

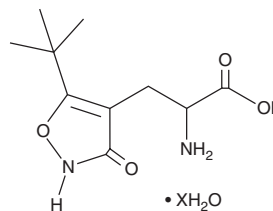
PRODUCT INFORMATION



ATPA (hydrate)

Item No. 30882

Formal Name: α -amino-5-(1,1-dimethylethyl)-2,3-dihydro-3-oxo-4-isoxazolepropanoic acid, hydrate
MF: $C_{10}H_{16}N_2O_4 \cdot XH_2O$
FW: 228.2
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

ATPA (hydrate) is supplied as a solid. A stock solution may be made by dissolving the ATPA (hydrate) in water. The solubility of ATPA (hydrate) in water is approximately 10 mM. We do not recommend storing the aqueous solution for more than one day.

Description

ATPA is an agonist of ionotropic glutamate receptor 5 (GluR5/GluK1; $K_i = 0.0043 \mu M$).¹ It is selective for GluR5 over GluR1-4 (K_i s = 6-14 μM), as well as GluR6 ($K_i = >1,000 \mu M$). ATPA induces inward currents in HEK293 cells expressing recombinant human GluR5 and isolated rat dorsal root ganglion (DRG) neurons (EC_{50} s = 2.1 and 0.6 μM , respectively).

Reference

1. Clarke, V.R., Ballyk, B.A., Hoo, K.H., *et al.* A hippocampal GluR5 kainate receptor regulating inhibitory synaptic transmission. *Nature* **389(6651)**, 599-603 (1997).

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, 11/02/2022

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