

# 6-Bnz-cAMP

## Membrane-permeant, site-selective PKA agonist

Site selective and membrane permeant activator of cAMP-dependent protein kinase. Increased metabolic stability against PDE compared to cAMP. Selects site A of both cAK isozymes and thus can be used for synergistic activation of PKA type II if combined with site B selective analogues such as Sp-8-CPT-cAMPS (Prod. No. [BLG-C012](#)) or 8-piperidino-cAMP (Prod. No. [BLG-P002](#)). For preferential stimulation of cAK type I a combination with the site B selective analogue 8-HA-cAMP (Prod. No. [BLG-H006](#)) or 8-AHA-cAMP (Prod. No. [BLG-A011](#)) can be used.

*BLG-B009-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

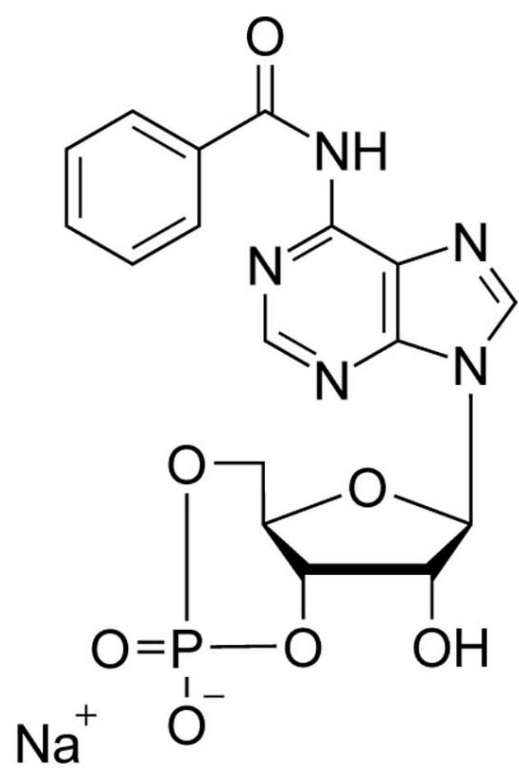
### Ordering Information

[Order Online »](#)

BLG-B009-50	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 N6-Benzoyladenine-3',5'-cyclic monophosphate . sodium salt

CAS 30275-80-0

Couple Type Activator, Modified nucleotides

Formula  $C_{17}H_{15}N_5O_7P \cdot Na$

MW 455.3

Purity > 98% HPLC

Quantity 10µmol≈4.6mg

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)