

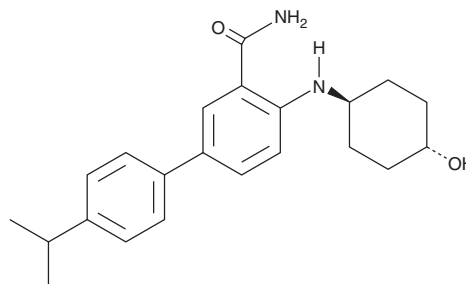
PRODUCT INFORMATION



GRP94 Inhibitor 1

Item No. 43140

CAS Registry No.: 2234897-35-7
Formal Name: 4-[(*trans*-4-hydroxycyclohexyl)amino]-4'-(1-methylethyl)-[1,1'-biphenyl]-3-carboxamide
MF: C₂₂H₂₈N₂O₂
FW: 352.5
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

GRP94 inhibitor 1 is supplied as a solid. A stock solution may be made by dissolving the GRP94 inhibitor 1 in the solvent of choice, which should be purged with an inert gas. GRP94 inhibitor 1 is soluble (≥10 mg/ml) in DMSO. GRP94 inhibitor 1 is sparingly soluble (1-10 mg/ml) in ethanol.

Description

GRP94 inhibitor 1 is a ligand of glucose-regulated protein 94 (GRP94; IC₅₀ = 2 nM in a fluorescence polarization assay).¹ It decreases the levels of integrin α2 and integrin αL in PANC-1 cells when used at concentrations of 2.5 or 5 μM. GRP94 inhibitor 1 (10 or 30 mg/kg twice per day) increases body weight, decreases disease severity, and reduces serum and colon cell levels of TNF-α and IL-6 in a mouse model of ulcerative colitis induced by dextran sodium sulfate (DSS).

Reference

1. Jiang, F., Guo, A.-P., Xu, J.-C., *et al.* Discovery of a potent Grp94 selective inhibitor with anti-inflammatory efficacy in a mouse model of ulcerative colitis. *J. Med. Chem.* **61**(21), 9513-9533 (2018).

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, 04/16/2025

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