

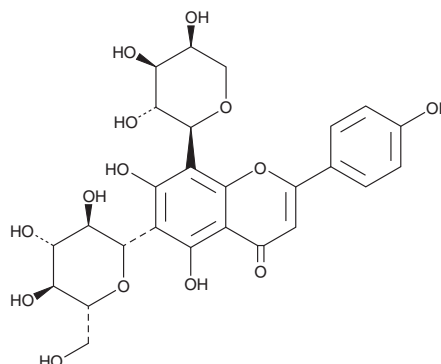
PRODUCT INFORMATION



Schaftoside

Item No. 27950

CAS Registry No.: 51938-32-0
Formal Name: 8- α -L-arabinopyranosyl-6- β -D-glucopyranosyl-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one
Synonym: Apigenin 6-C-glucoside-8-C-arabinoside
MF: C₂₆H₂₈O₁₄
FW: 564.5
Purity: \geq 98%
UV/Vis.: λ_{max} : 217, 274, 335 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years
Item Origin: Plant/*Desmodium styracifolium*



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

Laboratory Procedures

Schaftoside is supplied as a crystalline solid. A stock solution may be made by dissolving the schaftoside in the solvent of choice, which should be purged with an inert gas. Schaftoside is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of schaftoside in these solvents is approximately 30 and 20 mg/ml, respectively.

Schaftoside is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, schaftoside should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Schaftoside has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Schaftoside is a flavonoid C-glycoside that has been found in *C. annuum* and has antioxidant and anti-inflammatory activities.^{1,2} It inhibits xanthine oxidase with an IC₅₀ value of 51.4 μ M and has antioxidant activity in a β -carotene-linoleic acid bleaching assay (IC₅₀ = 46.3 μ M).¹ Schaftoside (400 μ g/kg) inhibits neutrophil influx into mouse lung by 62% in a model of airway inflammation induced by aerosol administration of LPS.²

References

1. Materska, M. Flavone C-glycosides from *Capsicum annuum* L.: Relationships between antioxidant activity and lipophilicity. *Eur. Food Res. Technol.* **240(3)**, 549-557 (2015).
2. De Melo, G.O., Muzitano, M.F., Legora-Machado, A., et al. C-Glycosylflavones from the aerial parts of *Eleusine indica* inhibit LPS-induced mouse lung inflammation. *Planta Med.* **71(4)**, 362-363 (2005).

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, 08/25/2023

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