

PRODUCT INFORMATION



Chaetoglobosin A

Item No. 19405

CAS Registry No.: 50335-03-0
Formal Name: (1E,4S,5E,7R,9E,11aR,14S,14aR,15S,15aR,16aS,16bR)-4,7,14,14a,15,15a,16a,16b-octahydro-7-hydroxy-14-(1H-indol-3-ylmethyl)-4,6,15,15a-tetramethyl-3H-cyclotridec[d]oxireno[f]isindole-8,11,12(13H)-trione

MF: C₃₂H₃₆N₂O₅

FW: 528.6

Purity: ≥98%

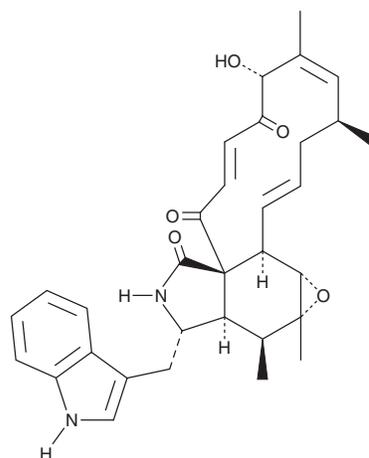
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Special Conditions: After reconstitution protect from light at -20°C

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



Laboratory Procedures

Chaetoglobosin A is supplied as a solid. A stock solution may be made by dissolving the chaetoglobosin A in the solvent of choice, which should be purged with an inert gas. Chaetoglobosin A is soluble in methanol.

Description

Chaetoglobosin A is a mycotoxic cytochalasin that was first isolated from the marine-derived endophytic fungus *C. globosum*.¹ It targets filamentous actin and demonstrates antibacterial and nematicidal effects as well as induces apoptosis in cancer cell lines.²⁻⁴

References

1. Cui, C.-M., Li, X.-M., Li, C.-S., *et al.* Cytoglobosins A-G, cytochalasins from a marine-derived endophytic fungus, *Chaetomium globosum* QEN-14. *J. Nat. Prod.* **73**(4), 729-733 (2010).
2. Yahara, I., Harada, F., Sekita, S., *et al.* Correlation between effects of 24 different cytochalasins on cellular structures and cellular events and those on actin *in vitro*. *J. Cell Biol.* **92**(1), 69-78 (1982).
3. Hu, Y., Zhang, W., Zhang, P., *et al.* Nematicidal activity of chaetoglobosin A produced by *Chaetomium globosum* NK102 against *Meloidogyne incognita*. *J. Agric. Food Chem.* **61**(1), 41-46 (2013).
4. Knudsen, P.B., Hanna, B., Ohl, S., *et al.* Chaetoglobosin A preferentially induces apoptosis in chronic lymphocytic leukemia cells by targeting the cytoskeleton. *Leukemia* **28**(6), 1289-1298 (2014)

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, 11/30/2022

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