

PRODUCT INFORMATION



5S(6R)-EET

Item No. 9004349

CAS Registry No.: 98103-47-0
Formal Name: (2S,3R)-3-(2Z,5Z,8Z)-2,5,8-tetradecatrien-1-yl-2-oxiranebutanoic acid

Synonyms: 5S(6R)-EpETrE,
5S(6R)-epoxy-8(Z),11(Z),14(Z)-
ETrE, 5S(6R)-epoxy-all-cis-8,11,14-
Eicosatrienoic Acid,
5S(6R)-Epoxyeicosatrienoic Acid,
FA 20:4;O

MF: C₂₀H₃₂O₃

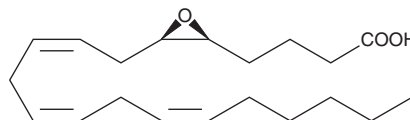
FW: 320.5

Purity: ≥95%

Supplied as: A 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

5S(6R)-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. 5S(6R)-EET is sparingly soluble (1-10 mg/ml) in acetonitrile.

Description

5S(6R)-EET is an oxylin. It depolarizes transepithelial voltage in isolated rabbit renal cortical collecting duct when applied to the luminal side at a concentration of 1 μM.¹ In solution, 5(6)-EET degrades into 5,6-DiHET and 5(6)-δ-lactone, which can be converted to 5(6)-DiHET and quantified by GC-MS.²

References

1. Sakairi, Y., Jacobson, H.R., Noland, T.D., *et al.* 5,6-EET inhibits ion transport in collecting duct by stimulating endogenous prostaglandin synthesis. *Am. J. Physiol.* **268(5 Pt 2)**, F931-F939 (1995).
2. Rashid, M., Manivet, P., Nishio, H., *et al.* Identification of the binding sites and selectivity of sarpogrelate, a novel 5-HT₂ antagonist, to human 5-HT_{2A}, 5-HT_{2B} and 5-HT_{2C} receptor subtypes by molecular modeling. *Life Sci.* **73(2)**, 193-207 (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, 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