

# 8-PIP-cAMP

## Site-selective PKA agonist

Site selective cyclic AMP analogue with high selectivity for site A of protein kinase A type I and for site B of type II. Acts synergistically together with analogues having opposite site-selectivity. Combination with e.g. 8-HA-cAMP (Prod. No. [BLG-H006](#)) preferentially activates type I, while e.g. with 6-MBC-cAMP (Prod. No. [BLG-M012](#)) selects type II. Increased metabolic stability and membrane permeability compared to cAMP.

*BLG-P002-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

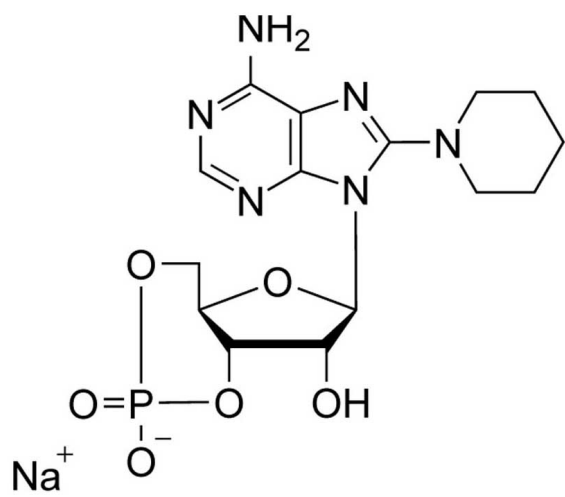
## Ordering Information

[Order Online »](#)

BLG-P002-10	10µmol
BLG-P002-50	5x10µmol

## 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-Piperidinoadenosine-3',5'-cyclic monophosphate . sodium salt

**CAS** 31357-06-9

**Couple Type** Modified nucleotides

**Formula**  $C_{15}H_{20}N_6O_6P \cdot Na$

**MW** 434.3

**Purity** > 98% HPLC

**Quantity** 10µmol≈4.3mg

**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)