

PRODUCT INFORMATION



Aerothionin

Item No. 18459

CAS Registry No.: 28714-26-3
Formal Name: (5S,5'S,10R,10'R)-N,N'-1,4-butanediylbis[7,9-dibromo-10-hydroxy-8-methoxy-1-oxa-2-azaspiro[4.5]deca-2,6,8-triene-3-carboxamide

MF: C₂₄H₂₆Br₄N₄O₈

FW: 818.1

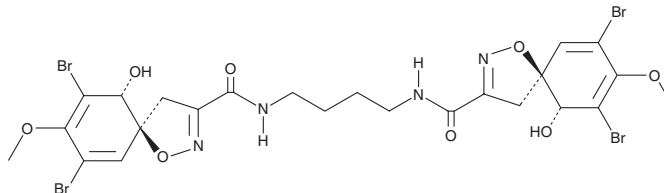
Purity: ≥97%

Supplied as: A film

Storage: -20°C

Stability: ≥4 years

Item Origin: Animal/*Suberea* sp.



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

Laboratory Procedures

Aerothionin is supplied as a film. A stock solution may be made by dissolving the aerothionin in the solvent of choice. Aerothionin is soluble in organic solvents such as ethanol and DMSO.

Description

Aerothionin is a dimeric spiroisoxazoline derived from marine sponges that demonstrates antimycobacterial activity.¹ It is cytotoxic to HeLa cells, inhibiting proliferation with an IC₅₀ value of 29 μM.¹

Reference

1. Shaala, L.A., Youssef, D.T.A., Badr, J.M., *et al.* Bioactive secondary metabolites from the red sea marine verongid sponge *Suberea* species. *Mar. Drugs* **13**, 1621-1631 (2015).

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, 12/07/2022

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