

PRODUCT INFORMATION



Cynaropicrin

Item No. 25099

CAS Registry No.: 35730-78-0
Formal Name: 2-(hydroxymethyl)-2-propenoic acid, (3aR,4S,6aR,8S,9aR,9bR)-dodecahydro-8-hydroxy-3,6,9-tris(methylene)-2-oxoazuleno[4,5-b]furan-4-yl ester

MF: C₁₉H₂₂O₆

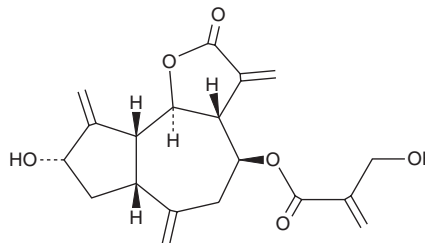
FW: 346.4

Purity: ≥98%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

Cynaropicrin is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. Cynaropicrin is miscible in these solvents.

Cynaropicrin is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of cynaropicrin should be diluted with the aqueous buffer of choice. Cynaropicrin has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Cynaropicrin is a sesquiterpene lactone originally isolated from artichoke (*C. scolymus*) that has diverse biological activities.¹⁻⁵ It inhibits the growth of SKOV3, LOX-IMVI, A549, MCF-7, HCT15, and PC-3 cancer cells (IC₅₀s = 1.1-8.7 μg/ml).¹ Cynaropicrin inhibits hepatitis C virus (HCV) replication in Huh7.5 cells with EC₅₀ values ranging from 0.4 to 1.4 μM for genotypes 1a, 1b, 2b, 3a, 4a, 5a, 6a, and 7a.² It inhibits release of TNF-α and nitric oxide (NO) from LPS-stimulated RAW264.7 cells (IC₅₀s = 8.24 and 1.1 μM, respectively) as well as LPS-induced lymphocyte proliferation (IC₅₀ = 0.9 μM).³ Cynaropicrin inhibits the growth of *T. cruzi* bloodstream trypomastigotes isolated from infected mice (EC₅₀ = 1 μg/ml).⁴ It also exhibits antifeedant activity against *S. granarius* beetles, *T. confusum* larvae, and *T. granarium* larvae.⁵

References

1. Elsebai, M.F., Mocan, A., and Atanasov, A.G. Cynaropicrin: A comprehensive research review and therapeutic potential as an anti-hepatitis C virus agent. *Front. Pharmacol.* **7:472**, (2016).
2. Elsebai, M.F., Koutsoudakis, G., Saludes, V., et al. Pan-genotypic hepatitis C virus inhibition by natural products derived from the wild egyptian artichoke. *J. Virol.* **90(4)**, 1918-1930 (2015).
3. Cho, J.Y., Baik, K.U., Jung, J.H., et al. *In vitro* anti-inflammatory effects of cynaropicrin, a sesquiterpene lactone, from *Saussurea lappa*. *Eur. J. Pharmacol.* **398(3)**, 399-407 (2000).
4. da Silva, C.F., Batista Dda, G., De Araújo, J.S., et al. Activities of psilostachyin A and cynaropicrin against *Trypanosoma cruzi* *in vitro* and *in vivo*. *Antimicrob. Agents Chemother.* **57(11)**, 5307-5314 (2013).
5. Cis, J., Nowark, G., and Kisiel, W. Antifeedant properties and chemotaxonomic implications of sesquiterpene lactones and syringin from *Rhaponticum pulchrum*. *Biochem. System. Ecol.* **34(12)**, 862-867 (2006).

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, 11/09/2018

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