

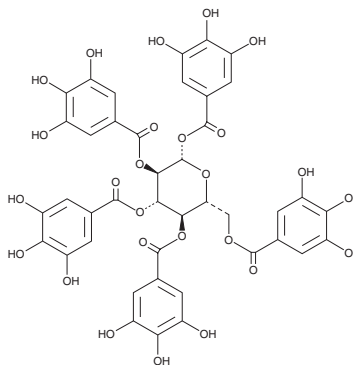
PRODUCT INFORMATION



1,2,3,4,6-Penta-O-galloyl-β-D-glucopyranose

Item No. 16007

CAS Registry No.: 14937-32-7
Formal Name: 1,2,3,4,6-pentakis(3,4,5-trihydroxybenzoate) β-D-glucopyranose
MF: C₄₁H₃₂O₂₆
FW: 940.7
Purity: ≥95%
UV/Vis.: λ_{max}: 218, 281 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

1,2,3,4,6-Penta-O-galloyl-β-D-glucopyranose is supplied as a crystalline solid. A stock solution may be made by dissolving the 1,2,3,4,6-penta-O-galloyl-β-D-glucopyranose in the solvent of choice, which should be purged with an inert gas. 1,2,3,4,6-Penta-O-galloyl-β-D-glucopyranose is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 1,2,3,4,6-penta-O-galloyl-β-D-glucopyranose in ethanol and DMSO is approximately 10 mg/ml and approximately 5 mg/ml in DMF.

1,2,3,4,6-Penta-O-galloyl-β-D-glucopyranose is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 1,2,3,4,6-penta-O-galloyl-β-D-glucopyranose should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. 1,2,3,4,6-Penta-O-galloyl-β-D-glucopyranose has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

1,2,3,4,6-Penta-O-galloyl-β-D-glucopyranose is a major component of the root of *P. suffruticosa*, a tree peony native to China and used in traditional mudanpi medicine practices.¹ At 30 μM, it has been reported to inhibit the growth of human hepatocellular carcinoma SK-HEP-1 cells by causing cell cycle arrest and suppressing the activation of NF-κB.¹

Reference

- Oh, G.S., Pae, H.O., Oh, H., *et al.* In vitro anti-proliferative effect of 1,2,3,4,6-penta-O-galloyl-beta-D-glucose on human hepatocellular carcinoma cell line, SK-HEP-1 cells. *Cancer Lett.* **174(1)**, 17-24 (2001).

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, 10/31/2022

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