

PRODUCT INFORMATION



SLM6031434 (hydrochloride)

Item No. 30434

CAS Registry No.: 1897379-34-8
Formal Name: (2S)-2-[3-[4-(octyloxy)-3-(trifluoromethyl)phenyl]-1,2,4-oxadiazol-5-yl]-1-pyrrolidinecarboximidamide, monohydrochloride

MF: C₂₂H₃₀F₃N₅O₂ • HCl
FW: 490

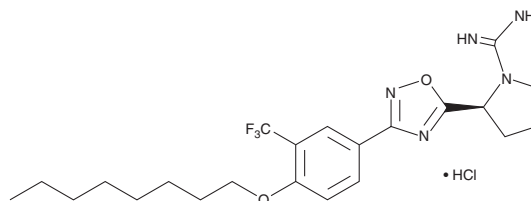
Purity: ≥98%

UV/Vis.: λ_{max}: 259 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 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SLM6031434 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SLM6031434 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SLM6031434 (hydrochloride) in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

SLM6031434 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SLM6031434 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SLM6031434 (hydrochloride) has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

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 ($K_i = >20 \mu\text{M}$). 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, 03/19/2020

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