

SLF1081851 . hydrochloride

Sphingosine-1-phosphate transporter inhibitor

SLF1081851 is a sphingosine-1-phosphate transporter (Spns2) inhibitor. It inhibited sphingosine-1-phosphate (S1P) release in HeLa cells. SLF1081851 is active *in vivo* and decreased lymphocyte counts and plasma S1P levels in mice and rats. Spns2 inhibitors represent potential alternative immunomodulatory agents to S1P receptor agonists as they are able to induce lymphopenia without drastic changes to systemic S1P levels or stimulating the S1P1-5 signaling pathway.

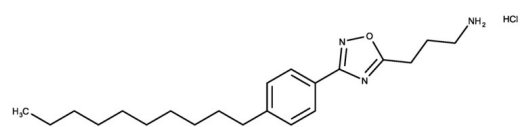
Ordering Information

[Order Online »](#)

ENZ-CHM475-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3-(3-(4-Decylphenyl)-1,2,4-oxadiazol-5-yl)propan-1-amine hydrochloride

Appearance White solid.

CAS 2999629-35-3

Couple Target SPNS2

Couple Type Inhibitor

Formula $C_{21}H_{33}N_3O \cdot HCl$

Identity Determined by NMR.

MW 379.79

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (30 mg/ml) or ethanol (7 mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)