

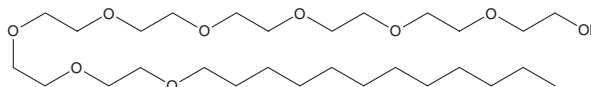
# PRODUCT INFORMATION



## Octaethylene Glycol monododecyl ether

Item No. 20938

CAS Registry No.: 3055-98-9  
Formal Name: 3,6,9,12,15,18,21,24-octaohexatriacontan-1-ol  
Synonym: C<sub>12</sub>E<sub>8</sub>  
MF: C<sub>28</sub>H<sub>58</sub>O<sub>9</sub>  
FW: 538.8  
Purity: ≥95%  
Supplied as: A low-melting 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

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

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of octaethylene glycol monododecyl ether can be prepared by directly dissolving the low-melting solid in aqueous buffers. The solubility of octaethylene glycol monododecyl ether in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Octaethylene glycol monododecyl ether is a nonionic surfactant formed by the ethoxylation of dodecanol, yielding a compound with eight repeated units of ethylene glycol. It can be used for solubilization of membrane-bound proteins.<sup>1</sup>

### Reference

1. Fogeron, M.L., Badillo, A., Jirasko, V., *et al.* Wheat germ cell-free expression: Two detergents with a low critical micelle concentration allow for production of soluble HCV membrane proteins. *Protein Expr. Purif.* **105**, 39-46 (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, 01/30/2025

#### 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