

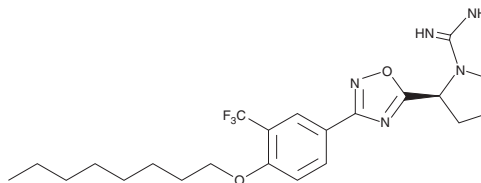
# PRODUCT INFORMATION



**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:** C<sub>22</sub>H<sub>30</sub>F<sub>3</sub>N<sub>5</sub>O<sub>2</sub>  
**FW:** 453.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 258 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.

## Description

SLM6031434 is an inhibitor of sphingosine kinase 2 (SPHK2) ( $K_i = 0.4 \mu\text{M}$  for the recombinant mouse enzyme).<sup>1</sup> 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*<sup>-/-</sup>, but not *Sphk2*<sup>-/-</sup>, 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, 02/07/2025

## 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](http://WWW.CAYMANCHEM.COM)