

(+)-Aphidicolin

Reversible inhibitor of DNA polymerase α and δ

Produced by *Cephalosporium aphidicola* and other fungi. A diterpene fungal metabolite that acts as reversible inhibitor of eukaryotic nuclear DNA replication. Useful for cell synchronization. Blocks the cell cycle at early S phase. Prolongs the half life of DNA methyltransferase. Specific inhibitor of DNA polymerase α and δ in eukaryotic cells and in some viruses of animal origin. Acts synergistically with vincristine and doxorubicin. Apoptosis inhibitor/inducer.

Citations: 21

[View Online »](#)

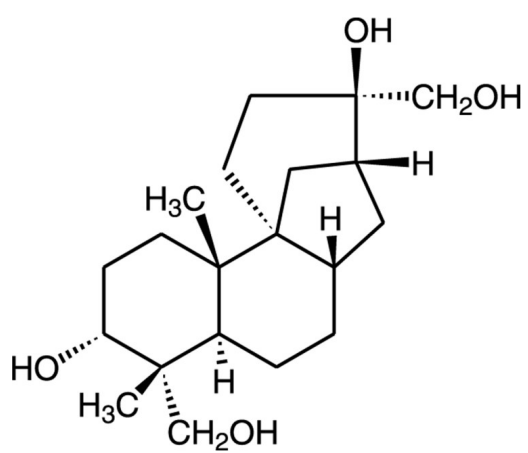
Ordering Information

[Order Online »](#)

BML-CC101-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	38966-21-1
Couple Target	DNA polymerase
Couple Type	Inhibitor
Formula	$C_{20}H_{34}O_4$
MI	14: 727
MW	338.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or ethanol (25mg/ml).

Last modified: May 29, 2024



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

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

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