

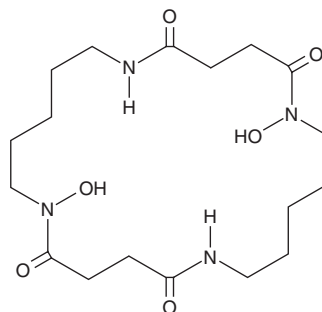
PRODUCT INFORMATION



Bisucaberin

Item No. 32890

CAS Registry No.: 112972-60-8
Formal Name: 1,12-dihydroxy-1,6,12,17-tetraazacyclodocosane-2,5,13,16-tetrone
Synonym: NSC 619796
MF: $C_{18}H_{32}N_4O_6$
FW: 400.5
Purity: $\geq 90\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Bacterium/*Alteromonas haloplankis*



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

Laboratory Procedures

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

Description

Bisucaberin is a siderophore and bacterial metabolite that has been found in *A. haloplanktis* and has anticancer activity.^{1,2} It inhibits the growth of L1210 and 1MC carcinoma cells (IC_{50} s = 9.7 and 12.7 μM , respectively) and sensitizes fibrosarcoma 1023 cells to macrophage-mediated cytolysis.

References

- Hou, Z., Raymond, K.N., O'Sullivan, B., *et al.* A preorganized siderophore: Thermodynamic and structural characterization of alcaligin and bisucaberin, microbial macrocyclic dihydroxamate chelating agents. *Inorg. Chem.* **37**(26), 6630-6637 (1998).
- Kameyama, T., Takahashi, A., Kurasawa, S., *et al.* Bisucaberin, a new siderophore, sensitizing tumor cells to macrophage-mediated cytolysis. I. Taxonomy of the producing organism, isolation and biological properties. *J. Antibiot. (Tokyo)* **40**(12), 1664-70 (1987).

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, 12/14/2022

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