

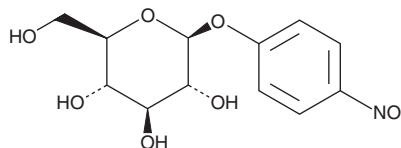
PRODUCT INFORMATION



4-Nitrophenyl β -D-Glucopyranoside

Item No. 35109

CAS Registry No.: 2492-87-7
Synonyms: *p*-Nitrophenyl β -D-glucopyranoside, PNPG
MF: C₁₂H₁₅NO₈
FW: 301.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 221, 297 nm
Supplied as: A 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-glucopyranoside is supplied as a solid. A stock solution may be made by dissolving the 4-nitrophenyl β -D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. 4-Nitrophenyl β -D-glucopyranoside is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 4-nitrophenyl β -D-glucopyranoside in ethanol is approximately 5 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 4-nitrophenyl β -D-glucopyranoside can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 4-nitrophenyl β -D-glucopyranoside in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

4-Nitrophenyl β -D-glucopyranoside is a chromogenic substrate for β -glucosidase.¹ Upon hydrolysis by β -glucosidase, 4-nitrophenol is released, which can be quantified by colorimetric detection at 405 nm as a measure of β -glucosidase activity.

Reference

1. Britton, J., Raston, C.L., and Weiss, G.A. Rapid protein immobilization for thin film continuous flow biocatalysis. *Chem. Commun. (Camb)* **52(66)**, 10159-10162 (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, 10/20/2022

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