

PRODUCT INFORMATION



O-7460

Item No. 14009

CAS Registry No.: 1572051-31-0
Formal Name: 2-[(fluoromethylphosphinyl)oxy]-1-
[(1-methylethoxy)methyl]ethyl ester,
9Z-octadecenoic acid

MF: C₂₅H₄₈FO₅P

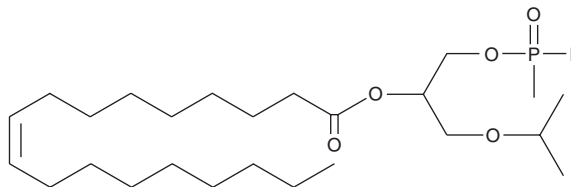
FW: 478.6

Purity: ≥95%

Supplied as: A 5 mg/ml solution in methyl acetate

Storage: -20°C

Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

In humans, two forms of diacylglycerol lipase, DAGL α and DAGL β , generate the endocannabinoid 2-arachidonoyl glycerol (2-AG; Item No. 62160) by attacking DAG at the *sn*-1 position. O-7460 is a selective inhibitor of 2-AG biosynthesis *via* DAGL α (IC₅₀ = 690 nM).¹ It demonstrates much weaker inhibition towards human monoacylglycerol lipase and rat brain fatty acid amide hydrolase (IC₅₀s > 10 μ M) and does not bind to CB₁ or CB₂ cannabinoid receptors (K_is > 10 μ M).¹ At 0-12 mg/kg, i.p. in mice, O-7460 was reported to dose-dependently inhibit high-fat diet intake and reduce body weight.¹

Reference

1. Bisogno, T., Mahadevan, A., Coccurello, R., *et al.* A novel fluorophosphonate inhibitor of the biosynthesis of the endocannabinoid 2-arachidonoylglycerol with potential anti-obesity effects. *Br. J. Pharmacol.* **169**(4), 784-793 (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, 12/08/2025

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