

PRODUCT INFORMATION



15(S)-HETE MaxSpec® Standard

Item No. 10007251

CAS Registry No.: 54845-95-3

Formal Name: 15S-hydroxy-5Z,8Z,11Z,13E-eicosatetraenoic acid

MF: C₂₀H₃₂O₃

FW: 320.5

Purity: ≥95%

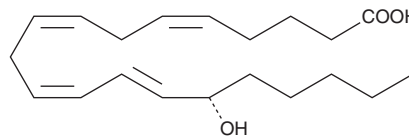
Supplied as: A solution in ethanol; in a deactivated glass ampule

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.



Description

15(S)-HETE is a major arachidonic acid metabolite from the 15-lipoxygenase pathway. In mammals, 15(S)-HETE is synthesized in the respiratory epithelium, leukocytes, and reticulocytes.¹ 15(S)-HETE is present in µg/ml concentrations in the nasal secretions of allergic rhinitis.²

15(S)-HETE MaxSpec® standard is a quantitative grade standard of 15(S)-HETE (Item No. 34720) 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 15(S)-HETE 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. Nadel, J.A., Conrad, D.J., Ueki, I.F., *et al.* Immunocytochemical localization of arachidonate 15-lipoxygenase in erythrocytes, leukocytes, and airway cells. *J. Clin. Invest.* **87(4)**, 1139-1145 (1991).
2. Ramis, I., Roselló-Catafau, J., Bulbena, O., *et al.* 15-Hydroxyeicosatetraenoic acid as a major eicosanoid in nasal secretions: Assay by high-performance liquid chromatographic-radioimmunoassay and gas chromatographic-mass spectrometric procedures. *J. Chromatogr.* **496(2)**, 416-422 (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, 05/19/2021

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