

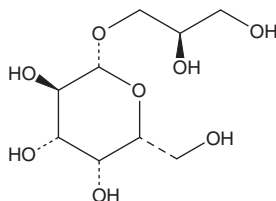
PRODUCT INFORMATION



(2R)-Glycerol-O-β-D-galactopyranoside

Item No. 17218

CAS Registry No.: 16232-91-0
Formal Name: 2R,3-dihydroxypropyl-β-D-galactopyranoside
MF: C₉H₁₈O₈
FW: 254.2
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Description

(2R)-Glycerol-O-β-D-galactopyranoside is a substrate for β-galactosidase, the lactose repressor, the galactose-binding protein, and the β-methylgalactoside transport system.^{1,2} It also serves as a substrate for the third *lac* operon encoded enzyme, thiogalactoside transacetylase.^{1,3}

References

1. Silhavy, T.J. and Boos, W. Selection procedure for mutants defective in the β-methylgalactoside transport system of *Escherichia coli* utilizing the compound 2R-glyceryl-β-D-galactopyranoside. *J. Bacteriol.* **120**(1), 424-432 (1974).
2. Silhavy, T.J. and Boos, W. A convenient synthesis of (2R)-glyceryl-β-D-galactopyranoside. A substrate for β-galactosidase, the lactose repressor, the galactose-binding protein, and the β-methylgalactoside transport system. *J. Biol. Chem.* **248**(18), 6571-6574 (1973).
3. Boos, W. and Wallenfels, K. Investigation on the induction of the Lac-enzyme. 2. The permeation of galactosylglycerol in *Escherichia coli*. *Eur. J. Biochem.* **3**(3), 360-363 (1968).

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, 11/28/2022

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