

PRODUCT INFORMATION



CP 346,086

Item No. 39156

CAS Registry No.: 186390-48-7
Formal Name: N-[1,2,3,4-tetrahydro-2-(1H-1,2,4-triazol-5-ylmethyl)-6-isoquinolinyl]-4'-(trifluoromethyl)-[1,1'-biphenyl]-2-carboxamide

MF: C₂₆H₂₂F₃N₅O

FW: 477.5

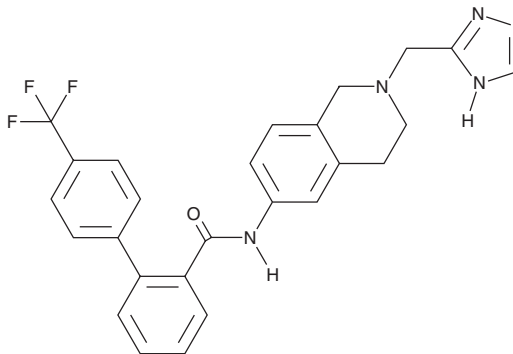
Purity: ≥98%

UV/Vis.: λ_{max}: 246 nm

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

CP 346,086 is supplied as a solid. A stock solution may be made by dissolving the CP 346,086 in the solvent of choice, which should be purged with an inert gas. CP 346,086 is soluble in DMSO.

Description

CP 346,086 is an inhibitor of microsomal triglyceride transfer protein (MTTP; IC₅₀ = 2 nM).¹ It inhibits apolipoprotein B (ApoB) and triglyceride secretion in HepG2 cells (IC₅₀ = 2.6 nM for both). CP 346,086 (10 mg/kg) lowers plasma triglyceride levels, as well as plasma total, VLDL, and LDL cholesterol levels in rats.

Reference

1. Chandler, C.E., Wilder, D.E., Pettini, J.L., *et al.* CP-346086: An MTP inhibitor that lowers plasma cholesterol and triglycerides in experimental animals and in humans. *J. Lipid Res.* **44**(10), 1887-1901 (2003).

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, 09/25/2023

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