

5'-AHC-2',3'-cAMP- Agarose

2',3'-cAMP immobilized on agarose gel

The signalling nucleotide 2',3'-cAMP (Biolog Cat. No. A 307) immobilized on agarose by an aminohexylcarbonyl spacer attached to the 5' ribose position of the ligand. The gel may be useful for affinity chromatography of 2',3'-cAMP-responsive proteins.

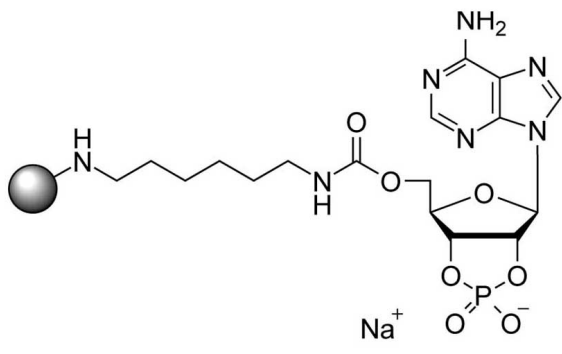
Ordering Information

[Order Online »](#)

BLG-A329-06	0.6ml
BLG-A329-25	2.5ml
BLG-A329-60	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 5'-O-(6-Aminohexylcarbamoyl)adenosine-2',3'-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

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