

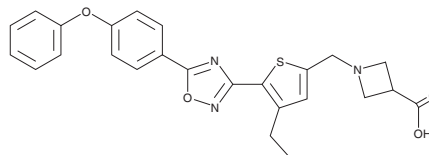
PRODUCT INFORMATION



CS 2100

Item No. 14396

CAS Registry No.: 913827-99-3
Formal Name: 1-[[4-ethyl-5-[5-(4-phenoxyphenyl)-1,2,4-oxadiazol-3-yl]-2-thienyl]methyl]-3-azetidincarboxylic acid
MF: C₂₅H₂₃N₃O₄S
FW: 461.5
Purity: ≥98%
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

CS 2100 is supplied as a solid. A stock solution may be made by dissolving the CS 2100 in the solvent of choice, which should be purged with an inert gas. CS 2100 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

CS 2100 is a sphingosine-1-phosphate receptor 1 (S1P₁) agonist.^{1,2} It selectively induces GTPγS binding in CHO-K1 cells expressing human or rat S1P₁ (EC₅₀s = 4 and 1.5 nM, respectively) over those expressing human or rat S1P₃ (EC₅₀s = >20 and 7.4 μM, respectively).² CS 2100 (0.1 and 1 mg/kg) transiently reduces blood lymphocyte levels in rats with the levels returning to baseline within 24-48 hours. It reduces hind paw volume in a rat model of adjuvant-induced arthritis (ID₅₀ = 0.44 mg/kg). CS 2100 (0.3 and 1 mg/kg) also decreases severity in a mouse model of experimental autoimmune encephalomyelitis (EAE).

References

1. Nakamura, T., Asano, M., Sekiguchi, Y., *et al.* Synthesis and evaluation of CS-2100, a potent, orally active and S1P₃- sparing S1P₁ agonist. *Eur. J. Med. Chem.* **51**, 92-98 (2012).
2. Nakamura, T., Asano, M., Sekiguchi, Y., *et al.* Discovery of CS-2100, a potent, orally active and S1P₃-sparing S1P₁ agonist. *Bioorg. Med. Chem. Lett.* **22**(4), 1788-1792 (2012).

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, 09/11/2024

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