

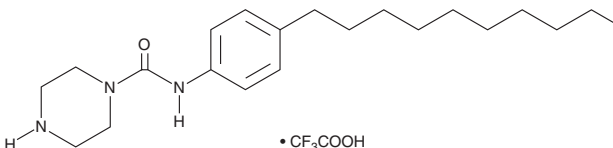
PRODUCT INFORMATION



SLF80821178 (trifluoroacetate salt)

Item No. 44781

Formal Name: N-(4-decylphenyl)-1-piperazinecarboxamide
MF: $C_{21}H_{35}N_3O \cdot CF_3COOH$
FW: 459.6
Purity: $\geq 98\%$
Supplied as: A 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

SLF80821178 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the SLF80821178 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SLF80821178 (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

SLF80821178 is an inhibitor of the sphingosine-1-phosphate transporter (SPNS2).¹ It reduces sphingosine-1-phosphate (S1P) secretion in HeLa cells with an IC_{50} value of 51 nM. Oral administration of SLF80821178 (100 mg/kg) decreases the number of circulating lymphocytes in mice, without reducing plasma S1P concentration. It reduces disease severity in a mouse model of experimental autoimmune encephalomyelitis (EAE) induced by MOG₃₅₋₅₅ (Item No. 35644) when administered at doses of 10 and 30 mg/kg per day.² Unlike the loss of Spns2 function in Spns2^{-/-} mice, inhibition of Spns2 by SLF80821178 does not induce neurosensory hearing defects.

References

1. Foster, D.J., Dunnivant, K., Shrader, C.W., *et al.* Discovery of potent, orally bioavailable sphingosine-1-phosphate transporter (Spns2) inhibitors. *J. Med. Chem.* **67**(13), 11273-11295 (2024).
2. Kharel, Y., Huang, T., Dunnivant, K., *et al.* Assessment of Spinster homologue 2 (Spns2)-dependent transport of sphingosine-1-phosphate as a therapeutic target. *Br. J. Pharmacol.* **182**(9), 2014-2030 (2025).

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, 02/09/2026

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