

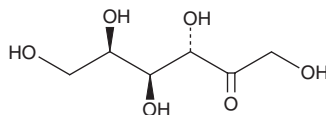
PRODUCT INFORMATION



D-(-)-Tagatose

Item No. 39233

CAS Registry No.: 87-81-0
Formal Name: D-tagatose
MF: C₆H₁₂O₆
FW: 180.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

D-(-)-Tagatose is supplied as a solid. A stock solution may be made by dissolving the D-(-)-tagatose in water. We do not recommend storing the aqueous solution for more than one day.

Description

D-(-)-Tagatose is a natural non-caloric sweetener and hexose monosaccharide that has been found in dairy products.¹ It is metabolized by gut bacteria in the colon of pigs.² Formulations containing D-(-)-tagatose have been used as a sweetening agents in foods and beverages.

References

1. Livesey, G. and Brown, J.C. D-tagatose is a bulk sweetener with zero energy determined in rats. *J. Nutr.* **126(6)**, 1601-1609 (1996).
2. Laerke, H.N. and Jensen, B.B. D-tagatose has low small intestinal digestibility but high large intestinal fermentability in pigs. *J. Nutr.* **129(5)**, 1002-1009 (1999).

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, 01/22/2024

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