

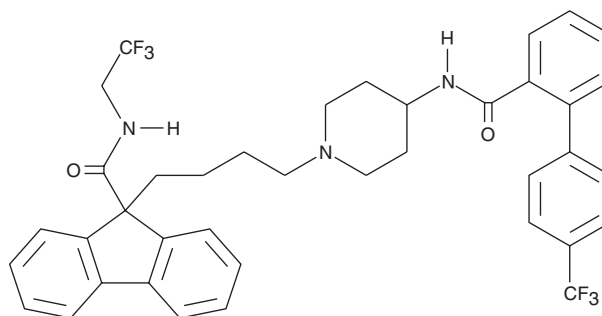
PRODUCT INFORMATION



Lomitapide

Item No. 10009610

CAS Registry No.: 182431-12-5
Formal Name: N-(2,2,2-trifluoroethyl)-9-[4-[4-[[[4'-(trifluoromethyl)[1,1'-biphenyl]-2-yl]carbonyl]amino]-1-piperidiny]butyl]-9H-fluorene-9-carboxamide
Synonyms: AEGR 773, BMS 201038
MF: C₃₉H₃₇F₆N₃O₂
FW: 693.7
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 259 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

Lomitapide is supplied as a crystalline solid. A stock solution may be made by dissolving the lomitapide in the solvent of choice, which should be purged with an inert gas. Lomitapide is soluble in organic solvents such as ethanol and DMSO. The solubility of lomitapide in these solvents is approximately 10 and 5 mg/ml, respectively.

Description

Lomitapide is an inhibitor of microsomal triglyceride transfer protein (MTTP; IC₅₀ = 0.5 nM in a triglyceride transfer assay).¹ It inhibits apolipoprotein B (ApoB) secretion in HepG2 cells (EC₅₀ = 0.8 nM). Lomitapide inhibits triglyceride secretion in fasted rats (ED₅₀ = 0.19 mg/kg, p.o.) and reduces total plasma cholesterol levels in hamsters (ED₅₀ = 2.4 mg/kg). It also decreases plasma cholesterol and triglyceride levels in the Watanabe-heritable hyperlipidemic (WHHL) rabbit model of homozygous familial hypercholesterolemia when administered at a dose of 10 mg/kg. Formulations containing lomitapide have been used as an adjunct to a low-fat diet and other lipid-lowering treatments in the treatment of homozygous familial hypercholesterolemia.

Reference

1. Wetterau, J.R., Gregg, R.E., Harrity, T.W., et al. An MTP inhibitor that normalizes atherogenic lipoprotein levels in WHHL rabbits. *Science* **282**(5389), 751-754 (1998).

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, 12/14/2022

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