7-Aminoactinomycin D

Fluorescent DNA binding agent

Fluorescent, intercalating DNA-binding agent useful for identifying apoptotic cells by flow cytometry. Capable of identifying populations representing live, apoptotic, and late apoptotic/dead cells in a quick, simple, reproducible, cost-effective fashion. Inhibitor of DNA-primed RNA polymerase and DNA polymerase. Cytochemical probe. Exhibits growth-inhibitory activity against certain leukemias and sarcomas. has antibacterial properties.

Citations: 17

View Online »

Ordering Information

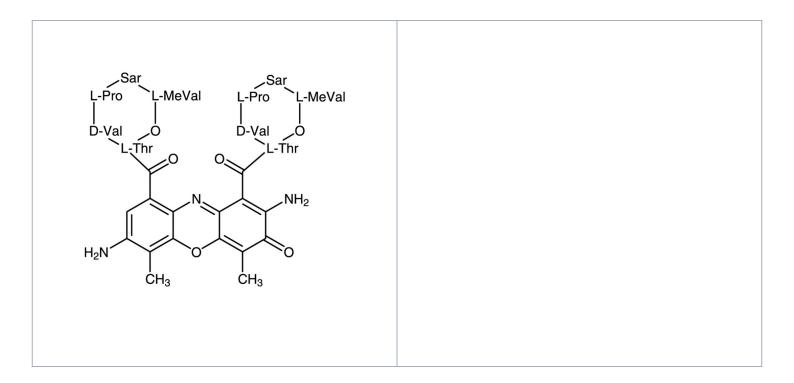
Order Online »

BML-AP400-0001

1mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

Handling Protect from light. Hygroscopic.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 7-AAD, 7AAD

Appearance Purple crystalline solid.

CAS 7240-37-1

Formula $C_{62}H_{87}N_{13}O_{16}$

MW 1270.4

Purity ≥97% (HPLC)

Solubility Soluble in DMSO (20mg/ml), methanol or 100% ethanol.

Slightly soluble in water.

Source Semisynthetic from actinomycin D (Prod. No. BML-GR300-

0005)

Last modified: May 29, 2024

