

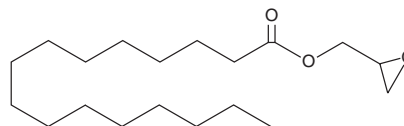
PRODUCT INFORMATION



Glycidyl Palmitate

Item No. 9003681

CAS Registry No.: 7501-44-2
Formal Name: hexadecanoic acid, 2-oxiranylmethyl ester
Synonyms: 18:0 GE, 18:0 Glycidyl ester, GE(18:0), NSC 406558, Palmitic Acid Glycidyl ester
MF: $C_{19}H_{36}O_3$
FW: 312.5
Purity: $\geq 98\%$
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

Glycidyl palmitate is supplied as a solid. A stock solution may be made by dissolving the glycidyl palmitate in the solvent of choice, which should be purged with an inert gas. Glycidyl palmitate is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

Glycidyl palmitate is a glycidyl fatty acid ester.¹ It has been found in palm and corn oils.

Reference

1. Blumhorst, M.R., Collison, M.W., Cantrill, R., et al. Collaborative study for the analysis of glycidyl fatty acid esters in edible oils using LC-MS. *J. Am. Oil Chem. Soc.* **90**(4), 493-500 (2013).

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, 11/10/2022

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