

PRODUCT INFORMATION



Astragaloside III

Item No. 22432

CAS Registry No.: 84687-42-3
Formal Name: (3 β ,6 α ,16 β ,20R,24S)-20,24-epoxy-6,16,25-trihydroxy-9,19-cyclolanostan-3-yl 2-O- β -D-glucopyranosyl- β -D-xylopyranoside

Synonyms: AS-III, AST-III

MF: C₄₁H₆₈O₁₄

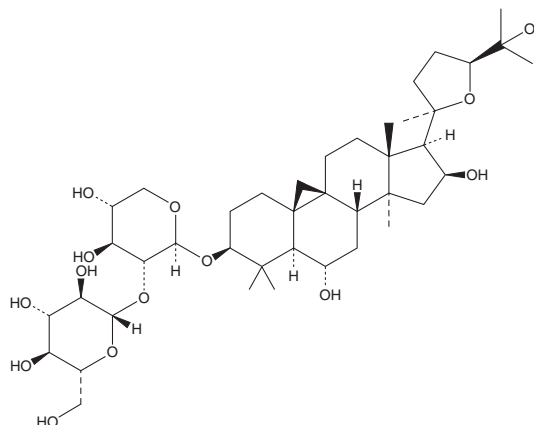
FW: 785.0

Purity: \geq 95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

Astragaloside III is supplied as a crystalline solid. A stock solution may be made by dissolving the astragaloside III in the solvent of choice, which should be purged with an inert gas. Astragaloside III is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of astragaloside III in these solvents is approximately 30 and 20 mg/ml, respectively.

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

Astragaloside III is a saponin first isolated from the dried plant roots of the genus *Astragalus*, which is used in traditional Chinese medicine.¹ It is bioavailable after oral administration, distributed mainly to the thymus and spleen, and has a half-life of approximately 2 hours.^{2,3}

References

1. Ren, S., Zhang, H., Mu, Y., *et al.* Pharmacological effects of Astragaloside IV: A literature review. *J. Tradit. Chin. Med.* **33**(3), 413-416 (2013).
2. Liu, X.-H., Guo, L., Yang, Y.-L., *et al.* Development and validation of a rapid and simple UPLC-ESI-MS method for pharmacokinetics and tissue distribution of astragaloside III in rats. *J. Chromatogr. Sci.* **54**(5), 811-818 (2016).
3. Zhai, Y., Li, P., Wang, M., *et al.* Determination of astragaloside III in rat plasma by liquid chromatography-tandem mass spectrometry and its application to a rat pharmacokinetic study. *Biomed. Chromatogr.* **30**(2), 105-110 (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, 12/15/2022

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