

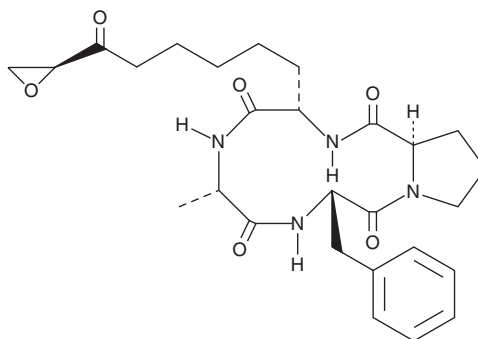
PRODUCT INFORMATION



1-Alaninechlamydocin

Item No. 28102

CAS Registry No.: 141446-96-0
Formal Name: cyclo[L-alanyl-L-phenylalanyl-D-prolyl-(α S,2S)- α -amino- η -oxo-2-oxiraneoctanoyl]
MF: $C_{27}H_{36}N_4O_6$
FW: 512.6
Purity: $\geq 95\%$
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

1-Alaninechlamydocin is supplied as a solid. A stock solution may be made by dissolving the 1-alaninechlamydocin in the solvent of choice. 1-Alaninechlamydocin is slightly soluble (0.1-1 mg/ml) in organic solvents such as acetonitrile.

Description

1-Alaninechlamydocin is a fungal metabolite originally isolated from a Great Lakes-derived *Tolypocladium* sp. and an inhibitor of histone deacetylases (HDACs).¹ It reduces total HDAC activity in HeLa cell lysates in a concentration-dependent manner. 1-Alaninechlamydocin reduces proliferation of MIA PaCa-2, PANC-1, and hTERT-HPNE cells (GI_{50} s = 5.3, 14, and 2.0 nM, respectively).

Reference

1. Du, L., Risinger, A.L., King, J.B., et al. A potent HDAC inhibitor, 1-alaninechlamydocin, from a *Tolypocladium* sp. induces G2/M cell cycle arrest and apoptosis in MIA PaCa-2 cells. *J. Nat. Prod.* **77**(7), 1753-1757 (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, 07/08/2024

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