

2'-AHC-c-diAMP- Agarose

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

The second messenger c-diAMP (Cat. No. BLG-C088) immobilized on agarose by an aminohexylcarbamoyl spacer attached to one of the two ribose 2' positions of the ligand. The gel can be used for affinity chromatography of various c-diAMP-responsive proteins. Normally all affinity gels are supplied in pre-packed columns. The free beads without a column are available upon request.

Ordering Information

[Order Online »](#)

BLG-A183-06	0.6ml
BLG-A183-25	2.5ml
BLG-A183-60	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 diadenosine monophosphate, immobilized on agarose

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

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