

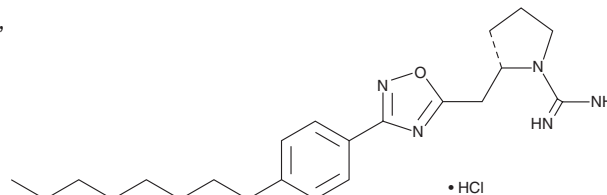
PRODUCT INFORMATION



SLP7111228 (hydrochloride)

Item No. 23290

CAS Registry No.: 1449768-48-2
Formal Name: (2S)-2-[[3-(4-octylphenyl)-1,2,4-oxadiazol-5-yl]methyl]-1-pyrrolidinecarboximidamide, monohydrochloride
MF: C₂₂H₃₃N₅O • HCl
FW: 420.0
Purity: ≥95%
UV/Vis.: λ_{max}: 249 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.

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, which should be purged with an inert gas. SLP7111228 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). 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_i values of 0.048 and >10 μM for SPHK1 and 2, respectively.¹ 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.

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