

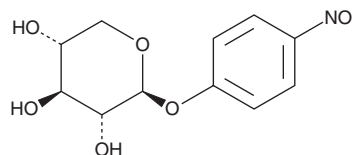
# PRODUCT INFORMATION



## 4-Nitrophenyl $\beta$ -D-Xylopyranoside

Item No. 38883

**CAS Registry No.:** 2001-96-9  
**Formal Name:** (2S,3R,4S,5R)-2-(4-nitrophenoxy)tetrahydro-2H-pyran-3,4,5-triol  
**Synonyms:** p-Nitrophenyl  $\beta$ -D-Xylopyranoside, NSC 371094, PNPX  
**MF:** C<sub>11</sub>H<sub>13</sub>NO<sub>7</sub>  
**FW:** 271.2  
**Purity:**  $\geq 98\%$   
**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  $\beta$ -D-xylopyranoside is supplied as a solid. A stock solution may be made by dissolving the 4-nitrophenyl  $\beta$ -D-xylopyranoside in the solvent of choice, which should be purged with an inert gas. 4-Nitrophenyl  $\beta$ -D-xylopyranoside is soluble in acetonitrile.

### Description

4-Nitrophenyl  $\beta$ -D-xylopyranoside is a chromogenic substrate for  $\beta$ -xylosidase.<sup>1</sup> Upon hydrolysis by  $\beta$ -xylosidase, 4-nitrophenol is released, which can be quantified by colorimetric detection at 405 nm as a measure of  $\beta$ -xylosidase activity.

### Reference

1. de Morais Souto, B., de Araújo, A.C.B., Hamann, P.R.V., *et al.* Functional screening of a Caatinga goat (*Capra hircus*) rumen metagenomic library reveals a novel GH3  $\beta$ -xylosidase. *PLoS One* **16**(1), e0245118 (2021).

#### 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, 12/04/2023

#### 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](http://WWW.CAYMANCHEM.COM)