

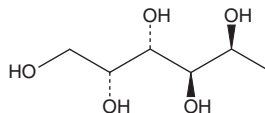
PRODUCT INFORMATION



L-Fucitol

Item No. 26540

CAS Registry No.: 13074-06-1
Formal Name: 1-deoxy-D-galactitol
Synonym: NSC 1957
MF: C₆H₁₄O₅
FW: 166.2
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

L-Fucitol is supplied as a crystalline solid. Aqueous solutions of L-fucitol can be prepared by directly dissolving the crystalline solid in aqueous buffers. L-Fucitol is slightly soluble in PBS, pH 7.2. It is also soluble in water. The solubility of L-fucitol in water is approximately 30 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-Fucitol is a reduced form of L(-)-fucose (Item No. 16479). It has been used to determine the structure of *E. coli* and *B. pallidus* L-fucose isomerase.^{1,2}

References

1. Seemann, J.E. and Schulz, G.E. Structure and mechanism of L-fucose isomerase from *Escherichia coli*. *J. Mol. Biol.* **273**(1), 256-268 (1997).
2. Takeda, K., Yoshida, H., Izumori, K., *et al.* X-ray structures of *Bacillus pallidus* D-arabinose isomerase and its complex with L-fucitol. *Biochim. Biophys. Acta.* **1804**(6), 1359-1368 (2010).

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, 10/24/2022

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