

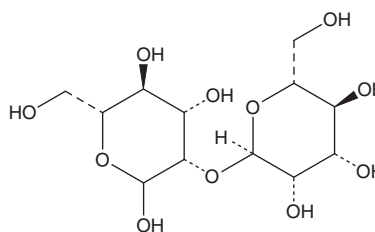
# PRODUCT INFORMATION



## 2α-Mannobiose

Item No. 21848

CAS Registry No.: 15548-39-7  
Formal Name: 2-O-α-D-mannopyranosyl-D-mannopyranose  
MF: C<sub>12</sub>H<sub>22</sub>O<sub>11</sub>  
FW: 342.3  
Purity: ≥95%  
UV/Vis.: λ<sub>max</sub>: 276 nm  
Supplied as: A crystalline 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

2α-Mannobiose is supplied as a crystalline solid. Aqueous solutions of 2α-mannobiose can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 2α-mannobiose in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

2α-Mannobiose is a disaccharide composed of two mannose molecules linked by a 1-2 glycosidic bond.<sup>1</sup> It has been immobilized on monolithic capillary columns for the affinity purification of mannose binding proteins.<sup>2</sup>

### References

1. Zhang, H., Brokman, S.M., Fang, N., *et al.* Linkage position and residue identification of disaccharides by tandem mass spectrometry and linear discriminant analysis. *Rapid Commun. Mass Spectrom.* **22**(10), 1579-1586 (2008).
2. Bedair, M. and El Rassi, Z. Affinity chromatography with monolithic capillary columns I. Polymethacrylate monoliths with immobilized mannan for the separation of mannose-binding proteins by capillary electrochromatography and nano-scale liquid chromatography. *J. Chromatogr. A* **1044**(1-2), 177-186 (2004).

#### 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, 12/01/2022

#### 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