

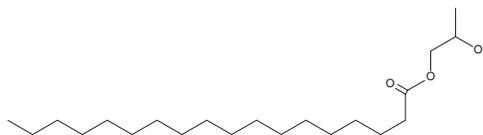
PRODUCT INFORMATION



2-Hydroxypropyl Octadecanoate

Item No. 41151

CAS Registry No.: 1323-39-3
Formal Name: octadecanoic acid, monoester with 1,2-propanediol
Synonyms: NSC 4841, 2-Hydroxypropyl Stearate, Propyleneglycol Monostearate
MF: C₂₁H₄₂O₃
FW: 342.6
Purity: ≥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

2-Hydroxypropyl octadecanoate is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxypropyl octadecanoate in the solvent of choice, which should be purged with an inert gas. 2-Hydroxypropyl octadecanoate is soluble in ethanol, ethyl ether, hexane, and chloroform.

Description

2-Hydroxypropyl octadecanoate is a lipid ester composed of a hydroxylated propylene glycol connected to the long-chain saturated fatty acid stearic acid (Item No. 10011298).¹ Formulations containing 2-hydroxypropyl octadecanoate have been used as emulsifiers in food-grade preparations and as topical pharmaceutical excipients.^{1,2}

References

1. Macdonald, I.A. Emulsifiers: Processing and quality control. *J. Am. Oil Chem. Soc.* **45**, 584A-586A (1968).
2. Searles, G.E., and DesGroseilliers, J.-P. Excipients in topical corticosteroid preparations in Canada. *CMAJ* **141**(5), 396-405 (1989).

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, 06/25/2024

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