

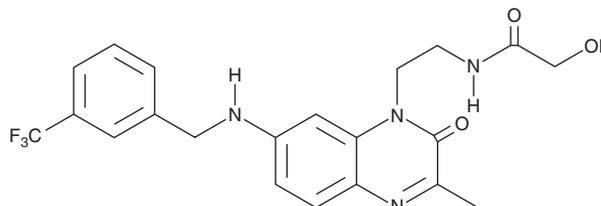
PRODUCT INFORMATION



CVT-12012

Item No. 27156

CAS Registry No.: 1018675-35-8
Formal Name: 2-hydroxy-N-[2-[3-methyl-2-oxo-7-[[[3-(trifluoromethyl)phenyl]methyl]amino]-1(2H)-quinoxaliny]ethyl]-acetamide
MF: C₂₁H₂₁F₃N₄O₃
FW: 434.4
Purity: ≥98%
UV/Vis.: λ_{max}: 228, 379 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

CVT-12012 is supplied as a crystalline solid. A stock solution may be made by dissolving the CVT-12012 in the solvent of choice, which should be purged with an inert gas. CVT-12012 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CVT-12012 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

CVT-12012 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CVT-12012 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CVT-12012 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CVT-12012 is an orally bioavailable inhibitor of stearoyl-CoA desaturase (SCD; IC₅₀ = 6.1 nM in HepG2 cells).¹ It dose-dependently reduces levels of the SCD desaturation products palmitoleic acid (Item No. 10009871) and vaccenic acid in the plasma and liver of rats fed a high-sucrose diet. CVT-12012 (20 mg/kg) also reduces hepatic triglyceride levels in sucrose-fed rats.

Reference

1. Koltun, D.O., Zilbershtein, T.M., Migulin, V.A., *et al.* Potent, orally bioavailable, liver-selective stearoyl-CoA desaturase (SCD) inhibitors. *Bioorg. Med. Chem. Lett.* **19(15)**, 4070-4074 (2009).

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, 10/19/2022

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