

5-Aza-2'- deoxycytidine

DNA methyltransferase inhibitor

Inhibitor of DNA methyltransferase. Restores mRNA and protein expression of caspase-8 and TRAIL (tumor necrosis factor-related apoptosis inducing ligand) sensitivity of resistant cell lines. Enhances apoptosis induced by HDAC (histone deacetylase) inhibitors. Reverses embryonic stem cell differentiation by demethylating differentiation-related genes.

Citations: 8

[View Online »](#)

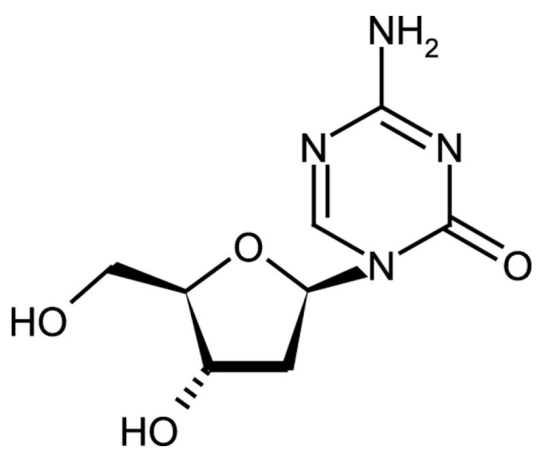
Ordering Information

[Order Online »](#)

ALX-480-096-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Packaged under inert gas. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-Aza-CdR, 5-Aza-dC, 5-Deoxy-2'-azacytidine
Appearance	White to off-white lyophilized solid.
CAS	2353-33-5
Couple Target	DNA methyltransferase
Couple Type	Inhibitor
Formula	$C_8H_{12}N_4O_4$
MW	228.2
Purity	≥98% (HPLC)
RTECS	XZ3012000
Solubility	25mg/ml soluble in DMSO or 50% acetic acid; 1mg/ml soluble in methanol. (Warming may be required to achieve complete solubilization.)

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)