

# Sp-2'-O-MB-cAMPS

Metabolically activatable form of the PDE-resistant PKA agonist Sp-cAMPS

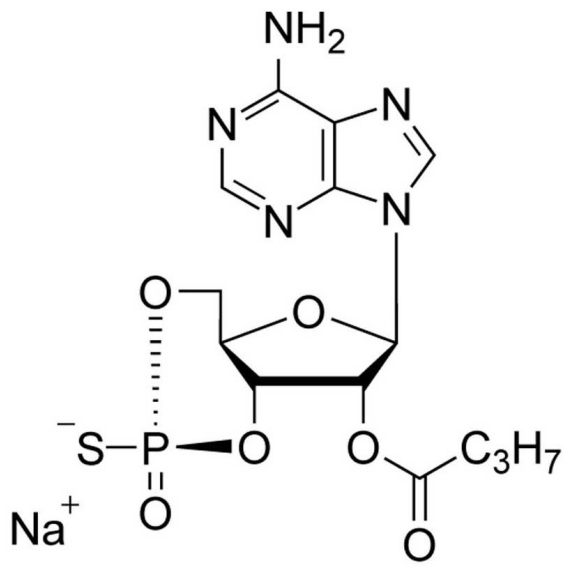
Sp-2'-O-MB-cAMPS is a lipophilic precursor of Sp-cAMPS, a PDE-resistant activator of protein kinase A. During metabolic activation by esterases Sp-cAMPS and butyrate are released. Significantly more lipophilic and membrane permeant compared to Sp-cAMPS.

*BLG-M005-25 (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-M005-05 | 5 $\mu$ mol   |
| BLG-M005-25 | 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** 2'-O-Monobutyryladenine-3',5'-cyclic monophosphorothioate . sodium salt, Sp-isomer

**CAS** 152218-23-0

**Couple Type** Modified nucleotides

**Formula**  $C_{14}H_{17}N_5O_6PS \cdot Na$

**MW** 437.4

**Purity** > 97% HPLC

**Quantity** 5µmol≈2.2mg

**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)