

# PRODUCT INFORMATION

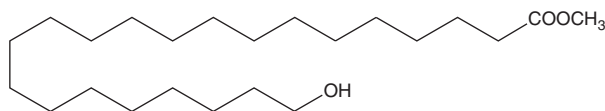


## 22-hydroxy Docosanoic Acid methyl ester

Item No. 24656

**CAS Registry No.:** 38646-51-4  
**Formal Name:** 22-hydroxy-docosanoic acid, methyl ester  
**Synonyms:**  $\omega$ -hydroxy Behenic Acid methyl ester,  
22-hydroxy DCA methyl ester,  
 $\omega$ -hydroxy Docosanoic Acid methyl ester,  
Methyl 22-hydroxydocosanoate, SFE 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

22-hydroxy Docosanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 22-hydroxy docosanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 22-hydroxy Docosanoic acid methyl ester is soluble in organic solvents such as chloroform, warm ethanol, and ethyl ether.

### Description

22-hydroxy Docosanoic acid methyl ester is a hydroxylated fatty acid methyl ester that has been found in *P. reticulata* cork extracts, peat samples from inland Florida, and sediment samples from the Harney River.<sup>1,2</sup>

### References

1. Mota, G.S., Sartori, C.J., Ferreira, J., *et al.* Cellular structure and chemical composition of cork from *Plathymenia reticulata* occurring in the Brazilian Cerrado. *Ind. Crops Prod.* **90**, 65-75 (2016).
2. Jaffé, R., Rushdi, A.I., Medeiros, P.M., *et al.* Natural product biomarkers as indicators of sources and transport of sedimentary organic matter in a subtropical river. *Chemosphere* **64(11)**, 1870-1884 (2006).

#### 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, 04/02/2024

#### 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