

AMI-1

PRMT inhibitor

Potent, cell permeable and specific inhibitor of protein arginine N-methyltransferases (PRMTs) (IC_{50} =8.8 μ M for human PRMT1 and 3.0 μ M for yeast Hmt1p). Does not compete for the AdoMet (S-adenosyl-L-methionine; SAM) binding site. Exhibits only minimal effects on lysine N-methyltransferases. Suppresses the transcriptional coactivator activity of PRMT1 and PRMT4. Inhibitor of HIV-1 RT polymerase (IC_{50} =5.0 μ M).

Citations: 3

[View Online »](#)

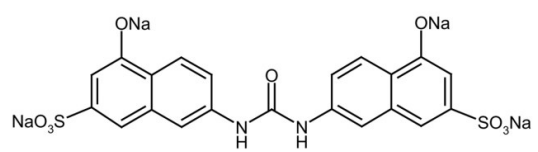
Ordering Information

[Order Online »](#)

ALX-270-440-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Hygroscopic. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Brown solid.
CAS	20324-87-2
Couple Target	PRMT
Couple Type	Inhibitor
Formula	$C_{21}H_{12}N_2O_9S_2Na_4$
Identity	Determined by NMR.
MW	592.4
Purity	≥98% (HPLC)
Solubility	Soluble in water, 100% ethanol or DMSO.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)