

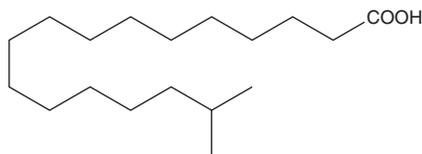
PRODUCT INFORMATION



Isostearic Acid

Item No. 43258

CAS Registry No.: 2724-58-5
Formal Name: 16-methyl-heptadecanoic acid
Synonym: Isooctadecanoic Acid
MF: C₁₈H₃₆O₂
FW: 284.5
Purity: ≥95% (mixture of isosterics)
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Isostearic acid is supplied as a neat oil. A stock solution may be made by dissolving the isostearic acid in the solvent of choice, which should be purged with an inert gas. Isostearic acid is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

Isostearic acid is a branched-chain fatty acid. Formulations containing isostearic acid have been used in skin creams and cosmetics as emulsifiers and texturing agents.

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, 08/13/2025

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