

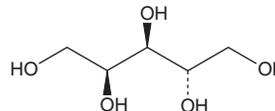
PRODUCT INFORMATION



L-Arabinitol

Item No. 38882

CAS Registry No.: 7643-75-6
Synonyms: L-arabino-Pentitol, (-)-Arabitol, L-Arabitol, L-(-)-Arabitol, L-Lyxitol, NSC 83253
MF: C₅H₁₂O₅
FW: 152.1
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-Arabinitol is supplied as a solid. A stock solution may be made by dissolving the L-arabinitol in water. We do not recommend storing the aqueous solution for more than one day.

Description

L-Arabinitol is an endogenous metabolite.^{1,2} The ratio of D-arabinitol (Item No. 33826) to L-arabinitol in urine has been used as a marker of *Candida* infections.

References

1. Stradomska, T.J., Bobula-Milewska, B., Bauer, A., *et al.* Urinary D-arabinitol/L-arabinitol levels in infants undergoing long-term antibiotic therapy. *J. Clin. Microbiol.* **43(10)**, 5351-5354 (2005).
2. Salonen, J.H., Rimpiläinen, M., Lehtonen, L., *et al.* Measurement of the D-arabinitol/L-arabinitol ratio in urine of neutropenic patients treated empirically with amphotericin B. *Eur. J. Clin. Microbiol. Infect. Dis.* **20(3)**, 179-184 (2001).

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, 11/06/2023

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