

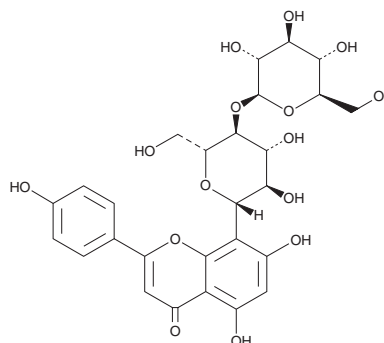
PRODUCT INFORMATION



Vitexin-4''-O-glucoside

Item No. 34923

CAS Registry No.: 178468-00-3
Formal Name: 8-(4-O-β-D-glucopyranosyl-β-D-glucopyranosyl)-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one
MF: C₂₇H₃₀O₁₅
FW: 594.5
Purity: ≥95%
UV/Vis.: λ_{max}: 216, 271, 335 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Crataegi fructus*



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

Laboratory Procedures

Vitexin-4''-O-glucoside is supplied as a solid. A stock solution may be made by dissolving the vitexin-4''-O-glucoside in the solvent of choice, which should be purged with an inert gas. Vitexin-4''-O-glucoside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of vitexin-4''-O-glucoside in these solvents is approximately 1 mg/ml.

Vitexin-4''-O-glucoside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vitexin-4''-O-glucoside should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Vitexin-4''-O-glucoside has a solubility of approximately 0.50 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Vitexin-4''-O-glucoside is a flavonoid glycoside and major component of *C. pinnatifida* that has antioxidant activity.¹ It reduces tertbutyl hydroperoxide-induced DNA damage, lipid peroxidation, and cell death in ECV304 cells when used at concentrations ranging from 16 to 128 μM.

Reference

1. Li, H.-B., Ying, X.X., and Lu, J. The mechanism of vitexin-4''-O-glucoside protecting ECV-304 cells against tertbutyl hydroperoxide induced injury. *Nat. Prod. Res.* **24(18)**, 1695-1703 (2010).

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, 12/02/2022

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