

Aclacinomycin A

Originally identified as an anti-tumor drug, aclacinomycin A is the first described non-peptidic inhibitor showing discrete specificity for the chymotrypsin-like activity of the 20S proteasome. Aclacinomycin A has no effect on cathepsin B, it stimulates trypsin activity but inhibits the activity of both chymotrypsin and calpain.

Citations: 8

[View Online »](#)

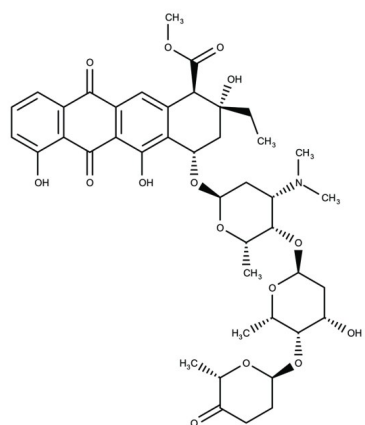
Ordering Information

[Order Online »](#)

BML-AW8655-0005	5mg
-----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Aclarubicin

Appearance Yellow powder with orange cast.

CAS 57576-44-0

Couple Target Calpain, Chymotrypsin, Proteasome, Trypsin

Couple Type Activator, Inhibitor

Formula $C_{42}H_{53}NO_{15}$

MW 811.9

Purity ≥95% (TLC)

Solubility Soluble in DMSO (25mg/ml) or dimethyl formamide (25mg/ml).



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-
be@enzolifesciences.com](mailto:info-be@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)