

Zebularine

Dnmt inhibitor

Zebularine is a DNA methylation inhibitor which acts by forming a covalent complex with DNMT. It displays antiangiogenic activity and markedly inhibits vessel formation in different tumor models. It reactivates silenced E-cadherin and potentiates the effects of other epigenetic therapeutics such as vorinostat. Induces differentiation of mesenchymal cells into cardiomyocytes. Inhibits cytidine deaminase.

Citations: 5

[View Online »](#)

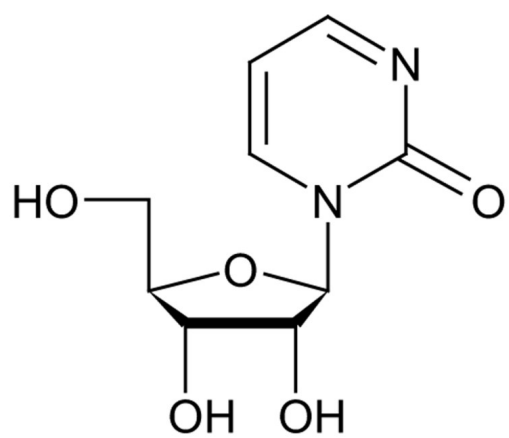
Ordering Information

[Order Online »](#)

BML-GR344-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	1-(β-D-Ribofuranosyl)-1,2-dihydropyrimidin-2-one, 2-Pyrimidone-1-β-D-ribose
Appearance	Off-white solid.
Application Notes	Can be used for <i>in vitro</i> or cell based studies.
CAS	3690-10-6
Couple Target	Dnmt
Couple Type	Inhibitor
Formula	$C_9H_{12}N_2O_5$
MW	228.2
Purity	≥98% (TLC)
Solubility	Soluble in water (>25 mg/ml) or DMSO (>25 mg/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-
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)