

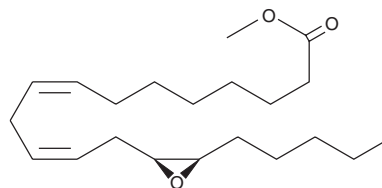
PRODUCT INFORMATION



(±)14(15)-EpEDE methyl ester

Item No. 10007528

Formal Name: (±)(8Z,11Z)-methyl 13-(3-pentylloxiran-2-yl)trideca-8,11-dienoate
Synonym: (±)14,15-Epoxyeicosadienoic Acid methyl ester
MF: C₂₁H₃₆O₃
FW: 336.5
Purity: ≥90%
Supplied as: A solution in hexane
Storage: -20°C
Stability: ≥2 years



NOTE: Relative stereochemistry shown in chemical structure

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

Laboratory Procedures

(±)14(15)-EpEDE methyl ester is supplied as a solution in hexane. To change the solvent, simply evaporate the hexane under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of (±)14(15)-EpEDE methyl ester in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

(±)14(15)-EpEDE methyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the hexane solution of (±)14(15)-EpEDE methyl ester should be diluted with the aqueous buffer of choice. (±)14(15)-EpEDE methyl ester has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(±)14(15)-EpEDE (Item No. 10007527) is an EpEDE formed from 8,11,14-eicosatrienoic acid, also known as dihomo-γ-linolenic acid (DGLA; Item No. 90230). (±)14(15)-EpEDE methyl ester is an analog of (±)14(15)-EpEDE with greater lipid solubility than the free base. The methyl ester may be more amenable for the formulation of fatty acid-containing diets and dietary supplements.

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, 02/15/2024

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