

PRODUCT INFORMATION



(±)11-HETE-d₈ Item No. 9002385

Formal Name: (±)11-hydroxy-5Z,8Z,12E,14Z-
eicosatetraenoic-5,6,8,9,11,12,14,15-d₈ acid

Synonym: (±)11-Hydroxyeicosatetraenoic Acid-d₈

MF: C₂₀H₂₄D₈O₃

FW: 328.5

Chemical Purity: ≥98% ((±)11-HETE)

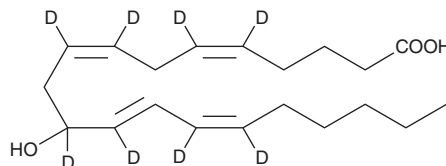
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀

UV/Vis.: λ_{max}: 234 nm

Supplied as: A 100 µg/ml solution in acetonitrile

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

(±)11-HETE-d₈ is intended for use as an internal standard for the quantification of 11-HETE 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-HETE-d₈ is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of (±)11-HETE-d₈ in ethanol is approximately 50 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Description

(±)11-HETE is formed non-enzymatically from arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607).^{1,2} 11(R)-HETE is formed from arachidonic acid by COX-1, COX-2, and aspirin-acetylated COX-2.^{3,4} 11(R)- and 11(S)-HETE are formed from arachidonic acid via cytochrome P450 (CYP), with 11(R)-HETE formed at a higher ratio than 11(S)-HETE.⁵

References

- Powell, W.S. and Rokach, J. Biosynthesis, biological effects, and receptors of hydroxyeicosatetraenoic acids (HETEs) and oxoeicosatetraenoic acids (oxo-ETEs) derived from arachidonic acid. *Biochim. Biophys. Acta* **1851(4)**, 340-355 (2014).
- Derogis, P.B.M.C., Chaves-Filho, A.B., and Miyamoto, S. Characterization of hydroxy and hydroperoxy polyunsaturated fatty acids by mass spectrometry. *Bioactive lipids in health and disease*. Trostchansky, A. and Rubbo, H., editors, *Springer* (2019).
- Thuresson, E.D., Lakkides, K.M., and Smith, W.L. Different catalytically competent arrangements of arachidonic acid within the cyclooxygenase active site of prostaglandin endoperoxide H synthase-1 lead to the formation of different oxygenated products. *J. Biol. Chem.* **275(12)**, 8501-8507 (2000).
- Xiao, G., Tsai, A.I., Palmer, G., et al. Analysis of hydroperoxide-induced tyrosyl radicals and lipoxygenase activity in aspirin-treated human prostaglandin H synthase-2. *Biochemistry* **36(7)**, 1836-1845 (1997).
- Capdevila, J., Yadagiri, P., Manna, S., et al. Absolute configuration of the hydroxyeicosatetraenoic acids (HETEs) formed during catalytic oxygenation of arachidonic acid by microsomal cytochrome P-450. *Biochem. Biophys. Res. Commun.* **141(3)**, 1007-1011 (1986).

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, 05/05/2026

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