

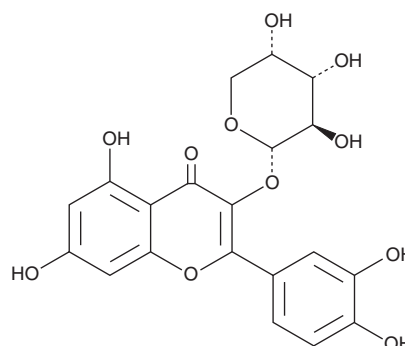
PRODUCT INFORMATION



Quercetin 3-O- α -L-arabinopyranoside

Item No. 30707

CAS Registry No.: 22255-13-6
Formal Name: 3-(α -L-arabinopyranosyloxy)-2-(3,4-dihydroxyphenyl)-5,7-dihydroxy-4H-1-benzopyran-4-one
Synonym: Guajaverin
MF: C₂₀H₁₈O₁₁
FW: 434.4
Purity: \geq 98%
UV/Vis.: λ_{max} : 256, 359 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 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- α -L-arabinopyranoside is supplied as a crystalline solid. A stock solution may be made by dissolving the quercetin 3-O- α -L-arabinopyranoside in the solvent of choice, which should be purged with an inert gas. Quercetin 3-O- α -L-arabinopyranoside is soluble in DMSO.

Description

Quercetin 3-O- α -L-arabinopyranoside is a phenol that has been found in *Byrsonima crassa* and has antioxidant activity.^{1,2} It scavenges 2,2-diphenyl-1-picrylhydrazyl (DPPH; Item No. 14805) radicals in a cell-free assay (IC₅₀ = 48.6 μ M).¹ Quercetin 3-O- α -L-arabinopyranoside inhibits *H. pylori*-induced oxidative burst of isolated rat polymorphonuclear (PMN) neutrophils (IC₅₀ = 75.3 μ M).² It reduces N-retinylidene-N-retinylethanolamine- and blue light-induced death of ARPE-19 retinal epithelial cells in a concentration-dependent manner.³ *In vivo*, quercetin 3-O- α -L-arabinopyranoside (25, 50, and 200 mg/kg) prevents blue light-induced retinal degeneration in a mouse model of macular degeneration.

References

1. Joshi, K.R., Devkota, H.P., Watanabe, T., *et al.* Phenolic compounds from the flowers of Nepalese medicinal plant *Aconogonon molle* and their DPPH free radical-scavenging activities. *Nat. Prod. Res.* **28(23)**, 2208-2210 (2014).
2. Bonacorsi, C., Raddi, M.S.G., da Fonseca, L.M., *et al.* Effect of *Byrsonima crassa* and phenolic constituents on *Helicobacter pylori*-induced neutrophils oxidative burst. *Int. J. Mol. Sci.* **13(1)**, 133-141 (2012).
3. Kim, J., Jin, H.L., Jang, D.S., *et al.* Quercetin-3-O- α -l-arabinopyranoside protects against retinal cell death via blue light-induced damage in human RPE cells and Balb-c mice. *Food Funct.* **9(4)**, 2171-2183 (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, 12/12/2022

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