

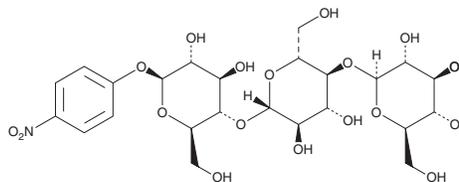
PRODUCT INFORMATION



4-Nitrophenyl β -D-Cellotrioside

Item No. 19523

CAS Registry No.: 106927-48-4
Formal Name: 4-nitrophenyl O- β -D-glucopyranosyl-(1 \rightarrow 4)-O- β -D-glucopyranosyl-(1 \rightarrow 4)- β -D-glucopyranoside
Synonym: *p*-Nitrophenyl β -D-Cellotrioside
MF: C₂₄H₃₅NO₁₈
FW: 625.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 221, 296 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

4-Nitrophenyl β -D-Cellotrioside is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-nitrophenyl β -D-cellotrioside in the solvent of choice. 4-Nitrophenyl β -D-Cellotrioside is slightly soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

4-Nitrophenyl β -D-cellotrioside is a chromogenic substrate for endoglucanases and cellobiohydrolases.¹ Hydrolysis of this substrate by these enzymes liberates 4-nitrophenol, which generates a yellow color that is measured by monitoring absorbance at 405 nm.

Reference

1. Hitomi, J., Hatada, Y., Kawaminami, S., *et al.* Amino acid sequence and stereoselective hydrolytic reaction of an endo-1,4- β -glucanase from a *Bacillus* strain. *Biosci. Biotechnol. Biochem.* **61(12)**, 2004-2009 (2016).

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, 11/03/2022

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