# PRODUCT INFORMATION



# SLP7111228 (hydrochloride)

Item No. 23290

CAS Registry No.: 1449768-48-2

(2S)-2-[[3-(4-octylphenyl)-1,2,4-oxadiazol-Formal Name:

5-yl]methyl]-1-pyrrolidinecarboximidamide,

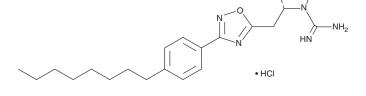
monohydrochloride

MF: C<sub>22</sub>H<sub>33</sub>N<sub>5</sub>O • HCl

420.0 FW: ≥95% **Purity:** UV/Vis.:  $\lambda_{\text{max}}$ : 249 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**

SLP7111228 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SLP7111228 (hydrochloride) in the solvent of choice. SLP7111228 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of SLP7111228 (hydrochloride) in ethanol and DMSO is approximately 30 mg/ml and approximately 12.5 mg/ml in DMF.

SLP7111228 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SLP7111228 (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SLP7111228 (hydrochloride) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

SLP7111228 is a potent and selective inhibitor of sphingosine kinase (SPHK) 1 with K<sub>1</sub> values of 0.048 and >10 μM for SPHK1 and 2, respectively. It induces a concentration-dependent decrease in sphingosine-1-phosphate (S1P; Item No. 62570) in U937 cells but has no effect on sphingosine levels. In vivo, SLP7111228 decreases blood levels of S1P in a dose-dependent manner in rats.

#### Reference

1. Patwardhan, N.N., Morris, E.A., Kharel, Y., et al. Structure-activity relationship studies and in vivo activity of guanidine-based sphingosine kinase inhibitors: Discovery of SphK1- and SphK2-selective inhibitors. J. Med. Chem. 58(4), 1879-1899 (2015).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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

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, 01/30/2018

## **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