

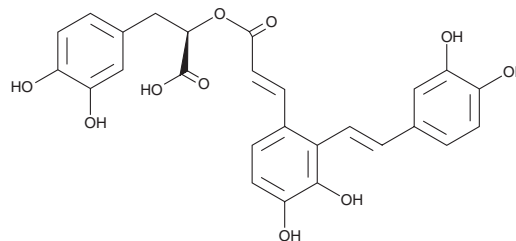
PRODUCT INFORMATION



Salvianolic Acid A

Item No. 17800

CAS Registry No.: 96574-01-5
Formal Name: αR-[[[(2E)-3-[2-[(1E)-2-(3,4-dihydroxyphenyl)ethenyl]-3,4-dihydroxyphenyl]-1-oxo-2-propen-1-yl]oxy]-3,4-dihydroxybenzenepropanoic acid
Synonyms: Dan Phenolic Acid A, (+)-Salvianolic Acid A
MF: C₂₆H₂₂O₁₀
FW: 494.5
Purity: ≥95%
UV/Vis.: λ_{max}: 289 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Salvianolic acid A is supplied as a crystalline solid. A stock solution may be made by dissolving the salvianolic acid A in the solvent of choice, which should be purged with an inert gas. Salvianolic acid A is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of salvianolic acid A in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of salvianolic acid A can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of salvianolic acid A in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Salvianolic acid A is an antioxidant and free radical scavenging compound extracted from *S. miltiorrhiza* that has been investigated for its cardioprotective and chemopreventative properties.^{1,2} It has been reported to reduce leukocyte-endothelial adherence, as well as inhibit inflammation and matrix metalloproteinases, suppress apoptosis, and reduce lipid peroxidation in damaged cardiac tissue.^{1,3}

References

1. Ho, J.H.C. and Hong, C.-Y. Salvianolic acids: Small compounds with multiple mechanisms for cardiovascular protection. *J. Biomed. Sci.* **18**, (2011).
2. Liu, G.-T., Zhang, T.-M., Wang, B., et al. Protective action of seven natural phenolic compounds against peroxidative damage to biomembranes. *Biochem. Pharmacol.* **43**, 147-152 (1992).
3. Zhao, Y., Guo, Y., and Gu, X. Salvianolic acid B, a potential chemopreventive agent, for head and neck squamous cell cancer. *J. Oncol.* **2011**, (2011).

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, 10/27/2022

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