

PRODUCT INFORMATION



14R(15S)-EET Item No. 9004351

CAS Registry No.: 98103-48-1
Formal Name: (5Z,8Z,11Z)-13-[(2R,3S)-3-pentyl-2-oxiranyl]-5,8,11-tridecatricarboxylic acid

Synonyms: 14R(15S)-EpETrE,
14R(15S)-epoxy-5(Z),8(Z),11(Z)-ETrE,
14R(15S)-epoxy-all-cis-5,8,11-Eicosatrienoic Acid, 14R(15S)-Epoxyeicosatrienoic Acid,
FA 20:4;O

MF: C₂₀H₃₂O₃

FW: 320.5

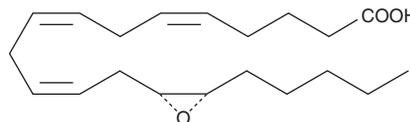
Purity: ≥95%

Supplied as: A 250 µg/ml solution in ethanol

Storage: -20°C

Stability: ≥1 year

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



Laboratory Procedures

14R(15S)-EET 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, which should be purged with an inert gas. 14R(15S)-EET is sparingly soluble (1-10 mg/ml) in acetonitrile.

Description

14R(15S)-EET is an oxylipin and a metabolite of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607).¹ It is formed *via* oxidation of arachidonic acid by the cytochrome P450 (CYP) isoforms CYP2C8 and CYP2C9. 14R(15S)-EET (10 µM) induces relaxation of precontracted isolated bovine coronary arteries.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. Falck, J.R., Krishna, U.M., Reddy, Y.K., *et al.* Comparison of vasodilatory properties of 14,15-EET analogs: Structural requirements for dilation. *Am. J. Physiol. Heart Circ. Physiol.* **284**(1), 337-349 (2003).

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, 11/24/2025

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