

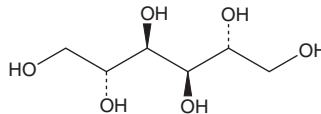
PRODUCT INFORMATION



D-Mannitol

Item No. 23372

CAS Registry No.: 69-65-8
MF: $C_6H_{14}O_6$
FW: 182.2
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 5 years



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

Description

D-Mannitol (Item No. 23372) is an analytical reference standard categorized as an alcohol sugar that is used as an adulterant.¹ It is commonly used as a cutting agent in heroin and cocaine. This product is intended for research and forensic applications.

Reference

1. Broséus, J., Gentile, N., and Esseiva, P. The cutting of cocaine and heroin: A critical review. *Forensic Sci. Int.* **262**, 73-83 (2016).

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, 03/01/2024

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