

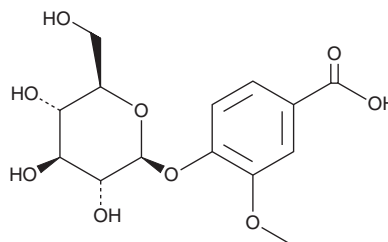
PRODUCT INFORMATION



Vanillic Acid 4-β-D-glucopyranoside

Item No. 19759

CAS Registry No.: 32142-31-7
Formal Name: 4-(β-D-glucopyranosyloxy)-3-methoxy-benzoic acid
MF: C₁₄H₁₈O₉
FW: 330.3
Purity: ≥98%
UV/Vis.: λ_{max}: 247, 288 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

Vanillic acid 4-β-D-glucopyranoside is supplied as a crystalline solid. A stock solution may be made by dissolving the vanillic acid 4-β-D-glucopyranoside in the solvent of choice, which should be purged with an inert gas. Vanillic acid 4-β-D-glucopyranoside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of vanillic acid 4-β-D-glucopyranoside in these solvents is approximately 20 and 10 mg/ml, respectively.

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 vanillic acid 4-β-D-glucopyranoside can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of vanillic acid 4-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

Vanillic Acid 4-β-D-glucopyranoside can be isolated from the fruits of *C. annuum* as well as the leaves of various additional plants. It belongs to a class of compounds known as hydrolyzable tannins and can be phytotoxic against different species.¹

Reference

1. Mkohtar, M., Soukup, J., Donato, P., *et al.* Determination of the polyphenolic content of a *C. annuum* L. extract by liquid chromatography coupled to photodiode array and mass spectrometry detection and evaluation of its biological activity. *J. Sep. Sci.* **38**(2), 171-178 (2015).

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