

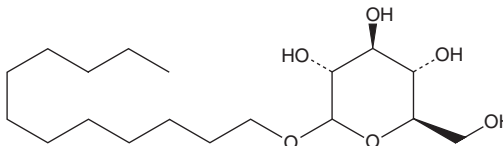
# PRODUCT INFORMATION



## Dodecyl-D-Glucopyranoside

Item No. 42721

CAS Registry No.: 27836-64-2  
Synonyms: Lauryl Glucoside, LG  
MF:  $C_{18}H_{36}O_6$   
FW: 348.5  
Purity:  $\geq 98\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Dodecyl-D-glucopyranoside is supplied as a solid. A stock solution may be made by dissolving the dodecyl-D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. Dodecyl-D-glucopyranoside is soluble ( $\geq 10$  mg/ml) in ethanol and DMSO.

### Description

Dodecyl-D-glucopyranoside is a nonionic detergent.<sup>1</sup> It has a critical micelle concentration (CMC) of 125  $\mu\text{M}$ . Formulations containing dodecyl-D-glucopyranoside have been used in personal care products, cosmetics, and laundry detergents.

### Reference

1. Qi, Y., Zhang, Z., Wu, Z., *et al.* Composite system of lauryl glucoside and lauryl glycoside sulfosuccinate: properties and applications in vitamin E nanoemulsions. *Chempluschem* **89**(8), e202400058 (2024).

#### 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, 03/24/2025

#### 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