

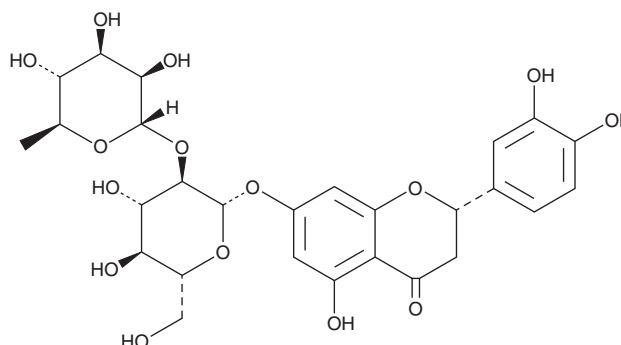
PRODUCT INFORMATION



Neoeriocitrin

Item No. 25839

CAS Registry No.: 13241-32-2
Formal Name: (2S)-7-[[2-O-(6-deoxy- α -L-mannopyranosyl)- β -D-glucopyranosyl]oxy]-2-(3,4-dihydroxyphenyl)-2,3-dihydro-5-hydroxy-4H-1-benzopyran-4-one
Synonym: Eriodictyol 7-O-neohesperidoside
MF: C₂₇H₃₂O₁₅
FW: 596.5
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 285 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/Aurantii fructus immaturus



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

Laboratory Procedures

Neoeriocitrin is supplied as a crystalline solid. A stock solution may be made by dissolving the neoeriocitrin in the solvent of choice. Neoeriocitrin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of neoeriocitrin in these solvents is approximately 1, 30, and 15 mg/ml, respectively.

Neoeriocitrin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, neoeriocitrin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Neoeriocitrin has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Neoeriocitrin is a flavonoid that has been found in *C. paradisi* and has antioxidative and osteogenic activities.^{1,2} It inhibits the production of 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals by 17.2% and the formation of superoxide radicals by 48.3% in cell-free assays, as well as decreases the rate of LDL oxidation *ex vivo* in isolated hamster plasma. Neoeriocitrin (2 μ g/ml) increases proliferation of MC3T3-E1 osteoblast precursor cells and reverses the antiproliferative effect of the MEK1 inhibitor PD 98059 (Item No. 10006726).² It also increases the expression of the osteogenic differentiation markers Runx2, Type 1 collagen, and osteocalcin.

References

1. Yu, J., Wang, L., Walzem, R.L., *et al.* Antioxidant activity of citrus limonoids, flavonoids, and coumarins. *J. Agric. Food Chem.* **53**(6), 2009-2014 (2005).
2. Li, L., Zeng, Z., and Cai, G. Comparison of neoeriocitrin and naringin on proliferation and osteogenic differentiation in MC3T3-E1. *Phytomedicine* **18**(11), 985-989 (2011).

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, 12/19/2022

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