

# Sp-8-OH-cAMPS

Polar analogue of the PDE-resistant PKA agonist Sp-cAMPS

Polar, membrane impermeant analogue of the protein kinase A activator Sp-cAMPS for the study of extracellular cAMP receptors and for prevention of membrane re-permeation during pipette application inside the cell. Highly resistant against mammalian cyclic nucleotide-dependent phosphodiesterases.

*BLG-H005-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

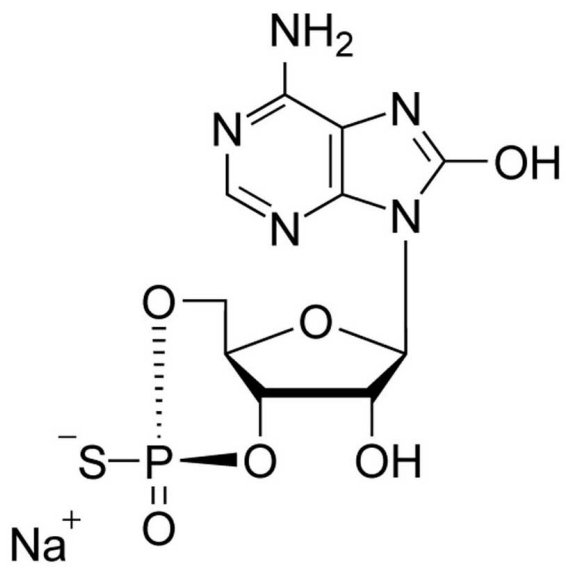
## Ordering Information

[Order Online »](#)

BLG-H005-05	5µmol
BLG-H005-25	5x5µmol

## 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-Hydroxyadenosine-3',5'-cyclic monophosphorothioate . sodium salt, Sp-isomer

**CAS** 749843-87-6

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{11}N_5O_6PS \cdot Na$

**MW** 383.3

**Purity** > 99% HPLC

**Quantity** 5µmol≈1.9mg

**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)