

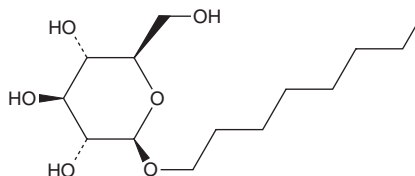
PRODUCT INFORMATION



n-Octylglucoside

Item No. 14327

CAS Registry No.: 29836-26-8
Formal Name: octyl-β-D-glucopyranoside
Synonym: BTB 11967
MF: C₁₄H₂₈O₆
FW: 292.4
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-Octylglucoside is supplied as a crystalline solid. A stock solution may be made by dissolving the n-Octylglucoside in the solvent of choice, which should be purged with an inert gas. n-Octylglucoside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of n-octylglucoside in ethanol is approximately 20 mg/ml and approximately 16 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-octylglucoside can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of n-octylglucoside 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-Octylglucoside is a mild, non-denaturing detergent that is used for the solubilization and reconstitution of membrane-bound proteins. The high critical micelle concentration (0.7%) of n-octylglucoside facilitates ready removal from final protein extracts by dialysis or gel filtration.^{1,2} n-Octylglucoside can be used in 2D electrophoresis and to improve selectivity of immunoprecipitation of phosphotyrosine modified proteins.^{1,3}

References

1. Zhang, G. and Neubert, T.A. Use of detergents to increase selectivity of immunoprecipitation of tyrosine phosphorylated peptides prior to identification by MALDI quadrupole-TOF MS. *Proteomics* **6**(2), 571-578 (2006).
2. Petri, W.A., Jr. and Wagner, R.R. Reconstitution into liposomes of the glycoprotein of vesicular stomatitis virus by detergent dialysis. *J. Biol. Chem.* **254**(11), 4313-4316 (1979).
3. Lambot, M., Frétier, S., Op De Beeck, A., et al. Reconstitution of hepatitis C virus envelope glycoproteins into liposomes as a surrogate model to study virus attachment. *J. Biol. Chem.* **277**(23), 20625-20630 (2002).

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