

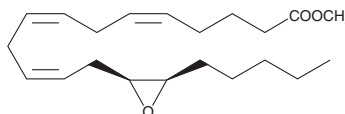
PRODUCT INFORMATION



14S(15R)-EET methyl ester

Item No. 28631

CAS Registry No.: 110901-52-5
Formal Name: (5Z,8Z,11Z)-13-[(2S,3R)-3-pentyloxiranyl]-5,8,11-tridecatricenoic acid, methyl ester
Synonym: 14S,15R-EpETrE
MF: C₂₁H₃₄O₃
FW: 334.5
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Description

14S(15R)-EET methyl ester is an esterified form of 14S(15R)-EET, which is an oxylipin and cytochrome P450 (CYP) metabolite of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607).¹ 14S(15R)-EET methyl ester is formed from 14S(15R)-EET to aid in the separation of unreacted arachidonic acid following the enzymatic epoxidation of arachidonic acid to 14S(15R)-EET.²

References

1. Daikh, B.E., Lasker, J.M., Raucy, J.L., *et al.* Regio- and stereoselective epoxidation of arachidonic acid by human cytochromes P450 2C8 and 2C91. *J. Pharmacol. Exp. Ther.* **271**(3), 1427-1433, 1994.
2. Woodman, J.W., Cinelli, M.A., Scharmen-Burgdorf, A., *et al.* Enzymatic synthesis of epoxidized metabolites of docosahexaenoic, eicosapentaenoic, and arachidonic acids. *J. Vis. Exp.* **148**, 10.3791/59770 (2019).

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, 10/18/2023

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