

# PRODUCT INFORMATION



## 19-O-Acetylchaetoglobosin A

Item No. 33164

CAS Registry No.: 50939-69-0  
Formal Name: (1E,4S,5E,7R,9E,11aR,14S,14aR,15S,15aR,16aS,16bR)

-7-(acetyloxy)-3,4,7,13,14,14a,15,15a,16a,16b-decahydro-14-(1H-indol-3-ylmethyl)-4,6,15,15a-tetramethyl-12H-cyclotridec[d]oxireno[f]isoindole-8,11,12-trione

Synonym: Chaetoglobosin A Acetate

MF: C<sub>34</sub>H<sub>38</sub>N<sub>2</sub>O<sub>6</sub>

FW: 570.7

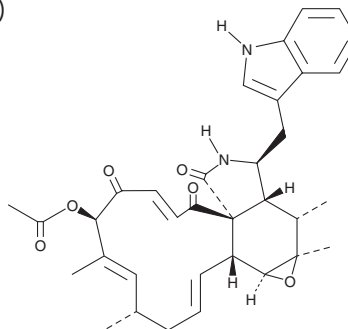
Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Fungus/*Acremonium* sp.



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

### Laboratory Procedures

19-O-Acetylchaetoglobosin A is supplied as a solid. A stock solution may be made by dissolving the 19-O-acetylchaetoglobosin A in the solvent of choice, which should be purged with an inert gas. 19-O-Acetylchaetoglobosin A is soluble in methanol and acetonitrile.

### Description

19-O-Acetylchaetoglobosin A is a fungal metabolite originally isolated from *C. globosum* that has actin polymerization inhibitory and cytotoxic activities.<sup>1,2</sup> It inhibits actin polymerization in a cell-free assay when used at a concentration of 2 μM.<sup>2</sup> 19-O-Acetylchaetoglobosin A (3.2, 10, and 32 μg/ml) is cytotoxic to HeLa cervical cancer cells.<sup>1</sup>

### References

1. Umeda, M., Ohtsubo, K., Saito, M., *et al.* Cytotoxicity of new cytochalasans from *Chaetomium globosum*. *Experientia* **31(4)**, 435-438 (1975).
2. Sekita, S., Yoshihira, K., Natori, S., *et al.* Structure-activity relationship of thirty-nine cytochalasans observed in the effects on cellular structures and cellular events and on actin polymerization *in vitro*. *J. Pharmacobiodyn.* **8(11)**, 906-916 (1985).

#### 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, 02/16/2024

#### 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