

# 2'-NH<sub>2</sub>-cAMP

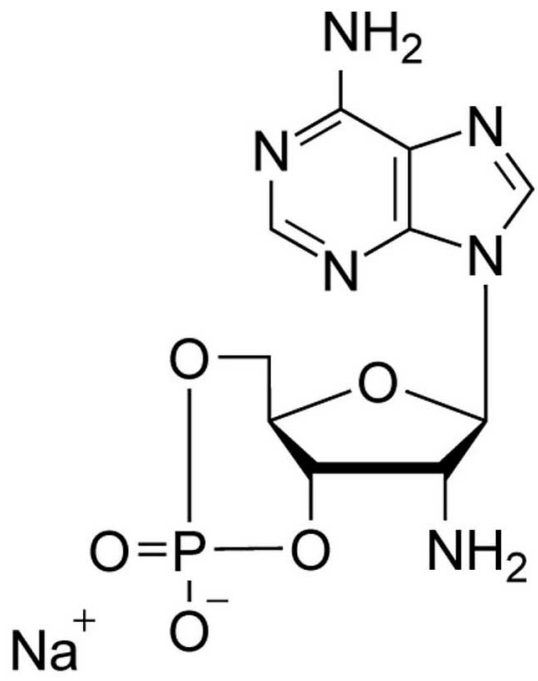
2'-NH<sub>2</sub>-cAMP has a reactive amino group suitable for modification with fluorophores and other markers or spacers and can be used as ligand for affinity chromatography of cAMP and cGMP binding proteins that do not require an intact 2'-OH group.

*BLG-A088-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

## Ordering Information [Order Online »](#)

BLG-A088-05	5µmol
BLG-A088-25	5x5µmol

## 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'-Amino-2'-deoxyadenosine-3',5'-cyclic monophosphate . sodium salt

**CAS** 85819-83-6

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{12}N_6O_5P \cdot Na$

**MW** 350.2

**Purity** > 98% HPLC

**Quantity** 5µmol≈1.8mg

**Solubility** Soluble in water.

**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)