

# P2X and P2Y Receptors

## Kenneth A. Jacobson

Molecular Recognition Section, Laboratory of Bioorganic Chemistry, National Institute of Diabetes and Digestive and Kidney Diseases, National Institutes of Health, Bethesda, Maryland 20892, USA. E-mail: kajacobs@helix.nih.gov

Kenneth Jacobson serves as Chief of the Laboratory of Bioorganic Chemistry and the Molecular Recognition Section at the National Institute of Diabetes and Digestive and Kidney Diseases, National Institutes of Health in Bethesda, Maryland, USA. Dr. Jacobson is a medicinal chemist with interests in the structure and pharmacology of G-protein-coupled receptors, in particular receptors for adenosine and for purine and pyrimidine nucleotides.

## Subtypes and Structures of P2 Receptor Families

The P2 receptors for extracellular nucleotides are widely distributed in the body and participate in regulation of nearly every physiological process.<sup>1,2</sup> Of particular interest are nucleotide receptors in the immune, inflammatory, cardiovascular, muscular, and central and peripheral nervous systems. The ubiquitous signaling properties of extracellular nucleotides acting at two distinct families of P2 receptors – fast P2X ion channels and P2Y receptors (G-protein-coupled receptors) – are now well recognized. These extracellular nucleotides are produced in response to tissue stress and cell damage and in the processes of neurotransmitter release and channel formation. Their concentrations can vary dramatically depending on circumstances. Thus, the state of activation of these receptors can be highly dependent on the stress conditions or disease states affecting a given organ. The P2 receptors respond to various extracellular mono- and dinucleotides (Table 1). The P2X receptors are more structurally restrictive than P2Y receptors in agonist selectivity. P2X receptors are activated principally by ATPs, while the P2Y receptors are activated by a group of five or more naturally occurring nucleotides. The P2X receptors are distributed throughout the nervous system (autonomic, central, enteric and sensory neurons, cochlear and retinal cells), vascular system (cardiomyocytes, endothelium and smooth muscle), the pulmonary and digestive systems (epithelium and visceral smooth muscle), skeletal muscle, bone, and hematopoietic cells. The P2X receptors consist of trimeric ligand-gated ion channels. The subunits are numbered P2X<sub>1</sub> through P2X<sub>7</sub>, and both heterotrimers and homotrimers occur. Activation of P2X receptors leads to influx of cations such as sodium and calcium, which depolarize excitable cells and activate cytosolic enzymes respectively. The P2X<sub>7</sub> receptor upon prolonged agonist exposure also opens a large pore, which can pass organic cations and dye molecules. The knowledge of P2X receptor structures was recently advanced with the X-ray crystallographic determination of the P2X<sub>4</sub> subunit.<sup>3,4</sup> However, this structure did not establish the precise ligand binding site within the protein. A major difficulty in designing new agonist and antagonist ligands for a given P2X receptor subtype is that the homotrimers and heterotrimers may have entirely different structural requirements.

The correspondence of the P2Y receptor subtypes with their native nucleotide ligands is shown in Table 1. The numbering of unique human P2Y receptors has some gaps – due to the fact that the assignment of numbers to certain putative P2Y receptors was later shown to be premature, with some of the previously designated sequences being P2Y species homologs and others being other types of receptors. Each of the native nucleotides may activate several P2Y receptor subtypes. The structures of representative adenine (Figure 1A, 1–10) and uracil (Figure 1B,

## Contents

Subtypes and Structures of P2 Receptor Families.....	1
Pharmacological Probes for P2X Receptors.....	3
Non-Selective P2X Ligands.....	3
P2X <sub>1</sub> and P2X <sub>2</sub> Receptors.....	3
P2X <sub>3</sub> Receptor.....	3
P2X <sub>4</sub> Receptor.....	3
P2X <sub>5</sub> Receptor.....	5
P2X <sub>7</sub> Receptor.....	5
Pharmacological Probes for P2Y Receptors.....	6
P2Y <sub>1</sub> Receptor.....	7
P2Y <sub>2</sub> and P2Y <sub>4</sub> Receptors.....	8
P2Y <sub>6</sub> Receptor.....	9
P2Y <sub>11</sub> Receptor.....	9
P2Y <sub>12</sub> Receptor.....	11
P2Y <sub>13</sub> Receptor.....	12
P2Y <sub>14</sub> Receptor.....	12
Conclusion.....	12
References.....	13
P2X and P2Y Receptor Compounds.....	15

**11–28**) nucleotides that activate P2 receptors are shown. The adenine nucleotide ATP is a full agonist at two human P2Y subtypes (P2Y<sub>2</sub> and P2Y<sub>11</sub> receptors), and the corresponding diphosphate ADP activates three different subtypes (P2Y<sub>1</sub>, P2Y<sub>12</sub>, and P2Y<sub>13</sub> receptors). The uracil nucleotide UTP activates two subtypes (P2Y<sub>2</sub> and P2Y<sub>4</sub> receptors), while UDP, previously thought to activate only a single subtype (P2Y<sub>6</sub> receptors), is now known to also activate P2Y<sub>14</sub> receptors along with the originally designated native agonist UDP-glucose.<sup>5</sup> The naturally occurring dinucleotide Ap<sub>4</sub>A and its homologs also activate various P2 receptors.

The structure, signaling, and regulation of P2Y receptors have been explored, and subfamilies of P2Y<sub>1</sub>-like and P2Y<sub>12</sub>-like receptors have been defined. These subfamilies constitute two pharmacologically distinct groups of P2Y receptors that also correlate with similarities in the function of key amino acid residues.<sup>6</sup> The preferential coupling of the first subfamily of P2Y<sub>1</sub>, P2Y<sub>2</sub>, P2Y<sub>4</sub>, P2Y<sub>6</sub>, and P2Y<sub>11</sub> receptors is to G<sub>q</sub>, leading to activation of phospholipase Cβ (PLCβ), and the second subfamily of P2Y<sub>12</sub>, P2Y<sub>13</sub>, and P2Y<sub>14</sub> receptors couple to G<sub>i</sub> resulting in the inhibition of adenylyl cyclase. P2Y<sub>11</sub> receptors also activate G<sub>s</sub> to stimulate adenylyl cyclase. Comparisons of the structural characteristics and functionally important amino acid residues within the family have been described. Specific cationic residues

and other residues in the TM region (e.g. Phe in TM3) and on the extracellular loops have conserved functions within the P2Y family. Molecular recognition in the P2Y<sub>1</sub>, P2Y<sub>2</sub>, P2Y<sub>11</sub>, and P2Y<sub>12</sub> receptors has been extensively explored using mutagenesis.<sup>7–10</sup>

P2Y receptor regulation has also been studied. In platelets, which express two ADP-responsive P2Y subtypes, the P2Y<sub>1</sub> receptor is more rapidly desensitized than the P2Y<sub>12</sub> receptor.<sup>11</sup> The P2Y<sub>1</sub> receptor is desensitized mainly through PKC-dependent processes, and the P2Y<sub>12</sub> receptor is a good substrate for the GPCR kinases (GRKs) leading to arrestin binding. Residues on the cytosolic C-terminal domain involved in the regulation of the P2Y<sub>1</sub> receptor have been probed. The internalization of the P2Y<sub>11</sub> receptor is dependent on coexpression of the rapidly desensitizing P2Y<sub>1</sub> receptor, suggesting the occurrence of receptor dimers.<sup>12</sup> Various heterodimers of P2Y receptors with other P2Y and non-P2Y GPCRs have been proposed. For example, a dimer of A<sub>1</sub> adenosine receptors and P2Y<sub>1</sub> receptors was characterized.<sup>13</sup>

Recurrent issues in the use of typical P2 receptor ligands include cross-reactivity with multiple P2 receptors and low bioavailability, due to polyanions, such as phosphates and sulfonates, present in the molecules. Another drawback of many of the currently used ligands is lability in biological systems. A large family of ectonucleotidase enzymes hydrolyzes the native nucleotides

**Table 1** | Subtypes of P2 receptors and their ligands (potency at the human homologs shown as *pEC*<sub>50</sub>, unless noted r = rat)

Receptor	Main Distribution	Agonists (native in bold, <i>pEC</i> <sub>50</sub> )	Antagonists
P2X <sub>1</sub>	Smooth muscle, platelets, cerebellum, dorsal horn spinal neurons	BzATP 8.7 > <b>ATP</b> 7.3, 2-MeSATP 7.3, α,β-MeATP 6.7 (rapid desensitization) >> CTP 4.4	NF 449 9.5 > IP <sub>5</sub> I 8.8 > TNP-ATP 8.2 > Ro 0437626 > NF 279 7.7
P2X <sub>2</sub>	Smooth muscle, brain, pancreas, retina, chromaffin cells, autonomic and sensory ganglia	<b>ATP</b> 5.9, 2-MeSATP 5.8 ≥ 2-MeSATP 5.8 >> α,β-MeATP <4	RB2 6.4 (r), iso PPADS 6.4 (r) > PPADS 5.4 (r) > Suramin 4.5 (r)
P2X <sub>3</sub>	Nociceptive sensory neurons, NTS, some sympathetic neurons	<b>ATP</b> 6.5, 2-MeSATP 6.5 ≥ Ap <sub>4</sub> A 6.3, α,β-MeATP 6.1 (rapid desensitization)	TNP-ATP 9.0, iso PPADS 7.1 > A317491 7.6 > NF 110 7.4 > PPADS 5.8
P2X <sub>4</sub>	Microglia, testis, colon, endothelial cells	<b>ATP</b> 6.3 >> α,β-MeATP 5.1 >> CTP 3.5, Ivermectin (potentiates) 6.6	5-BDBD >> TNP-ATP 4.8, PPADS 4.6 > BBG 3.9 (r)
P2X <sub>5</sub>	Proliferating cells in skin, gut, bladder, thymus, spinal cord, heart, adrenal medulla	ATP <sub>γ</sub> S 6.2 (r), <b>ATP</b> 6.0 >> α,β-MeATP <5.2	BBG 6.3 > PPADS 5.6, Suramin 5.4
P2X <sub>6</sub>	Brain, motor neurons in spinal cord	(no functional homomultimer)	–
P2X <sub>7</sub>	Macrophages, mast cells, microglia, pancreas, skin, endocrine organs	BzATP 5.3 > <b>ATP</b> 4.0 ≥ 2-MeSATP >4 >> α,β-MeATP >4	KN 62 7.5, BBG 8.0 (r)
P2Y <sub>1</sub>	Brain, epithelial and endothelial cells, platelets, immune cells, osteoclasts	MRS 2365 9.4 > 2-MeSADP 8.2 >> ADPβS 7.0 > <b>ADP</b> 5.1 > ATP	MRS 2500 9.0 > MRS 2279 7.3 > MRS 2179 6.5
P2Y <sub>2</sub>	Immune cells, epithelial and endothelial cells, kidney tubules, osteoblasts	<b>UTP</b> 8.1, MRS 2698 8.1 ≥ <b>ATP</b> 7.1, INS 365 7.0 > INS 37217 6.7, UTP <sub>γ</sub> S 6.6 > Ap <sub>4</sub> A 6.1 > MRS 2768 5.7	AR-C 126313 6 > Suramin 4.3 > RB2 >4
P2Y <sub>4</sub>	Endothelial cells, placenta	2'-azido-dUTP 7.1 > UTP <sub>γ</sub> S 5.8, <b>UTP</b> 5.6 ≥ <b>ATP</b> 5.7 (rat), > Ap <sub>4</sub> A 5.5 > CTP 5.2, ITP 5.1	ATP (human) 4.4 > RB2 >4 > Suramin >4
P2Y <sub>6</sub>	Airway and intestinal epithelial cells, spleen, placenta, T-cells, thymus	MRS 2693 7.8 > UDPβS 7.3, PBS 0474 7.2 > INS 48823 6.9, Up <sub>3</sub> U 6.6, <b>UDP</b> 6.5 > UTP >> ATP	MRS 2578 7.4 (noncompetitive) > RB2, PPADS
P2Y <sub>11</sub>	Spleen, intestine, granulocytes	ATP <sub>γ</sub> S 5.5 > ARC 67085MX 5.0 > BzATP 5.1 ≥ <b>ATP</b> 4.8	NF 157 7.4 > Suramin 4.8 > RB2 >4
P2Y <sub>12</sub>	Platelets, brain (glial cells), microglial cells	2-MeSADP 7.9 ≥ <b>ADP</b> 7.2	ARC 69931MX 8.4 > AZD 6140 7.9, INS 50589 7.8 > RB2 7.6 (r) > 2-MeSAMP 4.0
P2Y <sub>13</sub>	Spleen, brain, lymph nodes, bone marrow	<b>ADP</b> 7.9 = 2-MeSADP 7.9 > 2-MeSATP 7.1, ATP 6.6	ARC 69931MX 8.4 > ARC 67085 6.7 > MRS 2211 6.0
P2Y <sub>14</sub>	Placenta, mast cells, adipose tissue, stomach, intestine, discrete brain regions	MRS 2690 7.3 > <b>UDP</b> 6.8, <b>UDPglucose</b> 6.5 > UDP-galactose 6.2	–

leading to complications in interpretation of biological results. Adenine nucleotides are progressively converted enzymatically, in the last step by the action of CD73/5'-nucleotidase on AMP, to form adenosine, which activates its own family of four receptors. Selective inhibitors of ectonucleotidases which can serve as modulators of receptor function are being explored.<sup>14</sup> Moreover, many P2 receptor agonists and antagonists are known to inhibit ectonucleotidases at comparable concentrations. Known P2 antagonists often interact intracellularly with other signaling mediators, such as G proteins.

One reason for the relatively slow progress in identifying competitive antagonists of the P2 receptors is that there are few selective radioligands available for either the P2X or P2Y receptors. Previously, various radioactive nucleotides have been suggested to bind to particular P2 receptors, but many of these reports were later questioned, and currently only the P2Y<sub>1</sub>, P2Y<sub>12</sub>, P2X<sub>1</sub> and P2X<sub>3</sub> receptors have viable radioligands.<sup>15-17</sup> Thus, improved and more versatile affinity probes for the P<sub>2</sub> receptors are still needed. New selective agonists and antagonists have recently been identified for some of the eight mammalian subtypes of P2Y receptors and for a few of the seven mammalian subtypes of P2X receptors. A careful probing of the structure activity relationships (SARs) at relevant P2 receptors has resulted in subtype-selective nucleotide agonists. Selective antagonist ligands for P2 receptors have been reported as a result of library screening, conversion of agonists into antagonists, and the careful structural modification of known non-selective ligands. The structures of representative nucleotide (Figure 2A, 29-42) and non-nucleotide (Figure 2B, 43-73) antagonists of the P2 receptors are shown.

## Pharmacological Probes for P2X Receptors

The development of P2X receptor ligands for potential therapeutic application is underway. Selective P2X receptor antagonists are of interest in pain control, urinary incontinence, diabetic retinopathy, inflammatory diseases such as rheumatoid arthritis, and other conditions.

### Non-Selective P2X Ligands

ATP activates all subtypes of P2X receptors, but at different concentrations varying from the low nanomolar to the high micromolar.<sup>18</sup> ADP and AMP, when purified, are inactive at P2X receptors. 2-Methylthioadenosine 5'-triphosphate (2-MeSATP) is a potent agonist at multiple P2X receptors, for example, P2X<sub>1</sub> (EC<sub>50</sub> = 54 nM) and P2X<sub>3</sub> (EC<sub>50</sub> = 350 nM) receptors.  $\alpha,\beta$ -Methyleneadenosine 5'-triphosphate (Figure 1A 6) activates and desensitizes the P2X<sub>1</sub> receptor and is inactive at the P2X<sub>2</sub> receptor. In tritiated form it serves as a radioligand of P2X<sub>1</sub> and P2X<sub>3</sub> receptors.

Older, non-selective and weak P2X antagonists (Figures 2A and 2B), such as organic dyes 43 and 45, have been in use for decades. The antiparasitic drug polysulfonated Suramin and the pyridoxal phosphate derivatives PPADS and positional isomer iso-PPADS are relatively nonsubtype-selective P2X antagonists, that also block some P2Y subtypes.<sup>19</sup> The PPADS analog MRS 2159 is more potent than PPADS at the P2X<sub>1</sub> receptor and also antagonizes the P2X<sub>3</sub> receptor. The nucleotide derivative TNP-ATP

is a potent P2X antagonist that is selective for several subtypes.<sup>20</sup> It antagonizes P2X<sub>1</sub>, P2X<sub>3</sub> and heteromeric P2X<sub>2/3</sub> receptors with IC<sub>50</sub> values of 6, 0.9 and 7 nM respectively, and displays 1000-fold selectivity for P2X<sub>3</sub> over P2X<sub>2</sub>, P2X<sub>4</sub> and P2X<sub>7</sub> receptors. The polysulfonated biphenyl derivative Evans Blue acts as a P2X receptor antagonist, but it also affects other channels and amino acid binding sites.

### P2X<sub>1</sub> and P2X<sub>2</sub> Receptors

P2X<sub>1</sub> antagonists have been reported in several compound classes. For example, the Suramin derivative NF 157 is a P2X<sub>1</sub> antagonist that also blocks the P2Y<sub>11</sub> receptor.<sup>22</sup> Other Suramin derivatives that act as selective P2X<sub>1</sub> antagonists include: PPNSD, NF 279, and the more highly selective P2X<sub>1</sub> antagonist NF 449.<sup>23,24</sup> The earlier-reported Suramin analog NF 023 is a moderately selective, competitive P2X antagonist with IC<sub>50</sub> values of 0.21 and 28.9  $\mu$ M at human P2X<sub>1</sub> and P2X<sub>3</sub> receptors respectively, and is inactive at P2X<sub>2</sub> and P2X<sub>4</sub> receptors.<sup>25</sup> In a separate chemical series, the benzimidazole-2-carboxamide derivative Ro 0437626 was recently reported to be a selective P2X<sub>1</sub> antagonist (IC<sub>50</sub> = 3  $\mu$ M) that displays > 30-fold selectivity over P2X<sub>2</sub>, P2X<sub>3</sub> and P2X<sub>2/3</sub> receptors (IC<sub>50</sub> > 100  $\mu$ M).<sup>26</sup> The dinucleotide Ip<sub>3</sub>I was shown to antagonize the P2Y<sub>1</sub> receptor.<sup>42</sup> The pyridoxal phosphate derivative MRS 2219 is a weak potentiator of P2X<sub>1</sub>-mediated responses.<sup>27</sup>

There are no selective ligands for the P2X<sub>2</sub> receptor. The nonselective antagonists Suramin, TNP-ATP, RB-2, and isoPPADS have been used to study this receptor.

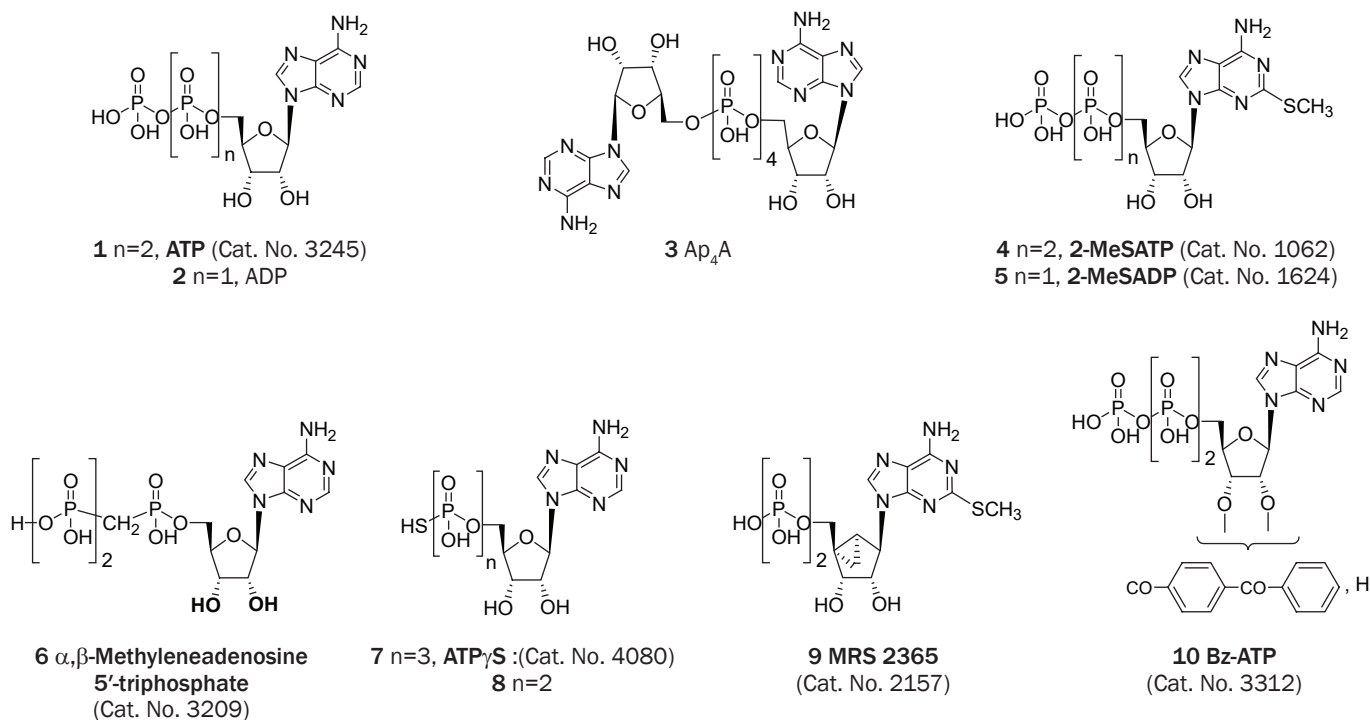
### P2X<sub>3</sub> Receptor

The P2X<sub>3</sub> receptor may exist as a homotrimer or as a heterotrimer in combination with P2X<sub>2</sub> subunits. The Suramin derivative NF 110 is a high affinity P2X<sub>3</sub> receptor antagonist (K<sub>i</sub> values are 36, 82 and 4140 nM for P2X<sub>3</sub>, P2X<sub>1</sub> and P2X<sub>2</sub> receptors respectively) that is inactive at P2Y<sub>1</sub>, P2Y<sub>2</sub> and P2Y<sub>11</sub> receptors (IC<sub>50</sub> > 10  $\mu$ M). A major advance was the introduction of the competitive antagonist by Abbott Laboratories, A 317491, which blocks P2X<sub>3</sub> (IC<sub>50</sub> = 22 nM) and P2X<sub>2/3</sub> (IC<sub>50</sub> = 92 nM) receptors and is roughly three orders of magnitude selective for P2X<sub>3</sub> in comparison to P2X<sub>1</sub> and P2X<sub>2</sub> receptors.<sup>17</sup> A 317491 is inactive at P2X<sub>4</sub> receptors and at all P2Y receptors. Due to the presence of three carboxylic acid groups, A 317491 is of limited bioavailability. Another potent P2X<sub>3</sub> antagonist is the pyrimidinediamine derivative RO-3, which is a selective antagonist of the homomeric P2X<sub>3</sub> and heteromeric P2X<sub>2/3</sub> receptors (pIC<sub>50</sub> values are 7.0 and 5.9 nM respectively) and is inactive at P2X<sub>1</sub>, P2X<sub>2</sub>, P2X<sub>4</sub>, P2X<sub>5</sub> and P2X<sub>7</sub> receptors (IC<sub>50</sub> > 10 M).<sup>29</sup> The endogenous heptapeptide Spinorphin (LVVYPWT) was found to be a very potent P2X<sub>3</sub> antagonist (IC<sub>50</sub> = 8.3  $\mu$ M).<sup>30</sup>

### P2X<sub>4</sub> Receptor

Few of the known P2X antagonists act at the P2X<sub>4</sub> receptor. The benzofurodiazepinone derivative 5-BDBD is an antagonist of P2X<sub>4</sub>-mediated currents (IC<sub>50</sub> = 0.50  $\mu$ M).<sup>31</sup> The bacteria-derived broad spectrum antiparasitic agent Ivermectin, which is a macrocyclic lactone, is a positive allosteric modulator for the P2X<sub>4</sub> receptor, but it also affects other ion channels, such as nicotinic acetylcholine receptors.<sup>32</sup>

**Figure 1A** | Adenine derivatives that have been useful as antagonists in the study of P2 receptors

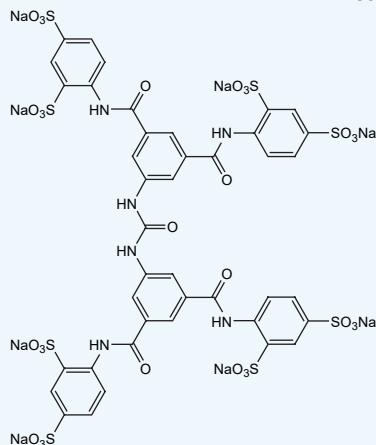


(**Bold text** denotes compounds available from Tocris at time of publication)

### P2X<sub>1</sub> Antagonist

**NF 449**

Cat. No. 1391



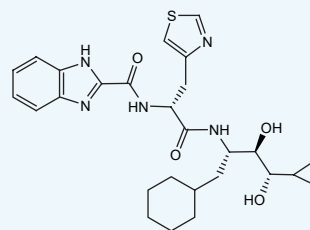
NF 449 is a potent purinergic receptor antagonist that displays high selectivity for P2X<sub>1</sub> (IC<sub>50</sub> values are 0.28, 0.69, 120, 1820, 47,000 and > 300,000 nM for rP2X<sub>1</sub>, rP2X<sub>1+5</sub>, rP2X<sub>2+3</sub>, rP2X<sub>3</sub>, rP2X<sub>2</sub> and P2X<sub>4</sub> receptors respectively). The compound provides antithrombotic protection *in vivo*. NF 449 also acts as a G<sub>sα</sub>-selective antagonist.

Hohenegger *et al.* (1998) Gsa-selective G protein antagonists. *Proc.Natl. Acad. Sci.* **95** 346. Hechler *et al.* (2005) Inhibition of platelet functions and thrombosis through selective or non-selective inhibition of the platelet P2 receptors with increasing doses of NF449 [4,4',4'',4'''-(carbonylbis(imino-5,1,3-benzenetriylbis-(carbonylimino)))tetrakis-benzene-1,3-disulfonic acid octasodium salt]. *J.Pharmacol.Exp.Ther.* **314** 232. Rettinger *et al.* (2005) Profiling at recombinant homomeric and heteromeric rat P2X receptors identifies the suramin analogue NF449 as a highly potent P2X<sub>1</sub> receptor antagonist. *Neuropharmacology.* **48** 461.

### P2X<sub>1</sub> Antagonist

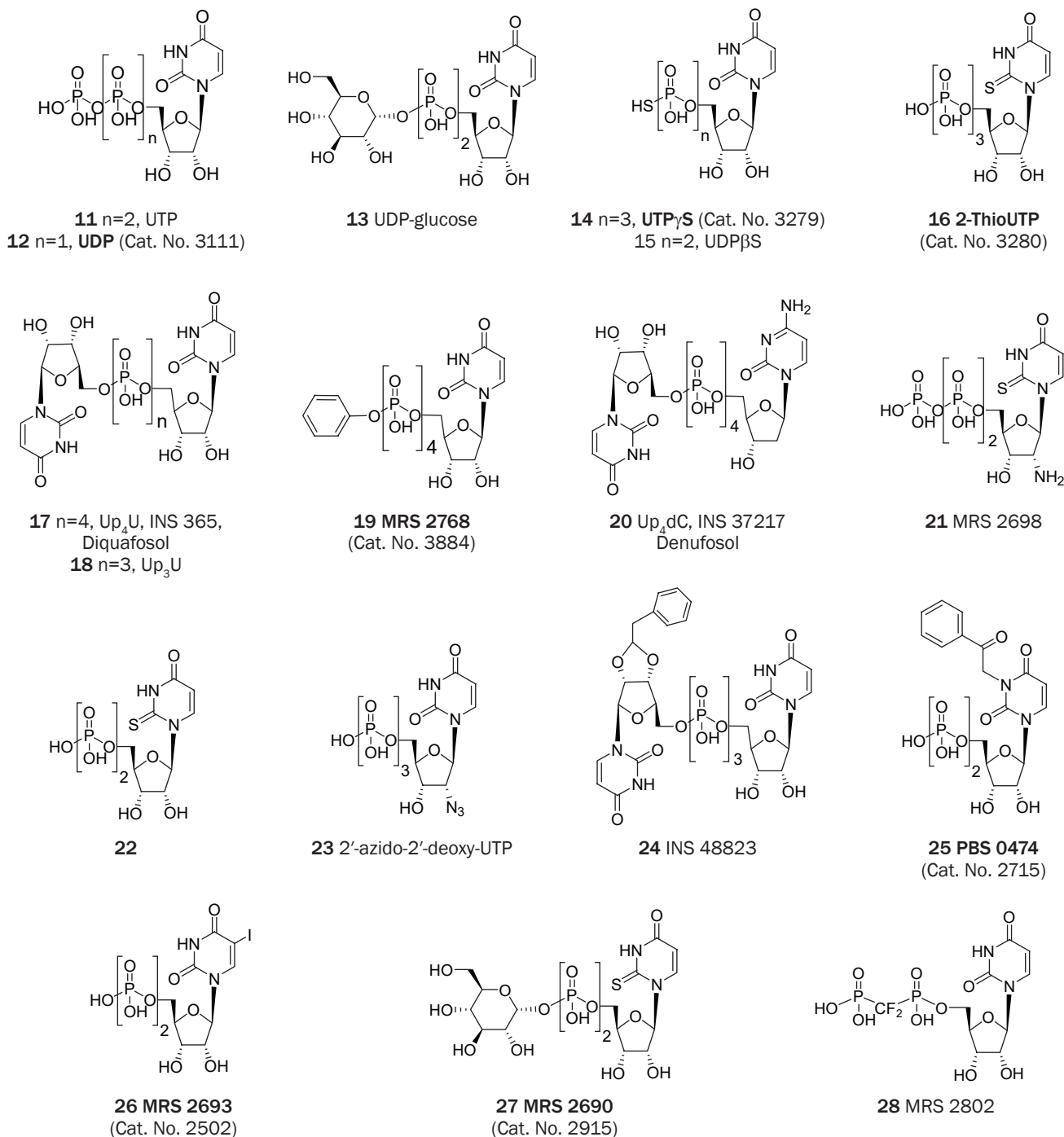
**Ro 0437626**

Cat. No. 2188



Ro 0437626 is a selective P2X<sub>1</sub> purinergic receptor antagonist (IC<sub>50</sub> = 3 μM) that displays > 30-fold selectivity over P2X<sub>2</sub>, P2X<sub>3</sub> and P2X<sub>2/3</sub> receptors (IC<sub>50</sub> > 100 μM).

Jaime-Figueroa *et al.* (2005) Discovery and synthesis of a novel and selective drug-like P2X<sub>1</sub> antagonist. *Bioorg.Med.Chem.Lett.* **15** 3292. King *et al.* (2004) Investigation of the effects of P2 purinoceptor ligands on the micturition reflex in female urethane-anaesthetized rats. *Br.J.Pharmacol.* **142** 519. Ford *et al.* (2006) Purinoceptors as therapeutic targets for lower urinary tract dysfunction. *Br.J.Pharmacol.* **147** S132.

**Figure 1B** | Uracil derivatives that have been useful as antagonists in the study of P2 receptors

(**Bold** text denotes compounds available from Tocris at time of publication)

## P2X<sub>5</sub> Receptor

There are no selective ligands for the P2X<sub>5</sub> and P2X<sub>6</sub> receptors. However, the dye Coomassie Brilliant blue G (BBG) has been used effectively to block P2X<sub>5</sub> receptor function ( $IC_{50} = 0.5 \mu\text{M}$ ),<sup>33</sup> but this dye also blocks P2X<sub>4</sub> receptors ( $IC_{50} = 3 \mu\text{M}$  at human receptors) and P2X<sub>7</sub> receptors ( $IC_{50}$  values are 10 nM and 267 nM at rat and human receptors respectively).

## P2X<sub>7</sub> Receptor

2'-(3')-O-(4-Benzoylbenzoyl)adenosine-5'-triphosphate (BzATP) is a

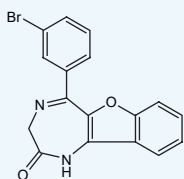
P2X<sub>7</sub> receptor agonist that exhibits an order of magnitude greater potency than ATP. It is also a partial agonist at P2X<sub>1</sub> ( $pEC_{50} = 8.7$ ) and P2Y<sub>1</sub> receptors.<sup>34</sup> One of the first antagonists of the P2X<sub>7</sub> receptor identified was the tyrosine and isoquinoline derivative KN-62, but it acts in a non-competitive fashion and is inactive at the rat P2X<sub>7</sub> homolog. KN-62 is also an inhibitor of CaM kinase II.<sup>35</sup> Oxidized-ATP (o-ATP) has also been used extensively to antagonize P2X<sub>7</sub> receptors.

It has been a challenge to identify antagonists that block the P2X<sub>7</sub> receptor in a species-independent manner. The

## P2X<sub>4</sub> Receptor Antagonist

5-BDBD

Cat. No. 3579



5-BDBD is a potent P2X<sub>4</sub> receptor antagonist. The compound blocks P2X<sub>4</sub>-mediated currents in Chinese hamster ovary cells (IC<sub>50</sub> = 0.50 μM).

Donnelly-Roberts *et al.* (2008) Painful purinergic receptors. *J.Pharmacol.Exp. Ther.* **324** 409.

tetrazolylmethylpyridine derivative A 438079 is a potent, selective, and competitive P2X<sub>7</sub> antagonist (pIC<sub>50</sub> = 6.9).<sup>36</sup> The quinolinamino derivative A 740003 is a potent and selective competitive P2X<sub>7</sub> receptor antagonist (IC<sub>50</sub> values are 40 and 18 nM for human and rat P2X<sub>7</sub> receptors respectively) that is also highly selective in comparison to various P2X and P2Y receptors.<sup>36,37</sup> The adamantyl derivative from AstraZeneca AZ 10606120 antagonizes the P2X<sub>7</sub> receptor with K<sub>D</sub> values of 1.4 and 19 nM at human and rat P2X<sub>7</sub> receptors respectively.<sup>38</sup> The biphenyl derivative AZ 11645373 potently antagonized the human P2X<sub>7</sub> receptor in a non-surmountable manner with K<sub>B</sub> values ranging from 5–20 nM and was inactive at the rat P2X<sub>7</sub> receptor and at all other P2X subtypes.<sup>39</sup> The substituted glycylic anilide derivative GW 791343 is a positive allosteric modulator of the rat P2X<sub>7</sub> receptor and a negative allosteric modulator of the human P2X<sub>7</sub> receptor (pIC<sub>50</sub> = 6.9–7.2).<sup>40</sup>

## Pharmacological Probes for P2Y Receptors

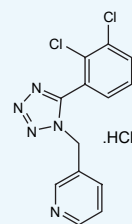
There has been progress in the development of selective agonist and antagonist ligands for P2Y receptors for preclinical development.<sup>2</sup> Until recently, the only P2Y receptor ligand in pharmaceutical use was the antithrombotic P2Y<sub>12</sub> receptor antagonist Clopidogrel (Plavix).<sup>41</sup> Therefore, there is much activity to identify new agents to act at the P2Y<sub>12</sub> receptor and at other P2Y receptors for pharmaceutical development. The rapidly accelerating progress in this field has already resulted in new drug candidates for pulmonary diseases, dry eye disease, and thrombosis.

Many selective ligand probes, both agonists and antagonists of the P2Y receptors, are now available. However, some subtypes, such as the P2Y<sub>4</sub> receptor, are entirely lacking such selective ligands. Detailed SAR analyses have been constructed for P2Y<sub>1</sub> and P2Y<sub>12</sub> receptors, which are both proaggregatory in platelets. Nucleotide agonists selective for P2Y<sub>1</sub>, P2Y<sub>2</sub>, P2Y<sub>6</sub>, and P2Y<sub>14</sub> receptors and nucleotide antagonists selective for P2Y<sub>1</sub> and P2Y<sub>12</sub> receptors have been described. Selective non-nucleotide antagonists are now sought to avoid issues of limited stability and bioavailability. Such antagonists have been reported for P2Y<sub>1</sub>, P2Y<sub>2</sub>, P2Y<sub>6</sub>, P2Y<sub>11</sub>, P2Y<sub>12</sub>, and P2Y<sub>13</sub> receptors. The screening of chemically diverse compound libraries has resulted in competitive P2Y<sub>12</sub> receptor antagonists that are being tested as potential antithrombotic agents.

## P2X<sub>7</sub> Receptor Antagonist

A 438079

Cat. No. 2972



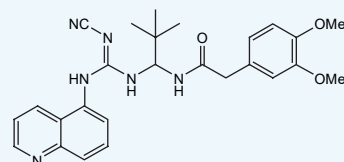
A 438079 is a competitive P2X<sub>7</sub> receptor antagonist (pIC<sub>50</sub> = 6.9 for the inhibition of Ca<sup>2+</sup> influx in a human recombinant cell line expressing P2X<sub>7</sub>). The compound is devoid of activity at other P2 receptors (IC<sub>50</sub> >> 10 μM). A 438079 possesses antinociceptive activity in models of neuropathic pain *in vivo*.

Nelson *et al.* (2006) Structure-activity relationship studies on a series of novel, substituted 1-benzyl-5-phenyltetrazole P2X<sub>7</sub> antagonists. *J.Med.Chem.* **49** 3659. Donnelly-Roberts and Jarvis (2007) Discovery of P2X<sub>7</sub> receptor-selective antagonists offers new insights into P2X<sub>7</sub> receptor function and indicates a role in chronic pain states. *Br.J.Pharmacol.* **151** 571. McGaraughty *et al.* (2007) P2X<sub>7</sub>-related modulation of pathological nociception in rats. *Neuroscience* **146** 1817.

## P2X<sub>7</sub> Receptor Antagonist

A 740003

Cat. No. 3701



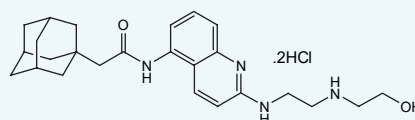
A 740003 is a potent, selective and competitive P2X<sub>7</sub> receptor antagonist (IC<sub>50</sub> values are 18 and 40 nM for rat and human receptors respectively). The compound displays selectivity over a variety of P2X and P2Y receptors up to a concentration of 100 μM. A 740003 reduces nociception in animal models of persistent neuropathic and inflammatory pain.

Honore *et al.* (2006) A-740003 [N-(1-((cyanoimino)(5-quinolinylamino)methyl)amino)-2,2-dimethylpropyl)-2-(3,4-dimethoxyphenyl)acetamide], a novel and selective P2X<sub>7</sub> receptor antagonist, dose-dependently reduces neuropathic pain in the rat. *J.Pharmacol.Exp.Ther.* **319** 1376. King (2007) Novel P2X<sub>7</sub> receptor antagonists ease the pain. *Br.J.Pharmacol.* **151** 565. Donnelly-Roberts *et al.* (2009) Mammalian P2X<sub>7</sub> receptor pharmacology: comparison of recombinant mouse, rat and human P2X<sub>7</sub> receptors. *Br.J.Pharmacol.* **157** 1203.

## P2X<sub>7</sub> Receptor Antagonist

AZ 10606120

Cat. No. 3323



AZ 10606120 is a potent P2X<sub>7</sub> receptor antagonist (K<sub>D</sub> values are 1.4 and 19 nM at human and rat P2X<sub>7</sub> receptors respectively). The compound binds in a positive cooperative manner to sites distinct from, but coupled to, the ATP binding site and acts as a negative allosteric modulator.

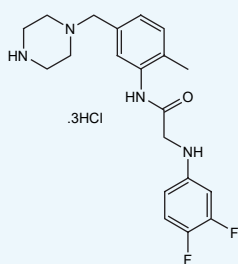
Michel *et al.* (2007) Direct labelling of the human P2X<sub>7</sub> receptor and identification of positive and negative cooperativity of binding. *Br.J.Pharmacol.* **151** 103. Michel and Fonfria (2007) Agonist potency at P2X<sub>7</sub> receptors is modulated by structurally diverse lipids. *Br.J.Pharmacol.* **152** 523. Michel *et al.* (2008) Negative and positive allosteric modulators of the P2X<sub>7</sub> receptor. *Br.J.Pharmacol.* **153** 737.

(Sold for research purposes under agreement from AstraZeneca)

P2X<sub>7</sub> Allosteric Modulator

GW 791343

Cat. No. 3385



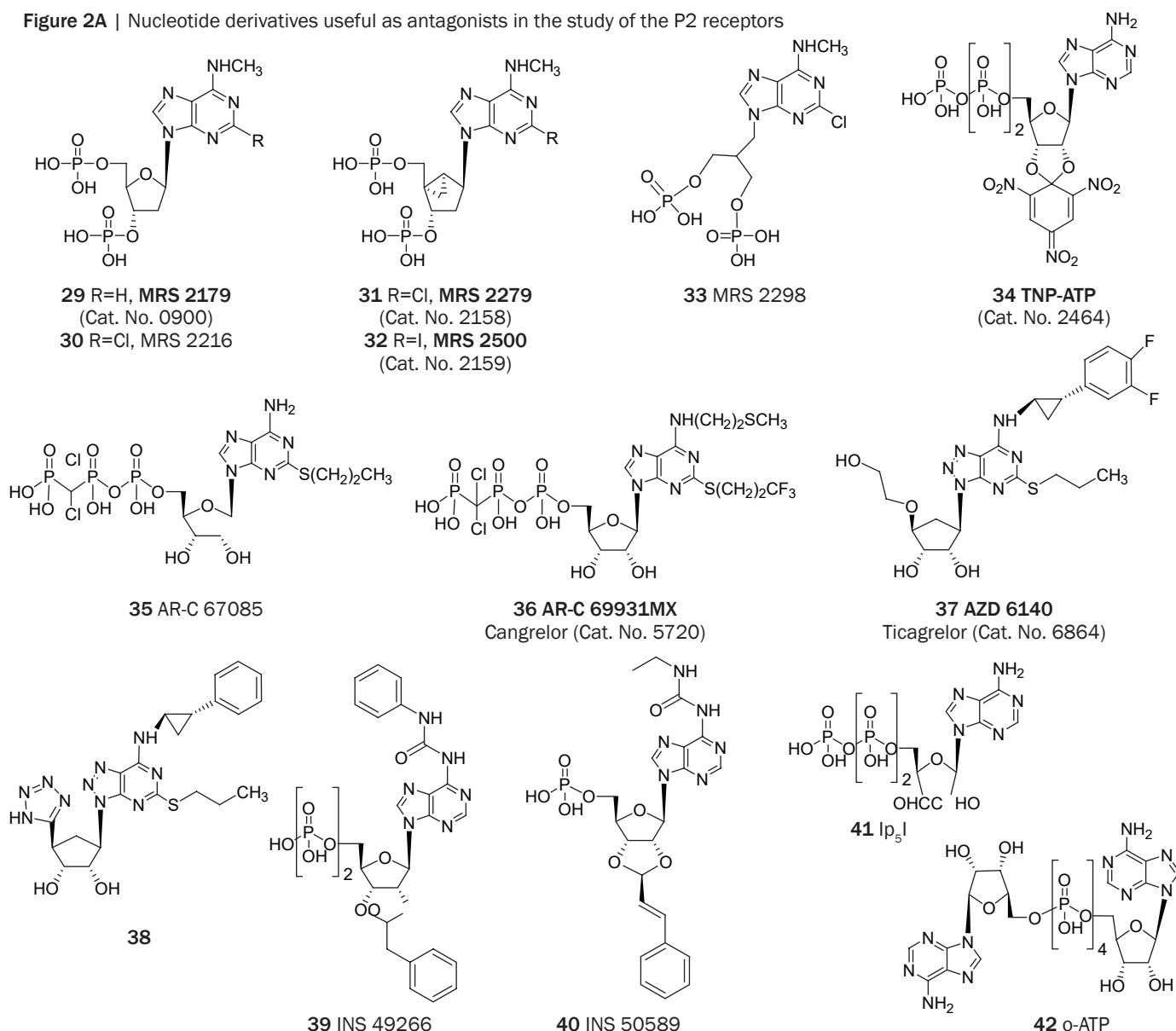
GW 791343 is a P2X<sub>7</sub> allosteric modulator. The compound exhibits species-specific activity and acts as a negative allosteric modulator of human P2X<sub>7</sub> (pIC<sub>50</sub> = 6.9–7.2) and a positive allosteric modulator of rat P2X<sub>7</sub>.

Michel *et al.* (2008) Negative and positive allosteric modulators of the P2X<sub>7</sub> receptor. *Br.J.Pharmacol.* **153** 737. Michel *et al.* (2008) Identification of regions of the P2X<sub>7</sub> receptor that contribute to human and rat species differences in antagonist effects. *Br.J.Pharmacol.* **155** 738.

P2Y<sub>1</sub> Receptor

2-MeSADP, like ADP, activates the P2Y<sub>1</sub>, P2Y<sub>12</sub> and P2Y<sub>13</sub> receptors. 2-MeSADP is a more potent agonist at the P2Y<sub>1</sub> receptor than 2-MeSATP. N<sup>6</sup>-methyl nucleotides are tolerated at the P2Y<sub>1</sub> receptor, consistent with a small hydrophobic pocket in the P2Y<sub>1</sub> receptor binding site surrounding the N<sup>6</sup>-position of adenine nucleotides. The favored ribose-ring conformation for each of the subtypes of the P2Y<sub>1</sub>-like subfamily has been established using conformationally-restricted (i.e. rigid) ribose equivalents, which also improve stability of the phosphate esters toward nucleotidases. In particular, the methanocarba ring system consisting of fused cyclopropane and cyclopentane has been useful in exploring the biologically active conformations of nucleoside and nucleotide derivatives. The North (N)-methanocarba analog of 2-MeSADP, MRS 2365 (EC<sub>50</sub> = 0.4 nM), is a selective, high affinity agonist of the P2Y<sub>1</sub> receptor.<sup>11</sup> Another means of improving hydrolytic stability is the introduction of a borano group within the phosphate moiety of P2Y receptor agonists.<sup>43</sup> Many nucleotide antagonists of the P2Y<sub>1</sub> receptor

**Figure 2A** | Nucleotide derivatives useful as antagonists in the study of the P2 receptors



(Bold text denotes compounds available from Tocris at time of publication)

have been introduced. Usually these are adenine nucleotides containing bisphosphate groups, for example a ribose 3',5'-bisphosphate moiety. A *N*<sup>6</sup>-methyl 2'-deoxyadenosine bisphosphate derivative MRS 2179 ( $pK_B = 6.99$ ) and its 2-chloro analog MRS 2216 are selective P2Y<sub>1</sub> antagonists.<sup>44</sup> The same (N)-conformational constraint of the ribose moiety that enhances agonist action also favors the potency and selectivity in nucleotide antagonists. For example, the ring-constrained (N)-methanocarba nucleotide bisphosphates MRS 2279 ( $pK_B = 8.10$ ) and MRS 2500 ( $pK_B = 9$ ) are selective, high affinity antagonists of the P2Y<sub>1</sub> receptor.<sup>45</sup> The antithrombotic action of MRS 2500 (by blocking the P2Y<sub>1</sub> receptor selectively) has been demonstrated *in vivo* in the mouse and other species.<sup>46,47</sup> Antagonists of the P2Y<sub>1</sub> receptor of moderate affinity may also be derived from acyclic nucleotides, such as the bisphosphates derivative MRS 2298.<sup>46</sup>

Non-nucleotide antagonists of the P2Y<sub>1</sub> receptor have been discovered through screening of structurally diverse chemical libraries. The first such compound to be reported was **63**, which is a selective and orally bioavailable antagonist of the human P2Y<sub>1</sub> receptor from Pfizer of novel chemotype with a  $K_i$  value of 90 nM.<sup>48</sup> Other structurally diverse antagonists of the P2Y<sub>1</sub> receptor have been reported. Pyridyl isatogen (PIT) is an allosteric modulator of the P2Y<sub>1</sub> receptor that displays mixed antagonism/potentialiation.<sup>49</sup>

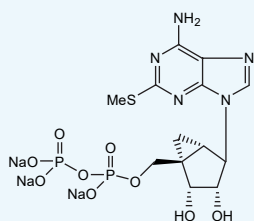
## P2Y<sub>2</sub> and P2Y<sub>4</sub> Receptors

Both P2Y<sub>2</sub> and P2Y<sub>4</sub> receptors are activated by UTP, and simple modifications enhance selectivity for the P2Y<sub>2</sub> receptor. UTP<sub>γ</sub>S is a more selective and stable agonist of the P2Y<sub>2</sub> receptor than UTP.<sup>50</sup> However, this compound is subject to chemical oxidation. 2-ThioUTP is also a selective agonist of the P2Y<sub>2</sub> receptor.<sup>51</sup> Combination of modifications of UTP in the selective P2Y<sub>2</sub> agonist MRS 2698 provided an  $EC_{50}$  of 8 nM and selectivity of 300-fold in comparison to the P2Y<sub>4</sub> receptor.<sup>52</sup> Certain dinucleoside tetraphosphates potently activate the P2Y<sub>2</sub> and P2Y<sub>4</sub> receptors. The uracil dinucleotides that have been in clinical trials are Up<sub>4</sub>U (INS 365, Diquafosol,  $EC_{50} = 0.1 \mu\text{M}$ ) and Up<sub>4</sub>dC (INS 37217, Denufosol,  $EC_{50} = 0.22 \mu\text{M}$ ).<sup>53</sup> Diquafosol was recently approved in Japan for use in treating dry eye. By virtue of being dinucleotides,

## P2Y<sub>1</sub> Agonist

MRS 2365

Cat. No. 2157



MRS 2365 is a highly potent, selective P2Y<sub>1</sub> receptor agonist ( $EC_{50} = 0.4 \text{ nM}$ ). The compound displays no activity at P2Y<sub>12</sub> receptors and only very low agonist activity at P2Y<sub>13</sub> receptors (at concentrations up to 1  $\mu\text{M}$ ).

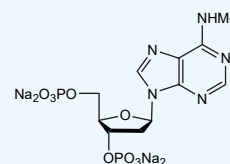
Ravi *et al.* (2002) Adenine nucleotide analogues locked in a northern methanocarba conformation: enhanced stability and potency as P2Y<sub>1</sub> receptor agonists. *J.Med.Chem.* **45** 2090. Chhatrivala *et al.* (2004) Induction of novel agonist selectivity for the ADP-activated P2Y<sub>1</sub> receptor versus the ADP-activated P2Y<sub>12</sub> and P2Y<sub>13</sub> receptors by conformational constraint of an ADP analog. *J.Pharmacol.Exp.Ther.* **311** 1038.

(Sold under license from the NIH. US Patent 10/169975)

## P2Y<sub>1</sub> Antagonist

MRS 2179

Cat. No. 0900



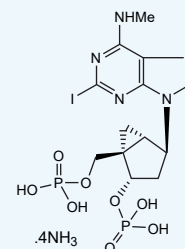
MRS 2179 is a competitive antagonist at P2Y<sub>1</sub> receptors ( $K_B = 100 \text{ nM}$ ). The compound is selective over P2X<sub>1</sub> ( $IC_{50} = 1.15 \mu\text{M}$ ), P2X<sub>3</sub> ( $IC_{50} = 12.9 \mu\text{M}$ ), P2X<sub>2</sub>, P2X<sub>4</sub>, P2Y<sub>2</sub>, P2Y<sub>4</sub> and P2Y<sub>6</sub> receptors.

Boyer *et al.* (1998) Competitive and selective antagonism of P2Y<sub>1</sub> receptors by *N*<sup>6</sup>-methyl 2'-deoxyadenosine 3',5'-bisphosphate. *Br.J.Pharmacol.* **124** 1. Moro *et al.* (1998) Human P2Y<sub>1</sub> receptor molecular modeling and site-directed mutagenesis as tools to identify agonist and antagonist recognition sites. *J.Med. Chem.* **41** 1456. Nandan *et al.* (2000) Synthesis, biological activity, and molecular modeling of ribose-modified deoxyadenosine bisphosphate analogues as P2Y<sub>1</sub> receptor ligands. *J.Med.Chem.* **43** 829. Brown *et al.* (2000) Activity of novel adenine nucleotide derivatives as agonists and antagonists at recombinant rat P2X receptors. *Drug Dev.Res.* **49** 253.

## P2Y<sub>1</sub> Antagonist

MRS 2500

Cat. No. 2159



MRS 2500 is a highly potent and selective antagonist of the platelet P2Y<sub>1</sub> receptor ( $K_i = 0.78 \text{ nM}$ ). The compound inhibits ADP-induced aggregation of human platelets with an  $IC_{50}$  value of 0.95 nM. MRS 2500 prevents thrombus formation *in vivo*.

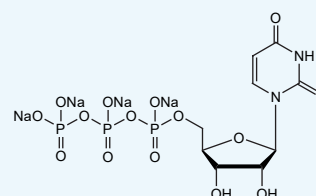
Kim *et al.* (2003) 2-Substitution of adenine nucleotide analogues containing a bicyclo[3.1.0]hexane ring system locked in a northern conformation: enhanced potency as P2Y<sub>1</sub> receptor antagonists. *J.Med.Chem.* **46** 4974. Cattaneo *et al.* (2004) Antiaggregatory activity in human platelets of potent antagonists of the P2Y<sub>1</sub> receptor. *Biochem.Pharmacol.* **68** 1995. Hechler *et al.* (2006) MRS2500 [2-iodo-*N*<sup>6</sup>-methyl-(*N*)-methanocarba-2'-deoxyadenosine-3',5'-bisphosphate], a potent, selective, and stable antagonist of the platelet P2Y<sub>1</sub> receptor with strong antithrombotic activity in mice. *J.Pharmacol.Exp.Ther.* **316** 556.

(Sold under license from the NIH, US Patent 60/029,855.)

## P2Y<sub>2</sub> Agonist

2-ThioUTP tetrasodium salt

Cat. No. 3280



2-ThioUTP is a potent and selective P2Y<sub>2</sub> agonist.  $EC_{50}$  values are 0.035, 0.35 and 1.5  $\mu\text{M}$  for hP2Y<sub>2</sub>, hP2Y<sub>4</sub> and hP2Y<sub>6</sub> receptors respectively.

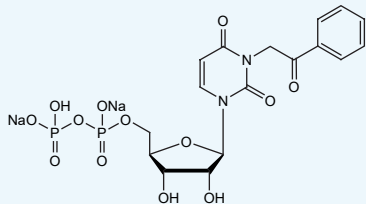
El-Tayeb *et al.* (2006) Synthesis and structure-activity relationships of uracil nucleotide derivatives and analogues as agonists at human P2Y<sub>2</sub>, P2Y<sub>4</sub> and P2Y<sub>6</sub> receptors. *J.Med.Chem.* **49** 7076. Ko *et al.* (2008) Synthesis and potency of novel uracil nucleotides and derivatives as P2Y<sub>2</sub> and P2Y<sub>6</sub> receptor agonists. *Bioorg.Med.Chem.* **16** 6319.



## P2Y<sub>6</sub> Agonist

PSB 0474

Cat. No. 2715



PSB 0474 is a potent and selective P2Y<sub>6</sub> receptor agonist. EC<sub>50</sub> values are 70, > 1000 and > 10,000 nM for P2Y<sub>6</sub>, P2Y<sub>2</sub> and P2Y<sub>4</sub> receptors respectively.

El-Tayeb et al. (2006) Synthesis and structure-activity relationships of uracil nucleotide derivatives and analogues as agonists at human P2Y<sub>2</sub>, P2Y<sub>4</sub> and P2Y<sub>6</sub> receptors. *J.Med.Chem.* **49** 7076.

they are more stable to enzymatic hydrolysis than nucleoside triphosphates, but these agonists are non-selective compared to the P2Y<sub>4</sub> receptor. The 2'-deoxycytidine (dC) moiety of 20 serves to enhance the *in vivo* stability toward ectonucleotidases. The agonist MRS 2768 (uridine tetraphosphate δ-phenyl ester) is selective for the P2Y<sub>2</sub> receptor with moderate potency (EC<sub>50</sub> = 1.89 μM).<sup>54</sup>

Definitive antagonists of the P2Y<sub>2</sub> receptor are not available. AR-C 126313 and its higher molecular weight analog AR-C 118925 were reported to selectively antagonize the P2Y<sub>2</sub> receptor, however it appears that these compounds are only micromolar in affinity (Figure 2B).<sup>55</sup> The large polyanionic molecules Reactive blue 2 (RB2, an anthraquinone dye) and Suramin are slightly selective antagonists of the P2Y<sub>2</sub> and P2Y<sub>4</sub> receptors, respectively. However, RB2 and Suramin also block various P2X receptors (Table 1).

There are no truly selective ligands for the P2Y<sub>4</sub> receptor. The agonist 2'-azido-2'-deoxy-UTP (Figure 1B 23) displayed 5-fold P2Y<sub>4</sub> selectivity.<sup>51</sup> Thus, new agonists and antagonists are needed to distinguish this subtype pharmacologically from the P2Y<sub>2</sub> receptor, which is also activated by UTP. The other native agonist of the P2Y<sub>2</sub> receptor, ATP, acts as an antagonist at the human, but not the rat, P2Y<sub>4</sub> receptor.

## P2Y<sub>6</sub> Receptors

UDP activates both the P2Y<sub>6</sub> and P2Y<sub>14</sub> receptors. It is worth noting, however, that extracellular UDP can serve as a substrate for the generation of UTP through the action of nucleoside diphosphokinase (NDPK), which may complicate pharmacological studies. The action of UDP at the P2Y<sub>14</sub> receptor has been controversial. UDP was initially described as inactive at the newly cloned P2Y<sub>14</sub> receptor, however later study found an antagonist action of UDP at the human but not rat P2Y<sub>14</sub> receptor. Finally the observed antagonist action of UDP was shown to occur in cells expressing an unnatural, engineered chimeric G protein. However, in HEK293 and other cells in which endogenous G<sub>i</sub> proteins mediate the functional response, UDP acts as a potent agonist.<sup>5</sup> Thus, inaccurate results might be obtained using UDP alone in pharmacological studies if multiple P2Y subtypes are present. UDPβS (15) and Up<sub>3</sub>U have been used as more stable activators of the P2Y<sub>6</sub> receptor subtype than UDP,<sup>50,53</sup> although UDPβS also activates the P2Y<sub>14</sub> receptor.

The SAR of nucleotide derivatives in activating the P2Y<sub>6</sub> receptor has been explored. Certain dinucleoside triphosphates have been explored as P2Y<sub>6</sub> receptor ligands, for example, INS 48823 (EC<sub>50</sub> = 125 nM) potently activates the receptor.<sup>56</sup> Other UDP derivatives, e.g. 3-phenacyl UDP (PSB 0474) and 5-iodo-UDP (MRS 2693), are selective P2Y<sub>6</sub> agonists with EC<sub>50</sub> values of 70 and 15 nM respectively.<sup>57,58</sup> Molecular modeling predicted that the South (S)-conformation of the ribose ring is the preferred conformation in receptor binding, which was then confirmed by synthesis of a conformationally constrained methanocarbonyl analog of UDP. The noncompetitive P2Y<sub>6</sub> receptor antagonist MRS 2578 is a diisothiocyanate derivative, which has low stability and solubility in aqueous medium.<sup>59</sup> Competitive antagonists of the P2Y<sub>6</sub> receptor have not yet been reported.

## P2Y<sub>11</sub> Receptors

ATPγS (Figure 1A) acts as a potent P2Y<sub>11</sub> receptor agonist. The P2Y<sub>12</sub> antagonist 2-propylthio-β,γ-dichloromethylene-ATP (AR-C 67085, Figure 2A) is also the most potent reported agonist of the P2Y<sub>11</sub> receptor (EC<sub>50</sub> = 8.9 μM).<sup>60</sup> Thus, it must be used with caution in pharmacological studies in which both P2Y subtypes might be present.

A potent antagonist NF 157, derived from nonselective P2 antagonist Suramin, has been reported to be a selective antagonist of the P2Y<sub>11</sub> receptor (pK<sub>i</sub> = 7.35).<sup>61</sup> However, this compound also antagonizes the P2X<sub>1</sub>, P2X<sub>2</sub>, and P2X<sub>3</sub> receptors.

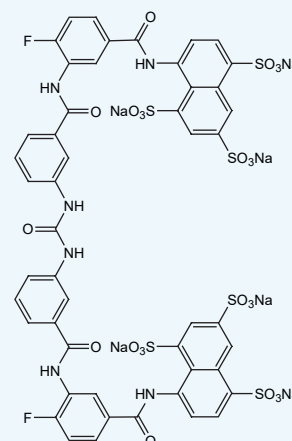
## P2Y<sub>12</sub> Receptors

The medicinal chemistry of the P2Y<sub>12</sub> receptor has been extensively explored. The thienopyridines, such as Clopidogrel (Figure 2B), act as liver-activated prodrugs that are irreversible inhibitors of the P2Y<sub>12</sub> receptor.<sup>62</sup> This thienopyridine P2Y<sub>12</sub>

## P2X<sub>1</sub>/P2YX<sub>11</sub> Antagonist

NF 157

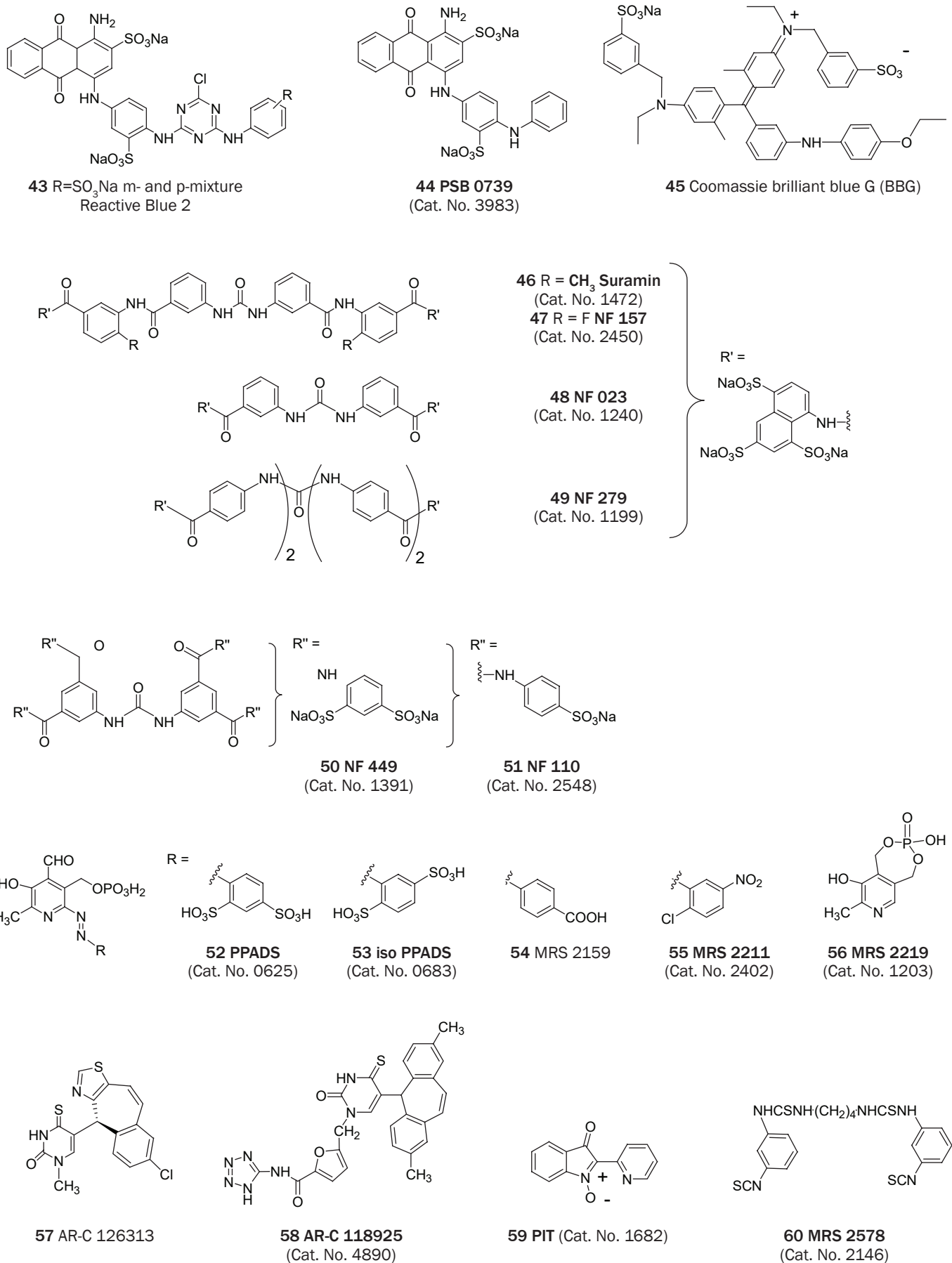
Cat. No. 2450



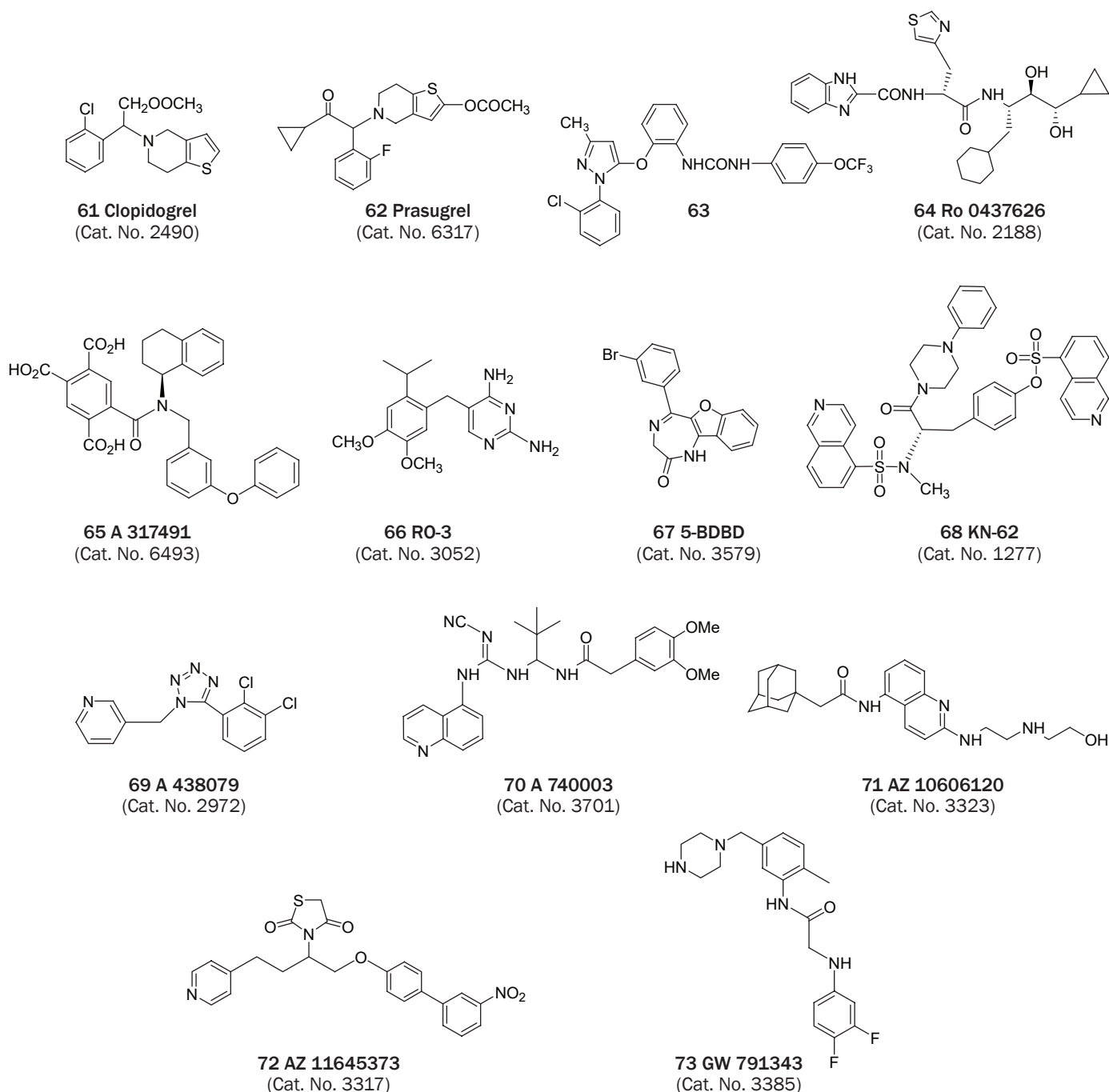
NF 157 is a purinergic receptor antagonist that potently inhibits P2Y<sub>11</sub> receptor activity (IC<sub>50</sub> = 463 nM). The compound displays selectivity for P2Y<sub>11</sub> and P2X<sub>1</sub> receptors over P2Y<sub>1</sub>, P2Y<sub>2</sub>, P2X<sub>2</sub>, P2X<sub>3</sub>, P2X<sub>4</sub> and P2X<sub>7</sub> receptors. NF 157 inhibits NAD<sup>+</sup>-induced activation of human granulocytes.

Ullmann et al. (2005) Synthesis and structure-activity relationships of suramin-derived P2Y<sub>11</sub> receptor antagonists with nanomolar potency. *J.Med.Chem.* **48** 7040. Moreschi et al. (2006) Extracellular NAD<sup>+</sup> is an agonist of the human P2Y<sub>11</sub> purinergic receptor in human granulocytes. *J.Biol.Chem.* **281** 31419.

**Figure 2B** | Non-nucleotides that have been useful antagonists in the study of P2 receptors



(Bold text denotes compounds available from Tocris at time of publication)

**Figure 2B** | Non-nucleotides that have been useful antagonists in the study of P2 receptors

(**Bold** text denotes compounds available from Tocris at time of publication)

receptor antagonist requires a two-step preactivation *in vivo* and therefore has a delayed onset of action and long reversal of the platelet effect after drug administration is stopped. Another thienopyridine antagonist that has been in clinical trials, Prasugrel is a more potent P2Y<sub>12</sub> antagonist, but displays a longer bleeding time. Prasugrel only requires one step of preactivation *in vivo*.<sup>63</sup>

Directly-acting P2Y<sub>12</sub> receptor antagonists have also been reported. The observation that ATP acts as an antagonist at this ADP-activated subtype has enabled the introduction of various 5'-triphosphate analogs as selective receptor probes and clinical

candidates. Thus, the antithrombotic nucleotide derivatives from AstraZeneca AR-C 67085 ( $EC_{50} = 30 \mu\text{M}$ ) and ARC 69931MX (Cangrelor,  $EC_{50} = 0.4 \text{ nM}$ ) have been tested clinically as antithrombotic agents.<sup>64</sup> A 5'-triphosphate group in adenine nucleotides is not strictly required for P2Y<sub>12</sub> receptor antagonism, as in the case of compound **38** and the potent antagonist and approved medicine ticagrelor (AZD 6140  $pIC_{50} = 7.9$ ).<sup>64,65</sup> Other nucleotide antagonists of the P2Y<sub>12</sub> receptor that have been reported are nucleotide derivatives from Inspire Pharmaceuticals, INS 49266 (an ADP derivative with  $EC_{50}$  of 52 nM) and INS 50589 (an AMP derivative with  $EC_{50}$  of 11 nM).<sup>66</sup> Library screening has aided greatly in the identification of novel chemotypes that act as

P2Y<sub>12</sub> receptor antagonists, and several of these compounds are being developed by the pharmaceutical industry. One very potent and selective competitive antagonist of the P2Y<sub>12</sub> receptor, PSB 0739, derived from RB2 was recently reported.

### P2Y<sub>13</sub> Receptors

ADP is also the preferred agonist ligand at the P2Y<sub>13</sub> receptor, and ATP is less potent. In the rat, ADP is 3-5-fold more potent than 2-MeSADP. A selective P2Y<sub>13</sub> receptor antagonist, MRS 2211, a derivative of PPADS, has a pK<sub>i</sub> of 6.0 at this receptor.<sup>67</sup>

### P2Y<sub>14</sub> Receptors

The SAR of analogs of UDP-glucose (EC<sub>50</sub> = 0.35 μM) and UDP at the P2Y<sub>14</sub> receptor was recently systematically explored.<sup>68</sup> Other naturally occurring UDP-sugars activate this receptor less potently. The 2-thio analog of UDP-glucose, MRS 2690, is a 6-fold more potent agonist for the P2Y<sub>14</sub> receptor and, unlike UDP-glucose, is inactive at the P2Y<sub>2</sub> receptor. The P2Y<sub>14</sub> receptor is structurally restrictive with respect to modification of the nucleobase, ribose, and phosphate moieties of agonist ligands. However, the glucose moiety may be deleted in UDP analogs, some of which still are very potent in receptor activation. For example, α,β-difluoromethylene-UDP, MRS 2802, is inactive at the P2Y<sub>6</sub> receptor and fully activates the human P2Y<sub>14</sub> receptor with an EC<sub>50</sub> of 63 nM.

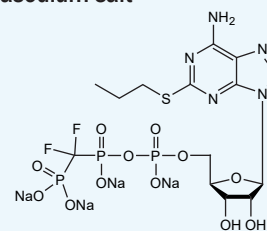
### Conclusion

Novel ligands for the P2X and P2Y receptor families are now available for use as tools in pharmacological studies. Selective nucleotide agonist ligands, although typically of low bioavailability and stability *in vivo*, have been designed. Recently, selective antagonist ligands for P2 receptors have been reported as a result of library screening, conversion of agonists into antagonist, and the careful structural modification of known non-selective ligands.

### P2Y<sub>12</sub> Antagonist

AR-C 66096 tetrasodium salt

Cat. No. 3321



AR-C 66096 is a potent and selective P2Y<sub>12</sub> receptor antagonist. The compound blocks ADP-induced inhibition of adenylyl cyclase *in vitro* (pK<sub>B</sub> = 7.6) and inhibits ADP-induced aggregation of washed human platelets (pIC<sub>50</sub> = 8.16).

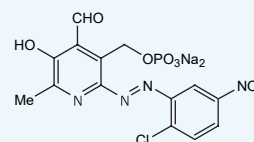
Humphries *et al.* (1994) FPL 66094: a novel, highly potent and selective antagonist at human platelet P<sub>2U</sub>-purinoceptors. *Br.J.Pharmacol.* **113** 1057. Ingall *et al.* (1999) Antagonists of the platelet P receptor: a novel approach to antithrombotic therapy. *J.Med.Chem.* **42** 213. Simon *et al.* (2001) Activity of adenosine diphosphates and triphosphates on a P2Y<sub>7</sub>-type receptor in brain capillary endothelial cells. *Br.J.Pharmacol.* **132** 173.

(Sold for research purposes under agreement from AstraZeneca)

### P2Y<sub>13</sub> Antagonist

MRS 2211

Cat. No. 2402



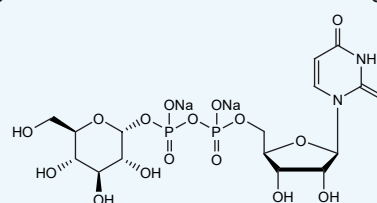
MRS 2211 is a competitive P2Y<sub>13</sub> receptor antagonist (pIC<sub>50</sub> = 5.97). The compound displays > 20-fold selectivity over P2Y<sub>1</sub> and P2Y<sub>12</sub> receptors.

Kim *et al.* (2005) Synthesis of pyridoxal phosphate derivatives with antagonist activity at the P2Y<sub>13</sub> receptor. *Biochem.Pharmacol.* **70** 266. von Kugelgen (2006) Pharmacological profiles of cloned mammalian P2Y-receptor subtypes. *Pharmacol.Ther.* **110** 415.

### P2Y<sub>14</sub> Agonist

MRS 2690

Cat. No. 2915



MRS 2690 is a potent P2Y<sub>14</sub> receptor agonist (EC<sub>50</sub> = 49 nM). The compound displays 7-fold higher potency than UDP-glucose.

Ko *et al.* (2007) Structure-activity relationship of uridine 5'-diphosphoglucose analogues as agonists of the human P2Y<sub>14</sub> receptor. *J.Med.Chem.* **50** 2030.

## References

1. Surprenant and North (2009) *Ann.Rev.Physiol.* **71** 333.
2. Abbracchio et al. (2006) *Pharmacol. Rev.* **58** 281.
3. Kawate et al. (2009) *Nature* **460** 592.
4. Young (2009) *Trends Biochem.Sci.* **35** 83.
5. Carter et al. (2009) *Mol.Pharmacol.* **76** 1341.
6. Ivanov et al. (2006) *J.Comput.Aided Mol.Des.* **20** 417.
7. Hoffmann et al. (1999) *J.Biol.Chem.* **274** 14639.
8. Hillmann et al. (2009) *J.Med.Chem.* **52** 2762.
9. Zylberg et al. (2007) *Biochem.J.* **405** 277.
10. Hoffmann et al. (2009) *J.Pharmacol.Exp.Ther.* **331** 648.
11. Bourdon et al. (2006) *J.Thromb.Haemost.* **4** 861.
12. Ecke et al. (2008) *Biochem.J.* **409** 107.
13. Nakata et al. (2003) *Drug Dev. Res.* **58** 340.
14. Baqi et al. (2010) *J.Med.Chem.* **53** 2076.
15. Houston et al. (2006) *Br.J.Pharmacol.* **147** 459.
16. El-Tayeb et al. (2005) *Bioorg.Med.Chem.Lett.* **15** 5450.
17. Jarvis et al. (2004) *J.Pharmacol.Exp.Ther.* **310** 407.
18. Jacobson et al. (2002) *J.Med.Chem.* **45** 4057.
19. Kim et al. (2001) *J.Med.Chem.* **44** 340.
20. Virginio et al. (1998) *Mol.Pharmacol.* **53** 969.
21. Whittenburg et al. (1996) *Naunyn-Schmiedeberg's.Arch.Pharmacol.* **354** 491.
22. Ullmann et al. (2005) *J.Med.Chem.* **48** 7040.
23. Rettinger et al. (2000) *Neuropharmacology* **39** 2044.
24. Rettinger et al. (2005) *Neuropharmacology* **48** 461.
25. Soto et al. (1999) *Neuropharmacology* **38** 141.
26. Jaime-Figueroa et al. (2005) *Bioorg.Med.Chem.Lett.* **15** 3292.
27. Jacobson et al. (1998) *J.Med.Chem.* **41** 2201.
28. Hausmann et al. (2006) *Mol.Pharmacol.* **69** 2058.
29. Carter et al. (2009) *Bioorg.Med.Chem.Lett.* **19** 1628.
30. Jung et al. (2007) *J.Med.Chem.* **50** 4543.
31. Donnelly-Roberts et al. (2008) *J.Pharmacol.Exp.Ther.* **324** 409.
32. Silberberg et al. (2007) *Neuron* **54** 263.
33. Bo et al. (2003) *Mol.Pharmacol.* **63** 1407.
34. Bianchi et al. (1999) *Eur.J.Pharmacol.* **376** 127.
35. Chessell et al. (1998) *Br.J.Pharmacol.* **124** 1314.
36. Donnelly-Roberts et al. (2009) *Br.J.Pharmacol.* **157** 1203.
37. Michel et al. (2008) *Br.J.Pharmacol.* **153** 737.
38. Stokes et al. (2006) *Br.J.Pharmacol.* **149** 880.
39. Michel et al. (2008) *Br.J.Pharmacol.* **155** 738.
40. Cattaneo (2010) *Circulation* **121** 171.
41. King et al. (1999) *Br.J.Pharmacol.* **128** 981.
42. Eliahu et al. (2009) *Eur.J.Med.Chem.* **44** 1525.
43. Nandanani et al. (1999) *J.Med.Chem.* **42** 1625.
44. Kim et al. (2003) *J.Med. Chem.* **46** 4974.
45. Cattaneo et al. (2004) *Biochem.Pharmacol.* **68** 1995.
46. Hechler et al. (2006) *J.Pharm.Exp.Therap.* **316** 556.
47. Pfefferkorn et al. (2008) *Bioorg.Med.Chem.Lett* **18** 3338.
48. Gao et al. (2004) *Biochem.Pharmacol.* **68** 231.
49. Hou et al. (2002) *Am.J.Phys.Heart and Circ.Phys.* **282** 784.
50. Jacobson et al. (2006) *Biochem.Pharmacol.* **71** 540.
51. Ivanov et al. (2007) *J.Med.Chem.* **50** 1166.
52. Shaver et al. (2005) *Purinerg. Sig.* **1** 183.
53. Ko et al. (2008) *Bioorg.Med.Chem.* **16** 6319.
54. Kindon et al. (1998) *World Pat.* **98** 54180.
55. Korcok et al. (2005) *J.Biol.Chem.* **280** 16909.
56. El-Tayeb et al. (2006) *J.Med.Chem.* **49** 7076.
57. Besada et al. (2006) *J.Med.Chem.* **49** 5532-5543.
58. Mamedova et al. (2004) *Biochem. Pharmacol.* **67** 1763.
59. Communi et al. (1999) *Br.J.Pharmacol.* **128** 1199.
60. Ullmann et al. (2005) *J.Med.Chem.* **48** 7040.
61. Savi et al. (2002) *Throm.Haemostas.* **84** 891.
62. Jakubowski et al. (2007) *Br.J.Clin.Pharmacol.* **63** 421.
63. Springthorpe et al. (2007) *Bioorg.Med.Chem.Lett.* **17** 6013.
64. Ye et al. (2008) *ChemMedChem* **3** 732.
65. Douglass et al. (2008) *J.Med.Chem.* **51** 1007.
66. Kim et al. (2005) *Biochem.Pharmacol.* **70** 266.
67. Das et al. (2010) *J.Med.Chem.* **53** 471.

## P2X and P2Y Receptor Compounds Available from Tocris

Cat. No.	Product Name	Primary Action
<b>Non-selective P2 Receptor</b>		
Agonists		
3245	<a href="#">ATP</a>	P2 agonist
4080	<a href="#">ATP<sub>γ</sub>S</a>	Non-selective P2 agonist; analog of ATP (Cat. No. 3245)
3312	<a href="#">BzATP</a>	P2X <sub>7</sub> agonist; also P2X <sub>1</sub> and P2Y <sub>1</sub> partial agonist; Photoaffinity label for ATPase
5157	<a href="#">DMNPE-caged ATP</a>	Caged ATP
3209	<a href="#">α,β-Methyleneadenosine 5'-triphosphate</a>	Non-selective P2 agonist
1062	<a href="#">2-Methylthioadenosine triphosphate (2-MeSATP)</a>	Non-selective P2 agonist
Antagonists		
2450	<a href="#">NF 157</a>	Selective P2Y <sub>11</sub> and P2X <sub>1</sub> antagonist
0625	<a href="#">PPADS tetrasodium salt</a>	Non-selective P2 antagonist
1472	<a href="#">Suramin hexasodium salt</a>	Non-selective P2 antagonist
<b>P2X Receptor</b>		
Antagonists		
6493	<a href="#">A 317491 sodium salt</a>	Selective, high affinity P2X <sub>3</sub> and P2X <sub>2/3</sub> antagonist; antinociceptive
2972	<a href="#">A 438079</a>	Competitive P2X <sub>7</sub> antagonist
3701	<a href="#">A 740003</a>	Potent and selective P2X <sub>7</sub> antagonist
4473	<a href="#">A 804598</a>	Potent and selective P2X <sub>7</sub> antagonist
4232	<a href="#">A 839977</a>	Potent P2X <sub>7</sub> antagonist
3323	<a href="#">AZ 10606120</a>	Potent P2X <sub>7</sub> antagonist
3317	<a href="#">AZ 11645373</a>	Potent and selective human P2X <sub>7</sub> antagonist
3579	<a href="#">5-BDBD</a>	Potent and selective P2X <sub>4</sub> antagonist
5545	<a href="#">BX 430</a>	Selective P2X <sub>4</sub> allosteric agonist
0845	<a href="#">Evans Blue tetrasodium salt</a>	Selective P2X antagonist; also non-NMDA iGluR antagonist
5299	<a href="#">JNJ 47965567</a>	Potent and selective P2X <sub>7</sub> antagonist; brain penetrant
1277	<a href="#">KN-62</a>	Non-competitive P2X <sub>7</sub> antagonist
1240	<a href="#">NF 023</a>	Selective, competitive P2X antagonist
2548	<a href="#">NF 110</a>	Potent P2X <sub>3</sub> antagonist
1199	<a href="#">NF 279</a>	Potent and selective P2X <sub>1</sub> antagonist
1391	<a href="#">NF 449</a>	Highly selective P2X <sub>1</sub> antagonist
0683	<a href="#">iso-PPADS tetrasodium salt</a>	P2X antagonist
1309	<a href="#">PPNDS</a>	Potent and selective P2X <sub>1</sub> antagonist
2188	<a href="#">Ro 0437626</a>	Selective P2X <sub>1</sub> antagonist
4391	<a href="#">Ro 51</a>	Potent P2X <sub>3</sub> and P2X <sub>2/3</sub> antagonist
3052	<a href="#">RO-3</a>	Selective P2X <sub>3</sub> and P2X <sub>2/3</sub> antagonist
2464	<a href="#">TNP-ATP</a>	Potent and selective P2X antagonist
Modulators		
3385	<a href="#">GW 791343</a>	P2X <sub>7</sub> allosteric modulator
1260	<a href="#">Ivermectin</a>	Positive allosteric modulator of P2X <sub>4</sub> receptor; also positive allosteric modulator of α7 nAChRs
5156	<a href="#">Ivermectin B1a-d<sub>2</sub></a>	Deuterated Ivermectin (Cat. No. 1260)
Other		
1203	<a href="#">MRS 2219</a>	Potentiates P2X <sub>1</sub> -mediated responses
<b>P2Y Receptors</b>		
Agonists		
1624	<a href="#">2-Methylthioadenosine diphosphate (2-MeSADP)</a>	Potent agonist at P2Y <sub>1</sub> , P2Y <sub>12</sub> and P2Y <sub>13</sub>
2157	<a href="#">MRS 2365</a>	Highly potent and selective P2Y <sub>1</sub> agonist
2915	<a href="#">MRS 2690</a>	Potent P2Y <sub>14</sub> agonist
2502	<a href="#">MRS 2693 trisodium salt</a>	Selective P2Y <sub>8</sub> agonist

Cat. No.	Product Name	Primary Action
3884	<a href="#">MRS 2768 tetrasodium salt</a>	Selective P2Y <sub>2</sub> agonist
4260	<a href="#">MRS 2957</a>	Potent and selective P2Y <sub>6</sub> agonist
4261	<a href="#">MRS 4062</a>	Selective P2Y <sub>4</sub> agonist
3892	<a href="#">NF 546</a>	Selective P2Y <sub>11</sub> agonist
2715	<a href="#">PSB 0474</a>	Potent and selective P2Y <sub>6</sub> agonist
4333	<a href="#">PSB 1114</a>	Potent and selective P2Y <sub>2</sub> agonist
3279	<a href="#">UTPyS trisodium salt</a>	Selective P2Y <sub>2/4</sub> agonist
<b>Antagonists</b>		
4890	<a href="#">AR-C 118925XX</a>	Selective and competitive P2Y <sub>2</sub> antagonist
3321	<a href="#">AR-C 66096 tetrasodium salt</a>	Potent and selective P2Y <sub>12</sub> antagonist
5720	<a href="#">AR-C 69931</a>	Highly potent P2Y <sub>12</sub> antagonist
6085	<a href="#">AZD 1283</a>	High affinity P2Y <sub>12</sub> antagonist
6078	<a href="#">BPTU</a>	P2Y <sub>1</sub> allosteric antagonist
1820	<a href="#">(+)-Clopidogrel hydrogen sulfate</a>	Selective P2Y <sub>12</sub> antagonist; active enantiomer of (±)-clopidogrel hydrochloride (Cat. No. 2490)
5316	<a href="#">Elinogrel</a>	P2Y <sub>12</sub> antagonist
0900	<a href="#">MRS 2179 tetrasodium salt</a>	Selective P2Y <sub>1</sub> antagonist
2402	<a href="#">MRS 2211</a>	Competitive P2Y <sub>13</sub> antagonist
2158	<a href="#">MRS 2279</a>	Selective, high affinity P2Y <sub>1</sub> antagonist
2159	<a href="#">MRS 2500</a>	Highly potent and selective P2Y <sub>1</sub> antagonist
2146	<a href="#">MRS 2578</a>	Selective P2Y <sub>6</sub> antagonist
4862	<a href="#">PPTN</a>	High affinity and selective P2Y <sub>14</sub> antagonist
6317	<a href="#">Prasugrel</a>	Irreversible P2Y <sub>12</sub> antagonist; antiplatelet and orally active
3983	<a href="#">PSB 0739</a>	Highly potent P2Y <sub>12</sub> antagonist
5650	<a href="#">SAR 216471</a>	Potent P2Y <sub>12</sub> antagonist; orally available
6864	<a href="#">Ticagrelor</a>	High affinity and reversible P2Y <sub>12</sub> antagonist; displays inverse agonism <i>in vitro</i> ; also inhibits ENT1
3931	<a href="#">Ticlopidine</a>	Selective P2Y <sub>12</sub> antagonist
3111	<a href="#">UDP disodium salt</a>	Competitive antagonist at P2Y <sub>14</sub> receptors; endogenous ligand
<b>Ectonucleotidases (NTPDase)</b>		
<b>Inhibitors</b>		
3633	<a href="#">Adenosine 5'-(<math>\alpha,\beta</math>-methylene)diphosphate sodium salt</a>	Ecto-5'-nucleotidase (CD73) inhibitor
1283	<a href="#">ARL 37156 trisodium salt</a>	NTPDase inhibitor
2689	<a href="#">POM 1</a>	Inhibitor of E-NTPDases
2574	<a href="#">PSB 06126</a>	NTPDase 3 inhibitor
2573	<a href="#">PSB 069</a>	Non-selective NTPDase inhibitor
<b>Nucleoside Transporters</b>		
<b>Inhibitors</b>		
1745	<a href="#">5-Iodotubercidin</a>	Nucleoside transporter inhibitor; also a broad spectrum kinase inhibitor
4588	<a href="#">8MDP</a>	Potent equilibrative nucleoside transporter (ENT) inhibitor
2924	<a href="#">NBMPR</a>	Equilibrative nucleoside transporter 1 (ENT1) inhibitor
5100	<a href="#">TC-T 6000</a>	Potent equilibrative nucleoside transporter 4 (ENT4) inhibitor

For a complete and up-to-date product listing please visit | [tocris.com](https://www.tocris.com)

**R&D** SYSTEMS

**NOVUS**  
BIOLOGICALS

**TOCRIS**

protein**simple**

**ACD**<sup>™</sup>

**bio**techne<sup>®</sup>

Global [info@bio-techne.com](mailto:info@bio-techne.com) [bio-techne.com/find-us/distributors](https://www.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](mailto:info.cn@bio-techne.com) TEL +86 (21) 52380373

[bio-techne.com](https://www.bio-techne.com)

