

# Sinefungin

## Methyltransferase inhibitor

Strong competitive inhibitor of methyltransferases which uses S-adenosyl-1-methionine as the methyl group donor to yield methylated products such as 5-methylcytosine or N<sup>6</sup>-methyl adenosine on DNA and RNA. In addition, sinefungin is involved in a number of physiological processes.

Citations: 11

[View Online »](#)

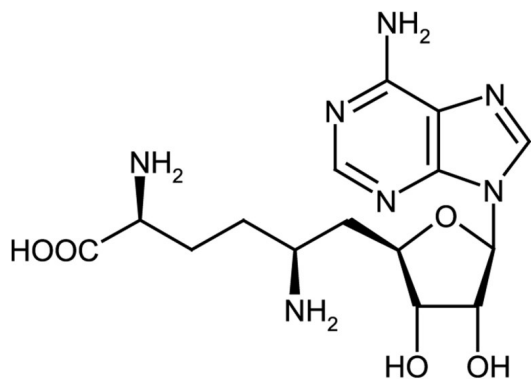
## Ordering Information

[Order Online »](#)

ALX-380-070-M001	1mg
ALX-380-070-M005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep under nitrogen. Packaged under nitrogen. Protect from light. Very hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	6,9-Diamino-1-(6-amino-9H-purin-9-yl)-1,5,6,7,8,9-hexadeoxy-D-glycero- $\alpha$ -L-talo-decafuranuronic acid, A-9145, Adenosylornithine
Appearance	White to yellow powder.
CAS	58944-73-3
Couple Target	Methyltransferase
Couple Type	Inhibitor
Formula	$C_{15}H_{23}N_7O_5$
MI	14: 8544
MW	381.4
Purity	≥95% (HPLC)
RTECS	HE3140000
Solubility	Soluble in water.

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