

Histamine Receptors

Iwan de Esch and Rob Leurs

Leiden/Amsterdam Center for Drug Research (LACDR), Division of Medicinal Chemistry, Faculty of Sciences, Vrije Universiteit Amsterdam, De Boelelaan 1083, 1081 HV, Amsterdam, The Netherlands

Dr. Iwan de Esch is an assistant professor and Prof. Rob Leurs is full professor and head of the Division of Medicinal Chemistry of the Leiden/Amsterdam Center of Drug Research (LACDR), VU University Amsterdam, The Netherlands. Since the seventies, histamine receptor research has been one of the traditional themes of the division. Molecular understanding of ligand-receptor interaction is obtained by combining pharmacology (signal transduction, proliferation), molecular biology, receptor modelling and the synthesis and identification of new ligands.

Introduction

Histamine is one of the aminergic neurotransmitters and plays an important role in the regulation of several (patho)physiological processes. In the mammalian brain histamine is synthesised in restricted populations of neurons that are located in the tuberomammillary nucleus of the posterior hypothalamus.¹ These neurons project diffusely to most cerebral areas and have been implicated in several brain functions (e.g. sleep/wakefulness, hormonal secretion, cardiovascular control, thermoregulation, food intake, and memory formation).² In peripheral tissues, histamine is stored in mast cells, eosinophils, basophils, enterochromaffin cells and probably also in some specific neurons. Mast cell histamine plays an important role in the pathogenesis of various allergic conditions. After mast cell degranulation, release of histamine leads to various well-known symptoms of allergic conditions in the skin and the airway system. In 1937, Bovet and Staub discovered compounds that antagonise the effect of histamine on these allergic reactions.³ Ever since, there has been intense research devoted towards finding novel ligands with (anti-) histaminergic activity. This research field has been fuelled by the consecutive discovery of four unique histamine receptor subtypes. Every receptor subtype has a very distinct (patho)physiological role and all of them belong to the superfamily of G-protein-coupled receptors (GPCRs). The histamine H₁ and H₂ receptor subtypes have proven to be excellent drug targets. Ligands for the histamine H₃ receptor subtype are currently entering clinical studies and the recently discovered histamine H₄ receptor subtype is subject of intense preclinical research.

The Histamine H₁ Receptor

Until the seventies, histamine research focused on the role of histamine in allergic diseases. This resulted in the development of several potent 'antihistamines' (e.g. mepyramine, see figure 1), which were useful in inhibiting pronounced symptoms of allergic conditions.⁴ However, the first generation of H₁ receptor antagonists that were developed for treating allergies revealed distinct side effects such as sedation. This particular physiological effect of the ligands was eliminated by structural modifications that prevent blood-brain-barrier penetration of the drugs. The first generation as well as the more recently developed antihistamines (originally termed antagonists but later reclassified as inverse agonists⁵) are clinically very successful and are widely available drugs. In 1991, the cDNA encoding a bovine H₁ receptor protein was cloned after an expression cloning strategy in *Xenopus oocytes*.⁶ The deduced amino acid sequence revealed a 491 amino acid protein of 56 kDa. Using the cDNA sequence encoding the bovine H₁ receptor, the cDNA sequences and intronless genes encoding the rat,^{3,7} guinea-pig,^{8,9} human^{10,11} and mouse¹² H₁ receptor proteins were cloned soon thereafter.

Contents

Introduction.....	1
The Histamine H ₁ Receptor	1
Ligands for H ₁ Receptors.....	2
The Histamine H ₂ Receptor	3
Ligands for H ₂ Receptors.....	3
The Histamine H ₃ Receptor	4
Ligands for H ₃ Receptors.....	4
The Histamine H ₄ Receptor	6
Ligands for H ₄ Receptors.....	7
Conclusion	7
References	8
Histamine Receptor Compounds.....	9

These receptor proteins are slightly different in length, highly homologous and do not show major pharmacological differences. Analysis of the 5'-flanking region of the human, rat and guinea-pig gene^{7,8,10} resulted in the identification of several DNA-binding motifs, including potential glucocorticoid responsive elements. The human H₁ receptor gene resides on chromosome 3.¹³ The H₁ receptor belongs to the large family of GPCRs. The receptor is associated with the phospholipase C-catalysed formation of inositol 1,4,5-triphosphate (IP₃) and 1,2-diacylglycerol (DAG). Histamine induces production of inositol phosphates in several tissues (including brain, airway, intestinal and vascular smooth muscle²⁴) via G_{αq} protein activation.¹⁴ In other tissues, activation of H₁ receptors can also stimulate adenylyl cyclase and formation of cAMP. Not all details of this signalling pathway are understood, for example questions remain about the G protein and the involvement of Ca²⁺.¹⁵ In any case, alternative signalling pathways can be mediated by the histamine H₁ receptor. Recent results indicate that the functional heterogeneity can be ligand-directed.¹⁶

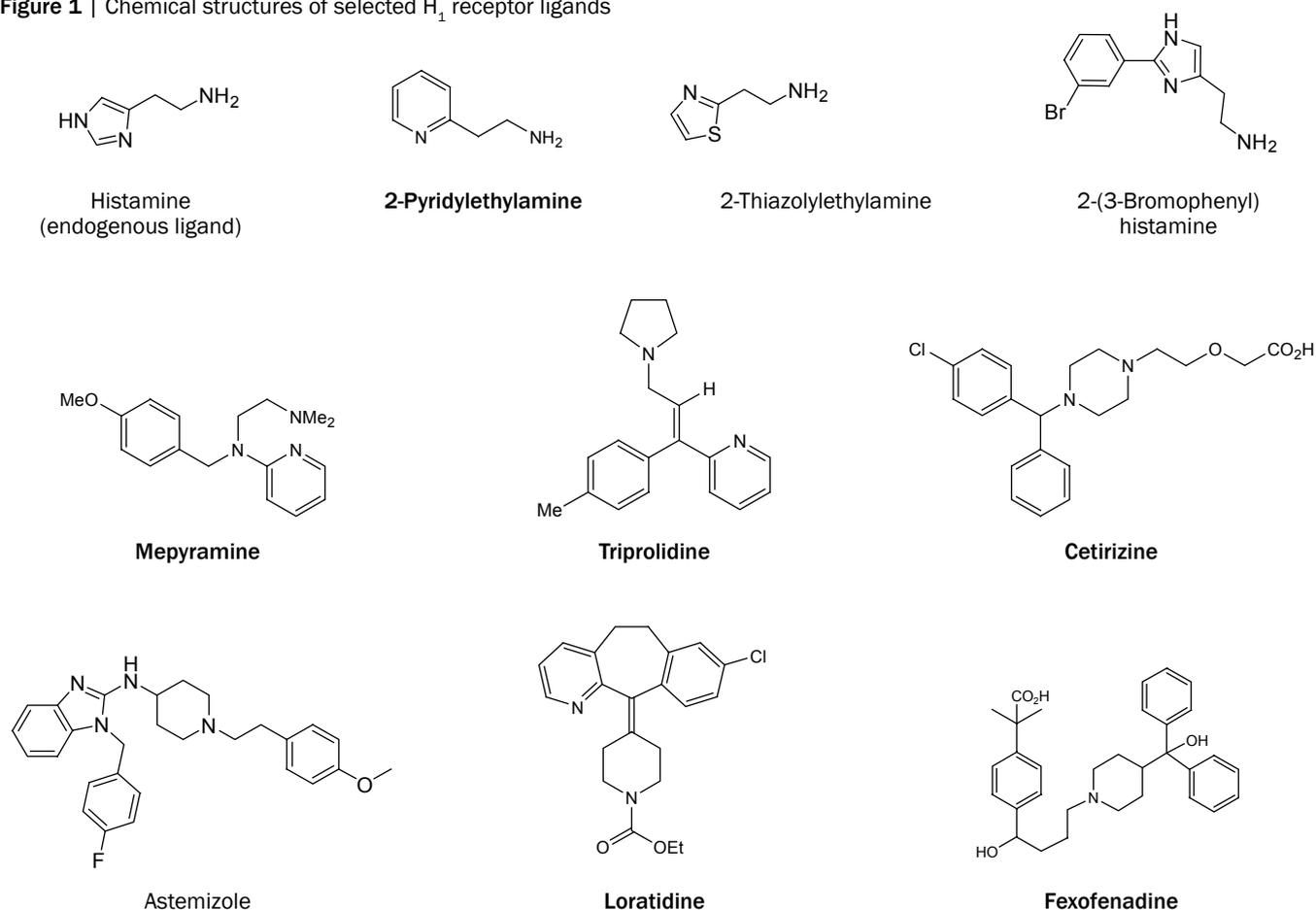
The histamine H₁ receptor is a well established drug target and has been thoroughly studied for decades. Nevertheless, H₁ receptor research continues to flourish as many new techniques and approaches are being developed by using this receptor as an archetypal GPCR target. Most notably, in the last few years, the histamine H₁ receptor has been the subject of various molecular biology studies (e.g., large-scale overproduction¹⁷ and GPCR

binding site elucidation¹⁸), biophysical approaches (such as solid-state NMR¹⁹) and investigations towards the general activation mechanisms of GPCRs.^{20,21}

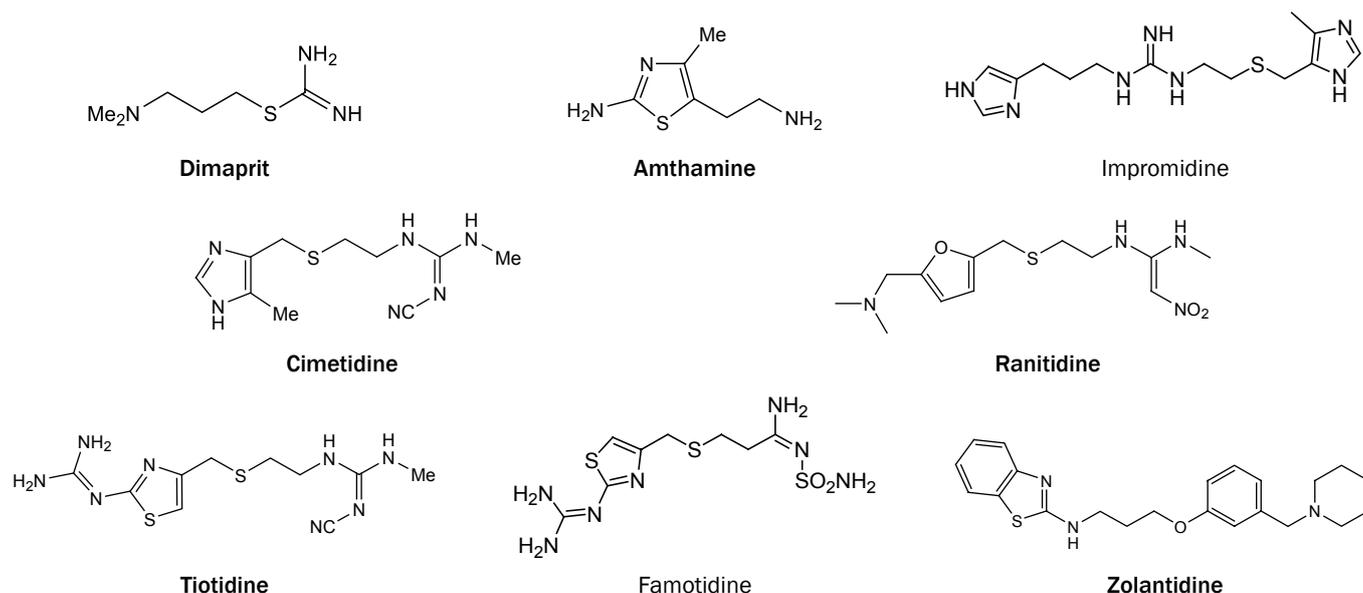
Ligands for H₁ Receptors

Modification of the imidazole moiety of histamine has been the most successful approach for obtaining selective H₁ agonists (figure 1). The presence of the tautomeric N⁻-N⁺ system of the imidazole ring is not obligatory, as reflected by the selective H₁ agonists 2-pyridylethylamine and 2-thiazolyethylamine. Substitution of the imidazole ring at the 2-position leads to relatively selective H₁ agonists. For example, 2-(*meta*-halogenated) phenylhistamines are relatively potent H₁ receptor agonists at the guinea-pig ileum;²² however, these compounds act as partial agonists in other systems.²³ A wide array of potent and selective H₁ antagonists are available.⁴ Compounds such as mepyramine (also called pyrilamine) and triprolidine are highly potent H₁ antagonists and very useful tools for pharmacological investigations. [³H]-mepyramine is, for example, successfully used as an H₁ receptor radioligand.²⁴ These so-called classical 'antihistamines' easily penetrate the brain and are therefore also useful in *in vivo* CNS studies.² Elimination of the blood-brain-barrier passage by some minor structural modifications has resulted in many new, non-sedating H₁ antagonists (e.g., cetirizine, astemizole, fexofenadine and loratidine).⁴

Figure 1 | Chemical structures of selected H₁ receptor ligands



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

Figure 2 | Chemical structures of selected H₂ receptor ligands

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

The Histamine H₂ Receptor

The observation that the classical ‘antihistamines’ (i.e. H₁ receptor inverse agonists) cannot antagonise all histamine-induced effects (e.g. at the stomach and the heart), led Ash and Schild in 1966 to propose two distinct subtypes of histamine receptors: H₁ and H₂.²⁵ This hypothesis became generally accepted when Black *et al*²⁶ succeeded in the synthesis of a series of new compounds (e.g. burimamide, cimetidine), which were able to block the effects of histamine on the stomach and the heart. These H₂ receptor antagonists proved to be very useful in the therapy of gastric ulcers. Gantz *et al*²⁷ were the first to clone a cDNA encoding a 359 amino acid H₂ receptor. Using degenerate primers based on the known sequence similarity of various GPCRs, the H₂ receptor sequence was obtained from canine gastric parietal cDNA by PCR. Soon thereafter, the intronless genes encoding the rat,²⁸ human,²⁹ guinea pig³⁰ and mouse³¹ H₂ receptor were cloned by means of homology screening. Identification of the promoter region of the human H₂ receptor gene revealed the existence of regulatory transcription sites and regions displaying stimulatory and inhibitory effects on gene expression monitored in a luciferase assay.³² Studies have indicated that the human H₂ receptor gene resides on chromosome 5.³⁰ Interestingly, several polymorphisms have been found in the human H₂ receptor gene³³ and one of the mutations has been linked to schizophrenia.³⁴ The histamine H₂ receptor is coupled to the adenylate cyclase system in a variety of tissues (e.g. brain, stomach, heart, gastric mucosa, lung).²⁴ Moreover, cell lines transfected with the cloned H₂ receptor genes showed an H₂ receptor-mediated increase of cAMP.^{35,36,37} In addition, alternative signalling pathways for the H₂ receptor have been identified. In differentiated HL-60 cells and CHO or HEPA cells transfected with the H₂ receptor cDNA, an H₂ receptor-mediated increase of the intracellular Ca²⁺ concentration and/or IP₃ levels

was observed.^{38,39} Moreover, in CHO cells expressing the rat H₂ receptor, activation of the H₂ receptor resulted in an inhibition of the release of arachidonic acid induced by either constitutive purinergic receptors or a Ca²⁺-ionophore,³⁶ as well as an increase in cAMP. These new signal transduction pathways are both regulated via unknown, cAMP-independent pathways.

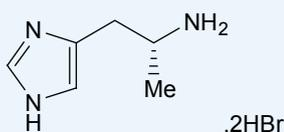
Ligands for H₂ Receptors

The first selective H₂ receptor agonist was dimaprit (figure 2). This compound is almost as active as histamine at the H₂ receptor but hardly displays any H₁ receptor activity.⁴⁰ Much later, it was shown that dimaprit is also a moderate H₃ receptor antagonist⁴¹ and a moderate H₄ receptor agonist.⁴² Amthamine can be considered a rigid dimaprit analogue.⁴³ This compound combines a high H₂ receptor selectivity with a potency which is slightly higher compared to histamine, both *in vitro* and *in vivo*.⁴⁴ An H₂ receptor agonist that is more potent than histamine is the guanidine derivative impromidine. This ligand actually combines a rather high H₂ receptor affinity with a reduced efficacy. Impromidine also shows moderate H₁- and potent H₃-receptor antagonistic activity^{45,46} as well as potent H₄ receptor partial agonistic activity.⁴²

The finding that N α -guanylhistamine acts as a partial H₂ agonist in a gastric acid secretion test did lead to the development of the relatively weak H₂ antagonist burimamide. Years later, it was shown that burimamide is also an H₃ and H₄ receptor partial agonist.⁴² Nevertheless, burimamide was a good lead for the development of selective and clinically useful H₂ receptor antagonists, such as cimetidine.^{45,47,48} The 4-methylimidazole moiety of cimetidine can easily be replaced by other heterocyclic groups (figure 2). Replacement by a substituted furan- (e.g. ranitidine) or thiazole ring (e.g. tiotidine and famotidine) leads to compounds that are usually more potent at the H₂ receptor than

(R)-(-)- α -Methylhistamine, High Affinity H₃ Agonist**(R)-(-)- α -Methylhistamine**

Cat. No. 0569



This potent and high affinity H₃ agonist displays >200-fold selectivity over H₄ receptors. The compound inhibits H₃ receptor-mediated histamine synthesis and release in the CNS and stimulates H₄ receptor-mediated eosinophil shape change (EC₅₀ = 66 nM).

Oishi *et al.* (1989) Effects of histamine H₃-agonist (R)- α -methylhistamine and the antagonist thioperamide on histamine modulation in the mouse and rat brain. *J.Neurochem.* **52** 1388. Hew *et al.* (1990) Characterization of histamine- H₃ receptors in guinea pig ileum with H₃-selective ligands. *Br.J.Pharmacol.* **101** 621. Schwartz *et al.* (1990) A third histamine receptor subtype - characterization, localization and functions of the H₃-receptor. *Agents Actions* **30** 13. Buckland *et al.* (2003) Histamine induces cytoskeletal changes in human eosinophils via the H₄ receptor. *Br.J.Pharmacol.* **140** 1117.

cimetidine. Moreover, the replacement of the imidazole moiety also eliminates the undesired inhibition of cytochrome P450.⁴⁸ The potent tritiated H₂ antagonist tiotidine and [¹²⁵I]-iodinated H₂ antagonist iodoaminopotentidine are successfully used as radioligands for the H₂ receptor.²⁴ The newly developed brain-penetrating H₂ antagonist zolantidine is an important tool for *in vivo* CNS studies.⁴⁹ The H₂ receptor was reported to be spontaneously active in transfected CHO cells.⁵⁰ Based on this concept, many H₂ antagonists were reclassified; cimetidine, ranitidine and famotidine are in fact inverse agonists, whereas burimamide acts in this model system as a neutral antagonist.⁵⁰

The Histamine H₃ Receptor

In the eighties, another physiological role of histamine became apparent, namely its role as a neurotransmitter. In 1983, Arrang and co-workers discovered that the inhibitory effect of histamine on its own release and synthesis was not mediated by the known H₁ and H₂ receptor subtypes as no correlation with either the H₁ or the H₂ receptor activity of known histaminergic compounds was observed.⁴⁶ Soon after, the H₃ receptor agonist (R)-(-)- α -methylhistamine and the antagonist thioperamide (see figure 3) were developed.⁵¹ It was confirmed that this receptor subtype indeed regulates the release and synthesis of histamine and in addition has a regulatory role in the release of other neurotransmitters, such as serotonin, noradrenalin and dopamine.⁵² Next to high expression in certain regions of the CNS (for example the basal ganglia, hippocampus and cortical areas, i.e. the parts of the brain that are associated with cognition) the H₃ receptor is present in the peripheral nervous system, e.g. in the gastrointestinal tract, the airways and the cardiovascular system.^{53,54} Initial efforts to identify the H₃ receptor gene, using the anticipated homology with the identified H₁ and H₂ receptor gene all failed. Eventually, the human H₃ receptor cDNA was cloned by Lovenberg and co-workers in 1999.⁵⁵ In search of novel GPCRs through homology searching of expressed sequence tag databases, a receptor with high similarity to the M₂ muscarinic acetylcholine receptor was identified. Expression of the gene and full characterisation established this protein as the histamine H₃ receptor. The cloning of the H₃ receptor of other species, including

rat,⁵⁶ guinea pig⁵⁷ and mouse⁵⁸ soon followed and it was revealed that major H₃ receptor species differences exist. The human H₃ receptor gene was assigned to the telomeric region of the q arm of chromosome 20 and contains three exons that are interrupted by two introns.⁵⁹ The H₃ receptor gene can undergo extensive alternative splicing, resulting in many H₃ receptor isoforms that have different signalling properties and expression profiles.^{59,60,61} It was shown that the H₃ receptor displays particularly high constitutive activity, again leading to a reclassification of existing ligands into agonists, neutral antagonists and inverse agonists.^{62,63} Gbahou and co-workers revealed that the compound proxyfan (figure 3) is a protean agonist both *in vitro* and *in vivo*, meaning that this remarkable compound behaves as agonist, neutral antagonist or inverse agonist, depending on the signalling pathway studied.⁶⁴

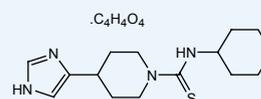
The H₃ receptor signals through G_{v/o} proteins as was shown by the pertussis toxin sensitive stimulation of [³⁵S]-GTP γ S binding in rat cortical membranes.⁶⁵ Through negative coupling to adenylyl cyclase, stimulation of the H₃ receptor results in lower levels of cAMP, thereby reducing downstream signalling events such as CREB-dependent gene transcription.⁵⁶ Alternative signalling pathways may be activated by the G_{v/o} proteins, including mitogen-activated protein kinase (MAPK)⁶⁰ and phosphatidylinositol 3-kinase (PI3K) pathways. G_{v/o} protein activation can also lead to the activation of phospholipase A₂ (PLA₂) to induce the release of arachidonic acid, the lowering of intracellular Ca²⁺ levels through voltage-gated ion channels⁶⁷ and the inhibition of the Na⁺/H⁺ exchanger (NHE).⁶⁸ With the recent progress that has been made in the characterisation of the H₃ receptor (as outlined above), many pharmaceutical companies have very active H₃ receptor drug development programs. As a result, several H₃ receptor ligands have entered clinical studies for a plethora of applications, including obesity, narcolepsy, dementia and migraine (among others).^{69,70}

Ligands for H₃ Receptors

At the H₃ receptor, histamine itself is a highly active agonist. Modification of the endogenous ligand by mono- or dimethylation of the amino function results in compounds that are more active

Thioperamide, a Brain-penetrant H₃ and H₄ Ligand**Thioperamide**

Cat. No. 0644



Thioperamide, the potent histamine H₃ and H₄ ligand, displays K_i values of 25 and 27 nM at recombinant H₃ and H₄ receptors respectively. The compound acts as an antagonist at H₃ receptors and displays inverse agonist activity at H₄ receptors. It freely crosses the blood-brain barrier.

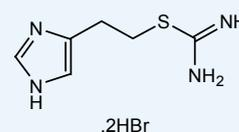
Hew *et al.* (1990) Characterisation of histamine H₃ receptors in guinea pig ileum with H₃ selective ligands. *Br.J.Pharmacol.* **101** 621. Liu *et al.* (2001) Cloning and pharmacological characterization of a fourth histamine receptor (H₄) expressed in bone marrow. *Mol. Pharmacol.* **59** 420. Ling *et al.* (2004) Histamine H₄ receptor mediates eosinophil chemotaxis with cell shape change and adhesion molecule upregulation. *Br.J.Pharmacol.* **142** 161. Gbahou *et al.* (2006) Compared pharmacology of human H₃ and H₄ receptors: structure-activity relationships of histamine derivatives. *Br.J.Pharmacol.* **147** 744.

and selective for the H₃ receptor. Methylation of the α-carbon atom of the ethylamine sidechain drastically increases the potency at the H₃ receptor. This increased activity resides completely in the R-isomer; the corresponding S-isomer is approximately 100-fold less potent. Since the methylation leads to highly reduced activity at both H₁ and H₂ receptors, (R)-(-)-α-methylhistamine (figure 3) was initially considered a selective agonist at the H₃ receptor. However, when the H₄ receptor was discovered it was shown that (R)-(-)-α-methylhistamine also has considerable affinity for this histamine receptor subtype.⁴² Nevertheless, in combination with its less active S-isomer, (R)-(-)-α-methylhistamine has proven to be highly useful for the pharmacological characterisation of H₃ receptor-mediated effects.⁴¹ Tritiated forms of Nα-methylhistamine and (R)-(-)-α-methylhistamine are available as radiolabelled agonists for the H₃ receptor. For potent H₃ agonism, the amine function of histamine can be replaced by an isothioureia group, as in imetit. This compound is also very active *in vitro* and *in vivo*,^{71,72,73} as is (R)-(-)-α-methylhistamine. The amine function can also be incorporated in ring structures. For example, immepip is a potent H₃ receptor agonist that is effective *in vitro* and *in vivo*.⁷⁴ Although the described first generation H₃ agonists were intensively used as reference ligands to study the H₃ receptor, all of them proved to also have considerable activity at the H₄ receptor. Therefore, a new generation of potent and selective H₃ receptor agonists has been developed, most notably immethridine⁷⁵ (pEC₅₀ = 9.74; displays 300-fold selectivity over the H₄ receptor) and

Imetit, High Affinity H₃/H₄ Agonist

Imetit

Cat. No. 0729



.2HBr

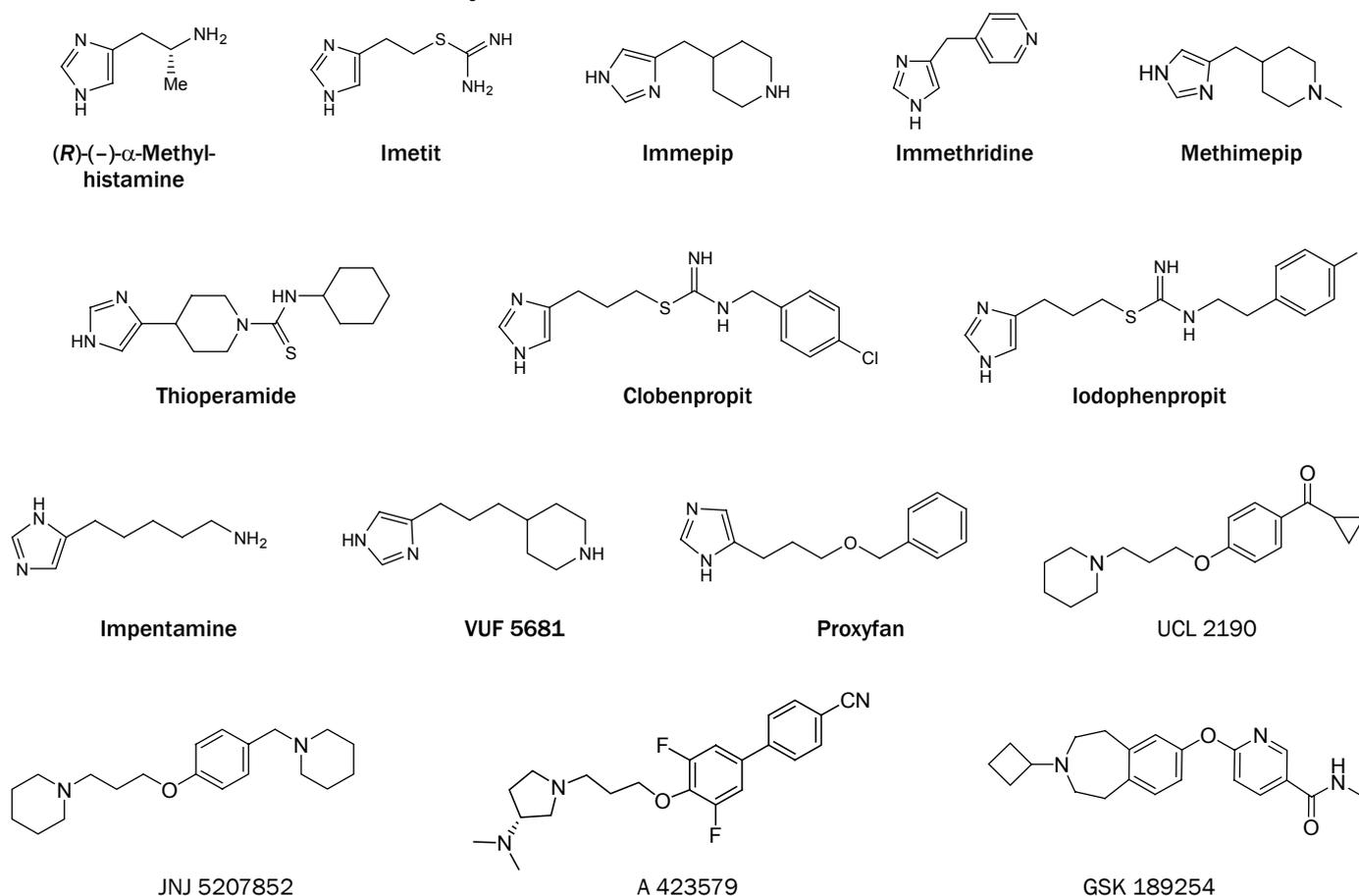
Imetit is an extremely high affinity, potent agonist at H₃ and H₄ receptors (K_i values are 0.3 and 2.7 nM respectively). The agonist induces shape change in eosinophils *in vitro* with an EC₅₀ of 25 nM and is centrally active following systemic administration *in vivo*.

Garbarg et al. (1992) S-[2-(4-Imidazolyl)ethyl]isothioureia, a highly specific and potent histamine H₃ receptor agonist. *J.Pharmacol.Exp.Ther.* **263** 304. **Farzin and Attarzadeh** (2000) Influence of different histamine receptor agonists and antagonists on apomorphine-induced licking behavior in rats. *Eur.J.Pharmacol.* **404** 169. **Ling et al.** (2004) Histamine H₄ receptor mediates eosinophil chemotaxis with cell shape change and adhesion molecule upregulation. *Br.J.Pharmacol.* **142** 161.

methimepip.⁷⁶ These latter compounds are devoid of high H₄ receptor activity.

As with the first generation H₃ receptor agonists, the first generation H₃ receptor antagonists (all of them possessing an imidazole heterocycle) turned out to have affinity for the H₄ receptor. The first potent H₃ receptor antagonist (later reclassified as an inverse agonist) that was devoid of H₁ receptor and H₂

Figure 3 | Chemical structures of selected H₃ receptor ligands



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

receptor activity was thioperamide (figure 3).⁵¹ This compound has been used in many H₃ receptor studies as a reference ligand and is active *in vitro* and *in vivo* (the compound is able to penetrate the CNS).⁷⁷ However, thioperamide displays some 5-HT₃ receptor antagonism⁷⁸ and is an inverse agonist at the H₄ receptor. The remarkable H₃ receptor species difference can be demonstrated with thioperamide as the compound has a 10-fold higher affinity for the rat H₃ receptor than for the human H₃ receptor.⁵⁶ Based on the H₃ receptor agonist imetit (*vide ante*), the highly potent H₃ receptor inverse agonist clobenpropit was developed (pA₂ = 9.9).⁷³ This compound has some 5-HT₃ receptor activity⁷⁸ and displays partial agonist activity at H₄ receptors. In addition, radioligands for the H₃ receptor have been described (e.g. [¹²⁵I]-iodophenpropit and [¹²⁵I]-iodoproxyfan).^{79,80} The moderately active H₂ receptor antagonist burimamide (pA₂ = 5.1, figure 2) also has good affinity for the H₃ (pK_i = 7.9) and the H₄ (pK_i = 7.4) receptor.⁴² Impentamine is a potent histamine H₃ receptor inverse agonist (pA₂ = 8.4). Like burimamide, this compound can act as a partial agonist in SK-N-MC cells expressing human H₃ receptors. It has also been shown that small structural modifications of impentamine, i.e. alkylation of the primary amine moiety of impentamine with e.g. methyl-, isopropyl- and p-chlorobenzyl- groups results in ligands that cover the complete spectrum of functional activity, i.e. agonism, neutral antagonism and inverse agonism.^{63,81} The compound VUF 5681 was reported as a neutral H₃ antagonist (or 'silent' antagonist),⁸² findings that were recently confirmed in studies investigating constitutive activity using rat brain cortex.⁸³ It has been shown that the functional activity of proxyfan depends on the system used, ranging from full agonist to inverse agonist. Thus, proxyfan can be classified as a protean agonist, both *in vitro* and *in vivo*.^{64,84}

The first non-imidazole H₃ receptor ligand was reported by Ganellin in 1998.⁸⁵ This elegant medicinal chemistry work did lead to the potent compound UCL 2190.⁸⁶ Following the cloning of the H₃ receptor in 1999, several pharmaceutical companies entered the H₃ research field and further explored this structural class. An example of these new ligands is JNJ 5207852, a compound that contains some typical structural features for H₃ receptor ligands, e.g. piperidiny propoxy side chain. This neutral antagonist is active in several models for cognition^{87,88} but does

Clobenpropit, Highly Potent H₃ Antagonist

Clobenpropit

Cat. No. 0752



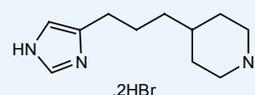
An extremely potent antagonist/inverse agonist at H₃ receptors (pA₂ = 9.93), clobenpropit also displays partial agonist activity at H₄ receptors. The ligand induces eosinophil shape change with an EC₅₀ of 3 nM.

Van der Goot et al. (1992) Isothiourea analogues of histamine as potent agonists or antagonists of the histamine H₃ receptor. *Eur.J.Med.Chem.* **27** 511.
Yokoyama et al. (1994) Clobenpropit (VUF-9153), a new histamine H₃ receptor antagonist, inhibits electrically induced convulsions in mice. *Eur.J.Pharmacol.* **260** 23.
Liu et al. (2001) Cloning and pharmacological characterization of a fourth histamine receptor (H₄) expressed in bone marrow. *Mol. Pharmacol.* **59** 420.
Buckland et al. (2003) Histamine induces cytoskeletal changes in human eosinophils via the H₄ receptor. *Br.J.Pharmacol.* **140** 1117.

VUF 5681, a Novel H₃ Antagonist

VUF 5681

Cat. No. 2493



VUF 5681 is a potent histamine H₃ receptor silent antagonist (pK_i = 8.35).

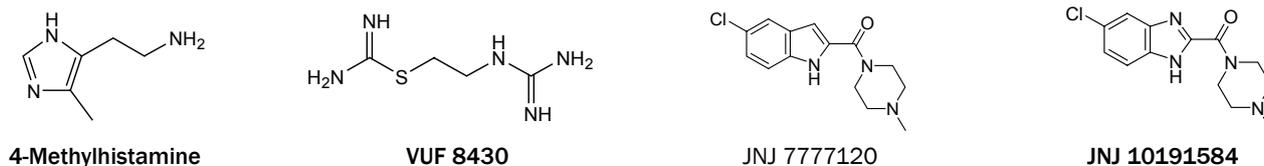
Kitbunnadaj et al. (2004) Identification of 4-(1*H*-imidazol-4(5*H*)-ylmethyl)pyridine (immethridine) as a novel, potent, and highly selective histamine H₃ receptor agonist. *J.Med.Chem.* **47** 2414.
Leurs et al. (2005) The histamine H₃ receptor: from gene cloning to H₃ receptor drugs. *Nat.Rev.Drug Discov.* **4** 107.
Moreno-Delgado et al. (2006) Constitutive activity of H₃ autoreceptors modulates histamine synthesis in rat brain through the cAMP/PKA pathway. *Neuropharmacology* **51** 517.

not act as an appetite suppressant and has no effect on food intake.⁸⁹ Other compounds, such as Abbott's A 423579 seem to have more efficacy in obesity models, while lacking clear cognitive effects.⁹⁰ At the time of writing the differences in efficacy for distinct clinical applications of the different classes of H₃ ligands is not understood and subject of intense research.^{70,91} Interestingly, GSK 189254 has been in trials for three different diseases: neuropathic pain, narcolepsy and dementia.⁹²

The Histamine H₄ Receptor

Immediately following the cloning of the H₃ receptor gene, several groups identified the homologous H₄ receptor sequence in the human genome databases.⁹³⁻⁹⁷ Indeed, the H₄ receptor has high sequence identity with the H₃ receptor (31% at the protein level, 54% in the transmembrane domains). The H₃ and H₄ receptor are also similar in gene structure. The human H₄ receptor gene is located on chromosome 18q11.2 in a single copy per haploid genome. The gene spans more than 21 kbp and contains three exons that are interrupted by two large introns.⁹⁸ To date, two H₄ receptor isoforms have been identified. Cloning of the genes that encode the mouse, rat, guinea-pig and pig H₄ receptors reveal only limited sequence homology with the human H₄ receptor. The H₄ receptor is mainly expressed in bone marrow and peripheral leukocytes, and mRNAs of the human H₄ receptor are detected in e.g. mast cells, dendritic cells, spleen and eosinophils.⁹³⁻⁹⁷ The receptor has a pronounced effect on the chemotaxis of several cell types that are associated with immune and inflammatory responses. The H₄ receptor is mainly coupled to G_{i/o} proteins, thereby leading to a decrease in the production of cAMP and subsequent downstream effects such as regulation of cAMP responsive element-binding protein (CREB) gene transcription. Furthermore, H₄ receptor stimulation affects the pertussis-toxin-sensitive activation of mitogen-activated protein (MAP) kinase pathways. Studying the increased levels of [³⁵S]GTPγS levels in H₄ transfected cells, it has been shown that the H₄ receptor is constitutively active. The Gβγ subunits of the G_{i/o} proteins activate phospholipase C, and thereby increase the Ca²⁺ concentrations. In mast cell and eosinophils, this Ca²⁺ response can be linked to cellular chemotaxis.

Considering the physiological role of the H₄ receptor, several applications are under preclinical investigation,^{99,100} including allergy and asthma,¹⁰¹ as well as chronic inflammations such as

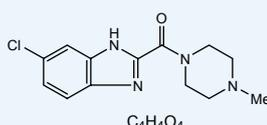
Figure 4 | The first selective H₄ receptor ligands reported in scientific literature

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

JNJ 10191584, H₄-Selective Antagonist

JNJ 10191584

Cat. No. 2441



JNJ 10191584 is a highly selective histamine H₄ receptor silent antagonist. It binds with high affinity to the human H₄ receptor ($K_i = 26$ nM) and is > 540-fold selective over the H₃ receptor ($K_i = 14.1$ μM). *In vitro* the antagonist inhibits mast cell and eosinophil chemotaxis with IC₅₀ values of 138 and 530 nM respectively. The antagonist is orally active *in vivo*.

Terzioglu *et al.* (2004) Synthesis and structure-activity relationships of indole and benzimidazole piperazines as histamine H₄ receptor antagonists. *Bioorg. Med. Chem. Lett.* **14** 5251. Venable *et al.* (2005) Preparation and biological evaluation of indole, benzimidazole, and thienopyrrole piperazine carboxamides: potent human histamine H₄ antagonists. *J. Med. Chem.* **48** 8289. Varga *et al.* (2005) Inhibitory effects of histamine H₄ receptor antagonists on experimental colitis in the rat. *Eur. J. Pharmacol.* **522** 130.

inflammatory bowel disease (IBD)¹⁰² and rheumatoid arthritis.¹⁰³ The H₄ receptor is also being associated with pruritus (itch)^{104,105} and is involved in the progression of colon cancer.¹⁰⁶

Ligands for H₄ Receptors

As was described above, most imidazole-containing, first generation H₃ receptor ligands have considerable affinity for the H₄ receptor as well. However, selective H₄ receptor ligands have been described. 4-Methylhistamine is a potent H₄ agonist while displaying more than a 100-fold selectivity over the other histamine receptor subtypes, including the H₂ receptor for which this ligand was originally developed.⁴² A slightly different and complimentary profile was reported for VUF 8430. This compound has a high H₄ receptor activity and affinity, minimal affinity for the H₁ and H₂ receptor and a 33-fold lower affinity for the H₃ receptor.¹⁰⁷

Potent and selective H₄ receptor antagonists are also emerging. The first reported neutral antagonist is JNJ 7777120, a compound that is frequently being used as H₄ receptor reference ligand. Currently, [³H]-JNJ 7777120 and [³H]-histamine are used as H₄ receptor radioligands. Unfortunately, JNJ 7777120 has a poor stability in human and rat liver microsomes and a limited half life of about two hours. The benzimidazole derivative JNJ 10191584 (VUF 6002) is also a neutral H₄ antagonist.¹⁰⁸ This compound is orally active *in vivo* and has an improved liver microsome stability but still a limited half life.^{102,109} More recently, 2-arylbenzimidazoles

have been described as ligands with low nanomolar affinity for the H₄ receptor.¹¹⁰ Considering the number of H₄ receptor-related patent applications that have recently been disclosed (as reviewed elsewhere¹⁰⁰), it can be anticipated that many new H₄ receptor ligands will be described in scientific literature in the near future.

Conclusions

The histamine receptor research field continues to blossom as both members of industry and academia find this family of receptors very rewarding, both scientifically and commercially. For more than seventy years, considerable efforts have been devoted to finding new ways to modulate the different physiological processes that are mediated by histamine. The consecutive discovery of new histamine receptor subtypes (distinction between H₁ and H₂ receptors in 1966, discovery of the H₃ receptor in 1983 and the H₄ receptor in 2000, note the interval of seventeen years) provides a complete and exciting research platform. The histamine receptor family also seems to hold the promise of GPCRs as excellent drug targets with two receptor subtypes (H₁ and H₂) addressed by blockbuster drugs, the third subtype (H₃) leading to frantic clinical studies and the latest addition to the family (H₄) leading to very interesting preclinical data. It can therefore be anticipated that histamine receptor research will continue to thrive in the years to come.

References

1. Schwartz et al. (1991) *Physiol.Rev.* **71** 1.
2. Yanai et al. (2007) *Pharmacol.Ther.* **113** 1.
3. Bovet et al. (1937) *C.R.Soc.Biol.Ses.Fil.* **123** 547.
4. Zhang et al. (1997) In: *Burger's Medicinal Chemistry and Drug Discovery*, Fifth Edition. Ed. Wolff, p 495. John Wiley & Sons, Inc.
5. Leurs et al. (2002) *Clin.Exp.Allergy* **32** 489.
6. Yamashita et al. (1991) *Proc.Natl.Acad.Sci.USA* **88** 11515.
7. Fujimoto et al. (1993) *Biochem.Biophys.Res.Comm.* **190** 294.
8. Horio et al. (1993) *J.Biol.Chem.* **114** 408.
9. Traiffort et al. (1994) *J.Neurochem.* **62** 507.
10. Fukui et al. (1994) *Biochem.Biophys.Res.Comm.* **201** 894.
11. De Backer et al. (1993) *Biochem.Biophys.Res.Comm.* **197** 1601.
12. Inoue et al. (1996) *Genomics* **36** 178.
13. Leconiat et al. (1994) *Hum.Genet.* **94** 186.
14. Hill et al. (1997) *Pharmacol.Rev.* **49** 253.
15. Leurs et al. (1994) *J.Neurochem.* **62** 519.
16. Noniri et al. (2004) *J.Pharmacol.Exp.Ther.* **311** 274.
17. Ratnala et al. (2004) *Eur.J.Biochem.* **271** 2636.
18. Bruysters et al. (2005) *Mol.Pharmacol.* **67** 1045.
19. Ratnala et al. (2007) *J.Am.Chem.Soc.* **129** 867.
20. Xie et al. (2006) *J.Pharmacol.Exp.Ther.* **317** 139.
21. Jongejan et al. (2005) *Nat.Chem.Biol.* **1** 371.
22. Zingel et al. (1995) In: *Progress in Drug Research*, Ed. Jucker, p 49, Birkhauser Verlag, Basel.
23. Leurs et al. (1994) *Biochem.Biophys.Res.Comm.* **201** 295.
24. Leurs et al. (1995) *Pharmacol.Ther.* **66** 413.
25. Ash and Schild (1966) *Br.J.Pharmacol.* **27** 427.
26. Black et al. (1972) *Nature* **236** 385.
27. Gantz et al. (1991) *Proc.Natl.Acad.Sci.USA* **88** 429.
28. Ruat et al. (1991) *Biochem.Biophys.Res.Comm.* **179** 1470.
29. Gantz et al. (1991) *Biochem.Biophys.Res.Comm.* **178** 1386.
30. Traiffort et al. (1995) *Biochem.Biophys.Res.Comm.* **211** 570.
31. Kobayashi et al. (1996) *Genomics* **37** 390.
32. Nishi et al. (1995) *Biochem.Biophys.Res.Comm.* **210** 616.
33. Orange et al. (1996) *Neuroreport* **7** 1293.
34. Orange et al. (1996) *Mol.Psychiatry* **1** 466.
35. Leurs et al. (1994) *Br.J.Pharmacol.* **112** 847.
36. Traiffort et al. (1992) *Proc.Natl.Acad.Sci.USA* **89** 2649.
37. Delvalle et al. (1992) *Am.J.Physiol.* **263** G967.
38. Mitsuhashi et al. (1989) *J.Biol.Chem.* **264** 18356.
39. Seifert et al. (1992) *Mol.Pharmacol.* **42** 227.
40. Parsons et al. (1977) *Agents Actions* **7** 31.
41. Leurs et al. (1995) In: *Progress in Drug Research* Ed. Jucker, p 107, Birkhauser Verlag, Basel.
42. Lim et al. (2005) *J.Pharmacol.Exp.Ther.* **314** 1310.
43. Eriks et al. (1992) *J.Med.Chem.* **35** 3239.
44. Coruzzi et al. (1996) *Arch.Pharmacol.* **353** 417.
45. Durant et al. (1978) *Nature* **276** 403.
46. Arrang et al. (1983) *Nature* **302** 832.
47. Ganellin (1992) In: *The Histamine Receptor*. Eds. Schwartz and Haas, p 1. Wiley-Liss, New York.
48. Van der Goot et al. (1991) In: *Handbook of Experimental Pharmacology*. Ed. Uvnas, p 573. Springer-Verlag, Berlin.
49. Young et al. (1988) *J.Med.Chem.* **31** 626.
50. Smit et al. (1996) *Proc.Natl.Acad.Sci.USA* **93** 6802.
51. Arrang et al. (1987) *Nature* **327** 117.
52. Schlicker et al. (1994) *Fund.Clin.Pharmacol.* **8** 128.
53. Martinez-Mir et al. (1990) *Brain Res.* **526** 322.
54. Bertaccini et al. (1998) In: *The histamine H3 receptor; a target for new drugs*. Eds. Leurs and Timmerman, p 59. Elsevier
55. Lovenberg et al. (1999) *Mol.Pharmacol.* **55** 1101.
56. Lovenberg et al. (2000) *J.Pharmacol.Exp.Ther.* **293** 771.
57. Tardivel-Lacombe et al. (2000) *Neuroreport* **11** 755.
58. Chen et al. (2003) *Eur.J.Pharmacol.* **467** 57.
59. Tardivel-Lacombe et al. (2001) *Neuroreport* **12** 321.
60. Drutel et al. (2001) *Mol.Pharmacol.* **59** 1.
61. Bakker et al. (2004) *Inflamm.Res.* **53** 509.
62. Morisset et al. (2000) *Nature* **408** 860.
63. Wieland et al. (2001) *J.Pharmacol.Exp.Ther.* **299** 908.
64. Gbahou et al. (2003) *Proc.Natl.Acad.Sci.USA* **100** 11086.
65. Clark and Hill (1996) *Eur.J.Pharmacol.* **296** 223.
66. Rouleau et al. (2002) *Br.J.Pharmacol.* **135** 383.
67. Silver et al. (2002) *Proc.Natl.Acad.Sci.USA* **99** 501.
68. Silver et al. (2001) *Proc.Natl.Acad.Sci.USA* **98** 2855.
69. Leurs et al. (2005) *Nat.Rev.Drug Discov.* **4** 107.
70. Celanire et al. (2005) *Drug Discov.Today.* **10** 1613.
71. Garbarg et al. (1992) *J.Pharmacol.Exp.Ther.* **263** 304.
72. Howson et al. (1992) *Bioorg.Med.Chem.Lett.* **2** 77.
73. Van der Goot et al. (1992) *Eur.J.Med.Chem.* **27** 511.
74. Vollinga et al. (1994) *J.Med.Chem.* **37** 332.
75. Kitbunnadaj et al. (2004) *J.Med.Chem.* **47** 2414.
76. Kitbunnadaj et al. (2005) *J.Med.Chem.* **48** 2100.
77. Mochizuki et al. (1996) *Brain.Res.* **743** 178.
78. Leurs et al. (1995) *Br.J.Pharmacol.* **116** 2315.
79. Ligneau et al. (1994) *J.Pharmacol.Exp.Ther.* **271** 452.
80. Jansen et al. (1994) *Br.J.Pharmacol.* **113** 355.
81. Govoni et al. (2006) *J.Med.Chem.* **49** 2549.
82. Kitbunnadaj et al. (2003) *J.Med.Chem.* **46** 5445.
83. Moreno-Delgado et al. (2006) *Neuropharmacology* **51** 517.
84. Krueger et al. (2005) *J.Pharmacol.Exp.Ther.* **314** 271.
85. Ganellin et al. (1998) *Arch.Pharmacol. (Weinheim)* **331** 395.
86. Schwartz et al. (2000) Patent W00006254.
87. Passani et al. (2004) *TIPS* **25** 618.
88. Hancock et al. (2004) *Exp.Opin.Invest.Drugs* **13** 1237.
89. Barbier et al. (2004) *Br.J.Pharmacol.* **143** 649.
90. Hancock et al. (2005) *Exp.Opin.Invest.Drugs* **14** 223.
91. Yoshimoto et al. (2006) *Proc.Natl.Acad.Sci.USA* **1003** 13866.
92. Woolley et al. (2006) Abstr. BAP Summer Meeting, Oxford, UK TE12
93. Oda et al. (2000) *J.Biol.Chem.* **275** 36781.
94. Liu et al. (2001) *Mol.Pharmacol.* **59** 420.
95. Nguyen et al. (2001) *Mol.Pharmacol.* **59** 427.
96. Morse et al. (2001) *J.Pharmacol.Exp.Ther.* **296** 1058.
97. Zhu et al. (2001) *Mol.Pharmacol.* **59** 434.
98. Coge et al. (2001) *Biochem.Biophys.Res.Comm.* **284** 301.
99. De Esch et al. (2005) *TIPS* **26** 462.
100. Lim et al. (2006) *Curr.Top.Med.Chem.* **6** 1365.
101. Fung-Leung et al. (2004) *Curr.Opin.Invest.Drugs* **5** 1174.
102. Varga et al. (2005) *Eur.J.Pharmacol.* **522** 130.
103. Ikawa et al. (2005) *Biol.Pharm.Bull.* **28** 2016.
104. O'Donoghue et al. (2007) *Dermatol.Ther.* **18** 333.
105. Dunford et al. (2007) *J.Allergy Clin.Immunol.* **119** 176.
106. Cianchi et al. (2005) *Clin.Cancer Res.* **11** 6807.
107. Lim et al. (2006) *J.Med.Chem.* **49** 6650.
108. Terzioglu et al. (2004) *Bioorg.Med.Chem.Lett.* **14** 5251.
109. Venable et al. (2005) *J.Med.Chem.* **48** 8289.
110. Lee-Dutra et al. (2006) *Bioorg.Med.Chem.Lett.* **16** 6043.

Histamine Receptor Compounds Available from Tocris

Catalog No.	Product Name	Primary Function
H₁ Receptors		
Agonists		
0646	HTMT	H ₁ /H ₂ agonist
2478	2-Pyridylethylamine	H ₁ receptor agonist
Antagonists		
2577	Cetirizine	Selective H ₁ antagonist
1453	Clemastine	H ₁ antagonist
1425	(S)-(+)-Dimethindene	H ₁ antagonist. Also M ₂ muscarinic antagonist
0508	Doxepin	Highly potent H ₁ antagonist. Also binds to H ₄ receptor
2429	Fexofenadine	H ₁ receptor antagonist; non-sedating antiallergic agent
0784	Ketotifen	H ₁ antagonist
1944	Loratidine	Peripheral H ₁ antagonist; anti-allergic agent
0660	Mepyramine	Selective H ₁ inverse agonist
2018	Mirtazepine	Potent H ₁ antagonist. Also 5-HT ₂ , 5-HT ₃ and α_2 -antagonist. Antidepressant
0662	<i>trans</i> -Triprolidine	Standard H ₁ antagonist, highly potent
H₂ Receptors		
Agonists		
0668	Amthamine	Highly selective standard H ₂ agonist
0506	Dimaprit	Standard H ₂ selective agonist
0646	HTMT	H ₂ /H ₁ agonist
Antagonists		
0721	Aminopotentidine	H ₂ antagonist
0902	Cimetidine	H ₂ antagonist, I ₁ agonist
0833	ICI 162,846	Potent H ₂ antagonist, active <i>in vivo</i>
1967	Ranitidine	Selective H ₂ antagonist
0826	Tiotidine	Potent, selective H ₂ antagonist
1070	Zolantidine	Potent, centrally active H ₂ antagonist
H₃ and H₄ Receptors		
Agonists		
0729	Imetit	Standard H ₃ and H ₄ agonist (H ₃ > H ₄)
0932	Immepip	Standard H ₃ agonist. Also H ₄ agonist
2315	Immethridine	Potent H ₃ agonist, highly selective over H ₄
0573	N ^α -Methylhistamine	Non-selective H ₃ agonist
0569	(R)-(-)- α -Methylhistamine	Potent, standard H ₃ agonist
0572	(S)-(+)- α -Methylhistamine	H ₃ agonist, less active enantiomer
2342	4-Methylhistamine	Selective, high affinity H ₄ agonist
2477	Proxyfan	High affinity H ₃ ligand
2494	VUF 8430	Potent, high affinity H ₄ agonist
Antagonists		
0752	Clobenpropit	Highly potent H ₃ antagonist/inverse agonist and H ₄ partial agonist
1858	Impentamine	Selective H ₃ antagonist
0779	Iodophenpropit	Potent, selective H ₃ antagonist
2441	JNJ 10191584	Selective H ₄ receptor antagonist; orally active
2477	Proxyfan	High affinity H ₃ ligand
2034	ROS 234	Potent H ₃ antagonist
0644	Thioperamide	H ₃ /H ₄ antagonist/inverse agonist. Active <i>in vivo</i>
2493	VUF 2681	Potent H ₃ receptor silent antagonist

R&D SYSTEMS

NOVUS
BIOLOGICALS

TOCRIS

protein simple

ACD™

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

