

## Mithramycin A

**Catalog No:** 14128, 14129

**Format:** 5 mg, 1 mg

**Chemical Properties:**

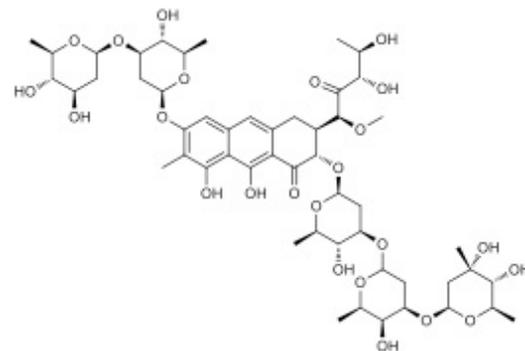
MW = 1085.2

C<sub>52</sub>H<sub>76</sub>O<sub>24</sub>

CAS 18378-89-7

Physical Properties: Yellow powder; MP = 177°C

**Names:** Mithramycin A, A 2371; Antibiotic LA 7017; Aureolic Acid; NSC 24559; PA 144; Plicamycin



**Pharmacology:** Inhibits DNA methyltransferase (ref 1). A selective Sp1 inhibitor, it binds to GC rich DNA sequences, displacing Sp1 transcription factor binding to oncogene promoters, inhibiting their expression (ref 2). Mithramycin A (at 10-200 nM) sensitizes tumor cells to TRAIL-induced apoptosis (ref 3).

**Solubilization:** May be dissolved in DMSO (20 mg/ml); or ethanol (10 mg/ml)

**Fluorescent Properties:** N/A

**Quality Control:**

>98% (HPLC); NMR (Conforms)

**References:**

1. RK Lin *et al.* *Anticancer Drugs* 2007, 18:157
2. Z Jia *et al.* *Cancer Res.* 2010, 70:1111
3. T-J Lee *et al.* *Mol. Cancer Ther.* 2006, 5:2737

**Storage and Guarantee:** Store desiccated as supplied at -20° for up to 2 years. Store solutions at -20°C for up to 1 month. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.