

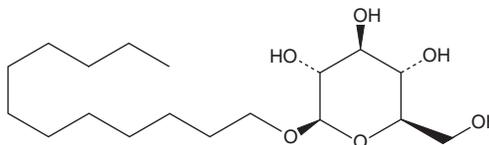
PRODUCT INFORMATION



n-Dodecyl-β-D-Glucopyranoside

Item No. 36030

CAS Registry No.: 59122-55-3
Synonyms: BDDG, Dodecyl β-D-glucoside, Dodecyl glucoside
MF: C₁₈H₃₆O₆
FW: 348.5
Purity: ≥95%
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-Dodecyl-β-D-glucopyranoside is supplied as a solid. A stock solution may be made by dissolving the n-dodecyl-β-D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. n-Dodecyl-β-D-glucopyranoside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of n-dodecyl-β-D-glucopyranoside in DMF is approximately 10 mg/ml and approximately 5 mg/ml in ethanol and DMSO.

Description

n-Dodecyl-β-D-glucopyranoside is a nonionic detergent.¹ It has a critical micelle concentration (CMC) of 190 μM.¹ It increases chloramphenicol acetyltransferase (CAT) activity when used at a concentration of 204 μM. n-Dodecyl-β-D-glucopyranoside inhibits the uptake of β-methyl glucoside in hamster small intestine *in vitro* (K_i = 200 μM).²

References

1. Lu, J. and Jiang, C. Inhibition kinetics of chloramphenicol acetyltransferase by selected detergents. *Biochem. Biophys. Res. Commun.* **196(1)**, 12-17 (1993).
2. Ramaswamy, K., Bhattacharyya, B.R., and Crane, R.K. Studies on the transport of aliphatic glucosides by hamster small intestine *in vitro*. *Biochim. Biophys. Acta* **433(1)**, 32-38 (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, 12/19/2022

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