

# 2-AHA-cAMP-Agarose

Cyclic AMP immobilized on agarose gel

Cyclic AMP immobilized on agarose by an aminohexylamino spacer attached to position 2 of the ligand. The gel can be used for affinity chromatography of various cyclic nucleotide-responsive proteins such as protein kinases, phosphodiesterases and others. The ligand is offered separately as well (Prod. No. [BLG-A053](#)).

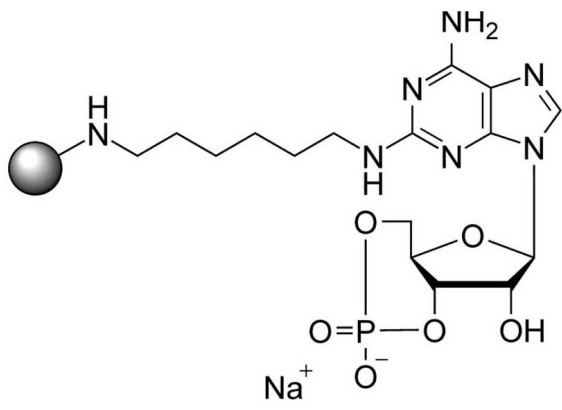
## Ordering Information

[Order Online »](#)

BLG-A054-06	0.6ml
BLG-A054-25	2.5ml
BLG-A054-60	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-Aminohexyl)aminoadenosine-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)