

PRODUCT INFORMATION

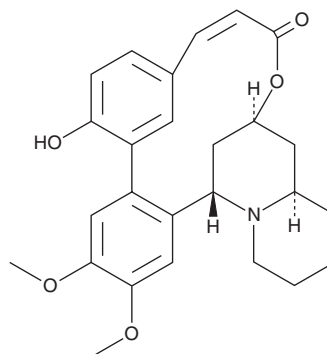


Cryogenine

Item No. 25287

CAS Registry No.: 10308-13-1
Formal Name: (4aR,6S,9Z,20S)-2,3,4,4a,5,6-hexahydro-14-hydroxy-17,18-dimethoxy-20H-6,20-methano-11,15-metheno-1H,8H-benzo[g]pyrido[2,1-d][1,5]oxaazacyclohexadecin-8-one

Synonym: NSC 272693
MF: C₂₆H₂₉NO₅
FW: 435.5
Purity: ≥98%
UV/Vis.: λ_{max}: 283 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years
Item Origin: Plant/*Heimia salicifolia*



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

Laboratory Procedures

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

Cryogenine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, cryogenine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Cryogenine has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cryogenine is an alkaloid originally isolated from *H. salicifolia* that has anti-inflammatory activity.¹ It inhibits prostaglandin synthetase (IC₅₀ = 424 μM). Cryogenine (100 mg/kg per day, p.o.) reduces paw edema and the mean arthritic index in a rat model of adjuvant-induced polyarthritis.²

References

1. Lema, W.J., Blankenship, J.W., and Malone, M.H. Prostaglandin synthetase inhibition by alkaloids of *Heimia salicifolia*. *J. Ethnopharmacol.* **15(2)**, 161-167 (1986).
2. Kosersky, D.S., Brown, J.K., and Malone, M.H. Effects of cryogenine on adjuvant-induced arthritis and serum turbidity in rats. *J. Pharm. Sci.* **62(12)**, 1965-1971 (1973).

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, 11/27/2019

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