

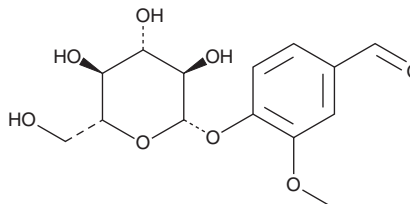
PRODUCT INFORMATION



Vanillin 4-O- β -D-Glucoside

Item No. 35908

CAS Registry No.: 494-08-6
Formal Name: 4-(β -D-glucopyranosyloxy)-3-methoxybenzaldehyde
Synonym: Glucovanillin
MF: C₁₄H₁₈O₈
FW: 314.3
Purity: \geq 98%
UV/Vis.: λ_{max} : 226, 271, 306 nm
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Synthetic



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

Laboratory Procedures

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

Description

Vanillin 4-O- β -D-glucoside is a vanillin precursor.¹ It is the main storage form of vanillin in green vanilla beans.

Reference

1. Paradío, V.T., Flores, A., López, K.M., *et al.* Effect of endogenous and exogenous enzymatic treatment of green vanilla beans on extraction of vanillin and main aromatic compounds. *J. Food Sci. Technol.* **55**(6), 2059-2067 (2018).

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