

PRODUCT INFORMATION



(±)14(15)-EET-d₁₁ methyl ester Item No. 29383

Formal Name: (±)14,15-epoxy-5Z,8Z,11Z-eicosatrienoic acid-d₁₁, methyl ester

Synonym: (±)14,15-EpETrE-d₁₁ methyl ester

MF: C₂₁H₂₃D₁₁O₃

FW: 345.6

Chemical Purity: ≥98% ((±)14(15)-EET methyl ester)

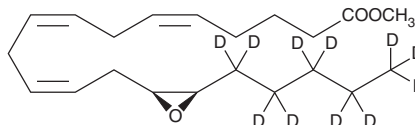
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₁₁); ≤1% d₀

Supplied as: A solution in ethanol

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)-EET-d₁₁ methyl ester is intended for use as an internal standard for the quantification of (±)14(15)-EET methyl ester (Item No. 50650) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

(±)14(15)-EET-d₁₁ methyl ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of (±)14(15)-EET-d₁₁ methyl ester in these solvents is approximately 50 mg/ml.

Description

(±)14(15)-EET methyl ester (Item No. 50650) is an esterified form of (±)14(15)-EET (Item No. 50651). 14(S),15(R)-EET and 14(R),15(S)-EET are formed via epoxidation of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) by a variety of cytochrome P450 (CYP) isoforms.¹ 14(S),15(R)-EET is produced at a higher proportion by CYP2C23, whereas 14(R),15(S)-EET is produced at a greater proportion by CYP2B2 and CYP2C24.

Reference

1. Capdevila, J.H., Falck, J.R., and Harris, R.C. Cytochrome P450 and arachidonic acid bioactivation: Molecular and functional properties of the arachidonate monooxygenase. *J. Lipid Res.* **41(2)**, 163-181 (2000).

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, 01/30/2024

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