

PRODUCT INFORMATION

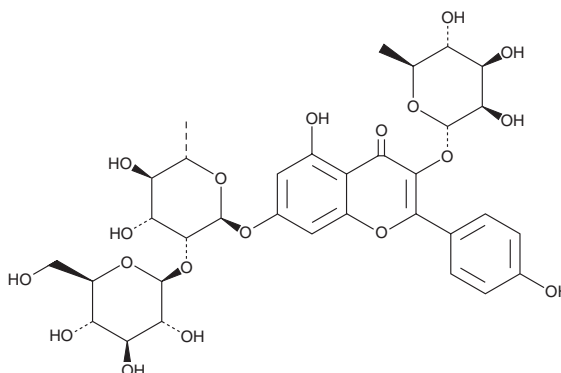


Grosvenorine

Item No. 34810

CAS Registry No.: 156980-60-8
Formal Name: 7-[[6-deoxy-2-O-β-D-glucopyranosyl-α-L-mannopyranosyl]oxy]-3-[[6-deoxy-α-L-mannopyranosyl]oxy]-5-hydroxy-2-(4-hydroxyphenyl)-4H-1-benzopyran-4-one

MF: C₃₃H₄₀O₁₉
FW: 740.7
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 266, 345 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Siraitia grosvenorii*



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

Laboratory Procedures

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

Description

Grosvenorine is a flavonol glycoside that has been found in *S. grosvenorii* and has antibacterial and antioxidant activities.¹ It is active against methicillin-sensitive and -resistant *S. aureus*, as well as *Enterococcus*, in disc diffusion assays when used at a concentration of 250 μg/ml. Grosvenorine scavenges ABTS (Item No. 27317) and DPPH (Item No. 14805) radicals in cell-free assays in a concentration-dependent manner.

Reference

1. Wang, M., Xing, S., Luu, T., *et al.* The gastrointestinal tract metabolism and pharmacological activities of grosvenorine, a major and characteristic flavonoid in the fruits of *Siraitia grosvenorii*. *Chem. Biodivers.* **12(11)**, 1652-1664 (2015).

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, 11/21/2022

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