

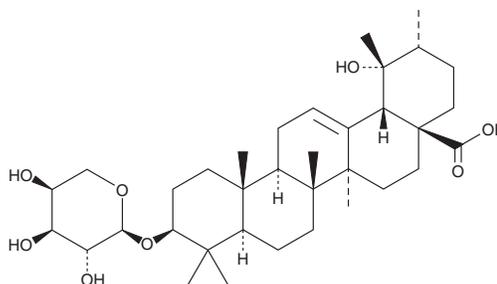
# PRODUCT INFORMATION



## Ziyuglycoside II

Item No. 40831

**CAS Registry No.:** 35286-59-0  
**Formal Name:** 3β-(α-L-arabinopyranosyloxy)-19-hydroxy-urs-12-en-28-oic acid  
**Synonym:** ZGS II  
**MF:** C<sub>35</sub>H<sub>56</sub>O<sub>8</sub>  
**FW:** 604.8  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Sanguisorba officinalis* L.



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

### Laboratory Procedures

Ziyuglycoside II is supplied as a solid. A stock solution may be made by dissolving the ziyuglycoside II in the solvent of choice, which should be purged with an inert gas. Ziyuglycoside II is slightly soluble (0.1-1 mg/ml) in ethanol and DMSO.

### Description

Ziyuglycoside II is a steroid glycoside that has been found in *S. officinalis* and has diverse biological activities.<sup>1-3</sup> It is cytotoxic to SMMC-7721, PC-9, HCT116, MCF-7, and MDA-MB-231 cancer cells when used at concentrations of 30 and 90 μg/ml.<sup>1</sup> Ziyuglycoside II (6, 12, and 18 mg/kg) increases the number of white blood cells in peripheral blood and the number of hematopoietic stem and progenitor cells (HSPCs) in a mouse model of leukopenia induced by the nitrogen mustard alkylating agent cyclophosphamide (Item No. 13849).<sup>2</sup> It also reduces small intestinal epithelial cell apoptosis and IL-1β, IL-6, and TNF-α levels, as well as improves diarrhea symptoms in a mouse model of rotavirus-induced diarrhea.<sup>3</sup>

### References

1. Liao, W., Fan, L., Zheng, Z., *et al.* Ziyuglycoside II exerts antiproliferative and antimetastasis effects on hepatocellular carcinoma cells. *Anticancer Drugs* **31(8)**, 819-827 (2020).
2. Fang, H., Xie, X., Liu, P., *et al.* Ziyuglycoside II alleviates cyclophosphamide-induced leukopenia in mice via regulation of HSPC proliferation and differentiation. *Biomed. Pharmacother.* **132:110862**, (2020).
3. Chen, X., Liu, L., Chen, W., *et al.* Ziyuglycoside II inhibits rotavirus induced diarrhea possibly via TLR4/NF-κB pathways. *Biol. Pharm. Bull.* **43(6)**, 932-937 (2020).

#### 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, 05/29/2024

#### 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