

PRODUCT INFORMATION



CBM 301940

Item No. 43420

CAS Registry No.: 902146-11-6

Formal Name: 3-[[[4,5-dihydro-5-[2,2,2-trifluoro-1-hydroxy-1-(trifluoromethyl)ethyl]-3-isoxazolyl]carbonyl]amino]-butanoic acid, 1,1-dimethylethyl ester

MF: C₁₅H₂₀F₆N₂O₅

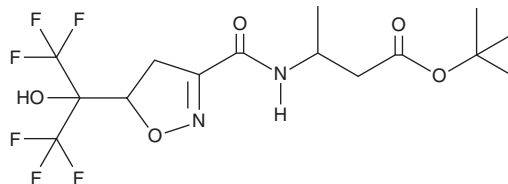
FW: 422.3

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

CBM 301940 is supplied as a solid. A stock solution may be made by dissolving the CBM 301940 in the solvent of choice, which should be purged with an inert gas. CBM 301940 is soluble (≥10 mg/ml) in acetonitrile.

Description

CBM 301940 is an inhibitor of malonyl-CoA decarboxylase (MCD; IC₅₀ = 23 nM).¹ It increases the rate of glucose oxidation in isolated and perfused rat hearts when used at a concentration of 1 μM.

Reference

1. Cheng, J.F., Huang, Y., Penuliar, R., *et al.* Discovery of potent and orally available malonyl-CoA decarboxylase inhibitors as cardioprotective agents. *J. Med. Chem.* **49**(14), 4055-4058 (2006).

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, 07/21/2025

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