

# 5'-AHC-2',3'-cAMP

Functionalized 2',3'-cAMP for immobilization or conjugation with markers and dyes

Functionalized analogue of 2',3'-cAMP (Biolog Cat. No. A 307) which is suitable as a ligand in affinity chromatography or for coupling of various labelling structures including fluorophores.

*BLG-A327-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

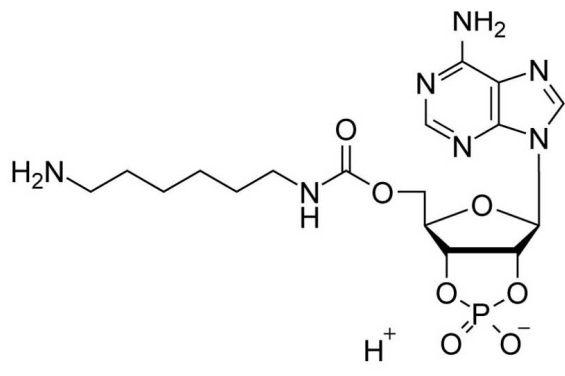
## Ordering Information

[Order Online »](#)

BLG-A327-05	5 $\mu$ mol
BLG-A327-25	5x5 $\mu$ 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** 5'-O-(6-Aminohexylcarbamoyl)adenosine-2',3'-cyclic monophosphate (free acid)

**Couple Type** Modified nucleotides

**Formula**  $C_{17}H_{26}N_7O_7P$

**MW** 471.4

**Purity** > 97% HPLC

**Quantity** 5µmol are equivalent to ~2.3 mg

**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-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)