

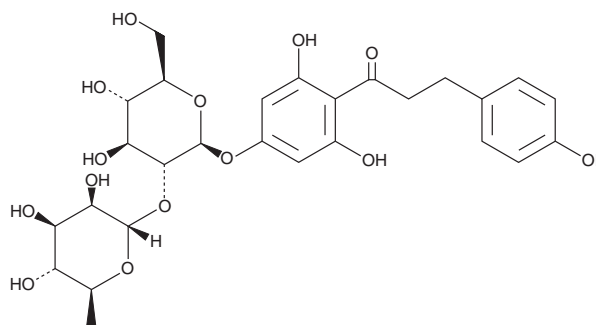
PRODUCT INFORMATION



Naringin Dihydrochalcone

Item No. 27791

CAS Registry No.: 18916-17-1
Formal Name: 1-[4-[[2-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranosyl]oxy]-2,6-dihydroxyphenyl]-3-(4-hydroxyphenyl)-1-propanone
MF: C₂₇H₃₄O₁₄
FW: 582.6
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 226, 284 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/Pomelo peel



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

Laboratory Procedures

Naringin dihydrochalcone is supplied as a crystalline solid. A stock solution may be made by dissolving the naringin dihydrochalcone in the solvent of choice, which should be purged with an inert gas. Naringin dihydrochalcone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of naringin dihydrochalcone in these solvents is approximately 30 mg/ml.

Description

Naringin dihydrochalcone is an artificial sweetener and flavonoid that has been isolated from various plants and has antioxidant activity.^{1,2} It scavenges ABTS (Item No. 27317), oxygen, and 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals in cell free assays (IC₅₀s = 24, 322.8, and 318.9 μ M, respectively).²

References

1. Gentili, B. and Horowitz, R.M. Chromatography of dihydrochalcone sweeteners and related compounds: A reagent for detecting dihydrochalcones *J. Chromatogr.* **63(5730)**, 467-469 (1971).
2. Li, X., Chen, B., Xie, H., *et al.* Antioxidant structure-activity relationship analysis of five dihydrochalcones. *Molecules* **23(5)**, E1162 (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, 11/09/2022

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