

# Decoyinine

## GMP synthase inhibitor

Decoyinine is a non-competitive and reversible inhibitor of GMP synthase. GMPS was shown to be a positive regulator of invasiveness in melanoma cells and decoyinine reduced invasion of melanoma cells by 30%. Decoyinine reduces intracellular levels of GMP, GDP and GTP. Numerous studies of decoyinine in bacteria have been performed showing that guanylates play key roles in bacterial growth, cell wall biosynthesis, sporulation, and pathogenicity.

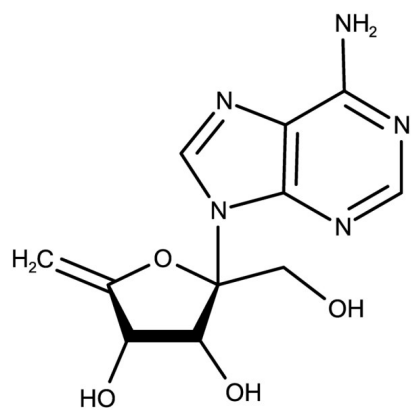
### Ordering Information

[Order Online »](#)

ENZ-CHM470-0005	5mg
-----------------	-----

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 4',5'-Didehydro-1'-C-(hydroxymethyl)-adenosine

**Appearance** Off-white solid.

**CAS** 2004-04-8

**Couple Target** GMP synthase

**Couple Type** Inhibitor

**Formula**  $C_{11}H_{13}N_5O_4$

**Identity** Determined by NMR.

**MW** 279.26

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO (50 mg/ml) or water (25 mg/ml).



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)