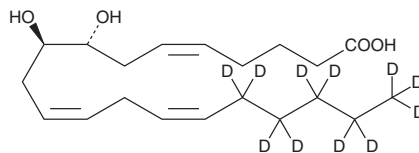


PRODUCT INFORMATION



(±)8(9)-DiHET-d₁₁ Item No. 10009998

Formal Name:	(±)8,9-dihydroxy-5Z,11Z,14Z-eicosatrienoic-16,16,17,17,18,18,19,19,20,20,20-d ₁₁ acid
Synonym:	(±)8,9-DiHETrE-d ₁₁
MF:	C ₂₀ H ₂₃ D ₁₁ O ₄
FW:	349.6
Chemical Purity:	≥95% (8(9)-DiHET)
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

(±)8(9)-DiHET-d₁₁ is intended for use as an internal standard for the quantification of 8(9)-DiHET 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).

(±)8(9)-DiHET-d₁₁ is supplied as a solution in ethanol. A stock solution may be made by dissolving the (±)8(9)-DiHET-d₁₁ in the solvent of choice, which should be purged with an inert gas. (±)8(9)-DiHET-d₁₁ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (±)8(9)-DiHET-d₁₁ in these solvents is approximately 50 mg/ml.

Description

8(S),9(S)-DiHET and 8(R),9(R)-DiHET are vicinal diols formed via enzymatic hydration of 8(9)-EET by cytosolic or soluble epoxide hydrolases.^{1,2} 8(S),9(S)-DiHET is produced at a greater proportion than 8(R),9(R)-DiHET by cytosolic epoxide hydrolase.¹

References

1. 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).
2. Zhang, J.Y., Prakash, C., Yamashita, K., *et al.* Regiospecific and enantioselective metabolism of 8,9-epoxyeicosatrienoic acid by cyclooxygenase. *Biochem. Biophys. Res. Commun.* **183**, 138-143 (1992).

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