

mGlu Antagonists Kit ab120322

概述

产品名称

mGlu Antagonists试剂盒

特异性

A collection of 6 antagonists for mGlu receptor subtypes. Products are freeze-dried to an exact weight so that addition of 1 ml or 0.5 ml of solvent provides a stock solution of 5 mM or 10 mM, respectively. The kit contains LY 341495 sodium salt ([ab120400](#)), (R,S)-MCPG sodium salt ([ab120252](#)), LY 367385 sodium salt ([ab120418](#)), YM 298198 hydrochloride ([ab120015](#)), MTEP hydrochloride ([ab120035](#)) and MMPIP ([ab120245](#)).

产品概述

A collection of 6 antagonists for mGlu receptor subtypes. Products are freeze-dried to an exact weight so that addition of 1 ml or 0.5 ml of solvent provides a stock solution of 5 mM or 10 mM, respectively. The kit contains LY 341495 sodium salt ([ab120400](#)), (R,S)-MCPG sodium salt ([ab120252](#)), LY 367385 sodium salt ([ab120418](#)), YM 298198 hydrochloride ([ab120015](#)), MTEP hydrochloride ([ab120035](#)) and MMPIP ([ab120245](#)).

Receptor

Description

All mGlu
(mGlu_{1,2,3,4,5,6,7,8})

LY 341495 sodium salt:

Very selective nanomolar potent antagonist for group II mGlu receptors. Also a relatively potent antagonist for group III mGlu receptors at high nanomolar to low micromolar concentrations. K_d values are 1.67, 0.75, 31.6, 72.7, 14 nM at human mGlu₂, mGlu₃, mGlu₆, mGlu_{7a}, and human mGlu_{8a} receptors, respectively. IC_{50} values are 7.8, 8.2 and 22 μ M for mGlu_{1a}, mGlu_{5a} and mGlu₄ receptors, respectively. Readily brain penetrant and active *in vivo*.

Suggested Solvent: Water

Group I & II

(R,S)-MCPG sodium salt:

Group I / group II metabotropic glutamate receptor antagonist. Water soluble form.

Suggested Solvent: Water

mGlu₁

LY 367385 sodium salt:

Potent, selective mGlu_{1a} antagonist. Neuroprotective and anticonvulsant *in vivo*.

Suggested Solvent: Water

mGlu₁

YM298198 hydrochloride:

Potent, selective, water-soluble, non-competitive mGlu₁ antagonist (IC₅₀ = 16 nM) Selective over mGlu_{2,3,4,5,6} and 7 and ionotropic receptors (IC₅₀ > 10 μM). Orally active analgesic *in vivo*.

Suggested Solvent: Water

mGlu₅

MTEP hydrochloride:

Brain penetrant, potent, selective and non-competitive mGlu₅ receptor antagonist (IC₅₀ = 5 nM in Ca²⁺-flux assay; K_i = 16 nM). *In vivo* anxiolytic, devoid of side effects seen with MPEP and benzodiazepines.

Suggested Solvent: Water

mGlu₇

MMPIP:

Potent, selective allosteric antagonist at mGlu₇ receptors. Inhibits agonist-induced intracellular calcium mobilisation and cAMP accumulation (IC₅₀ values are 26 and 610 nM). Also displays intrinsic activity; increasing forskolin-induced cAMP accumulation in the absence of an agonist. Exhibits no significant activity at mGluR1, mGluR2, mGluR3, mGluR4, mGluR5, and mGluR8.

Suggested Solvent: DMSO

说明

Providing storage is as stated on the product vial and the vial is kept tightly sealed, the product can be stored for up to 6 months.

Wherever possible, you should prepare and use solutions on the same day. However, if you need to make up stock solutions in advance, we recommend that you store the solution as aliquots in tightly sealed vials at -20°C. Generally, these will be useable for up to one week. Before use, and prior to opening the vial we recommend that you allow your product to equilibrate to room temperature for at least 1 hour.

KIT

性能

存放说明

Store at -20°C. Please refer to protocols.

组件	1 kit
ab120252 - (R,S)-MCPG sodium salt	1 vial
ab120400 - LY 341495 sodium salt	1 vial
ab120418 - LY 367385 sodium salt	1 vial
ab120245 - MMPIP	1 vial
ab120035 - MTEP hydrochloride	1 vial
ab120015 - YM298198 hydrochloride	1 vial

相关性

Metabotropic Glutamate Receptor contains a Periplasmic Binding Protein Type 1 region, which belongs to Type 1 periplasmic binding fold superfamily. This model and hierarchy represent the ligand binding domains of the LacI family of transcriptional regulators, periplasmic binding proteins of the ABC-type transport systems.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours

- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors