

PRODUCT INFORMATION



5(S),6(R)-DiHETE MaxSpec® Standard

Item No. 10007252

CAS Registry No.: 82948-88-7

Formal Name: 5S,6R-dihydroxy-7E,9E,11Z,14Z-eicosatetraenoic acid

MF: C₂₀H₃₂O₄

FW: 336.5

Purity: ≥95%

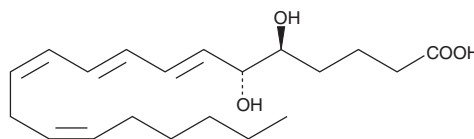
Supplied as: A solution in ethanol; in a deactivated glass ampule

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

Storage: -20°C

Stability: ≥2 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.



Description

5(S),6(R)-DiHETE is a dihydroxy polyunsaturated fatty acid and a nonenzymatic hydrolysis product of leukotriene A₄ (LTA₄). Mouse liver cytosolic epoxide hydrolase catalyzes the conversion of LTA₄ to 5(S),6(R)-DiHETE.¹ 5(S),6(R)-DiHETE is a weak LTD₄ (Item No. 20310) receptor agonist in guinea pig lung membranes.² It induces guinea pig ileum contraction with an ED₅₀ value of 1.3 µM.²

5(S),6(R)-DiHETE MaxSpec® standard is a quantitative grade standard of 5(S),6(R)-DiHETE (Item No. 35200) 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 5(S),6(R)-DiHETE 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. Haeggström, J., Wetterholm, A., Hamberg, M., et al. Enzymatic formation of 5,6-dihydroxy-7,9,11,14-eicosatetraenoic acid: Kinetics of the reaction and stereochemistry of the product. *Biochim. Biophys. Acta.* **958(3)**, 469-476 (1988).
2. Muller, A., Rechenq, E., Kugel, C., et al. Comparative biological activities of the four synthetic (5,6)-diHETE isomers. *Prostaglandins* **38(6)**, 635-644 (1989).

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, 01/07/2019

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