

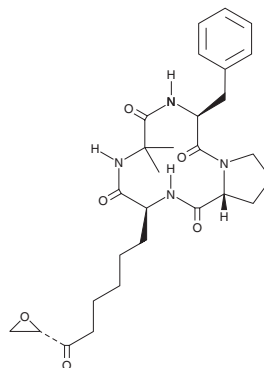
PRODUCT INFORMATION



Chlamydocin

Item No. 28468

CAS Registry No.: 53342-16-8
Formal Name: cyclo[2-methylalanyl-L-phenylalanyl-D-prolyl-(αS,2S)-α-amino-η-oxo-2-oxiraneoctanoyl]
MF: C₂₈H₃₈N₄O₆
FW: 526.6
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

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

Description

Chlamydocin is a histone deacetylase (HDAC) inhibitor that was originally isolated from *D. chlamydosporia* and has anticancer properties.¹ It is selective for HDAC1 over HDAC6 (IC₅₀ = 0.15 and 1,100 nM, respectively).² Chlamydocin increases acetylation of histone H3 and histone H4 in A2780 cells when used at concentrations ranging from 1 to 1,000 nM.¹ It inhibits growth of A2780, Malme-3M, MCF-7, HT-29, and HeLa cancer cells (IC₅₀s = 0.36, 45, 5.3, 4.3, and 14 nM, respectively). Chlamydocin increases lifespan by 10% in a P185 mouse allograft model when administered at doses ranging from 20 to 160 mg/kg.³

References

1. De Schepper, S., Bruwiere, H., Verhulst, T., *et al.* Inhibition of histone deacetylases by chlamydocin induces apoptosis and proteasome-mediated degradation of survivin. *J. Pharmacol. Exp. Ther.* **304**(2), 881-888 (2003).
2. Furumai, R., Komatsu, Y., Nishino, N., *et al.* Potent histone deacetylase inhibitors built from trichostatin A and cyclic tetrapeptide antibiotics including trapoxin. *Proc. Natl. Acad. Sci. U.S.A.* **98**(1), 87-92 (2001).
3. Stähelin, H. and Trippmacher, A. Cytostatic activity of chlamydocin, a rapidly inactivated cyclic tetrapeptide. *Eur. J. Cancer* **10**(12), 801-808 (1974).

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, 08/16/2023

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