

PRODUCT INFORMATION



Octacosanoic Acid

Item No. 31735

CAS Registry No.: 506-48-9

Synonyms: C28:0, FA 28:0,
Montanic Acid,
NSC 407311

MF: $C_{28}H_{56}O_2$

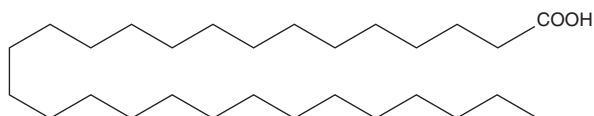
FW: 424.7

Purity: $\geq 95\%$

Supplied as: A crystalline 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

Octacosanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the octacosanoic acid in the solvent of choice, which should be purged with an inert gas. Octacosanoic acid is slightly soluble in ethanol and chloroform.

Description

Octacosanoic acid is a very long-chain saturated fatty acid. It is the major component of D-003, a mixture of very long-chain aliphatic acids purified from sugar cane wax that has antiplatelet and cholesterol-lowering activities in animal models.¹

Reference

1. Gámez, R., Rodeiro, I., Fernández, I., *et al.* Preliminary evaluation of the cytotoxic and genotoxic potential of D-003: Mixture of very long chain fatty acids. *Teratog. Carcinoq. Mutagen.* **22**(3), 175-181 (2002).

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, 02/20/2024

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