

PRODUCT INFORMATION



Quercitrin

Item No. 19866

CAS Registry No.: 522-12-3
Formal Name: 3-[[6-deoxy- α -L-mannopyranosyl)oxy]-2-(3,4-dihydroxyphenyl)-5,7-dihydroxy-4H-1-benzopyran-4-one
Synonyms: C.I. 75720, NSC 9221,

Quercetin 3-rhamnoside,
Quercetin 3-L-rhamnoside

MF: C₂₁H₂₀O₁₁

FW: 448.4

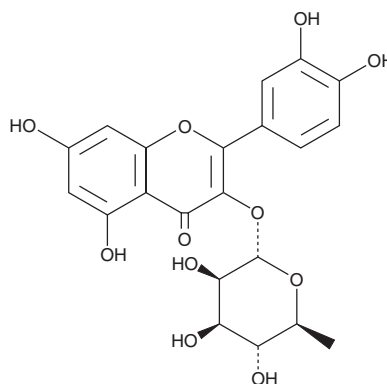
Purity: \geq 98%

UV/Vis.: λ_{max} : 254, 351 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

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

Quercitrin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, quercitrin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Quercitrin 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

Quercitrin is a glycoside formed from the flavonoid quercetin (Item No. 10005169) and the deoxy sugar rhamnose. It can be found in a wide range of medicinal plants and has been reported to have antioxidant, antiviral, and anti-inflammatory properties.^{1,2,3}

References

1. Li, X., Jiang, Q., Wang, T., *et al.* Comparison of the antioxidant effects of quercitrin and isoquercitrin: Understanding the role of the 6''-OH group. *Molecules* **21(9)**, (2016).
2. Chiow, K.H., Phoon, M. C., Putti, T., *et al.* Evaluation of antiviral activities of *Houttuynia cordata* Thunb. extract, quercetin, quercitrin and cinanserin on murine coronavirus and dengue virus infection. *Asian Pac. J. Trop. Med.* **9(1)**, 1-7 (2016).
3. Jurikova, T., Rop, J. Mlcek, O., *et al.* Phenolic profile of edible honeysuckle berries (genus *Lonicera*) and their biological effects. *Molecules* **17(1)**, 61-79 (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, 12/22/2022

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