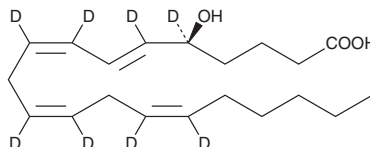


PRODUCT INFORMATION

5(S)-HETE-d₈ Item No. 334230

CAS Registry No.: 330796-62-8
Formal Name: 5S-hydroxy-6E,8Z,11Z,14Z-eicosatetraenoic-5,6,8,9,11,12,14,15-d₈ acid
Synonym: 5(S)-Hydroxyeicosatetraenoic Acid-d₈
MF: C₂₀H₂₄D₈O₃
FW: 328.5
Chemical Purity: ≥95% (5-HETE)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₈); ≤1% d₀
UV/Vis.: λ_{max}: 236 nm
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

5(S)-HETE-d₈ is intended for use as an internal standard for the quantification of 5-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).

5(S)-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 purged with an inert gas can be used. 5(S)-HETE-d₈ is miscible in these solvents. The solubility of 5(S)-HETE-d₈ in 0.1 M Na₂CO₃ is approximately 2 mg/ml.

Description

(±)5-HETE is formed *via* non-enzymatic oxidation of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607).¹ 5(S)- and 5(R)-HETE are formed by lipoxygenase-mediated oxidation of arachidonic acid.^{2,3}

References

1. Astarita, G., Kendall, A.C., Dennis, E.A., *et al.* Targeted lipidomics strategies for oxygenated metabolites of polyunsaturated fatty acid. *Biochim. Biophys. Acta* **1851**(4), 456-468 (2015).
2. Dodge, W. and Thomas, M. The effect of 5-hydroxyeicosatetraenoic acid on the proliferation of granulocyte progenitors and embryonic fibroblasts of the chick. *Biochem. Biophys. Res. Commun.* **131**(2), 731-735 (1985).
3. Hada, T., Swift, L.L., and Brash, A.R. Discovery of 5R-lipoxygenase activity in oocytes of the surf clam, *Spisula solidissima*. *Biochim. Biophys. Acta* **1346**(2), 109-119 (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, 11/10/2022

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