

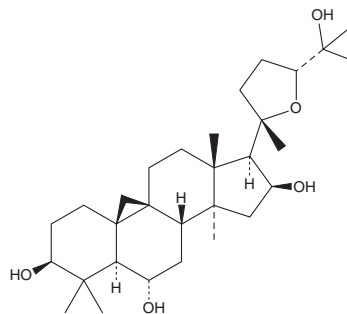
PRODUCT INFORMATION



Cycloastragenol

Item No. 11722

CAS Registry No.: 84605-18-5
Formal Name: 20,24R-epoxy-9,19-cyclolanostane-3 β ,6 α ,16 β ,25-tetrol
Synonyms: Astramembrangenin, Cyclogalegigenin
MF: C₃₀H₅₀O₅
FW: 490.7
Purity: \geq 98%
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

Cycloastragenol is supplied as a crystalline solid. A stock solution may be made by dissolving the cycloastragenol in the solvent of choice, which should be purged with an inert gas. Cycloastragenol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of cycloastragenol in these solvents is approximately 10, 3, and 25 mg/ml, respectively.

Cycloastragenol is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cycloastragenol should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Cycloastragenol has a solubility of approximately 0.25 mg/ml in a 1:8 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cycloastragenol, the aglycone derivative of astragaloside A (Item No. 16677), can be found in various species of *Astragalus* and is reported to activate telomerase.¹ It has been shown to slow telomere loss, increase replicative capacity, and improve immune function in CD8⁺ T lymphocytes from HIV-1-infected patients.²

References

1. Zhu, J., Lee, S., Ho, M.K.C., *et al.* *In vitro* intestinal absorption and first-pass intestinal and hepatic metabolism of cycloastragenol, a potent small molecule telomerase activator. *Drug Metab. Pharmacokinet.* **25**(5), 477-486 (2010).
2. Fauce, S.R., Jamieson, B.D., Chin, A.C., *et al.* Telomerase-based pharmacologic enhancement of antiviral function of human CD8⁺ T lymphocytes. *J. Immunol.* **181**(10), 7400-7406 (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, 09/30/2022

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