

# Sp-6-AH-cAMPS- Agarose

PDE-resistant Sp-cAMPS immobilized on agarose gel

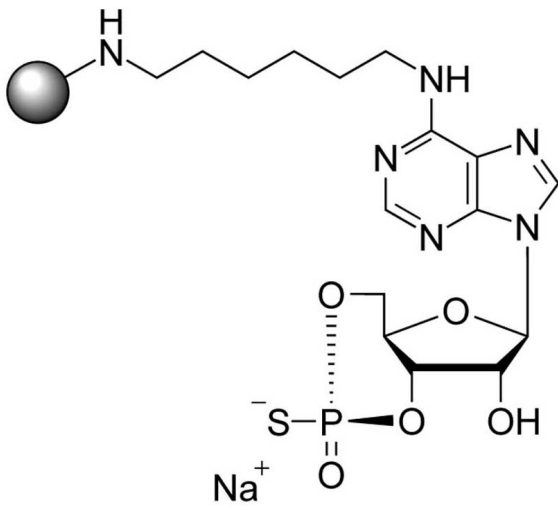
## Ordering Information

[Order Online »](#)

BLG-A108-06	0.6ml
BLG-A108-25	2.5ml
BLG-A108-60	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** N6-(6-Aminohexyl)adenosine-3',5'-cyclic monophosphorothioate, Sp-isomer; immobilized on agarose gel

**Couple Type** Modified nucleotides



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)