

# PRODUCT INFORMATION

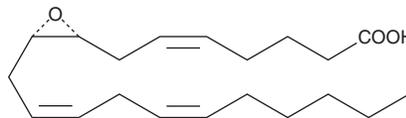


## 8R(9S)-EET

Item No. 9004347

**CAS Registry No.:** 129172-62-9  
**Formal Name:** (5Z)-7-[(2R,3S)-3-(2Z,5Z)-2,5-undecadien-

**Synonyms:** 8R(9S)-epoxy-all-cis-5,11,14-Eicosatrienoic Acid, 8R(9S)-EpETrE, 8R(9S)-Epoxyeicosatrienoic Acid, 8R(9S)-epoxy-5(Z),11(Z),14(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:** ≥2 years

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

### Laboratory Procedures

8R(9S)-EET is supplied as a solution in ethanol. A stock solution may be made by dissolving the 8R(9S)-EET in the solvent of choice, which should be purged with an inert gas. 8R(9S)-EET is sparingly soluble (1-10 mg/ml) in ethanol and acetonitrile.

### Description

8R(9S)-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) isoform CYP2C11. 8R(9S)-EET induces relaxation of precontracted isolated dog epicardial arterioles (EC<sub>50</sub> = 15 pM).<sup>2</sup>

### References

1. Capdevila, J.H., Karara, A., Waxman, D.J., *et al.* Cytochrome P-450 enzyme-specific control of the regio- and enantiofacial selectivity of the microsomal arachidonic acid epoxygenase. *J. Biol. Chem.* **265**(19), 10865-10871 (1990).
2. Zhang, Y., Oltman, C.L., Lu, T., *et al.* EET homologs potently dilate coronary microvessels and activate BK<sub>Ca</sub> channels. *Am. J. Physiol. Heart Circ. Physiol.* **280**(6), H2430-H2440 (2001).

#### 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