

# 2'-AHC-c-diAMP

## Functionalized c-diAMP for labelling or immobilization

Analogue of the bacterial second messenger c-diAMP (Cat. No. BLG-C088) which is suitable as a ligand in affinity chromatography or for coupling of various labelling structures including fluorophores.

*BLG-A182-005 (5 x 0.1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

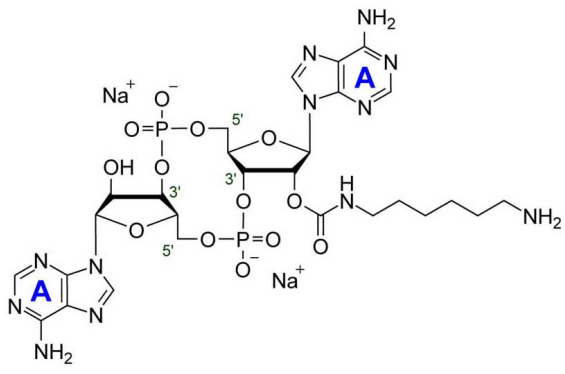
### Ordering Information

[Order Online »](#)

|              |                 |
|--------------|-----------------|
| BLG-A182-001 | 0.1 $\mu$ mol   |
| BLG-A182-005 | 5x0.1 $\mu$ mol |

### 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-Aminohexylcarbamoyl)-cyclic diadenosine monophosphate . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{27}H_{38}N_{12}O_{13}P_2$

**MW** 800.6 (free acid)

**Purity** > 95% HPLC

**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)