

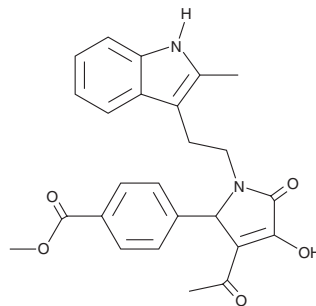
PRODUCT INFORMATION



PYD-106

Item No. 28837

CAS Registry No.: 1560894-05-4
Formal Name: 4-[3-acetyl-2,5-dihydro-4-hydroxy-1-[2-(2-methyl-1H-indol-3-yl)ethyl]-5-oxo-1H-pyrrol-2-yl]-benzoic acid, methyl ester
MF: C₂₅H₂₄N₂O₅
FW: 432.5
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 248 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

PYD-106 is supplied as a solid. A stock solution may be made by dissolving the PYD-106 in the solvent of choice, which should be purged with an inert gas. PYD-106 is soluble in the organic solvent DMSO.

Description

PYD-106 is a positive allosteric modulator of NR2C subunit-containing NMDA receptors (NMDARs).¹ It enhances glutamate-induced maximal current responses in HEK293 cells expressing NR1/NR2C NMDA receptors (EC₅₀ = 13 μM). PYD-106 is selective for NR2C subunit-containing NMDA receptors over NRA-, NR2B-, and NR2D-subunit containing NMDA receptors in *Xenopus* oocytes expressing NR1/NR2 heteromeric NMDA receptors at 30 μM.

Reference

1. Khatri, A., Burger, P.B., Swanger, S.A., *et al.* Structural determinants and mechanism of action of a GluN2C-selective NMDA receptor positive allosteric modulator. *Mol. Pharmacol.* **86**(5), 548-560 (2014).

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, 11/21/2022

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