

# MANT-5'-dAMP

Fluorescent analogue of 2'-deoxy-5'-AMP

*BLG-D081-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

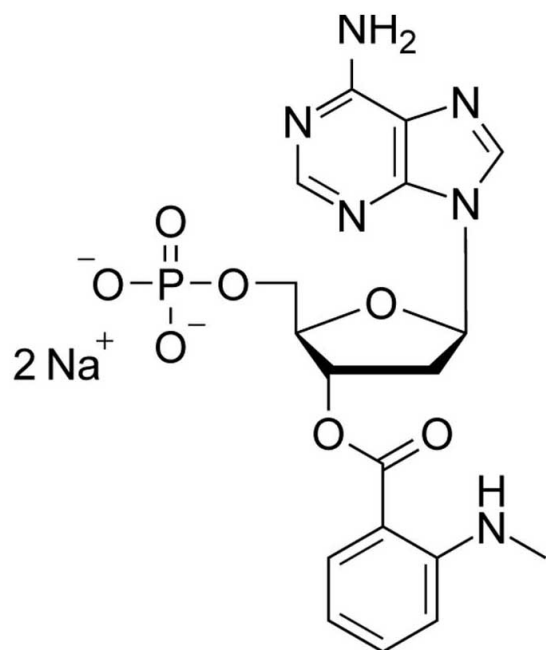
## Ordering Information

[Order Online »](#)

BLG-D081-05	5 $\mu$ mol
BLG-D081-25	5x5 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	2'-Deoxy-3'-O-(N'-methylantraniloyl)adenosine-5'-O-monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{18}H_{21}N_6O_7P$ (free acid)
MW	464.4
Purity	> 97% HPLC
Quantity	5µmol≈2.3mg
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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