

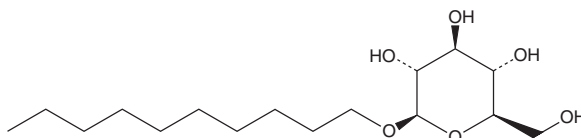
PRODUCT INFORMATION



n-Decyl-β-D-Glucopyranoside

Item No. 42722

CAS Registry No.: 58846-77-8
Formal Name: decyl, β-D-glucopyranoside
Synonyms: C10βGlc, Decyl Glucoside
MF: C₁₆H₃₂O₆
FW: 320.4
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

n-Decyl-β-D-glucopyranoside is supplied as a solid. A stock solution may be made by dissolving the n-decyl-β-D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. n-Decyl-β-D-glucopyranoside is soluble (≥10 mg/ml) in ethanol and 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 n-decyl-β-D-glucopyranoside can be prepared by directly dissolving the solid in aqueous buffers. n-Decyl-β-D-glucopyranoside is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

n-Decyl-β-D-glucopyranoside is a nonionic alkyl glucoside chiral surfactant with a critical micelle concentration (CMC) value of 0.8 mM.¹ It has been used to solubilize membrane proteins.² Formulations containing n-decyl-β-D-glucopyranoside are associated with allergic contact dermatitis.

References

1. Matsumura, S., Imai, K., Yoshikawa, S., *et al.* Surface activities, biodegradability and antimicrobial properties of n-alkyl glucosides, mannosides and galactosides. *J. Am. Oil Chem. Soc.* **67**, 996-1001 (1990).
2. Stubbs, G.W., Smith, H.G., Jr., and Litman, B.J. Alkyl glucosides as effective solubilizing agents for bovine rhodopsin. A comparison with several commonly used detergents. *Biochim. Biophys. Acta* **426(1)**, 46-56 (1976).

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, 02/24/2025

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