
Chaetocin

(Catalog No. M64001)

Alternative Names

Description

Chaetocin, an antitumor antibiotic, is a thiodioxopiperazine natural product produced by Chaetomium species. Specific inhibitor of the lysine-specific methyltransferase SU (VAR)3-9 both in vitro and in vivo. It displays potent antimyeloma activity in IL-6-dependent myeloma cell lines. Its antimyeloma activity appears to be due to induction of oxidative stress and consequent apoptosis.

CAS Number

28097-03-2

Chemical Formula

$C_{30}H_{28}N_6O_6S_4$

Molecular Weight

696.9

Size

200 µg

Purity

≥ 98%

Solubility

Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml), dimethyl formamide or methanol. Poorly soluble in water.

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
Chaetocin	200 µg	M64001 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.