

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 »](#)

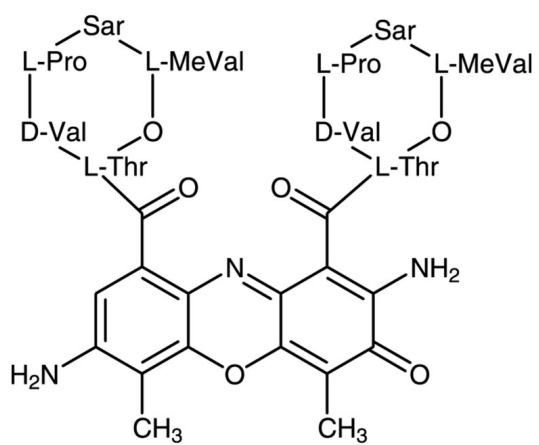
Ordering Information

[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)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)