

PRODUCT INFORMATION



Mauritianin

Item No. 34326

CAS Registry No.: 109008-28-8

Formal Name: 3-[[O-6-deoxy- α -L-mannopyranosyl-(1 $\rightarrow$ 2)-O-[6-deoxy- α -L-mannopyranosyl-(1 $\rightarrow$ 6)]]- β -D-galactopyranosyl]oxy]-5,7-dihydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one

Synonyms: Kaempferol 3-O-(2'',6''-di-O- α -rhamnopyranosyl)- β -galactopyranoside, Kaempferol 3-O-{ α -L-rhamnopyranosyl(1 $\rightarrow$ 6)-[α -L-rhamnopyranosyl-(1 $\rightarrow$ 2)]}- β -D-galactopyranoside

MF: C₃₃H₄₀O₁₉

FW: 740.7

Purity: $\geq 95\%$

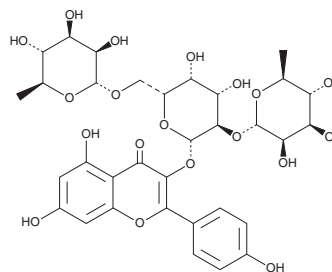
UV/Vis.: $\lambda_{\max}$: 267, 347 nm

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

Item Origin: Plant/*Hypericum japonicum*



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

Laboratory Procedures

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

Description

Mauritianin is a flavonoid glycoside that has been found in *R. anguifera* and is a DNA topoisomerase I inhibitor.¹ It stabilizes the covalent complex between DNA and recombinant human topoisomerase I when used at a concentration of 100 μ M.

Reference

1. Ma, J., Jones, S.H., Marshall, R.A., et al. DNA topoisomerase I inhibitors from *Rinorea anguifera*. *Bioorg. Med. Chem. Lett.* **15**(3), 813-816 (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, 10/29/2024

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