

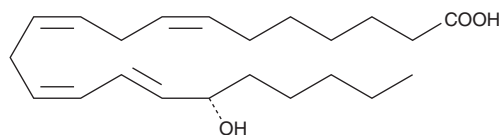
PRODUCT INFORMATION



17(S)-HDoTE

Item No. 34745

CAS Registry No.: 105835-39-0
Formal Name: 17S-hydroxy-7Z,10Z,13Z,15E-docosatetraenoic acid
Synonyms: 17(S)-HDTA, 17(S)-hydroxy Docosatetraenoic Acid
MF: C₂₂H₃₆O₃
FW: 348.5
Purity: ≥98%
UV/Vis.: λ_{max}: 236 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.

Description

17(S)-HDoTE is a metabolite of adrenic acid (Item No. 90300).¹ It is formed from adrenic acid by 15-lipoxygenase (15-LO) via a 17-HpDoTE intermediate.

Reference

1. Lee, I.-G., An, J.-U., Ko, Y.-J., *et al.* Enzymatic synthesis of new hepoxilins and trioxilins from polyunsaturated fatty acids. *Green Chem.* **21(11)**, 3172-3181 (2019).

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, 09/14/2021

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