

# PRODUCT INFORMATION



## SLF1081851 (hydrochloride)

Item No. 37150

**Formal Name:** 3-(4-decylphenyl)-1,2,4-oxadiazole-5-propanamine, hydrochloride

**MF:** C<sub>21</sub>H<sub>33</sub>N<sub>3</sub>O • HCl

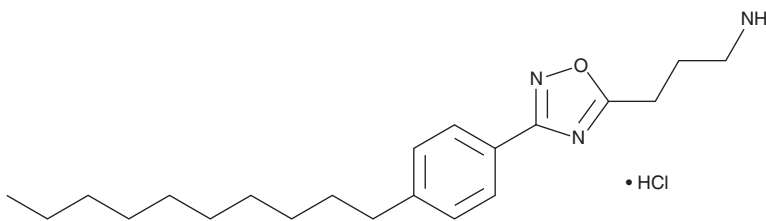
**FW:** 380.0

**Purity:** ≥95%

**Supplied as:** A 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

SLF1081851 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the SLF1081851 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SLF1081851 (hydrochloride) is slightly soluble in acetonitrile.

SLF1081851 (hydrochloride) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

SLF1081851 is an inhibitor of sphingosine-1-phosphate (S1P) transporter (SPNS2).<sup>1</sup> It inhibits mouse Spns2 by 67% when used at a concentration of 2 μM. SLF1081851 inhibits S1P export from HeLa cells expressing mouse Spns2 (IC<sub>50</sub> = 1.93 μM). *In vivo*, SLF1081851 (20 mg/kg) decreases blood lymphocyte numbers and plasma S1P levels in mice.

### Reference

1. Fritzscheier, R., Foster, D., Peralta, A., *et al.* Discovery of *in vivo* active sphingosine-1-phosphate transporter (Spns2) inhibitors. *J. Med. Chem.* **65(11)**, 7656-7681 (2022).

#### 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, 01/16/2024

#### 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