

PRODUCT INFORMATION



(±)10,11-EDT

Item No. 38935

CAS Registry No.: 355016-20-5

Formal Name: (7Z)-rel-9-[(2R,3S)-3-(2Z,5Z)-2,5-undecadien-1-yl-2-oxiranyl]-7-nonenic acid

Synonyms: (±)-Dihomo-10,11-EET, (±)10,11-EpDoTrE, (±)10,11-Epoxydocosatrienoic Acid

MF: C₂₂H₃₆O₃

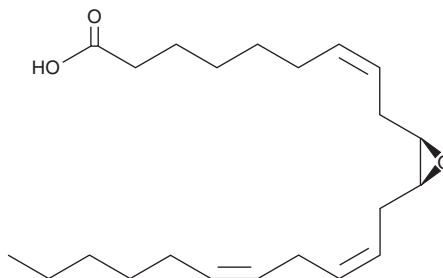
FW: 348.5

Purity: ≥98%

Supplied as: A 100 µg/ml solution in ethanol

Storage: -20°C

Stability: ≥4 years



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

Description

(±)10,11-EDT is an oxylipin and a metabolite of adrenic acid (Item No. 90300) formed via the cytochrome P450 (CYP) pathway.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, 07/21/2026

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