

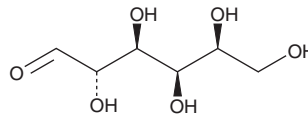
PRODUCT INFORMATION



L-Gulose

Item No. 34049

CAS Registry No.: 6027-89-0
MF: C₆H₁₂O₆
FW: 180.2
Purity: ≥95%
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

L-Gulose is supplied as a solid. A stock solution may be made by dissolving the L-gulose in the solvent of choice, which should be purged with an inert gas. L-Gulose is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of L-gulose in these solvents is approximately 10 and 5 mg/ml, respectively.

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 L-gulose can be prepared by directly dissolving the solid in aqueous buffers. The solubility of L-gulose in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Gulose is a carbohydrate starting material.^{1,2} It has been used as a starting material in the synthesis of L-nucleoside-based anti-HIV agents.

References

1. Woodyer, R.D., Christ, T.N., and Deweese, K.A. Single-step bioconversion for the preparation of L-gulose and L-galactose. *Carbohydr. Res.* **345(3)**, 363-368 (2010).
2. Jeong, L.S., Schinazi, R.F., Beach, J.W., *et al.* Asymmetric synthesis and biological evaluation of β-L-(2R,5S)- and α-L-(2R,5R)-1,3-oxathiolane-pyrimidine and -purine nucleosides as potential anti-HIV agents. *J. Med. Chem.* **36(2)**, 181-195 (1993).

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, 11/17/2022

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