

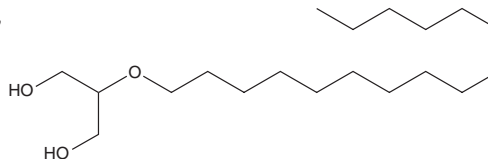
PRODUCT INFORMATION



2-O-Hexadecyl Glycerol

Item No. 11924

CAS Registry No.: 1931-78-8
Formal Name: 2-(hexadecyloxy)-1,3-propanediol
Synonyms: 16 OGE, β -Chimyl Alcohol, 2-Chimyl Alcohol, 2-Palmityl Glycerol
MF: C₁₉H₄₀O₃
FW: 316.5
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

2-O-Hexadecyl glycerol is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-O-hexadecyl glycerol in the solvent of choice, which should be purged with an inert gas. 2-O-Hexadecyl glycerol is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of 2-O-hexadecyl glycerol in these solvents is approximately 15 mg/ml.

Description

2-O-Hexadecyl glycerol is a β -monoglyceride of palmitic acid (Item No. 10006627). It is an intermediate in the metabolism of triacylglycerols or diacylglycerols and an analog of 2-palmitoyl glycerol (Item No. 17882), an endocannabinoid that fails to bind to cannabinoid CB₁ or CB₂ receptors, but rather potentiates the biological effects 2-arachidonoyl glycerol (Item No. 62160).¹

Reference

1. Ben-Shabat, S., Frider, E., Sheskin, T., *et al.* An entourage effect: Inactive endogenous fatty acid glycerol esters enhance 2-arachidonoyl-glycerol cannabinoid activity. *Eur. J. Pharmacol.* **353**, 23-31 (1998).

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, 11/17/2022

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