

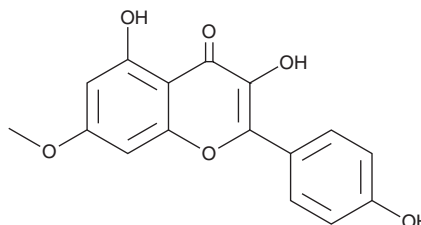
PRODUCT INFORMATION



Rhamnocitrin

Item No. 36272

CAS Registry No.: 569-92-6
Formal Name: 3,5-dihydroxy-2-(4-hydroxyphenyl)-7-methoxy-4H-1-benzopyran-4-one
Synonyms: C.I. 75650, Kaempferol 7-O-methyl ether
MF: $C_{16}H_{12}O_6$
FW: 300.3
Purity: $\geq 95\%$
UV/Vis.: $\lambda_{\max}$: 267, 366 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Belamcanda chinensis*



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

Laboratory Procedures

Rhamnocitrin is supplied as a solid. A stock solution may be made by dissolving the rhamnocitrin in the solvent of choice, which should be purged with an inert gas. Rhamnocitrin is soluble in the organic solvent methanol.

Description

Rhamnocitrin is a flavonoid that has been found in *C. erythrophyllum* and has diverse biological activities.^{1,2} It is active against *S. sonei*, *S. aureus*, and *M. luteus* when used at concentrations ranging from 25 to 100 $\mu\text{g/ml}$.¹ Rhamnocitrin scavenges DPPH (Item No. 14805) radicals in a cell-free assay ($\text{IC}_{50} = \sim 25 \mu\text{M}$).² It also inhibits LPS- and IFN- γ -induced production of nitric oxide (NO), TNF- α , and IL-6 in isolated mouse peritoneal macrophages.

References

1. Martini, N.D., Katerere, D.R.P., and Eloff, J.N. Biological activity of five antibacterial flavonoids from *Combretum erythrophyllum* (Combretaceae). *J. Ethnopharmacol.* **93**(2-3), 207-212 (2004).
2. Fang, S.-H., Rao, Y.K., and Tzeng, Y.-M. Anti-oxidant and inflammatory mediator's growth inhibitory effects of compounds isolated from *Phyllanthus urinaria*. *J. Ethnopharmacol.* **116**(2), 333-340 (2008).

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, 01/03/2023

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