

# PRODUCT INFORMATION



## (±)11(12)-EET MaxSpec® Standard

Item No. 10007262

**Formal Name:** (±)11,12-epoxy-5Z,8Z,14Z-eicosatrienoic acid

**Synonym:** (±)11,12-EpETrE

**MF:** C<sub>20</sub>H<sub>32</sub>O<sub>3</sub>

**FW:** 320.5

**Purity:** ≥95%

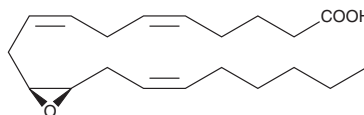
**Supplied as:** A solution in ethanol; in a deactivated glass ampule

**Concentration:** 100 µg/ml (nominal); see certificate of analysis for verified concentration

**Storage:** -20°C

**Stability:** ≥5 years; *Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and product expiry date will be updated upon completion of testing.*

**Special Conditions:** Store upright and unopened at -20°C. Warm to room temperature prior to opening. Light sensitive.



NOTE: Relative stereochemistry shown in chemical structure

### Description

(±)11(12)-EET is biosynthesized in rat and rabbit liver microsomes by CYP450.<sup>1,2</sup> (±)11(12)-EET has been shown, along with (±)8(9)-EET (Item No. 50351), to play a role in the recovery of depleted Ca<sup>2+</sup> pools in cultured smooth muscle cells.<sup>3</sup>

(±)11(12)-EET MaxSpec® standard is a quantitative grade standard of (±)11(12)-EET (Item No. 50511) that has been prepared specifically for mass spectrometry or any application where quantitative reproducibility is required. The solution has been prepared gravimetrically and is supplied in a deactivated glass ampule sealed under argon. The concentration was verified by comparison to an independently prepared calibration standard. This (±)11(12)-EET MaxSpec® standard is guaranteed to meet identity, purity, stability, and concentration specifications and is provided with a batch-specific certificate of analysis. Ongoing stability testing is performed to ensure the concentration remains accurate throughout the shelf life of the product.

**Note:** *The amount of solution added to the vial is in excess of the listed amount. Therefore, it is necessary to accurately measure volumes for preparation of calibration standards. Follow recommended storage and handling conditions to maintain product quality.*

### References

1. Chacos, N., Falck, J.R., Wixtrom, C., *et al.* Novel epoxides formed during the liver cytochrome P-450 oxidation of arachidonic acid. *Biochem. Biophys. Res. Commun.* **104(3)**, 916-922 (1982).
2. Oliw, E.H., Guengerich, F.P., and Oates, J.A. Oxygenation of arachidonic acid by hepatic monooxygenases. Isolation and metabolism of four epoxide intermediates. *J. Biol. Chem.* **257(7)**, 3771-3781 (1982).
3. Graber, M.N., Alfonso, A., and Gill, D.L. Recovery of Ca<sup>2+</sup> pools and growth in Ca<sup>2+</sup> pool-depleted cells is mediated by specific epoxyeicosatrienoic acids derived from arachidonic acid. *J. Biol. Chem.* **272(47)**, 29546-29553 (1997).

#### 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, 03/09/2021

#### 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