

PRODUCT INFORMATION

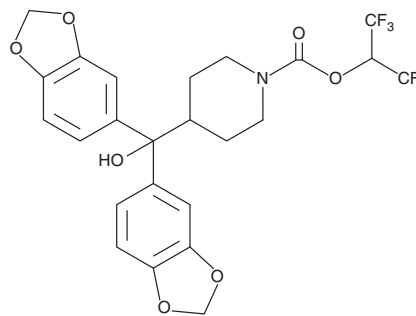


KML29

Item No. 11777

Sold under license from The Scripps Research Institute

CAS Registry No.: 1380424-42-9
Formal Name: 1,1,1,3,3,3-hexafluoropropan-2-yl
4-(bis(benzo[d][1,3]dioxol-5-yl)(hydroxy)
methyl)piperidine-1-carboxylate
MF: C₂₄H₂₁F₆NO₇
FW: 549.4
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 287 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

KML29 is supplied as a crystalline solid. A stock solution may be made by dissolving the KML29 in the solvent of choice, which should be purged with an inert gas. KML29 is soluble in the organic solvent DMSO at a concentration of approximately 2 mg/ml.

Description

Endocannabinoids such as 2-arachidonoyl glycerol (2-AG) and arachidonoyl ethanolamide are biologically active lipids that are involved in a number of synaptic processes including activation of cannabinoid receptors. Monoacylglycerol lipase (MAGL) is a serine hydrolase responsible for the hydrolysis of 2-AG to arachidonic acid and glycerol, thus terminating its biological function. KML29 is an O-hexafluoroisopropyl carbamate analog of JZL 184 (Item No. 13158) that potently and selectively inhibits MAGL (IC₅₀s = 15, 43, and 5.9 nM in mouse, rat, and human brain proteomes, respectively) over FAAH (IC₅₀s > 50 μM).¹ At 5-20 mg/kg, KML29 dose-dependently blocks mouse brain MAGL activity *in vivo*, without any measurable effect on FAAH activity.¹ As a second generation MAGL inhibitor, KML29 supersedes the low-level cross reactivity that JZL 184 displays for FAAH yet still maintains comparable potency to its parent compound.

Reference

1. Chang, J.W., Niphakis, M.J., Lum, K.M., *et al.* Highly selective inhibitors of monoacylglycerol lipase bearing a reactive group that is bioisosteric with endocannabinoid substrates. *Chem. Biol.* **19**(5), 579-588 (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, 11/16/2022

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