

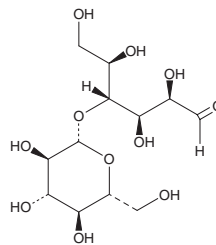
# PRODUCT INFORMATION



## D-Cellobiose

Item No. 20364

<b>CAS Registry No.:</b>	528-50-7
<b>Formal Name:</b>	4-O-β-D-glucopyranosyl-D-glucose
<b>Synonyms:</b>	Cellobiose 90, D-(+)-Cellobiose
<b>MF:</b>	C <sub>12</sub> H <sub>22</sub> O <sub>11</sub>
<b>FW:</b>	342.3
<b>Purity:</b>	≥95%
<b>Supplied as:</b>	A crystalline solid
<b>Storage:</b>	Room temperature
<b>Stability:</b>	As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



### Laboratory Procedures

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

### Description

D-Cellobiose is a disaccharide that consists of two β-glucose molecules linked by a β-1,4 glycosidic bond. It can be used as a substrate of β-glucosidase.<sup>1</sup>

### Reference

1. Lawley, B., *et al.* Whole-transcriptome shotgun sequencing (RNA-seq) screen reveals upregulation of cellobiose and motility operons of *Lactobacillus ruminis* L5 during growth on tetrasaccharides derived from barley β-glucan. *Appl. Environ. Microbiol.* **79(18)**, 5661-5669 (2013).

#### 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, 02/23/2017

#### 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