

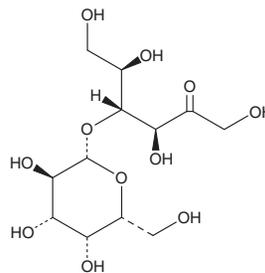
PRODUCT INFORMATION



Lactulose

Item No. 23795

CAS Registry No.: 4618-18-2
Formal Name: 4-O-β-D-galactopyranosyl-D-fructose
MF: C₁₂H₂₂O₁₁
FW: 342.3
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

Lactulose is supplied as a solid. A stock solution may be made by dissolving the lactulose in water. We do not recommend storing the aqueous solution for more than one day.

Description

Lactulose is a synthetic and non-digestible sugar.¹ It reduces colonic pH to 3.5 and stimulates circular muscle activity and propulsion of fluid in isolated guinea pig ileum and colon. Formulations containing lactulose have been used to treat constipation.

Reference

1. Bennett, A. and Eley, K.G. Intestinal pH and propulsion: An explanation of diarrhoea in lactase deficiency and laxation by lactulose. *J. Pharm. Pharmacol.* **28(3)**, 192-195 (1976).

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, 11/23/2022

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