

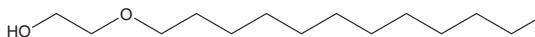
PRODUCT INFORMATION



Ethylene Glycol monododecyl ether

Item No. 36857

CAS Registry No.: 4536-30-5
Formal Name: 2-(dodecyloxy)-ethanol
Synonyms: C₁₂E₁, EGDE
MF: C₁₄H₃₀O₂
FW: 230.4
Purity: ≥98%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Ethylene glycol monododecyl ether is supplied as a neat oil. A stock solution may be made by dissolving the ethylene glycol monododecyl ether in the solvent of choice, which should be purged with an inert gas. Ethylene glycol monododecyl ether is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ethylene glycol monododecyl ether in ethanol is approximately 15 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Ethylene glycol monododecyl ether is a nonionic surfactant with a critical micelle concentration (CMC) of 27.5 μM at 25°C.¹ It enhances the transdermal permeation of the antiemetic ondansetron across isolated porcine skin when used at a concentration of 160 mM, however, it decreases the viability of human epidermal keratinocytes and human dermal fibroblasts when used at a concentration of 160 μM.²

References

1. Islam, M.N. and Kato, T. Temperature dependence of the surface phase behavior and micelle formation of some nonionic surfactants. *J. Phys. Chem. B* **107**(4), 965-971 (2003).
2. Silva, S.M.C., Hu, L., Sousa, J.J.S., *et al.* A combination of nonionic surfactants and iontophoresis to enhance the transdermal drug delivery of ondansetron HCl and diltiazem HCl. *Eur. J. Pharm. Biopharm.* **80**(3), 663-673 (2012).

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, 12/21/2022

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