

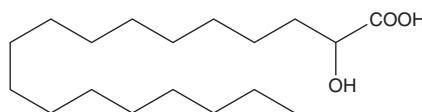
PRODUCT INFORMATION



2-hydroxy Stearic Acid

Item No. 24590

CAS Registry No.: 629-22-1
Formal Name: 2-hydroxy-octadecanoic acid
Synonyms: NSC 907, (±)- α -hydroxy Stearic Acid, (±)- α -hydroxy Stearic Acid
MF: C₁₈H₃₆O₃
FW: 300.5
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 Stearic acid is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxy stearic acid in the solvent of choice. 2-hydroxy Stearic acid is soluble in a 5:1 solution of chloroform:methanol.

Description

2-hydroxy Stearic acid is a naturally occurring hydroxylated fatty acid that has been found in humans and *U. armoricana*.¹⁻³ It reduces growth of Ehrlich ascites tumor (EAT) cells *in vitro* when used at a concentration of 100 μ M.⁴ Levels of 2-hydroxy stearic acid are elevated in human joint synovial fluid after ankle fracture and in patients with Buruli ulcer disease.^{2,3}

References

1. Kendel, M., Wielgosz-Collin, G., Bertrand, S., *et al.* Lipid composition, fatty acids and sterols in the seaweeds *Ulva armoricana*, and *Solieria chordalis* from Brittany (France): An analysis from nutritional, chemotaxonomic, and antiproliferative activity perspectives. *Mar. Drugs* **13(9)**, 5606-5628 (2015).
2. Leimer, E.M., Pappan, K.L., Nettles, D.L., *et al.* Lipid profile of human synovial fluid following intra-articular ankle fracture. *J. Orthop. Res.* **35(3)**, 657-666 (2017).
3. Niang, F., Sarfo, F.S., Frimpong, M., *et al.* Metabolomic profiles delineate mycolactone signature in Buruli ulcer disease. *Sci. Rep.* **5:17693**, (2015).
4. Kusmoto, K., Kageyama, K., Matsuda, T., *et al.* Effects of ethyl-esterization, chain-lengths, unsaturation degrees, and hyperthermia on carcinostatic effect of omega-hydroxylated fatty acids. *Exp. Oncol.* **29(2)**, 106-110 (2007).

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