

PRODUCT INFORMATION

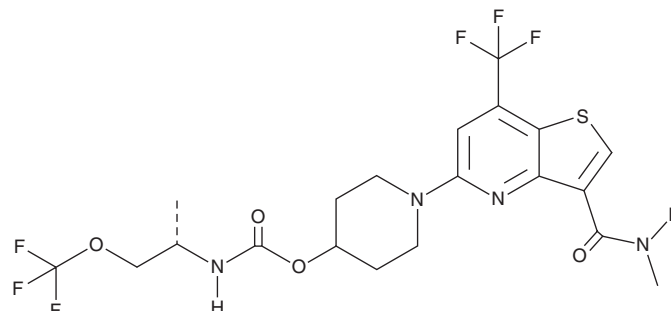


UGT8 Inhibitor 19

Item No. 32723

CAS Registry No.: 2414349-93-0
Formal Name: N-[(1S)-1-methyl-2-(trifluoromethoxy)ethyl]-carbamic acid, 1-[3-[(methylamino)carbonyl]-7-(trifluoromethyl)thieno[3,2-b]pyridin-5-yl]-4-piperidinyl ester
Synonyms: Ceramide Galactosyltransferase Inhibitor 19, CGT Inhibitor 19, UDP-galactose Ceramide Galactosyltransferase Inhibitor 19

MF: C₂₀H₂₂F₆N₄O₄S
FW: 528.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 252 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

UGT8 inhibitor 19 is supplied as a crystalline solid. A stock solution may be made by dissolving the UGT8 inhibitor 19 in the solvent of choice, which should be purged with an inert gas. UGT8 inhibitor 19 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of UGT8 inhibitor 19 in ethanol is approximately 25 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

UGT8 inhibitor 19 is an inhibitor of UDP-galactose ceramide galactosyltransferase (UGT8), also known as ceramide galactosyltransferase (CGT; IC₅₀ = 0.2 nM).¹ It is selective for UGT8 over UGT1A1, -1A6, -2B7, -2B15, and -2B17 (IC₅₀s = >10 μM for all). *In vivo*, UGT8 inhibitor 19 inhibits the synthesis of galactosylceramide (GalCer) and sulfatide (SFT) in juvenile mouse brain (ED₅₀s = <3 mg/kg for both) and adult mouse kidney when administered at doses of 3, 10, and 30 mg/kg.

Reference

1. Thurairatnam, S., Lim, S., Barker, R.H., Jr., *et al.* Brain penetrable inhibitors of ceramide galactosyltransferase for the treatment of lysosomal storage disorders. *ACS Med. Chem. Lett.* **11(10)**, 2010-2016 (2020).

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, 11/09/2022

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