

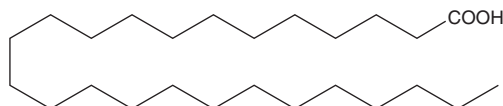
PRODUCT INFORMATION



Pentacosanoic Acid

Item No. 15197

CAS Registry No.: 506-38-7
Formal Name: pentacosanoic acid
Synonyms: FA 25:0, Hyenic Acid, NSC 89289
MF: C₂₅H₅₀O₂
FW: 382.7
Purity: ≥98%
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

Pentacosanoic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the pentacosanoic acid in the solvent of choice, which should be purged with an inert gas. Pentacosanoic acid is soluble in the organic solvent chloroform at a concentration of approximately 2.5 mg/ml.

Description

Pentacosanoic acid is a 25-carbon, fully saturated long-chain fatty acid found in plant lipids.

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, 02/07/2024

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