

# 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.<sup>1</sup> 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).<sup>2</sup> 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.<sup>3</sup> 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 H<sub>1</sub> and H<sub>2</sub> receptor subtypes have proven to be excellent drug targets. Ligands for the histamine H<sub>3</sub> receptor subtype are currently entering clinical studies and the recently discovered histamine H<sub>4</sub> receptor subtype is subject of intense preclinical research.

## The Histamine H<sub>1</sub> 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.<sup>4</sup> However, the first generation of H<sub>1</sub> 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 agonists<sup>5</sup>) are clinically very successful and are widely available drugs. In 1991, the cDNA encoding a bovine H<sub>1</sub> receptor protein was cloned after an expression cloning strategy in *Xenopus oocytes*.<sup>6</sup> The deduced amino acid sequence revealed a 491 amino acid protein of 56 kDa. Using the cDNA sequence encoding the bovine H<sub>1</sub> receptor, the cDNA sequences and intronless genes encoding the rat,<sup>3,7</sup> guinea-pig,<sup>8,9</sup> human<sup>10,11</sup> and mouse<sup>12</sup> H<sub>1</sub> receptor proteins were cloned soon thereafter.

### Contents

Introduction.....	1
The Histamine H <sub>1</sub> Receptor .....	1
Ligands for H <sub>1</sub> Receptors .....	2
The Histamine H <sub>2</sub> Receptor .....	3
Ligands for H <sub>2</sub> Receptors .....	3
The Histamine H <sub>3</sub> Receptor .....	4
Ligands for H <sub>3</sub> Receptors .....	4
The Histamine H <sub>4</sub> Receptor .....	6
Ligands for H <sub>4</sub> 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<sup>7,8,10</sup> resulted in the identification of several DNA-binding motifs, including potential glucocorticoid responsive elements. The human H<sub>1</sub> receptor gene resides on chromosome 3.<sup>13</sup> The H<sub>1</sub> receptor belongs to the large family of GPCRs. The receptor is associated with the phospholipase C-catalysed formation of inositol 1,4,5-triphosphate (IP<sub>3</sub>) and 1,2-diacylglycerol (DAG). Histamine induces production of inositol phosphates in several tissues (including brain, airway, intestinal and vascular smooth muscle<sup>24</sup>) via G<sub>αq</sub> protein activation.<sup>14</sup> In other tissues, activation of H<sub>1</sub> 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 Ca<sup>2+</sup>.<sup>15</sup> In any case, alternative signalling pathways can be mediated by the histamine H<sub>1</sub> receptor. Recent results indicate that the functional heterogeneity can be ligand-directed.<sup>16</sup>

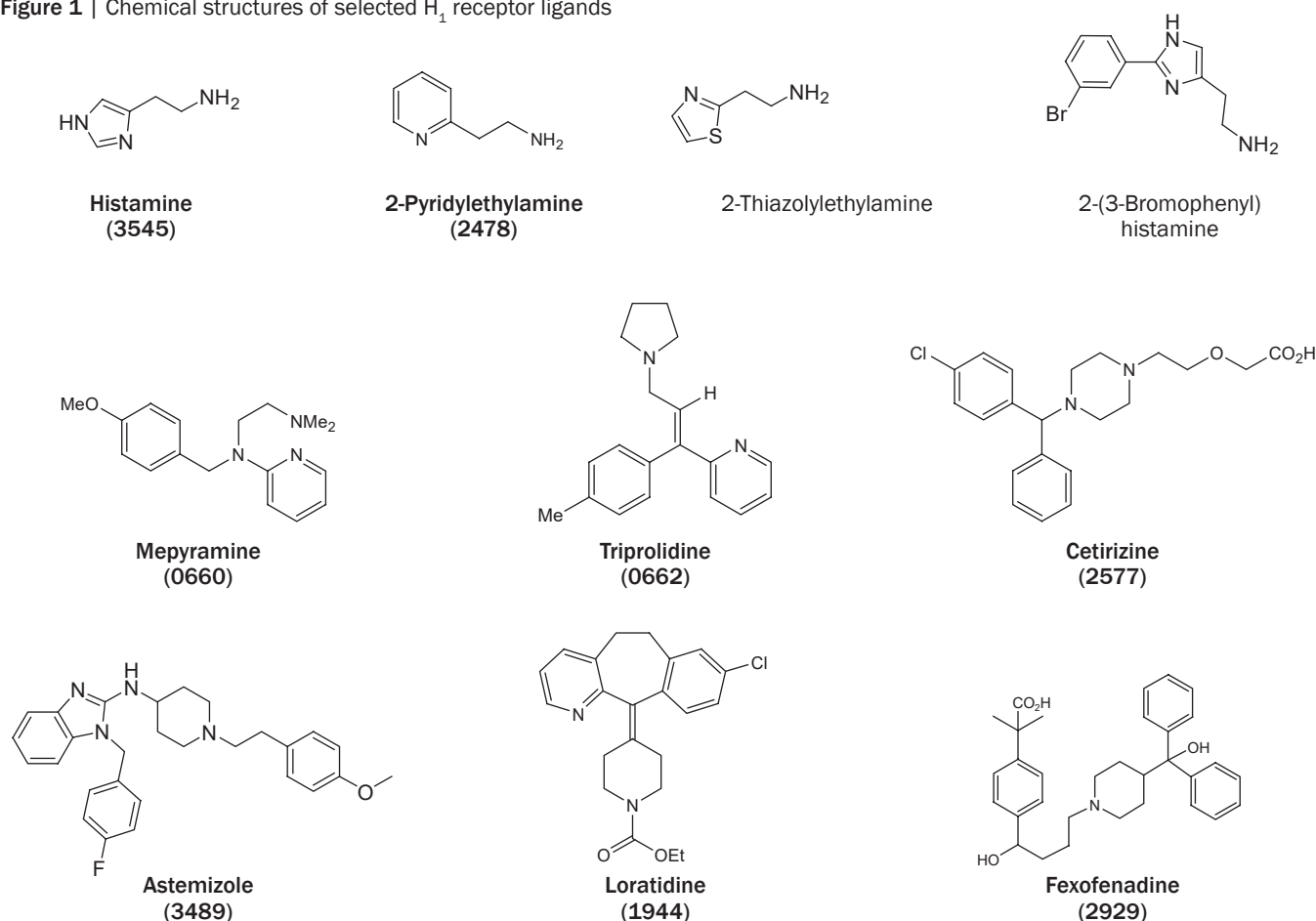
The histamine H<sub>1</sub> receptor is a well established drug target and has been thoroughly studied for decades. Nevertheless, H<sub>1</sub> receptor research continues to flourish as many new techniques and approaches are being developed by using this receptor as an archetypal GPCR target. Most notably, in the last few years, the histamine H<sub>1</sub> receptor has been the subject of various molecular biology studies (e.g., large-scale overproduction<sup>17</sup> and GPCR

binding site elucidation<sup>18</sup>), biophysical approaches (such as solid-state NMR<sup>19</sup>) and investigations towards the general activation mechanisms of GPCRs.<sup>20,21</sup>

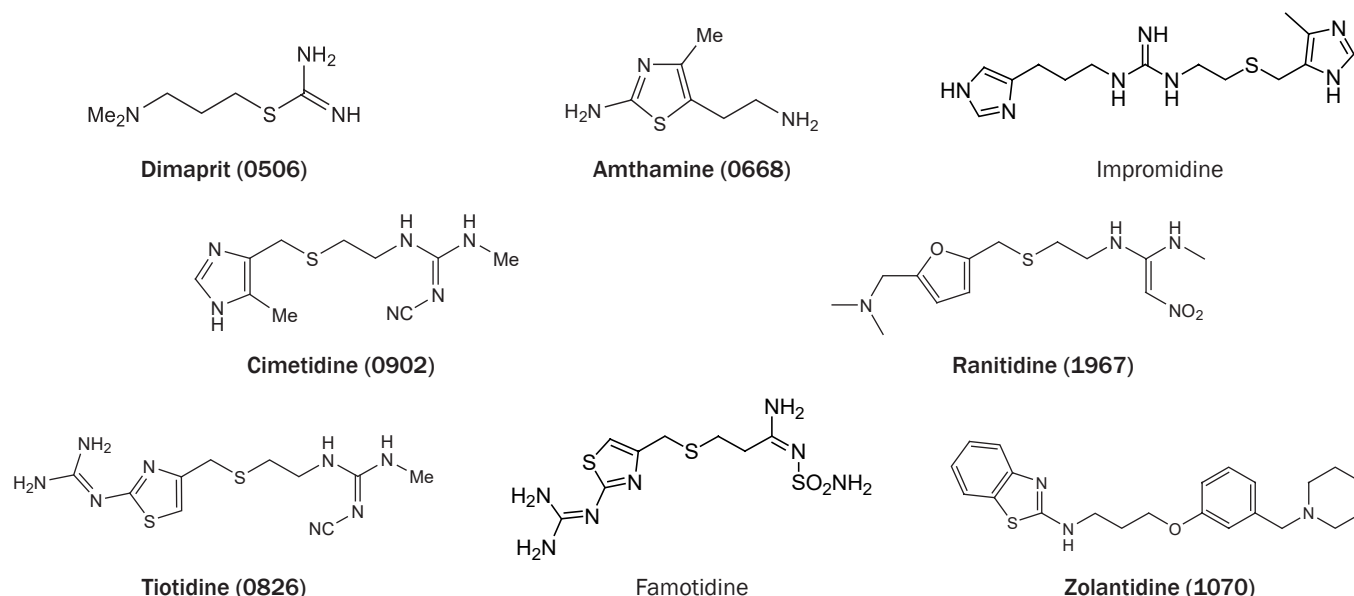
## Ligands for H<sub>1</sub> Receptors

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

**Figure 1** | Chemical structures of selected H<sub>1</sub> receptor ligands



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

**Figure 2** | Chemical structures of selected H<sub>2</sub> receptor ligands

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

## The Histamine H<sub>2</sub> Receptor

The observation that the classical ‘antihistamines’ (i.e. H<sub>1</sub> 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: H<sub>1</sub> and H<sub>2</sub>.<sup>25</sup> This hypothesis became generally accepted when Black *et al*<sup>26</sup> 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 H<sub>2</sub> receptor antagonists proved to be very useful in the therapy of gastric ulcers. Gantz *et al*<sup>27</sup> were the first to clone a cDNA encoding a 359 amino acid H<sub>2</sub> receptor. Using degenerate primers based on the known sequence similarity of various GPCRs, the H<sub>2</sub> receptor sequence was obtained from canine gastric parietal cDNA by PCR. Soon thereafter, the intronless genes encoding the rat,<sup>28</sup> human,<sup>29</sup> guinea pig<sup>30</sup> and mouse<sup>31</sup> H<sub>2</sub> receptor were cloned by means of homology screening. Identification of the promoter region of the human H<sub>2</sub> receptor gene revealed the existence of regulatory transcription sites and regions displaying stimulatory and inhibitory effects on gene expression monitored in a luciferase assay.<sup>32</sup> Studies have indicated that the human H<sub>2</sub> receptor gene resides on chromosome 5.<sup>30</sup> Interestingly, several polymorphisms have been found in the human H<sub>2</sub> receptor gene<sup>33</sup> and one of the mutations has been linked to schizophrenia.<sup>34</sup> The histamine H<sub>2</sub> receptor is coupled to the adenylate cyclase system in a variety of tissues (e.g. brain, stomach, heart, gastric mucosa, lung).<sup>24</sup> Moreover, cell lines transfected with the cloned H<sub>2</sub> receptor genes showed an H<sub>2</sub> receptor-mediated increase of cAMP.<sup>35,36,37</sup> In addition, alternative signalling pathways for the H<sub>2</sub> receptor have been identified. In differentiated HL-60 cells and CHO or HEPA cells transfected with the H<sub>2</sub> receptor cDNA, an H<sub>2</sub> receptor-mediated increase of the intracellular Ca<sup>2+</sup> concentration and/or IP<sub>3</sub> levels

was observed.<sup>38,39</sup> Moreover, in CHO cells expressing the rat H<sub>2</sub> receptor, activation of the H<sub>2</sub> receptor resulted in an inhibition of the release of arachidonic acid induced by either constitutive purinergic receptors or a Ca<sup>2+</sup>-ionophore,<sup>36</sup> as well as an increase in cAMP. These new signal transduction pathways are both regulated via unknown, cAMP-independent pathways.

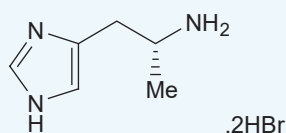
## Ligands for H<sub>2</sub> Receptors

The first selective H<sub>2</sub> receptor agonist was dimaprit (figure 2). This compound is almost as active as histamine at the H<sub>2</sub> receptor but hardly displays any H<sub>1</sub> receptor activity.<sup>40</sup> Much later, it was shown that dimaprit is also a moderate H<sub>3</sub> receptor antagonist<sup>41</sup> and a moderate H<sub>4</sub> receptor agonist.<sup>42</sup> Amthamine can be considered a rigid dimaprit analogue.<sup>43</sup> This compound combines a high H<sub>2</sub> receptor selectivity with a potency which is slightly higher compared to histamine, both *in vitro* and *in vivo*.<sup>44</sup> An H<sub>2</sub> receptor agonist that is more potent than histamine is the guanidine derivative impromidine. This ligand actually combines a rather high H<sub>2</sub> receptor affinity with a reduced efficacy. Impromidine also shows moderate H<sub>1</sub>- and potent H<sub>3</sub>-receptor antagonistic activity<sup>45,46</sup> as well as potent H<sub>4</sub> receptor partial agonistic activity.<sup>42</sup>

The finding that N $\alpha$ -guanylhistamine acts as a partial H<sub>2</sub> agonist in a gastric acid secretion test did lead to the development of the relatively weak H<sub>2</sub> antagonist burimamide. Years later, it was shown that burimamide is also an H<sub>3</sub> and H<sub>4</sub> receptor partial agonist.<sup>42</sup> Nevertheless, burimamide was a good lead for the development of selective and clinically useful H<sub>2</sub> receptor antagonists, such as cimetidine.<sup>45,47,48</sup> 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 H<sub>2</sub> receptor than

**(R)-(-)- $\alpha$ -Methylhistamine, High Affinity  $H_3$  Agonist****(R)-(-)- $\alpha$ -Methylhistamine**

Cat. No. 0569



This potent and high affinity  $H_3$  agonist displays >200-fold selectivity over  $H_4$  receptors. The compound inhibits  $H_3$  receptor-mediated histamine synthesis and release in the CNS and stimulates  $H_4$  receptor-mediated eosinophil shape change ( $EC_{50} = 66$  nM).

**Oishi et al.** (1989) Effects of histamine  $H_3$ -agonist (R)- $\alpha$ -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-  $H_3$  receptors in guinea pig ileum with  $H_3$ -selective ligands. *Br.J.Pharmacol.* **101** 621. **Schwartz et al.** (1990) A third histamine receptor subtype - characterization, localization and functions of the  $H_3$ -receptor. *Agents Actions* **30** 13. **Buckland et al.** (2003) Histamine induces cytoskeletal changes in human eosinophils via the  $H_4$  receptor. *Br.J.Pharmacol.* **140** 1117.

cimetidine. Moreover, the replacement of the imidazole moiety also eliminates the undesired inhibition of cytochrome P450.<sup>48</sup> The potent tritiated  $H_2$  antagonist tiotidine and [<sup>125</sup>I]-iodinated  $H_2$  antagonist iodoaminopotentidine are successfully used as radioligands for the  $H_2$  receptor.<sup>24</sup> The newly developed brain-penetrating  $H_2$  antagonist zolantidine is an important tool for *in vivo* CNS studies.<sup>49</sup> The  $H_2$  receptor was reported to be spontaneously active in transfected CHO cells.<sup>50</sup> Based on this concept, many  $H_2$  antagonists were reclassified; cimetidine, ranitidine and famotidine are in fact inverse agonists, whereas burimamide acts in this model system as a neutral antagonist.<sup>50</sup>

## The Histamine $H_3$ 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  $H_1$  and  $H_2$  receptor subtypes as no correlation with either the  $H_1$  or the  $H_2$  receptor activity of known histaminergic compounds was observed.<sup>46</sup> Soon after, the  $H_3$  receptor agonist (R)-(-)- $\alpha$ -methylhistamine and the antagonist thioperamide (see figure 3) were developed.<sup>51</sup> 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.<sup>52</sup> 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  $H_3$  receptor is present in the peripheral nervous system, e.g. in the gastrointestinal tract, the airways and the cardiovascular system.<sup>53,54</sup> Initial efforts to identify the  $H_3$  receptor gene, using the anticipated homology with the identified  $H_1$  and  $H_2$  receptor gene all failed. Eventually, the human  $H_3$  receptor cDNA was cloned by Lovenberg and co-workers in 1999.<sup>55</sup> In search of novel GPCRs through homology searching of expressed sequence tag databases, a receptor with high similarity to the  $M_2$  muscarinic acetylcholine receptor was identified. Expression of the gene and full characterisation established this protein as the histamine  $H_3$  receptor. The cloning of the  $H_3$  receptor of other species, including

rat,<sup>56</sup> guinea pig<sup>57</sup> and mouse<sup>58</sup> soon followed and it was revealed that major  $H_3$  receptor species differences exist. The human  $H_3$  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.<sup>59</sup> The  $H_3$  receptor gene can undergo extensive alternative splicing, resulting in many  $H_3$  receptor isoforms that have different signalling properties and expression profiles.<sup>59,60,61</sup> It was shown that the  $H_3$  receptor displays particularly high constitutive activity, again leading to a reclassification of existing ligands into agonists, neutral antagonists and inverse agonists.<sup>62,63</sup> 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.<sup>64</sup>

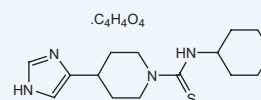
The  $H_3$  receptor signals through  $G_{i/o}$  proteins as was shown by the pertussis toxin sensitive stimulation of [<sup>35</sup>S]-GTP $\gamma$ S binding in rat cortical membranes.<sup>65</sup> Through negative coupling to adenylyl cyclase, stimulation of the  $H_3$  receptor results in lower levels of cAMP, thereby reducing downstream signalling events such as CREB-dependent gene transcription.<sup>56</sup> Alternative signalling pathways may be activated by the  $G_{i/o}$  proteins, including mitogen-activated protein kinase (MAPK)<sup>60</sup> and phosphatidylinositol 3-kinase (PI3K) pathways.  $G_{i/o}$  protein activation can also lead to the activation of phospholipase  $A_2$  ( $PLA_2$ ) to induce the release of arachidonic acid, the lowering of intracellular  $Ca^{2+}$  levels through voltage-gated ion channels<sup>67</sup> and the inhibition of the  $Na^+/H^+$  exchanger (NHE).<sup>68</sup> With the recent progress that has been made in the characterisation of the  $H_3$  receptor (as outlined above), many pharmaceutical companies have very active  $H_3$  receptor drug development programs. As a result, several  $H_3$  receptor ligands have entered clinical studies for a plethora of applications, including obesity, narcolepsy, dementia and migraine (among others).<sup>69,70</sup>

## Ligands for $H_3$ Receptors

At the  $H_3$  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  $H_3$  and  $H_4$  Ligand****Thioperamide**

Cat. No. 0644



Thioperamide, the potent histamine  $H_3$  and  $H_4$  ligand, displays  $K_i$  values of 25 and 27 nM at recombinant  $H_3$  and  $H_4$  receptors respectively. The compound acts as an antagonist at  $H_3$  receptors and displays inverse agonist activity at  $H_4$  receptors. It freely crosses the blood-brain barrier.

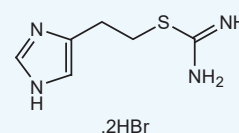
**Hew et al.** (1990) Characterisation of histamine  $H_3$  receptors in guinea pig ileum with  $H_3$  selective ligands. *Br.J.Pharmacol.* **101** 621. **Liu et al.** (2001) Cloning and pharmacological characterization of a fourth histamine receptor ( $H_4$ ) expressed in bone marrow. *Mol. Pharmacol.* **59** 420. **Ling et al.** (2004) Histamine  $H_4$  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  $H_3$  and  $H_4$  receptors: structure-activity relationships of histamine derivatives. *Br.J.Pharmacol.* **147** 744.

and selective for the H<sub>3</sub> receptor. Methylation of the α-carbon atom of the ethylamine sidechain drastically increases the potency at the H<sub>3</sub> 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 H<sub>1</sub> and H<sub>2</sub> receptors, (R)-(-)-α-methylhistamine (figure 3) was initially considered a selective agonist at the H<sub>3</sub> receptor. However, when the H<sub>4</sub> receptor was discovered it was shown that (R)-(-)-α-methylhistamine also has considerable affinity for this histamine receptor subtype.<sup>42</sup> Nevertheless, in combination with its less active S-isomer, (R)-(-)-α-methylhistamine has proven to be highly useful for the pharmacological characterisation of H<sub>3</sub> receptor-mediated effects.<sup>41</sup> Tritiated forms of α-methylhistamine and (R)-(-)-α-methylhistamine are available as radiolabelled agonists for the H<sub>3</sub> receptor. For potent H<sub>3</sub> 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*,<sup>71,72,73</sup> as is (R)-(-)-α-methylhistamine. The amine function can also be incorporated in ring structures. For example, immepip is a potent H<sub>3</sub> receptor agonist that is effective *in vitro* and *in vivo*.<sup>74</sup> Although the described first generation H<sub>3</sub> agonists were intensively used as reference ligands to study the H<sub>3</sub> receptor, all of them proved to also have considerable activity at the H<sub>4</sub> receptor. Therefore, a new generation of potent and selective H<sub>3</sub> receptor agonists has been developed, most notably immethridine<sup>75</sup> (pEC<sub>50</sub> = 9.74; displays 300-fold selectivity over the H<sub>4</sub> receptor) and

### Imetit, High Affinity H<sub>3</sub>/H<sub>4</sub> Agonist

Imetit

Cat. No. 0729



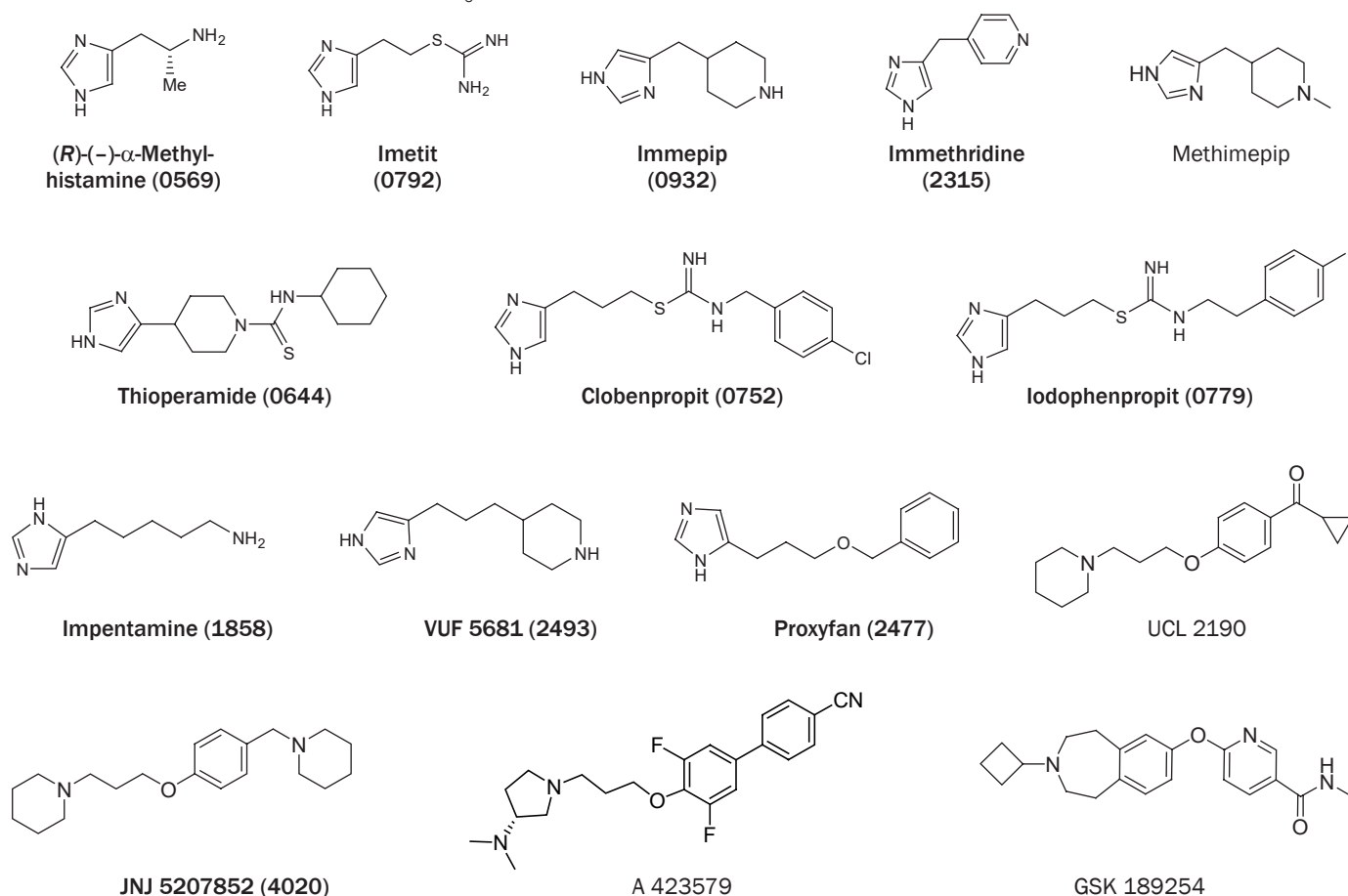
Imetit is an extremely high affinity, potent agonist at H<sub>3</sub> and H<sub>4</sub> receptors (K<sub>i</sub> values are 0.3 and 2.7 nM respectively). The agonist induces shape change in eosinophils *in vitro* with an EC<sub>50</sub> of 25 nM and is centrally active following systemic administration *in vivo*.

Garbarg *et al.* (1992) S-[2-(4-Imidazolyl)ethyl]isothioureia, a highly specific and potent histamine H<sub>3</sub> 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 H<sub>4</sub> receptor mediates eosinophil chemotaxis with cell shape change and adhesion molecule upregulation. *Br.J.Pharmacol.* **142** 161.

methimepip.<sup>76</sup> These latter compounds are devoid of high H<sub>4</sub> receptor activity.

As with the first generation H<sub>3</sub> receptor agonists, the first generation H<sub>3</sub> receptor antagonists (all of them possessing an imidazole heterocycle) turned out to have affinity for the H<sub>4</sub> receptor. The first potent H<sub>3</sub> receptor antagonist (later reclassified as an inverse agonist) that was devoid of H<sub>1</sub> receptor and H<sub>2</sub>

**Figure 3** | Chemical structures of selected H<sub>3</sub> receptor ligands



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



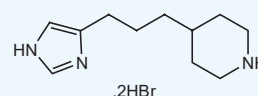
receptor activity was thioperamide (figure 3).<sup>51</sup> This compound has been used in many H<sub>3</sub> receptor studies as a reference ligand and is active *in vitro* and *in vivo* (the compound is able to penetrate the CNS).<sup>77</sup> However, thioperamide displays some 5-HT<sub>3</sub> receptor antagonism<sup>78</sup> and is an inverse agonist at the H<sub>4</sub> receptor. The remarkable H<sub>3</sub> receptor species difference can be demonstrated with thioperamide as the compound has a 10-fold higher affinity for the rat H<sub>3</sub> receptor than for the human H<sub>3</sub> receptor.<sup>56</sup> Based on the H<sub>3</sub> receptor agonist imetit (*vide ante*), the highly potent H<sub>3</sub> receptor inverse agonist clobenpropit was developed (pA<sub>2</sub> = 9.9).<sup>73</sup> This compound has some 5-HT<sub>3</sub> receptor activity<sup>78</sup> and displays partial agonist activity at H<sub>4</sub> receptors. In addition, radioligands for the H<sub>3</sub> receptor have been described (e.g. [<sup>125</sup>I]-iodophenpropit and [<sup>125</sup>I]-iodoproxyfan).<sup>79,80</sup> The moderately active H<sub>2</sub> receptor antagonist burimamide (pA<sub>2</sub> = 5.1, figure 2) also has good affinity for the H<sub>3</sub> (pK<sub>i</sub> = 7.9) and the H<sub>4</sub> (pK<sub>i</sub> = 7.4) receptor.<sup>42</sup> Impentamine is a potent histamine H<sub>3</sub> receptor inverse agonist (pA<sub>2</sub> = 8.4). Like burimamide, this compound can act as a partial agonist in SK-N-MC cells expressing human H<sub>3</sub> 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.<sup>63,81</sup> The compound VUF 5681 was reported as a neutral H<sub>3</sub> antagonist (or 'silent' antagonist),<sup>82</sup> findings that were recently confirmed in studies investigating constitutive activity using rat brain cortex.<sup>83</sup> 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*.<sup>64,84</sup>

The first non-imidazole H<sub>3</sub> receptor ligand was reported by Ganellin in 1998.<sup>85</sup> This elegant medicinal chemistry work did lead to the potent compound UCL 2190.<sup>86</sup> Following the cloning of the H<sub>3</sub> receptor in 1999, several pharmaceutical companies entered the H<sub>3</sub> 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 H<sub>3</sub> receptor ligands, e.g. piperidiny propyloxy side chain. This neutral antagonist is active in several models for cognition<sup>87,88</sup> but does

### VUF 5681, a Novel H<sub>3</sub> Antagonist

VUF 5681

Cat. No. 2493



VUF 5681 is a potent histamine H<sub>3</sub> receptor silent antagonist (pK<sub>i</sub> = 8.35).

**Kitbunnadaj et al.** (2004) Identification of 4-(1H-imidazol-4(5)-ylmethyl)pyridine (immethridine) as a novel, potent, and highly selective histamine H<sub>3</sub> receptor agonist. *J. Med. Chem.* **47** 2414. **Leurs et al.** (2005) The histamine H<sub>3</sub> receptor: from gene cloning to H<sub>3</sub> receptor drugs. *Nat. Rev. Drug Discov.* **4** 107. **Moreno-Delgado et al.** (2006) Constitutive activity of H<sub>3</sub> 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.<sup>89</sup> Other compounds, such as Abbott's A 423579 seem to have more efficacy in obesity models, while lacking clear cognitive effects.<sup>90</sup> At the time of writing the differences in efficacy for distinct clinical applications of the different classes of H<sub>3</sub> ligands is not understood and subject of intense research.<sup>70,91</sup> Interestingly, GSK 189254 has been in trials for three different diseases: neuropathic pain, narcolepsy and dementia.<sup>92</sup>

### The Histamine H<sub>4</sub> Receptor

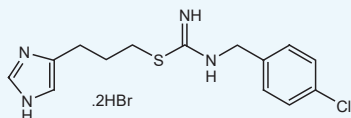
Immediately following the cloning of the H<sub>3</sub> receptor gene, several groups identified the homologous H<sub>4</sub> receptor sequence in the human genome databases.<sup>93–97</sup> Indeed, the H<sub>4</sub> receptor has high sequence identity with the H<sub>3</sub> receptor (31% at the protein level, 54% in the transmembrane domains). The H<sub>3</sub> and H<sub>4</sub> receptor are also similar in gene structure. The human H<sub>4</sub> 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.<sup>98</sup> To date, two H<sub>4</sub> receptor isoforms have been identified. Cloning of the genes that encode the mouse, rat, guinea-pig and pig H<sub>4</sub> receptors reveal only limited sequence homology with the human H<sub>4</sub> receptor. The H<sub>4</sub> receptor is mainly expressed in bone marrow and peripheral leukocytes, and mRNAs of the human H<sub>4</sub> receptor are detected in e.g. mast cells, dendritic cells, spleen and eosinophils.<sup>93–97</sup> The receptor has a pronounced effect on the chemotaxis of several cell types that are associated with immune and inflammatory responses. The H<sub>4</sub> receptor is mainly coupled to G<sub>i/o</sub> 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, H<sub>4</sub> receptor stimulation affects the pertussis-toxin-sensitive activation of mitogen-activated protein (MAP) kinase pathways. Studying the increased levels of [<sup>35</sup>S] GTPγS levels in H<sub>4</sub> transfected cells, it has been shown that the H<sub>4</sub> receptor is constitutively active. The Gβγ subunits of the G<sub>i/o</sub> proteins activate phospholipase C, and thereby increase the Ca<sup>2+</sup> concentrations. In mast cell and eosinophils, this Ca<sup>2+</sup> response can be linked to cellular chemotaxis.

Considering the physiological role of the H<sub>4</sub> receptor, several applications are under preclinical investigation,<sup>99,100</sup> including allergy and asthma,<sup>101</sup> as well as chronic inflammations such as

### Clobenpropit, Highly Potent H<sub>3</sub> Antagonist

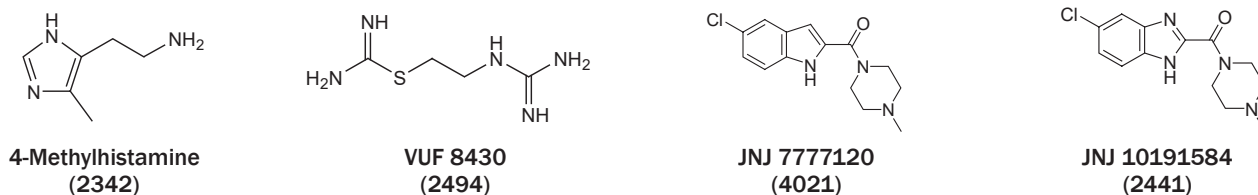
Clobenpropit

Cat. No. 0752



An extremely potent antagonist/inverse agonist at H<sub>3</sub> receptors (pA<sub>2</sub> = 9.93), clobenpropit also displays partial agonist activity at H<sub>4</sub> receptors. The ligand induces eosinophil shape change with an EC<sub>50</sub> of 3 nM.

**Van der Goot et al.** (1992) Isothiourea analogues of histamine as potent agonists or antagonists of the histamine H<sub>3</sub> receptor. *Eur. J. Med. Chem.* **27** 511. **Yokoyama et al.** (1994) Clobenpropit (VUF-9153), a new histamine H<sub>3</sub> 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 (H<sub>4</sub>) expressed in bone marrow. *Mol. Pharmacol.* **59** 420. **Buckland et al.** (2003) Histamine induces cytoskeletal changes in human eosinophils via the H<sub>4</sub> receptor. *Br. J. Pharmacol.* **140** 1117.

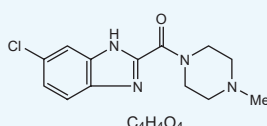
**Figure 4** | The first selective H<sub>4</sub> receptor ligands reported in scientific literature

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

### **JNJ 10191584, H<sub>4</sub>-Selective Antagonist**

**JNJ 10191584**

**Cat. No. 2441**



JNJ 10191584 is a highly selective histamine H<sub>4</sub> receptor silent antagonist. It binds with high affinity to the human H<sub>4</sub> receptor ( $K_i = 26$  nM) and is > 540-fold selective over the H<sub>3</sub> receptor ( $K_i = 14.1$  μM). *In vitro* the antagonist inhibits mast cell and eosinophil chemotaxis with IC<sub>50</sub> 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 H<sub>4</sub> 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 H<sub>4</sub> antagonists. *J. Med. Chem.* **48** 8289. Varga *et al.* (2005) Inhibitory effects of histamine H<sub>4</sub> receptor antagonists on experimental colitis in the rat. *Eur. J. Pharmacol.* **522** 130.

inflammatory bowel disease (IBD)<sup>102</sup> and rheumatoid arthritis.<sup>103</sup> The H<sub>4</sub> receptor is also being associated with pruritus (itch)<sup>104,105</sup> and is involved in the progression of colon cancer.<sup>106</sup>

### Ligands for H<sub>4</sub> Receptors

As was described above, most imidazole-containing, first generation H<sub>3</sub> receptor ligands have considerable affinity for the H<sub>4</sub> receptor as well. However, selective H<sub>4</sub> receptor ligands have been described. 4-Methylhistamine is a potent H<sub>4</sub> agonist while displaying more than a 100-fold selectivity over the other histamine receptor subtypes, including the H<sub>2</sub> receptor for which this ligand was originally developed.<sup>42</sup> A slightly different and complimentary profile was reported for VUF 8430. This compound has a high H<sub>4</sub> receptor activity and affinity, minimal affinity for the H<sub>1</sub> and H<sub>2</sub> receptor and a 33-fold lower affinity for the H<sub>3</sub> receptor.<sup>107</sup>

Potent and selective H<sub>4</sub> receptor antagonists are also emerging. The first reported neutral antagonist is JNJ 777120, a compound that is frequently being used as H<sub>4</sub> receptor reference ligand. Currently, [<sup>3</sup>H]-JNJ 777120 and [<sup>3</sup>H]-histamine are used as H<sub>4</sub> receptor radioligands. Unfortunately, JNJ 777120 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 H<sub>4</sub> antagonist.<sup>108</sup> This compound is orally active *in vivo* and has an improved liver microsome stability but still a limited half life.<sup>102,109</sup> More recently, 2-arylbenzimidazoles

have been described as ligands with low nanomolar affinity for the H<sub>4</sub> receptor.<sup>110</sup> Considering the number of H<sub>4</sub> receptor-related patent applications that have recently been disclosed (as reviewed elsewhere<sup>100</sup>), it can be anticipated that many new H<sub>4</sub> 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 H<sub>1</sub> and H<sub>2</sub> receptors in 1966, discovery of the H<sub>3</sub> receptor in 1983 and the H<sub>4</sub> 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 (H<sub>1</sub> and H<sub>2</sub>) addressed by blockbuster drugs, the third subtype (H<sub>3</sub>) leading to frantic clinical studies and the latest addition to the family (H<sub>4</sub>) 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

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



## Histamine Receptor Compounds Available from Tocris

Catalog No.	Product Name	Primary Function
<b>H<sub>1</sub> Receptors</b>		
<b>Agonists</b>		
2478	<a href="#">2-Pyridylethylamine</a>	H <sub>1</sub> agonist
<b>Inverse Agonists</b>		
0660	<a href="#">Mepyramine</a>	Selective H <sub>1</sub> inverse agonist
<b>Antagonists</b>		
3489	<a href="#">Astemizole</a>	Orally active, potent H <sub>1</sub> antagonist; also K <sub>v</sub> 11.1 (hERG) channel blocker
2577	<a href="#">Cetirizine</a>	Selective H <sub>1</sub> antagonist
1453	<a href="#">Clemastine</a>	H <sub>1</sub> antagonist
5371	<a href="#">Clemizole</a>	H <sub>1</sub> antagonist; also TRPC5 blocker
5958	<a href="#">Desloratadine</a>	High affinity H <sub>1</sub> antagonist; anti-inflammatory
1425	<a href="#">(S)-(+)-Dimethindene</a>	H <sub>1</sub> antagonist; also M <sub>2</sub> muscarinic antagonist
3072	<a href="#">Diphenhydramine</a>	H <sub>1</sub> antagonist
0508	<a href="#">Doxepin</a>	Highly potent H <sub>1</sub> antagonist; also binds to H <sub>4</sub> receptor
2429	<a href="#">Fexofenadine</a>	H <sub>1</sub> receptor antagonist; non-sedating antiallergic agent
0784	<a href="#">Ketotifen</a>	H <sub>1</sub> antagonist
1944	<a href="#">Loratadine</a>	Peripheral H <sub>1</sub> antagonist; anti-allergic agent
4245	<a href="#">Meclizine</a>	H <sub>1</sub> antagonist; also human pregnane X receptor agonist
2018	<a href="#">Mirtazepine</a>	Potent H <sub>1</sub> antagonist; also 5-HT <sub>2</sub> , 5-HT <sub>3</sub> and $\alpha_2$ -antagonist; antidepressant
4241	<a href="#">Olopatadine</a>	H <sub>1</sub> antagonist; ocular anti-allergic agent
3948	<a href="#">Terfenadine</a>	H <sub>1</sub> antagonist; also K <sub>v</sub> 11.1 (hERG) and K <sub>v</sub> 6 (KATP) channel blocker
0662	<a href="#">trans-Tripolidine</a>	Highly potent H <sub>1</sub> antagonist
3996	<a href="#">Zotepine</a>	H <sub>1</sub> antagonist; also 5-HT <sub>2A</sub> and D <sub>2</sub> antagonist
<b>H<sub>2</sub> Receptors</b>		
<b>Agonists</b>		
0668	<a href="#">Amthamine</a>	Highly selective standard H <sub>2</sub> agonist
0506	<a href="#">Dimaprit</a>	Standard H <sub>2</sub> selective agonist
<b>Antagonists</b>		
0721	<a href="#">Aminopotentidine</a>	H <sub>2</sub> antagonist
0902	<a href="#">Cimetidine</a>	H <sub>2</sub> antagonist; also I <sub>1</sub> agonist
0833	<a href="#">ICI 162,846</a>	Potent H <sub>2</sub> antagonist; active <i>in vivo</i>
1967	<a href="#">Ranitidine</a>	Selective H <sub>2</sub> antagonist
0826	<a href="#">Tiotidine</a>	Potent and selective H <sub>2</sub> antagonist
1070	<a href="#">Zolantidine</a>	Potent, centrally active H <sub>2</sub> antagonist
<b>H<sub>3</sub> Receptors</b>		
<b>Agonists</b>		
2315	<a href="#">Immethridine</a>	Potent H <sub>3</sub> agonist; highly selective over H <sub>4</sub>
0573	<a href="#">N<sup>α</sup>-Methylhistamine</a>	Non-selective H <sub>3</sub> agonist
0569	<a href="#">(R)-(-)-<math>\alpha</math>-Methylhistamine</a>	Potent standard H <sub>3</sub> agonist
0572	<a href="#">(S)-(+)-<math>\alpha</math>-Methylhistamine</a>	H <sub>3</sub> agonist; less active enantiomer
2477	<a href="#">Proxyfan</a>	High affinity H <sub>3</sub> ligand
<b>Inverse Agonists</b>		
3743	<a href="#">BF 2649</a>	Potent and selective H <sub>3</sub> inverse agonist
<b>Antagonists</b>		
4697	<a href="#">A 331440</a>	Selective H <sub>3</sub> antagonist
2211	<a href="#">Carcinine</a>	Highly selective H <sub>3</sub> antagonist
2419	<a href="#">GT 2016</a>	Selective H <sub>3</sub> antagonist
1858	<a href="#">Impentamine</a>	Selective H <sub>3</sub> antagonist
0779	<a href="#">Iodophenpropit</a>	Potent and selective H <sub>3</sub> antagonist

Catalog No.	Product Name	Primary Function
4019	<a href="#">JNJ 10181457</a>	H <sub>3</sub> antagonist
4020	<a href="#">JNJ 5207852</a>	High affinity H <sub>3</sub> antagonist
2034	<a href="#">ROS 234</a>	Potent H <sub>3</sub> antagonist
4441	<a href="#">SEN 12333</a>	H <sub>3</sub> antagonist; also $\alpha 7$ nAChR agonist
2493	<a href="#">VUF 5681</a>	Potent H <sub>3</sub> receptor silent antagonist
<b>H<sub>4</sub> Receptors</b>		
<b>Agonists</b>		
2342	<a href="#">4-Methylhistamine</a>	Selective, high affinity H <sub>4</sub> agonist
4769	<a href="#">VUF 10460</a>	Selective H <sub>4</sub> agonist
2494	<a href="#">VUF 8430</a>	Potent, high affinity H <sub>4</sub> agonist
<b>Antagonists</b>		
3753	<a href="#">A 943931</a>	Potent and selective H <sub>4</sub> antagonist
3640	<a href="#">A 987306</a>	Potent and selective H <sub>4</sub> antagonist
2441	<a href="#">JNJ 10191584</a>	Selective H <sub>4</sub> antagonist; orally active
6279	<a href="#">JNJ 39758979</a>	High affinity and selective H <sub>4</sub> antagonist; orally bioavailable
4021	<a href="#">JNJ 7777120</a>	Selective H <sub>4</sub> antagonist
<b>Non-selective Histamine Compounds</b>		
<b>Agonists</b>		
3545	<a href="#">Histamine</a>	Endogenous histamine agonist
0646	<a href="#">HTMT</a>	H <sub>1</sub> and H <sub>2</sub> agonist
0729	<a href="#">Imetit</a>	Standard H <sub>3</sub> and H <sub>4</sub> agonist (H <sub>3</sub> > H <sub>4</sub> )
0932	<a href="#">Immepip</a>	H <sub>3</sub> and H <sub>4</sub> agonist
<b>Antagonists</b>		
0752	<a href="#">Clobenpropit</a>	Highly potent H <sub>3</sub> antagonist; also H <sub>4</sub> partial agonist
0644	<a href="#">Thioperamide</a>	H <sub>3</sub> antagonist; also H <sub>4</sub> inverse agonist
<b>Histaminergic-related Compounds</b>		
4857	<a href="#">Amlexanox</a>	Inhibits histamine release from rat mast cells; anti-allergic agent
3201	<a href="#">Dimebon</a>	Non-selective antihistamine; displays neuroprotectant and cognitive enhancing abilities
0743	<a href="#">DPPE</a>	Inhibitor of histamine binding at the intracellular binding site
3848	<a href="#">Sinomenine</a>	Anti-inflammatory; causes mast cell degranulation and histamine release
0512	<a href="#">SKF 91488</a>	Histamine N-methyltransferase inhibitor
6424	<a href="#">Sodium Cromoglicate</a>	Inhibits histamine release and mast cell degranulation

## Life Science Literature from Tocris

Tocris provides a wide range of scientific literature, including the following titles:

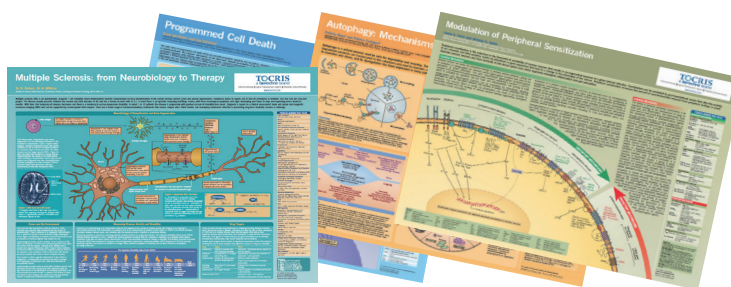
### Product Guides & Listings

- Neurodegeneration
- G Protein-Coupled Receptors
- Pain Research
- Ion Channels



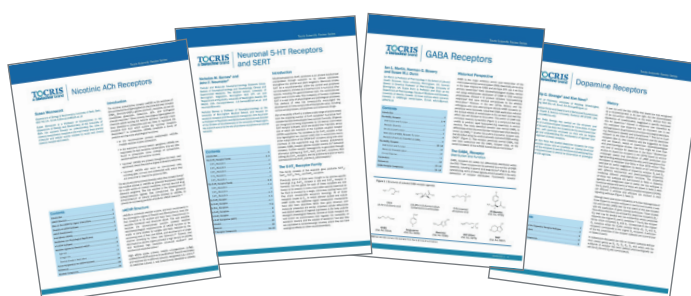
### Life Science Posters

- Multiple Sclerosis
- Pain
- Parkinson's Disease
- Schizophrenia



### Scientific Reviews

- Dopamine Receptors
- GABA Receptors
- mGlutamate Receptors
- Nicotinic ACh Receptors



For a complete selection of Tocris literature please visit | [tocris.com/literature](https://tocris.com/literature)

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



For research use or manufacturing purposes only. Trademarks and registered trademarks are the property of their respective owners.

TLT19-1505