

# 6-AE-8-N<sub>3</sub>-cAMP

Precursor analogue for photoaffinity labeling

*BLG-A101-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada.*

*Please [contact us](#) for available options.*

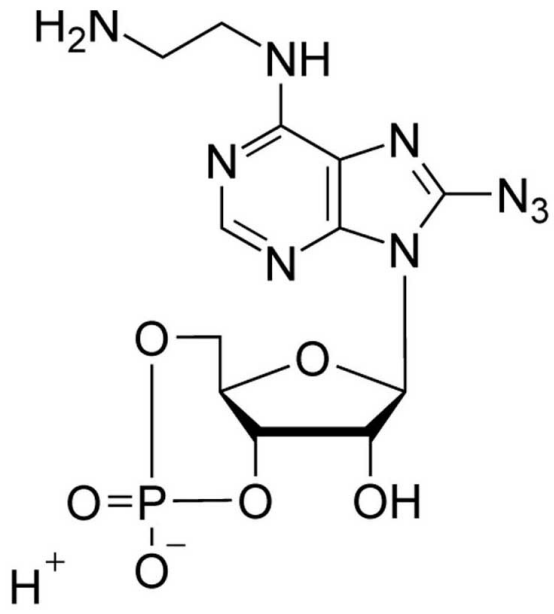
## Ordering Information

[Order Online »](#)

BLG-A101-01	1µmol
BLG-A101-05	5x1µmol

## 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** N6-(2-Aminoethyl)-8-azidoadenosine-3',5'-cyclic monophosphate

**Couple Type** Modified nucleotides

**Formula**  $C_{12}H_{16}N_9O_6P$

**MW** 413.3

**Purity** > 97% HPLC

**Quantity** 1µmol≈0.4mg

**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)