

PRODUCT INFORMATION



HPI-1 (hydrate)

Item No. 15221

Formal Name: 1,4,5,6,7,8-hexahydro-4-(3-hydroxyphenyl)-7-(2-methoxyphenyl)-2-methyl-5-oxo-3-quinolinecarboxylic acid-2-methoxyethyl ester, hydrate

MF: C₂₇H₂₉NO₆ • XH₂O

FW: 463.5

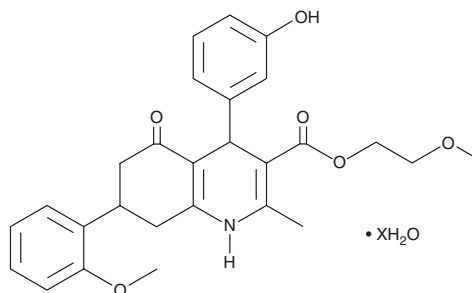
Purity: ≥95%

UV/Vis.: λ_{max}: 218, 365 nm

Supplied as: A crystalline 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

HPI-1 (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the HPI-1 (hydrate) in the solvent of choice. HPI-1 (hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of HPI-1 (hydrate) in these solvents is approximately 20, 3, and 10 mg/ml, respectively.

HPI-1 (hydrate) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

HPI-1 (hydrate) is a Hedgehog (Hh) pathway inhibitor that suppresses signaling through Sonic Hh (IC₅₀ = 1.5 μM) without significantly affecting Wnt signaling (IC₅₀ ≥ 30 μM).¹ HPI-1 suppresses Hh activation induced by loss of Suppressor of Fused or by Gli overexpression, suggesting action at posttranslational modification of Gli protein or at the interaction of Gli with a co-factor.¹ HPI-1 (hydrate) also inhibits signaling through the oncogenic Smoothed (Smo) mutant SmoM2 in neuron precursors, preventing cell proliferation.¹

Reference

1. Hyman, J.M., Firestone, A.J., Heine, V.M., *et al.* Small-molecule inhibitors reveal multiple strategies for Hedgehog pathway blockade. *Proc. Natl. Acad. Sci. USA* **106**(33), 14132-14137 (2009).

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