

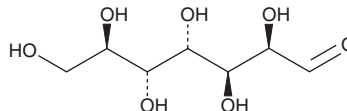
PRODUCT INFORMATION



D-glycero-D-galacto-Heptose

Item No. 46208

CAS Registry No.: 5328-64-3
Formal Name: 2R,3S,4S,5R,6R,7-hexahydroxyheptanal
Synonym: NSC 1977
MF: $C_7H_{14}O_7$
FW: 210.2
Purity: $\geq 98\%$
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

D-glycero-D-galacto-Heptose is supplied as a solid. A stock solution may be made by dissolving the D-glycero-D-galacto-heptose in the solvent of choice, which should be purged with an inert gas. D-glycero-D-galacto-Heptose is soluble (≥ 10 mg/ml) in DMSO.

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 D-glycero-D-galacto-heptose can be prepared by directly dissolving the solid in aqueous buffers. D-glycero-D-galacto-Heptose is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

D-glycero-D-galacto-Heptose is a seven-carbon monosaccharide and the antigenic component of a polysaccharide from *E. saburreum*, a bacteria found in the human oral cavity.¹ It has also been found in avocado and cactus fruits.^{2,3}

References

1. Nakazawa, F., Ito, T., Sato, N., et al. Chemical composition and immunological characterization of polysaccharide antigen from *Eubacterium saburreum* T18. *Infect. Immun.* **55**(4), 871-876 (1987).
2. Sephton, H.H. and Richtmyer, N.K. The isolation of a second octulose and of a heptose from the avocado: D-Glycero-L-galacto-octulose and D-glycero-D-galacto-heptose. *J. Org. Chem.* **28**(6), 1691-1694 (1963).
3. Alsaad, A.J.A. and Sameer, L.M. Study of the antioxidant activity of cactus (*Opuntia delillenii*) fruits (pulp and peels) and characterisation of their bioactive compounds by GC-MS. *Basrah J. Agric. Sci.* **34**(2), 204-219 (2021).

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, 08/03/2026

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