

PRODUCT INFORMATION

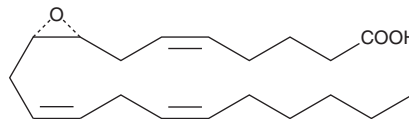


8R(9S)-EET

Item No. 9004347

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

Synonyms: 1-yl-2-oxiranyl]-5-heptenoic acid
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: C₂₀H₃₂O₃
FW: 320.5
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.

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).¹ 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 (EC₅₀ = 15 pM).²

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_{Ca} 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, 04/09/2024

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