

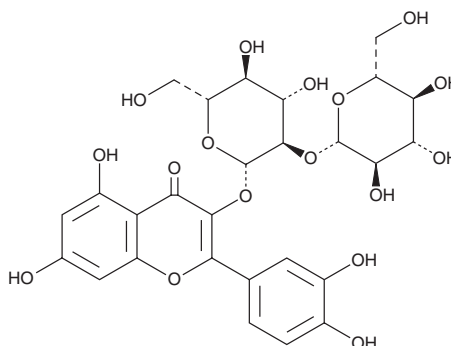
PRODUCT INFORMATION



Quercetin 3-O-sophoroside

Item No. 28592

CAS Registry No.:	18609-17-1
Formal Name:	2-(3,4-dihydroxyphenyl)-3-[(2-O-β-D-glucopyranosyl-β-D-glucopyranosyl)oxy]-5,7-dihydroxy-4H-1-benzopyran-4-one
Synonyms:	Baimaside, Quercetin 3-sophoroside
MF:	C ₂₇ H ₃₀ O ₁₇
FW:	626.5
Purity:	≥98%
UV/Vis.:	λ _{max} : 207, 257, 355 nm
Supplied as:	A crystalline solid
Storage:	-20°C
Stability:	≥4 years
Item Origin:	Plant/ <i>A. venetum</i> L.



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

Laboratory Procedures

Quercetin 3-O-sophoroside is supplied as a crystalline solid. A stock solution may be made by dissolving the quercetin 3-O-sophoroside in the solvent of choice, which should be purged with an inert gas. Quercetin 3-O-sophoroside is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of quercetin 3-O-sophoroside in these solvents is approximately 10 and 3 mg/ml, respectively.

Quercetin 3-O-sophoroside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, quercetin 3-O-sophoroside should first be dissolved in DMSO or DMF and then diluted with the aqueous buffer of choice. Quercetin 3-O-sophoroside has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) and approximately approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Quercetin 3-O-sophoroside is a flavonoid glycoside that has been found in *B. napus* (rapeseed) and has antioxidant activity.¹⁻³ Quercetin 3-O-sophoroside has a relative antioxidant capacity of 1.45 compared to Trolox (Item No. 10011659) in an ABTS (Item No. 27317) assay.³ It also inhibits lipid peroxidation in a cell-free assay using phospholipid liposomes with an IC₅₀ value of 9.2 μM.

References

1. Ferreres, F., Taveira, M., Pereira, D.M., *et al.* Tomato (*Lycopersicon esculentum*) seeds: new flavonols and cytotoxic effect. *J. Agric. Food Chem.* **58(5)**, 2854-2861 (2010).
2. Wang, Y., Meng, G., Chen, S., *et al.* Correlation analysis of phenolic contents and antioxidation in yellow- and black-seeded *Brassica napus*. *Molecules* **23(7)**, e1815 (2018).
3. Plumb, G.W., Price, K.R., Modes, M.J.C., *et al.* Antioxidant properties of the major polyphenolic compounds in broccoli. *Free Radic. Res.* **27(4)**, 429-435 (1997).

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