

PRODUCT INFORMATION

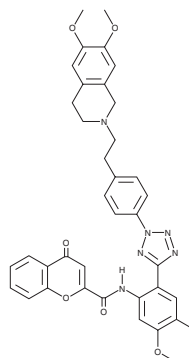


HM30181

Item No. 32873

CAS Registry No.: 849675-66-7
Formal Name: N-[2-[2-[4-[2-(3,4-dihydro-6,7-dimethoxy-2(1H)-isoquinolinyl)ethyl]phenyl]-2H-tetrazol-5-yl]-4,5-dimethoxyphenyl]-4-oxo-4H-1-benzopyran-2-carboxamide

Synonym: Encequidar
MF: C₃₈H₃₆N₆O₇
FW: 688.7
Purity: ≥98%
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

HM30181 is supplied as a crystalline solid. A stock solution may be made by dissolving the HM30181 in the solvent of choice, which should be purged with an inert gas. HM30181 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of HM30181 in these solvents is approximately 30 mg/ml.

Description

HM30181 is an inhibitor of P-glycoprotein (P-gp), also known as multidrug resistance protein 1 (MDR1; IC₅₀ = 0.63 nM).¹ It is selective for P-gp over breast cancer resistance protein (BCRP; IC₅₀ = 3,700 nM), as well as multidrug resistance-associated protein 1 (MRP1), MRP2, and MRP3 at 100 μM. It inhibits transepithelial basal-to-apical transport of paclitaxel in MDCK monolayers overexpressing P-gp (IC₅₀ = 35.4 nM). HM30181 inhibits tumor growth and induces tumor regression in an HT-29 mouse xenograft model when administered at doses of 10 and 20 mg/kg, respectively, in combination with paclitaxel (Item No. 10461).

Reference

1. Kwak, J.-O., Lee, S.H., Lee, G.S., *et al.* Selective inhibition of MDR1 (ABCB1) by HM30181 increases oral bioavailability and therapeutic efficacy of paclitaxel. *Eur. J. Pharmacol.* **627**(1-3), 92-98 (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, 10/27/2022

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