

2'-AHC-c-diGMP- Agarose

2'-AHC-c-diGMP immobilized on agarose gel

2'-AHC-c-diGMP-Agarose is suitable for affinity chromatography of c-diGMP-responsive proteins. c-diGMP (Prod. No. [BLG-C057](#)) was found to be a near ubiquitous signalling molecule in bacteria involved in a wide variety of physiological processes such as biofilm formation.

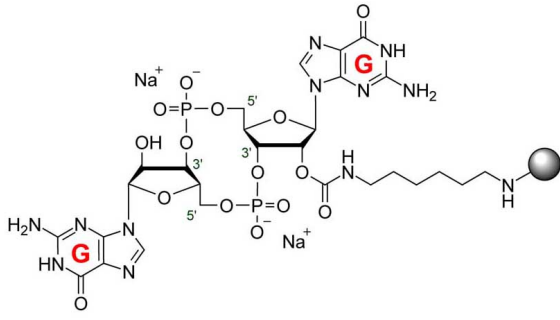
Ordering Information

[Order Online »](#)

BLG-A153-06	0.6ml
BLG-A153-25	2.5ml
BLG-A153-60	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'-O-(6-Aminohexylcarbamoyl)-cyclic diguanosine monophosphate, immobilized on agarose

Couple Type Modified nucleotides

Formulation Suspension in 30mM NaH₂PO₄ buffer (pH 7), containing 1 % sodium azide as preservative. Ligand density: approximately 1 µmol/ml of settled gel. UV: λ_{max} 253 nm/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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com