

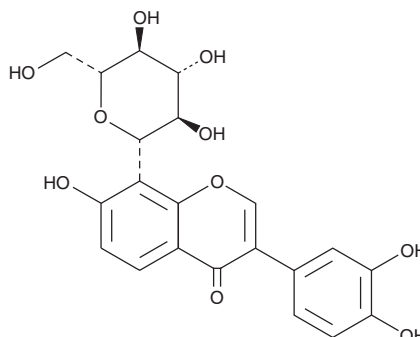
# PRODUCT INFORMATION



## 3'-hydroxy Puerarin

Item No. 36771

**CAS Registry No.:** 117060-54-5  
**Formal Name:** 3-(3,4-dihydroxyphenyl)-8-β-D-glucopyranosyl-7-hydroxy-4H-1-benzopyran-4-one  
**Synonym:** NPI 031H  
**MF:** C<sub>21</sub>H<sub>20</sub>O<sub>10</sub>  
**FW:** 432.4  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Pueraria lobata*



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

### Laboratory Procedures

3'-hydroxy Puerarin is supplied as a solid. A stock solution may be made by dissolving the 3'-hydroxy puerarin in the solvent of choice, which should be purged with an inert gas. 3'-hydroxy Puerarin is soluble in the organic solvent DMSO.

### Description

3'-hydroxy Puerarin is an isoflavonoid that has been found in *P. lobata* and has antioxidant activity.<sup>1</sup> It scavenges peroxy nitrite, nitric oxide radicals, and total reactive oxygen species (ROS) with IC<sub>50</sub> values of 1.36, 1.13, and 6.51 μM, respectively.

### Reference

1. Jin, S.E., Son, Y.K., Min, B.-S., *et al.* Anti-inflammatory and antioxidant activities of constituents isolated from *Pueraria lobata* roots. *Arch. Pharm. Res.* **35**(5), 823-837 (2012).

#### 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, 11/14/2022

#### 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