

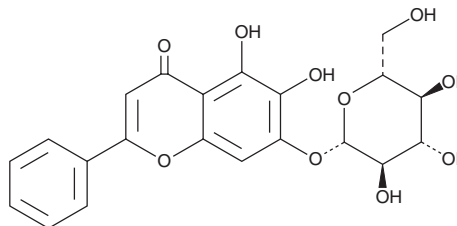
PRODUCT INFORMATION



Oroxin A

Item No. 34924

CAS Registry No.: 57396-78-8
Formal Name: 7-(β-D-glucopyranosyloxy)-5,6-dihydroxy-2-phenyl-4H-1-benzopyran-4-one
Synonym: Baicalein-7-O-Glucoside
MF: C₂₁H₂₀O₁₀
FW: 432.4
Purity: ≥98%
UV/Vis.: λ_{max}: 216, 246, 279, 315 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Plant/*Oroxylum indicum*



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

Laboratory Procedures

Oroxin A is supplied as a solid. A stock solution may be made by dissolving the oroxin A in the solvent of choice, which should be purged with an inert gas. Oroxin A is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of oroxin A in these solvents is approximately 1 mg/ml.

Oroxin A is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, oroxin A should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Oroxin A has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Oroxin A is a flavonoid glycoside that has been found in *O. indicum*.¹

Reference

1. Chen, L.-J., Games, D.E., and Jones, J. Isolation and identification of four flavonoid constituents from the seeds of *Oroxylum indicum* by high-speed counter-current chromatography. *J. Chromatogr. A* **988**(1), 95-105 (2003).

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, 10/19/2022

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