

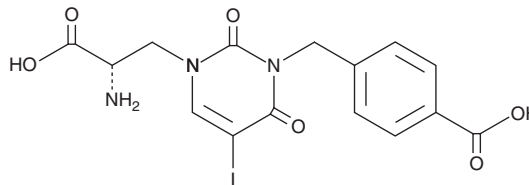
PRODUCT INFORMATION



UBP 301

Item No. 17597

CAS Registry No.: 569371-10-4
Formal Name: (αS)-α-amino-3-[(4-carboxyphenyl)methyl]-3,4-dihydro-5-iodo-2,4-dioxo-1(2H)-pyrimidinepropanoic acid
MF: C₁₅H₁₄IN₃O₆
FW: 459.2
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 280 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

UBP 301 is supplied as a crystalline solid. A stock solution may be made by dissolving the UBP 301 in the solvent of choice, which should be purged with an inert gas. UBP 301 is soluble in the organic solvent DMSO (warmed) at a concentration of approximately 0.5 mg/ml.

Description

UBP 301 is an antagonist of the kainate receptor, an ionotropic glutamate receptor (apparent $K_d = 5.94 \mu\text{M}$).¹ It displays ~30-fold selectivity over AMPA receptors.¹ UBP 301 is used to assess the roles of kainate receptors in neuronal signaling, particularly in reference to neuropathic pain.^{2,3}

References

1. More, J.C.A., Troop, H.M., Dolman, N.P., *et al.* Structural requirements for novel willardiine derivatives acting as AMPA and kainate receptor antagonists. *Br. J. Pharmacol.* **138(6)**, 1093-1100 (2003).
2. Doolen, S., Blake, C.B., Smith, B.N., *et al.* Peripheral nerve injury increases glutamate-evoked calcium mobilization in adult spinal cord neurons. *Mol. Pain* **8:56**, (2012).
3. Osgood, D.B., Harrington, W.F., Kenney, E.V., *et al.* The utility of ionotropic glutamate receptor antagonists in the treatment of nociception induced by epidural glutamate infusion in rats. *Surg. Neurol. Int.* **4**, 106 (2013).

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, 02/20/2023

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