

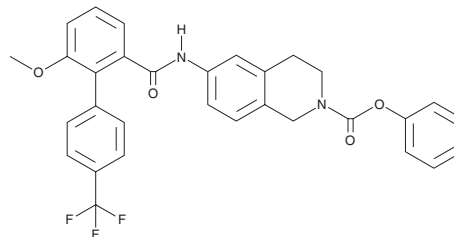
PRODUCT INFORMATION



SLX-4090

Item No. 41979

CAS Registry No.: 913541-47-6
Formal Name: 3,4-dihydro-6-[[[6-methoxy-4'-(trifluoromethyl)[1,1'-biphenyl]-2-yl]carbonyl]amino]-2(1H)-isoquinolinecarboxylic acid, phenyl ester
MF: C₃₁H₂₅F₃N₂O₄
FW: 546.5
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

SLX-4090 is supplied as a solid. A stock solution may be made by dissolving the SLX-4090 in the solvent of choice, which should be purged with an inert gas. SLX-4090 is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

SLX-4090 is an inhibitor of microsomal triglyceride transfer protein (MTTP; IC₅₀ = 8 nM).¹ It selectively inhibits MTTP-mediated secretion of apolipoprotein B (ApoB) over ApoA1 in primary human hepatocytes (IC₅₀s = 81 and >30,000 nM, respectively). *In vivo*, SLX-4090 reduces postprandial increases in serum triglyceride levels in rats (ED₅₀ = 7 mg/kg). It reduces serum triglyceride levels, weight gain, and total cholesterol levels in ApoE^{-/-} mice fed a high-fat diet when administered at a dose of 5.8 mg/kg per day.

Reference

1. Kim, E., Campbell, S., Schueller, O., *et al.* A small-molecule inhibitor of enterocytic microsomal triglyceride transfer protein, SLx-4090: Biochemical, pharmacodynamic, pharmacokinetic, and safety profile. *J. Pharmacol. Exp. Ther.* **337**(3), 775-785 (2011).

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, 01/13/2025

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