

Metabotropic Glutamate Receptors

Molecular Pharmacology

TOCRIS
bioscience

Francine C Acher

Laboratoire de Chimie et Biochimie Pharmacologiques et Toxicologiques, UMR8601-CNRS, Université René Descartes-Paris V, 45 rue des Saints-Pères, 75270 Paris Cedex 06, France. E-mail: Francine.Acher@univ-paris5.fr

Francine C Acher is currently a CNRS Research Director within the Biomedical Institute of University Paris-V (France). Her research focuses on structure/function studies and drug discovery using chemical tools (synthetic chemistry, molecular modeling), molecular biology and pharmacology within interdisciplinary collaborations.

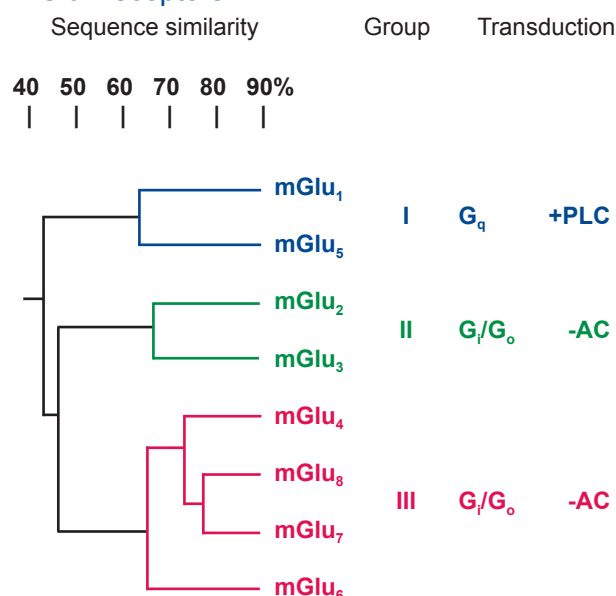
Introduction

Glutamate is the major excitatory amino acid transmitter in the brain. It is released from presynaptic vesicles and activates postsynaptic ligand-gated ion channel receptors (NMDA, AMPA and kainate receptors) to secure fast synaptic transmission.¹ Glutamate also activates metabotropic glutamate (mGlu) receptors, which modulate its release, postsynaptic response, as well as the activity of other synapses.^{2,3} Glutamate has been shown to be involved in many neuropathologies such as anxiety, pain, ischemia, Parkinson's disease, epilepsy and schizophrenia. Thus, because of their modulating properties, mGlu receptors are recognized as promising therapeutic targets.⁴ It is expected that drugs acting at mGlu receptors will regulate the glutamatergic system without affecting the normal synaptic transmission.

mGlu receptors are G-protein-coupled receptors (GPCRs). Eight subtypes have been identified and classified into three groups (I-III) based upon sequence homology, transduction mechanism and pharmacological profile (see Figure 1). Group I includes mGlu₁ and mGlu₅ receptors, which couple to G_q and activate phospholipase C (PLC). Group II (mGlu₂, mGlu₃) and group III (mGlu₄, mGlu₆, mGlu₇ and mGlu₈) receptors couple to G_i/G_o and inhibit adenylyl cyclase (AC). Group I receptors are mostly located postsynaptically, thus their activation increases excitability. On the other hand, group II/III receptors are generally presynaptic and their activation reduces glutamate release. Specific ligands have been found for each group and some of the subtypes, as described hereafter.^{5,6}

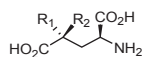
mGlu receptors belong to family 3 of the GPCR superfamily.⁷ Similar to all GPCRs, mGlu receptors contain a heptahelical domain (HD) in the membrane region. In addition, like all members of family 3, mGlu receptors are characterized by a large extracellular amino terminal domain (ATD) where

Figure 1 | Classification of the 8 Subtypes of mGlu Receptors

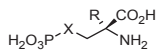


the glutamate binding site is found (see Figure 3). This domain adopts a bilobate structure similar to LIVBP (Leucine Isoleucine Valine Binding Protein), a bacterial periplasmic protein involved in the transport of hydrophobic amino acids;⁸⁻¹⁰ these amino acids bind to an open conformation of the protein, which closes subsequently to trap them in between the two lobes. A similar binding mode has been proposed for glutamate and competitive agonists in the LIVBP domain (LIVBPD) of mGlu receptors. Moreover, it was shown that the closed conformation of this domain is required for receptor activation.¹¹ Examination of the glutamate binding site in the eight mGlu receptor subtype crystal structures (mGlu₁)¹⁰ or homology models¹²⁻¹⁶ reveals a common binding motif for the α -amino and α -carboxylic functions of glutamate,¹⁷ while residues that bind the distal γ -carboxylate vary

Figure 2 | Competitive mGlu Receptor Ligand Structures

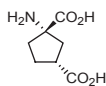


$R_1, R_2 = H$ **L-Glutamic acid**
 $R_1 = H, R_2 = CH_2CH(Ph)_2$ **ADED**
 (LY 310225)

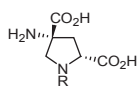


$X = CH_2, R = H$ **(S)-AP4**
 $X = O, R = H$ **(S)-SOP**
 $X = CH_2, R = Me$ **MAP4**

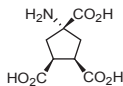
Ac-Asp-Glu
NAAG



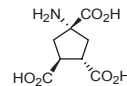
(1S,3R)-ACPD



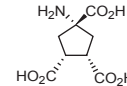
$R = H$ **(2R,4R)-APDC**
 $R = CH_2Ph$ **BnAPDC**
 $R = CH_2Naphthyl$ **NM-APDC**
 $R = CO_2H$ **APTC**



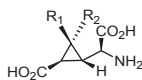
ACPT-I



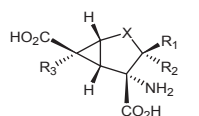
(+)-ACPT-III



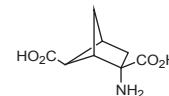
ACPT-II



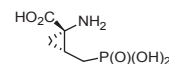
$R_1, R_2 = H$ **L-CCG-I**
 $R_1 = H, R_2 = 3'-CH_3$ **3'Me-CCG**
 $R_1 = H, R_2 = 3'-CH_2OH$ **3'HM-CCG**
 $R_1 = H, R_2 = 3'-CH_2SH$ **3'SM-CCG**
 $R_1 = H, R_2 = CO_2H$ **DCG IV**
 $R_1 = H, R_2 = 9'$ xanthylethyl **XE-CCG-I**



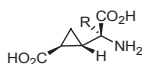
$X = CH_2, R_1, R_2, R_3 = H$ **LY 354740**
 $X = O, R_1, R_2, R_3 = H$ **LY 379268**
 $X = S, R_1, R_2, R_3 = H$ **LY 389795**
 $X = CH_2, R_1, R_3 = H, R_2 = F$ **MGS0008**
 $X = CH_2, R_1, R_2 = H, R_3 = F$ **LY 354740-6F**
 $X = CO, R_1, R_2 = H, R_3 = F$ **MGS0028**
 $X = CH_2, R_2, R_3 = H, R_1 = OH$ **HYDIA**
 $X = CH_2, R_1 = 3,4-Cl_2PhCH_2O$
 $R_2 = H, R_3 = F$ **MGS0039**



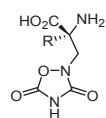
ABHxD-I



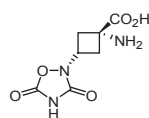
(1S,2R)-APCPr



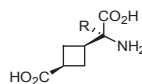
$R = Me$ **MCCG-I**
 $R = 9'$ -xanthylmethyl **LY 341495**
 $R = (3-ClPh)_2Et$ **mCD-CCG**



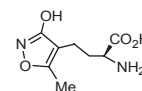
$R = H$ **Quis**
 $R = benzyl$ **BnQuis**



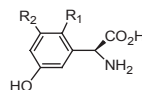
Z-CBQA



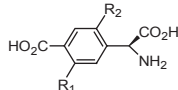
$R = 9'$ -thioxanthylmethyl
LY 393675



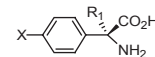
(S)-HomoAMPA



$R_1 = H, R_2 = OH$ **(S)-3,5-DHPG**
 $R_1 = Cl, R_2 = H$ **CHPG**



$R_1, R_2 = H$ **(S)-4CPG**
 $R_1 = CO_2H, R_2 = H$ **(S)-3,4-DCPG**
 $R_1 = H, R_2 = Me$ **4C2MPG**
(LY 367385)
 $R_1 = OH, R_2 = Me$ **4C3H2MPG**
(LY 339840)



$X = CO_2H, R_1 = Me$ **(S)-MCPG**
 $X = CO_2H, R_1 = 9'$ -thioxanthylmethyl
LY 367366
 $X = PO_3H_2, R_1 = H$ **(S)-PPG**
 $X = PO_3H_2, R_1 = Me$ **MPPG**
 $X = PO_3H_2, R_1 = cyclopropyl$ **CPPG**

Agonists are shown in turquoise
 Antagonists are shown in dark blue

(Bold Text Denotes Compounds Available From Toctris)

from one subtype to another.¹⁴ Thus, not surprisingly, all competitive agonists are α -amino acids, bearing various selective functional groups on their side chain⁶ (see Figure 2). The first generation of orthosteric ligands was followed by a second generation of allosteric modulators that bind in the HD.¹⁸ The first molecule described as a non-competitive mGlu antagonist was CPCCOEt in the late nineties.¹⁹ Since then, numerous allosteric modulators have been discovered by high-throughput screening (HTS) in pharmaceutical companies.²⁰⁻²²

The purpose of the present article is to review our actual knowledge of the pharmacology of mGlu receptors. Several detailed reviews^{2,3,5,6} have been published; thus only the most potent and selective known ligands will be presented and emphasis will be placed on compounds that were recently disclosed.

Competitive Ligands

An α -amino acid moiety can be found in all mGlu receptor competitive ligands (agonists and antagonists) and most of the side chains hold an acidic function. In the ligand active conformations, the spatial disposition of these functional groups is that of glutamate in an extended conformation, as predicted by pharmacophore²³ and homology models.¹⁴ For many years these compounds have

are already quite hydrophilic²⁵ and few side effects are predicted. Other glutamate analogs were also shown to be systemically active: (2*R*,4*R*)-APDC, (*S*)-DCPG, 3'-Me-CCG, 3'-HM-CCG and ACPT-I (Figure 2). Desensitization was also feared with continuous activation in the case of group II/III receptors, yet little was observed after several days of agonist activation. Altogether these results promote a renewed interest in mGlu receptor competitive ligands.

Agonists (Table 1 and Figure 2)

The first agonist that was able to discriminate between ionotropic and metabotropic glutamate receptors was *trans*-ACPD (1*S*,3*R* isomer).²⁶ The ligand contributed considerably to the study of metabotropic glutamate receptors despite its lack of subtype selectivity.^{2,3,5} A limited number of molecules possess agonist activity across all mGlu receptors. The endogenous agonist L-glutamate, L-CCG-I and ABHxD-I are the most potent.^{2,3,5} It can be noted that L-CCG-I and ABHxD-I are conformationally constrained and mimic the bioactive extended glutamate conformation.²³ Selectivity can be gained by adding new chemical groups onto these structures.

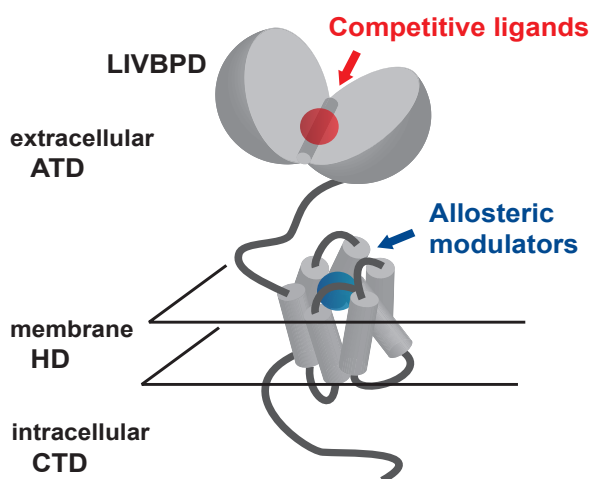
Group I

Quisqualate (Quis) is the most potent group I agonist. However, it also activates AMPA receptors, thus its use is restricted. The most popular group I selective agonist is (*S*)-3,5-DHPG, yet it exhibits only moderate potency.^{2,3,5} CHPG²⁷ and *Z/E*-CBQA²⁸ have been claimed to specifically activate mGlu₅ receptors, although the affinity of the former is quite low. To date, no specific mGlu₁ competitive agonists have been disclosed.

Group II

LY 354740 was the first mGlu agonist reported to exhibit a nanomolar affinity.²⁴ It is group II selective, as are its oxy (LY 379268) and thia (LY 389795) derivatives.²⁹ The introduction of a fluorine atom at position 3 (MGS0008) or 6 (LY 354740-6F) retained the potent activity, which was even enhanced when a carbonyl group was added, as in the case of MGS0028.³⁰ This series of bicyclic glutamate analogs was derived from the general agonist L-CCG-I, where increased potency and group II selectivity was gained through the second hydrocarbon ring. However, it was recently shown that a methyl or hydroxymethyl substituent in the 3' position (3'-Me-CCG and 3'-HM-CCG) provided agonists with similar potency.^{31,32} Replacement of the hydroxyl functionality at C3' of 3'-HM-CCG, by a sulfhydryl results in decreased affinity at mGlu_{2/3}. Interestingly, this analog (3'-SM-CCG) remains an mGlu₂ agonist but is a full antagonist at mGlu₃.³³ A similar selectivity was also reported for the C4 β -methyl-substituted analog of LY 354740³⁴. These two compounds selectively activate mGlu₂, while NAAG is the only

Figure 3 | Schematic Representation of an mGlu Receptor: the Two Orthosteric and Allosteric Binding Sites are Indicated



been considered as valuable research tools, but not as drug candidates, because of their poor LogP, related to their highly polar chemical structures. Jim Monn and colleagues from Eli Lilly were the first to show that such a glutamate analog, LY 354740, was able to pass the biological barriers and thus be orally active as an anticonvulsant and anxiolytic.²⁴ Moreover, such drugs are poorly metabolized as they

reported mGlu₃ competitive agonist to discriminate between the two group II subtypes. Other group II selective agonists have been described with submicromolar affinity: (2*R*,4*R*)-APDC and DCG IV.

Group III

Most of the potent group III selective agonists bear a diacidic side chain, which can interact with the highly

basic distal binding pocket.^{14,16} (S)-AP4 (L-AP4), (S)-SOP(L-SOP) and (1*S*,2*R*)-APCPr^{35,36} are the most potent, displaying submicromolar affinities at cloned receptors, except for mGlu₇, to which all bind with weak affinity. (S)-PPG,^{37,38} (S)-3,4-DCPG,³⁹ ACPT-I and (+)-ACPT-III⁴⁰ have also been described as micromolar agonists. Interestingly, a CCG derivative bearing a hydroxymethyl group in the 3' position

Table 1 | Potencies of Selective and Non-selective mGlu Receptor Agonists^a

Receptor		Group I		Group II		Group III			
		mGlu ₁	mGlu ₅	mGlu ₂	mGlu ₃	mGlu ₄	mGlu ₆	mGlu ₇	mGlu ₈
Non-selective agonists	L-Glu^{c,d}	1-13	3-11	0.3-12	2-9	3-17	5-38	2300	8-10
	L-CCG-I^{c,d}	2	3	0.5	0.4	9	6	230	3
	ABHxD-I ^{c,d}	2	0.7	0.3	2	23	5	–	–
	(1<i>S</i>,3<i>R</i>)-ACPD^{c,d}	5-80	5-40	7-18	6-17	100-1000	300	n.e.	45-166
Group I subtype-selective agonists	Quis^{c,d}	0.03-3	0.02-0.3	100-1000	40-220	100-1000	n.e.	n.e.	720
	(S)-3,5-DHPG^{c,d}	6	2	n.e.	n.e.	n.e.	–	n.e.	n.e.
	CHPG^c	> 10000	750	–	–	–	–	–	–
	Z-CBQA ^c	> 1000	11	> 100	–	> 100	–	–	–
Group II subtype-selective agonists	LY 354740^{b,c}	> 100	> 100	0.01	0.04	> 100	3	> 100	12
	LY 379268^{b,c}	> 100	> 100	0.003	0.005	21	0.4	> 100	2
	LY 389795 ^{b,c}	> 100	> 100	0.004	0.008	> 100	2	> 100	7
	MGS0008 ^e	> 100	> 100	0.029	0.049	> 100	> 100	> 100	–
	LY 354740-6F ^e	> 100	> 100	0.017	0.081	> 100	> 100	> 100	–
	MGS0028 ^e	> 100	> 100	0.0006	0.0021	> 100	> 100	> 100	–
	3'Me-CCG ^f	> 100	> 100	0.008	0.038	> 100	1.198	> 100	1.32
	(+)-3'HM-CCG ^g	> 100	> 100	0.004	0.007	1.8	0.147	> 100	0.010
	2<i>R</i>,4<i>R</i>-APDC^{b,c}	> 100	> 100	0.4	0.4	> 300	110	> 300	> 100
	DCG IV^{c,d}	ant.	ant.	0.1-0.4	0.1-0.2	ant.	ant.	ant.	ant.
	NAAG^{c,d}	> 300	> 300	134-1000	10-65	> 300	> 300	–	–
Group III subtype-selective agonists	(S)-AP4^{c,d}	> 1000	> 1000	> 1000	> 1000	0.2-1.2	0.6-0.9	160-500	0.06-0.9
	(S)-SOP^{c,d}	n.e.	n.e.	ant.	ant.	1-4	3	160-1200	2
	(1 <i>S</i> ,2 <i>R</i>)-APCPr ^h	–	–	–	–	0.6	1.9	602	0.3
	ACPT-I^{c,d,i}	ant.	–	n.e.	–	6.5	18.4	–	10.1
	(+)-ACPT-III ^{c,d,i}	–	–	–	–	8.8	19.2	–	7.0
	(S)-PPG ^{b,j}	> 500	> 500	> 300	> 200	3.2 (5.2)	(4.7)	48 (185)	(0.21)
	(S)-HomoAMPA ^c	> 1000	> 1000	> 1000	–	> 1000	58	> 5000	–
	BnAPDC ^c	> 1000	ant.	ant.	> 100	> 300	20	–	> 300
	(S)-3,4-DCPG^{b,k}	ant.	> 100	> 100	> 100	8.8	3.6	> 100	0.031

(Bold Text Denotes Compounds Available From Toctris)

^aEC₅₀ or K_d values (μM) measured with rat or human (when indicated^b) cloned receptors. ant. = antagonist; n.e. = no effect. References for agonist potencies that have been cited in reviews ⁵ and/or ⁶ are referred to as such. ^bEC₅₀ or K_d values (μM) obtained with human mGlu receptors.

^cSchoepp *et al* (1999)⁵ ^dPin *et al* (1999)⁶ ^eNakazato *et al* (2000)³⁰ ^fCollado *et al* (2002)³¹ ^gCollado *et al* (2004)³² ^hKroona *et al* (1991)³⁵ ⁱSibille *et al* (unpublished results) ^jBessis *et al* (unpublished results) ^kGasparini *et al* (1999)³⁷ and (2000);³⁸ data in parentheses refer to (±)-PPG³⁷ ^lThomas *et al* (2001)³⁹

(3'HM-CCG) displays similar affinity for mGlu₈ and mGlu_{2/3} receptors.³² Again, very few compounds are subtype-selective: *N*-benzyl-APDC (BnAPDC)⁴¹ and (*S*)-homoAMPA⁴² at mGlu₆; (*S*)-3,4-DCPG at mGlu₈, with an EC₅₀ over two orders of magnitude lower than at other group III receptors.³⁹ Very recently Faust Pharmaceuticals discovered a selective mGlu₄ agonist (FP429) in the APTC family.⁴³

Antagonists (Table 2 and Figure 2)

Most of the competitive antagonists prevent the complete closing of the two lobes of the LIVBP domain. Substitution of the α -proton of glutamate analogs by a methyl group as in MCCG, MCPG and MAP4, or a bulkier group as in LY 341495, turns the corresponding agonists (4CPG, AP4 and L-CCG-I) into antagonists. However, agonist properties can be recovered when residues responsible for the hindrance are mutated.¹¹ Closing can also be disturbed by ionic repulsion as in the case of ACPT-II.¹¹

Group I

The first generation of group I mGlu antagonists was composed of 4-carboxyphenylglycine derivatives, such as (*S*)-MCPG, which has been widely used. Its

affinity was improved when the α -methyl group was changed to α -thioxanthylmethyl, as in LY 367366, but this derivative is also able to antagonize group II/III receptor activation.⁵ The highest potency was then found with α -substituted 3-carboxycyclobutylglycines such as LY 393675 (*cis* isomer) and its *trans* isomer⁵ or a *cis/trans* mixture (LY 393053).⁴⁴ This latter mixture was shown to be systemically active and inhibit both mGlu₁ and mGlu₅, as well as activate other group II/III mGlu receptors.⁴⁴ Although slightly less potent, LY 367385 (4C2MPG) and LY 339840 (4C3H2MPG) display subtype 1 selectivity.⁴⁵ However, LY 367385 was also shown to inhibit the cystine/glutamate exchanger.⁴⁶ No mGlu₅ selective and competitive antagonists have been described.

Group II

As most potent group II agonists are derived from L-CCG-I, the most potent group II antagonists are obtained when aryl substituents are introduced in specific positions of that glutamate analog. Thus

Table 2 | Potencies of Selective and Non-selective mGlu Receptor Competitive Antagonists^a

Receptor		Group I		Group II		Group III			
		mGlu ₁	mGlu ₅	mGlu ₂	mGlu ₃	mGlu ₄	mGlu ₆	mGlu ₇	mGlu ₈
Non-selective antagonists	LY 341495^{b,c,i}	6.8-9.7	8.2	0.021	0.014	2.6-22	1.1-1.8	0.99	0.17
	LY 393053 ^{b,e}	1.0	1.6	3.0	–	> 100	–	20	3.0
	ACPT-II^d	115	–	88	–	77	–	–	123
Group I subtype-selective antagonists	LY 367385^{b,f}	8.8	> 300	> 300	–	> 300	–	–	–
	LY 367366 ^{b,c}	6.6	5.6	–	–	–	–	–	–
	LY 339840 ^{b,f}	7.5	140	> 300	–	> 300	–	–	–
	(S)-MCPG^{c,d}	40-320	195-460	15-340	300-1000	> 1000	> 100	> 1000	> 300
Group II subtype-selective antagonists	ADED ^{b,c}	> 300	> 300	18	6.1	> 300	–	> 300	> 300
	(<i>S</i>)-BnQuis ^{b,c}	300	300	7.1	–	n.e.	n.e.	–	–
	mCD-CCG ^g	43	49	0.007	0.010	–	–	–	1.8
	HYDIA ^h	> 100	> 100	0.105	0.102	22	–	–	15
	MSG0039 ⁱ	> 100	–	0.020	0.024	1.7	2.1	–	–
	NMAPDC ^{b,c}	> 300	> 300	20	8.6	> 300	–	–	> 300
	XE-CCG ^{b,j}	–	–	0.20	0.075	–	–	–	–
Group III subtype-selective antagonists	DCG IV^d	390	630	ago.	ago.	22	40	25-40	15-32
	MAP4^{c,d}	n.e.	–	500	–	90-190	–	–	25-105
	CPPG^{b,c,k}	–	–	–	–	12	4	17	11
	MPPG^{c,d}	> 1000	n.e.	11-320	–	54-110	480	300	20-50

(Bold Text Denotes Compounds Available From Tocris)

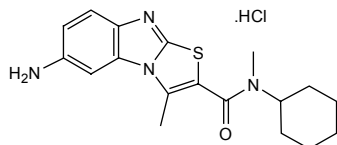
^aIC₅₀ or K_b values (μ M) measured with rat or human (when indicated^b) cloned receptors. ago. = agonist; n.e. = no effect. References for antagonist potencies that have been cited in reviews ⁵ and/or ⁶ are referred to as such.

^bIC₅₀ or K_b values (μ M) obtained with human mGlu receptors.

^cSchoepp *et al* (1999)⁵ ^dPin *et al* (1999)⁶ ^eChen *et al* (2000)⁴⁴ ^fKingston *et al* (2002)⁴⁵ ^gSørensen *et al* (2003)⁴⁸ ^hAdam *et al* (1999)⁵¹ ⁱChaki *et al* (2004)⁴⁹ ^jPellicciari *et al* (2001)⁴⁷ ^kConway *et al* (2001);⁵² Naples and Hampson (2001);¹¹⁷ Wright *et al* (2000)⁵⁴

YM 298198, Selective mGlu₁ Antagonist

YM 298198
Cat. No. 2448



YM 298198 is a newly characterized, non-competitive antagonist with high affinity and selectivity for mGlu₁ receptors ($K_i = 19$ nM). The compound is inactive at other mGlu receptor subtypes (mGlu₂₋₇), ionotropic receptors and glutamate transporters at concentrations up to 10 μ M. YM 298198 inhibits glutamate-induced IP production more potently than CPCCOEt (IC_{50} values are 16 nM and 6.3 μ M respectively), and is orally active *in vivo*, demonstrating an antinociceptive effect in hyperalgesic mice.

Kohara et al (2005) Radioligand binding properties and pharmacological characterization of 6-Amino-N-cyclohexyl-N,3-dimethyl-thiazolo[3,2-a]-benzimidazole-2-carboxamide (YM-298198), a high-affinity, selective and noncompetitive antagonist of metabotropic glutamate receptor type 1. *J.Pharmacol.Exp.Ther.* **315** 163.

(Sold with the permission of Astellas Pharma Inc.)

LY 341495⁵ and XE-CCG⁴⁷, containing a 9'-xanthylmethyl or 9'-xanthylethyl moiety in the α - or 3'-position, display nanomolar affinities. Potency is retained when the α -xanthyl moiety is replaced by two substituted phenyl groups (e.g. mCD-CCG).⁴⁸ It was recently reported that MGS0039 also exhibited high competitive group II antagonist activity.^{49,50} The addition of a dichlorobenzyl group to a close analog of HYDIA⁵¹ notably increases its affinity. MGS0039 and HYDIA are derivatives of the well-known agonist LY 354740. Indeed, it was previously demonstrated that substitution at the 3-position of the bicyclohexane is critical for agonist/antagonist properties.¹⁴ Systemic and antidepressant-like effects were observed with both LY 341495 and MGS0039.⁴⁹ Other arylalkyl-substituted glutamate analogs such as ADED (LY 310225), (S)-BnQuis and NM-APDC display group II selectivity with IC_{50} values in the micromolar range.⁵

Group III

Highly potent and group III-selective competitive antagonists have not yet been reported. The best agonist (S)-AP4 becomes a moderate antagonist when its α -proton is substituted by a methyl group, in MAP4. MCPG, a weak group I/II antagonist, becomes a moderate group III antagonist when the 4-carboxylate is replaced by a phosphonate, in MPPG. Addition of a substituent in the 3-position leads to a similar group III antagonist activity but increases selectivity for group III over group II.⁵² CPPG, the analog of MPPG bearing an α -cyclopropyl group, exhibits slightly increased potency^{5,52} in the same range as DCG IV, which is also a group II agonist.⁵³ Thus, the best activity is found with the non-selective antagonist LY 341495.⁵⁴

Allosteric Modulators

Allosteric modulators are non-competitive ligands that bind in the transmembrane heptahelical domain (HD). Negative and positive modulators have been identified.²⁰⁻²² Negative modulators inhibit receptor activation without affecting agonist binding, while positive modulators enhance agonist activation but do not activate receptors alone. Among the numerous mGlu receptor modulators that have been described (mostly in patents), only those for which biological activities are available will be presented here. These compounds are generally highly potent and subtype-selective, which is not the case for most competitive ligands.

Group I (Figure 4)

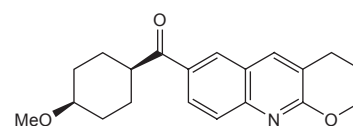
Both non-competitive inhibitors and enhancers have been disclosed for group I receptors.

mGlu₁ Antagonists

Detailed studies have been devoted to CPCCOEt, the first negative mGlu receptor modulator.^{19,55,56} In particular, specific residues of the HD that bind to CPCCOEt were identified by a group from Novartis.¹⁹ Following this, other compounds with higher affinities were discovered by HTS and subsequent optimization, in various companies: NPS2390^{57,58} (NPS Pharma Inc.), BAY 36-7620⁵⁹ (Bayer AG), LY 456066^{60,61} (Eli Lilly), R214127⁵⁸/JNJ 16259685^{62,63} (Johnson & Johnson), 3,5-dimethyl-pyrrole-2,4-dicarboxylic acid diesters of which DM-PPP is the most potent derivative⁶⁴ (GlaxoSmithKline), several analogs of

JNJ 16259685, Highly Potent mGlu₁ Antagonist

JNJ 16259685
Cat. No. 2333



JNJ 16259685 is a non-competitive mGlu₁ antagonist ($K_i = 0.34$ nM) that exhibits low nanomolar potency. It inhibits glutamate-induced Ca^{2+} mobilisation at the human mGlu₁ receptor with an IC_{50} value of 1.21 nM and is approximately 6000 times more potent than CPCCOEt and 50 times more potent than BAY 36-7620. The antagonist is selective over mGlu₅ (> 400-fold) and displays no activity at mGlu₂, mGlu₃, mGlu₄, mGlu₆, AMPA or NMDA receptors ($IC_{50} > 10$ μ M). JNJ 16259685 is centrally active following systemic administration.

Lavreysen et al (2004) JNJ16259685, a highly potent, selective and systemically active mGlu1 receptor antagonist. *Neuropharmacology* **47** 961.
Mabire et al (2005) Synthesis, structure-activity relationship, and receptor pharmacology of a new series of quinoline derivatives acting as selective, noncompetitive mGlu1 antagonists. *J.Med.Chem.* **48** 2134.

(Sold under license from Johnson & Johnson
Pharmaceutical Research & Development, a division
of Janssen Pharmaceutica NV)

EM-TBPC^{65,66} (Hoffmann-La Roche) and recently YM 298198⁶⁷ (Yamanouchi Pharma) and triazafluorenones⁶⁸ (Abbott Laboratories). A homology model of the mGlu₁ allosteric binding site has been generated and a binding mode proposed for EM-TBPC, which has been validated by mutagenesis and functional assays.⁶⁵ Additionally, it was shown that several inhibitors (R214127, CPCCOEt, NPS2390 and BAY 36-7620) bind to this same site.⁵⁸ Promising anxiolytic and analgesic effects have been reported with allosteric mGlu₁ receptor antagonists. However, cognition impairment was also found with JNJ 16259685.⁶⁹

mGlu₁ Positive Modulators

The first allosteric potentiators of rat mGlu₁ receptors to be disclosed were Ro 01-6128, Ro 67-4853⁷⁰ and Ro 67-7476.^{71,72} Chimeric and mutated receptors were constructed to confirm the transmembrane localization of the binding site of these ligands, which are subtype 1 selective.⁷¹ Interestingly, Ro 67-7476 and Ro 01-6128 have little or no effect on human mGlu₁ receptor activation while Ro 67-4853 produces a pronounced enhancement.⁷¹

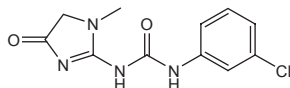
mGlu₅ Antagonists

SIB 1757 and SIB 1893⁷³ were initially found and optimized into MPEP,⁷⁴ which has been widely used⁷⁵ to explore the physiological roles of mGlu₅ receptors as a potential therapeutic target. Further investigations led to a methoxy derivative M-MPEP that can easily be radiolabeled.⁷⁶ More recently MTEP, a pyridine derivative of MPEP with improved aqueous solubility, was described with similar high mGlu₅ affinity⁷⁷ as well as its radiolabeled methoxymethyl derivative MM-MTEP,^{78,79} M-PEPy⁷⁸ and bipyridyl derivative MTEB.⁸⁰ The Merck group also disclosed new families

Fenobam, Selective mGlu₅ Antagonist

Fenobam

Cat. No. 2386



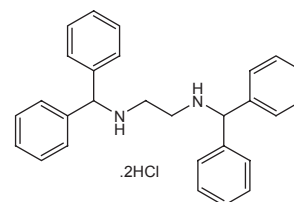
Fenobam is a potent and selective non-competitive mGlu₅ antagonist that displays inverse agonist activity. It blocks mGlu₅ constitutive activity *in vitro* with an IC₅₀ value of 87 nM. Fenobam acts at an allosteric modulatory site shared with MPEP and binds the receptor with K_d values of 54 and 31 nM for rat and human receptors respectively. The compound displays anxiolytic activity following oral administration *in vivo*.

Porter *et al* (2005) Description of a clinically validated anxiolytic with mGlu₅ antagonist properties. *Neuropharmacology* **49** (Suppl. 1) 267. Porter *et al* (2005) Fenobam: a clinically validated nonbenzodiazepine anxiolytic is a potent, selective, and noncompetitive mGlu₅ receptor antagonist with inverse agonist activity. *J.Pharmacol.Exp.Ther.* **315** 711.

AMN 082, The First mGlu₇-Selective Agonist

AMN 082

Cat. No. 2385



AMN 082 is the first selective mGlu₇ agonist. It potently inhibits cAMP accumulation and stimulates GTPγS binding in recombinant cells and on membranes expressing mGlu₇ (EC₅₀ = 64 - 290 nM). AMN 082 is selective over other mGluR subtypes and selected ionotropic glutamate receptors up to 10 μM. The agonist acts via a novel allosteric site and is orally active and brain penetrant.

Flor *et al* (2005) AMN082, the first selective mGluR7 agonist: activation of receptor signaling via an allosteric site in the transmembrane domain modulates stress parameters *in vivo*. *Neuropharmacology* **49** (Suppl. 1) 244. Mitsukawa *et al* (2005) A selective metabotropic glutamate receptor 7 agonist: Activation of receptor signaling via an allosteric site modulates stress parameters *in vivo*. *Proc.Natl.Acad.Sci.USA* **102** 18712.

of potent mGlu₅ antagonists: aryl benzoxazoles⁸¹ (illustrated by BOMA), heteroarylazoles⁸² (such as tetrazole PTeB⁸² and its derivatives,^{83,84} dipyrindyl amides⁸⁵ and dipyrindyl amines⁸⁶). Recently it was demonstrated that the known anxiolytic drug fenobam was, in fact, a potent non-competitive mGlu₅ antagonist.⁸⁷ MPEP and derivatives were shown to exhibit anxiolytic effects in animal models.^{75,82} Additionally, molecular determinants of the high affinity binding site of MPEP have been defined,⁸⁸ a striking similarity with critical residues of the mGlu₁ binding site was observed.

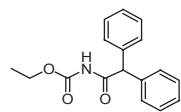
mGlu₅ Positive Modulators

Two mGlu₅ potentiators, DFB⁸⁹ and CPPHA,⁹⁰ were first identified by the Merck group. It should be noted that replacing the fluorine atoms of DFB with methoxy substituents turns this ligand into an antagonist (DMeOB),⁸⁹ while dichlorobenzaldazine (DCB) is a neutral modulator that attenuates the potentiation conferred by DFB.⁸⁹ Similar modulations were found with close analogs of MPEP; moving the methyl substituent of the MPEP pyridyl ring to the neighbouring carbon turns this analog (5MPEP) into a neutral modulator.⁹¹ Thus, it is suggested that the positive allosteric binding site overlaps the negative one.

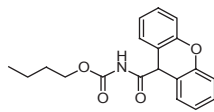
New series of mGlu₅ positive modulators (diphenylpyrazolebenzamide CDPPB^{92,93} and oxadiazol ADX47273^{94,95}) demonstrated anti-psychotic effects in animal models of schizophrenia.

Figure 4 | Group I Allosteric Modulator Structures and Potencies

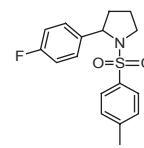
A. mGlu₁ Receptor Potentiators



Ro 01-6128
EC₅₀ = 0.2 μM

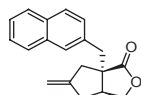


Ro 67-4853
EC₅₀ = 0.07 μM

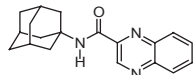


Ro 67-7476
EC₅₀ = 0.2 μM

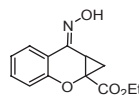
B. mGlu₁ Receptor Antagonists



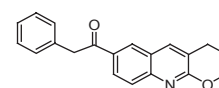
BAY 36-7620
IC₅₀ = 160 nM
K_i = 11.2 nM



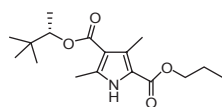
NPS2390
IC₅₀ = 5.2 nM (chimera CaSR/mGlu₁)
K_i = 1.4 nM



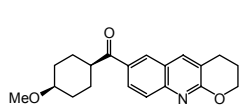
CPCCOEt
IC₅₀ = 10.3 μM
K_i = 4.9 μM
(- isomer active)



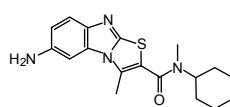
R214127
IC₅₀ = 21.6 nM
K_i = 1.4 nM



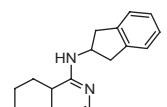
DM-PPP
IC₅₀ = 16 nM



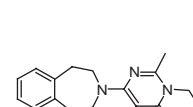
JNJ 16259685
IC₅₀ = 3.24 nM
K_i = 0.34 nM



YM 298198
IC₅₀ = 16 nM
K_i = 19 nM

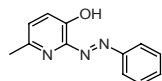


LY 456066
IC₅₀ = 12 nM (hmGlu₁)

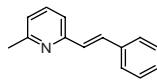


EM-TBPC
K_i = 11 nM

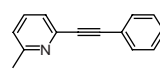
C. mGlu₅ Receptor Antagonists



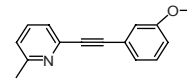
SIB 1757
IC₅₀ = 0.37 μM (hmGlu₅)



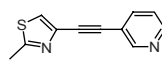
SIB 1893
IC₅₀ = 0.29 μM (hmGlu₅)



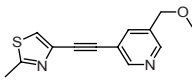
MPEP
IC₅₀ = 36 nM (hmGlu₅)



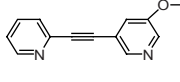
M-MPEP
IC₅₀ = 10 nM (hmGlu₅)



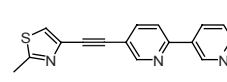
MTEP
IC₅₀ = 5 nM (hmGlu₅)



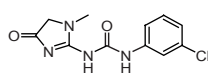
MM-MTEP
IC₅₀ = 7 nM (hmGlu₅)



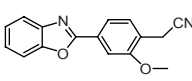
M-PEPy
IC₅₀ = 1 nM (hmGlu₅)



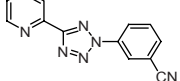
MTEB
IC₅₀ = 2 nM (hmGlu₅)



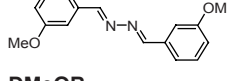
Fenobam
IC₅₀ = 58 nM



BOMA
IC₅₀ = 3 nM

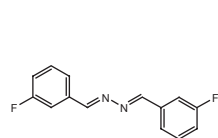


PTeB
IC₅₀ = 0.07 μM (hmGlu₅)

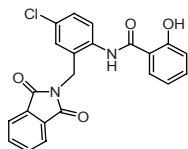


DMeOB
IC₅₀ = 3 μM (mGlu₅)
IC₅₀ = 35 μM (mGlu₄)
IC₅₀ = 50 μM (mGlu₃)

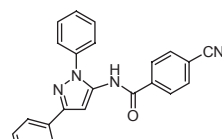
D. mGlu₅ Receptor Potentiators



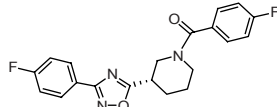
DFB
EC₅₀ = 2.4 μM (mGlu₅)



CPPHA
EC₅₀ = 0.14 μM (mGlu₅)

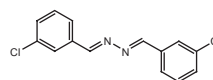


CDPBB
EC₅₀ = 27 nM (hmGlu₅)

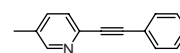


ADX47273

E. mGlu₅ Receptor Neutral Modulators



DCB
IC₅₀ = 2.6 μM for DFB potentiation attenuation



5MPEP
EC₅₀ = 2.32 μM for blocking MPEP inhibition
IC₅₀ = 1.71 μM for DFB and CDPBB potentiation attenuation

(Bold Text Denotes Compounds Available From Toctris)

Group II (Figures 5A and 5B)**mGlu₂ Antagonists**

Heterocyclic enol ethers such as Ro 64-5229 were reported as the first selective non-competitive mGlu₂ receptor antagonists.^{96,97} A series of dihydrobenzo[b][1,4]diazepin-2-one derivatives, typified by CH-DBO, was later claimed also to contain mGlu₂ receptor antagonists with nanomolar affinities.⁹⁸

mGlu₂ Positive Modulators

LY 487379, a pyridylmethylsulfonamide, has been reported to potentiate the activity of glutamate at mGlu₂ receptors with an EC₅₀ of 0.3 μM and be highly selective for this subtype.⁹⁹ It was also demonstrated that LY 487379 binds to a pocket in the transmembrane domain that is different from the orthosteric site in the ATD.⁹⁹ Further SAR studies led to the discovery of the 1-methylbutoxy analog (2,2,2-TEMPS) with improved potency (EC₅₀ = 14 nM) and selectivity.^{100,101} Recently, a new chemical series of phenyl-tetrazolyl acetophenones (e.g. PTBE) was disclosed as selective mGlu₂ potentiators,¹⁰² followed by extensive SAR studies.¹⁰³⁻¹⁰⁷ Both types of potentiators showed *in vivo* activity.^{106,108}

mGlu₃

No specific allosteric modulators have yet been reported in the literature for this mGlu receptor subtype. Yet, a benzotriazole derivative of PTBE was noted to potentiate both mGlu₂ and mGlu₃ receptors.¹⁰⁶

Group III (Figures 5C, 5D and 5E)

To date, very few group III modulators have been disclosed. These include mGlu₄, mGlu₇ and mGlu₈ potentiators. PHCCC was initially described as an mGlu₁ receptor antagonist.⁵⁵ However, it was

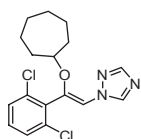
recently found that its (-) enantiomer potentiates mGlu₄ receptor activation.^{109,110} Two other mGlu₅ antagonists SIB 1893 and MPEP were reported to enhance agonist potency and efficacy at human mGlu₄ at higher concentrations.¹¹¹ Recently, potentiators of mGlu₇ (AMN 082 from Novartis)¹¹² and mGlu₈ (Thiomethylanilide A and B from NPS Pharmaceuticals)¹¹³ were disclosed.

Radiolabeled Ligands and PET Radiotracers

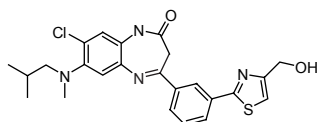
Several of the potent mGlu ligands have been radiolabeled: Quis, DCG-IV, LY 354740, (S)-AP4, (S)-DCPG, LY 341495, CPPG, EM-TBPC, R214127, LY 456066, MPEP and MMTEP. Some mGlu receptor ligands have been labeled with carbon-11 or fluorine-18 and used for PET imaging.¹¹⁴⁻¹¹⁶

Conclusion

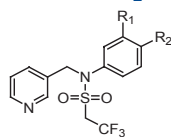
In the last few years some interesting new competitive ligands have been discovered, such as the selective mGlu₈ receptor agonist (S)-3,4-DCPG. However, the largest advances in mGlu receptor pharmacology have been made with allosteric modulators. These compounds are generally highly potent and selective. Moreover, many of them display *in vivo* activity and open the way to new therapeutic agents. Although some further subtype-selective compounds are still awaited, particularly for group III mGlu receptors, the panel of available mGlu receptor ligands is now rather broad and is allowing investigators to shed important new light on the physiological and pathological roles of the various mGlu receptor subtypes in the normal and diseased brain. This is currently ongoing in many laboratories and we anticipate watching the results unfold with great interest.

Figure 5 | Group II and Group III Allosteric Modulator Structures and Potencies**A. mGlu₂ Receptor Antagonists**

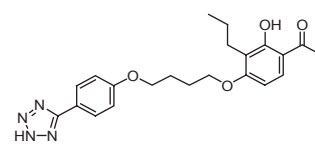
Ro 64-5229
IC₅₀ = 109 nM



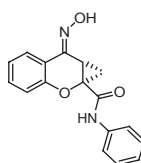
CH-DBO
K_i = 3 nM

B. mGlu₂ Receptor Potentiators

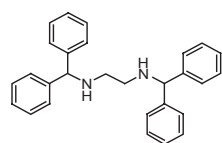
R₁ = H R₂ = 2-methoxyphenoxy
LY487379 (EC₅₀ = 0.3 μM)
R₁ = 1-methylbutoxy R₂ = H
2,2,2-TEMPS (EC₅₀ = 14 nM)



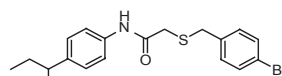
PTBE
EC₅₀ = 0.43 μM

C. mGlu₄ Receptor Potentiator

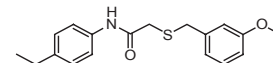
(-)-PHCCC
EC₅₀ = 3 μM

D. mGlu₇ Receptor Potentiator

AMN 082
EC₅₀ = 0.06-0.29 μM

E. mGlu₈ Receptor Potentiators

Thiomethylanilide A
EC₅₀ = 1.0 μM



Thiomethylanilide B
EC₅₀ = 5.4 μM

(Bold Text Denotes Compounds Available From Tocris)

List of Acronyms

ABHxD	2-aminobicyclo[2.1.1]hexane-2,5-dicarboxylic acid
ACPD	1-aminocyclopentane-1,3-dicarboxylic acid
ACPT-I	(1S,3R,4S)-1-aminocyclopentane-1,3,4-tricarboxylic acid
ACPT-II	(1R,3R,4S)-1-aminocyclopentane-1,3,4-tricarboxylic acid
(+)-ACPT-III	(3S,4S)-1-aminocyclopentane-1,3,4-tricarboxylic acid
ADED	(2S,4S)-2-amino-4-(2,2-diphenylethyl)pentane-1,5-dioic acid
ADX47273	(S)-[4-fluoro-phenyl]-[3-[3-(4-fluoro-phenyl)-[1,2,4]oxadiazol-5-yl]-piperidin-1-yl]-methanone
AMN 082	N,N'-dibenzhydrylethane-1,2-diamine
AMPA	2-amino-3-(3-hydroxy-5-methylisoxazol-4-yl)propionic acid
HomoAMPA	2-amino-4-(3-hydroxy-5-methylisoxazol-4-yl)butyric acid
AP4	2-amino-4-phosphono-butyric acid
APCPr	1-amino-2-phosphonomethylcyclopropane carboxylic acid
APDC	4-aminopyrrolidine-2,4-dicarboxylic acid
APTc	4-aminopyrrolidine-1,2,4-tricarboxylic acid
BAY 36-7620	(3aS,6aS)-6a-naphthalen-2-ylmethyl-5-methyliden-hexahydro-cyclopental[c]furan-1-on
BnAPDC	N-benzyl-4-aminopyrrolidine-2,4-dicarboxylic acid
BnQuis	α -benzylquisqualic acid
BOMA	2-[4-(1,3-benzoxazol-2-yl)-2-methoxyphenyl]acetone
L-CCG-I	(2S,1'S,2'S)-2-(carboxycyclopropyl)glycine
3'Me-CCG	(2S,1'S,2'S,3'R)-2-(3'-methyl-2'-carboxycyclopropyl)glycine
3'HM-CCG	(2S,1'S,2'R,3'R)-2-(3'-hydroxymethyl-2'-carboxycyclopropyl)glycine
3'SM-CCG	(2S,1'R,2'R,3'R)-2-(3'-mercaptopmethyl-2'-carboxycyclopropyl)glycine
mCD-CCG	2-(di-meta-chlorophenylethyl)-2-(2'-carboxycyclopropyl)glycine
XE-CCG	(2S,1'S,2'S,3'R)-2-(3'-xanthenylethyl-2'-carboxycyclopropyl)glycine
CBQA	1-amino-3-[3',5'-dioxo-1',2',4'-oxadiazolidinyl]cyclobutane-1-carboxylic acid
CDPPB	3-cyano-N-(1,3-diphenyl-1H-pyrazol-5-yl)benzamide
CH-DBO	8-chloro-4-[3-(4-hydroxymethyl-thiazol-2-yl)-phenyl]-7-isobutyl-methylamino)-1,3-dihydro-benzo[b][1,4]diazepin-2-one
CHPG	2-chloro-5-hydroxyphenylglycine
4C3H2MPG	4-carboxy-3-hydroxy-2-methylphenylglycine
4C2MPG	(+)-4-carboxy-2-methylphenylglycine
4CPG	4-carboxyphenylglycine
(-)-CPCCOEt	(1aS,7aS)-(2-hydroxyimino-1a,2-dihydro-1H-7-oxacyclopropa[b]naphthalene-7a-carboxylic acid ethyl ester
CPPG	α -cyclopropyl-4-phosphonophenylglycine
CPPHA	N-(4-chloro-2-[[1,3-dioxo-1,3-dihydro-2H-isoindol-2-yl)methyl]phenyl)-2-hydroxybenzamide
DCG IV	(2S,1'R,2'R)-2-(2',3'-dicarboxycyclopropyl)glycine
3,4-DCPG	3,4-dicarboxyphenylglycine
3,5-DHPG	3,5-dihydroxyphenylglycine
DCB	3,3'-dichlorobenzaldazine
DFB	3,3'-difluorobenzaldazine
DMeOB	3,3'-dimethoxybenzaldazine
DM-PPP	3,5-dimethyl-pyrrole-2,4-dicarboxylic acid 2-propylester 4-[(S)-1,2,2-trimethylpropyl]ester
EM-TBPC	1-ethyl-2-methyl-6-oxo-4-(1,2,4,5-tetrahydro-benzo[d]azepin-3-yl)-1,6-dihydro-pyrimidine-5-carbonitrile
L-Glu	L-glutamate
GPCR	G-protein-coupled receptor
HYDIA	(1S,2R,3R,5R,6S)-3-hydroxy-2-aminobicyclo[3.1.0]hexane-2,6-dicarboxylic acid
JNJ 16259685	(3,4-dihydro-2H-pyrano[2,3-b]quinolin-7-yl)-(cis-4-methoxycyclohexyl)-methanone
LY 339840 (4C3H2MPG)	(RS)-4-carboxy-3-hydroxy-2-methylphenylglycine
LY 341495	(2S,1'S,2'S)-2-(9-xanthylmethyl)-2-(2'-carboxycyclopropyl)glycine
LY 354740	(1S,2S,5R,6S)-2-aminobicyclo[3.1.0]hexane-2,6-dicarboxylic acid
LY 354740-6F	(1R,2S,5R,6R)-2-amino-6-fluorobicyclo[3.1.0]hexane-2,6-dicarboxylic acid
LY 367366	α -thioxanthylmethyl-4-carboxyphenylglycine
LY 367385 (4C2MPG)	(+)-4-carboxy-2-methylphenylglycine
LY 379268	2-oxa-4-aminobicyclo[3.1.0]hexane-4,6-dicarboxylic acid
LY 389795	2-thia-4-aminobicyclo[3.1.0]hexane-4,6-dicarboxylic acid
LY 393053	(\pm)-2-amino-2-(3-cis and trans-carboxycyclobutyl)-3-(9-thioxanthyl)propionic acid
LY 393675	(S)-cis- α -thioxanthylmethyl-3-carboxycyclobutylglycine
LY 397366	α -thioxanthylmethyl-4-carboxyphenylglycine
LY 456066	2-[4-(indan-2-ylamino)-5,6,7,8-tetrahydroquinazolin-2-yl]sulfanyl]-ethanol
LY 487379	N-(4-(2-methoxyphenoxy)phenyl)-N-(2,2,2-trifluoroethylsulfonyl)pyrid-3-ylmethylamine
MAP4	2-methyl-2-amino-4-phosphono-butyric acid
MCCG	(2S,3S,4S)-2-methyl-2-(carboxycyclopropyl)glycine
MCPG	α -methyl-4-carboxyphenylglycine
MGS0008	(1S,2S,3S,5R,6S)-2-amino-3-fluorobicyclo[3.1.0]hexane-2,6-dicarboxylic acid
MGS0028	(1R,2S,5S,6S)-2-amino-6-fluoro-4-oxobicyclo[3.1.0]hexane-2,6-dicarboxylic acid
MGS0039	(1R,2R,3R,5R,6R)-2-amino-3-(3,4-dichlorobenzyloxy)-6-fluorobicyclo[3.1.0]hexane-2,6-dicarboxylic acid
MPPG	α -methyl-4-phosphonophenylglycine
MPEP	2-methyl-6-(phenylethynyl)pyridine
5MPEP	5-methyl-2-(phenylethynyl)pyridine
M-MPEP	2-[(3-methoxyphenyl)ethynyl]-6-methylpyridine
MTEB	5-[(2-methyl-1,3-thiazol-4-yl)ethynyl]-2,3'-bipyridine
MTEP	3-[(2-methyl-1,3-thiazol-4-yl)ethynyl]pyridine
MM-MTEP	3-(methoxymethyl)-5-[(2-methyl-1,3-thiazol-4-yl)-ethynyl]pyridine
NAAG	N-acetyl-L-aspartyl-L-glutamate
NM-APDC	(2R,4R)-4-amino-1-(1-naphthylmethyl)pyrrolidine-2,4-dicarboxylic acid
NMDA	N-methyl-D-aspartate
NPS2390	N-(1-adamantyl)-2-quinoxaline-carboxamide
PPG	4-phosphonophenylglycine
PHCCC	N-phenyl-7-(hydroxylimino)cyclopropa[b]chromen-1a-carboxamide
3,5-dimethyl PPP	3,5-dimethyl-pyrrole-2,4-dicarboxylic acid 2-propylester 4-[(S)-1,2,2-trimethyl-propyl]ester
PTBE	1-(2-hydroxy-3-propyl-4-4-[4-(2H-tetrazol-5-yl)phenoxy]butoxyphenyl)ethanone
PTeB	3-(5-pyridin-2-yl-2H-tetrazol-2-yl)benzotrile
Quis	quisqualate
R214127	1-(3,4-dihydro-2H-pyrano[2,3-b]quinolin-7-yl)-2-phenyl-1-ethanone
Ro 01-6128	diphenylacetyl-carbamic acid ethyl ester
Ro 64-5229	1-Z-[2-cycloheptyloxy-2-(2,6-dichlorophenyl)viny]-[1,2,4-triazole)
Ro 67-4853	(9H-xanthen-9-carbonyl)-carbamic acid butyl ester
Ro 67-7476	(S)-2-(4-fluoro-phenyl)-1-(toluene-4-sulfonyl)-pyrrolidine
SIB 1757	6-methyl-2-(phenylazo)-3-pyridinol
SIB 1893	(E)-2-methyl-6-(2-phenylethynyl)pyridine
SOP	Serine-O-phosphate
2,2,2-TEMPS	2,2,2-trifluoroethyl-N-[3-(1-methylbutoxy)phenyl]-N-(3-pyridinylmethyl)sulfonamide
Thiomethylaniide A	2-[(4-bromobenzyl)thio]-N-(4-isobutylphenyl)acetamide
Thiomethylaniide B	2-[(4-methoxybenzyl)thio]-N-(4-ethylphenyl)acetamide
YM 298198	6-Amino-N-cyclohexyl-N,3-dimethyl-thiazolo[3,2-a]benzimidazole-2-carboxamide

References

- Dingledine *et al* (1999) *Pharmacol. Rev.* **51** 7.
- Conn and Pin (1997) *Ann. Rev. Pharmacol. Toxicol.* **37** 205.
- Pin and Acher (2002) *Curr. Drug Targets - CNS & Neur. Dis.* **1** 297.
- Marek (2004) *Curr. Opin. Pharmacol.* **4** 18.
- Schoepp *et al* (1999) *Neuropharmacology* **38** 1431.
- Pin *et al* (1999) *Eur. J. Pharmacol.* **375** 277.
- Pin *et al* (2003) *Pharmacol. Ther.* **98** 325.
- O'Hara *et al* (1993) *Neuron* **11** 41.
- Bessis *et al* (2000) *Protein Sci.* **9** 2200.
- Kunishima *et al* (2000) *Nature* **407** 971.
- Bessis *et al* (2002) *Proc. Natl. Acad. Sci. USA* **99** 11097.
- Hampson *et al* (1999) *J. Biol. Chem.* **274** 33488.
- Malherbe *et al* (2001) *Mol. Pharmacol.* **60** 944.
- Bertrand *et al* (2002) *J. Med. Chem.* **45** 3171.
- Rosemond *et al* (2002) *J. Biol. Chem.* **277** 7333.
- Rosemond *et al* (2004) *Mol. Pharmacol.* **66** 834.
- Acher and Bertrand (2005) *Biopolymers* **80** 357.
- May and Christopoulos (2003) *Curr. Opin. Pharmacol.* **3** 551.
- Litschig *et al* (1999) *Mol. Pharmacol.* **55** 453.
- Augelli-Szafran and Schwarz (2003) *Ann. Report Med. Chem.* **38** 21.
- Kew (2004) *Pharmacol. Ther.* **104** 233.
- Sabbatini and Micheli (2004) *Exp. Opin. Ther. Patents* **14** 1593.
- Bessis *et al* (1999) *Neuropharmacology* **38** 1543.
- Monn *et al* (1997) *J. Med. Chem.* **40** 528.
- Johnson *et al* (2002) *Drug Metab. Disp.* **30** 27.
- Palmer *et al* (1989) *Eur. J. Pharmacol.* **166** 585.
- Doherty *et al* (1997) *Neuropharmacology* **36** 265.
- Littman *et al* (1999) *J. Med. Chem.* **42** 1639.
- Monn *et al* (1999) *J. Med. Chem.* **42** 1027.
- Nakazato *et al* (2000) *J. Med. Chem.* **43** 4893.
- Collado *et al* (2002) *J. Med. Chem.* **45** 3619.
- Collado *et al* (2004) *J. Med. Chem.* **47** 456.
- Gonzalez *et al* (2005) *Bioorg. Med. Chem.* **13** 6556.
- Dominguez *et al* (2005) *J. Med. Chem.* **48** 3605.
- Kroona *et al* (1991) *J. Med. Chem.* **34** 1692.
- Sibille *et al* (2002) *Drugs Fut.* **27** (Suppl. A) 450.
- Gasparini *et al* (1999) *J. Pharmacol. Exp. Ther.* **290** 1678.
- Gasparini *et al* (2000) *Bioorg. Med. Chem. Lett.* **10** 1241.
- Thomas *et al* (2001) *Neuropharmacology* **40** 311.
- Acher *et al* (1997) *J. Med. Chem.* **40** 3119.
- Tückmantel *et al* (1997) *Bioorg. Med. Chem. Lett.* **7** 601.
- Ahmadian *et al* (1997) *J. Med. Chem.* **40** 3700.
- Schann *et al* (2005) *Neuropharmacology* **49** (Suppl. 1) 272.
- Chen *et al* (2000) *Neuroscience* **95** 787.
- Kingston *et al* (2002) *Neurosci. Lett.* **330** 127.
- Melendez *et al* (2005) *J. Pharmacol. Exp. Ther.* **314** 139.
- Pellicciari *et al* (2001) *Bioorg. Med. Chem. Lett.* **11** 3179.
- Sørensen *et al* (2003) *Bioorg. Med. Chem.* **11** 197.
- Chaki *et al* (2004) *Neuropharmacology* **46** 457.
- Nakazato *et al* (2004) *J. Med. Chem.* **47** 4570.
- Adam *et al* (1999) *Neuropharmacology* **38** A1 abstract 3.
- Conway *et al* (2001) *Bioorg. Med. Chem. Lett.* **11** 777.
- Brabet *et al* (1998) *Neuropharmacology* **37** 1043.
- Wright *et al* (2000) *Naunyn-Schmied. Arch. Pharmacol.* **362** 546.
- Annoura *et al* (1996) *Bioorg. Med. Chem. Lett.* **6** 763.
- Ott *et al* (2000) *J. Med. Chem.* **43** 4428.
- Van Wagenen *et al* (1998) *Soc. Neurosci. Abstract* **24**.
- Lavreysen *et al* (2003) *Mol. Pharmacol.* **63** 1082.
- Carroll *et al* (2001) *Mol. Pharmacol.* **59** 965.
- Li *et al* (2002) *Neuropharmacology* **43** A79.
- Ambler *et al* (2001) *Patent WO 01/32632*.
- Lavreysen *et al* (2004) *Neuropharmacology* **47** 961.
- Mabire *et al* (2005) *J. Med. Chem.* **48** 2134.
- Micheli *et al* (2003) *Bioorg. Med. Chem.* **11** 171.
- Malherbe *et al* (2003) *J. Biol. Chem.* **278** 8340.
- Binggeli *et al* (2002) *Patent WO 02051418*.
- Kohara *et al* (2005) *J. Pharmacol. Exp. Ther.* **315** 163.
- Zheng *et al* (2005) *J. Med. Chem.* **48** 7374.
- Steckler *et al* (2005) *Behav. Brain Res.* **164** 52.
- Bleicher *et al* (2000) *Patent WO 0063166*.
- Knoflach *et al* (2001) *Proc. Natl. Acad. Sci. USA* **98** 13402.
- Wichmann *et al* (2002) *Farmacol.* **57** 989.
- Varney *et al* (1999) *J. Pharmacol. Exp. Ther.* **290** 170.
- Gasparini *et al* (1999) *Neuropharmacology* **38** 1493.
- Spooren *et al* (2001) *TIPS* **22** 331.
- Gasparini *et al* (2002) *Bioorg. Med. Chem. Lett.* **12** 407.
- Cosford *et al* (2003) *J. Med. Chem.* **46** 204.
- Cosford *et al* (2003) *Bioorg. Med. Chem. Lett.* **13** 351.
- Anderson *et al* (2002) *J. Pharmacol. Exp. Ther.* **303** 1044.
- Roppe *et al* (2004) *Bioorg. Med. Chem. Lett.* **14** 3993.
- Wang *et al* (2004) *Bioorg. Med. Chem.* **12** 17.
- Roppe *et al* (2004) *J. Med. Chem.* **47** 4645.
- Poon *et al* (2004) *Bioorg. Med. Chem. Lett.* **14** 5477.
- Tehrani *et al* (2005) *Bioorg. Med. Chem. Lett.* **15** 5061.
- Bonnefous *et al* (2005) *Bioorg. Med. Chem. Lett.* **15** 1197.
- Kamenecka *et al* (2005) *Bioorg. Med. Chem. Lett.* **15** 4350.
- Porter *et al* (2005) *J. Pharmacol. Exp. Ther.* **315** 711.
- Malherbe *et al* (2003) *Mol. Pharmacol.* **64** 823.
- O'Brien *et al* (2003) *Mol. Pharmacol.* **64** 731.
- O'Brien *et al* (2004) *J. Pharmacol. Exp. Ther.* **309** 568.
- Rodríguez *et al* (2005) *Mol. Pharmacol.* **68** 1793.
- Lindsley *et al* (2004) *J. Med. Chem.* **47** 5825.
- Kinney *et al* (2005) *J. Pharmacol. Exp. Ther.* **313** 199.
- Bessis *et al* (2005) *Patent WO2005044797*.
- Le Poul *et al* (2005) *Neuropharmacology* **49** (Suppl. 1) 252.
- Kolczewski *et al* (1999) *Bioorg. Med. Chem. Lett.* **9** 2173.
- Adam *et al* (1999) *Neuropharmacology* **38** A1 abstract 2.
- Adam *et al* (2002) *Patent US2002/0193367*.
- Schaffhauser *et al* (2003) *Mol. Pharmacol.* **64** 798.
- Barda *et al* (2004) *Bioorg. Med. Chem. Lett.* **14** 3099.
- Hu *et al* (2004) *Bioorg. Med. Chem. Lett.* **14** 5071.
- Pinkerton *et al* (2004) *J. Med. Chem.* **47** 4595.
- Pinkerton *et al* (2004) *Bioorg. Med. Chem. Lett.* **14** 5867.
- Pinkerton *et al* (2005) *Bioorg. Med. Chem. Lett.* **15** 1565.
- Cube *et al* (2005) *Bioorg. Med. Chem. Lett.* **15** 2389.
- Govek *et al* (2005) *Bioorg. Med. Chem. Lett.* **15** 4068.
- Bonnefous *et al* (2005) *Bioorg. Med. Chem. Lett.* **15** 4354.
- Johnson *et al* (2005) *Psychopharmacology* **179** 271.
- Maj *et al* (2003) *Neuropharmacology* **45** 895.
- Marino *et al* (2003) *Proc. Natl. Acad. Sci. USA* **100** 13668.
- Mathiesen *et al* (2003) *Br. J. Pharmacol.* **138** 1026.
- Mitsukawa *et al* (2005) *Proc. Natl. Acad. Sci. USA* **102** 18712.
- Wilson *et al* (2005) *Neuropharmacology* **49** (Suppl. 1) 278.
- Hamill *et al* (2005) *Synapse* **56** 205.
- Huang *et al* (2005) *J. Med. Chem.* **48** 5096.
- Yu *et al* (2005) *Nucl. Med. Biol.* **32** 631.
- Naples and Hampson (2001) *Neuropharmacology* **40** 170.

Metabotropic Glutamate Receptor Compounds Available from Tocris

Group I Selective Metabotropic Glutamate Receptor Ligands

Agonists

- 0187** (\pm)-*trans*-ACPD
Group I/group II mGlu agonist
- 0284** (1*S*,3*R*)-ACPD
Group I/group II mGlu agonist
- 0860** *t*ADA
Group I agonist; some mGlu₅ selectivity
- 1049** CHPG
mGlu₅ selective agonist
- 0342** (RS)-3,5-DHPG
Selective group I mGlu agonist
- 0805** (S)-3,5-DHPG
Active enantiomer of (0342)
- 1058** L-3'-F₂CCG-I
Potent group I/group II agonist
- 0218** L-Glutamic acid
Endogenous, non-selective agonist
- 1826** Group I mGlu Receptor Tocriset
Selection of 5 group I mGlu receptor ligands
- 0324** (RS)-3-Hydroxyphenylglycine
Group I mGlu agonist
- 0326** (S)-3-Hydroxyphenylglycine
Group I mGlu agonist, active isomer
- 1829** Mixed mGlu Receptor Tocriset
Selection of 5 mixed mGlu receptor ligands
- 1490** MNI-caged-L-glutamate
Stable photoreleaser of L-glutamate
- 0188** L-Quisqualic acid
Very potent group I mGlu agonist
- 0162** S-Sulfo-L-cysteine
Group I agonist

Antagonists

- 2254** ACDPP
Selective mGlu₅ receptor antagonist
- 0904** AIDA
Potent, selective group I mGlu antagonist
- 0125** DL-AP3
Group I mGlu antagonist
- 0329** (S)-3-Carboxy-4-hydroxyphenylglycine
Group I antagonist/group II agonist

- 0320** (S)-4-Carboxy-3-hydroxyphenylglycine
Group I antagonist/group II agonist
- 0323** (S)-4-Carboxyphenylglycine
Competitive group I mGlu antagonist/weak group II agonist
- 1028** CPCCOEt
Selective, non-competitive mGlu₁ receptor antagonist
- 1009** E4CPG
Group I/group II mGlu antagonist
- 2386** Fenobam
Potent and selective mGlu₅ antagonist
- 1826** Group I mGlu Receptor Tocriset
Selection of 5 group I mGlu receptor ligands
- 1749** HexylHIBO
Group I mGlu antagonist
- 1750** (S)-HexylHIBO
Group I mGlu antagonist
- 2333** JNJ 16259685
Extremely potent, mGlu₁-selective non-competitive antagonist
- 1237** LY 367385
Selective mGlu_{1a} antagonist
- 2196** 3-MATIDA
Potent, selective mGlu₁ antagonist
- 0336** (RS)-MCPG
Non-selective mGlu antagonist
- 0337** (S)-MCPG
Active isomer of (0336)
- 1829** Mixed mGlu Receptor Tocriset
Selection of 5 mixed mGlu receptor ligands
- 1212** MPEP
mGlu₅ subtype-selective antagonist
- 2390** MPMQ
Selective mGlu₁ antagonist
- 1027** PHCCC
Potent group I mGlu antagonist. Also mGlu₄ potentiator
- 1215** SIB 1757
Highly selective mGlu₅ antagonist
- 1214** SIB 1893
Highly selective mGlu₅ antagonist
- 2448** YM 298198
Highly potent, selective non-competitive mGlu₁ antagonist
- 2447** Desmethyl-YM 298198
Derivative of Cat. No. 2448

Other

- 1952 DCB**
Neutral allosteric modulator at mGlu₅
- 1625 DFB**
Allosteric potentiator at mGlu₅
- 1953 DMeOB**
Negative allosteric modulator at mGlu₅
- 2028 Anti-mGlu₁**
Antibody recognising rat mGlu₁ receptors
- 2032 Anti-mGlu₅**
Antibody recognising rat mGlu₅ receptors

Group II Selective Metabotropic Glutamate Receptor Ligands

Agonists

- 0187 (±)-trans-ACPD**
Group II/group I mGlu agonist
- 0284 (1S,3R)-ACPD**
Group II/group I mGlu agonist
- 1208 (2R,4R)-APDC**
Highly selective group II agonist
- 0329 (S)-3-Carboxy-4-hydroxyphenylglycine**
Selective group II mGlu agonist, also group I mGlu antagonist
- 0320 (S)-4-Carboxy-3-hydroxyphenylglycine**
Group II agonist/group I antagonist
- 0323 (S)-4-Carboxyphenylglycine**
Group I mGlu antagonist/weak group II agonist
- 0333 L-CCG-I**
Potent group II mGlu agonist
- 0975 DCG IV**
Potent group II mGlu agonist. Also group III mGlu antagonist and NMDA agonist
- 1058 L-3'-F₂-CCG-I**
Potent group II/group I agonist
- 0218 L-Glutamic acid**
Endogenous, non-selective agonist
- 1827 Group II mGlu Receptor Tocriset**
Selection of 5 group II mGlu receptor ligands
- 1829 Mixed mGlu Receptor Tocriset**
Selection of 5 mixed mGlu receptor ligands
- 1490 MNI-caged-L-glutamate**
Stable photoreleaser of L-glutamate
- 0391 Spaglumic acid**
Selective mGlu₃ agonist

Antagonists

- 1073 (RS)-APICA**
Selective group II antagonist
- 1009 E4CPG**
Group II/group I mGlu antagonist
- 0971 EGLU**
Highly selective group II mGlu antagonist
- 1827 Group II mGlu Receptor Tocriset**
Selection of 5 group II mGlu receptor ligands
- 1209 LY 341495**
Highly potent, selective group II antagonist
- 0336 (RS)-MCPG**
Non-selective mGlu antagonist
- 0337 (S)-MCPG**
Active isomer of (0336)
- 1829 Mixed mGlu Receptor Tocriset**
Selection of 5 mixed mGlu receptor ligands
- 0854 MSPG**
Group II/group III mGlu antagonist
- 0855 MTPG**
Group II/group III mGlu antagonist. More selective for group II than group III

Other

- 2027 Anti-mGlu₂**
Antibody recognising rat mGlu₂ receptors
- 2029 Anti-mGlu_{2/3}**
Antibody recognising human mGlu₂ and mGlu₃ receptors

Group III Selective Metabotropic Glutamate Receptor Ligands

Agonists

- 1111 ACPT-I**
Group III mGlu agonist
- 1113 (±)-ACPT-III**
Selective group III agonist
- 2385 AMN 082**
The first selective mGlu₇ agonist

- 0103 L-AP4**
Selective group III mGlu agonist
- 1394 (RS)-3,4-DCPG**
Potent systemically active anticonvulsant. Racemate of (1302)
- 1302 (S)-3,4-DCPG**
Potent, selective mGlu_{8a} agonist
- 0218 L-Glutamic acid**
Endogenous, non-selective agonist
- 1828 Group III mGlu Receptor Tocriset**
Selection of 5 group III mGlu receptor ligands
- 1026 HomoAMPA**
Potent, highly selective mGlu₆ agonist
- 1829 Mixed mGlu Receptor Tocriset**
Selection of 5 mixed mGlu receptor ligands
- 1490 MNI-caged-L-glutamate**
Stable photoreleaser of L-glutamate
- 0238 O-Phospho-L-serine**
Group III mGlu agonist
- 1220 (RS)-PPG**
Potent group III mGlu agonist

Antagonists

- 0972 CPPG**
Very potent group III mGlu antagonist
- 0975 DCG IV**
Group III antagonist/group II mGlu agonist
- 1828 Group III mGlu Receptor Tocriset**
Selection of 5 group III mGlu receptor ligands
- 1209 LY 341495**
Group II/III antagonist
- 0711 MAP4**
Selective group III mGlu antagonist
- 1829 Mixed mGlu Receptor Tocriset**
Selection of 5 mixed mGlu receptor ligands
- 0853 MPPG**
Group III/group II mGlu antagonist. More selective for group III than group II
- 0803 MSOP**
Specific group III mGlu antagonist
- 0854 MSPG**
Group III/group II mGlu antagonist
- 1369 UBP1112**
Group III mGlu antagonist

Other

- 2031 Anti-mGlu₇**
Antibody recognising human mGlu₇ receptors

Quisqualate-Sensitised AP6 Site

- 0101 DL-AP4**
Agonist/broad spectrum EAA antagonist
- 0102 D-AP4**
Less potent enantiomer
- 0103 L-AP4**
Agonist; also group III mGlu agonist
- 0105 DL-AP5**
Agonist; also NMDA antagonist
- 0106 D-AP5**
Less potent enantiomer. Also NMDA antagonist
- 0107 L-AP5**
Agonist; also NMDA antagonist
- 0341 L-AP6**
Selective agonist, highly potent
- 0188 L-Quisqualic acid**
Sensitiser

Miscellaneous Metabotropic Glutamate Receptor Compounds

- 1112 ACPT-II**
Competitive mGlu receptor antagonist
- 0216 L-Cysteinesulfinic acid**
Agonist at PLD-coupled mGlu receptor, also NMDA agonist
- 0285 Ibotenic acid**
Non-selective mGlu agonist, also NMDA agonist
- 1611 Lamotrigine**
Inhibits glutamate release. Anticonvulsant
- 2289 Lamotrigine isethionate**
Water-soluble salt of Cat. No. 1611
- 1721 Tocriscreen Glutamate Ligands**
Collection of glutamate receptor and related compounds