

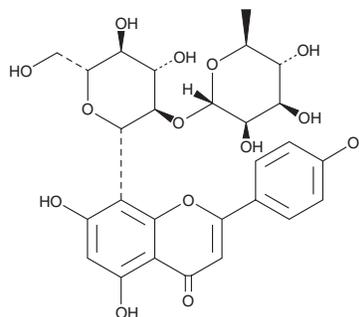
PRODUCT INFORMATION



Vitexin 2-O-rhamnoside

Item No. 26816

CAS Registry No.: 64820-99-1
Formal Name: 8-[2-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranosyl]-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one
MF: C₂₇H₃₀O₁₄
FW: 578.5
Purity: \geq 95%
UV/Vis.: λ_{max} : 215, 271, 332 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Crataegus* sp.



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

Laboratory Procedures

Vitexin 2-O-rhamnoside is supplied as a crystalline solid. A stock solution may be made by dissolving the vitexin 2-O-rhamnoside in the solvent of choice, which should be purged with an inert gas. Vitexin 2-O-rhamnoside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of vitexin 2-O-rhamnoside in these solvents is approximately 15 and 10 mg/ml, respectively.

Vitexin 2-O-rhamnoside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vitexin 2-O-rhamnoside should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Vitexin 2-O-rhamnoside has a solubility of approximately 0.5 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 2-O-rhamnoside is a flavonoid glycoside that has been found in *B. vulgaris* species.¹ It inhibits DNA synthesis in MCF-7 breast cancer cells (IC₅₀ = 17.5 μ M).

Reference

1. Ninfali, P., Bacchiocca, M., Antonelli, A., *et al.* Characterization and biological activity of the main flavonoids from Swiss Chard (*Beta vulgaris* subspecies *cycla*). *Phytomedicine* **14**(2-3), 216-221 (2007).

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, 12/07/2022

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