

PRODUCT INFORMATION

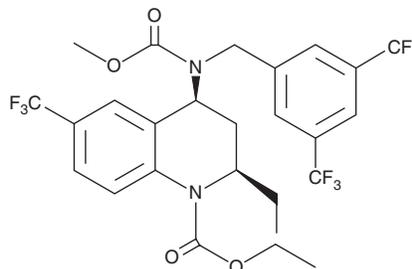


Torcetrapib

Item No. 15028

CAS Registry No.: 262352-17-0
Formal Name: 4S-[[[3,5-bis(trifluoromethyl)phenyl]methyl](methoxycarbonyl)amino]-2R-ethyl-3,4-dihydro-6-(trifluoromethyl)-1(2H)-quinolinecarboxylic acid, ethyl ester
Synonyms: CP 529,414, (-)-Torcetrapib

MF: C₂₆H₂₅F₉N₂O₄
FW: 600.5
Purity: ≥98%
UV/Vis.: λ_{max}: 251 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

Torcetrapib is supplied as a crystalline solid. A stock solution may be made by dissolving the torcetrapib in the solvent of choice, which should be purged with an inert gas. Torcetrapib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of torcetrapib in these solvents is approximately 2, 1, and 0.5 mg/ml, respectively.

Description

Torcetrapib is an inhibitor of cholesteryl ester transfer protein (CETP; IC₅₀s = 47 and 51 nM for cholesteryl and triglyceride transfer, respectively).¹ It is selective for CETP over phospholipid transfer protein (PLTP), lecithin:cholesterol acyltransferase (LCAT), and hepatic and lipoprotein lipases at 10 μM. Torcetrapib (30 mg/kg) inhibits CETP activity and increases HDL cholesterol levels in serum from transgenic mice expressing human CETP and ApoA1.² It increases mean arterial blood pressure in Zucker *fa/fa* diabetic rats when administered at doses of 60 and 200 mg/kg.

References

1. Clark, R.W., Ruggeri, R.B., Cunningham, D., *et al.* Description of the torcetrapib series of cholesteryl ester transfer protein inhibitors, including mechanism of action. *J. Lipid Res.* **47(3)**, 537-552 (2006).
2. Cao, G., Beyer, T.P., Zhang, Y., *et al.* Evacetrapib is a novel, potent, and selective inhibitor of cholesteryl ester transfer protein that elevates HDL cholesterol without inducing aldosterone or increasing blood pressure. *J. Lipid Res.* **52(12)**, 2169-2176 (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, 10/21/2022

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