

# Sp-8-AHA-cAMP

**PDE-resistant Sp-cAMPS functionalized for immobilization or conjugation with markers and dyes**

PDE-stable cAMP analog suitable for immobilization as affinity ligand (e.g. for purification of phosphodiesterase) or for coupling of various labeling structures including fluorophores.

*BLG-A071-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

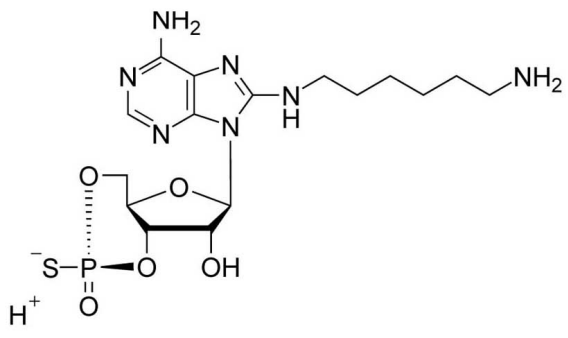
## Ordering Information

[Order Online »](#)

BLG-A071-05	5 $\mu$ mol
BLG-A071-25	5x5 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 8-(6-Aminohexyl)aminoadenosine-3',5'-cyclic monophosphate, Sp-isomer

**CAS** 214272-03-4

**Couple Type** Modified nucleotides

**Formula**  $C_{16}H_{26}N_7O_5PS$

**MW** 459.5

**Purity** > 98% HPLC

**Quantity** 5µmol≈2.3mg

**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)