

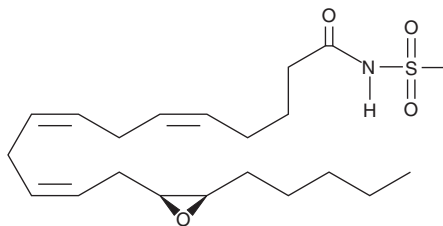
Product Information



(±)14(15)-EET-SI

Catalog No. 10009286

CAS Registry No.: 218461-97-3
Formal Name: N-(methylsulfonyl)-13-(3-pentylxiranyl)-5Z,8Z,11Z-tridecatrienamide
Synonyms: (±)14,15-EpETrE-SI, (±)14(15)-EET-sulfonimide
MF: C₂₁H₃₅NO₄S
FW: 397.6
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that (±)14(15)-EET-SI be stored as supplied at -20°C. It should be stable for at least one year.

(±)14(15)-EET-SI is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the (±)14(15)-EET-SI under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of (±)14(15)-EET-SI in these solvents is approximately 10 mg/ml.

(±)14(15)-EET-SI is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of (±)14(15)-EET-SI should be diluted with the aqueous buffer of choice. (±)14(15)-EET-SI has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Arachidonic acid is metabolized in the vascular endothelium to epoxytrienoic acids (EETs; EETs) by cytochrome P450 enzymes. The EETs are released in response to acetylcholine, bradykinin, arachidonic acid, or cyclic stretch.¹ (±)14(15)-EET-SI is the methyl sulfonamide analog of (±)14(15)-EET. This substitution results in a metabolically more stable compound because it is not sensitive to β-oxidation or membrane esterification. (±)14(15)-EET-SI is equipotent to (±)14(15)-EET in vascular agonist activity as measured by relaxation of precontracted bovine coronary arteries.¹ In addition, 14(15)-EET and the methyl sulfonamide analog both stimulate tyrosine phosphorylation and induce mitogenesis in renal epithelial cells.²

References

- Gauthier, K.M., Falck, J.R., Reddy, L.M., *et al.* 14,15-EET analogs: Characterization of structural requirements for agonist and antagonist activity in bovine coronary arteries. *Pharmacol. Res.* **49**, 515-524 (2004).
- Chen, J.-K., Falck, J.R., Reddy, K.M., *et al.* Epoxyicosatrienoic acids and their sulfonimide derivatives stimulate tyrosine phosphorylation and induce mitogenesis in renal epithelial cells. *J. Biol. Chem.* **273(44)**, 29254-29261 (1998).

Related Products

(±)14(15)-EET methyl ester - Cat. No. 50650 • (±)14(15)-EET - Cat. No. 50651

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 11/17/2009

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com