

PRODUCT INFORMATION



2-Hydroxypropyl Hexadecanoate

Item No. 41150

CAS Registry No.: 29013-28-3
Formal Name: hexadecanoic acid, monoester with 1,2-propanediol
Synonyms: 2-Hydroxypropyl Palmitate, Propyleneglycol Monopalmitate
MF: C₁₉H₃₈O₃
FW: 314.5
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 hexadecanoate is supplied as a solid. A stock solution may be made by dissolving the 2-hydroxypropyl hexadecanoate in the solvent of choice, which should be purged with an inert gas. 2-Hydroxypropyl hexadecanoate is soluble in ethanol, ethyl ether, chloroform, and hexane.

Description

2-Hydroxypropyl hexadecanoate is a medium-chain fatty acid monoester.¹ Formulations containing 2-hydroxypropyl hexadecanoate have been used as emulsifying agents in food preparations.²

References

1. Bunchongprasert, K. and Shao, J. Effect of fatty acid ester structure on cytotoxicity of self-emulsified nanoemulsion and transport of nanoemulsion droplets. *Colloids Surf. B Biointerfaces* **194**, 111220 (2020).
2. Uematsu, Y., Hirokado, M., Hirata, K., *et al.* Determination of monoglycerides and propylene glycol esters of fatty acids in food additive preparations by GC without derivatization. *Shokuhin Eiseigaku Zasshi / Journal of the Food Hygienic Society of Japan* **38(4)**, 248-258 (1997).

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, 06/26/2024

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