

# PRODUCT INFORMATION



## Myricetin 3-galactoside

Item No. 36467

**CAS Registry No.:** 15648-86-9  
**Formal Name:** 3-(β-D-galactopyranosyloxy)-5,7-dihydroxy-2-(3,4,5-trihydroxyphenyl)-4H-1-benzopyran-4-one  
**Synonyms:** Myricetin 3-O-β-galactopyranoside, Myricetin 3-O-β-D-galactopyranoside, Myricetin 3-O-galactoside, Myricetin 3-β-D-galactoside

**MF:** C<sub>21</sub>H<sub>20</sub>O<sub>13</sub>

**FW:** 480.4

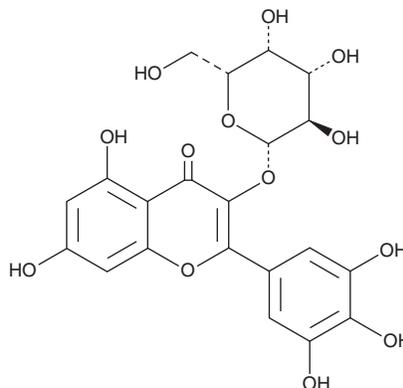
**Purity:** ≥98%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Plant/*Myrica rubra*



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

### Laboratory Procedures

Myricetin 3-galactoside is supplied as a solid. A stock solution may be made by dissolving the myricetin 3-galactoside in the solvent of choice, which should be purged with an inert gas. Myricetin 3-galactoside is soluble in acetone, chloroform, dichloromethane, ethyl acetate and DMSO.

### Description

Myricetin 3-galactoside is a flavonoid that has been found in *M. communis* and has diverse biological activities.<sup>1,2</sup> It inhibits xanthine oxidase in a cell-free assay when used at a concentration of 100 µg/ml.<sup>1</sup> Myricetin 3-galactoside is cytotoxic to K562 cells (IC<sub>50</sub> = 220 µg/ml) and inhibits hydrogen peroxide-induced malondialdehyde (MDA) formation in the same cells when used at concentrations of 110, 220, and 440 µg/ml.<sup>1</sup> *In vivo*, myricetin 3-galactoside (0.26 and 0.78 mg/kg) reduces carrageenan-induced paw edema in wild-type but not *nos2*<sup>-/-</sup> mice.<sup>2</sup> It also reduces formalin-induced paw licking in mice.

### References

- Hayder, N., Bouhlel, I., Skandrani, I., *et al.* In vitro antioxidant and antigenotoxic potentials of myricetin-3-o-galactoside and myricetin-3-o-rhamnoside from *Myrtus communis*: modulation of expression of genes involved in cell defence system using cDNA microarray. *Toxicol. In Vitro* **22(3)**, 567-581 (2008).
- de Oliveira Azevedo, A., Campos, J.J., de Souza, G.G., *et al.* Antinociceptive and anti-inflammatory effects of myricetin 3-O-β-galactoside isolated from *Davilla elliptica*: Involvement of the nitrenergic system. *J. Nat. Med.* **69(4)**, 487-493 (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, 12/19/2022

#### 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