

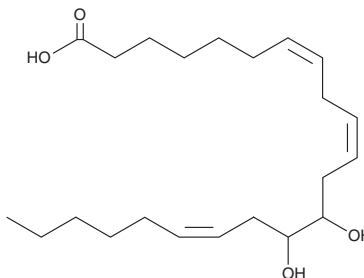
PRODUCT INFORMATION



(±)13,14-DiHDoTrE

Item No. 38938

CAS Registry No.: 944908-94-5
Formal Name: 13,14-dihydroxy-7Z,10Z,16Z-docosatrienoic acid
Synonyms: (±)13,14-DHDT,
(±)13,14-Dihydroxydocosatrienoic Acid
MF: C₂₂H₃₈O₄
FW: 366.5
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

(±)13,14-DiHDoTrE is an oxylipin and a racemic mixture of the (R)/(S) enantiomeric forms biosynthesized from the adrenic acid metabolite (±)13,14-EDT (Item No. 38937) by soluble epoxide hydrolase (sEH).¹

Reference

1. Singh, N., Barnych, B., Wagner, K.M., *et al.* Adrenic acid-derived epoxy fatty acids are naturally occurring lipids and their methyl ester prodrug reduces endoplasmic reticulum stress and inflammatory pain. *ACS Omega* **6**(10), 7165-7174 (2021).

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, 10/04/2023

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