

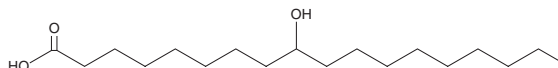
# PRODUCT INFORMATION



## 9-hydroxy Stearic Acid

Item No. 18016

**CAS Registry No.:** 3384-24-5  
**Formal Name:** 9-hydroxy-octadecanoic acid  
**Synonyms:** 9-HSA, 9-hydroxy Octadecanoic Acid  
**MF:** C<sub>18</sub>H<sub>36</sub>O<sub>3</sub>  
**FW:** 300.5  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

9-hydroxy Stearic acid is supplied as a solid. A stock solution may be made by dissolving the 9-hydroxy stearic acid in the solvent of choice, which should be purged with an inert gas. 9-hydroxy Stearic acid is soluble in the organic solvent chloroform.

### Description

9-hydroxy Stearic acid is a hydroxy fatty acid and an active metabolite of 9-PAHSA (Item No. 17037).<sup>1,2</sup> It is formed from 9-PAHSA by liver and pancreatic carboxyl ester lipases.<sup>2</sup> 9-hydroxy Stearic acid (5 μM) inhibits histone deacetylase 1 (HDAC1) in HT-29 colon cancer cell lysates.<sup>3</sup> It inhibits the proliferation of, and induces cell cycle arrest at the G<sub>0</sub>/G<sub>1</sub> phase in, HT-29 cells when used at a concentration of 100 μM.<sup>1</sup>

### References

1. Calonghi, N., Cappadone, C., Pagnotta, E., *et al.* 9-Hydroxystearic acid upregulates p21<sup>WAF1</sup> in HT29 cancer cells. *Biochem. Biophys. Res. Commun.* **314(1)**, 138-142 (2004).
2. Kolar, M.J., Kamat, S.S., Parsons, W.H., *et al.* Branched fatty acid esters of hydroxy fatty acids are preferred substrates of the MODY8 protein carboxyl ester lipase. *Biochemistry* **55(33)**, 4636-4641 (2016).
3. Calonghi, N., Cappadone, C., Pagnotta, E., *et al.* Histone deacetylase 1: A target of 9-hydroxystearic acid in the inhibition of cell growth in human colon cancer. *J. Lipid Res.* **46(8)**, 1596-1603 (2005).

#### 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, 11/03/2022

#### 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