

Ian L. Martin, Norman G. Bowery and Susan M.J. Dunn

Ian Martin is Professor of Pharmacology in the School of Life and Health Sciences, Aston University, Birmingham, UK. Norman Bowery is Emeritus Professor of Pharmacology, University of Birmingham, UK. Susan Dunn is Professor and Chair at the Department of Pharmacology, Faculty of Medicine and Dentistry, University of Alberta, Canada. All three authors share common interests in GABAergic transmission. E-mail: sdunn@pmcol.ualberta.ca

Contents

Introduction.....	1
The GABA_A Receptor.....	1-9
Distribution and Function.....	1
Receptor Diversity.....	2
Structure and Function.....	3
Modulators of GABA _A Receptor Function.....	6
Receptors Previously Classified as GABA _C	9
The GABA_B Receptor.....	10-12
Distribution and Function.....	10
Molecular Characterization.....	10
Clinical Potential.....	11
Conclusions.....	12
References.....	12-13
GABA Receptor Compounds.....	14-15

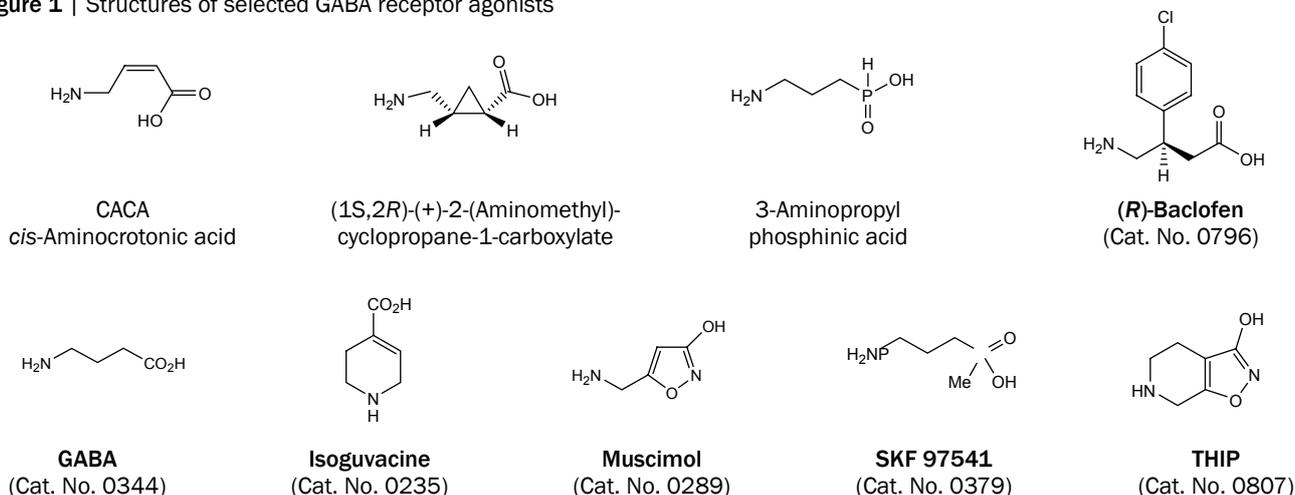
Historical Perspective

GABA is the major inhibitory amino acid transmitter of the mammalian central nervous system (CNS). Essentially all neurons in the brain respond to GABA and perhaps 20% use it as their primary transmitter.¹ Early electrophysiological studies, carried out using iontophoretic application of GABA to CNS neuronal preparations, showed it to produce inhibitory hyperpolarizing responses² that were blocked competitively by the alkaloid bicuculline.³ However, in the late 1970s, Bowery and his colleagues, who were attempting to identify GABA receptors on peripheral nerve terminals, noted that GABA application reduced the evoked release of noradrenalin in the rat heart and that this effect was not blocked by bicuculline. This action of GABA was mimicked, however, by baclofen (Figure 1), a compound that was unable to produce rapid hyperpolarizing responses in central neurons. This newly identified receptor was named GABA_B to differentiate it from the more familiar receptor type which became known as GABA_A.^{4,5} Another bicuculline-insensitive receptor was first identified using the conformationally restricted GABA analog, CACA^{6,7} (Figure 1). This receptor, previously termed GABA_C, has now been subsumed into the GABA_A receptor class, on the recommendation of the IUPHAR Nomenclature Committee.⁸

The GABA_A Receptor Distribution and Function

GABA_A receptors are widely but differentially distributed within the CNS.⁹ These receptors can be activated by a number of GABA isosteres, including muscimol and isoguvacine¹⁰ (Figure 1). After radiolabeling, some of these ligands proved valuable in the early delineation of receptor distribution. Functionally, receptor

Figure 1 | Structures of selected GABA receptor agonists



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

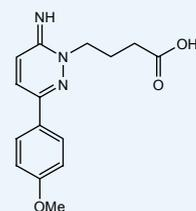
activation results in an increased membrane chloride conductance,^{11,12} usually causing an influx of Cl⁻ and membrane hyperpolarization. In general, concentration-response curves exhibit positive cooperativity, which is consistent with the presence of at least two agonist binding sites on each receptor molecule.¹³⁻¹⁵ On continued exposure to high agonist concentrations, the agonist-induced current decreases as a consequence of receptor desensitization.¹⁶⁻¹⁸ Biophysical characterization of the receptor, carried out initially using noise analysis of neurons in primary culture, provided the first estimates of mean single channel conductance and average channel open times,¹⁹ the latter of which varied with the nature of the activating agonist.²⁰ Development of single channel recording techniques²¹ provided further detail on the nature of single channel events with the demonstration of multiple single channel conductances: 44, 30, 19 and 12pS,²² the 30pS conductance being the most prevalent. Both channel opening times and opening frequency are dependent on agonist concentration and the competitive antagonist, bicuculline, reduces the conductance by modulating both of these parameters.^{23,24} Other competitive antagonists include the pyridazinyl GABA derivative, SR 95531 (Figure 2). The receptor can also be blocked non-competitively by picrotoxin and a number of bicyclophosphates.²⁵ In addition, penicillin decreases channel open probability in a manner that is compatible with open channel block.²⁶

Receptor Diversity

Purification of the bovine brain receptor in the early 1980s revealed two major subunits of the GABA_A receptor, which were named α and β . Elucidation of partial amino acid sequences of these subunits allowed subunit-specific monoclonal antibodies to be raised, thus providing the opportunity to explore the fine anatomical detail of receptor distribution.²⁷ The sequence data

SR 95531 hydrobromide, a selective, competitive GABA_A antagonist

Cat. No. 1262



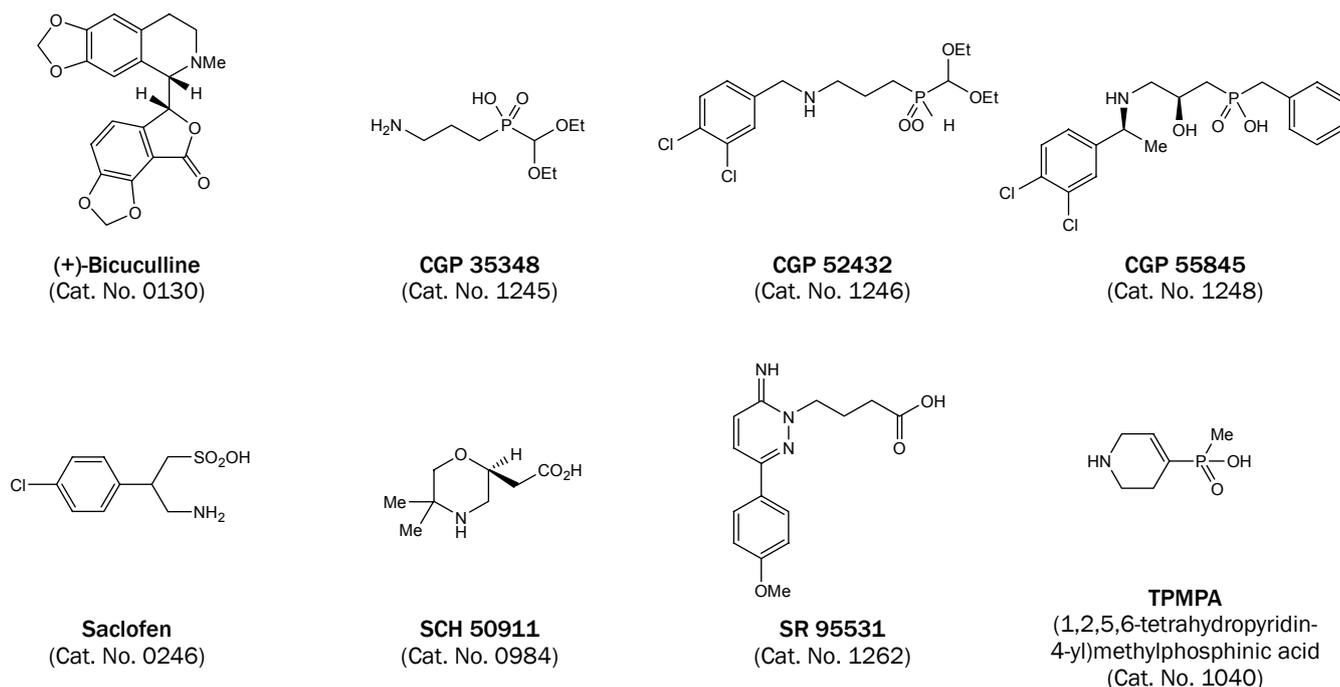
SR 95531 is a selective, competitive GABA_A receptor antagonist that displaces [³H]-GABA from rat brain membranes with a K_i value of 150 nM. Unlike bicuculline, SR 95531 selectively antagonizes GABA-induced Cl⁻ currents with little action on pentobarbitone-induced currents. The compound also acts as a low affinity glycine receptor antagonist.

Heaulme *et al.* (1986) *Brain Res.* **384** 224. Uchida *et al.* (1996) *Eur.J.Pharmacol.* **307** 89. Beato *et al.* (2007) *J.Physiol.* **580** 171.

also facilitated the cloning of the first two GABA_A receptor subunit isoforms.^{28,29}

Subsequent molecular studies revealed a multiplicity of protein subunits that have now been divided into seven classes, based on the extent of similarities in their deduced amino acid sequences. Within these classes there are further subdivisions into subunit isoforms, some of which exhibit alternate splice variants. In man, six α -, three β -, three γ - and three ρ - subunit isoforms are presently known, together with single representatives of the δ , ϵ , π and θ classes. Within a single subunit class, the sequence homology is about 70% but between classes this falls to around 30%. Additional isoforms of some of these classes are known in other species.³⁰ In the earlier receptor nomenclature,

Figure 2 | Structures of selected GABA receptor agonists



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

the three ρ -subunits were considered to define the GABA_c receptor.⁸

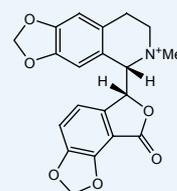
Deduced amino acid sequences from each of the subunits reveal homologies and a common structural organization which places them firmly within the so-called Cys-loop ligand-gated ion channel (LGIC) family. These receptors are pentamers of homologous subunits that assemble to form a central ion channel traversing the cell membrane. The archetypal member of the family is the peripheral nicotinic acetylcholine receptor (nAChR) with other members of the family including glycine and 5-HT₃ receptors. Each subunit has a long amino terminal domain of more than 200 amino acids which carries the signature cys-cys loop. This extracellular domain is followed by four hydrophobic segments, each of which is about 20 amino acids long. These four segments, termed TM1–TM4, were predicted to form transmembrane domains with TM2 contributing to formation of the ion channel lining. Between TM3 and TM4 there is a large intracellular loop, which is the most divergent part of the sequence within the GABA_A receptor subfamily.

Despite the plethora of receptor subunits, current evidence suggests that only a limited number of GABA_A receptor subunit combinations are expressed *in vivo*.³¹ Each subunit is encoded by a separate gene and a combination of *in situ* hybridization and immunohistochemical studies has revealed a distinct distribution of the various gene products in the CNS.^{32,33} This is consistent with the idea that each receptor subtype, made up of different combinations of subunits, serves defined physiological roles. In turn, this provides valuable information for development of subtype-selective pharmaceutical agents. However, an added complexity is that the expression patterns of individual subunits are not immutable. These can change during development, in response to normal physiological cycles and also as a consequence of pharmacological intervention.^{34–37}

Most receptors in the mammalian CNS comprise α -, β - and γ -subunits, with the most ubiquitous receptor subtype containing the $\alpha 1$, $\beta 2$ and $\gamma 2$ isoforms.³⁸ The recognition and functional

(-)-Bicuculline methochloride, a water-soluble GABA_A antagonist

Cat. No. 0131



(-)-Bicuculline methochloride is a water soluble and more stable salt of (+)-bicuculline (Cat. No. 0130) that acts as a competitive GABA_A receptor antagonist. The compound blocks inhibitory hyperpolarizing responses and reduces Cl⁻ conductance by modulating channel opening time and frequency.

Kemp et al. (1986) *Br.J.Pharmacol.* **87** 677. MacDonald et al. (1989) *J.Physiol.* **410** 479. Seutin and Johnson (1999) *TIPS* **20** 268.

characteristics of individual GABA_A receptor subtypes have been explored extensively using recombinant receptors expressed in mammalian cells or *Xenopus* oocytes. Many mutagenesis studies have been carried out to determine the roles of individual subunits, peptide segments and specific amino acids in receptor function. It is clear that in order to interpret mutagenic results effectively at the molecular level, it is essential to have an accurate view of the overall structure of the receptor.

Structure and Function

When the sequence homologies of many subunits of the Cys-loop LGIC family were first revealed, it seemed reasonable to predict that all members would share the same structural organization as the *Torpedo* nAChR. This is the best characterized member of the family and it has been elegantly imaged using cryoelectron microscopy, most recently at a resolution of 4Å.³⁹ It is a pentamer of homologous subunits that are arranged pseudosymmetrically around an integral ion channel. Using electron microscopy to

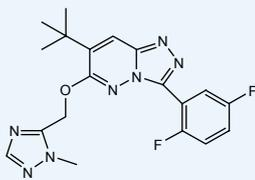
Table 1 | Comparative pharmacology of GABA receptors

Compound	GABA _A	GABA _B	GABA _{A-p} (formerly GABA _C)	Reference
GABA	Agonist	Agonist	Agonist	-
Muscimol	Agonist	Inactive	Partial agonist	5, 7, 158
Isoguvacine	Agonist	Inactive	Antagonist	5, 7
THIP	Agonist	Inactive	Antagonist	5, 7
P4S	Agonist	Inactive	Antagonist	5, 7
TACA	Agonist	Inactive	Agonist	7
CACA	Inactive	Inactive	Partial agonist	7
(R)-Baclofen	Inactive	Agonist	Inactive	5, 7
Bicuculline	Antagonist	Inactive	Inactive	5, 7
Picrotoxin	Antagonist	Inactive	Antagonist	5, 7
CGP 35348	Inactive	Antagonist	Inactive	159
CGP 54626	Inactive	Antagonist	Inactive	159
CGP 64213	Inactive	Antagonist	Inactive	159
SCH 50911	Inactive	Antagonist	Inactive	159
TPMPA	Inactive	Inactive	Antagonist	7, 160, 161

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

L-838,417, a subtype-selective GABA_A partial agonist

Cat. No. 3250



L-838,417 is a subtype-selective GABA_A receptor partial agonist. It selectively binds to $\alpha 1$, $\alpha 2$, $\alpha 3$ and $\alpha 5$ subunits (K_i values are 0.79, 0.67, 0.67 and 2.25 nM respectively) but displays no efficacy at $\alpha 1$ ($\alpha 1$ -sparing). The compound exhibits non-sedative anxiolytic, antinociceptive and anti-inflammatory activity *in vivo*.

McCabe et al. (2004) *Neuropharmacology* 46 171. McMahon and France (2006) *Br.J.Pharmacol.* 147 260. Knabi et al. (2008) *Nature* 451 330.

image the purified porcine GABA_A receptor by negative staining, a pentameric structure of similar diameter (about 8 nm across the pentamer) was revealed.⁴⁰ It is now believed that the most abundant $\alpha 1\beta 2\gamma 2$ GABA_A receptor subtype comprises two copies each of the $\alpha 1$ - and $\beta 2$ -subunits together with a single $\gamma 2$ -subunit.⁴¹ The arrangement of the subunits within the pentamer was first studied by concatenation,⁴² an approach that involves physical linking of the cDNAs encoding two or more subunits prior to their ectopic expression with other subunits. Such studies have demonstrated a pentameric subunit arrangement of β - α - β - α - γ lying in an anticlockwise direction when viewed from the outside of the cell.⁴³ With this information, it has been possible to use the 4Å structure of the *Torpedo* nAChR as a template to construct *in silico* models of this most common GABA_A receptor subtype^{44,45} (Figures 3a and 3b). These models provide a means to explore the similarities and differences in the structure and function of different members of the GABA_A receptor subfamily.

The exploration of ligand recognition in the extracellular domains of the Cys-loop family of receptors has been continuing for almost four decades. In 2001, rejuvenation of interest in this area came from a somewhat unexpected source. The structure of a water-soluble acetylcholine binding protein (AChBP) from *Lymnaea stagnalis* was determined at 2.7Å resolution,⁴⁶ a structure that soon proved to be a valuable homolog of the extracellular segment of the nAChR and other members of the family. This protein was the first to be used as a template to model the extracellular domain of the $\alpha 1\beta 2\gamma 2$ GABA_A receptor.⁴⁷ Together with the plethora of mutagenic data available in the literature at the time, this model furnished the first direct structural evidence that was compatible with the long-standing idea that GABA recognition sites were located at the β - α interfaces. In addition, it rationalized a great deal of experimental data which suggested that an allosteric site for the classical benzodiazepines lies in a similar position at the adjacent α - γ interface. In the case of the GABA activation sites, the current consensus is that the primary determinants of agonist recognition are found within at least six non-contiguous stretches ('loops') of amino acids in the extracellular domains of each subunit, loops A-C being contributed by the 'principal' subunit (β) and loops D-F by the neighboring 'subordinate' subunit (α , Figure 3c). Sequence

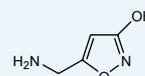
comparisons of the 'recognition loops' in different subunits of the receptor superfamily reveals some homology. However, it is the structural divergence within these loops that provides the exquisite acuity of ligand recognition which differentiates the family members.

The value of this *in silico* approach has proved significant. Not only does it allow visualization of the disposition of the amino acids involved in ligand recognition, but also, using theoretical ligand docking approaches, it becomes feasible to address receptor subtype-selective ligand design, an area that is of undoubted commercial interest.⁴⁸ Since binding sites for both neurotransmitters and allosteric modulators occur at subunit-subunit interfaces, we must again consider the importance of the subunit arrangement within each pentameric receptor. As discussed above, there is considerable theoretical and experimental evidence to assume that the subunit arrangement of the $\alpha 1\beta 2\gamma 2$ receptor is secure. However, there is no *a priori* reason to assume that less abundant receptors should adopt a similar pentameric architecture. What, for example, is the arrangement of subunits in receptor subtypes comprising $\alpha\beta\delta$ -subunits? To address this question, atomic force microscopy (AFM) was recently used to investigate the $\alpha 4\beta 3\delta$ subtype.⁴⁹ The subunits were C-terminally tagged with different epitopes and, after ectopic expression and decoration with the appropriate antibodies, the receptor-antibody complexes were visualized by AFM. The results suggested a similar arrangement to the $\alpha 1\beta 2\gamma 2$ subtype with the δ -subunit simply replacing the γ -subunit within the pentamer. However, the possibility of heterogeneity in receptor assembly cannot be excluded and, for example, results from concatenation studies suggest that the δ -subunit may be a little more promiscuous than first suggested.⁵⁰

Comparison of the two receptor subtypes described above ($\alpha 1\beta 2\gamma 2$ and $\alpha 4\beta 3\delta$) is important from both a physiological and pharmacological perspective. During the last several years, it has become clear that there are two major types of GABA_A receptor-mediated inhibitory responses i.e. phasic and tonic.⁵¹ Phasic inhibition results from activation of GABA_A receptors that are localized primarily to the synapse, such as the abundant $\alpha 1\beta 2\gamma 2$ subtype. Tonic transmission is mediated by less abundant extrasynaptic receptors, including the $\alpha 4\beta 3\delta$ subtype, that are thought to be activated by the low concentrations of the natural agonist which escape the efficient re-uptake machinery found in both neurons and glia. There is currently considerable interest in developing drugs that have differential effects on these two forms of inhibitory neurotransmission. Although the natural

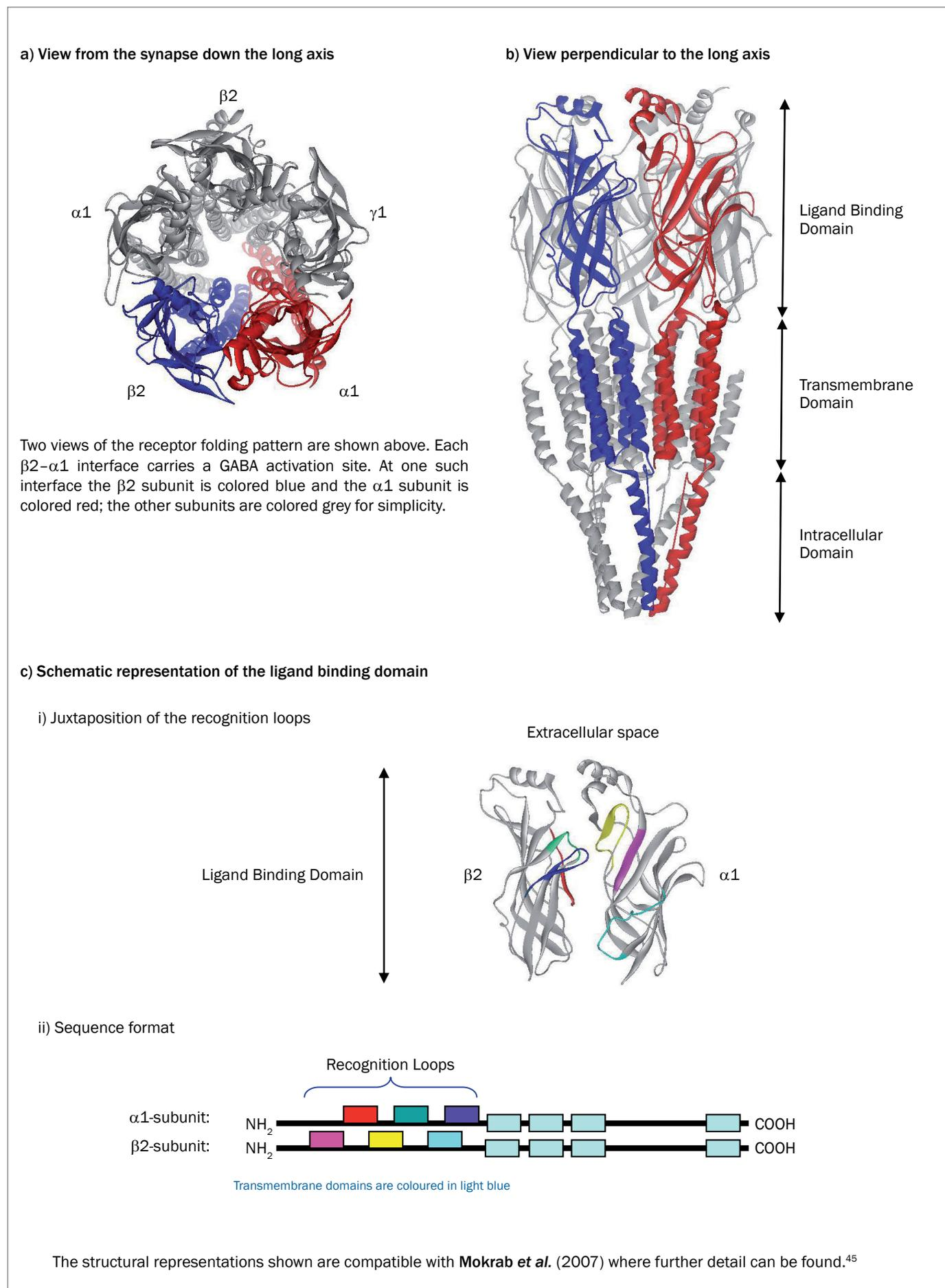
Muscimol, a potent GABA_A agonist

Cat. No. 0289



Muscimol is a potent GABA_A receptor agonist and partial GABA_C receptor agonist. The compound inhibits memory retention via central GABA_A receptors and attenuates airway constriction via peripheral GABA_A receptors.

Johnstone et al. (1996) *TIPS* 17 319. Gleason et al. (2009) *J.Appl.Physiol.* 106 1257. Jafari-Sabet and Jannat-Dastjerdi (2009) *Behav.Brain Res* 202 5.

Figure 3 | Model of the $\alpha 1\beta 2\gamma 2$ GABA_A receptor structure

agonist, GABA, appears to be a full agonist at the $\alpha 1\beta 2\gamma 2$ receptor, its conformationally restricted analog, THIP (also known as gaboxadol, Figure 1), is a partial agonist. In contrast, THIP is a full agonist at the $\alpha 4\beta 3\delta$ receptor where GABA acts as a partial agonist. Interestingly THIP exhibits hypnotic properties⁵² which are functionally quite distinct from those seen with the most widely used hypnotics, namely zopiclone and the $\alpha 1$ -selective agents, zolpidem and zaleplon. The latter compounds facilitate phasic inhibition by interacting with classical benzodiazepine-sensitive receptors at the synapse. THIP appears to produce its effects by modulating tonic inhibition mediated by extrasynaptic receptors,⁵³ which may also be selective targets for general anesthetics.

Modulators of GABA_A Receptor Function

The Benzodiazepines

The therapeutic importance of the benzodiazepines has been a significant impetus to GABA_A receptor research. Classical benzodiazepines potentiate agonist-mediated activation of the GABA_A receptor by causing a parallel leftward shift of the GABA concentration-response curve. In 1976, the discovery of saturable, high affinity binding sites for [³H]-diazepam in the brain^{54,55} provided an important experimental tool for their study. All of the overt effects of the benzodiazepines: sedative, anxiolytic, anticonvulsant, muscle relaxant and amnestic, are mediated by GABA_A receptors. However, not all the GABA_A receptors recognize the benzodiazepines. The particular α -subunit isoform present within an individual GABA_A receptor subtype is the primary determinant of benzodiazepine recognition (Table 2). If the $\alpha 1$ -subunit of the most common GABA_A receptor ($\alpha 1\beta 2\gamma 2$) is replaced by $\alpha 4$ or $\alpha 6$ the receptor fails to recognize the classical benzodiazepines. It is now clear from both biochemical and mutational analysis that this insensitivity can be attributed to a single amino acid substitution in the extracellular N-terminal domain: a histidine (H101) in the $\alpha 1$ -, $\alpha 2$ -, $\alpha 3$ - and $\alpha 5$ -subunits is replaced by an arginine residue in $\alpha 4$ and $\alpha 6$.^{56,57} When receptors containing the former subunits are expressed with a β - and $\gamma 2$ -subunit, all are recognized by the classical benzodiazepines. However, several agents differentiate between the subtypes on the basis of the particular α -subunit isoform present in the pentamer. The first of these compounds to be identified was the triazolopyridazine CL 21887258 (Figure 4),

which is related to the recently introduced hypnotic, zaleplon. Similarly β -carboline-3-carboxylic acid esters also show a preference for certain α -subunit-containing receptors.⁵⁹ Zolpidem (Figure 4), currently the most widely prescribed hypnotic in the USA, has been shown to have high affinity for $\alpha 1$ -containing receptors, lower affinity for receptors carrying $\alpha 2$ or $\alpha 3$ very low affinity for those containing $\alpha 5$ ^{60,61} (Table 2) and no observable interaction with receptors which contain the $\alpha 4$ - or $\alpha 6$ -subunits.

Using knockin (KI) technology, the importance of the α -subunit histidine-arginine substitution has been turned into an advantage. The exquisite specificity of this switch dictates that, by replacing the $\alpha 1$ histidine with an arginine (H101R) in the germ line, the KI adult animals will differ from their wild-type counterparts only in the ability of their $\alpha 1$ -containing receptors to recognize the benzodiazepines. Thus, it was expected that characterization of the knockin mouse phenotype would allow the complex pharmacological effects of benzodiazepines to be dissected based on their interactions with specific GABA_A receptor subtypes. Extensions of this approach have proved particularly valuable; it is now clear that the $\alpha 1$ -subunit is responsible for the sedative, anterograde amnestic and some of the anticonvulsant effects of the benzodiazepines,^{62,63} whereas the $\alpha 2$ -subunit has been associated with their anxiolytic actions.⁶⁴ Not all of the results are clear cut; for example, the pharmacodynamic profile of the $\alpha 3$ -selective ligand, TP003, suggests a contribution of this subunit to both anxiolytic and anticonvulsant effects.^{65,66} Also, receptors containing the $\alpha 5$ -subunit have been implicated in learning and memory processes.⁶⁷ This approach has been significantly advanced recently using conditional knockin studies, which have revealed selective changes in the ability of GABA_A receptors within particular cell groups to recognize the hypnotic, zolpidem.⁶⁸ Unfortunately, attempts to delineate the functional importance of individual GABA_A receptor subunits using gene knockout technology have proved frustrating. It is clear that ablation of subunit expression frequently results in compensatory changes in the expression of other subunits, providing significant challenges in assigning specific responsibility for the resulting phenotype.³⁸

Perhaps one of the most interesting phenomenological observations to arise from studies of benzodiazepine interactions with the GABA_A receptors has been the development of the inverse agonist concept. Non-benzodiazepine ligands were

Table 2 | Affinity of benzodiazepine site ligands for GABA_A receptor subtypes

Compound	$\alpha 1\beta 2\gamma 2$	$\alpha 2\beta 2\gamma 2$	$\alpha 3\beta 2\gamma 2$	$\alpha 4\beta 2\gamma 2$	$\alpha 5\beta 2\gamma 2$	$\alpha 6\beta 2\gamma 2$
Diazepam	16.1	16.9	17.0	>10,000	14.9	>10,000
Clonazepam	1.3	1.7	2.0	–	–	>10,000
Triazolam	1.8	1.2	3.0	–	1.2	–
Bretazenil	1.2	1.2	1.3	–	2.4	–
Flumazenil	1.0	1.1	1.5	107	0.4	90
Ro 15-4513	10.4	5.5	7.8	5.0	0.5	5.1
CL 218872	130	1820	1530	>10,000	490	>10,000
β -CCM	1.7	6.5	4.1	–	27	2050
Zolpidem	17	291	357	–	>15,000	–

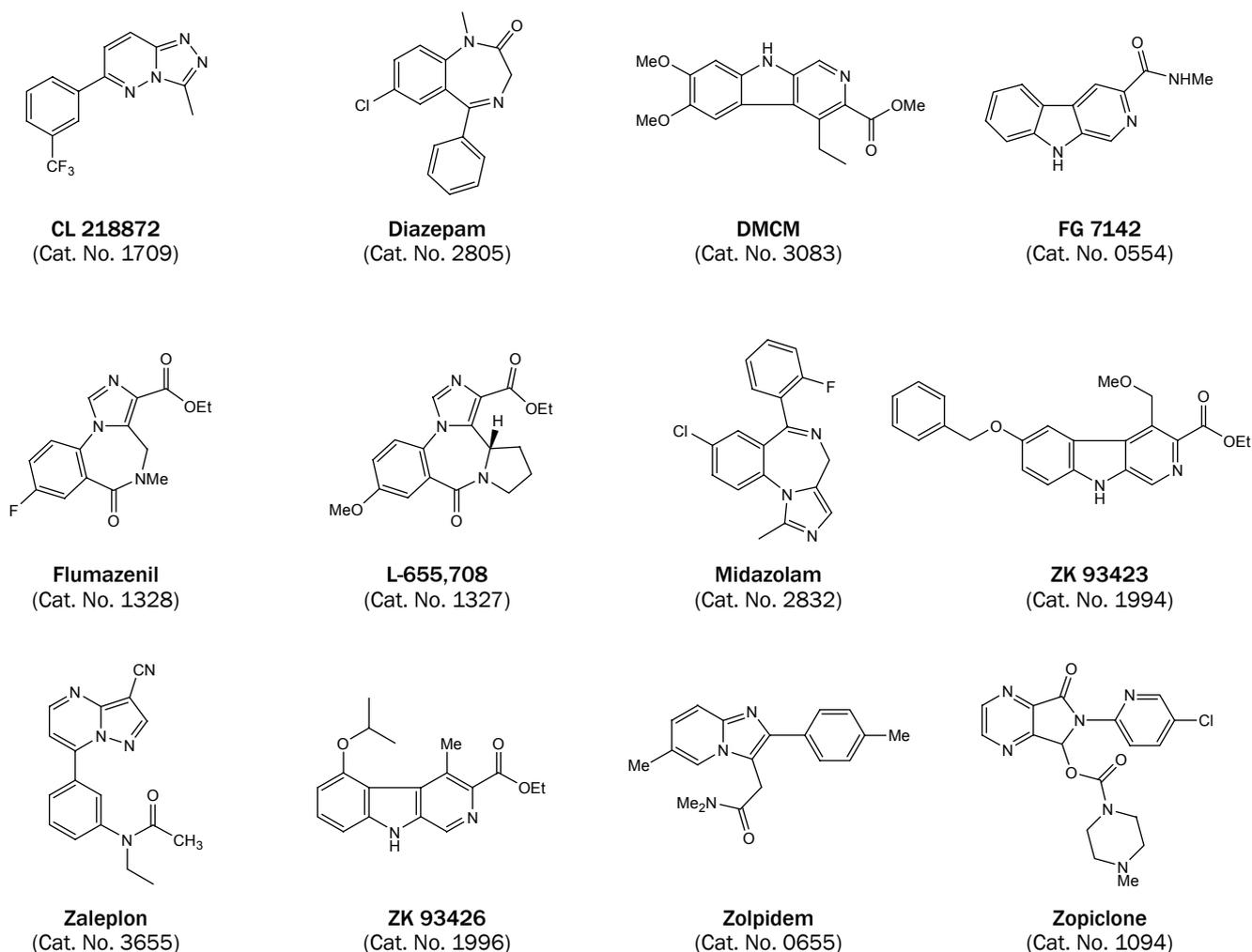
(Bold text denotes compounds available from Tocris at time of publication)

discovered that were able to displace a radiolabeled benzodiazepine from its binding sites. One of the first of these was ethyl β -carboline-3-carboxylate (β -CCE) which was shown to have effects that were diametrically opposed to those of the classical benzodiazepines e.g. it is proconvulsant. This led to a new terminology; β -CCE became known as an inverse agonist with the classical benzodiazepines then being classified as agonists.^{69,70} *In vitro* electrophysiological experiments using inverse agonists show that they shift the GABA concentration-response curve to the right, decreasing the potency of the natural transmitter. Thus, while the agonist benzodiazepine site ligands increase channel opening frequency, the inverse agonists decrease it.⁷¹ The full efficacy spectrum is found within the β -carboline series: the ethyl ester is proconvulsant and thus acts as a partial inverse agonist, the propyl ester is essentially devoid of efficacy leading it to be termed an antagonist,⁷² while aromatic substitution in the A ring produces agonists with similar properties to the classical benzodiazepines⁷³ (Figure 4). The therapeutic potential afforded by the inverse agonist concept has not escaped the attention of the pharmaceutical industry with the development of partial inverse agonists selective for $\alpha 5$ -containing receptor subtypes as cognition enhancers.⁷⁴

Steroids

The observation that 5 α -pregnan-3 α -ol-11,20-dione (alphaxalone; Figure 5), a synthetic steroidal anesthetic, was able to enhance stimulation-evoked inhibition produced by GABA_A receptor agonists in rat cuneate nucleus slices,⁷⁵ was the first evidence for allosteric steroid sites on these receptors. Subsequent voltage clamp studies conducted on both neurons and adrenomedullary chromaffin cells^{76,77} confirmed the stereoselective activity of the progesterone metabolites 5 α -pregnan-3 α -ol-20-one (allopregnanolone), 5 β -pregnan-3 α -ol-20-one (pregnanolone) and 5 α -pregnan-3 α ,21-diol-20-one (allotetrahydrodeoxycorticosterone). Mechanistically, the action of these compounds appeared to be similar to that of the barbiturates which, at low concentrations, potentiate the effects of GABA by increasing channel open times and, at higher concentrations, directly activate the receptor.^{78–82} Later studies revealed that the sites of barbiturate and steroid action are distinct.⁸³ Conserved residues within the α - and β -subunit membrane spanning domains of the $\alpha 1\beta 2\gamma 2$ receptor, which are important for both steroid facilitation and direct activation, have been identified.⁸⁴

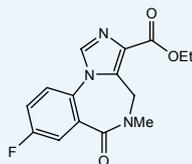
Figure 4 | Structures of selected benzodiazepine site ligands



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

Flumazenil, a benzodiazepine antagonist

Cat. No. 1328



Flumazenil is a benzodiazepine antagonist that is nonselective for $\alpha 1$, $\alpha 2$, $\alpha 3$ or $\alpha 5$ -containing GABA_A receptors. The compound reverses benzodiazepine sedation and is centrally active upon systemic administration *in vivo*.

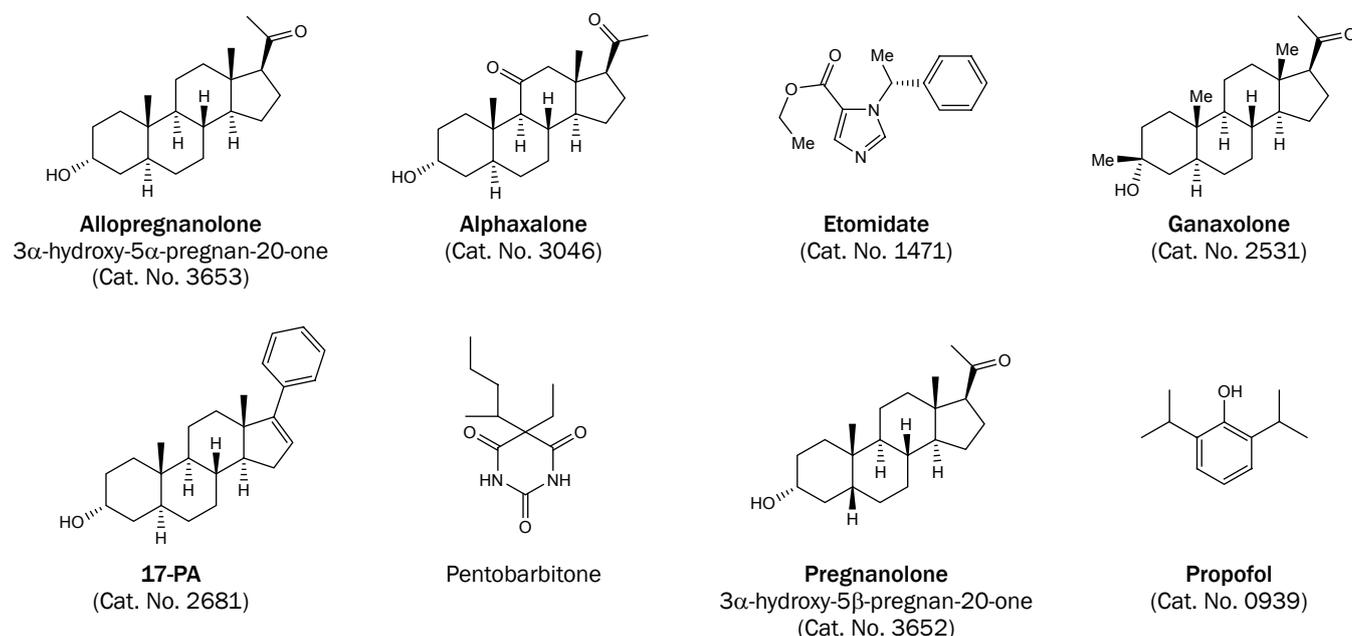
Polc et al. (1981) *Naunyn-Schmied.Arch.Pharmacol.* **316** 317. **Atack et al.** (1999) *Neuropsychopharmacology* **20** 255. **Doble** (1999) *J.Psychopharmacol.* **13** S11.

Studies with ectopically expressed receptors comprising $\alpha\beta\gamma$ -subunits demonstrated limited impact of subunit composition on the functional effects of the steroids.⁸⁵ At putative extrasynaptic receptors, where δ replaces the γ -subunit, there is evidence for increased steroid potency.⁸⁶ However, these observations may be explained, at least in part, by the reduced efficacy of the endogenous neurotransmitter, GABA, at these receptor subtypes.⁸⁷ There have been many literature reports to demonstrate that the potency of steroids varies in different brain regions and there is also evidence to suggest that the observed effects may be influenced by changes in receptor phosphorylation and modulation of enzymatic activity in the steroid metabolic pathways.⁸⁸ This functional complexity is amplified further by normal physiological fluctuations in steroid levels associated with, for example, pregnancy and the ovarian cycle. These can lead to altered patterns of subunit expression that may contribute to the mood swings that are associated with these events.^{89,37}

General Anesthetics

It is now clear that GABA_A receptors play a significant role in general anesthesia. Many of the receptor subtypes are sensitive to clinically relevant concentrations of general anesthetics and exhibit the appropriate stereospecificity. The characteristics of these agents are diffuse; they exhibit sedative, hypnotic, analgesic and amnesic properties in addition to producing a loss of mobility.⁹⁰ This multiplicity of effects, together with their structural diversity, has meant that it is very difficult to dissect the actions of general anesthetics at the molecular level. One agent, ketamine, mainly affects glutamatergic excitatory responses mediated by NMDA receptors and there is no evidence that the older anesthetics, nitrous oxide and xenon, modulate GABAergic inhibition. The evidence for interactions of other inhalational and intravenous agents with the GABA_A receptors continues to grow. Since general anesthetics are hydrophobic and need to access the CNS, it is perhaps not surprising that they target hydrophobic pockets within the transmembrane domains of the receptor. Initial evidence suggested that the inhalational anesthetics favored the α -subunits⁹¹ while *in vitro* and *in vivo* evidence has accumulated to suggest that intravenous anesthetics interact with the β -subunits.^{92,93} Over the past decade it has become increasingly clear that significant effects of the general anesthetics occur not by their ability to potentiate the fast phasic inhibition mediated by synaptically located receptors but as a result of their effects on receptors that are located extrasynaptically. The extrasynaptic $\alpha 5\beta 3\gamma 2$ receptor in the hippocampus is probably associated with the amnesic actions of many of these agents,⁹⁴ while those receptors containing δ -subunits in the ventrobasal thalamic nucleus provide the intriguing link between the reversible loss of consciousness in man, a sleep-related phenomenon that is a primary characteristic induced by the general anesthetics.⁹⁵ It is clear that the diversity

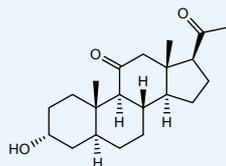
Figure 5 | Structures of selected compounds active at allosteric sites of GABA_A receptors



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

Alphaxalone, a direct activator and potentiator of GABA_A

Cat. No. 3046



Alphaxalone is a neurosteroid anesthetic that directly activates and potentiates the GABA_A receptor-activated membrane current (I_{GABA}). Efficacy, but not potency, of this compound is determined by the alpha subunit of the receptor (EC_{50} values are 1.4, 1.8, 2.1, 2.4 and 2.5 μ M for $\alpha 1\beta 1\gamma 3$, $\alpha 1\beta 1\gamma 1$, $\beta 1\gamma 1$, $\alpha 2\beta 1\gamma 2L$ and $\alpha 1\beta 1\gamma 2L$ isoforms respectively).

Maitra et al. (1999) *Brain Res.* **819** 75. Wegner et al. (2007) *Neuropharmacology* **52** 672. Zecharia et al. (2009) *J. Neurosci.* **29** 2177.

of agents and targets provide valuable clues that must be addressed systematically to optimize the potential for development of novel anesthetic agents.⁹⁶

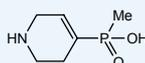
Alcohol

The receptors responsible for the pharmacological effects of ethanol have been the subject of much speculation but there are only a small number of putative targets that are responsive to low concentrations of ethanol (< 20 mM). Initial observations of altered ethanol-induced behaviors in δ -subunit knockout mice⁹⁷ were followed by *in vitro* studies of recombinant GABA_A receptors. There was significant excitement when it was reported that agonist activation of δ -containing GABA_A receptors could be facilitated by 10 mM ethanol;^{98,99} however, replication of this response has not proved possible and these discrepancies remain unexplained.¹⁰⁰ It has been suggested that many of the *in vivo* effects of ethanol may be attributed to indirect effects arising from its ability to increase levels of several endogenous steroids which, in turn, can potentiate GABA_A receptor-mediated responses. Evidence in favor of this idea comes from observations that, not only do the consequential effects of ethanol correlate well with an increase in steroid levels, but they are also inhibited by blockers of steroid synthesis.¹⁰¹

Phosphorylation may again play a significant role since it has been noted that in PKC δ knockout mice, the pharmacological

TPMPA, a selective GABA_{A-p} antagonist

Cat. No. 1040



TPMPA is a selective, competitive GABA_{A-p} antagonist which exhibits only minimal effects on GABA_A and GABA_B receptors (K_b values are 2.1 μ M (antagonist), 320 μ M (antagonist) and $EC_{50} \sim 500$ μ M (weak agonist) respectively). It displays 8-fold selectivity for human recombinant $\rho 1$ receptors over $\rho 2$ receptors and blocks the paired-pulse depression component of inhibitory post-synaptic currents *in vitro*.

Murata et al. (1996) *Bioorg. Med. Chem. Lett.* **6** 2073. Chebib et al. (1998) *Eur. J. Pharmacol.* **357** 227. Xu et al. (2009) *Exp. Neurol.* **216** 243.

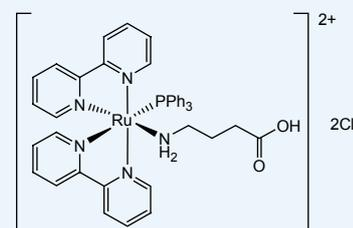
effects of ethanol are reduced as are the ataxic responses to both pentobarbital and pregnanolone. Since the flunitrazepam response remained intact in these animals, it was suggested that the overt effects were mediated by benzodiazepine insensitive GABA_A receptors. Supplementary studies showed that the PKC δ -dependent effects of ethanol could be observed in ectopically expressed $\alpha 4\beta 3\delta$ receptors.¹⁰² Thus, assignment of the effects of ethanol to specific GABA_A receptors remains enigmatic.

GABA_{A-p} Receptors (Previously Classified as GABA_C)

Although the receptors that were originally designated as GABA_C are now considered to be members of the GABA_A family,⁸ it is useful to highlight their distinguishing features. These receptors were originally classified on the basis of their unique pharmacology. The natural agonist, GABA, was reported to be about an order of magnitude more potent at this subclass than at other GABA_A receptors and, although CACA activated this receptor, this agent was not recognized by either the GABA_A or GABA_B classes (Figure 1). GABA_C receptor responses were not inhibited by bicuculline but, like the GABA_A receptors, they were blocked by picrotoxin. A selective GABA_C receptor antagonist, (1,2,5,6-tetrahydropyridine-4-yl)methylphosphonic acid7 (TPMPA, Figure 2) was later identified. Additional pharmacological differences from the GABA_A receptors included its lack of modulation by the benzodiazepines, barbiturates or neuroactive steroids. Receptors displaying these characteristics were shown to have a restricted distribution, initially being found in the spinal cord and subsequently in the retina,^{6,103} the source from which the first ρ -subunit was cloned.¹⁰⁴ Three homologous ρ -subunits, $\rho 1$ to $\rho 3$, have now been identified. These can be expressed as either homomers or heteromers^{105,106} and the ectopically expressed receptors exhibit the pharmacological characteristics of the elusive GABA_C receptors. There is only limited evidence that the ρ -subunits co-assemble with any of the other GABA_A receptor subunits.¹⁰⁷ The genes encoding the $\rho 1$ - and $\rho 2$ -subunits are found on chromosome 6 in humans, and are thus distinct from the clusters of GABA_A receptor subunit genes which are

RuBi-GABA, excitable by visible wavelength

Cat. No. 1040

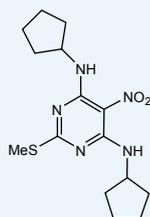


RuBi-GABA (Ruthenium-bipyridine-triphenylphosphine caged GABA) is excited by visible wavelengths and has two-photon uncaging capabilities. It provides greater tissue penetration, less phototoxicity, faster photorelease kinetics and better spatial resolution than UV light-sensitive caged compounds. Following photolysis, the compound produces GABA receptor-mediated currents in pyramidal neurons *in vitro* and displays no effect on endogenous GABAergic or glutamatergic transmission at concentrations effective for uncaging.

Zayat et al. (2003) *J. Am. Chem. Soc.* **125** 882. Nikolenko et al. (2005) *Chem. Comm.* **7** 1752. Rial Verde et al. (2008) *Front. Neural Circuits* **2** 2.

GS 39783, a positive modulator at GABA_B receptors

Cat. No. 2001



GS 39783 is a positive allosteric modulator of GABA_B receptor function. It potentiates the effects of GABA on [³⁵S]GTPγS binding at recombinant and native GABA_B receptors (EC₅₀ values are 2.1 and 3.1 μM respectively). The compound decreases cocaine self-administration, blocks the rewarding properties of nicotine and produces anxiolytic-like activity without the side effects associated with baclofen or benzodiazepines *in vivo*.

Urwyler *et al.* (2003) *J.Pharmacol.Exp.Ther.* **307** 322. Cryan *et al.* (2004) *J.Pharmacol.Exp.Ther.* **310** 952. Mombereau *et al.* (2007) *J.Pharmacol.Exp.Ther.* **321** 172.

found on chromosomes 4, 5, 15 and X with the exception of δ, which is found on chromosome 1. The ρ-subunit sequences display between 30 and 38% homology to the GABA_A receptor subunits at the amino acid level but, interestingly, in the important TM2 region of the sequence, they show greater homology to the glycine α-subunits than to any of the GABA_A receptor subunits.

The GABA_B Receptor

The other major class of GABA receptors is the metabotropic (G-protein-coupled) GABA_B receptor. These exhibit a distinct ligand recognition profile to the GABA_A receptor family,⁴ and are differentially distributed within the mammalian CNS.¹⁰⁸ Functionally they inhibit adenylyl cyclase activity¹⁰⁹ and presynaptic calcium channels, decreasing transmitter release,¹¹⁰ and activate postsynaptic potassium channels, producing the late inhibitory postsynaptic potential.¹¹¹

Distribution and Function

The initial observation that GABA inhibited the release of noradrenalin from rat atria *in vitro*, an effect not blocked by bicuculline methobromide but mimicked by (*R*)-baclofen, provided the seminal evidence to distinguish the GABA_B receptor from more familiar members of the GABA_A receptor family.⁴ Subsequent studies, using both functional and radioligand binding techniques, have further refined the structure-activity profile at GABA_B receptors¹¹² (Figures 1 and 2). Although the receptor is widely distributed within the mammalian CNS it is generally found at lower densities than GABA_A receptors, and exhibits a distinct distribution: the highest concentrations being found in the molecular layer of the cerebellum, the frontal cortex and certain thalamic nuclei.¹⁰⁸ The receptor is also found in the periphery where its activation modulates autonomic control of the intestine and decreases esophageal reflux.^{112,113} The receptor is coupled to adenylyl cyclase via G_i and G_o proteins. While the consequences remain poorly defined, activation of presynaptic GABA_B receptors also leads to the inhibition of high voltage-

activated Ca²⁺ channels, an effect that is mediated by the G protein βγ subunits. This results in decreased transmitter release and possibly also limits synaptic vesicle recruitment to the active zone.¹¹⁴

GABA activation of postsynaptic GABA_B receptors produces hyperpolarization via the modulation of inwardly rectifying K_{IR,3} type K⁺ channels¹¹⁵ that mediate the late phase of the inhibitory postsynaptic potential.

Molecular Characterization

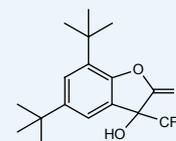
The molecular characterization of the GABA_B receptor was achieved in 1997 when the availability of specific high affinity antagonists allowed the expression cloning of the GABA_{B1} subunit.¹¹⁶ Subsequent studies demonstrated that while this protein showed many of the expected characteristics, when expressed ectopically, it coupled poorly to its effector machinery, and exhibited a remarkably low affinity for agonists compared to the native receptor; it appeared to be retained within the endoplasmic reticulum.¹¹⁷ Subsequent studies revealed the identity of an additional subunit, GABA_{B2},¹¹⁸⁻¹²⁰ which interacts with the GABA_{B1} subunit C-terminus, masking the ER retention signal of the GABA_{B1} subunit¹²¹ and facilitating the trafficking of the GABA_{B1} subunit to the cell surface. This provided the first secure evidence of receptor dimerization.

The GABA_B receptor belongs to the class C of the GPCRs, together with the metabotropic glutamate receptors mGlu1-8 and the calcium-sensing receptor.¹²² Each subunit comprises a large N-terminal extracellular domain exhibiting the venus fly-trap motif, followed by 7-transmembrane helices and an intracellular C-terminus. Two splice variants for the GABA_{B1} subunit are known; they are encoded by the same gene and arise by alternate promoter usage to produce GABA_{B1a} and GABA_{B1b}.¹²³ These differ only in their N-terminal domains, GABA_{B1a} contains a repeat of a conserved protein binding motif, so-called 'sushi domains', that are lacking in GABA_{B1b}; the first 147 amino acids of GABA_{B1a} are replaced by 18 amino acids in GABA_{B1b}.¹²⁴

While both subunits within the GABA_B heterodimer exhibit the venus fly-trap motif at the extracellular N-terminus, it is the GABA_{B1} subunit that is responsible for both agonist and antagonist recognition; the residues responsible are not conserved in the GABA_{B2} subunit.¹²⁵ Within this recognition domain there is also a serine residue that appears to be responsible for the ability of the

Rac BHFF, a potent, selective GABA_B positive allosteric modulator

Cat. No. 3313



Rac BHFF is a potent and selective GABA_B receptor positive allosteric modulator that increases the potency and efficacy of GABA (> 15-fold and > 149% respectively). The compound exhibits anxiolytic activity *in vivo* and is orally active.

Maiherbe *et al.* (2008) *Br.J.Pharmacol.* **154** 797.

receptor to sense Ca^{2+} concentrations.¹²⁶ While the $\text{GABA}_{\text{B}2}$ subunit is not primarily responsible for agonist recognition, its presence markedly increases the agonist affinity of the $\text{GABA}_{\text{B}1}$ subunit.¹²⁷ The $\text{GABA}_{\text{B}2}$ subunit mediates G protein-coupling, the second intracellular loop being particularly important,^{128,129} although it is clear that $\text{GABA}_{\text{B}1}$ is important in facilitating this process. It is the $\text{GABA}_{\text{B}2}$ subunit that appears to be the interaction site for an increasing family of positive allosteric modulators¹³⁰ (Figure 6), where binding occurs within the transmembrane domain¹³¹ to augment agonist tone while exhibiting no direct agonist activity.¹³²

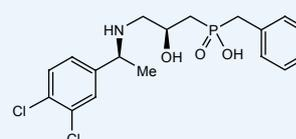
The restricted molecular heterogeneity found in the GABA_{B} receptor population has proved a significant frustration, since ectopic expression studies have failed to provide support for the varied functional responses ascribed to these receptors *in vivo*.¹³³ Knockout studies targeting $\text{GABA}_{\text{B}1}$ or $\text{GABA}_{\text{B}2}$ have not relieved these difficulties, both deletions producing similar phenotypes, although each compromised the expression of the conjugate subunit.^{134,135} Functional distinctions between the $\text{GABA}_{\text{B}1}$ subunit isoforms have started to emerge suggesting that the $\text{GABA}_{\text{B}1a}$ isoform is primarily associated with the heteroreceptors controlling glutamate release.¹³⁶⁻¹³⁹ It has been suggested that this differential cellular localization may be associated with the presence of the sushi repeats, present in $\text{GABA}_{\text{B}1a}$ but not in $\text{GABA}_{\text{B}1b}$, that are known to be important in protein-protein interactions in other environments.¹⁴⁰ Interestingly it has recently been shown that a soluble truncated form of the $\text{GABA}_{\text{B}1a}$ subunit, named $\text{GABA}_{\text{B}1j}$, exhibits nanomolar affinity for neuronal membranes. It is identical to the first 157 amino acids of the $\text{GABA}_{\text{B}1a}$ subunit and contains the sushi repeats together with a 72 amino acid C-terminal extension with no homology to other known proteins. In its presence both basal and stimulated glutamate release are decreased, but GABA_{B} receptor function at presynaptic autoreceptors or postsynaptic receptors remains unaffected.¹⁴¹

Clinical Potential

Baclofen remains the only clinically available agent that targets the GABA_{B} receptor. It was introduced into clinical practice in 1972 long before the discovery of the GABA_{B} receptor and remains the intervention of choice in spasticity associated with multiple sclerosis and cerebral palsy. Baclofen exhibits a challenging side effect profile on systemic administration, producing drowsiness, nausea, muscle weakness and mental

CGP 55845, a potent, selective GABA_{B} antagonist

Cat. No. 1248



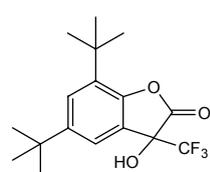
CGP 55845 is a potent, selective GABA_{B} receptor antagonist ($\text{IC}_{50} = 5 \text{ nM}$) that prevents agonist binding ($\text{pK}_i = 8.35$) and inhibits GABA and glutamate release (pEC_{50} values are 8.08 and 7.85 respectively). The compound inhibits GABA_{B} receptor responses to baclofen ($\text{IC}_{50} = 130 \text{ nM}$ in an isoproterenol assay) and potentiates the hypoglycemic response to glucose *in vitro*.

Waldmeier *et al.* (1994) *Br.J.Pharmacol.* **113** 1515. Cunningham and Enna (1996) *Brain Res.* **720** 220. Zhang *et al.* (2009) *J.Physiol.* **578** 735.

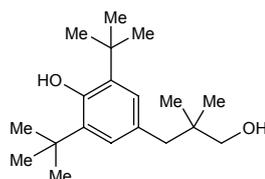
confusion largely due to poor brain penetration necessitating the use of high oral doses.¹⁴² The muscle relaxant effects mediated within the spinal cord can be secured by intrathecal administration, allowing a marked reduction in dose, thus limiting the systemic effects and the development of tolerance.¹⁴³ Baclofen also exhibits antinociceptive properties at the spinal level, which again allows local administration,¹⁴⁴ however significant analgesia is also mediated within the ventrobasal nucleus of the thalamus,¹⁴⁵ a site which requires systemic administration. The analgesic effects of baclofen currently have limited application in humans.¹⁴⁶

The early specific GABA_{B} receptor antagonists suffered from a limited potency, with phaclofen, for example, displaying an affinity of only $100 \mu\text{M}$. A number of selective, high affinity and systemically active antagonists are now available (Figure 2), that may have significant clinical potential in absence epilepsy.¹⁴² Mice overexpressing the $\text{GABA}_{\text{B}1a}$ isoform exhibit characteristics associated with atypical absence epilepsy.¹⁴⁷ In contrast, recent reports suggest that impaired GABA_{B} receptor function may contribute to repetitive firing in human temporal lobe epilepsy tissue.¹⁴⁸ The first exploration of GABA_{B} receptor antagonists clinically was an open trial with SGS 742 (CGP 36742) in mild cognitive deficit in man.¹⁴⁹ While the initial results appeared promising further clinical reports have not reached the literature. Recent studies suggest that mechanistically phospho-protein

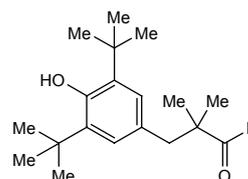
Figure 6 | Structures of selected allosteric modulators of the GABA_{B} receptor



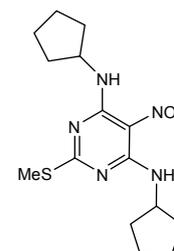
rac BFFF
(Cat. No. 3313)



CGP 7930
(Cat. No. 1513)



CGP 13501
(Cat. No. 1514)

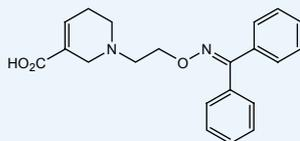


GS 39783
(Cat. No. 2001)

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

NNC 711, a selective inhibitor of GAT-1

Cat. No. 1779



NNC 711 is a potent and selective inhibitor of GABA uptake by GAT-1 (IC_{50} values are 0.04, 171, 1700 and 622 μ M for hGAT-1, rGAT-2, hGAT-3 and hBGT-1 respectively). The inhibitor displays anticonvulsant activity following systemic administration *in vivo*.

Suzdak *et al.* (1992) *Eur.J.Pharmacol.* **223** 189. Borden *et al.* (1994) *Eur. J.Pharmacol.* **269** 219. O'Connell *et al.* (2001) *Eur.J.Pharmacol.* **424** 37.

kinase A (pPKA) plays a significant role in the effects of this antagonist in the Morris water maze.¹⁵⁰

It has been known for some time that GABA_B receptor activation effectively reduces the craving for addictive drugs, first demonstrated as a reduction in cocaine self-administration in rats,¹⁵¹ and similar findings have emerged with other drugs of abuse.¹⁴² Positive allosteric modulators at the receptor may prove to be a more attractive means of control. These agents could reasonably be expected to facilitate GABA_B receptor mediated tone circumventing the side effect profile associated with the use of systemic agonists. Indeed, recent studies suggest that compounds of this type significantly reduce cocaine self-administration in rats^{152,153} (Figure 5), with similar approaches providing some support for their potential as anxiolytics.¹⁵⁴

The GABA_B receptors remain somewhat enigmatic, promissory notes of significant therapeutic potential have not thus far materialized and it has been argued that the lack of readily differentiable receptor subtypes has limited the opportunity for discrete drug targeting. Evidence for their functional importance continues to expand with recent studies highlighting their impact on both the tegmental pedunculopontine nucleus, important in the acute rewarding effects of the opiates¹⁵⁵ and orexin neurons, associated with sleep/ wakefulness cycles.¹⁵⁶ Clinical applications remain elusive; perhaps the allosteric modulators may yet prove as valuable in the modulation of tonic activity at the GABA_B receptors as the benzodiazepines within the GABA_A receptor family. Recent development of novel GABA_B receptor agonists, the structure of which effectively restricts them to the peripheral compartment, are currently under investigation for intervention in gastroesophageal reflux in man, after proving efficacious and importantly devoid of the baclofen-associated central side effects in preclinical studies.¹⁵⁷

Conclusions

Despite the overwhelming representation of the GPCRs in the human genome, it is the ionotropic receptor for the major inhibitory neurotransmitter GABA that has achieved the most visibility to date. Its serendipitous exploitation within the clinical arena has stimulated a plethora of intriguing insights into the mechanisms by which communication within the nervous system is achieved and the nuances of modulation which provide opportunities for further refinement of pharmacological

intervention. The paucity of molecular heterogeneity exhibited by the GABA_B receptors has proved problematic for specific drug targeting. The functional characterization of this receptor in the mammalian system was predicted to lead to significant clinical developments. Although this goal has not yet been achieved, recent research provides novel opportunities for therapeutic intervention. The next decade will undoubtedly prove exciting.

References

1. Somogyi *et al.* (1998) *Brain Res.Rev.* **26** 113.
2. Krnjevic and Schwartz (1967) *Exp.Brain Res.* **3** 320.
3. Curtis *et al.* (1970) *Nature* **226** 1222.
4. Bowery *et al.* (1980) *Nature* **283** 92.
5. Hill and Bowery (1981) *Nature* **290** 149.
6. Johnston *et al.* (1975) *J.Neurochem.* **24** 157.
7. Johnston (1996) *TIPS* **17** 319.
8. Barnard *et al.* (1998) *Pharmacol.Rev.* **65** 68-76.
9. Palacios *et al.* (1981) *Brain Res.* **222** 285.
10. Ebert *et al.* (1997) *Mol.Pharmacol.* **52** 1150.
11. Curtis *et al.* (1968) *Exp.Brain Res.* **5** 1.
12. Kelly *et al.* (1969) *Exp.Brain Res.* **7** 11.
13. Dichter (1980) *Brain.Res.* **190** 111.
14. Sakmann *et al.* (1983) *J.Neural Transm.* **18** 83.
15. Bormann and Clapham (1985) *Proc.Natl.Acad.Sci.USA* **82** 2168.
16. Adams and Brown (1975) *J.Physiol.* **250** 85.
17. Akaike *et al.* (1987) *J.Physiol.* **391** 219.
18. Mathers (1987) *Synapse* **1** 96.
19. Barker *et al.* (1982) *J.Physiol.* **322** 365.
20. Barker and Mathers (1981) *Science* **212** 358.
21. Hamill *et al.* (1981) *Pflugers Arch.* **391** 85.
22. Bormann *et al.* (1987) *J.Physiol.* **385** 243.
23. Macdonald *et al.* (1989) *J.Physiol.* **410** 479.
24. Twyman *et al.* (1990) *J.Physiol.* **423** 193.
25. Bowery *et al.* (1976) *Br.J.Pharmacol.* **57** 435.
26. Twyman *et al.* (1992) *J.Physiol.* **445** 97.
27. Richards *et al.* (1987) *J.Neurosci.* **7** 1866.
28. Stephenson and Barnard (1986) in Benzodiazepine/GABA chloride channels: structural and functional properties. p261 Alan R Liss.
29. Schofield *et al.* (1987) *Nature* **328** 221.
30. Hevers and Luddens (1998) *Mol.Neurobiol.* **18** 35.
31. Olsen and Sieghart (2008) *Pharmacol.Rev.* **60** 243.
32. Wisden *et al.* (1992) *J.Neurosci.* **12** 1040.
33. Pirker *et al.* (2000) *Neuroscience* **101** 815.
34. Laurie (1992) *J.Neurosci.* **12** 4151.
35. Holt *et al.* (1996) *Neuropharmacology* **35** 1457.
36. Smith *et al.* (1998) *J.Neurosci.* **18** 5275.
37. Griffiths and Lovick (2005) *Neuroscience* **136** 457.
38. Rudolph and Mohler (2004) *Ann.Rev.Pharmacol.Toxicol.* **44** 475.
39. Unwin (2005) *J.Mol.Biol.* **346** 967.
40. Nayeem *et al.* (1994) *J.Neurochem.* **62** 815.
41. Farrar *et al.* (1999) *J.Biol.Chem.* **274** 10100.
42. Im (1995) *J.Biol.Chem.* **270** 26063.
43. Baumann (2002) *J.Biol.Chem.* **277** 46020.
44. Ernst *et al.* (2005) *Mol.Pharmacol.* **68** 1291.
45. Mokrab *et al.* (2007) *J.Mol.Graph Model.* **26** 760.
46. Brejc *et al.* (2001) *Nature* **411** 269.
47. Cromer *et al.* (2002) *TIBS* **27** 280.
48. Clayton *et al.* (2007) *Curr.Med.Chem.* **14** 2755.
49. Barrera *et al.* (2008) *Mol.Pharmacol.* **73** 960.
50. Kaur (2009) *J.Biol.Chem.* **284** 7889.
51. Farrant and Nusser (2005) *Nature Neurosci.Rev.* **6** 215.
52. Krosggaard-Larson *et al.* (2004) *Biochem.Pharmacol.* **68** 1573.

53. Drasbek and Jensen (2006) *Cereb.Cortex* **16** 1134.
54. Squires and Braestrup (1976) *Nature* **266** 732.
55. Mohler and Okada (1977) *Science* **198** 849.
56. Wieland et al. (1992) *J.Biol.Chem.* **267** 1426.
57. Duncalfe et al. (1996) *J.Biol.Chem.* **271** 9209.
58. Squires et al. (1979) *Pharmacol.Biochem.Behav.* **10** 825.
59. Nielsen and Braestrup (1980) *Nature* **286** 606.
60. Hadingham et al. (1993) *Mol.Pharmacol.* **43** 970.
61. Sanger et al. (1994) *Neurosci.Behav.Rev.* **18** 355.
62. Rudolph et al. (1999) *Nature* **401** 796.
63. McKernan et al. (2000) *Nature Neurosci.* **3** 587.
64. Low et al. (2000) *Science* **290** 131.
65. Dias et al. (2005) *J.Neurosci.* **25** 10682.
66. Fradley et al. (2007) *J.Psychopharmacol.* **21** 384.
67. Collinson et al. (2000) *Eur.J.Neurosci.* **12** 171.
68. Wulff et al. (2007) *Nature Neurosci.* **10** 923.
69. Tenen and Hirsch (1980) *Nature* **288** 609.
70. Cowan et al. (1981) *Nature* **290** 54.
71. Barker et al. (1984) in *Actions and interactions of GABA and the benzodiazepines*. p203 Publ Raven Press.
72. Smiechen et al. (1993) in *Anxiolytic β -carbolines: from molecular biology to the clinic*. p7 Publ Springer Verlag.
73. Stephens et al. (1987) *Brain Res.Bull.* **19** 309.
74. Sternfield et al. (2004) *J.Med.Chem.* **47** 2176.
75. Harrison and Simmonds (1984) *Brain.Res.* **323** 287.
76. Cottrell et al. (1987) *Br.J.Pharmacol.* **90** 491.
77. Harrison et al. (1987) *J.Neurosci.* **7** 604.
78. Study and Barker (1981) *Proc.Natl.Acad.Sci.USA* **78** 7180.
79. Twyman et al. (1989) *Ann.Neurol.* **25** 213.
80. Callachan et al. (1987) *Neurosci.Letts.* **29** S21.
81. Nicoll et al. (1975) *Nature* **258** 625.
82. Simmonds (1981) *Br.J.Pharmacol* **73** 739.
83. Peters et al. (1988) *Br.J.Pharmacol.* **94** 1257.
84. Hosie et al. 2006 *Nature* **444** 486.
85. Belelli et al. (2006) *Neuroscience* **138** 821.
86. Brown et al. (2002) *Br.J.Pharmacol.* **136** 965.
87. Bianchi and MacDonald (2003) *J.Neurosci.* **23** 10934.
88. Herd et al. (2007) *Pharmacol.Ther.* **116** 20.
89. Koksmas et al. (2003) *J.Neurosci.* **23** 788.
90. Bonin and Orser (2008) *Pharmacol.Biochem.Behav.* **90** 105.
91. Nishikawa et al. (2002) *Neuropharmacology* **42** 337 .
92. Belelli et al. (1997) *Proc.Natl.Acad.Sci.USA* **94** 11031.
93. Jurd et al. (2003) *FASEB J.* **17** 250.
94. Cheng et al. (2006) *J.Neurosci.* **26** 3713.
95. Franks (2008) *Nature Neurosci.Rev.* **9** 370.
96. Zeller et al. (2007) *BMC Pharmacol.* **7** 2.
97. Mihalek et al. (2001) *Alcohol Clin.Exp.Res.* **25** 1708.
98. Sundstrom-Poromaa (2002) *Nature Neurosci.* **5** 721.
99. Wallner et al. (2003) *Proc.Natl.Acad.Sci.USA* **100** 15218.
100. Borghese et al. (2006) *J.Pharmacol.Exp.Ther.* **316** 1360.
101. Morrow et al. (2001) *Brain Res.Rev.* **37** 98.
102. Choi et al. (2008) *J.Neurosci.* **28** 11890.
103. Polenzani et al. (1991) *Proc.Natl.Acad.Sci.USA* **88** 4318.
104. Cutting et al. (1991) *Proc.Natl.Acad.Sci.USA* **88** 2673.
105. Enz and Cutting (1998) *Vision Res.* **38** 1431.
106. Enz and Cutting (1999) *Eur.J.Neurosci.* **11** 41.
107. Sieghart & Ernst (2005) *Curr.Med.Chem.Cent.Nerv.Sys. Agents* **5** 217.
108. Bowery et al. (1987) *Neuroscience* **20** 365.
109. Hill (1985) *Br.J.Pharmacol.* **84** 249.
110. Dunlap and Fishbach (1981) *J.Physiol.* **317** 519.
111. Newberry and Nicoll (1985) *J.Physiol.* **360** 161.
112. Bowery et al. (2002) *Pharmacol.Rev.* **54** 247.
113. Ong and Kerr (1990) *Life Sci.* **46** 1489.
114. Sakaba and Neher (2003) *Nature* **424** 775.
115. Luscher et al. (1997) *Neuron* **19** 687.
116. Kaupmann et al. (1997) *Nature* **368** 239.
117. Couve et (1998) *J.Biol.Chem.* **273** 26361.
118. Jones et al. (1998) *Nature* **396** 674.
119. Kaupmann et al. (1998) *Nature* **396** 683.
120. White et al. (1998) *Nature* **396** 679.
121. Bettler et al. (2004) *Physiol.Rev.* **84** 835.
122. Pin et al. (2007) *Pharmacol.Rev.* **59** 5.
123. Steiger et al. (2004) *J.Neurosci.* **24** 6115.
124. Blein (2004) *J.Biol.Chem.* **279** 48292.
125. Kniazeff et al. (2002) *J.Neurosci.* **22** 7352.
126. Galvez et al. (2000) *Mol.Pharmacol.* **57** 419.
127. Galvez et al. (2001) *EMBO J.* **20** 2152.
128. Robbins et al. (2001) *J.Neurosci.* **21** 8043.
129. Duthey et al. (2002) *J.Biol.Chem.* **277** 3236.
130. Urwyler et al. (2001) *Mol.Pharmacol.* **60** 963.
131. Dupuis et al. (2006) *Mol.Pharmacol.* **70** 2027.
132. Pin et al. (2005) *FEBS J.* **272** 2947.
133. Ullrich and Bettler (2007) *Curr.Opin.Neurobiol.* **17** 298.
134. Schuler et al. (2001) *Neuron* **31** 47.
135. Gassmann et al. (2004) *J.Neurosci.* **24** 6086.
136. Vigot et al. (2006) *Neuron* **50** 589.
137. Perez-Garci et al. (2006) *Neuron* **50** 603.
138. Waldmeier et al. (2008) *J.Neural Transm.* **115** 1401.
139. Guetg et al. (2009) *J.Neurosci.* **29** 1414.
140. Grace et al. (2004) *Proc.Natl.Acad.Sci.USA* **101** 12836.
141. Tiao et al. (2008) *J.Biol.Chem.* **283** 31005.
142. Bowery (2006) *Curr.Opin.Pharmacol.* **6** 37.
143. Penn (1989) *N.Engl.J.Med.* **320** 1517.
144. Ataka et al. (2000) *Pain* **86** 273.
145. Potes et al. (2006) *J.Neurosci.Res.* **83** 515.
146. Pan et al. (2008) *Pharmacol.Ther.* **117** 141.
147. Wu et al. (2007) *Neurobiol.Dis.* **26** 439.
148. Teichgräber et al. (2009) *Epilepsia* 2009 **50** 1697.
149. Froestl et al. (2004) *Biochem.Pharmacol.* **68** 1479.
150. Sunyer et al. (2009) *Hippocampus* **19** 90.
151. Roberts and Andrews (1997) *Psychopharmacology* **131** 271.
152. Smith et al. (2004) *Psychopharmacology* **173** 105.
153. Slattery et al. (2005) *Neuropsychopharmacology* **30** 2065.
154. Malherbe et al. (2008) *Br.J.Pharmacol.* **154** 797.
155. Heinmiller et al. (2009) *Behav.Neurosci.* **123** 145.
156. Matsuki et al. (2009) *Proc.Natl.Acad.Sci.USA* **106** 4459.
157. Alstermark et al. (2008) *J.Med.Chem.* **51** 4315.
158. Zhang et al. (2001) *TIPS* **22** 121.
159. Bowery and Enna (2000) *J.Pharmacol.Exp.Ther.* **292** 2.
160. Murata et al. (1996) *Bioorg.Med.Chem.Lett.* **6** 2071.
161. Ragozzino et al. (1996) *Mol.Pharmacol.* **50** 1024.
162. Sieghart (1995) *Pharmacol.Rev.* **47** 181.

GABA Receptor Compounds Available from Tocris

Cat. No.	Product Name	Primary Action
GABA_A Receptors		
Agonists		
0344	GABA	Endogenous agonist
0235	Isoguvacine hydrochloride	Selective GABA _A agonist
3250	L-838,417	Subtype-selective GABA _A partial agonist
0289	Muscimol	Potent GABA _A agonist
0181	TACA	GABA _A agonist. Also GABA-T substrate and GABA uptake inhibitor
0807	THIP hydrochloride	GABA _A agonist
0180	ZAPA sulfate	Agonist at 'low affinity' GABA _A receptor. More potent than GABA/muscimol
Antagonists		
0130	(+)-Bicuculline	Potent GABA _A antagonist
2503	(-)-Bicuculline methiodide	Water-soluble GABA _A antagonist
0109	(-)-Bicuculline methobromide	Water-soluble GABA _A antagonist
0131	(-)-Bicuculline methochloride	Water-soluble GABA _A antagonist
3109	Furosemide	GABA _A receptor antagonist. Also Na ⁺ /2Cl ⁻ /K ⁺ cotransporter blocker
1128	Picrotoxin	GABA _A receptor antagonist
2143	SCS	Selective GABA _A receptor antagonist; β1-subunit-selective
1262	SR 95531 hydrobromide	Selective, competitive GABA _A receptor antagonist
2905	TB 21007	α5-selective GABA _A inverse agonist
2745	U 93631	GABA _A receptor antagonist
Benzodiazepines		
3568	Bretazenil	Benzodiazepine partial agonist
0405	β-CCB	Benzodiazepine inverse agonist, putative endogenous ligand
2467	CGS 20625	Selective central benzodiazepine receptor partial agonist
0456	Chlormezanone	Skeletal muscle relaxant
1709	CL 218872	Benzodiazepine agonist
2805	Diazepam	Acts at the benzodiazepine modulatory site
0505	Dihydroergotoxine mesylate	Binds to GABA _A receptor Cl ⁻ channel; allosteric modulator of benzodiazepine site
3083	DMCM hydrochloride	Benzodiazepine inverse agonist
0554	FG 7142	Benzodiazepine inverse agonist
0658	FGIN-1-27	Potent, specific ligand for mitochondrial DBI receptor (TSPO)
0659	FGIN-1-43	Potent, specific ligand for mitochondrial DBI receptor
1328	Flumazenil	Benzodiazepine antagonist
0770	GBLD 345	High affinity benzodiazepine agonist
2174	Hispidulin	Partial positive allosteric modulator at the benzodiazepine receptor

Cat. No.	Product Name	Primary Action
1327	L-655,708	Selective for α5-containing GABA _A receptors
3087	Lorazepam	Acts at the benzodiazepine modulatory site
2832	Midazolam hydrochloride	Benzodiazepine agonist
0670	PK 11195	Antagonist at peripheral benzodiazepine receptors
1997	Ro 15-4513	Benzodiazepine partial inverse agonist
1995	Ro 19-4603	Benzodiazepine inverse agonist
3655	Zaleplon	Benzodiazepine agonist
1994	ZK 93423 hydrochloride	Potent benzodiazepine agonist
1996	ZK 93426 hydrochloride	Potent, competitive benzodiazepine antagonist
0655	Zolpidem	Benzodiazepine agonist
1094	Zopiclone	Benzodiazepine agonist
Modulators		
3653	Allopregnanolone	GABA _A receptor positive allosteric modulator
3046	Alphaxalone	Direct activator and potentiator of GABA _A
0881	Chlormethiazole hydrochloride	Potentiates GABA _A receptor function
1471	Etomidate	GABA mimetic and GABA modulatory agent
2531	Ganaxalone	Potent, positive allosteric modulator of GABA _A receptors
3597	Indiplon	Subtype-selective GABA _A receptor positive allosteric modulator
1295	Loreclezole hydrochloride	Subtype-selective GABA _A receptor modulator
2738	Org 20599	Positive allosteric modulator and direct agonist of GABA _A receptors
2681	7-PA	Antagonist of neurosteroid potentiation and direct gating of GABA _A
3652	Pregnanolone	GABA _A receptor positive allosteric modulator
0830	Primidone	Potentiates GABA _A receptor function
1512	SB 205384	GABA _A receptor modulator
3620	Topiramate	GABA _A receptor positive allosteric modulator. Also GluR5 antagonist
1558	Tracazolate hydrochloride	Subtype-selective GABA _A allosteric modulator
2734	U 89843A	Positive allosteric modulator of GABA _A receptors
2733	U 90042	GABA _A receptor ligand
3048	Valerenic acid	Positive allosteric modulator of GABA _A receptors
GABA_B Receptors		
Agonists		
0417	(RS)-Baclofen	Selective GABA _B agonist
0796	(R)-Baclofen	Active enantiomer of Cat. No. 0417
0344	GABA	Endogenous agonist
0379	SKF 97541	Extremely potent GABA _B agonist

Cat. No.	Product Name	Primary Action
Antagonists		
1245	CGP 35348	Brain penetrant, selective GABA _B antagonist
1247	CGP 46381	Brain penetrant, selective GABA _B antagonist
1246	CGP 52432	Potent, selective GABA _B antagonist
1088	CGP 54626 hydrochloride	Potent, selective GABA _B antagonist
1248	CGP 55845	Potent, selective GABA _B antagonist
0245	2-Hydroxysaclofen	Selective GABA _B antagonist, more potent than Cat. No. 0246
0178	Phaclofen	Selective GABA _B antagonist
0246	Saclofen	Selective GABA _B antagonist
0984	SCH 50911	Selective, competitive, orally active GABA _B
Modulators		
3313	rac BHFf	Potent, selective GABA _B positive allosteric modulator
1513	CGP 7930	Positive modulator at GABA _B receptors
1514	CGP 13501	Positive modulator at GABA _B receptors
2001	GS 39783	Positive modulator at GABA _B receptors
GABA_A (GABA_C) Receptors		
Agonists		
0344	GABA	Endogenous agonist
0289	Muscimol	Partial GABA _C agonist
0181	TACA	GABA _C agonist
Antagonists		
0379	SKF 97541	GABA _C antagonist. Also potent GABA _B agonist
0807	THIP hydrochloride	GABA _C antagonist
1040	TPMPA	Selective GABA _C antagonist
0180	ZAPA sulfate	GABA _C antagonist
GABA Transporters		
0206	β-Alanine	GABA uptake inhibitor (GAT-2 and -3). Also glycine receptor agonist
1296	CI 966	Selective inhibitor of GAT-1
0234	Guvacine	Specific GABA uptake inhibitor
0236	(±)-Nipecotic acid	GABA uptake inhibitor
2747	NNC 05-2090	GABA uptake inhibitor; moderately BGT-1 selective
1779	NNC 711	Selective inhibitor of GAT-1
0768	Riluzole	GABA uptake inhibitor. Also glutamate release inhibitor
1081	SKF 89976A	Potent GABA uptake inhibitor. Penetrates blood brain barrier
1561	(S)-SNAP 5114	GABA uptake inhibitor
0181	TACA	GABA uptake inhibitor. Also GABA _A agonist and substrate for GABA-T
Miscellaneous GABA		
1814	N-ArachidonylGABA	Inhibits pain <i>in vivo</i>
0806	Gabapentin	Anticonvulsant. Increases brain GABA

Cat. No.	Product Name	Primary Action
0538	<i>trans</i> -4-Hydroxycrotonic acid	GHB receptor ligand
0386	3-Methyl-GABA	Activator of GABA amino-transferase
1811	Modafinil	Psychostimulant
0780	NCS-382	Antagonist of γ-hydroxybutyric acid
2687	Pentylentetrazole	CNS stimulant
0939	Propofol	Potentiates GABA _A receptor-mediated inhibition and inhibits glutamate receptor-mediated excitation
3400	RuBi-GABA	Caged GABA; excitable by visible wavelength
2815	Valproic acid, sodium salt	Increases GABA levels; anticonvulsant
0808	Vigabatrin	GABA-T inhibitor
2625	Zonisamide	Anticonvulsant, modulates GABA neurotransmission

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

R&D SYSTEMS

NOVUS
BIOLOGICALS

TOCRIS

protein **simple**

ACD

biotechne®

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

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

