610	Potent CXCR2 antagonist	1 mg 10 mg 50 mg
<b>Cholesterol Regulation</b>				
Inhibitors	3776	Atorvastatin	HMG-CoA reductase inhibitor	10 mg 50 mg
	1639	AY 9944	Inhibitor of hedgehog (Hh) signaling; inhibits $\Delta^7$ -dehydrocholesterol reductase	10 mg
	2227	CI 976	Acyl-CoA:cholesterol acyltransferase (ACAT) inhibitor	10 mg 50 mg
	3540	Orlistat	Pancreatic, gastric and carboxylester lipase inhibitor; antiobesity and antihypercholesterolemic activity	10 mg 50 mg
	2775	Probucol	Antioxidant, anti-inflammatory and hypocholesterolemic agent	100 mg
	1965	Simvastatin	HMG-CoA reductase inhibitor	50 mg
	4184	Torcetrapib	Inhibitor of cholesteryl ester transfer protein (CETP)	10 mg 50 mg
	1638	U 18666A	Inhibitor of hedgehog (Hh) signaling; inhibits cholesterol synthesis	10 mg
	3039	YM 750	Acyl-CoA:cholesterol acyltransferase (ACAT) inhibitor	10 mg 50 mg
<b>Cyclic GMP</b>				
Activators	1089	8-Bromo-cGMP	cGMP analog; activates PKG	10 mg 50 mg

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>Cyclooxygenases</b>				
<i>Inhibitors</i>	3786	Celecoxib	Selective cyclooxygenase-2 (COX-2) inhibitor	10 mg 50 mg
	4454	Diclofenac	Cyclooxygenase inhibitor; NSAID	50 mg
	1418	Resveratrol	Cyclooxygenase inhibitor	100 mg
	4206	Valdecoxib	Selective and potent COX-2 inhibitor	10 mg 50 mg
<b>Cytokine Receptors</b>				
<i>Antagonists</i>	1793	AF 12198	Potent, selective human type I IL-1 receptor antagonist	1 mg
	2265	Lyn peptide inhibitor	Inhibits Lyn-dependent activities of IL-5 receptor; cell-permeable	1 mg
<b>Elastases</b>				
<i>Inhibitors</i>	3535	Sivelestat	Selective leukocyte elastase inhibitor	10 mg 50 mg
	2506	SSR 69071	Potent, orally active human leukocyte elastase inhibitor	10 mg
<b>Endothelin Receptors</b>				
<i>Agonists</i>	1189	BQ-3020	Selective ET <sub>B</sub> agonist	500 µg
	1160	Endothelin 1 (human, porcine)	Potent vasoconstrictor peptide	100 µg
	1899	Sarafotoxin S6a	Endothelin receptor agonist	100 µg
<i>Antagonists</i>	1441	BMS 182874	Highly selective, orally active non-peptide ET <sub>A</sub> antagonist	10 mg 50 mg
	1500	BQ 788	Selective ET <sub>B</sub> antagonist	1 mg
	1188	BQ-123	Selective ET <sub>A</sub> antagonist	500 µg
	1210	FR 139317	Highly potent, selective ET <sub>A</sub> antagonist	10 mg 50 mg
<b>Fatty Acid Oxidation</b>				
<i>Other</i>	4539	(R)-(+)-Etomoxir	Carnitine palmitoyltransferase I (CPT1) inhibitor	10 mg 50 mg
	0548	(±)-Lauroylcarnitine	Intermediate in lipid metabolism	50 mg
	0567	(±)-Myristoylcarnitine	Intermediate in lipid metabolism	50 mg
	0605	(±)-Octanoylcarnitine	Intermediate in lipid metabolism	50 mg
	0611	(±)-Propionylcarnitine	Intermediate in lipid metabolism	50 mg
	1484	Oleylethanolamide	GPR55 agonist; also PPARα agonist	10 mg 50 mg
	3118	Ranolazine	Antianginal; activates pyruvate dehydrogenase	50 mg
<b>GRK2</b>				
<i>Antagonists</i>	3594	GRK2i	GRK2 inhibitory polypeptide; G <sub>βγ</sub> antagonist	1 mg
<b>IGF-1 Receptor</b>				
<i>Inhibitors</i>	2768	PQ 401	IGF1R inhibitor	10 mg 50 mg
	2956	Picropodophyllotoxin	Selective IGF1R inhibitor	10 mg
<b>Lipoxygenases</b>				
<i>Inhibitors</i>	3541	BAY-X 1005	Orally active 5-lipoxygenase activating protein (FLAP) inhibitor	1 mg
	1311	MK 886	Inhibitor of 5-lipoxygenase-activating protein (FLAP)	10 mg 50 mg
	2850	PD 146176	Selective 15-lipoxygenase inhibitor	10 mg 50 mg
	0645	2-TEDC	5-, 12-, 15-Lipoxygenase inhibitor	10 mg 50 mg
	3308	Zileuton	Orally active 5-LOX inhibitor	10 mg 50 mg



## Cardiovascular Research Compounds – continued

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>Matrix Metalloproteinases</b>				
<i>Inhibitors</i>	2961	Batimastat	Potent, broad spectrum MMP inhibitor	1 mg 10 mg
	2631	Marimastat	Broad spectrum MMP inhibitor	1 mg 10 mg
	2628	ONO 4817	Broad spectrum MMP inhibitor	10 mg
	2916	Ro 32-3555	Potent, collagenase-selective MMP inhibitor	10 mg
	4187	UK 356618	Potent and selective MMP-3 inhibitor	10 mg
	2633	WAY 170523	Potent and selective inhibitor of MMP-13	1 mg 10 mg
<b>Mitochondrial Calcium Uniporter</b>				
<i>Inhibitors</i>	3603	Kaempferol	Mitochondrial Ca <sup>2+</sup> uniporter (MCU) activator; proapoptotic	50 mg
	1244	KB-R7943	MCU inhibitor; also inhibits Na <sup>+</sup> /Ca <sup>2+</sup> exchange	10 mg 50 mg
<b>Mitochondrial Permeability Transition Pore</b>				
<i>Inhibitors</i>	1101	Cyclosporin A	Calcineurin inhibitor	100 mg
	4110	Oligomycin A	Inhibitor of mitochondrial ATPase	5 mg
	2906	TRO 19622	Binds voltage-dependent anion channel (VDAC)	10 mg 50 mg
<b>Muscarinic Receptors (mAChRs)</b>				
<i>Antagonists</i>	1105	AF-DX 116	Selective M <sub>2</sub> antagonist	10 mg 50 mg
	1345	AF-DX 384	Potent M <sub>2</sub> /M <sub>4</sub> antagonist	10 mg 50 mg
	0482	4-DAMP	Muscarinic M <sub>3</sub> antagonist	50 mg
	2096	DAU 5884	M <sub>3</sub> receptor antagonist	10 mg 50 mg
	2507	J 104129	Potent, selective M <sub>3</sub> antagonist	10 mg
<b>Na<sup>+</sup>/Ca<sup>2+</sup> Exchanger</b>				
<i>Inhibitors</i>	4117	Bepidil	Nonselective calcium channel blocker	50 mg
	1114	CGP 37157	Antagonist of mitochondrial Na <sup>+</sup> /Ca <sup>2+</sup> exchange	10 mg 50 mg
	1244	KB-R7943	Na <sup>+</sup> /Ca <sup>2+</sup> exchange inhibitor (reverse mode)	10 mg 50 mg
	2184	SN-6	Selective Na <sup>+</sup> /Ca <sup>2+</sup> exchange inhibitor (reverse mode)	10 mg 50 mg
<b>Na<sup>+</sup>/H<sup>+</sup> Exchanger</b>				
<i>Inhibitors</i>	0890	Amiloride	Na <sup>+</sup> channel blocker; also I <sub>2</sub> imidazoline ligand	100 mg
	2727	Zoniporide	Selective NHE1 inhibitor	10 mg 50 mg
<b>Na<sup>+</sup>/K<sup>+</sup> ATPase</b>				
<i>Inhibitors</i>	4583	Digoxin	Na <sup>+</sup> /K <sup>+</sup> ATPase inhibitor	50 mg
	1076	Ouabain	Na <sup>+</sup> /K <sup>+</sup> ATPase inhibitor	100 mg
<b>Natriuretic Peptide Receptors</b>				
<i>Agonists</i>	1912	Atrial natriuretic factor (1-28) (rat)	Endogenous peptide regulating blood pressure	1 mg
	1906	Atrial natriuretic factor (1-28) (human, porcine)	Endogenous peptide regulating blood pressure	1 mg
	3520	C-type natriuretic factor peptide (1-22) (human, rat, swine)	Endogenous peptide agonist at NPR2	500 µg

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>Nitric Oxide</b>				
<i>Donors</i>	2147	Nicorandil	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel opener and NO donor	50 mg
	0756	SIN-1	Water-soluble NO donor	50 mg
	0603	SNOG	NO carrier, breaks down to release NO	10 mg 50 mg
	1135	Spermine NONOate	Slow NO releasing agent	10 mg 50 mg
<i>Inhibitors</i>	0546	L-NIO	Potent eNOS inhibitor	10 mg 50 mg
<i>Other</i>	0598	SNAP	A stable analog of endogenous S-nitroso compounds	10 mg 50 mg
<i>Substrates</i>	0722	<i>N</i> -Acetyl- <i>N</i> -acetoxy-4-chlorobenzenesulfonamide	Nitroxyl precursor	10 mg 50 mg
	0663	L-Arginine	Endogenous substrate for NOS	100 mg
<b>NKCC Cotransporter</b>				
<i>Inhibitors</i>	3108	Bumetanide	Na <sup>+</sup> /2Cl <sup>-</sup> /K <sup>+</sup> (NKCC) cotransporter inhibitor	50 mg
	3109	Furosemide	Na <sup>+</sup> /2Cl <sup>-</sup> /K <sup>+</sup> (NKCC) cotransporter inhibitor; also antagonizes GABA <sub>A</sub>	50 mg
<b>Oxidative Phosphorylation</b>				
<i>Inhibitors</i>	0452	CCCP	Oxidative phosphorylation uncoupler	500 mg
	3612	Enterostatin	Binds to β-subunit of F <sub>1</sub> -ATPase; anorexigenic peptide	1 mg
	0453	FCCP	Oxidative phosphorylation uncoupler	10 mg 50 mg
	3616	Rotenone	Inhibits complex I of the mitochondrial electron transport chain	50 mg
<b>PDGF Receptors</b>				
<i>Inhibitors</i>	4274	AP 24534	Potent multi-kinase and pan-BCR-ABL inhibitor	10 mg 50 mg
	1222	DMPQ	Potent, selective inhibitor of PDGFRβ	10 mg 50 mg
	3785	PD 166285	Potent Src inhibitor; also inhibits FGFR1, PDGFRβ and Wee1	1 mg 10 mg
	3304	SU 16f	Potent and selective PDGFRβ inhibitor	10 mg 50 mg
	3335	SU 6668	PDGFR, VEGFR and FGFR inhibitor	10 mg 50 mg
	3768	Sunitinib	Potent VEGFR, PDGFRβ and KIT inhibitor	10 mg 50 mg
<b>Phosphodiesterases</b>				
<i>Inhibitors</i>	0691	Dipyridamole	PDE inhibitor; coronary vasodilator	500 mg
	3053	Mesopram	Orally active PDE4 inhibitor	10 mg 50 mg
	1349	( <i>R</i> )-(-)-Rolipram	PDE4 inhibitor; more active enantiomer of rolipram (Cat. No. 0905)	10 mg 50 mg
	3784	Sildenafil	Orally active, potent PDE5 inhibitor	10 mg 50 mg
	1676	T 0156	Highly potent, selective PDE5 inhibitor	10 mg 50 mg
	1046	Zardaverine	PDE3/4 inhibitor	10 mg 50 mg

## Cardiovascular Research Compounds – continued

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>Phospholipases</b>				
<i>Inhibitors</i>	1462	AACOCF <sub>3</sub>	Phospholipase A <sub>2</sub> inhibitor	5 mg 25 mg
	1437	D609	Selective PC-PLC inhibitor	10 mg 50 mg
	3022	Edelfosine	Selective PI-PLC inhibitor; also PAF receptor agonist	10 mg
	1941	<i>m</i> -3M3FBS	Phospholipase C activator	10 mg
	1942	<i>o</i> -3M3FBS	Inactive analog of <i>m</i> -3M3FBS (Cat. No. 1941)	10 mg
	0606	OBAA	Phospholipase A <sub>2</sub> inhibitor	10 mg 50 mg
	1268	U 73122	Phospholipase C inhibitor	10 mg 50 mg
	4133	U 73343	Inactive analog of U 73122 (Cat. No. 1268)	10 mg 50 mg
<b>PI 3-Kinase</b>				
<i>Activators</i>	1983	740 Y-P	Cell-permeable PI 3-kinase activator	1 mg
<i>Inhibitors</i>	3578	AS 605240	Potent and selective PI 3-kinase $\gamma$ (PI3K $\gamma$ ) inhibitor	10 mg 50 mg
	3606	BAG 956	Dual PI 3-kinase and PDK1 inhibitor	10 mg 50 mg
	2814	PI 828	PI 3-kinase inhibitor	1 mg 10 mg 50 mg
<b>Potassium Channels</b>				
<i>Activators</i>	1377	Cromakalim	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel opener	10 mg 50 mg
	1378	Levcromakalim	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel opener; active enantiomer of cromakalim (Cat. No. 1377)	10 mg 50 mg
	0583	Minoxidil	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel opener	100 mg
	4519	ML 213	K <sub>v</sub> 7.2 and K <sub>v</sub> 7.4 channel opener	10 mg 50 mg
	2147	Nicorandil	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel opener and NO donor	50 mg
	4462	NS 3623	K <sub>v</sub> 11.1 (hERG) channel activator; antiarrhythmic	10 mg 50 mg
	1355	P1075	Potent K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel opener	10 mg 50 mg
<i>Blockers</i>	2533	DPO-1	Blocker of K <sub>v</sub> 1.5 channels; prevents atrial arrhythmia	10 mg 50 mg
	1808	E-4031	K <sub>v</sub> 11.1 (hERG) channel blocker; class III antiarrhythmic agent	10 mg 50 mg
	0911	Glibenclamide	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel blocker	100 mg
	2396	Glimepiride	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel blocker	10 mg 50 mg
	3899	JNJ 303	Potent and selective I <sub>Ks</sub> blocker	10 mg 50 mg
	4231	Nateglinide	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) blocker; displays high affinity for SUR1/K <sub>ir</sub> 6.2 channels	10 mg 50 mg
	2095	PNU 37883	Vascular K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel blocker	10 mg 50 mg
	3805	Repaglinide	K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel blocker	50 mg
	3948	Terfenadine	K <sub>v</sub> 11.1 (hERG) and K <sub>ir</sub> 6 (K <sub>ATP</sub> ) channel blocker; also H <sub>1</sub> receptor antagonist	50 mg

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>PPAR</b>				
<i>Agonists</i>	1307	Ciglitazone	Selective PPAR $\gamma$ agonist	10 mg 50 mg
	2229	GW 0742	Highly selective, potent PPAR $\delta$ agonist	10 mg 50 mg
	1664	GW 1929	Selective PPAR $\gamma$ agonist; orally active	10 mg 50 mg
	1677	GW 7647	Highly selective, potent PPAR $\alpha$ agonist; orally active	10 mg 50 mg
	4124	Pioglitazone	Selective PPAR $\gamma$ agonist; antidiabetic agent	10 mg 50 mg
	3114	Troglitazone	Selective PPAR $\gamma$ agonist; antidiabetic agent	10 mg 50 mg
<i>Antagonists</i>	3961	GSK 3787	Potent and selective PPAR $\delta$ antagonist	10 mg 50 mg
	4618	GW 6471	PPAR $\alpha$ antagonist	10 mg 50 mg
<b>Prostanoid Receptors</b>				
<i>Agonists</i>	1442	BMV 45778	Non-prostanoid prostacyclin IP receptor partial agonist	10 mg 50 mg
	2989	Epoprostenol	Endogenous IP receptor agonist	10 mg
	1932	U 46619	Potent, stable thromboxane A <sub>2</sub> (TP) receptor agonist	1 mg
<i>Antagonists</i>	0671	AH 6809	EP <sub>1</sub> and EP <sub>2</sub> receptor antagonist	10 mg 50 mg
	2514	L-161,982	Selective EP <sub>4</sub> receptor antagonist	10 mg
	3342	L-798,106	Potent and highly selective EP <sub>3</sub> antagonist	10 mg 50 mg
<i>Other</i>	1620	Alprostadil	Prostaglandin; vasodilator and antiplatelet agent <i>in vivo</i>	10 mg 50 mg
	2296	Prostaglandin E <sub>2</sub>	Major endogenous prostanoid	10 mg
	4214	Prostaglandin F <sub>2<math>\alpha</math></sub>	Naturally-occurring prostanoid; potent vasoconstrictor	10 mg
<b>Protease-activated Receptors</b>				
<i>Agonists</i>	1464	TFLLR-NH <sub>2</sub>	PAR <sub>1</sub> -activating peptide	1 mg
	3497	TRAP-6	PAR <sub>1</sub> peptide fragment (residues 42-47); acts as a PAR <sub>1</sub> agonist	5 mg
<i>Antagonists</i>	3643	FR 171113	PAR <sub>1</sub> antagonist	10 mg 50 mg
	2614	RWJ 56110	Selective PAR <sub>1</sub> antagonist	1 mg
	1592	SCH 79797	Potent, selective non-peptide PAR <sub>1</sub> antagonist	10 mg 50 mg
	1488	tcY-NH <sub>2</sub>	Selective PAR <sub>4</sub> antagonist	1 mg
<i>Other</i>	3393	RLLFT-NH <sub>2</sub>	Control peptide for TFLLR-NH <sub>2</sub> (Cat. No. 1464)	1 mg
	1185	Thrombin Receptor Agonist Peptide	Causes platelet aggregation and secretion	1 mg
<b>Purinergic P2 Receptors</b>				
<i>Agonists</i>	3312	BzATP	P2X <sub>7</sub> agonist; also P2X <sub>1</sub> and P2Y <sub>1</sub> partial agonist	1 mg
	2157	MRS 2365	Highly potent and selective P2Y <sub>1</sub> agonist	1 mg
	2915	MRS 2690	Potent P2Y <sub>14</sub> agonist	1 mg
<i>Antagonists</i>	3321	AR-C 66096	Potent and selective P2Y <sub>12</sub> antagonist	1 mg
	2490	( $\pm$ )-Clopidogrel	Selective P2Y <sub>12</sub> antagonist	10 mg 50 mg
	0900	MRS 2179	Selective P2Y <sub>1</sub> antagonist	10 mg 50 mg
	2159	MRS 2500	Extremely potent and selective P2Y <sub>1</sub> antagonist	1 mg
	1240	NF 023	Selective, competitive P2X <sub>1</sub> antagonist	10 mg 50 mg



## Cardiovascular Research Compounds – continued

Class	Cat. No.	Product Name	Primary Action	Unit Size
	3983	PSB 0739	Highly potent P2Y <sub>12</sub> receptor antagonist	10 mg 50 mg
	1472	Suramin	Non-selective P2 antagonist	100 mg
	3931	Ticlopidine	Selective P2Y <sub>12</sub> antagonist	50 mg
	2464	TNP-ATP	Potent, selective P2X antagonist	5 mg
<b>Rho-kinase</b>				
<i>Inhibitors</i>	0541	Fasudil	Inhibitor of cyclic nucleotide dependent- and Rho-kinases	10 mg 50 mg
	4009	GSK 269962	Potent and selective ROCK inhibitor	10 mg 50 mg
	2414	H 1152	Selective Rho-kinase (ROCK) inhibitor	1 mg
	4118	SB 772077B	Potent Rho-kinase inhibitor; vasodilator	10 mg 50 mg
	1254	Y-27632	Selective p160ROCK inhibitor	1 mg 10 mg 50 mg
<b>SERCA</b>				
<i>Inhibitors</i>	1235	Cyclopiazonic acid	Inhibitor of SERCA	10 mg 50 mg
	2006	Paxilline	SERCA blocker; also potent BK <sub>Ca</sub> channel blocker	10 mg
	1138	Thapsigargin	Potent inhibitor of SERCA	1 mg
<b>Sodium Channels</b>				
<i>Activators</i>	2918	Veratridine	Voltage-gated Na <sup>+</sup> channel opener	10 mg 50 mg
<i>Blockers</i>	1470	Flecainide	Cardiac Na <sup>+</sup> channel blocker; antiarrhythmic	10 mg 50 mg
	3251	KC 12291	Orally active atypical Na <sup>+</sup> blocker; cardioprotective	10 mg 50 mg
	1014	QX 314	Na <sup>+</sup> channel blocker	100 mg
	2313	QX 314 chloride	Na <sup>+</sup> channel blocker	50 mg
	4435	TC-N 1752	Selective Na <sub>v</sub> 1.7 blocker	10 mg 50 mg
	1078	Tetrodotoxin	Na <sup>+</sup> channel blocker	1 mg
	1069	Tetrodotoxin citrate	Na <sup>+</sup> channel blocker; citrate salt of tetrodotoxin (Cat. No. 1078)	1 mg
<b>Soluble Guanylyl Cyclase</b>				
<i>Activators</i>	2753	A 350619	Soluble guanylyl cyclase (sGC) activator	10 mg 50 mg
	4430	BAY 41-2272	Soluble guanylyl cyclase (sGC) activator	10 mg 50 mg
<i>Inhibitors</i>	4517	NS 2028	Potent soluble guanylyl cyclase (sGC) inhibitor	10 mg 50 mg
	0880	ODQ	Selective inhibitor of NO-sensitive guanylyl cyclase	10 mg 50 mg
<b>Stem Cells</b>				
<i>Other</i>	3842	5-Azacytidine	DNA methyltransferase inhibitor; induces cardiomyogenesis in MSCs	50 mg
	3851	Cardiogenol C	Induces cardiomyogenesis in ESCs	10 mg 50 mg
	3748	XAV 939	Inhibits Wnt signaling; promotes cardiomyogenesis	10 mg 50 mg
	2293	Zebularine	DNA methyltransferase inhibitor; induces cardiomyogenesis in MSCs	10 mg

Class	Cat. No.	Product Name	Primary Action	Unit Size
<b>TGF-<math>\beta</math> Receptors</b>				
<i>Inhibitors</i>	2939	A 83-01	Selective inhibitor of TGF- $\beta$ RI, ALK4 and ALK7	10 mg 50 mg
	3264	GW 788388	Selective inhibitor of TGF- $\beta$ RI	10 mg 50 mg
	2718	LY 364947	Selective inhibitor of TGF- $\beta$ RI	1 mg 10 mg
	1614	SB 431542	Potent, selective inhibitor of TGF- $\beta$ RI, ALK4 and ALK7	1 mg 10 mg
	3263	SB 505124	Selective inhibitor of TGF- $\beta$ RI, ALK4 and ALK7	10 mg 50 mg
	3211	SB 525334	Selective inhibitor of TGF- $\beta$ RI	10 mg 50 mg
	3269	SD 208	Potent ATP-competitive TGF- $\beta$ RI inhibitor	10 mg 50 mg
	3742	SJN 2511	Selective inhibitor of TGF- $\beta$ RI	10 mg 50 mg
<b>Urokinase</b>				
<i>Inhibitors</i>	4372	BC 11	Selective urokinase (uPA) inhibitor	10 mg 50 mg
	0442	4-Chlorophenylguanidine	Urokinase inhibitor	100mg
<b>Urotensin II</b>				
<i>Agonists</i>	2484	( $\pm$ )-AC 7954	Non-peptide UT receptor agonist	10 mg 50 mg
	1642	Urotensin II (human)	Endogenous vasoactive agonist for the UT receptor	1 mg
<i>Antagonists</i>	3162	[Orn <sup>5</sup> ]-URP	Urotensin-II (UT) receptor antagonist	1 mg
	1839	BIM 23127	NMB receptor antagonist; also UT receptor antagonist	1 mg
	3571	SB 657510	Selective urotensin-II (UT) receptor antagonist	10 mg 50 mg

# Further Reading

Please refer to the list of recommended papers for more information.

## Hypertension

- Harvey and Belevych** (2003) Muscarinic regulation of cardiac ion channels. *Br.J.Pharmacol.* **139** 1074
- Humbert et al** (2004) Cellular and molecular pathobiology of pulmonary arterial hypertension. *J.Am.Coll.Cardiol.* **43** 13S
- Rabinovitch** (2008) Molecular pathogenesis of pulmonary arterial hypertension. *J.Clin.Invest.* **118** 2372
- Unger et al** (2011) Therapeutic perspectives in hypertension: novel means for renin-angiotensin-aldosterone system modulation and emerging device-based approaches. *Eur.Heart J.* **32** 2739
- Castrillo and Tontonoz** (2004) PPARs in atherosclerosis: the clot thickens. *J.Clin.Invest.* **114** 1538

## Atherosclerosis

- Glass and Witztum** (2001) Atherosclerosis: the road ahead. *Cell* **104** 503
- Shiraishi et al** (2008) Chronic urotensin II infusion enhances macrophage foam cell formation and atherosclerosis in apolipoprotein E-knockout mice. *J.Hypertens.* **26** 1955
- Zernecke et al** (2008) Chemokines in atherosclerosis: an update. *Arterioscler.Thromb.Vasc.Biol.* **28** 1897

## Thrombosis and Hemostasis

- Born and Patrono** (2006) Antiplatelet drugs. *Br.J.Pharmacol.* **147** S241
- Michelson** (2011) Advances in antiplatelet therapy. *Hematology Am.Soc.Hematol.Educ.Program* **1** 62
- Clemetson and Clemetson** (1995) Platelet GPIb-V-IX complex. Structure, function, physiology, and pathology. *Semin.Thromb.Hemost.* **21** 130
- Jennings** (2009) Mechanisms of platelet activation: need for new strategies to protect against platelet-mediated atherothrombosis. *Thromb.Haemost.* **102** 248

## Myocardial Infarction

- Rakhit and Marber** (2001) Nitric oxide: an emerging role in cardioprotection? *Heart* **86** 368
- Mubagwa and Flameng** (2001) Adenosine, adenosine receptors and myocardial protection: an updated overview. *Cardiovasc.Res.* **52** 25

## Arrhythmia

- Grant** (2009) Cardiac ion channels. *Circ.Arrhythm.Electrophysiol.* **2** 185
- Sanguinetti and Tristani-Firouzi** (2006) hERG potassium channels and cardiac arrhythmia. *Nature* **440** 463

## Myocardial I/R Injury

- Halestrap et al** (2004) Mitochondrial permeability transition pore opening during myocardial reperfusion – a target for cardioprotection. *Cardiovasc.Res.* **61** 372
- Headrick and Lasley** (2009) Adenosine receptors and reperfusion injury of the heart. *Handb.Exp.Pharmacol.* **193** 189
- Murphy and Steenbergen** (2008) Mechanisms underlying acute protection from cardiac ischemia-reperfusion injury. *Physiol.Rev.* **88** 581
- Zorov et al** (2009) Regulation and pharmacology of the mitochondrial permeability transition pore. *Cardiovasc.Res.* **83** 213
- Auchampach and Bolli** (1999) Adenosine receptor subtypes in the heart: therapeutic opportunities and challenges. *Am.J.Physiol.* **276** H1113

## Heart Failure

- Lymeropoulos** (2011) GRK2 and  $\beta$ -arrestins in cardiovascular disease: something old, something new. *Am.J.Cardiovasc.Dis.* **1** 126
- Creemers et al** (2001) Matrix metalloproteinase inhibition after myocardial infarction. A new approach to prevent heart failure? *Circ.Res.* **89** 201

**R&D** SYSTEMS

**NOVUS**  
BIOLOGICALS

**TOCRIS**

protein**simple**

**biotechne**

Global info@bio-techne.com bio-techne.com/find-us/distributors TEL +1 612 379 2956  
North America TEL 800 343 7475 Europe | Middle East | Africa TEL +44 (0)1235 529449  
China info.cn@bio-techne.com TEL +86 (21) 52380373

bio-techne.com

