

# Sp-8-Br-2'-O-Me-cAMPS

**PDE-resistant, membrane-permeant Epac agonist**

Sp-8-Br-2'-O-Me-cAMPS is a hydrolysis-resistant form of 8-Br-2'-O-Me-cAMP (Prod. No. BLG-B022) which is a potent membrane-permeant stimulator of the exchange factors directly activated by cAMP (Epac or cAMP-GEF).

Since a free 2'-ribose hydroxyl group in cyclic AMP is essential for stimulation of cAMP-dependent protein kinase (PKA), the methylated structure of Sp-8-Br-2'-O-Me-cAMPS is an extremely poor PKA activator and allows for specific discrimination between both signalling pathways. On the other hand, potent activators of PKA carrying a modified 6 position at the adenine nucleobase can be used as Epac-negative controls. N<sup>6</sup>-modified cyclic AMP analogues such as N<sup>6</sup>-Benzoyl-cAMP (Prod. No. BLG-B009) or N<sup>6</sup>-Phenyl-cAMP (Prod. No. BLG-P006) are specific PKA agonists, but show only neglectable agonistic properties on Epac.

*BLG-B031-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

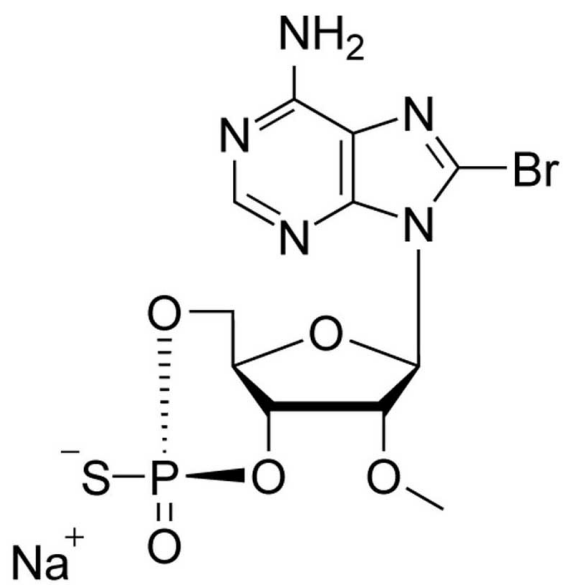
## Ordering Information

[Order Online »](#)

BLG-B031-01	1µmol
BLG-B031-05	5x1µmol

## 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-Bromo-2'-O-methyladenosine-3',5'-cyclic monophosphorothioate, Sp-isomer . sodium salt

**CAS** 634208-34-7

**Couple Type** Modified nucleotides

**Formula**  $C_{11}H_{12}BrN_5O_5PS \cdot Na$

**MW** 460.2

**Purity** > 98% HPLC

**Quantity** 1µmol≈0.5 mg

**Solubility** Soluble in water (≥10mM).

**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)