

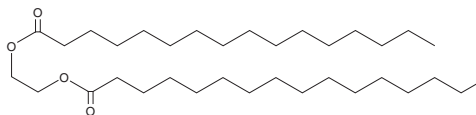
PRODUCT INFORMATION



1,2-Dipalmitoyl Ethylene Glycol

Item No. 25590

CAS Registry No.: 624-03-3
Formal Name: hexadecanoic acid,
1,1'-(1,2-ethanediyl) ester
Synonyms: 16:0 Ethylene Glycol, Ethylene
Glycol Dihexadecanoate, Ethylene
Glycol Dipalmitate, NSC 6819
MF: $C_{34}H_{66}O_4$
FW: 538.9
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

1,2-Dipalmitoyl ethylene glycol is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2-dipalmitoyl ethylene glycol in the solvent of choice, which should be purged with an inert gas. 1,2-Dipalmitoyl ethylene glycol is soluble in the organic solvent chloroform at a concentration of approximately 30 mg/ml.

Description

1,2-Dipalmitoyl ethylene glycol is a fatty acid ester. It has been used in the manufacture of plastics, moldable flame-retardant resin compositions, and as a component of antifouling paint.¹⁻³ 1,2-Dipalmitoyl ethylene glycol can migrate out of plastics but is not toxic to rats during prenatal development when administered at a dose of 900 mg/kg per day.³ Formulations containing 1,2-dipalmitoyl ethylene glycol have been used in a variety of cosmetic and personal care products.

References

1. Hanabusa, K. Flame-retardant polybutylene terephthalate resin compositions and mold parts formed therefrom. *Polyplastics Co., Ltd.* **US005258434A** (1993).
2. Hamade, R., Yamamori, N., and Okahata, Y. Glucoside derivatives for enzyme modification, lipid-coated enzymes, method of producing such enzymes and antifouling paint composition. *Nippon Paint Co., Ltd.* **US5770188A** (1998).
3. European Food Safety Authority. Scientific opinion on the safety assessment of the substance ethylene glycol dipalmitate, CAS No 624-03-3, for use in food contact materials. *EFSA J.* **13(2)**, 4019 (2015).

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, 11/28/2022

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