

# Rp-8-Br-cAMPS

Lipophilic analogue of the cyclic AMP antagonist Rp-cAMPS

Rp-8-Br-cAMPS is a novel combination of the well accepted protein kinase inhibitor Rp-cAMPS with the widely used 8-bromo cyclic AMP resulting in a membrane permeant cyclic AMP antagonist which is not metabolized by mammalian cyclic nucleotide phosphodiesterases.

Rp-8-Br-cAMPS is about 1.5 times more lipophilic compared to 8-Br-cAMP, and 2 times and 3 times more compared to Rp-cAMPS and cAMP respectively.

In contrast to common ATP-site inhibitors or to its parent compound Rp-cAMPS, Rp-8-Br-cAMPS discriminates between both isozymes of protein kinase A, preferring type I<sup>13</sup> and thus provides additional selectivity. By occupying cAMP binding sites Rp-8-Br-cAMPS prevents the kinase holoenzyme from dissociation and thus from activation.

Due to this working principle preincubation of the inhibitor prior to the activation step is necessary for optimal results.

*BLG-B001-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

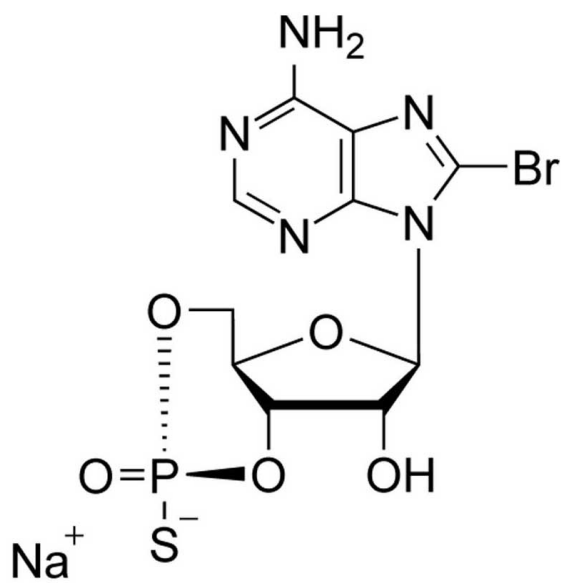
## Ordering Information

[Order Online »](#)

BLG-B001-05	5µmol
BLG-B001-25	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-Bromoadenosine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

**Appearance** Lyophilized or crystallized sodium salt.

**CAS** 129735-00-8

**Couple Type** Inhibitor, Modified nucleotides

**Formula**  $C_{10}H_{10}BrN_5O_5PS \cdot Na$

**MW** 446.2

**Purity** > 99% HPLC

**Quantity** 5µmol≈2.2mg

**Reconstitution** Rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

**Solubility** Soluble in water or buffer.

**Technical Info / Product** For the Original Manufacturer's data sheet please [click here](#).

### Notes



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)