

PRODUCT INFORMATION

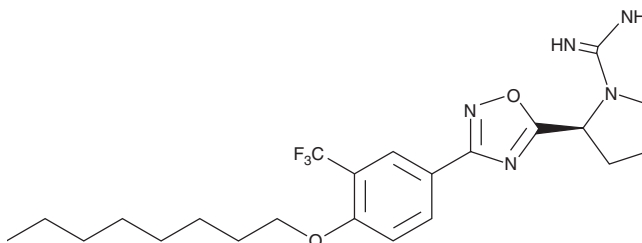


SLM6031434

Item No. 23289

CAS Registry No.: 1897379-33-7
Formal Name: (2S)-2-[3-[4-(octyloxy)-3-(trifluoromethyl)phenyl]-1,2,4-oxadiazol-5-yl]-1-pyrrolidinecarboximidamide

MF: C₂₂H₃₀F₃N₅O₂
FW: 453.5
Purity: ≥98%
UV/Vis.: λ_{max}: 258 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

SLM6031434 is supplied as a crystalline solid. A stock solution may be made by dissolving the SLM6031434 in the solvent of choice, which should be purged with an inert gas. SLM6031434 is slightly soluble in ethanol, methanol, DMSO, and dimethyl formamide.

Description

SLM6031434 is an inhibitor of sphingosine kinase 2 (SPHK2) ($K_i = 0.4 \mu\text{M}$ for the recombinant mouse enzyme).¹ It is selective for SPHK2 over SPHK1 ($K_{iS} > 20 \mu\text{M}$). It decreases sphingosine-1-phosphate (S1P; Item No. 62570) and increases sphingosine levels in U937 monocytic leukemia cells in a concentration-dependent manner. SLM6031434 (5 mg/kg) reduces blood S1P levels in *Sphk1*^{-/-}, but not *Sphk2*^{-/-}, mice. SLM6031434 also increases blood S1P levels in wild-type mice and rats.

Reference

1. Kharel, Y., Morris, E.A., Congdon, M.D., *et al.* Sphingosine kinase 2 inhibition and blood sphingosine 1-phosphate levels. *J. Pharmacol. Exp. Ther.* **355**(1), 23-31 (2015).

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, 03/12/2020

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