

PRODUCT INFORMATION



(±)16(17)-EpDPA MaxSpec® Standard

Item No. 26240

CAS Registry No.: 155073-46-4

Formal Name: (±)16,17-epoxy-4Z,7Z,10Z,13Z,19Z-docosapentaenoic acid

Synonyms: (±)16,17 EDP, (±)16,17-epoxy Docosapentaenoic Acid, (±)16,17-epoxy DPA, (±)16,17-EpDPE

MF: C₂₂H₃₂O₃

FW: 344.5

Purity: ≥95%

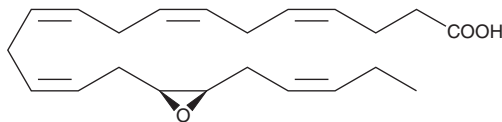
Supplied as: A solution in ethanol; in a deactivated glass ampule

Concentration: 100 µg/ml (nominal); see certificate of analysis for verified concentration

Storage: -20°C

Stability: ≥5 years; Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and product expiry date will be updated upon completion of testing.

Special Conditions: Store upright and unopened at -20°C. Warm to room temperature prior to opening. Light sensitive.



NOTE: Relative stereochemistry shown in chemical structure

Description

EDHF (endothelium-derived hyperpolarizing factor) is an unidentified mediator released from vascular endothelial cells in response to acetylcholine and bradykinin which is distinct from the NOS- (nitric oxide) and COX-derived (prostacyclin) vasodilators.^{1,2} Cytochrome P450 (CYP450) metabolism of polyunsaturated fatty acids produces epoxides such as (±)14(15)-EET (Item No. 50651) which are prime candidates for the actual active mediator.³ However, the CYP450 metabolites of eicosapentaenoic acid (EPA; Item No. 90110) and docosahexaenoic acid (DHA; Item No. 90310) have been little studied relative to arachidonate epoxigenase metabolites. (±)16(17)-EpDPA is the DHA homolog of (±)14(15)-EpETE, derived via epoxidation of the 16,17-double bond of DHA. The EDHF activity of (±)16(17)-EpDPA has not yet been determined. The epoxigenase metabolites of DHA have also been detected in a mouse inflammation model.⁴

(±)16(17)-EpDPA MaxSpec® standard is a quantitative grade standard of (±)16(17)-EpDPA (Item No. 10174) that has been prepared specifically for mass spectrometry or any application where quantitative reproducibility is required. The solution has been prepared gravimetrically and is supplied in a deactivated glass ampule sealed under argon. The concentration was verified by comparison to an independently prepared calibration standard. This (±)16(17)-EpDPA MaxSpec® standard is guaranteed to meet identity, purity, stability, and concentration specifications and is provided with a batch-specific certificate of analysis. Ongoing stability testing is performed to ensure the concentration remains accurate throughout the shelf life of the product. **Note:** The amount of solution added to the vial is in excess of the listed amount. Therefore, it is necessary to accurately measure volumes for preparation of calibration standards. Follow recommended storage and handling conditions to maintain product quality.

References

1. Chataigneau, T., Félétou, M., Duhault, J., et al. *Br. J. Pharmacol.* **123**(3), 574-580 (1998).
2. Fisslthaler, B., Popp, R., Kiss, L., et al. *Nature* **401**(6752), 493-497 (1999).
3. Baron, A., Frieden, M., and Bény, J.-L. *J. Physiol.* **504**(Pt 3), 537-543 (1997).
4. Serhan, C.N., Hong, S., Gronert, K., et al. *J. Exp. Med.* **196**(8), 1025-1037 (2002).

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, 03/23/2021

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