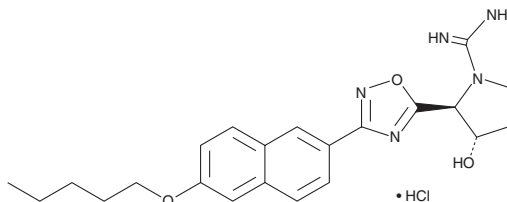


PRODUCT INFORMATION

SLC5111312 (hydrochloride)

Item No. 23288

CAS Registry No.: 1870811-01-0
Formal Name: (2S,3S)-3-hydroxy-2-[3-[6-(pentyloxy)-2-naphthalenyl]-1,2,4-oxadiazol-5-yl]-1-pyrrolidinecarboximidamide, monohydrochloride
MF: $C_{22}H_{27}N_5O_3 \cdot HCl$
FW: 445.9
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 237 nm
Supplied as: A crystalline solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SLC5111312 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the SLC5111312 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SLC5111312 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SLC5111312 (hydrochloride) in these solvents is approximately 14, 3, and 2 mg/ml, respectively.

SLC5111312 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SLC5111312 (hydrochloride) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SLC5111312 (hydrochloride) has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SLC5111312 is a dual inhibitor of sphingosine kinase 1 (SPHK1) and SPHK2 (K_i s = 0.73 and 0.9 μM , respectively, using human recombinant kinases).¹ SLC5111312 binding is selective for Sphk2 over Sphk1 in mice (K_i s = 1 and 20 μM , respectively) but not rat (K_i s = 1.1 and 0.8 μM , respectively).² It decreases cell-associated sphingosine-1-phosphate (S1P) and increases sphingosine in U937 monocytic leukemia cells at concentrations of 0.1 and 0.3 μM . SLC5111312 (10 mg/kg) also decreases S1P *in vivo* in *Sphk1*^{-/-} but not in *Sphk2*^{-/-} mice and in rat, showing a functional selectivity for Sphk2 in mice.

References

1. Congdon, M.D., Kharel, Y., Brown, A.M., *et al.* Structure-activity relationship studies and molecular modeling of naphthalene-based sphingosine kinase 2 inhibitors. *ACS Med. Chem. Lett.* **7**(3), 229-234 (2016).
2. 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