

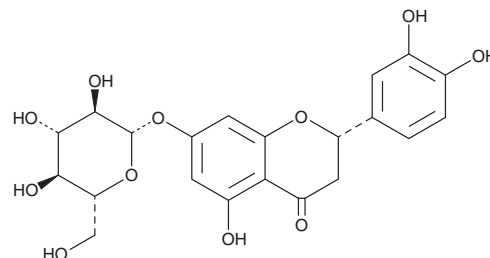
# PRODUCT INFORMATION



## Eriodictyol 7-O-glucoside

Item No. 36003

**CAS Registry No.:** 38965-51-4  
**Formal Name:** (2S)-2-(3,4-dihydroxyphenyl)-7-(β-D-glucopyranosyloxy)-2,3-dihydro-5-hydroxy-4H-1-benzopyran-4-one  
**Synonyms:** 7-O-β-D-Glucopyranosyl-eriodictyol, (2S)-Eriodictyol 7-O-β-D-glucopyranoside  
**MF:** C<sub>21</sub>H<sub>22</sub>O<sub>11</sub>  
**FW:** 450.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 283 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Viscum articulatum*



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

### Laboratory Procedures

Eriodictyol 7-O-glucoside is supplied as a solid. A stock solution may be made by dissolving the eriodictyol 7-O-glucoside in the solvent of choice, which should be purged with an inert gas. Eriodictyol 7-O-glucoside is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

### Description

Eriodictyol 7-O-glucoside is a flavanone glucoside that has been found in *V. coloratum* and has diverse biological activities.<sup>1-3</sup> It scavenges hydroxyl and superoxide anion radicals in cell-free assays (IC<sub>50</sub>s = 0.28 and 0.3 mM, respectively).<sup>1</sup> Eriodictyol 7-O-glucoside activates nuclear factor E2-related factor 2 (Nrf2) in a reporter assay.<sup>2</sup> It protects against cytotoxicity induced by cisplatin (Item No. 13119) in human renal mesangial cells (HRMCs), but not A549 lung or MDA-MB-231 breast cancer cells, when used at a concentration of 80 μM. Eriodictyol 7-O-glucoside (30 mg/kg) attenuates neurological deficits and reduces infarct volume in a rat model of cerebral ischemia induced by middle cerebral artery occlusion (MCAO).<sup>3</sup>

### References

1. Yao, H., Liao, Z.-X., Wu, Q., et al. Antioxidative flavanone glycosides from the branches and leaves of *Viscum coloratum*. *Chem. Pharm. Bull. (Tokyo)* **54**(1), 133-135 (2006).
2. Hu, Q., Zhang, D.D., Wang, L., et al. Eriodictyol-7-O-glucoside, a novel Nrf2 activator, confers protection against cisplatin-induced toxicity. *Food Chem. Toxicol.* **50**(6), 1927-1932 (2012).
3. Jing, X., Ren, D., Wei, X., et al. Eriodictyol-7-O-glucoside activates Nrf2 and protects against cerebral ischemic injury. *Toxicol. Appl. Pharmacol.* **273**(3), 672-679 (2013).

#### 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, 11/03/2022

#### 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