

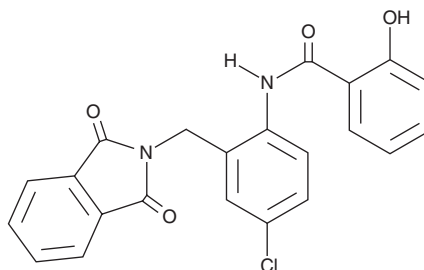
PRODUCT INFORMATION



CPPHA

Item No. 21275

CAS Registry No.: 693288-97-0
Formal Name: N-[4-chloro-2-[(1,3-dihydro-1,3-dioxo-2H-isindol-2-yl)methyl]phenyl]-2-hydroxy-benzamide
MF: C₂₂H₁₅ClN₂O₄
FW: 406.8
Purity: ≥98%
UV/Vis.: λ_{max}: 219 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

CPPHA is supplied as a solid. A stock solution may be made by dissolving the CPPHA in the solvent of choice, which should be purged with an inert gas. CPPHA is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

CPPHA is a positive allosteric modulator of metabotropic glutamate receptor 5 (mGluR5) that potentiates human and rat mGluR5 activation by glutamate, 3,5-DHPG (Item No. 14411), and quisqualate (EC₅₀s = 0.316-0.5 and 0.634-1.16 μM for human and rat mGluR5, respectively).¹ It is selective for mGluR5 over other mGluRs but exhibits submicromolar activity at PDE6, δ- and κ-opioid receptors, and the ether-a-go-go related gene (ERG) potassium channel (K_is = 500, 352, 474, and 813 nM, respectively). CPPHA potentiates 3,5-DHPG-induced NMDA currents in CA1 pyramidal cells and the depolarization of rat subthalamic nucleus (STN) neurons.

Reference

1. O'Brien, J.A., Lemaire, W., Wittmann, M., *et al.* A novel selective allosteric modulator potentiates the activity of native metabotropic glutamate receptor subtype 5 in rat forebrain. *J. Pharm. Exp. Ther.* **309**(2), 568-577 (2004).

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, 07/01/2025

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