

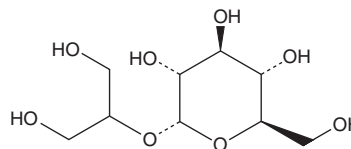
PRODUCT INFORMATION



2-O-(α -D-Glucopyranosyl)glycerol

Item No. 34492

CAS Registry No.: 22160-26-5
Formal Name: α -D-glucopyranoside, 2-hydroxy-1-(hydroxymethyl)ethyl
Synonyms: Glucosylglycerol, Glycooin
MF: $C_9H_{18}O_8$
FW: 254.2
Purity: $\geq 95\%$
Supplied as: A solution in water
Storage: -20°C
Stability: ≥ 2 years
Item Origin: Natural sucrose and glycerine



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

Laboratory Procedures

2-O-(α -D-Glucopyranosyl)glycerol is supplied as a solution in water. To change the solvent, simply evaporate the water under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 2-O-(α -D-glucopyranosyl)glycerol in these solvents is approximately 30 mg/ml.

Description

2-O-(α -D-Glucopyranosyl)glycerol is a compatible solute that has been found in various bacteria.^{1,2} It prevents drying-induced membrane fusion and carboxyfluorescein (CF) leakage in liposomes.² 2-O-(α -D-Glucopyranosyl)glycerol levels are increased in response to osmotic stress in bacteria.¹

References

1. Pocard, J.A., Smith, L.T., Smith, G.M., *et al.* A prominent role for glucosylglycerol in the adaptation of *Pseudomonas mendocina* SKB70 to osmotic stress. *J. Bacteriol.* **176(22)**, 6877-6884 (1994).
2. Hinch, D.K. and Hagemann, M. Stabilization of model membranes during drying by compatible solutes involved in the stress tolerance of plants and microorganisms. *Biochem. J.* **383(Pt 2)**, 277-283 (2004).

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, 11/23/2022

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