

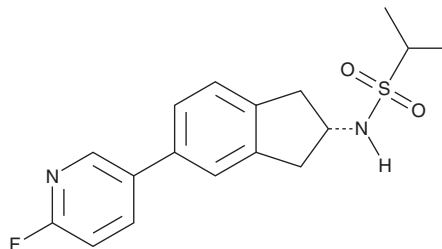
PRODUCT INFORMATION



UoS 12258

Item No. 36571

CAS Registry No.: 875927-64-3
Formal Name: N-[(2S)-5-(6-fluoro-3-pyridinyl)-2,3-dihydro-1H-inden-2-yl]-2-propanesulfonamide
MF: C₁₇H₁₉FN₂O₂S
FW: 334.4
Purity: ≥98%
UV/Vis.: λ_{max}: 250 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

UoS 12258 is supplied as a solid. A stock solution may be made by dissolving the UoS 12258 in the solvent of choice, which should be purged with an inert gas. UoS 12258 is soluble in organic solvents such as ethanol and DMSO. The solubility of UoS 12258 in these solvents is approximately 1 and 2 mg/ml, respectively.

Description

UoS 12258 is a positive allosteric modulator of AMPA receptors.¹ It potentiates glutamate-induced currents in HEK293 cells expressing human ionotropic glutamate receptor 2 (GluR2) homomeric AMPA receptors (EC₅₀ = 2.51 μM), as well as HEK293 cells expressing GluR1, GluR3, or GluR4 homomeric AMPA receptors. UoS 12258 (0.3 and 1 mg/kg) increases novel object exploration in a novel object recognition test in rats. It also reduces escape latencies in the Morris water maze in aged rats.

Reference

1. Ward, S.E., Beswick, P., Calcinaghi, N., *et al.* Pharmacological characterization of N-[(2S)-5-(6-fluoro-3-pyridinyl)-2, 3-dihydro-1H-inden-2-yl]-2-propanesulfonamide: a novel, clinical AMPA receptor positive allosteric modulator. *Br. J. Pharmacol.* **174**(5), 370-385 (2017).

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, 12/20/2022

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