

PRODUCT INFORMATION

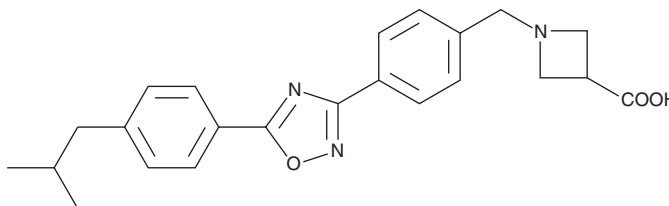


CAY10734

Item No. 10007833

CAS Registry No.: 635701-59-6
Formal Name: 1-[[4-[5-[4-(2-methylpropyl)phenyl]-1,2,4-oxadiazol-3-yl]phenyl]methyl]-3-azetidincarboxylic acid

MF: C₂₃H₂₅N₃O₃
FW: 391.5
Purity: ≥98%
UV/Vis.: λ_{max}: 251 nm
Supplied as: A crystalline 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

CAY10734 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10734 in the solvent of choice, which should be purged with an inert gas. CAY10734 is soluble in the organic solvent dimethyl formamide (DMF) at a concentration of approximately 5 mg/ml. CAY10734 is also slightly soluble in ethanol.

CAY10734 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10734 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10734 has a solubility of approximately 0.1 mg/ml in a 1:10 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10734 is an agonist of sphingosine-1-phosphate receptor 1 (S1P₁; IC₅₀ = 0.6 nM).¹ It selectively binds S1P₁ over S1P₂, S1P₃, and S1P₄ receptors (IC₅₀s = >10,000, 12,000, and 70 nM, respectively) but does also bind S1P₅ receptors (IC₅₀ = 1 nM) in radioligand binding assays. CAY10734 reduces the number of peripheral blood lymphocytes in mice, rats, and dogs, with the maximal response achieved when administered at doses of 10, 0.5, and 0.5 mg/kg, respectively. It increases graft survival in a rat skin allograft model when administered at a dose of 5 mg/kg per day in combination with cyclosporin A (CsA; Item No. 12088).

Reference

1. Li, Z., Chen, W., Hale, J.J., *et al.* Discovery of potent 3,5-diphenyl-1,2,4-oxadiazole sphingosine-1-phosphate-1 (S1P₁) receptor agonists with exceptional selectivity against S1P₂ and S1P₃. *J. Med. Chem.* **48**(20), 6169-6173 (2005).

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, 11/09/2022

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