

# 2'-AEC-cAMP-Agarose

cyclic AMP immobilized on agarose gel

2'-AEC-cAMP-Agarose is a gel for affinity chromatography of various cyclic nucleotide-responsive proteins, especially those which tolerate modification of the ribose 2'-hydroxy group such as some phosphodiesterases. This type of gel is also offered with a longer spacer (2'-AHC-cAMP-Agarose, Prod No. BLG-A058) and as a PDE-resistant version (Sp-2'-AEC-cAMPS-Agarose, Prod. No. BLG-A122). The free ligand 2'-AEC-cAMP/2'-EDA-cAMP (Prod. No. BLG-A098) is available as well. For immobilized cGMP or cAMP with other spacer types, attachment positions, different immobilization techniques or on other gels please inquire.

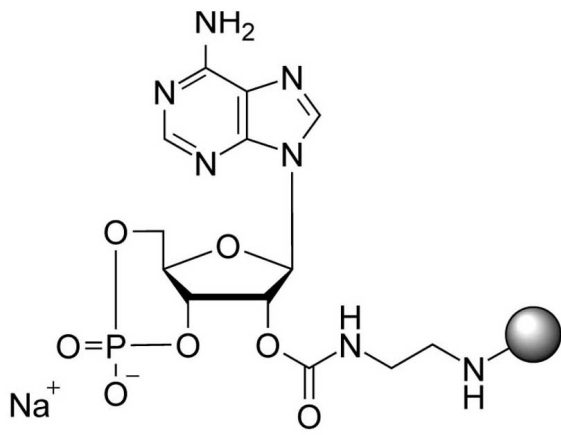
## Ordering Information

[Order Online »](#)

<b>BLG-A112-25</b>	2.5ml
<b>BLG-A112-60</b>	6ml

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 2'-O-(6-Aminoethylcarbamoyl)adenosine-3',5'-cyclic monophosphate; immobilized on agarose gel

**Couple Type** Modified nucleotides

**Formulation** Suspension in 30mM Na<sub>2</sub>HPO<sub>4</sub> buffer (pH 7), containing 0.1 % sodium azide as preservative. Ligand density: approximately 6 µmol/ml of settled gel. UV: λ<sub>max</sub> 258nm/suspension in glycol.

**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)