

# 2'-Biotin-16-c-diAMP

Biotinylated analogue of the bacterial second messenger c-diAMP

Analogue of c-diAMP with biotin conjugate most probably suitable as tracer in immunoassays.

*BLG-B106-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

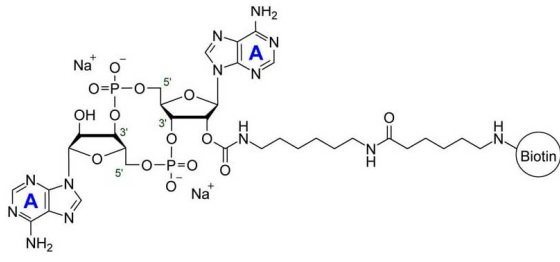
## Ordering Information

[Order Online »](#)

BLG-B106-001	0.1µmol
BLG-B106-005	5x0.1µmol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 2'-O-(6-[Biotinyl]aminohexylcarbamoyl)cyclic diadenosine monophosphate . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{43}H_{63}N_{15}O_{16}P_2S$  (free acid)

**MW** 1140.1 (free acid)

**Purity** > 95% HPLC

**Quantity** 0.1µmol≈~0.1mg

**Solubility** Soluble in water or in aqueous buffers (>1.8mM).

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)