

# PRODUCT INFORMATION

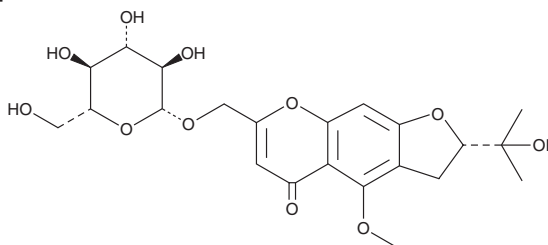


## Prim-O-Glucosylcimifugin

Item No. 34921

**CAS Registry No.:** 80681-45-4  
**Formal Name:** (2S)-7-[[β-D-glucopyranosyloxy)methyl]-2,3-dihydro-2-(1-hydroxy-1-methylethyl)-4-methoxy-5H-furo[3,2-g][1]benzopyran-5-one  
**Synonyms:** 1''-O-Glucosylcimifugin, PGCN

**MF:** C<sub>22</sub>H<sub>28</sub>O<sub>11</sub>  
**FW:** 468.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 215, 292 nm  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Cimicifugae rhizoma*



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

### Laboratory Procedures

Prim-O-glucosylcimifugin is supplied as a solid. A stock solution may be made by dissolving the prim-O-glucosylcimifugin in the solvent of choice, which should be purged with an inert gas. Prim-O-glucosylcimifugin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of prim-O-glucosylcimifugin in these solvents is approximately 15 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of prim-O-glucosylcimifugin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of prim-O-glucosylcimifugin in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Prim-O-glucosylcimifugin is a chromone that has been found in *Cimicifuga dahurica* and has diverse biological activities.<sup>1-4</sup> It inhibits the proliferation of EGF-stimulated MCF-7 breast cancer cells when used at a concentration of 10 μM.<sup>1</sup> Prim-O-glucosylcimifugin (15, 50, and 100 μg/ml) reduces LPS-induced increases in the production of nitric oxide (NO), TNF-α, IL-6, and IL-1β in RAW 264.7 macrophages.<sup>2</sup> It decreases pulmonary edema in a mouse model of LPS-induced acute lung injury when administered at doses ranging from 2.5 to 10 mg/kg.<sup>3</sup> Prim-O-glucosylcimifugin inhibits formalin-induced tonic pain in rats (ED<sub>50</sub> = 1.6 mg/animal).<sup>4</sup>

### References

1. Huyan, C.T.T., Luyen, B.T.T., Khan, G.J., *et al.* Chemical constituents from *Cimicifuga dahurica* and their anti-proliferative effects on MCF-7 breast cancer cells. *Molecules* **23**(5), 1083 (2018).
2. Zhou, J., Sun, Y.-Y., Sun, M.-Y., *et al.* Prim-O-glucosylcimifugin attenuates lipopolysaccharide-induced inflammatory response in RAW 264.7 macrophages. *Pharmacogn. Mag.* **13**(51), 378-384 (2017).
3. Chen, N., Wu, Q., Chi, G., *et al.* Prime-O-glucosylcimifugin attenuates lipopolysaccharide-induced acute lung injury in mice. *Int. Immunopharmacol.* **16**(2), 139-147 (2013).
4. Wu, L.-Q., Li, Y., Li, Y.-Y., *et al.* Antinociceptive effects of prim-O-glucosylcimifugin in inflammatory nociception via reducing spinal COX-2. *Biomol. Ther. (Seoul)* **24**(4), 418-425 (2016).

#### 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