

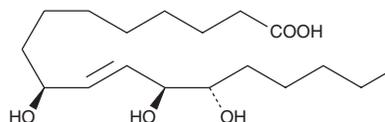
# PRODUCT INFORMATION



## 9(S),12(S),13(S)-TriHOME

Item No. 10005143

**CAS Registry No.:** 97134-11-7  
**Formal Name:** 9S,12S,13S-trihydroxy-10E-octadecenoic acid  
**Synonyms:** (-)-Pinellic Acid, 9S,12S,13S-Pinellic Acid  
**MF:** C<sub>18</sub>H<sub>34</sub>O<sub>5</sub>  
**FW:** 330.5  
**Purity:** ≥98%  
**Supplied as:** A 200 µg/ml 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

9(S),12(S),13(S)-TriHOME is a linoleic acid-derived oxylipin that has diverse biological activities.<sup>1-4</sup> It has been found in various plants and is produced in human eosinophils in a 15-lipoxygenase-dependent, soluble epoxide hydrolase-independent manner.<sup>1,5</sup> 9(S),12(S),13(S)-TriHOME inhibits antigen-induced β-hexosaminidase release from RBL-2H3 mast cells (IC<sub>50</sub> = 28.7 µg/ml).<sup>2</sup> It inhibits LPS-induced nitric oxide (NO) production in BV-2 microglia (IC<sub>50</sub> = 40.95 µM).<sup>3</sup> *In vivo*, 9(S),12(S),13(S)-TriHOME (1 g/animal) enhances the antiviral IgA and IgG antibody responses induced by a nasal influenza hemagglutinin (HA) vaccine by 5.2- and 2-fold, respectively, in mice.<sup>4</sup>

### References

1. Hamberg, M. and Hamberg, G. Peroxygenase-catalyzed fatty acid epoxidation in cereal seeds: Sequential oxidation of linoleic acid into 9(S),12(S),13(S)-trihydroxy-10(E)-octadecenoic acid. *Plant Physiol.* **110(3)**, 807-815 (1996).
2. Hong, S.S. and Oh, J.S. Inhibitors of antigen-induced degranulation of RBL-2H3 cells isolated from wheat bran. *J. Korean Soc. Appl. Biol. Chem.* **55**, 69-74 (2012).
3. Kim, C.S., Kwon, O.W., Kim, S.Y., *et al.* Five new oxylipins from *Chaenomeles sinensis*. *Lipids* **49(11)**, 1151-1159 (2014).
4. Shirahata, T., Sunazuka, T., Yoshida, K., *et al.* Total synthesis, elucidation of absolute stereochemistry, and adjuvant activity of trihydroxy fatty acids. *Tetrahedron* **62(40)**, 9483-9496 (2006).
5. Fuchs, D., Tang, X., Johnsson, A.-K., *et al.* Eosinophils synthesize trihydroxyoctadecenoic acids (TriHOMEs) via a 15-lipoxygenase dependent process. *Biochim. Biophys. Acta Mol. Cell Biol. Lipids* **1865(4)**, 158611 (2020).

#### 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, 05/21/2026

#### 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