

# 2'-AHC-cGMP- Agarose

Cyclic GMP immobilized on agarose gel

The protein kinase G activator cGMP immobilized on agarose by an aminohexylcarbamoyl spacer attached to the ribose 2' position of the ligand. The gel can be used for affinity chromatography of various cyclic nucleotide-responsive proteins, especially those which tolerate modification of the 2' ribose hydroxyl group, such as some phosphodiesterases. This type of gel is also available with immobilized PDE-resistant Rp-cGMPS (Rp-2'-AHC-cGMPS-Agarose (Prod. No. [BLG-A051](#))) and Sp-cGMPS (Sp-2'-AHC-cGMPS-Agarose (Prod. No. [BLG-A052](#))).

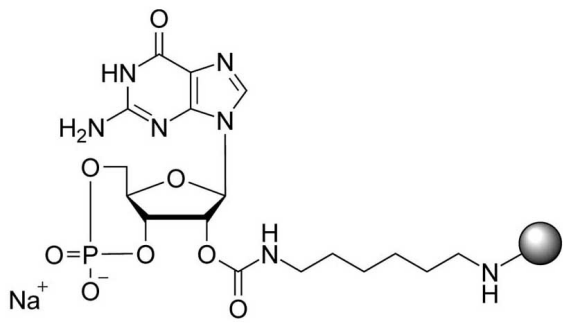
## Ordering Information

[Order Online »](#)

BLG-A059-25	2.5ml
BLG-A059-60	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'-(6-Aminohexylcarbamoyl)guanosine-3',5'-cyclic monophosphate; immobilized on agarose gel

**Couple Type** Modified nucleotides

**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)