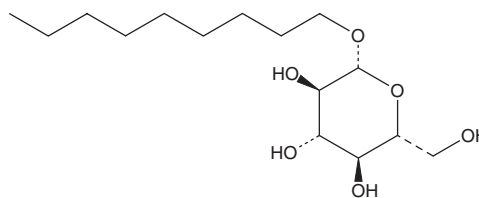


PRODUCT INFORMATION



n-Nonyl-β-D-glucopyranoside Item No. 25705

CAS Registry No.: 69984-73-2
Formal Name: nonyl, β-D-glucopyranoside
MF: C₁₅H₃₀O₆
FW: 306.4
Purity: ≥95%
Supplied as: A crystalline 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-Nonyl-β-D-glucopyranoside is supplied as a crystalline solid. A stock solution may be made by dissolving the n-nonyl-β-D-glucopyranoside in the solvent of choice. n-Nonyl-β-D-glucopyranoside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of n-nonyl-β-D-glucopyranoside in ethanol is approximately 20 mg/ml and approximately 15 mg/ml in DMSO and DMF.

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-nonyl-β-D-glucopyranoside can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of n-nonyl-β-D-glucopyranoside in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

n-Nonyl-β-D-glucopyranoside is an anionic alkylglucoside chiral surfactant that is commonly used for the solubilization and crystallization of biological membrane proteins.¹ It has also been used in the separation of drug enantiomers by micellar electrokinetic chromatography.²

References

1. Zhang, R., Marone, P.A., Thiyagarajan, P., *et al.* Structure and molecular fluctuations of n-alkyl-β-D-glucopyranoside micelles determined by x-ray and neutron scattering. *Langmuir* **15(22)**, 7510-7519 (1999).
2. Otsuka, K. and Terabe, S. Enantiomer separation of drugs by micellar electrokinetic chromatography using chiral surfactants. *J. Chromatogr. A.* **875(1-2)**, 163-178 (2000).

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, 12/07/2022

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