

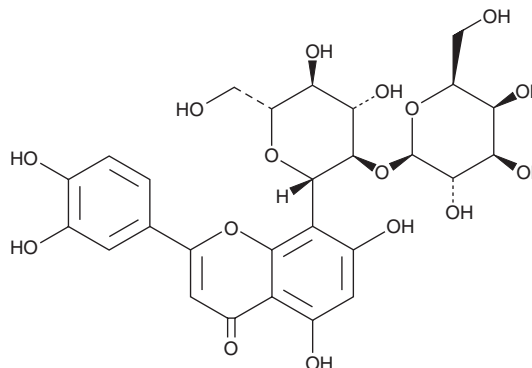
PRODUCT INFORMATION



2''-O-β-L-Galactopyranosylorientin

Item No. 36156

CAS Registry No.: 861691-37-4
Formal Name: 2-(3,4-dihydroxyphenyl)-8-(2-O-β-L-galactopyranosyl-β-D-glucopyranosyl)-5,7-dihydroxy-4H-1-benzopyran-4-one
Synonym: Orientin 2''-O-β-L-galactoside
MF: C₂₇H₃₀O₁₆
FW: 610.5
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Lophatherum gracile*



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

Laboratory Procedures

2''-O-β-L-Galactopyranosylorientin is supplied as a solid. A stock solution may be made by dissolving the 2''-O-β-L-Galactopyranosylorientin in the solvent of choice, which should be purged with an inert gas. 2''-O-β-L-Galactopyranosylorientin is soluble in DMSO.

Description

2''-O-β-L-Galactopyranosylorientin is a flavonoid glycoside that has been found in *T. chinensis* and has anti-inflammatory activity.¹ It decreases LPS-induced production of nitric oxide (NO), IL-6, and TNF-α in RAW 264.7 cells when used at concentrations ranging from 100 to 200 μM.

Reference

1. Liu, L.-J., Hu, X.-H., Guo, L.-N., *et al.* Anti-inflammatory effect of the compounds from the flowers of *Trollius chinensis*. *Pak. J. Pharm. Sci.* **31(5)**, 1951-1957 (2018).

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, 11/09/2022

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