

PRODUCT INFORMATION

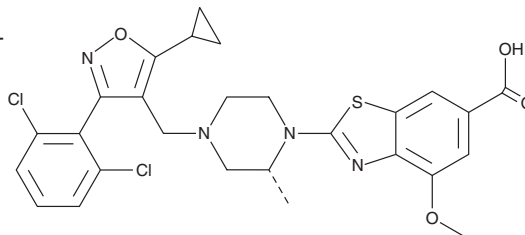


HPG1860

Item No. 39481

CAS Registry No.: 2226133-29-3
Formal Name: 2-[(2R)-4-[[5-cyclopropyl-3-(2,6-dichlorophenyl)-4-isoxazolyl]methyl]-2-methyl-1-piperazinyl]-4-methoxy-6-benzothiazolecarboxylic acid

MF: C₂₇H₂₆Cl₂N₄O₄S
FW: 573.5
Purity: ≥98%
Supplied as: A neat 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

HPG1860 is supplied as a neat solid. A stock solution may be made by dissolving the HPG1860 in the solvent of choice, which should be purged with an inert gas. HPG1860 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

HPG1860 is a farnesoid X receptor (FXR) agonist.¹ It induces luminescence in a reporter assay using HEK293T cells expressing human FXR (EC₅₀ = 18 nM). *In vivo*, HPG1860 (1, 3, or 10 mg/kg per day) decreases the serum levels of alanine transaminase (ALT) and total cholesterol, as well as hepatic inflammation, fat levels, and fibrosis, in a mouse model of non-alcoholic steatohepatitis (NASH) induced by a high-fat diet and carbon tetrachloride (CCl₄).

Reference

1. Mo, C., Xu, X., Zhang, P., *et al.* Discovery of HPG1860, a structurally novel nonbile acid FXR agonist currently in clinical development for the treatment of nonalcoholic steatohepatitis. *J. Med. Chem.* **66**(14), 9363-9375 (2023).

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, 08/12/2024

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