

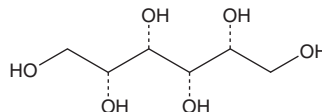
PRODUCT INFORMATION



D-Iditol

Item No. 35597

CAS Registry No.: 25878-23-3
Synonym: NSC 227898
MF: $C_6H_{14}O_6$
FW: 182.2
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

D-Iditol is supplied as a solid. A stock solution may be made by dissolving the D-Iditol in the solvent of choice, which should be purged with an inert gas. D-Iditol is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of D-Iditol in these solvents is approximately 2 and 1 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 D-Iditol can be prepared by directly dissolving the D-Iditol in aqueous buffers. The solubility of D-Iditol in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

D-Iditol is a polyol.¹ It inhibits glucosidase I, but not glucosidase II, when used at a concentration of ~ 1 mM.²

References

1. Sasahara, H. and Izumori, K. Production of D-iditol from D-sorbose by *Rhodotolura rubra* RY10 isolated from miso paste. *J. Biosci. Bioeng.* **87**(4), 548-550 (1999).
2. Muniruzzaman, S., Pan, Y.T., Zeng, Y., *et al.* Inhibition of glycoprotein processing by L-fructose and L-xylulose. *Glycobiology* **6**(8), 795-803 (1996).

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, 11/09/2022

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