

mGlu Agonists Kit ab120321

製品の概要

製品名	mGlu Agonists Kit
特異性	<p>A collection of 7 agonists and modulators for a range of mGlu receptor subtypes. Products are freeze-dried as an exact weight so that addition of 1 ml or 0.5 ml of water provides a stock solution of 5 mM or 10 mM*, respectively.</p> <p>The kit contains (S)-3,5-DHPG (ab120007), (R,S)-CHPG sodium salt (ab120221), LY 379268 (ab120196), L-AP4 (ab120002), VU0155041 sodium salt (ab120248), AMN082 dihydrochloride (ab120011) and (S)-3,4-DCPG (ab120006). *Please note that AMN082 dihydrochloride (ab120011) is only soluble to 5 mM in water.</p>

製品の概要	<p>A collection of 7 agonists and modulators for a range of mGlu receptor subtypes. Products are freeze-dried as an exact weight so that addition of 1 ml or 0.5 ml of water provides a stock solution of 5 mM or 10 mM*, respectively.</p> <p>The kit contains (S)-3,5-DHPG (ab120007), (R,S)-CHPG sodium salt (ab120221), LY 379268 (ab120196), L-AP4 (ab120002), VU0155041 sodium salt (ab120248), AMN082 dihydrochloride (ab120011) and (S)-3,4-DCPG (ab120006).</p> <p>*Please note that AMN082 dihydrochloride (ab120011) is only soluble to 5 mM in water.</p> <p>Receptor</p> <p>Description.</p> <p>mGlu_{1,5}</p> <p>(S)-3,5-DHPG:</p> <p>Selective group I mGlu agonist.</p> <p>Suggested Solvent: Water</p> <p>mGlu₅</p> <p>(R,S)-CHPG sodium salt:</p> <p>Water soluble form of (R,S)-CHPG (ab120039), a selective mGlu₅ agonist.</p> <p>Suggested Solvent: Water</p> <p>mGlu_{2,3}</p> <p>LY379268:</p> <p>Highly potent and systemically active mGlu₂ and mGlu₃ agonist. EC₅₀ values are 2.69 and 4.48 nM for hmGlu₂ and hmGlu₃, respectively and displays > 80-fold selectivity over group I and group III receptors. Effective in several animal models of stroke, epilepsy, drug abuse, schizophrenia, and pain.</p>
-------	--

Suggested Solvent: Water

mGlu_{4,6,7,8} (group III)

L-AP4:

Selective group III mGlu receptor agonist.

Suggested Solvent: Water

mGlu₄

VU0155041 sodium salt:

Water soluble, positive allosteric selective modulator at mGluR₄ (EC₅₀ values are 798 nM and 693 nM at human and rat mGluR4 receptors, respectively). It has advantages over PHCCC ([ab120043](#)) as it is water soluble and more potent. Active in behavioural models of Parkinsons Disease.

Suggested Solvent: Water

mGlu₇

AMN082 dihydrochloride:

Potent, selective, orally active and brain penetrant mGlu₇ receptor agonist. Potently inhibits cAMP accumulation and stimulates GTP binding (EC₅₀ = 64-290 nM) in mGluR7 transfected cells. Selective over mGlu_{1, 2, 3, 4, 5, 6} and ₈ and ionotropic receptors (up to 10 µM). Elevates plasma stress hormones *in vivo*. Induces mGlu₇ receptor internalisation in dissociated hippocampal neurones.

Suggested Solvent: Water (5 mM) / DMSO (10 mM)

mGlu₈

(S)-3,4-DCPG:

Potent, selective mGlu_{8a} agonist (EC₅₀ = 31 nM). Displays > 100-fold selectivity over mGlu₇. Systemically active *in vivo*.

Suggested Solvent: Water

特記事項

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

製品の特性

保存方法

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

内容	1 kit
ab120221 - (RS)-CHPG sodium salt	1 vial

内容	1 kit
<u>ab120006 - (S)-3,4-DCPG</u>	1 vial
<u>ab120007 - (S)-3,5-DHPG</u>	1 vial
<u>ab120011 - AMN082 dihydrochloride</u>	1 vial
<u>ab120002 - L-AP4</u>	1 vial
<u>ab120196 - LY 379268</u>	1 vial
<u>ab120248 - VU0155041 sodium salt</u>	1 vial

関連性 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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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