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 in induce lymphopenia without drastic changes to systemic S1P levels or stimulating the S1P1-5 signaling pathway.

Ordering Information

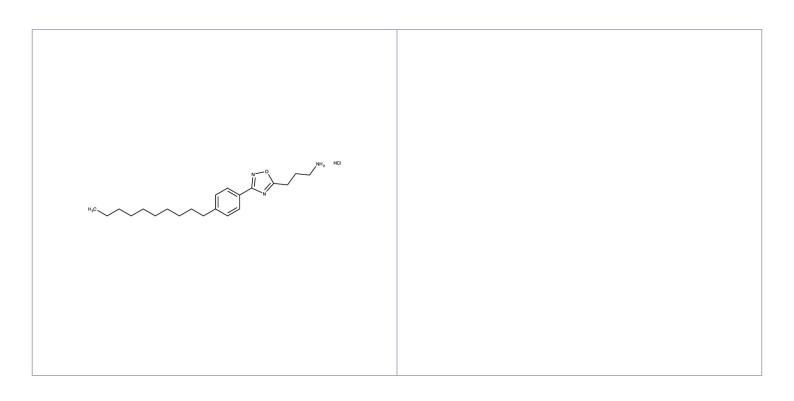
Order Online »

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 $C_{21}H_{33}N_3O$. HCl

Identity Determined by NMR.

MW 379.79

Purity ≥98% (HPLC)

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

Last modified: February 27, 2025



European Sales Office