

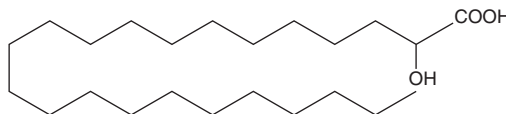
# PRODUCT INFORMATION



## 2-hydroxy Tricosanoic Acid

Item No. 24597

CAS Registry No.: 2718-37-8  
Formal Name: 2-hydroxy-tricosanoic acid  
Synonym: FA 23:0;O  
MF:  $C_{23}H_{46}O_3$   
FW: 370.6  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

2-hydroxy Tricosanoic acid is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy tricosanoic acid in the solvent of choice, which should be purged with an inert gas. 2-hydroxy Tricosanoic acid is soluble in a 5:1 solution of chloroform:methanol.

### Description

2-hydroxy Tricosanoic acid is a hydroxylated fatty acid that has been found in the inner bark of *E. globulus*, the leaves of wild, kiyomi, and zweigeltrepe grape vines, *C. elegans*, and *Pseudosuberites* and *S. massa* sea sponges.<sup>1-4</sup>

### References

- Freire, C.S.R., Silvestre, A.J.D., Neto, C.P., *et al.* Lipophilic extractives of the inner and outer barks of *Eucalyptus globulus*. *Holzforchung* **56(4)**, 372-379 (2002).
- Barnathan, G., Kornprobst, J.-M., Doumenq, P., *et al.* Sponge fatty acids, 5. Characterization of complete series of 2-hydroxy long-chain fatty acids in phospholipids of two Senegalese marine sponges from the family suberitidae: *Pseudosuberites* sp. and *Suberites massa*. *J. Nat. Prod.* **56(12)**, 2104-2113 (2004).
- Kawaguchi, M., Imai, H., Naoe, M., *et al.* Cerebrosides in grapevine leaves: Distinct composition of sphingoid bases among the grapevine species having different tolerances to freezing temperature. *Biosci. Biotechnol. Biochem.* **64(6)**, 1271-1273 (2000).
- Gerdt, S., Lochnit, G., Dennis, R.D., *et al.* Isolation and structural analysis of three neutral glycosphingolipids from a mixed population of *Caenorhabditis elegans* (Nematoda: Rhabditida). *Glycobiology* **7(2)**, 265-275 (1997).

#### 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, 02/23/2024

#### 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