

## Chaetocin

**Catalog No:** 14051

**Format:** 200 µg

**Chemical Properties:**

MW = 696.9

$C_{30}H_{28}N_6O_6S_4$

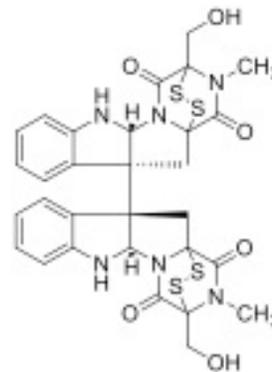
CAS 28097-03-2

Physical Properties: lyophilized solid

**Pharmacology:** Inhibits the lysine methyltransferase SU(VAR)3-9 (ref 1).

Displays anti-myeloma activity in IL-6 dependent myeloma cell lines (ref 2).

Cell permeable.



**Solubilization:** May be dissolved in DMSO (25 mg/ml); Ethanol (25 mg/ml)

**Fluorescent Properties:** N/A

**Quality Control:**

>99% (TLC); NMR (Conforms)

**References:**

1. D Greiner *et al. Nat. Chem. Biol.* 2005, 1:143
2. CR Isham *et al. Blood* 2007, 109:2579

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of arrival.