

# 2'-AHC-cGMP

Functionalized cGMP for immobilization or conjugation with markers and dyes

Cyclic GMP analogue suitable for immobilization as affinity ligand (e.g. for phosphodiesterases ) or for coupling of various labeling structures including fluorophores.

This structure is also offered as ligand immobilized to agarose (Prod. No. [BLG-A059](#)).

*BLG-A048-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

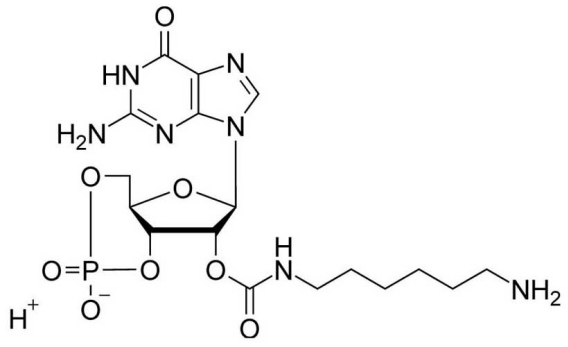
## Ordering Information

[Order Online »](#)

BLG-A048-05	5 $\mu$ mol
BLG-A048-25	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** 2'-(6-Aminohexylcarbamoyl)guanosine-3',5'-cyclic monophosphate (free acid)

**CAS** 1262749-60-9

**Couple Type** Modified nucleotides

**Formula**  $C_{17}H_{26}N_7O_8P$

**MW** 487.4

**Purity** > 98% HPLC

**Quantity** 5µmol≈2.4mg

**Solubility** Soluble in water or buffer.

**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)