

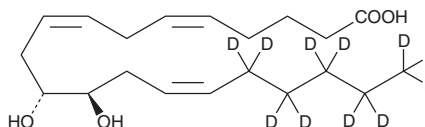
PRODUCT INFORMATION



(±)11(12)-DiHET-d₁₁

Item No. 10007975

Formal Name:	(±)11,12-dihydroxy-5Z,11Z,14Z-eicosatrienoic-16,16,17,17,18,18,19,19,20,20,20-d ₁₁ acid
Synonyms:	(±)11(12)-DHET-d ₁₁ , (±)11,12-DiHETrE-d ₁₁
MF:	C ₂₀ H ₂₃ D ₁₁ O ₄
FW:	349.6
Chemical Purity:	≥95%
Deuterium Incorporation:	≥99% deuterated forms (d ₁ -d ₁₁); ≤1% d ₀
Supplied as:	A solution in ethanol
Storage:	-20°C
Stability:	≥1 year



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

It is intended for use as an internal standard for the quantification of (±)11(12)-DiHET (Item No. 51511) 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).

(±)11(12)-DiHET-d₁₁ 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. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of (±)11(12)-DiHET-d₁₁ in these solvents is approximately 50 mg/ml.

Description

11(S),12(S)-DiHET and 11(R),12(R)-DiHET are vicinal diols formed via enzymatic hydration of 11(12)-EET by cytosolic or soluble epoxide hydrolases in a non-stereoselective manner.¹⁻³

References

1. Fang, X., Kaduce, T.L., Weintraub, N.L., *et al.* Functional implications of a newly characterized pathway of 11,12-epoxyeicosatrienoic acid metabolism in arterial smooth muscle. *Circ. Res.* **79**, 784-793 (1996).
2. Zeldin, D.C., Kobayashi, J., Falck, J.R., *et al.* Regio- and enantiofacial selectivity of epoxyeicosatrienoic acid hydration by cytosolic epoxide hydrolase. *J. Biol. Chem.* **268**(9), 6402-6407 (1993).
3. Zhang, G., Kodani, S., and Hammock, B.D. Stabilized epoxygenated fatty acids regulate inflammation, pain, angiogenesis and cancer. *Prog. Lipid Res.* **53**, 108-123 (2014).

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, 05/18/2022

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