

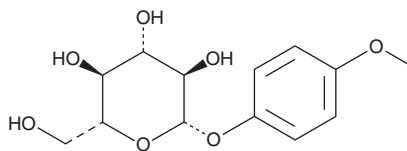
PRODUCT INFORMATION



4-Methoxyphenyl-β-D-Glucopyranoside

Item No. 35818

CAS Registry No.: 6032-32-2
Synonyms: Methylarbutin, *p*-Methoxyphenyl-β-D-Glucopyranoside
MF: C₁₃H₁₈O₇
FW: 286.3
Purity: ≥98%
UV/Vis.: λ_{max}: 224, 246 252, 280 nm
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

4-Methoxyphenyl-β-D-glucopyranoside is supplied as a solid. A stock solution may be made by dissolving the 4-methoxyphenyl-β-D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. 4-Methoxyphenyl-β-D-glucopyranoside is soluble in DMSO.

Description

4-Methoxyphenyl-β-D-glucopyranoside is an aryl glycoside that has been found in *Pterospermum heterophyllum* roots and has ice recrystallization inhibitory activity.^{1,2} It reduces mean grain size in a cell-free ice recrystallization assay and increases the percentage of intact isolated red blood cells (RBCs) in PBS post-freeze-thaw when used at a concentration of 100 mM.² 4-Methoxyphenyl-β-D-glucopyranoside has also been used as an intermediate in the synthesis of glucan oligosaccharides.³

References

1. Wang, M.-M., Li, S., Luo, G.-M., *et al.* Chemical constituents from roots of *Pterospermum heterophyllum*. *Zhongcaoyao* **43(9)**, 1699-1703 (2012).
2. Briard, J.G., Poisson, J.S., Turner, T.R., *et al.* Small molecule ice recrystallization inhibitors mitigate red blood cell lysis during freezing, transient warming and thawing. *Sci. Rep.* **6(1)**, 23619 (2016).
3. Montel, E., Hrmova, M., Fincher, G.B., *et al.* A chemoenzymatic route to conjugatable β(1→3)-glucan oligosaccharides. *Aus. J. Chem.* **62(6)**, 575-584 (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, 12/15/2022

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