

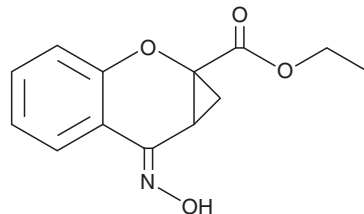
PRODUCT INFORMATION



CPCCOEt

Item No. 21486

CAS Registry No.: 179067-99-3
Formal Name: 7,7a-dihydro-7-(hydroxyimino)-benzo[b]cyclopropa[e]pyran-1a(1H)-carboxylic acid, ethyl ester
MF: C₁₃H₁₃NO₄
FW: 247.3
Purity: ≥98%
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

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

Description

CPCCOEt is a low affinity, selective, non-competitive antagonist of the metabotropic glutamate receptor subtype mGluR1b. It has been shown to inhibit glutamate-induced increases in intracellular calcium at human mGluR1b with an IC₅₀ value of 6.5 μM while having no agonist or antagonist activity at mGluR2, -4a, -5a, -7b, and -8a at concentrations up to 100 μM.¹

Reference

1. Litschig, S., Gasparini, F., Rueegg, D., *et al.* CPCCOEt, a noncompetitive metabotropic glutamate receptor 1 antagonist, inhibits receptor signaling without affecting glutamate binding. *Mol. Pharmacol.* **55(3)**, 453-461 (1999).

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, 11/08/2022

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