

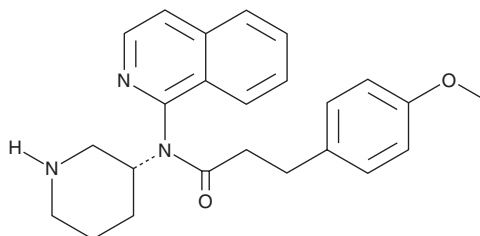
PRODUCT INFORMATION



(R)-IMPP

Item No. 20861

CAS Registry No.: 2133832-83-2
Formal Name: N-1-isoquinolinyl-4-methoxy-N-(3R)-3-piperidinyl-benzenepropanamide
MF: C₂₄H₂₇N₃O₂
FW: 389.5
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 275, 323 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

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

Description

(R)-IMPP is a cell-permeable inhibitor of the secretion of proprotein convertase subtilisin/kexin type 9 (PCSK9; IC₅₀ = 4.8 μM).¹ It promotes the uptake of LDL cholesterol in hepatoma cells by increasing LDL receptor levels.¹ (R)-IMPP apparently impairs PCSK9 secretion by causing transcript-dependent inhibition of PCSK9 translation.¹

Reference

1. Petersen, D.N., Hawkins, J., Ruangsiriluk, W., *et al.* A small-molecule anti-secretagogue of PCSK9 targets the 80S ribosome to inhibit PCSK9 protein translation. *Cell. Chem. Biol.* **23(11)**, 1362-1371 (2016).

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, 02/06/2025

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