

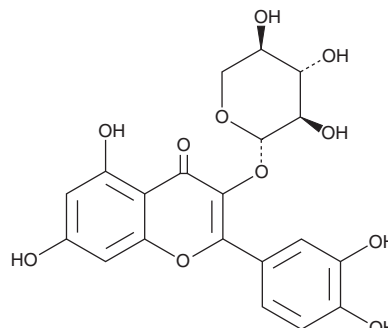
PRODUCT INFORMATION



Quercetin 3-O-β-D-xylopyranoside

Item No. 35141

CAS Registry No.: 549-32-6
Formal Name: 2-(3,4-dihydroxyphenyl)-5,7-dihydroxy-3-(β-D-xylopyranosyloxy)-4H-1-benzopyran-4-one
Synonyms: Reinutrin, Reynoutrin
MF: C₂₀H₁₈O₁₁
FW: 434.4
Purity: ≥98%
UV/Vis.: λ_{max}: 258, 361 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Psidium guajava*



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

Laboratory Procedures

Quercetin 3-O-β-D-xylopyranoside is supplied as a solid. A stock solution may be made by dissolving the Quercetin 3-O-β-D-xylopyranoside in the solvent of choice, which should be purged with an inert gas. Quercetin 3-O-β-D-xylopyranoside is soluble in DMSO.

Description

Quercetin 3-O-β-D-xylopyranoside is a flavonoid glycoside that has been found in *N. candida* and has diverse biological activities.¹⁻³ It reduces malondialdehyde (MDA) levels in isolated rat brain when used at a concentration of 10 μg/ml.¹ Quercetin 3-O-β-D-xylopyranoside (10, 50, and 100 μg/ml) reduces LPS- and cerulein-induced increases in reactive oxygen species (ROS) production and levels of GRP78, a marker of endoplasmic reticulum stress, in an AR42J rat pancreatic acinar cell model of pancreatitis.² *In vivo*, quercetin 3-O-β-D-xylopyranoside (12.5, 25, and 50 mg/kg) decreases serum TNF-α and IL-6 levels, myocardial MDA levels, and myocardial fibrosis, as well as improves cardiac function, in a rat model of ischemic heart failure.³

References

1. Liu, R.-N., Wang, W., Ding, Y., *et al.* A new flavonol glycoside and activity of compounds from the flower of *Nymphaea candida*. *J. Asian Nat. Prod. Res.* **9**(3-5), 333-338 (2007).
2. Seo, J.Y., Pandey, R.P., Lee, J., *et al.* Quercetin 3-O-xyloside ameliorates acute pancreatitis *in vitro* via the reduction of ER stress and enhancement of apoptosis. *Phytomedicine* **55**, 40-49 (2019).
3. Yang, W., Tu, H., Tang, K., *et al.* Reynoutrin improves ischemic heart failure in rats via targeting S100A1. *Front. Pharmacol.* **12**, 703962 (2021).

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, 11/11/2022

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