

PRODUCT INFORMATION



19S(20R)-EpDPA

Item No. 28627

CAS Registry No.: 2460392-58-7
Formal Name: 18-[(2S,3R)-3-ethyl-2-oxiranyl]-4Z,7Z,10Z,13Z,16Z-octadecapentaenoic acid
Synonyms: 19S,20R-EDP, 19S,20R-EpDPE, 19S,20R-epoxy DPA, 19S,20R-epoxy Docosapentaenoic Acid

MF: C₂₂H₃₂O₃

FW: 344.5

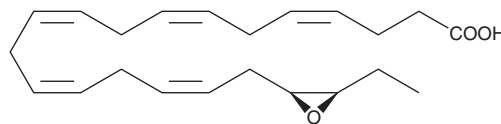
Purity: ≥90%

Supplied as: A 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

19S(20R)-EpDPA is an oxylipin and a metabolite of docosahexaenoic acid (DHA; Item No. 90310).^{1,2} It is formed from DHA by various cytochrome P450 (CYP) isoforms in a stereoselective manner.² Unlike 19R(20S)-EpDPA (Item No. 28628), 19S(20R)-EpDPA does not prevent calcium-induced increases in the spontaneous beating of isolated neonatal rat cardiomyocytes (NRCMs).³

References

1. Cinelli, M.A., Yang, J., Scharmen, A., *et al.* Enzymatic synthesis and chemical inversion provide both enantiomers of bioactive epoxydocosapentaenoic acids. *J. Lipid Res.* **59(11)**, 2237-2252 (2018).
2. Lucas, D., Goulitquer, S., Marienhagen, J., *et al.* Stereoselective epoxidation of the last double bond of polyunsaturated fatty acids by human cytochromes P450. *J. Lipid Res.* **51(5)**, 1125-1133 (2010).
3. Burke, J.E. and Dennis, E.A. Phospholipase A₂ structure/function, mechanism, and signaling. *J. Lipid Res.* **50(Suppl)**, S237-S242 (2009).

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, 01/22/2026

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