

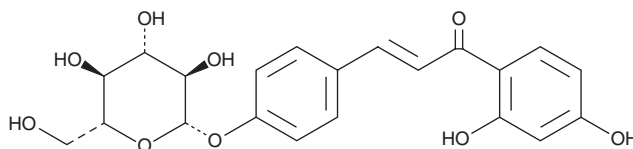
# PRODUCT INFORMATION



## Isoliquiritin

Item No. 30270

**CAS Registry No.:** 5041-81-6  
**Formal Name:** (2E)-1-(2,4-dihydroxyphenyl)-3-[4-(β-D-glucopyranosyloxy)phenyl]-2-propen-1-one  
**Synonym:** Isoliquiritoside  
**MF:** C<sub>21</sub>H<sub>22</sub>O<sub>9</sub>  
**FW:** 418.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 229, 362 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Glycyrrhiza glabra*



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

### Laboratory Procedures

Isoliquiritin is supplied as a crystalline solid. A stock solution may be made by dissolving the isoliquiritin in the solvent of choice, which should be purged with an inert gas. Isoliquiritin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of isoliquiritin in these solvents is approximately 20 mg/ml.

Isoliquiritin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, isoliquiritin should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Isoliquiritin 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

Isoliquiritin is a flavonoid that has been found in *G. uralensis* roots and has diverse biological activities.<sup>1,2,3</sup> It inhibits LPS-induced nitrite and prostaglandin E<sub>2</sub> (PGE<sub>2</sub>; Item No. 14010) production in RAW 264.7 cells when used at a concentration of 1.6 μM.<sup>1</sup> Isoliquiritin (5-500 μg/ml) inhibits tube formation by isolated rat aortic endothelial cells.<sup>2</sup> *In vivo*, isoliquiritin reduces carmine content, a marker of blood vessel formation, and pouch fluid weight in a mouse model of adjuvant-induced pouch granuloma formation (ED<sub>50</sub>s = 1.46 and 0.771 mg/kg, respectively). It increases cortical, hippocampal, and hypothalamic serotonin (5-HT) and norepinephrine levels and decreases immobility time in the forced swim and tail suspension tests in mice when administered at doses ranging from 10 to 40 mg/kg.<sup>3</sup>

### References

1. Kim, J.Y., Park, S.J., Yun, K.J., *et al.* Isoliquiritigenin isolated from the roots of *Glycyrrhiza uralensis* inhibits LPS-induced iNOS and COX-2 expression via the attenuation of NF-κB in RAW 264.7 macrophages. *Eur. J. Pharmacol.* **584**(1), 175-184 (2008).
2. Kobayashi, S., Miyamoto, T., Kimura, I., *et al.* Inhibitory effect of isoliquiritin, a compound in licorice root, on angiogenesis *in vivo* and tube formation *in vitro*. *Biol. Pharm. Bull.* **18**(10), 1382-1386 (1995).
3. Wang, W., Hu, X., Zhao, Z., *et al.* Antidepressant-like effects of liquiritin and isoliquiritin from *Glycyrrhiza uralensis* in the forced swimming test and tail suspension test in mice. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **32**(5), 1179-1184 (2008).

#### 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, 10/03/2022

#### 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