

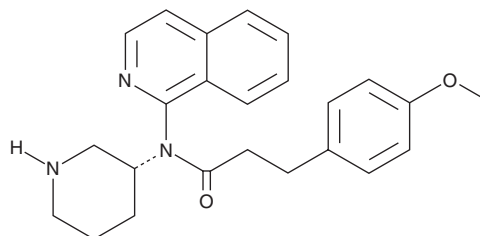
# PRODUCT INFORMATION



## (R)-IMPP

Item No. 20861

**CAS Registry No.:** 2133832-83-2  
**Formal Name:** N-1-isoquinolinyl-4-methoxy-N-(3R)-3-piperidinyl-benzenepropanamide  
**MF:** C<sub>24</sub>H<sub>27</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 389.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 220, 275, 323 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 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. (R)-IMPP is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. 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; IC<sub>50</sub> = 4.8 μM).<sup>1</sup> It promotes the uptake of LDL cholesterol in hepatoma cells by increasing LDL receptor levels.<sup>1</sup> (R)-IMPP apparently impairs PCSK9 secretion by causing transcript-dependent inhibition of PCSK9 translation.<sup>1</sup>

### 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**, 1-10 (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, 09/06/2022

#### 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