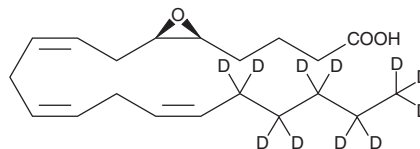


# PRODUCT INFORMATION



(±)5(6)-EET-d<sub>11</sub>  
Item No. 10009984

**Formal Name:** (±)5(6)-epoxy-8Z,11Z,14Z-eicosatrienoic-16,16,17,17,18,18,19,19,20,20,20-d<sub>11</sub> acid  
**Synonym:** (±)5,6-EET-d<sub>11</sub>  
**MF:** C<sub>20</sub>H<sub>21</sub>D<sub>11</sub>O<sub>3</sub>  
**FW:** 331.5  
**Chemical Purity:** ≥95% (mixture of free acid and lactone)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>11</sub>); ≤1% d<sub>0</sub>  
**Supplied as:** A solution in methyl acetate  
**Storage:** -80°C  
**Stability:** ≥6 months



NOTE: Relative stereochemistry shown in chemical structure

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

## Laboratory Procedures

(±)5(6)-EET-d<sub>11</sub> is intended for use as an internal standard for the quantification of 5(6)-EET by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

(±)5(6)-EET-d<sub>11</sub> is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate 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 (±)5(6)-EET-d<sub>11</sub> in these solvents is approximately 50 mg/ml.

## Description

5(S),6(R)-EET and 5(R),6(S)-EET are formed *via* epoxidation of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) by a variety of cytochrome P450 (CYP) isoforms, including CYP2B1, CYP2B2, CYP2C11, CYP2C23, and CYP2C24.<sup>1,2</sup>

## References

1. Spector, A.A. and Norris, A.W. Action of epoxyeicosatrienoic acids on cellular function. *Am. J. Physiol. Cell Physiol.* **292**(3), C996-C1012 (2007).
2. Capdevila, J.H., Falck, J.R., and Harris, R.C. Cytochrome P450 and arachidonic acid bioactivation: Molecular and functional properties of the arachidonate monooxygenase. *J. Lipid Res.* **41**(2), 163-181 (2000).

### 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, 02/25/2021

## 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