

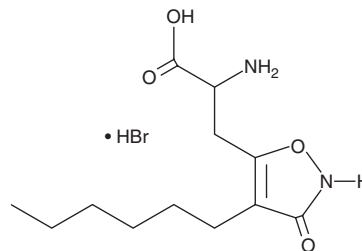
PRODUCT INFORMATION



HexylHIBO (hydrobromide)

Item No. 29164

Formal Name: α -amino-4-hexyl-2,3-dihydro-3-oxo-5-isoxazolepropanoic acid, monohydrobromide
Synonym: Hexylhomoibotenic Acid
MF: $C_{12}H_{20}N_2O_4 \cdot HBr$
FW: 337.2
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

HexylHIBO (hydrobromide) is supplied as a solid. A stock solution may be made by dissolving the HexylHIBO (hydrobromide) in the solvent of choice, which should be purged with an inert gas. HexylHIBO (hydrobromide) is slightly soluble (0.1-1 mg/ml) in methanol.

Description

HexylHIBO is an antagonist of group I metabotropic glutamate receptors (mGluRs; K_b s = 140 and 110 μM for mGluR1 α and mGluR5 α , respectively).¹ It is selective for group I mGluRs over mGluR2 (group II) and mGluR4 α (group III) (K_b s = $>1,000 \mu M$ for both), as well as AMPA, NMDA, and kainate receptors (IC_{50} s = $>100 \mu M$ for all). HexylHIBO (6.3 and 12.5 mg/kg, i.c.v.) inhibits NMDA-induced seizures in mice. It also potentiates increases in plasma adrenocorticotrophic hormone (ACTH) levels in a rat model of restraint stress.²

References

1. Madsen, U., Bräuner-Osborne, H., Frydenvang, K., *et al.* Synthesis and pharmacology of 3-isoxazolol amino acids as selective antagonists at group I metabotropic glutamic acid receptors. *J. Med. Chem.* **44**(7), 1051-1059 (2001).
2. Evanson, N.K. and Herman, J.P. Metabotropic glutamate receptor-mediated signaling dampens the HPA axis response to restraint stress. *Physiol. Behav.* **150**, 2-7 (2015).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/29/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM