

# Sp-8-AHA-cAMPS- Agarose

## Immobilized PDE-resistant PKA agonist

The protein kinase A activator Sp-cAMPS immobilized on agarose by an aminohexylamino spacer attached to position 8 of the ligand. Free Sp-cAMPS is only very slowly metabolized by corresponding phosphodiesterases and thus represents an extraordinary stable ligand. In addition, its reduced binding affinity towards protein kinase A compared to cAMP should result in milder desorption conditions in affinity chromatography. The gel can be used for affinity chromatography of various cyclic nucleotide-responsive proteins such as protein kinases, phosphodiesterases and others. This ligand is also available with a shorter spacer arm (Sp-8-AEA-cAMPS-Agarose (Prod. No. [BLG-A008](#))).

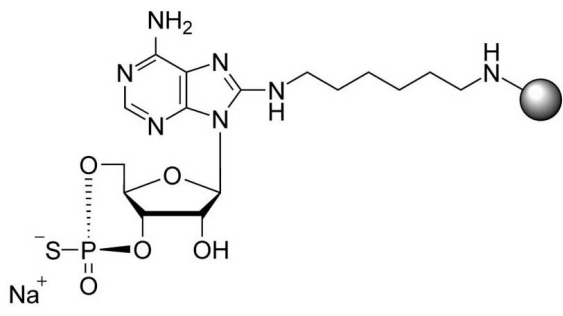
### Ordering Information

[Order Online »](#)

<b>BLG-A013-06</b>	0.6ml
<b>BLG-A013-25</b>	2.5ml
<b>BLG-A013-60</b>	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** 8-(6-Aminohexyl)aminoadenosine-3',5'-cyclic monophosphorothioate, Sp-isomer; immobilized on agarose gel

**Couple Type** Modified nucleotides

**Formulation** Liquid. 0.05M Na<sub>2</sub>HPO<sub>4</sub> containing 0.1% sodium azide.

**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)