

# PRODUCT INFORMATION



## Xylotetraose

Item No. 35387

CAS Registry No.: 22416-58-6

Formal Name: O-β-D-xylopyranosyl-(1→4)-O-β-D-xylopyranosyl-(1→4)-O-β-D-xylopyranosyl-(1→4)-D-xylose

MF: C<sub>20</sub>H<sub>34</sub>O<sub>17</sub>

FW: 546.5

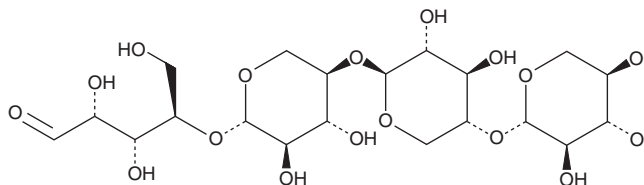
Purity: ≥98%

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

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

### Description

Xylotetraose is a xylooligosaccharide that has been found in corn and various plant-based agricultural wastes.<sup>1,2</sup>

### References

1. Chapla, D., Pandit, P., and Shah, A. Production of xylooligosaccharides from corncob xylan by fungal xylanase and their utilization by probiotics. *Bioresour. Technol.* **115**, 215-221 (2012).
2. Akpınar, O., Erdogan, K., and Bostanci, S. Enzymatic production of xylooligosaccharide from selected agricultural wastes. *Food Bioprod. Process.* **87(2)**, 145-151 (2009).

#### 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, 12/12/2022

#### 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