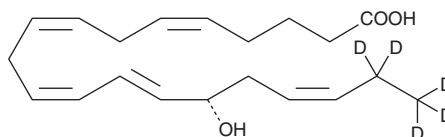


# PRODUCT INFORMATION

## 15(S)-HEPE-d<sub>5</sub> Item No. 31221

**CAS Registry No.:** 2750534-74-6  
**Formal Name:** 15(S)-hydroxy-5Z,8Z,11Z,13E,17Z-eicosapentaenoic-19,19,20,20,20-d<sub>5</sub> acid  
**Synonyms:** 15(S)-hydroxy Eicosapentaenoic Acid-d<sub>5</sub>, 15(S)-hydroxy EPA-d<sub>5</sub>  
**MF:** C<sub>20</sub>H<sub>25</sub>D<sub>5</sub>O<sub>3</sub>  
**FW:** 323.5  
**Chemical Purity:** ≥98% (15(S)-HEPE)  
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>5</sub>); ≤1% d<sub>0</sub>  
**UV/Vis.:** λ<sub>max</sub>: 237 nm  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

15(S)-HEPE-d<sub>5</sub> is intended for use as an internal standard for the quantification of 15-HEPE 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).

15(S)-HEPE-d<sub>5</sub> 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 DMSO and dimethyl formamide purged with an inert gas can be used.

### Description

(±)-15-HEPE and 15(S)-HEPE are formed *via* non-enzymatic and 15-lipoxygenase-mediated oxidation of eicosapentaenoic acid (Item Nos. 90110 | 90110.1 | 21908), respectively.<sup>1,2</sup>

### References

1. Miller, C., Yamaguchi, R.Y., and Ziboh, V.A. Guinea pig epidermis generates putative anti-inflammatory metabolites from fish oil polyunsaturated fatty acids. *Lipids* **24**(12), 998-1003 (1989).
2. Tan, L., Xin, X., Zhai, L., *et al.* *Drosophila* fed ARA and EPA yields eicosanoids, 15S-hydroxy-5Z,8Z, 11Z, 13E-eicosatetraenoic acid, and 15S-hydroxy-5Z,8Z,11Z,13E,17Z-eicosapentaenoic acid. *Lipids* **51**(4), 435-449 (2016).

#### 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, 01/25/2024

#### 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