

PRODUCT INFORMATION



9(S),12(S),13(S)-TriHODE

Item No. 38950

CAS Registry No.: 95341-44-9

Formal Name: 9S,12S,13S-trihydroxy-10E,15Z-octadecadienoic acid

Synonyms: Corchorifatty Acid F, FA 18:2;O3, Fulgidic Acid, 9(S),12(S),13(S)-trihydroxy-10(E),15(Z)-ODE, 9(S),12(S),13(S)-Trihydroxyoctadecadienoic Acid, 9(S),12(S),13(S)-trihydroxy-*trans*-10,*cis*-15-Octadecadienoic Acid

MF: C₁₈H₃₂O₅

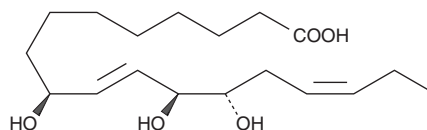
FW: 328.4

Purity: ≥95%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Description

9(S),12(S),13(S)-TriHODE is an α -linolenic acid-derived oxylipin that has been found in *C. olitorius*.¹⁻³ It is formed from α -linolenic acid (Item Nos. 90210 | 21910) via a 9(S)-HpOTrE (Item No. 45120) intermediate.¹ Tumor levels of 9(S),12(S),13(S)-TriHODE are decreased compared with adjacent healthy tissue in dogs with mammary cancer.⁴

References

1. Hamberg, M., Sanz, A., Rodriguez, M.J., *et al.* Activation of the fatty acid α -dioxygenase pathway during bacterial infection of tobacco leaves. Formation of oxylipins protecting against cell death. *J. Biol. Chem.* **278**(51), 51796-51805 (2003).
2. Yoshikawa, M., Murakami, T., Shimada, H., *et al.* Medicinal foodstuffs. XIV. On the bioactive constituents of moroheiya. (2): New fatty acids, corchorifatty acids A, B, C, D, E, and F, from the leaves of *Corchorus olitorius* L. (Tiliaceae): Structures and inhibitory effect on NO production in mouse peritoneal macrophages. *Chem. Pharm. Bull. (Tokyo)* **46**(6), 1008-1014 (1998).
3. Prost, I., Dhondt, S., Rothe, G., *et al.* Evaluation of the antimicrobial activities of plant oxylipins supports their involvement in defense against pathogens. *Plant Physiol.* **139**(4), 1902-1913 (2005).
4. Yu, C., Zheng, H.-H., Zhang, Y.-Z., *et al.* Identification of canine mammary tumor-associated metabolites using untargeted metabolomics. *Theriogenology* **211**, 84-96 (2023).

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, 03/18/2024

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