

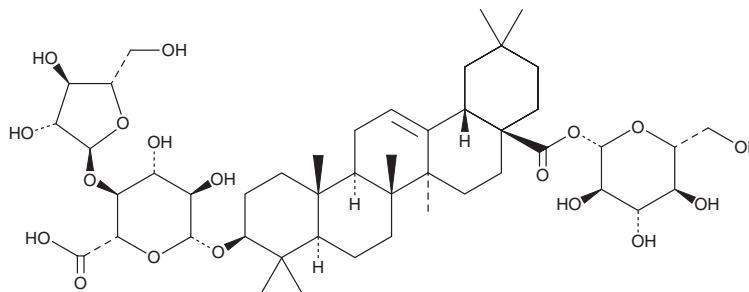
# PRODUCT INFORMATION



## Araloside A

Item No. 30219

**CAS Registry No.:** 7518-22-1  
**Formal Name:** (3β)-28-(β-D-glucopyranosyloxy)-28-oxoolean-12-en-3-yl 4-O-α-L-arabinofuranosyl-β-D-glucopyranosiduronic acid  
**Synonyms:** Chikusetsusaponin 4, Chikusetsusaponin IV, Oleanoside E  
**MF:** C<sub>47</sub>H<sub>74</sub>O<sub>18</sub>  
**FW:** 927.1  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Plant/*Aralia chinensis*



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

### Laboratory Procedures

Araloside A is supplied as a crystalline solid. Aqueous solutions of araloside A can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of araloside A in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Araloside A is a triterpenoid saponin that has been found in *A. taibaiensis* and has anti-inflammatory and gastroprotective effects.<sup>1,2</sup> It increases apoptosis of and decreases production of IL-1β and IL-6 in human MH7A rheumatoid arthritis fibroblast-like synoviocytes when used at concentrations ranging from 4 to 16 μM.<sup>1</sup> Araloside A (50 and 100 mg/kg) reduces gastric acid secretion and lesion size in rat models of hydrochloric acid and ethanol-, aspirin-, water immersion stress-, or pyloric ligation-induced ulcer formation.<sup>2</sup>

### References

- Ding, Y., Zhao, Q., and Wang, L. Pro-apoptotic and anti-Inflammatory effects of araloside A on human rheumatoid arthritis fibroblast-like synoviocytes. *Chem. Biol. Interact.* **306**, 131-137 (2019).
- Lee, E.B., Kim, O.J., Kang, S.S., et al. Araloside A, an antiulcer constituent from the root bark of *Aralia elata*. *Biol. Pharm. Bull.* **28(3)**, 523-526 (2005).

#### 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, 11/29/2022

#### 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