

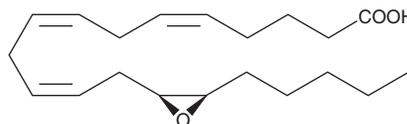
# PRODUCT INFORMATION



## 14S(15R)-EET

Item No. 9004352

**CAS Registry No.:** 105304-92-5  
**Formal Name:** 13-[(2S,3R)-3-pentyl-2-oxiranyl]-5Z,8Z,11Z-tridecatrienoic acid  
**Synonyms:** 14S(15R)-EpETrE, 14S(15R)-epoxy-all-cis-5,8,11 Eicosatrienoic Acid, 14S(15R)-Epoxyeicosatrienoic Acid, 14S(15R)-epoxy-5(Z),8(Z),11(Z)-ETrE, FA 20:4;O  
**MF:** C<sub>20</sub>H<sub>32</sub>O<sub>3</sub>  
**FW:** 320.5  
**Purity:** ≥95%  
**Supplied as:** A 250 µg/ml solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 years



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

### Laboratory Procedures

14S(15R)-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. 14S(15R)-EET is sparingly soluble (1-10 mg/ml) in acetonitrile

### Description

14S(15R)-EET is an oxylipin and a metabolite of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607).<sup>1</sup> It is formed *via* oxidation of arachidonic acid by the cytochrome P450 (CYP) isoforms CYP2C8 and CYP2C9. 14S(15R)-EET (1 µM) induces relaxation in precontracted isolated bovine coronary arteries.<sup>2</sup>

### 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