

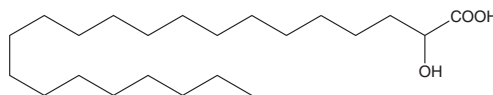
# PRODUCT INFORMATION



## 2-hydroxy Docosanoic Acid

Item No. 24594

**CAS Registry No.:** 13980-14-8  
**Formal Name:** 2-hydroxy-docosanoic acid  
**Synonyms:** 2-hydroxy Behenic Acid,  
α-hydroxy Behenic Acid,  
2-hydroxy DCA, α-hydroxy DCA,  
(±)-2-hydroxy DCA, (±)-α-hydroxy DCA,  
(±)-2-hydroxy Docosanoic Acid,  
(±)-α-hydroxy Docosanoic Acid



**MF:** C<sub>22</sub>H<sub>44</sub>O<sub>3</sub>  
**FW:** 356.6  
**Purity:** ≥98%  
**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

2-hydroxy Docosanoic acid is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy docosanoic acid in the solvent of choice. 2-hydroxy Docosanoic acid is soluble in a 5:1 solution of chloroform:methanol.

### Description

2-hydroxy Docosanoic acid is a hydroxylated fatty acid that has been found in the mature black epidermis of the Antarctic minke whale, as the N-acyl chain of galactosylceramides in mouse brain, and the inner bark of *E. globulus*.<sup>1-3</sup> It is upregulated in prostate carcinoma tissue as compared to normal prostate epithelium.<sup>4</sup>

### References

1. Yunoki, K., Ishikawa, H., Fukui, Y., *et al.* Chemical properties of epidermal lipids, especially sphingolipids, of the Antarctic minke whale. *Lipids* **43(2)**, 151-159 (2008).
2. Alderson, N.L., Maldonado, E.N., Kern, M.J., *et al.* FA2H-dependent fatty acid 2-hydroxylation in postnatal mouse brain. *J. Lipid Res.* **47(12)**, 2772-2780 (2006).
3. Freire, C.S.R., Silvestre, A.J.D., Neto, C.P., *et al.* Lipophilic extractives of the inner and outer barks of *Eucalyptus globulus*. *Holzforschung* **56(4)**, 372-379 (2002).
4. Jung, K., Reszka, R., Kamlage, B., *et al.* Tissue metabolite profiling identifies differentiating and prognostic biomarkers for prostate carcinoma. *Int. J. Cancer* **133(12)**, 2914-2924 (2013).

#### 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, 12/19/2022

#### 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