

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.

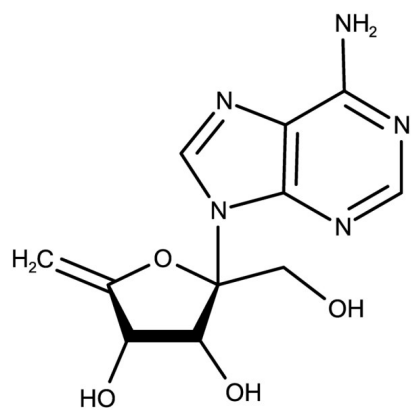
Ordering Information

[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).

Last modified: February 27, 2025



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)