

PRODUCT INFORMATION



3-hydroxy Docosanoic Acid

Item No. 22666

CAS Registry No.: 89946-08-7
Formal Name: 3-hydroxy-docosanoic acid
Synonyms: 3-hydroxy Behenic Acid,
3-hydroxy DCA, β -hydroxy DCA,
 β -hydroxy Docosanoic Acid

MF: C₂₂H₄₄O₃

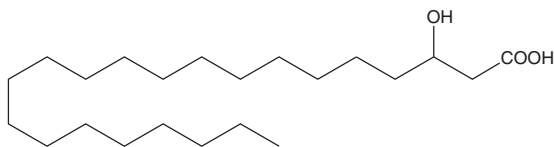
FW: 356.6

Purity: $\geq 95\%$

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥ 2 years



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

Laboratory Procedures

3-hydroxy Docosanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-hydroxy docosanoic acid in the solvent of choice. 3-hydroxy Docosanoic acid is soluble in the organic solvent chloroform, which should be purged with an inert gas, at a concentration of approximately 0.3 mg/ml.

Description

3-hydroxy Docosanoic acid is a hydroxylated form of the 22-carbon saturated docosanoic acid (Item No. 9000338). It is found in lipopolysaccharides from *C. trachomatis* and *C. psittaci* and in *S. synnaedendra*.^{1,2} 3-hydroxy Docosanoic acid, in the form of an acyl-CoA metabolite, is an intermediate in fatty acid chain elongation from arachidic acid (C20:0; Item No. 9000339) to docosanoic acid (C22:0).³

References

1. Heine, H., Müller-Loennies, S., Brade, L., et al. Endotoxic activity and chemical structure of lipopolysaccharides from *Chlamydia trachomatis* serotypes E and L2 and *Chlamydomphila psittaci* 6BC. *Eur. J. Biochem.* **270**(3), 440-450 (2003).
2. Sebolai, O.M., Kock, J.L., Pohl, C.H., et al. Report on the discovery of a novel 3-hydroxy oxylipin cascade in the yeast *Saccharomycopsis synnaedendra*. *Prostaglandins Other Lipid Mediat.* **74**(1-4), 139-146 (2004).
3. Guillou, H., Zdravec, D., Martin, P.G., et al. The key roles of elongases and desaturases in mammalian fatty acid metabolism: Insights from transgenic mice. *Prog. Lipid Res.* **49**(2), 186-199 (2010).

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, 05/01/2018

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