

Metabotropic Glutamate Receptors

Molecular Pharmacology

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.

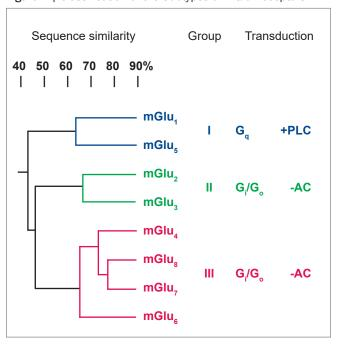
Contents Introduction ______1 Competitive Ligands 2-11 Agonists2 Group I......2 Antagonists......5 Group I11 mGlu₁ Positive Modulators.....11 mGlu₂ Positive Modulators......14 mGlu_{2/3} Antagonists......14 Metabotropic Glutamate Receptor Compounds21-23

Introduction

Glutamate is the major excitatory neurotransmitter 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. Glutamate also activates metabotropic glutamate (mGlu) receptors which modulate its release and postsynaptic response as well as the activity of other synapses. 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. Hi 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 their sequence homology, transduction mechanism and pharmacological profile (see Figure 1). Group I includes mGlu $_{\rm 1}$ and mGlu $_{\rm 5}$ receptors which couple to G $_{\rm q}$ and activate phospholipase C (PLC). Group II (mGlu $_{\rm 2}$, mGlu $_{\rm 3}$) and group III (mGlu $_{\rm 4}$, mGlu $_{\rm 6}$, mGlu $_{\rm 7}$, mGlu $_{\rm 8}$) receptors couple G $_{\rm 7}/{\rm G}_{\rm 9}$ and inhibit adenylyl cyclase (AC). Group I receptors are mostly located postsynaptically, thus their activation increases excitability. Conversely, group II/III receptors are generally presynaptic and their activation reduces glutamate release. Selective ligands have been found for each group and some of the subtypes, as described hereafter. $^{2.5-8}$

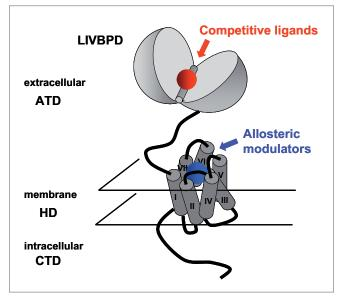
Figure 1 | Classification of the Subtypes of mGlu Receptors



mGlu receptors belong to class C of the GPCR superfamily.9 Like all GPCRs they hold a heptahelical domain (HD) in the membrane region. In addition, similar to all members of class C, they are characterized by a large extracellular amino terminal domain (ATD) where the glutamate binding site is found (see Figure 2). 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. 10-13 These amino acids bind to an open conformation of the protein, which subsequently closes 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 (see Figure 2). Moreover, it was shown that the closed conformation of this domain is required for receptor activation.¹⁴ Examination of the glutamate binding site in the eight mGlu receptor subtype crystal structures (mGlu₁, mGlu₂, mGlu₂)^{12,13} or homology models¹⁵⁻¹⁹ reveals a common binding motif for the α -amino and α -carboxylic functions of glutamate, ²⁰ while residues that bind the distal γ -carboxylate vary from one subtype to another.¹⁷ Thus, not surprisingly, all competitive agonists are α -amino acids, bearing various selective functional groups on their side chain⁶ including virtual screening hits and derivatives (see Figures 3A and 3B). The first generation of orthosteric ligands was followed by a second generation of allosteric modulators which bind in the HD.21 The first molecule described as a non-competitive mGlu receptor antagonist was CPCCOEt in the late nineties.²² Since then, numerous allosteric modulators have been discovered by high-throughput screening (HTS) in pharmaceutical companies.7,8,23

The purpose of the present article is to review our current knowledge of the pharmacology of mGlu receptors. Several detailed reviews^{2,4–6,8,24} have been published, therefore only the most potent and selective known ligands will be presented and emphasis will be placed on compounds that were more recently disclosed.

Figure 2 | Schematic Representation of an mGlu Receptor: the Two Orthosteric and Allosteric binding Sites are Indicated



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 pharmacophore²⁵ and homology models¹⁷ and found in X-ray structures. 12,13 For many years these compounds have been considered as valuable research tools but not as drug candidates because of their poor partition coefficient (LogP) related to their highly polar chemical structures and their lack of selectivity. Researchers at Eli Lilly were the first to show that such a glutamate analog, LY 354740, was able to pass the blood brain barrier and that its peptidyl prodrug, LY 544344, was orally active as an anticonvulsant and anxiolytic.26,27 A sulfonyl analog, LY 404039, orally administered as a methionine amide prodrug LY 2140023, being developed for the treatment of schizophrenia^{28,29} reached phase III clinical trials. Another advantage is that such drugs are barely metabolized since they are already quite hydrophilic³⁰ and few side effects are predicted. Other glutamate analogs were also shown to be systemically active such as (2R,4R)-APDC, 3'Me-CCG, 3'HM-CCG, (S)-DCPG, ACPT-I and LSP1-2111 (Figure 3A). Desensitization was also feared with continuous activation in the case of group II/III receptors, yet they are resistant to agonist-induced desensitization.31 Altogether these results promote a renewed interest in mGlu receptor competitive ligands.

Agonists (Table 1)

The first agonist that was able to discriminate between ionotropic and metabotropic glutamate receptors was *trans*-ACPD (1S,3*R* isomer).³² This ligand contributed considerably to the study of mGlu receptors despite its lack of subtype selectivity.^{2,5,24} 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.^{2,5,24} It can be noted that L-CCG-I and ABHxD-I are conformationally constrained and mimic the bioactive extended glutamate conformation common to all mGlu receptors.²⁵ When adding new chemical groups onto these structures, selectivity can be gained (Figure 3A).

Group I

Quisqualate (Quis) is the most potent group I agonist; however it also activates AMPA receptors, therefore its use is restricted. The most widely used group I selective agonist is (S)-3,5-DHPG, yet it exhibits only moderate potency. 2,5,24 CHPG 33 and Z/E-CBQA 34 have been claimed to specifically activate mGlu $_{\scriptscriptstyle 5}$ receptors although the affinity of the former is quite low. No specific mGlu $_{\scriptscriptstyle 1}$ competitive agonists have been disclosed to date.

Group I

LY 354740 was the first mGlu agonist reported to exhibit a nanomolar affinity.²⁶ It is group II selective, as are its oxy (LY 379268), thia (LY 389795)³⁵ and sulfonyl derivatives (LY 404039).³⁶ Introducing a fluorine atom at position 3 (MGS0008) or 6 (MGS0022) retained the potent activity which was further enhanced when a carbonyl group was added, as in the case of MGS0028.³⁷ This series of bicyclic glutamate analogs derives from the general agonist L-CCG-I where increased potency and

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

Pecentors		Group I		Group II		Group III			
Receptors		mGlu ₁	mGlu ₅	mGlu ₂	mGlu ₃	mGlu ₄	mGlu ₆	mGlu ₇	mGlu
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	-	-
	Quis ^{c,d}	0.03-3	0.02-0.3	100-1000	40-220	100-1000	n.e.	n.e.	720
Group I	(S)-3,5-DHPG ^{c,d}	6	2	n.e.	n.e.	n.e.	_	n.e.	n.e.
subtype-selective agonists	CHPG°	> 10000	750	_	-	_	-	_	_
	Z-CBQA°	> 1000	11	> 100	-	>100	_	_	_
	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
	MGS0008e	> 100	> 100	0.029	0.049	> 100	> 100	> 100	-
	MGS0022e	> 100	> 100	0.017	0.081	> 100	> 100	> 100	_
Group II	MGS0028 ^e	> 100	> 100	0.0006	0.0021	> 100	> 100	> 100	-
subtype-selective agonists	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	0.01
	LY 541850	n.e. ^t	n.e.t	0.16	ant.	n.e.t	_	n.e.t	n.e.
	2R,4R-APDCb,c	> 100	> 100	0.4	0.4	>300	110	>300	> 100
	DCG IVb,c	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	-	-
	(S)-AP4 ^{c,d}	> 1000	> 1000	> 1000	> 1000	0.2-1.2	0.6-0.9	160-500	0.06-0
	(S)-thioAP4 ^e	n.e. ^t	n.e. ^t	n.e. ^t	n.e.t	0.04	0.7	200	0.05
	(S)-SOP ^{c,d}	n.e.	n.e.	ant.	ant.	1-4	3	160-1200	2
	(1S,2R)-APCPr ^j	-	-	-	-	0.6	1.9	602	0.3
	LSP1-3081 ^k	n.e. ^t	n.e.t	n.e. ^t	n.e.t	0.16	3.3	419	0.51
	LSP1-2111	n.e. ^t	n.e.t	n.e.t	n.e.t	2.2	1.7	53	66
	LSP4-2022 ^m	n.e. ^t	n.e. ^t	n.e. ^t	n.e. ^t	0.11	4.2	12	29
Group III	ACPT-I ^{c,d,n}	ant.	> 1000	> 1000	-	7.2	18.4	-	10.1
subtype-selective agonists	(+)-ACPT-III ^{c,d,n}	ant.	-	ant.	-	8.8	19.2	_	7.0
	FP429 ^{n,o}	>5000	>5000	>5000	>5000	48	380	-	56 ^u
	PCG-1 ^p	> 1000	> 1000	> 1000	-	9.4	13	700	63 ^v
	(S)-PBPG ^q	> 1000	> 1000	310	-	4.2	66	> 1000	4.4v
	(S)-PPG ^{b,r}	>500	>500	>300	>200	3.2 (5.2)	(4.7)	48 (185)	(0.21
	(S)-HomoAMPA°	> 1000	> 1000	> 1000	-	> 1000	58	>5000	_
	BnAPDC°	> 1000	ant.	ant.	> 100	>300	20	_	>300
	(S)-3,4-DCPG ^{b,s}	ant.	> 100	> 100	> 100	8.8	3.6	> 100	0.032

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

 $^{^{}a}$ EC $_{50}$ or K $_{b}$ values (μ M) measured with rat or human (when indicated b) cloned receptors. ant. = antagonist; n. e. = no effect. References for agonist potencies which have been cited in reviews 5 and/or 6 are referred as such.

^b EC₅₀ or K_b values obtained with human mGlu receptors °Schoepp et al. $(1999)^5$ d Pin et al. $(1999)^6$ e Selvam et al. $(2007)^{45}$ f Nakazato et al. $(2000)^{37}$ g Collado et al. $(2002)^{38}$ h Collado et al. $(2004)^{39}$ i Dominguez et al. $(2005)^{42}$ j Kroona et al. $(1991)^{45}$, Sibille et al. $(2007)^{47}$ k Cuomo et al. $(2009)^{56}$ Beurrier et al. $(2009)^{57}$ m Selvam et al. $(2011)^{56}$ n Schann et al. $(2006)^{60}$ o Frauli et al. $(2007)^{61}$ p Amori et al. $(2006)^{52}$ q Filosa et al. $(2006)^{51}$

Gasparini et al. (1999)⁴⁷ and (2000)⁴⁸; data in parentheses refer to (±)-PPG⁴⁷ * Thomas et al. (2001)⁴⁹ † n.e. = no effect at 100 µM

^u partial agonist 36% Glu max⁶¹ ^v K, value

Figure 3A | Competitive mGlu Receptor Ligand Structures

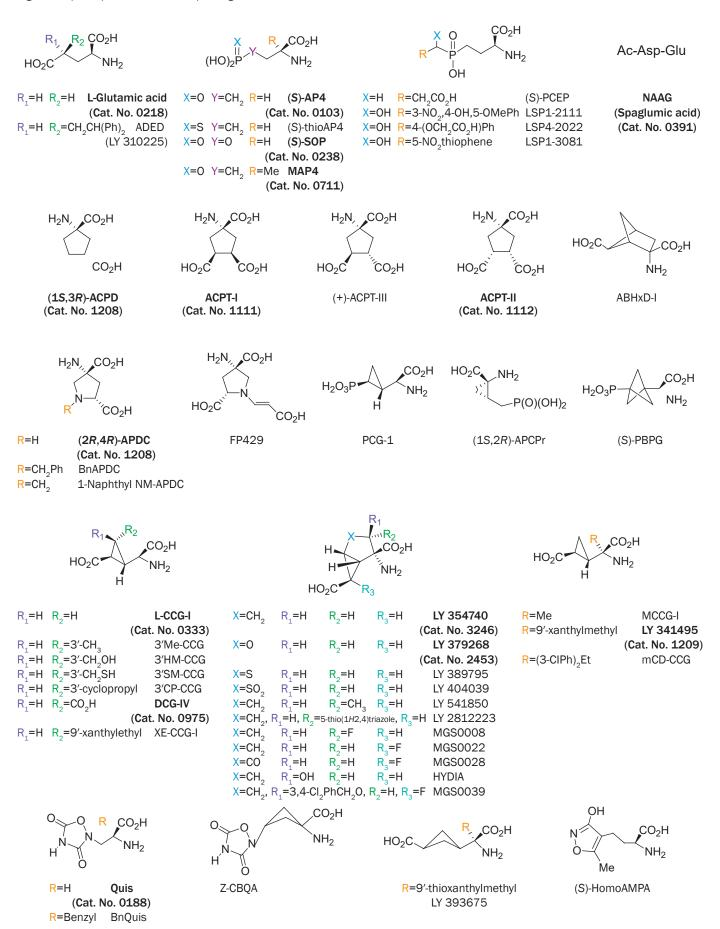


Figure 3B | Competitive mGlu Receptor Ligand Structures

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

group II selectivity were gained through the second hydrocarbon ring. However, it was shown that a methyl, hydroxymethyl or cyclopropyl substituent in the 3' position (3'Me-CCG, 3'HM-CCG and 3'CP-CCG respectively) provided agonists with similar potency.38-40 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 mGlu, agonist but is a full antagonist at mGlu₃.41 A similar selectivity was also reported for the C4β-methyl-substituted analog of LY 354740 (LY 541850). 42 Substitution by a thiotriazole group at this same position (LY 2812233) confers different pharmacological activity at the two subtypes.43 These three compounds selectively activate mGlu, while NAAG was the only reported mGlu, competitive agonist expected to discriminate between the two group II subtypes, however this was recently proved untrue.44 Other group II selective agonists have been described with submicromolar affinity, these include (2R,4R)-APDC and DCG-IV.

Group III

Most potent group III selective agonists bear a diacidic side chain that can interact with the highly basic distal binding pocket. 17,19 (S)-AP4 (L-AP4), (S)-thioAP4, 45 (S)-SOP (L-SOP) and (1S,2R)-APCPr46, 47 are the most potent, exhibiting submicromolar affinities at cloned receptors except for mGlu $_{7}$, to which all binding affinities are weak. (S)-PPG, 48,49 (S)-3,4-DCPG, 50 ACPT-I and (+)-ACPT-III, 51 (S)-PBPG and PCG-1 have also been described as micromolar agonists. Interestingly, a CCG derivative bearing a hydroxymethyl group in the 3′ position (3′HMCCG)

displays similar affinity for ${\rm mGlu_8}$ and ${\rm mGlu_{2/3}}$ receptors.³⁹ A new series of agonists deriving from a virtual screening hit, PCEP, was recently disclosed. 54-56 Among these is LSP1-308157 which displays potency close to L-AP4 and LSP1-211158 and LSP4-2022⁵⁶ that show a preference or selectivity for the mGlu₄ receptor respectively.56 In addition these agonists alleviate Parkinson's disease and anxiolytic symptoms following systemic injection in animal models.58-60 Nevertheless, very few group III mGlu receptor agonists are subtype-selective. FP429 is a full mGlu₄ and partial mGlu₈ agonist, 61,62 N-benzyl- APDC (BnAPDC)62 and (S)-homoAMPA $^{\rm 64}$ act at $\rm mGlu_{\rm 6}$ and (S)-3,4-DCPG at $\rm mGlu_{\rm 8}$ with an EC₅₀ over 2 orders of magnitude lower than at other group III receptors. 50 Interestingly, cinnabarinic acid, an endogenous metabolite of the kynurenine pathway, was demonstrated to be a weak mGlu, agonist, the first orthosteric agonist with non-αamino acid structure.65

Antagonists (Table 2)

Most competitive antagonists prevent the complete closing of the two lobes of the LIVBPD. Substitution of the α -proton of glutamate analogs by a methyl group, as in the case of MCCG, MCPG and MAP4, or a bulkier group as seen in LY 341495, turns the corresponding agonists (4CPG, AP4 and L-CCG-I) into antagonists. However, agonist properties can be recovered when the residues responsible for the hindrance are mutated. 14 Closing can also be disturbed by ionic repulsion, as seen with ACPT-II. 14

Figure 4 | Group I Allosteric Modulator Structures and Potencies

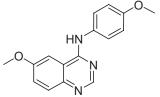
A. mGlu, Receptor Antagonists

Bay 36-7620 (Cat. No. 2501)

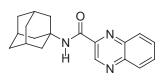
IC₅₀=160 nM K_i=11.2 nM DM-PPP IC₅₀=16 nM EM-TBPC K_i=11 nM JNJ 16259685 (Cat. No. 2333)

IC₅₀=0.55 nM K_i=0.34 nM

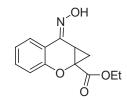
LY 456066 $IC_{50} = 12 \text{ nM (hmGlu}_{1})$



LY 456236 (Cat. No. 2390) $IC_{50} = 140 \text{ nM}$



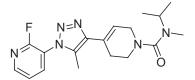
NPS 2390 (Cat. No. 4134) IC_{50} =5.2 nM chimera CaSR/mGlu₁ K_{i} = 1.4 nM



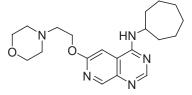
CPCCOEt (Cat. No. 1028) $IC_{50} = 10.3 \mu M$

 $K_i = 4.9 \mu M$ (-) isomer active

pyrazine-2-carboxamide PChPC K_i=9 nM



FTIDC (Cat. No. 5000)

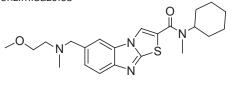


azaquinazole CMPPA K_i=6 nM

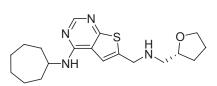
thiazolobenzimidazoles

$$H_2N$$

YM 298198 (Cat. No. 2448) IC_{50} =16 nM



YM 202074 (Cat. No. 3413) IC₅₀=8.6 nM



YM 230888 (Cat. No. 2986) IC₅₀=16 nM

$$\bigcup_{N=N}^{O} \bigvee_{N=N}^{N} \bigvee_{F}$$

 $\begin{array}{c} \rm isoindolone \\ \rm CFMTI \\ \rm IC_{50} = 2.6 \ nM \ (hmGlu_1) \end{array}$

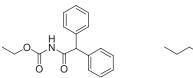
adamantyl methanone AdPyM K_i=24 nM

triazafluorenone A 841720 IC_{50} =10 nM (hmGlu₁)

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

Figure 4 | Group I Allosteric Modulator Structures and Potencies

B. mGlu, Receptor Potentiators



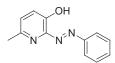
Ro 01-6128 (Cat. No. 4348) $EC_{50} = 0.2 \ \mu M$

Ro 67-4853 (Cat. No. 4347) $EC_{50} = 0.07 \mu M$

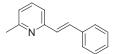
VU 71 EC₅₀= 2.4 μM

Ro 67-7476 (Cat. No. 4376) EC_{50} = 0.2 μ M

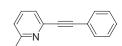
C. mGlu_s Receptor Antagonists; Alkyne Series



SIB 1757 (Cat. No 1215) $IC_{50} = 0.37 \ \mu M \ (hmGlu_5)$



SIB 1893 (Cat. No. 1214) $IC_{50} = 0.29 \mu M \text{ (hmGlu}_5)$



MPEP (Cat. No. 1212) IC₅₀ = 36 nM (hmGlu₅)

 $\begin{array}{c} \text{M-MPEP} \\ \text{IC}_{50} \text{= 10 nM (hmGlu}_{5}) \end{array}$

R=H MTEP (Cat. No. 2921) R=CH₂OMe MM-MTEP $IC_{50} = 5 \text{ nM}$ (hmGlu₅) $IC_{50} = 7 \text{ nM}$ (hmGlu₅)

M-PEP γ IC₅₀=1 nM (hmGlu₅)

 $\mathbb{S}_{N}^{\text{opt}} = \mathbb{Q}_{N}^{\text{opt}}$

MTEBP IC₅₀=2 nM (hmGlu₅)

S CN

R=H F-MTEB IC₅₀=0.08 nM R=F SP203 IC₅₀=0.036 nM

OMe H₂N

N N

AFQ056

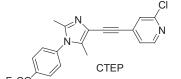
ADX10059 (Cat. No. 4416)

ADX48621

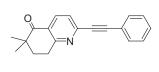
ethynylbenzamide IC₅₀=8 nM

ethynylpyrrolopyrazine $IC_{50} = 0.63$

5-(phenylethynyl)pyrimidine R=3-Me antago IC_{50} =7.5 nM R=4-Me ago IC_{50} =3.3 μ M



CTEP $IC_{50} = 6.4 \text{ nM (hmGlu}_{5})$



MRZ 8676 $IC_{50} = 23 \text{ nM (hmGlu}_5)$

Group I

The first generation of group I mGlu receptor 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 seen in LY 367366, but this derivative is also able to antagonize group II/III receptor activation. The highest potency was then found with α -substituted 3-carboxycyclobutylglycines such as LY 393675 (cis isomer) and its trans isomer, or a cis/trans mixture like LY 393053. This latter mixture was shown to be systemically active and to inhibit both mGlu $_1$ and mGlu $_5$ as well as other group II/III mGlu receptors. Although slightly less potent, LY 367385 (4C2MPG) and LY 339840 (4C3H2MPG) display subtype I selectivity; however, LY 367385 was also shown to inhibit the cystine/glutamate exchanger. No mGlu $_5$ selective and competitive antagonists have been described to date.

Group II

As most potent group II agonists derive 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 LY 341495,⁵ a fluorinated derivative⁶⁹ and XECCG69 holding

a 9'-xanthylmethyl or 9'-xanthylethyl moiety in the α - or 3'-position, display nanomolar affinities. The α -xanthyl moiety can be replaced by two substituted phenyl groups while retaining potency (e.g. mCD-CCG). As reported previously, stereospecific substitution at the 3-position of the agonist LY 354740 is critical for agonist/antagonist property. HYDIA71, and several O-benzyl derivatives such as MGS0039 exhibit high competitive group II antagonist activity. As Systemic and antidepressant-like effects were observed with both LY 341495 and MGS0039. Other arylalkylsubstituted glutamate and glutamate analogs such as ADED (LY 310225), (S)-BnQuis and NM-APDC display group II selectivity with ICs0 values in the micromolar range.

Group III

No highly potent and group III-selective competitive antagonists have been reported to date. 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, as in the case of MPPG. Addition of a substituent in the 3-position leads to similar group III antagonist activity but increases selectivity for group III over group II. 76 CPPG, the analog of MPPG bearing an α -cyclopropyl

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

Popontors		Group II Group II		up II	Group III				
Receptors		mGlu ₁	mGlu₅	mGlu ₂	mGlu ₃	m G lu ₄	mGlu ₆	mGlu ₇	mGlu ₈
Non-selective agonists	LY 341495b,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
	LY 367385b,f	8.8	>300	>300	-	>300	-	-	-
Group I	LY 367366 ^{b,c}	6.6	5.6	-	-	-	-	-	-
subtype-selective agonists	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
	ADED ^{b,c}	>300	>300	18	6.1	>300	-	>300	>300
	(S)-BnQuis ^{b,c}	300	300	7.1	-	n.e.	n.e.	-	-
Group II	mCD-CCG ^g	43	49	0.007	0.010	-	_	-	1.8
subtype-selective	HYDIA ^h	>100	> 100	0.10	0.11	22	-	-	15 (ago)
agonists	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 agonists	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

($\operatorname{\textbf{Bold}}$ text denotes compounds available from Tocris at time of publication)

^a IC₅₀ or K_b values (μM) measured with rat or human (when indicated^b) cloned receptors. ago. = agonist; n.e. = no effect. References for antagonist potencies which have been cited in reviews ⁵ and/or ⁶ are referred as such.

^b IC₅₀ or K_b values obtained with human mGlu receptors ^c Schoepp et al. (1999)⁵ ^d Pin et al. (1999)⁶ ^e Chen et al. (2000)⁶⁶ ^f Kingston et al. (2002)⁶⁷

⁵⁰ Sorensen (2003)⁷¹ h Adam (1999)^{72,216} Chaki et al. (2004)⁷⁴ Pellicciari et. al. (2001)⁷⁰ k Conway (2001)⁷⁷; Naples (2001)²¹⁷; Wright (2000)⁷⁹

Figure 4 | Group I Allosteric Modulator Structures and Potencies

D. mGlu_s Receptor Antagonists; Alkyne Biostere Series

tetrazole $IC_{50} = 4 \text{ nM (hmGlu}_5)$

oxadiazole VU 0285683 IC_{50} = 24 nM

7-arylquinoline $IC_{50} = 0.8 \text{ nM}$

ACDPP (Cat. No. 2254) IC₅₀= 134 nM

oxazolo-azepine IC_{50} = 16 nM

arylcarboxamide $IC_{50} = 14 \text{ nM}$

E. mGlu₅ Receptor Antagonists; Other Series

arylbenzoxazole BOMA IC₅₀= 3 nM

phenyltetrazole IC₅₀= 213 nM

thiazolotriazole GSK 2210875 IC_{50} = 40 nM (hmGlu₅)

carbamoyloxime $IC_{50} = 15 \text{ nM}$

pyrrolidinylpyridine $IC_{50} = 17 \text{ nM (hmGlu}_{5})$

piperidylamide IC₅₀= 32 nM

$$O = N N N N C$$

Fenobam (Cat. No. 2386) IC_{50} = 58 nM

anilinoquinazoline IC_{50} = 96 nM

benzimidazole IC₅₀= 24 nM

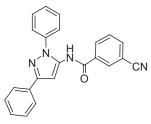
 $\begin{array}{lll} \text{DMeOB (Cat. No. 1953)} \\ \text{mGIu}_{5} & \text{IC}_{50} \! = \! 3 \; \mu\text{M} \\ \text{mGIu}_{4} & \text{IC}_{50} \! = \! 35 \; \mu\text{M} \\ \text{mGIu}_{8} & \text{IC}_{50} \! = \! 50 \; \mu\text{M} \end{array}$

Figure 4 | Group | Allosteric Modulator Structures and Potencies

F. mGlu₅ Receptor Potentiators

DFB (Cat. No. 1625) $EC_{50} = 2.4 \, \mu M \, (mGlu_5)$

CPPHA (Cat. No. 4787) $EC_{50} = 0.14 \, \mu M \, (mGlu_5)$



CDPPB (Cat. No. 3235) $EC_{50} = 20 \text{ nM (hmGlu}_5)$

ADX47273 $EC_{50} = 170 \text{ nM (hmGlu}_5)$

VU 1545 (Cat. No. 3325) EC₅₀=9.6 nM

EC₅₀=38 nM EC₅₀=5.9 nM X=CH MRZ 3573 X=NH cyclopentyl

$$R^2$$
 N N N X

VU 0360172 (Cat. No. 4323) $EC_{50} = 16 \text{ nM}$

2-aminomethyl-pyrimidine $EC_{50} = 14 \text{ nM}$

pyrido-oxazine $EC_{50} = 50 \text{ nM}$

VU 0357121 (Cat. No. 4437) $EC_{50} = 33 \text{ nM}$

G. mGlu₅ Receptor Neutral Modulators

DCB (Cat. No. 1952) IC_{50} =2.6 μ M for DFB potentiation attenuation

VU 0365396

group, exhibits slightly increased potency 5,77 in the same range as DCG-IV, which is also a group II agonist. 78 The best activity is found with the nonselective antagonist LY 341495. 79

Allosteric Modulators

Allosteric modulators are non-competitive ligands which bind in the transmembrane heptahelical domain. Both negative (NAMs) and positive modulators (PAMs) have been identified. RAMs inhibit receptor activation without affecting agonist binding while PAMs 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.

mGlu, Antagonists

Detailed studies have been devoted to CPCCOEt the first negative mGlu receptor modulator.^{22,81,82} In particular, specific residues of the HD that bind CPCCOEt were identified by a group from Novartis.²² Following this, other compounds with higher affinities were discovered by HTS and subsequent optimization, in various companies.83 These include: NPS 239084,85 (NPS Pharmaceuticals Inc.), Bay 36-762086 (Bayer AG), LY 45606687,88 and LY 45623689 (Eli Lilly), R21412785/JNJ 1625968590,91 (Johnson & Johnson), 3,5-dimethyl-pyrrole-2,4-dicarboxylic acid diesters (of which DM-PPP is the most potent derivative 92,93) (GlaxoSmithKline), several analogs of EM-TBPC94,95 (Hoffmann-La Roche), thiazolobenzimidazoles YM 298198,96 YM 20207497 and thienopyrimidine YM 230888 (Yamanouchi Pharma), triazafluorenones such as A 84172098 and more selective tetracyclic derivatives99 (Abbott Laboratories, Schering-Plough), CFMTI^{100,101} (Banyu Pharmaceutical Co.), pyrazines-2-carboxamides $PChPC^{102}$ and azaquinazolines such as CMPPA103 (Pfizer) and adamantyl methanone AdPyM¹⁰⁴ (Merz Pharmaceuticals). A homology model of the mGlu, allosteric binding site has been generated and a binding mode proposed for EM-TBPC which was validated by mutagenesis and functional assays.94 Additionally, it was shown that several inhibitors (R214127, CPCCOEt, NPS 2390, Bay 36-7620) bind to this same site.85 Promising anxiolytic and analgesic effects have been reported with allosteric mGlu, receptor antagonists; however potential side effects such as locomotor and cognition impairment were also discovered, impeding their development.83,105

mGlu₁ Positive Modulators

The first allosteric potentiators of rat ${\rm mGlu}_1$ receptors to be disclosed were Ro 01-6128, Ro 67-4853105, 107 and Ro 67-7476. 108,109 Chimeric and mutated receptors were constructed to confirm the transmembrane localization of the binding site of these ligands, which are subtype I selective. 108 Interestingly, Ro 67-7476 and Ro 01-6128 have little or no effect on human ${\rm mGlu}_1$ receptor activation whereas Ro 67-4853 produces a pronounced enhancement. 108 While CDPPB was known as an ${\rm mGlu}_5$ selective potentiator (see Figure 4F), VU 71 – which has the phenyl

substituent of the pyrazole core in the 4 rather than the 3 position – was discovered to be a selective mGlu1 potentiator, interacting with a site distinct from that of NAMs. 110

mGlu₅ Antagonists

SIB 1757 and SIB 1893¹¹¹ were initially found and optimized into MPEP¹¹² which has been widely used to explore the physiological roles of mGlu_s receptors as a potential therapeutic target. 113 Further investigations at Novartis led to a methoxy derivative M-MPEP, that can be easily tritium-labeled 114 and lead optimization resulted in AFQ056 (Mavoglurant) which has been in clinical trials for the symptomatic treatment of Parkinson's disease levodopa-induced dyskinesia (PD-LID) and Fragile X Syndrome. The therapeutic potential of mGlu_e antagonists prompted numerous groups to search for new ligands. 115-117 Early series contained an alkyne core while more recently extensive efforts focused on alternative chemotypes. MTEP, a thiazol derivative of MPEP with improved aqueous solubility, was described with similar high mGlu_s affinity¹¹⁸ as well as its tritiumlabeled methoxymethyl derivative MM-MTEP, 119,120 M-PEPy119 bipyridyl derivative MTEBP121 and fluorine derivatives for PET imaging (F-MTEB and SP203).122,123 Since these initial MPEP/ MTEP derivatives, 124 numerous disubstituted alkyne compounds have been described which include: ADX10059 (efficient for migraine and gastroesophageal reflux but also led to liver function abnormalities in patients); ADX48621¹¹⁷ (Dipraglurant), in phase II clinical trials for PD-LID; ethynylbenzamides (efficient in anxiety models);¹²⁵ ethynylpyrrolopyrazines;¹²⁶ MRZ 8676127 and CTEP, which displays high oral bioavailability and a long halflife of 18h.128 However, during development of the ethynyl series, it soon became apparent that minor structural changes unexpectedly modulated the pharmacology (a "molecular switch"), turning full NAMs into partial antagonists, PAMs or silent/neutral allosteric modulators (SAMs). 129,130 This is exemplified with 5-MPEP, where moving the methyl substituent of the MPEP pyridyl ring to the neighboring carbon turns this analog into a neutral modulator, 129 or with the 5-(phenylethynyl) pyrimidine series where the 3-methylphenyl derivative is a potent antagonist and the 4-methyl isomer a potentiator. 130

Several biosteric replacements of the alkyne core have been proposed: carboxamides, 132 arylquinolines, 133,134 heterocycles (e.g. tetrazole), 335 oxazolo-azepine or oxadiazole (VU 0285683).137 In parallel, HTS campaigns provided new scaffolds that were modulated into a plethora of chemical structures 116 for example aryl benzoxazoles138 (illustrated by BOMA), dipyridyl amides (ACDPP),139 phenyloxadiazoles and phenyltetrazoles,140 carbamoyloximes,141 thiazolotriazoles (such as GSK 2210875),142 pyrrolidinylpyridines, 143 piperidylamides, 144 benzimidazoles, 145 and anilinoquinazolines.146 In one of the HTS campaigns, it was found that the known anxiolytic drug fenobam was in fact a potent non-competitive mGlu₅ antagonist. 147 Based on this discovery, new derivatives were also developed.148 Potential therapeutic application of mGlu_r antagonists have been detailed in several reviews. 3,115,124,149 Additionally, molecular determinants of the high affinity binding site of MPEP have been defined 150 and a striking similarity with critical residues of the mGlu, binding site was observed.151

Figure 5 | Group II and Group III Allosteric Modulator Structures and Potencies

A. mGlu, Receptor Potentiators

$$O$$
 N
 O
 S
 CF_3

phenyl-tetrazolyl acetophenone PTBE EC_{50} =0.43 μM

oxazolidinone-1 EC₅₀=30 nM

oxazolidinone-2 EC₅₀=82 nM

BINA (Cat. No. 4048) EC₅₀=111 nM

TBPCOB EC₅₀=29 nM

benzimidazole-1 EC₅₀=30 nM

 $\begin{array}{ccc} & \text{benzimidazole-2} \\ \text{GSK 1331268} & \text{EC}_{50}\text{=}126 \text{ nM} \end{array}$

isoquinolone EC₅₀=250 nM

imidazopyridine EC_{50} =186 nM

pyridone (Cid et al 2010)¹⁸⁷ EC_{50} =525 nM

Imidazomethylpiperidine EC_{50} =35 nM

 $\begin{array}{c} \text{THIIC} \\ \text{EC}_{50} \text{=23 nM (hmGlu}_2) \end{array}$

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

Figure 5 | Group II and Group III Allosteric Modulator Structures and Potencies

B. $mGlu_{2/3}$ Receptor Antagonists

Ro 64-5229 (Cat. No. 2913) benzodiazepinone
$$IC_{50} = 109 \text{ nM}$$
 $IC_{50} = 34 \text{ nM}$ $IC_{50} = 34 \text{ nM}$ $IC_{50} = 2 \text{ nM}$ $IC_{50} = 2 \text{ nM}$

Ro 5488608

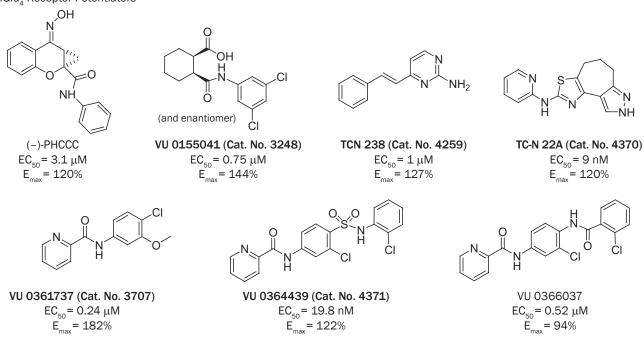
 $IC_{50} = 2.5 \text{ nM}$

Ro 4988546 IC₅₀= 8.7 nM

C. mGlu₂ Receptor Antagonists - mGlu₃ Receptor Agonists

	hmGlu ₂ K _i (μM)	hmGlu ₂ IC ₅₀ (μM)	hmGlu ₃ EC ₅₀ (μM)
R=H PHCCC (Cat. No. 1672)	N.E.	N.E.	N.E.
R=F	6.6	N.E. (SAM)	N.E. (SAM)
R=CI	1.0	0.8 (NAM)	13.4 (PAM)
R=Me	0.6	1.5 (NAM)	8.9 (PAM)
R=OMe	0.8	1.0 (NAM)	10.4 (PAM)

D. mGlu₄ Receptor Potentiators



phenylpicolinamide series

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

Figure 5 continued | Group II and Group III Allosteric Modulator Structures and Potencies

E. mGlu., Receptor Allosteric Agonists

AMN 082 (Cat. No. 2385) EC₅₀= 64 nM F. mGlu., Receptor Allosteric Antagonists

MMPIP (Cat. No. 2963)

G. mGlu Receptor PAM

AZ 12216052 (Cat. No. 4832) $EC_{50} = 11.0 \ \mu\text{M} \\ E_{max} = 77\%$

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

mGlu₅ Positive Modulators

Significant reports based on the first PAMs supported the development of mGlu_e potentiators as promising novel antipsychotics. 152 Consequently, they stimulated numerous research programs which have been conducted over the last few years. 152 The first mGlu_s PAMs to be identified were DFB, 153 CPPHA, 154,155 CDPPB 156,157 and ADX47273.158,159 As observed with mGlu_s NAMs, substituent modifications of mGlu_s PAMs led to a molecular switch: replacing the fluorine atoms of DFB by methoxy groups turns this ligand into an antagonist, while dichlorobenzaldazine (DCB) is a neutral modulator which attenuates the potentiation conferred by DFB.153 Similar modulations were found with close analogs of MPEP, as described above. 130,160 Acetylenic mGlu, PAM development led to MRZ 3573,¹⁵¹ VU 0360172¹³⁷ and to the 2-aminomethyl-pyrimidine phenylethynyl derivative that is a pure PAM, unlike several other mGlu_s PAMs that are 'ago-potentiators'. 131 Efforts to improve the metabolic stability of these PAMs resulted in the N-aryl piperazine (VU 0364289) 161,162 and piperidine amide series 162 and the phenoxymethyl pyridooxazines that are devoid of phenylacetylene and carbonyl functionalities. 163 Molecular switching 164 was also observed when building SAR around the ADX47273 structure¹⁶⁵ but not around CDPPB, 166 although moving a phenyl substituent changes the selectivity of VU 1545 (an mGlu, PAM) to VU 71 (an mGlu, PAM). A benzamide scaffold was also identified by HTS, chemical modulation led to the discovery of VU 0357121 but also to neutral (silent) modulators (such as VU 0365396).167 CPPHA and analogs appear to bind to a different site than MPEP while ethynyl PAM and NAM binding sites overlap.

Group II (Figure 5)

mGlu₂ Positive Modulators

The possible treatment of psychiatric diseases with mGlu $_2$ potentiators led to the launch of numerous research programs which resulted in the discovery of multiple modulators. 168,169 LY 487379, a pyridylmethylsulfonamide, was the first reported to potentiate the activity of glutamate at mGlu $_2$ receptors with an EC $_{50}$ value of 0.3 μM and to be highly selective for this subtype. 170 It was also demonstrated that LY 487379 binds to a pocket in the transmembrane domain which is different from the orthosteric site in the ATD. 170 Further SAR studies led to the discovery of

1-methylbutoxy analog (2,2,2-TEMPS) with improved potency (EC $_{50}$ = 14 nM) and selectivity. 171,172 Soon after, a new chemical series of phenyl-tetrazolyl acetophenones (e.g. PTBE) was disclosed as selective mGlu $_2$ potentiators, 173 followed by extensive SAR studies. $^{174-177}$ New chemotypes were later disclosed as a result of additional HTS hits and SAR studies. 178 Compounds presented here are mostly those selected among the series for *in vivo* assays and provide the best compromise between potency and metabolic stability: biphenylindanone (BINA), 179 recently optimized into benzizothiazolone, 180 benzimidazole- 181 and benzimidazole-2 (GSK 1331268), 182 oxazolidinone- 183 and oxazolidinone- 2184 optimized into oxazolobenzimidazoles (TBPCOB), 185 imidazopyridine, 186 1,5-disubstituted pyridine, 187 imidazole carboxamide (THIIC), 188 isoquinolones, 189 and imidazomethylpiperidine. 190

mGlu_{2/3} Antagonists

To date, only ${\rm mGlu_{2/3}}$ NAMs have been disclosed, mostly by researchers at Hoffmann-La Roche. Heterocyclic enol ethers such as Ro 64-5229 were reported as first selective noncompetitive ${\rm mGlu_2}$ receptor antagonists. A series of dihydrobenzo[b][1,4] diazepin-2-one derivatives was later disclosed, which exhibited nanomolar inhibition of receptor activation by LY 354740. Phis series was further improved in several derivatives, such as Ro 4491533 that was tested in vivo. More recently two novel antagonists, Ro 4988546 (from a new pyrazolo[1,5-a]pyrimidine scaffold) and Ro 5488608, were disclosed and used to investigate the structural determinant at the mGlu, NAM binding site. Phis such as Ro Hotelosed and site.

mGlu₂

A recent screening campaign provided specific mGlu₃ PAMs and NAMs, however the chemical structures were yet to be disclosed at the time of writing.¹⁹⁶ Interestingly, it was found that varying a substituent on the PHCCC structure resulted in a mGlu_{2/3} SAM or conferred dual mGlu₂ NAM - mGlu₃ PAM properties.¹⁹⁹

Group III

Group III modulators were the latest to be identified, mostly including mGlu₄ potentiators. PHCCC, which was initially described as an mGlu₄ receptor antagonist, 81 was the first mGlu₄

receptor PAM to be found as its (-) enantiomer.200,201 Two other mGlu_s antagonists, SIB 1893 and MPEP, were reported to enhance agonist potency and efficacy at human mGlu, at higher concentrations.202 Several mGlu, PAMs were subsequently discovered by HTS and hit optimization: VU 0155041;203 a series of phenylpicolinamides VU 0361737,²⁰⁴ VU 0364439,²⁰⁵ VU 366037;²⁰⁶ styryl aminopyrimidine;²⁰⁷ and thiazolopyrazole.^{208,209} Several of these ligands showed good brain penetration and benefits in motor dysfunction models but may possess intrinsic agonist activity as in the case of VU 0155041, and are therefore named ago-potentiators. 203 AMN 082 was described as an mGlu, allosteric agonist²¹⁰ however a recent study revealed a fast metabolism.²¹¹ Isoxazolpyridones such as MMPIP were determined as mGlu, antagonists^{212,213} but this effect may be context dependent.214 AZ 12216052, an mGlu PAM, was found to be systemically active in an animal model of anxiety.²¹⁵

Conclusion

In the early years, mGlu receptor molecular pharmacology efforts provided group selective competitive ligands. Although it now seems possible to discover subtype-selective orthosteric ligands, most of the recent advances 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 enabling investigators to shed 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.

List of Acronyms

A-841720 9-(Dimethylamino)-3-(hexahydro-1*H*-azepin-1-yl)pyrido[3',2':4,5]thieno[3,2-d]pyrimidin-4(3*H*)-one

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

ACDPP 3-Amino-6-chloro-5-dimethylamino-N-2-pyridinylpyrazinecarboxamide hydrochloride

AdPyM Adamantan-1-yl-[2-(6-morpholin-4-yl-2-pyridin-3-yl)-cyclopropyl]-methanone

ADX10059 2-((3-Fluorophenyl)ethynyl)-4,6-dimethylpyridin-3-amine

ADX48621 6-Fluoro-2-[4-(pyridin-2-yl)but-3-yn-1-yl]imidazo[1,2-a]pyridine (dipraglurant)

ADX47273 (S)-(4-Fluorophenyl)-[3-[3-(4-fluorophenyl)-[1,2,4]oxadiazol-5-yl]-piperidin-1-yl}-methanone AFQ056 (3aR,4S,7aR)-Methyl 4-hydroxy-4-(m-tolylethynyl)octahydro-1H-indole-1-carboxylate

AMN 082 *N,N'*-Bis(diphenylmethyl)-1,2-ethanediamine

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-phosphonobutyric acid

APCPr 1-Amino-2-(phosphonomethyl)cyclopropane carboxylic acid

APDC 4-Aminopyrrolidine-2,4-dicarboxylic acid

AZ 12216052 2-(4-Bromobenzylthio)-N-(4-sec-butylphenyl)acetamide

Bay 36-7620 (3aS,6aS)-6a-Naphtalen-2-ylmethyl-5-methyliden-hexahydro-cyclopental[c]furan-1-one

BINA 3'-((2-Cyclopentyl-6,7-dimethyl-1-oxo-2,3-dihydro-1*H*-inden-5-yloxy)methyl)biphenyl-4-carboxylic acid

BnAPDC N-Benzyl-(2R,4R)-4-aminopyrrolidine-2,4-dicarboxylic acid

BnQuis α -Benzylquisqualic acid

BOMA 2-[4-(1,3-Benzoxazol-2-yl)-2-methoxyphenyl]acetonitrile

L-CCG-I (2S, 1'S, 2'S)-2-(Carboxycyclopropyl)glycine

3'Me-CCG (2S,1'S,2'S,3'R)-2-(2'-Carboxy-3'-methylcyclopropyl)glycine

3'HM-CCG (2S,1'S,2'R,3'R)-2-(2'-Carboxy-3'-hydroxymethylcyclopropyl)glycine mCD-CCG 2-[2',2'-di(3-Chlorophenyl)ethyl]-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-1*H*-pyrazol-5-yl)benzamide

CHPG 2-Chloro-5-hydroxyphenylglycine

4C3H2MPG 4-Carboxy-3-hydroxy-2-methylphenylglycine
4C2MPG (+)-4-Carboxy-2-methylphenylglycine

4CPG 4-Carboxyphenylglycine

 $\hbox{CFMTI} \qquad \hbox{2-Cyclopropyl-5-[1-(2-fluoro-3-pyridinyl)-5-methyl-1$H-1,2,3-triazol-4-yl]-2,3-dihydro-1$H-isoindol-1-one } \\$

CMPPA N-Cycloheptyl-6-(2-morpholinoethoxy)pyrido[3,4-d]pyrimidin-4-amine

(-)-CPCCOEt (1aS,7aS)-(2-Hydroxyimino-1a,2-dihydro-1H-7-oxacyclopropa[b]naphthalene-7a-carboxylic acid ethyl ester

 ${\it CPPG} \hspace{1cm} \alpha\hbox{-cyclopropyl-4-phosphonophenylglycine}$

CPPHA N-[5-Chloro-2-[(1,3-dioxoisoindolin-2-yl)methyl]phenyl]-2-hydroxybenzamide CPPZ 1-(4-(2-Chloro-4-fluorophenyl)piperazin-1-yl)-2-(pyridin-4-ylmethoxy)ethanone

 $\hbox{CTEP} \qquad \hbox{2-Chloro-4-((2,5-dimethyl-1-(4-(trifluoromethoxy)phenyl)-1$$H$-imidazol-4-yl)ethynyl) pyridined for the properties of the p$

DCG-IV (2S, 1'R, 2'R)-2-(2',3'-Dicarboxycyclopropyl)glycine

3,4-DCPG3,4-Dicarboxyphenylglycine3,5-DHPG3,5-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-trimethyl-propyl)ester

EM-TBPC 1-Ethyl-2methyl-6-oxo-4-(1,2,4,5-tetrahydro-benzo[d]azepin-3-yl)-1,6-dihydro-pyrimidine-5-carbonitrile

FP429 (2S,4S)-4-Amino-1-[(E)-3-carboxyacryloyl]pyrrolidine-2,4-dicarboxylic acid

L-Glu L-Glutamate

GSK 1331268 2-((4-(5-Chloropyridin-2-yl)piperazin-1-yl)methyl)-1-methyl-1*H*-benzo[d]imidazole

GSK 2210875 (R)-1-(6-Methylthiazolo[3,2-b][1,2,4]triazol-5-yl)ethyl phenylcarbamate

HYDIA (1S,2R,3R,5R,6S)-3-Hydroxy-2-aminobicyclo[3.1.0]hexane-2,6-dicarboxylic acid
JNJ 16259685 (3,4-Dihydro-2*H*-pyrano[2,3-*b*]quinolin-7-yl)-(*cis*-4-methoxycyclohexyl)-methanone

LSP1-2111 [((3S)-3-Amino-3-carboxy)propyl][(4-hydroxy-5-methoxy-3-nitrophenyl)hydroxymethyl]phosphinic acid

LSP1-3081 [(3S)-3-(3-Amino-3-carboxypropyl(hydroxy)phosphinyl)-hydroxymethyl]-5-nitrothiophene

LSP4-2022 [((3S)-3-Amino-3-carboxy)propyl)][(4-(carboxymethoxy)phenyl)hydroxymethyl]phosphinic acid

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 367385 (4C2MPG) (+)-4-Carboxy-2-methylphenylglycine

LY 379268 2-0xa-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 \ (+/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(9-thioxanthyl) propionic \ acid \ (-/-)-2-Amino-2-(3-cis \ and \ trans-carboxycyclobutyl-3-(3-cis \ and \ trans-carboxycyclobutyl-3$

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-tetrahydro-quinazolin-2-ylsulfanyl]-ethanol

LY 456236 6-Methoxy-N-(4-methoxyphenyl)-4-quinazolinamine

LY 487379 N-(4-(2-Methoxyphenoxy)phenyl)-N-(2,2,2-trifluoroethylsulfonyl)pyrid-3-ylmethylamine

LY 541850 (1S,2S,3S,5R,6S)-2-Amino-3-methylbicyclo[3.1.0]hexane-2,6-dicarboxylic acid

LY 2812223 (1S,2S,3S,5R,6S)-3-(1*H*-1,2,4-Triazol-3-ylthio)-2-aminobicyclo[3.1.0]hexane-2,6-dicarboxylic acid

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 MGS0022 (1R,2S,5R,6R)-2-Amino-6-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

MPPG α -Methyl-4-phosphonophenylglycine MPEP 2-Methyl-6-(phenylethynyl)pyridine

M-MPEP 2-[(3-Methoxyphenyl)ethynyl]-6-methylpyridine

MTEB 5-[(2-Methyl-1,3-thiazol-4-yl)ethynyl]-benzonitrile

F-MTEB 3-Fluoro-5-[(2-methyl-1,3-thiazole-4-yl)ethynyl]-benzonitrile

MTEBP 5-[(2-Methyl-1,3-thiazol-4-yl)ethynyl]-2,3'-bipyridine

MTEP 3-[(2-Methyl-1,3-thiazol-4-yl)ethynyl]pyridine

 $\label{eq:mm-map} \mbox{MM-MTEP} \qquad \qquad \mbox{3-(Methoxymethyl)-5-[(2-methyl-1,3-thiazol-4-yl)-ethynyl]pyridine}$

MMPIP 6-(4-Methoxyphenyl)-5-methyl-3-(pyridin-4-yl)isoxazolo[4,5-c]pyridin-4(5H)-one

M-PEPy 3-Methoxy-5-(pyridin-2-ylethynyl)pyridine

MRZ 3573 2-(Phenylethynyl)-7,8-dihydroguinolin-5(6H)-one

MRZ 8676 6,6-Dimethyl-2-(phenylethynyl)-7,8-dihydroquinolin-5(6*H*)-one

Tocris Scientific Review Series

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
PBPG (2S)-2-(3'-Phosphonobicyclo[1.1.1]pentyl)glycine

PCEP 3-Amino-3-carboxypropyl-2'-carboxyethyl phosphinic acid PCG-1 trans-(2S,1'R,2'S)-2-(2'-Phosphonocyclopropyl) glycine

PChPC 5-(4-(Hydroxymethyl)piperidin-1-yl)-N-(trans-4-methylcyclohexyl)pyrazine-2-carboxamide

PPG 4-Phosphonophenylglycine

PHCCC N-Phenyl-7-(hydroxyimino)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-(2*H*-tetrazol-5-yl)phenoxy]butoxyphenyl)ethanone

Quis Quisqualate

R214127 1-(3,4-Dihydro-2*H*-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)vinyl]-(1,2,4-triazole)

Ro 67-4853 (9*H*-Xanthene-9-carbonyl)-carbamic acid butyl ester
Ro 67-7476 (S)-2-(4-Fluoro-phenyl)-1-(toluene-4-sulfonyl)-pyrrolidine

Ro 4988546 5-[7-Trifluoromethyl-5-(4-trifluoromethyl-phenyl)-pyrazolo[1,5-a]pyrimidin-3-ylethynyl]-pyridine-3-sulphonic acid Ro 5488608 3'-(8-Methyl-4-oxo-7-trifluoromethyl-4,5-dihydro-3*H*-benzo[b][1,4]diazepin-2-yl)-biphenyl-3-sulphonic acid

Ro 645229 (Z)-1-(2-(Cycloheptyloxy)-2-(2,6-dichlorophenyl)vinyl)-1*H*-1,2,4-triazole

SIB1757 6-Methyl-2-(phenylazo)-3-pyridinol

SIB1893 (E)-2-Methyl-6-(2-phenylethenyl)pyridine

SOP Serine-O-phosphate

SP203 3-Fluoro-5-[[2-(fluoromethyl)thiazol-4-yl]ethynyl]-benzonitrile

TBPCOB (S)-2-((6-tert-Butylpyridin-3-yloxy)methyl)-2,3-dihydrobenzo[d]oxazolo[3,2-a]imidazole-7-carbonitrile

TC-N 22A 4,5,6,8-Tetrahydro-*N*-2-pyridinylpyrazolo[3',4':6,7]cyclohepta[1,2]thiazol-2-amine

TCN 238 (E)-4-(2-Phenylethenyl)-2-pyrimidinamine

2,2,2-TEMPS 2,2,2-Trifluoro-*N*-(4-(4-(pentan-2-yl)phenoxy)phenyl)-*N*-(pyridin-3-ylmethyl)ethanesulfonamide

 $\label{eq:hamiltonian} THIIC \qquad \qquad N-(4-((2-(Trifluoromethyl)-3-hydroxy-4-(isobutyryl)phenoxy)methyl)benzyl)-1-methyl-1<math>H$ -imidazole-4-carboxamide

VU 71 4-Nitro-*N*-(1,4-diphenyl-1*H*-pyrazol-5-yl)benzamide

VU 1545 4-Nitro-N-(1-(2-fluorophenyl)-3-phenyl-1*H*-pyrazol-5-yl)benzamide
VU 0155041 *cis*-2-(3,5-Dichlorphenylcarbamoyl)cyclohexanecarboxylic acid
VU 0285683 3-Fluoro-5-(3-(pyridine-2-yl)-1,2,4-oxadiazol-5-yl)benzonitrile

VU 0357121 4-Butoxy-N-(2,4-difluorophenyl)benzamide

VU 0360172 N-Cyclobutyl-6-((3-fluorophenyl)ethynyl)nicotinamide

VU 0361737 N-(4-Chloro-3-methoxyphenyl)picolinamide

VU 0364439 N-(3-Chloro-4-(N-(2-chlorophenyl)sulfamoyl)phenyl)picolinamide

VU 366037 N-(3-Chloro-4-(2-chlorobenzamido)phenyl)picolinamide
VU 0364289 2-(4-(2-(Benzyloxy)acetyl)piperazin-1-yl)benzonitrile

VU 0365396 4-Butoxy-N-(2,6-difluorophenyl)benzamide

YM 202074 N-Cyclohexyl-6-{[(2-methoxyethyl)(methyl)amino]methyl}-N-methylthiazolo[3,2-a]benzimidazole-2-carboxamide

 $YM\ 230888 \qquad \qquad (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methyl]amino]methyl) thieno[2,3-d]pyrimidin-4-ylamine (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methyl]amino]methyl (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methyl]amino[2,3-d]pyrimidin-4-ylamine (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methyl]amino[2,3-d]pyrimidin-4-ylamine (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methyl]amino[2,3-d]pyrimidin-4-ylamine (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methyl]amino[2,3-d]pyrimidin-4-ylamine (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methyl]amino[2,3-d]pyrimidin-4-ylamine (R)-N-Cycloheptyl-6-([(tetrahydro-2-furyl)methy$

References

- 1. Dingledine et al. (1999) Pharmacol. Rev. 51 7.
- 2. Pin and Acher (2002) Curr.Drug Targets: CNS Neurol.Disord. 1 297.
- Niswender and Conn (2010) Annu. Rev. Pharmacol. Toxicol. 50 295.
- Recasens et al. (2007) Curr.Drug Targets 8 651.
- 5. Schoepp et al. (1999) Neuropharmacology 38 1431.
- 6. Pin et al. (1999) Eur.J.Pharmacol. 375 277.
- 7. Gasparini and Spooren (2007) Curr.Neuropharmacol. 5 187.
- 8. Urwyler (2011) Pharmacol. Rev. 63 59.
- 9. Pin et al. (2003) Pharmacol. Ther. 98 325.
- 10. O'Hara et al. (2003) Neuron 11 41.
- 11. Bessis et al. (2000) Protein Sci. 9 2200.
- 12. Kunishima et al. (2000) Nature 407 971.
- 13. Muto et al. (2007) Proc.Natl.Acad.Sci.U.S.A. 104 3759.
- 14. Bessis et al. (2002) Proc.Natl.Acad.Sci.U.S.A. 99 11097.
- 15. Hampson et al. (1999) J.Biol.Chem. 274 33488.
- 16. Malherbe et al. (2001) Mol. Pharmacol. 60 944.
- 17. Bertrand et al. (2002) J.Med.Chem. 45 3171.
- 18. Rosemond et al. (2002) J.Biol.Chem. 277 7333.
- 19. Rosemond et al. (2004) Mol. Pharmacol. 66 834.
- 20. Acher and Bertrand (2005). Biopolymers 80 357.
- 21. May and Christopoulos (2003) Curr.Opin.Pharmacol. 3 551.
- 22. Litschig et al. (1999) Mol. Pharmacol. 55 453.
- 23. Kew (2004) Pharm.Ther. 104 233.
- Conn and Pin (1997) Ann. Rev. Pharmacol. Toxicol. 37 205.
- Bessis et al. (1999) Neuropharmacology 38 1543.
- 26. Monn et al. (1997) J.Med.Chem. 40 528.
- 27. Dunayevich et al. (2008) Neuropsychopharmacology 33 1603.
- 28. Patil et al. (2007) Nat. Med. 13 1102.
- 29. Mezler et al. (2010) Curr.Opin.Investig.Drugs 11 833.
- 30. Johnson et al. (2002) Drug Metab.Disp. 30 27.
- 31. Lennon et al. (2010) Eur.J.Pharmacol. 649 29.
- 32. Palmer et al. (1989) Eur.J.Pharmacol. 166 585.
- 33. Doherty et al. (1997) Neuropharmacology 36 265.
- 34. Littman et al. (1999) J.Med.Chem. 42 1639.
- 35. Monn et al. (1999) J.Med.Chem. 42 1027.
- 36. Monn et al. (2007) J.Med.Chem. 50 233.
- 37. Nakazato et al. (2000) J.Med.Chem. 43 4893.
- 38. Collado et al. (2002) J.Med.Chem. 45 3619.
- 39. Collado et al. (2004) J.Med.Chem. 47 456.
- 40. Stanley et al. (2010) Bioorg. Med. Chem. 18 6089.
- 41. Gonzalez et al. (2005) Bioorg. Med. Chem. 13 6556.
- 42. Dominguez et al. (2005) J.Med.Chem. 48 3605.
- 43. Monn et al. (2011) Curr. Neuropharmacol. 9 Suppl. 1, 44.
- Chopra et al. (2009) J.Pharmacol.Exp.Ther. 330 212.
- Selvam et al. (2007) J.Med.Chem. 50 4656.
- 46. Kroona et al. (1991) J. Med. Chem. 34 1692.
- 47. Sibille et al. (2007) J.Med.Chem. 50 3585.
- 48. Gasparini et al. (1999) J.Pharm.Exp.Ther. 290 1678.
- 49. Gasparini et al. (2000) Bioorg.Med.Chem.Lett. 10 1241.
- 50. Thomas et al. (2001) Neuropharmacology 40 311.
- 51. Acher et al. (1997) J.Med.Chem. 40 3119.
- 52. Filosa et al. (2006) Bioorg. Med. Chem. 14 3811.
- 53. Amori et al. (2006) Bioorg. Med. Chem. Lett. 16 196.
- 54. Selvam et al. (2010) J.Med.Chem. 53 2797.
- 55. Triballeau et al. (2005) J.Med.Chem. 48 2534.
- 56. Selvam et al. (2011) submitted.
- 57. Cuomo et al. (2009) J.Neurochem. 109 1096.
- 58. Beurrier et al. (2009) FASEB J. 23 3619.
- 59. Wieronska et al. (2010) Neuropharmacology 59 627.
- 60. Goudet et al. (2011) submitted.
- 61. Schann et al. (2006) Bioorg. Med. Chem. Lett. 16 4856.

- 62. Frauli et al. (2007) Mol. Pharmacol. 71 704.
- 63. Tückmantel et al. (1997) Bioorg. Med. Chem. Lett. 7 601.
- Ahmadian et al. (1997) J.Med.Chem. 40 3700.
- Fazio et al. (2012) Mol. Pharmacol. 81 643.
- Chen et al. (2000) Neurosci. 95 787.
- 67. Kingston et al. (2002) Neurosci.Lett. 330 127.
- Melendez et al. (2005) J.Pharmacol.Exp.Ther. 314 139. 68.
- 69. Sakagami et al. (2008) Bioorg. Med. Chem. 16 4359.
- 70. Pellicciari et al. (2001) Bioorg. Med. Chem. Lett. 11 3179.
- 71. Sørensen et al. (2003) Bioorg. Med. Chem. 11 197.
- 72. Woltering et al. (2008) ChemMedChem 3 323.
- 73. Lundstrom et al. (2009) ChemMedChem 4 1086.
- 74. Chaki et al. (2004) Neuropharmacology 46, 457.
- 75. Nakazato et al. (2004) J.Med.Chem. 47 4570.
- 76. Yasuhara et al. (2006) Bioorg. Med. Chem. 14 3405.
- 77. Conway et al. (2001) Bioorg. Med. Chem. Lett. 11 777.
- 78. Brabet et al. (1998) Neuropharmacology 37 1043.
- 79. Wright et al. (2000) Naunyn Schmiedebergs Arch. Pharmacol. 362 546.
- 80. Sabbatini and Micheli (2004) Expert Opin. Ther. Patents 14 1593.
- 81. Annoura et al. (1996) Bioog. Med. Chem. Lett. 6 763.
- 82. Ott et al. (2000) J.Med.Chem. 43 4428.
- 83. Owen (2011) ACS.Chem.Neurosci. 2 394.
- 84. Van Wagenen et al. (1998) Society for Neuroscience Abstract 24 576.
- Lavreysen et al. (2003) Mol. Pharmacol. 63 1082.
- Caroll et al. (2001) Mol. Pharmacol. 59 965.
- 87. Li et al. (2002) Neuropharmacology 43 A79.
- Ambler and Baker (2001) Patent WO 01/32632. 88.
- Shannon et al. (2005) Neuropharmacology 49 188. 89.
- 90. Lavrevsen et al. (2004) Neuropharmacology 47 961.
- 91. Mabire et al. (2005) J.Med.Chem. 48 2134.
- 92. Micheli et al. (2003) Bioorg. Med. Chem. 11 171.
- 93. Di Fabio et al. (2007) Bioorg. Med. Chem. Lett. 17 2254.
- 94. Malherbe et al. (2003) J.Biol.Chem. 278 8340.
- 95. Binggeli et al. (2002) Patent W002051418.
- 96. Kohara et al. (2005) J.Pharmacol.Exp.Ther. 315 163.
- 97. Kohara et al. (2008) Brain Res. 1191 168.
- 98. Zheng et al. (2005) J.Med.Chem. 48 7374.
- 99. Wu et al. (2007) J.Med.Chem. 50 5550. 100. Ito et al. (2009) Bioorg. Med. Chem. Lett 19 5310.
- 101. Satow et al. (2009) J.Pharmacol.Exp.Ther. 330 179.
- 102. Owen et al. (2007) Bioorg. Med. Chem. Lett. 17 486.
- 103. Mantell et al. (2009) Bioorg. Med. Chem. Lett. 19 2190.
- 104. Noeske et al. (2009) Bioorg. Med. Chem. 17 5708.
- 105. Lesage and Steckler (2010) Eur.J.Pharmacol. 639 2.
- 106. Bleicher et al. (2000) Patent W00063166.
- 107. Vieira et al. (2009) Bioorg. Med. Chem. Lett. 19 1666.
- 108. Knoflach et al. (2001) Proc.Nat.Acad.Sci. 98 13402.
- 109. Wichmann et al. (2002) Farmaco. 57 989.
- 110. Hemstapat et al. (2006) Mol. Pharmacol. 70 616.
- 111. Varney et al. (1999) J.Pharmacol.Exp.Ther. 290 170. 112. Gasparini et al. (1999) Neuropharmacology 38 1493.
- 113. Spooren et al. (2001) T.I.P.S. 22 331.
- 114. Gasparini et al. (2002) Bioorg. Med. Chem. Lett. 12 407.
- 115. Jaeschke et al. (2008) Expert. Opin. Ther. Pat. 18 123.
- 116. Emmitte (2011) ACS.Chem.Neurosci. 2 411.
- 117. Rocher et al. (2011) Curr.Top Med.Chem. 11 680.
- 118. Cosford et al. (2003) J.Med.Chem. 46 204.
- 119. Cosford et al. (2003) Bioorg. Med. Chem. Lett. 13 351.
- 120. Anderson et al. (2002) J.Pharmacol.Exp.Ther. 303 1044. 121. Roppe et al. (2004) Bioorg. Med. Chem. Lett. 14 3993.
- 122. Simeon et al. (2007) J.Med.Chem. 50 3256.

Tocris Scientific Review Series

- 123. Simeon et al. (2011) J.Med.Chem. 54 901.
- 124. Slassi et al. (2005) Curr.Top.Med.Chem. 5 897.
- 125. Gilbert et al. (2011) Bioorg. Med. Chem. Lett. 21 195.
- 126. Micheli et al. (2008) Bioorg. Med. Chem. Lett. 18 1804.
- 127. Dekundy et al. (2010) J.Neural Transm. 118 1703.
- 128. Lindemann et al. (2011) J.Pharmacol.Exp.Ther. 339 474.
- 129. Rodriguez et al. (2005) Mol. Pharmacol. 68 1793.
- 130. Wood et al. (2011) Biochemistry 50 2403.
- 131. Sharma et al. (2009) J.Med.Chem. 52 4103.
- 132. Kulkarni et al. (2009) J.Med.Chem. 52 3563.
- 133. Milbank et al. (2007) Bioorg. Med. Chem. Lett. 17 4415.
- 134. Zhang et al. (2010) Bioorg.Med.Chem. 18 3026.
- 135. Roppe et al. (2004) J.Med.Chem. 47 4645.
- 136. Burdi et al. (2010) J.Med.Chem. 53 7107.
- 137. Rodriguez et al. (2010) Mol. Pharmacol. 78 1105.
- 138. Wang et al. (2004) Bioorg. Med. Chem. 12 17.
- 139. Bonnefous et al. (2005) Bioorg.Med.Chem.Lett. 15 1197.
- 140. Wagner et al. (2010) Bioorg. Med. Chem. Lett. 20 3737.
- 141. Galambos et al. (2010) Bioorg. Med. Chem Lett. 20 4371.
- 142. Pilla et al. (2010) Bioorg. Med. Chem. Lett. 20 7521.
- 143. Weiss et al. (2011) Bioorg.Med.Chem.Lett. 21 4891.
- 144. Spanka et al. (2010) Bioorg. Med. Chem. Lett. 20 184.
- 145. Carcache et al. (2011) ACS.Med.Chem.Lett. 2 58.
- 146. Felts et al. (2009) Bioorg. Med. Chem. Lett. 19 6623.
- 147. Porter et al. (2005) J.Pharmacol.Exp.Ther. 315 711.
- 148. Jaeschke et al. (2007) Bioorg. Med. Chem. Lett. 17 1307.
- 149. Gasparini et al. (2008) Curr.Opin.Drug Discov.Devel. 11 655.
- 150. Malherbe et al. (2003) Mol. Pharmacol. 64 823.
- 151. Vanejevs et al. (2008) J.Med.Chem. 51 634.
- 152. Stauffer et al. (2011) ACS.Chem.Neurosci. 2 450.
- 153. O'Brien et al. (2003) Mol. Pharmacol. 64 731.
- 154. O'Brien et al. (2004) J.Pharmacol.Exp.Ther. 309 568.
- 155. **Zhao** et al. (2007) Bioorg.Med.Chem.Lett. **17** 1386.
- 156. Lindsley et al. (2004) J.Med.Chem. 47 5825.
- 157. Kinney et al. (2005) J.Pharmacol.Exp.Ther. 313 199.
- 158. Le Poul et al. (2005) Neuropharmacology 49 252.
- 159. Liu et al. (2008) J.Pharmacol.Exp.Ther. 327 827.
- 160. Sams et al. (2011) Bioorg. Med. Chem. Lett. 21 3407.
- 161. **Zhou** et al. (2010) ACS.Chem.Neurosci. **1** 433.
- 162. Xiong et al. (2010) Bioorg. Med. Chem. Lett. 20 7381.
- 163. Varnes et al. (2011) Bioorg.Med.Chem.Lett. 21 1402.
- 164. Lamb et al. (2011) Bioorg.Med.Chem.Lett. 21 2711.
- 165. Engers et al. (2009) ChemMedChem 4 505.
- 166. de Paulis et al. (2006) J.Med.Chem. 49 3332.
- 167. Hammond et al. (2010) ACS.Chem.Neurosci. 1 702.
- 168. Trabanco et al. (2011) Curr.Med.Chem. 18 47.
- 169. Sheffler et al. (2011) ACS.Chem.Neurosci. 2 282.
- 170. Schaffhauser et al. (2003) Mol.Pharmacol. 64 798.
- 171. Barda et al. (2004) Bioorg.Med.Chem.Lett. 14 3099.
- 172. Hu et al. (2004) Bioorg. Med. Chem. Lett. 14 5071.
- 173. Pinkerton et al. (2004) J.Med.Chem. 47 4595.
- 174. Pinkerton et al. (2004) Bioorg. Med. Chem. Lett. 14 5867.
- 175. Pinkerton et al. (2005) Bioorg.Med.Chem.Lett. 15 1565.
- 176. **Cube** et al. (2005) Bioorg.Med.Chem.Lett. **15** 2389.
- Govek et al. (2005) Bioorg.Med.Chem.Lett. 15 4068.
 Fraley (2009) Expert Opin.Ther.Pat. 19 1259.
- 179. Bonnefous et al. (2005) Bioorg Med Chem Lett. 15 4354.
- 180. Dhanya et al. (2011) J.Med.Chem. 54 342.
- 181. **Zhang** et al. (2008) Bioorg.Med.Chem.Lett. **18** 5493.
- 182. **D'Alessandro** et al. (2010) Bioorg.Med.Chem.Lett. **20** 759.
- Duplantier et al. (2009) Bioorg.Med.Chem.Lett. 19 2524.
 Brnardic et al. (2010) Bioorg.Med.Chem.Lett. 20 3129.
- 185. Garbaccio et al. (2010) ACS.Med.Chem.Lett. 1 406.

- 186. Tresadern et al. (2010) Bioorg. Med. Chem. Lett. 20 175.
- 187. Cid et al. (2010) ACS.Chem.Neurosci. 1 788.
- 188. Fell et al. (2011) J.Pharmacol.Exp.Ther. 336 165.
- 189. **Trabanco** et al. (2011) Bioorg.Med.Chem.Lett. **21** 971.
- 190. Zhang et al. (2011) J.Med.Chem. 54 1724.
- 191. Kolczewski et al. (1999) Bioorg. Med. Chem. Lett. 9 2173.
- 192. Woltering et al. (2007) Bioorg. Med. Chem. Lett. 17 6811.
- 193. Woltering et al. (2008) Bioorg.Med.Chem.Lett. 18 1091.
- 194. Woltering et al. (2008) Bioorg.Med.Chem.Lett. 18 2725.
- 195. Woltering et al. (2010) Bioorg.Med.Chem.Lett. 20 6969.
- 196. Hemstapat et al. (2007) J.Pharmacol.Exp.Ther. 322 254.
- 197. Lundstrom et al. (2011) Br.J.Pharmacol.
- 198. Pratt et al. (2011) Comb.Chem.High Throughput Screen 14 631.
- 199. Schann et al. (2010) J.Med.Chem. 53 8775.
- 200. Maj et al. (2003) Neuropharmacology 45 895.
- 201. Marino et al. (2003) Proc.Natl.Acad.Sci.U.S.A. 100 13668.
- 202. Mathiesen et al. (2003) Br.J.Pharmacol. 138 1026.
- 203. Niswender et al. (2008) Mol. Pharmacol. 74 1345.
- 204. Engers et al. (2009) J.Med.Chem. 52 4115.
- 205. Engers et al. (2010) Bioorg. Med. Chem. Lett. 20 5175.
- 206. Engers et al. (2011) J.Med.Chem. **54** 1106.
- 207. East et al. (2010) Bioorg. Med. Chem. Lett. 20 4901
- 208. East and Gerlach (2010) Expert Opin. Ther. Pat. 20 441.
- 209. Hong et al. (2011) J.Med.Chem. 54 5070.
- 210. Mitsukawa et al. (2005) Proc.Natl.Acad.Sci.U.S.A. **102** 18712.
- 211. Sukoff Rizzo et al. (2011) J.Pharmacol.Exp.Ther. 338 345.
- 212. **Suzuki** et al. (2007) J.Pharmacol.Exp.Ther. **323** 147.
- 213. Nakamura et al. (2010) Bioorg.Med.Chem.Lett. **20** 726.
- 214. **Niswender** et al. (2010) Mol.Pharmacol. **77** 459.
- 215. **Duvoisin** et al. (2010) Behav.Brain Res. **212** 168.
- 216. Adam et al. (1999) Neuropharmacology 38 A1 abstract 3.
- 217. Naples and Hampson (2001) Neuropharmacology 40 170.

Metabotropic Glutamate Receptor Compounds Available from Tocris

Cat. No.	Product Name	Primary Action
Group I mGI	u Receptors	
Agonists		
0284	(1S,3R)-ACPD	Group I and group II mGlu agonist; active isomer of (±)-trans-ACPD (Cat. No. 0187)
0187	(±)-trans-ACPD	Group I and II mGlu agonist
1049	CHPG	mGlu ₅ selective agonist
3695	CHPG sodium salt	Selective mGlu ₅ agonist; sodium salt of CHPG (Cat. No. 1049)
0342	(RS)-3,5-DHPG	Selective group I mGlu agonist
0805	(S)-3,5-DHPG	Selective group I mGlu agonist; active enantiomer of (RS)-3,5-DHPG (Cat. No. 0342)
0188	L-Quisqualic acid	Group I mGlu agonist; also AMPA agonist
0162	S-Sulfo-L-cysteine sodium salt	Group I agonist
Inverse Agonis	sts	
2501	Bay 36-7620	mGlu ₁ inverse agonist
Antagonists		
5229	ABP 688	High affinity human mGlu _s antagonist
0904	AIDA	Potent and selective group I mGlu antagonist
5614	AZD 2066	mGlu ₅ antagonist; orally bioavailable and brain penetrant
5613	AZD 9272	Potent and selective mGlu ₅ antagonist; brain penetrant
0125	DL-AP3	Group I mGlu antagonist
0329	(S)-3-Carboxy-4-hydroxyphenylglycine	Group I mGlu antagonist; also group II mGlu agonist
0323	(S)-4-Carboxyphenylglycine	Competitive group I mGlu antagonist; also weak group II agonist
1028	CPCCOEt	Selective non-competitive mGlu, antagonist
2333	JNJ 16259685	Highly potent, mGlu ₁ -selective non-competitive antagonist
1237	LY 367385	Selective mGlu _{1a} antagonist
2390	LY 456236	Selective mGlu, antagonist
2196	3-MATIDA	Potent and selective mGlu ₁ antagonist
1212	MPEP	Potent mGlu ₅ antagonist; also positive allosteric modulator at mGlu ₄ receptors
2921	MTEP	Potent and selective mGlu ₅ antagonist
4134	NPS 2390	Group I mGlu antagonist
1027	PHCCC	Potent group I mGlu antagonist
1215	SIB 1757	Highly selective mGlu _s antagonist
2986	YM 230888	Selective mGlu ₁ antagonist
2448	YM 298198	Highly potent, selective non-competitive mGlu ₁ antagonist
Modulators		
3235	CDPPB	Positive allosteric modulator of mGlu _s receptors
1952	DCB	Allosteric potentiator of mGlu ₅ receptors
1625	DFB	Positive allosteric modulator of mGlu _s receptors
5000	FTIDC	Potent and selective negative allosteric modulator of ${\rm mGlu}_{\scriptscriptstyle 1}$ receptors; also ${\rm mGlu}_{\scriptscriptstyle 1}$ inverse agonist
5275	LSN 2463359	Potent and selective positive allosteric modulator of $\mathrm{mGlu}_{\scriptscriptstyle{5}}$ receptors
4348	Ro 01-6128	Positive allosteric modulator of mGlu ₁ receptors
4347	Ro 67-4853	Positive allosteric modulator of group I mGlu receptors
4346	Ro 67-7476	Positive allosteric modulator of mGlu ₁ receptors
4323	VU 0360172	Positive allosteric modulator of mGlu ₅ receptors
5693	VU 0409551	Selective positive allosteric modulator of $\mathrm{mGlu}_{\scriptscriptstyle{5}}$ receptors; brain penetrant and orally bioavailable
	V/U 04606E0	Potent and selective negative allosteric modulator of mGlu, receptors
5379	VU 0469650	Total and selective negative anosteric modulator of main, receptors
5379	VU 0483605	Positive allosteric modulator of mGlu ₁ receptors

Group II mGli	u Recentors	
Activators	a recopiors	
4120	Xanthurenic acid	Selectively activates group II mGlu receptors
Agonists	7.6	Colocatory destructed group it make recorption
0187	(±)-trans-ACPD	Group I and II 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
0333	L-CCG-I	Potent group II mGlu agonist
0975	DCG IV	Highly potent group II mGlu agonist; also NMDA agonist
3246	LY 354740	Potent and highly selective group II mGlu agonist
2453	LY 379268	Highly selective group II mGlu agonist
5064	LY 379268 disodium salt	Selective group II mGlu agonist; sodium salt of LY 379268 (Cat. No. 2453)
0391	Spaglumic acid	Selective mGlu ₃ agonist
Antagonists		Concerno mana ₃ agomet
1073	(RS)-APICA	Selective group II mGlu antagonist
0971	EGLU	Highly selective group II mGlu antagonist
1209	LY 341495	Highly potent and selective group II mGlu antagonist
4062	LY 341495 disodium salt	Potent and selective group II mGlu antagonist; disodium salt of LY 341495 (Cat. No. 1209)
2913	Ro 64-5229	Selective, non-competitive mGlu ₂ antagonist
Modulators		2
4048	BINA	Selective positive allosteric modulator of mGlu ₂ receptors
3949	CBiPES	Positive allosteric modulator of mGlu, receptors
3283	LY 487379	Selective positive allosteric modulator of mGlu, receptors
4388	MNI 137	Selective negative allosteric modulator of group II mGlu receptors
5362	TASP 0433864	Selective positive allosteric modulator of mGlu ₂ receptors
Group III mGI	lu Receptors	
Agonists	-	
1111	ACPT-I	Group III mGlu agonist
2385	AMN 082	Selective mGlu, agonist
0103	L-AP4	Selective group III mGlu agonist
4119	Cinnabarinic acid	Selective mGlu ₄ agonist
1302	(S)-3,4-DCPG	Potent and selective mGlu _{sa} agonist
0238	O-Phospho-L-serine	Group III mGlu agonist
1220	(RS)-PPG	Potent and selective mGlu _s agonist
Antagonists		Ü
0972	CPPG	Potent group III mGlu antagonist
0853	MPPG	Group III and group II mGlu antagonist; more selective for group III than group II
0803	MSOP	Selective group III mGlu antagonist
0854	MSPG	Group II and group II mGlu antagonist
1369	UBP1112	Group III mGlu antagonist
5248	XAP 044	Potent and selective mGlu ₇ antagonist
Modulators		
5715	(±)-ADX 71743	Negative allosteric modulator of mGlu, receptors; brain penetrant
4832	AZ 12216052	Positive allosteric modulator of mGlu ₈ receptors
2963	MMPIP	Potent and selective negative allosteric modulator of mGlu ₇ receptors
3248	VU 0155041	Potent positive allosteric modulator of mGlu ₄ receptors
3311	VU 0155041 sodium salt	Potent positive allosteric modulator of mGlu ₄ receptors; sodium salt of VU 0155041 (Cat. No. 3248)
3707	VU 0361737	Selective positive allosteric modulator of mGlu ₄ receptors
5378	VU 0422288	Selective positive allosteric modulator of group III mGlu receptors

Non-selectiv	ve mGlu Receptors	
Agonists		
0218	L-Glutamic acid	Endogenous, non-selective glutamate receptor agonist
0285	Ibotenic acid	Non-selective mGlu agonist, also NMDA agonist
Antagonists		
0112	γDGG	Broad spectrum glutamatergic antagonist
0101	DL-AP4	Broad spectrum glutamatergic antagemet Broad spectrum glutamatergic antagonist
0223	Kynurenic acid	Broad spectrum glutamatergic antagemet Broad spectrum glutamatergic antagonist
0336	(RS)-MCPG	Non-selective mGlu antagonist
3696	(RS)-MCPG disodium salt	Non-selective mGlu antagonist; disodium salt of (RS)-MCPG (Cat. No. 0336)
0337	(S)-MCPG	Non-selective mGlu antagonist; active isomer of (RS)-MCPG (Cat. No. 0336)
	ptor Ligand Sets	ion concern make an agency accine isomore in (i.e., ii.e. a (cast ii.e. cocc)
1826	Group I mGlu Receptor Tocriset™	Selection of 5 group I mGlu receptor ligands (Cat. Nos. 0805, 0188, 1237, 1212 and 0337)
1827	Group II mGlu Receptor Tocriset™	Selection of 5 group II mGlu receptor ligands (Cat. Nos. 1208, 0975, 1209, 0971 and 0337)
1828	Group II mGlu Receptor Tocriset™	Selection of 5 group III mGlu receptor ligands (Cat. Nos. 0103, 1220, 0972, 0803 and 1209)
1829	Mixed mGlu Receptor Tocriset™	Selection of 5 mixed mGlu receptor ligands (Cat. Nos. 0805, 0975, 0103, 0337 and 1209)
Caged Gluta	amate Compounds	
6553	JF-NP-26	Caged Raseglurant (Cat. No. 4416)
5785	MDNI-caged-L-gluatmate	Stable photoreleaser of L-glutamate
1490	MNI-caged-L-glutamate	Stable photoreleaser of L-glutamate
3332	NPEC-caged-LY 379268	Caged version of LY 379268 (Cat. No. 2453)
3574	RuBi-Glutamate	Caged glutamate; excited by visible wavelengths
Miscellaneo	ous Glutamate	
3618	Acamprosate	Glutamatergic modulator and GABA agonist
1611	Lamotrigine	Inhibits glutamate release; anticonvulsant
2289	Lamotrigine isethionate	Inhibits glutamate release; water-soluble salt of Lamotrigine (Cat. No. 1611)
2538	L-BMAA	Glutamate agonist; neurotoxic amino acid
0768	Riluzole	Glutamate release inhibitor; also inhibits GABA uptake and blocks Na, channels
2625	Zonisamide	Anticonvulsant, modulates glutamate neurotransmission
Carboxypep	otidase	
Inhibitors		
5033	2-MPPA	Selective glutamate carboxypeptidase II (GCP II) inhibitor; orally bioavailable
1380	PMPA (NAALADase inhibitor)	Highly potent, selective NAALADase (GCP II) inhibitor
2675	ZJ 43	Glutamate carboxypeptidase II and III (NAALADase, NAAG peptidase) inhibitor
Glutamate 1	Transporters (EAATs)	
Inhibitors	,	
0237	7-Chlorokynurenic acid	Potent competitive inhibitor of L-glutamate uptake
3697	7-Chlorokynurenic acid sodium salt	Potent competitive inhibitor of L-glutamate uptake; sodium salt of 7-Chlorokynurenic acid (Cat. No. 0237)
0111	Dihydrokainic acid	Non-transportable inhibitor of EAAT2 (GLT-1)
1223	DL-TBOA	Selective non-transportable inhibitor of EAATs
2532	TFB-TBOA	High affinity EAAT1 and EAAT2 blocker
0183	L-(-)-threo-3-Hydroxyaspartic acid	Transportable EAAT1-4 inhibitor and non-transportable EAAT5 inhibitor
0811	(±)-threo-3-Hydroxyaspartic acid	EAAT2 and EAAT4 blocker
2652	WAY 213613	Potent, non-substrate EAAT2 inhibitor
Modulators		
6578	GT 949	Potent and selective positive allosteric modulator of EAAT2
Other		
	Coftriovono dipadium polt	Increases EAAT2 expression and activity; neuroprotective
3732	Ceftriaxone disodium salt	increases LAATZ expression and activity, neuroprotective

















