

# PRODUCT INFORMATION



## 27-hydroxy Heptacosanoic Acid methyl ester

Item No. 24657

CAS Registry No.: 369635-50-7

Formal Name: 27-hydroxy-heptacosanoic acid,  
methyl ester

Synonyms: Methyl 27-hydroxyheptacosanoate,  
SFE 28:0;O

MF:  $C_{28}H_{56}O_3$

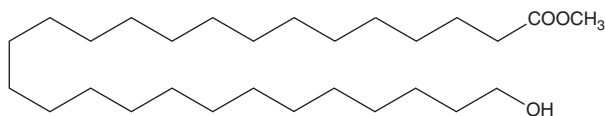
FW: 440.7

Purity:  $\geq 97\%$

Supplied as: A solid

Storage: -20

Stability:  $\geq 4$  years



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

### Laboratory Procedures

27-hydroxy Heptacosanoic acid methyl ester is supplied as a solid. A stock solution may be made by dissolving the 27-hydroxy heptacosanoic acid methyl ester in the solvent of choice, which should be purged with an inert gas. 27-hydroxy Heptacosanoic acid methyl ester is soluble in the organic solvent chloroform.

### Description

27-hydroxy Heptacosanoic acid methyl ester is a hydroxylated fatty acid methyl ester form of heptacosanoic acid.

#### 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, 04/05/2024

#### 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