

# 6-Bnz-cAMP-AM

The acetoxymethyl group of 6-Bnz-cAMP-AM masks the charged polar phosphate and thus makes the molecule highly membrane-permeant. Inside the cell esterases release the more polar 6-Bnz-cAMP which is a potent, site-selective activator of cAMP-dependent protein kinase. 6-Bnz-cAMP does not activate Epac and thus can be used as an Epac-negative control.

Phosphate tris(acetoxymethyl)ester,  $\text{PO}_4\text{-AM}_3$  (Prod. No. [BLG-P030](#)), is recommended as control reagent in 6-Bnz-cAMP-AM applications to test for side effects of enzymatically released acetic acid and formaldehyde, two metabolites with potential biological functions.

*BLG-B079-05 (5 x 1  $\mu\text{mol}$  pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

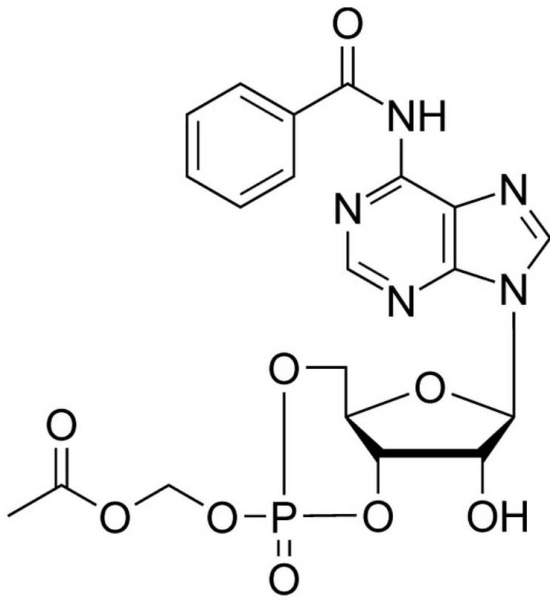
## Ordering Information

[Order Online »](#)

BLG-B079-01	1 $\mu\text{mol}$
BLG-B079-05	5x1 $\mu\text{mol}$

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** Aqueous solutions are rather labile and should be freshly prepared immediately before use. Stock solutions in anhydrous DMSO or DMF should be relatively stable when stored frozen at -20°C to -80°C.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N6-Benzoyladenine-3',5'-cyclic monophosphate, acetoxymethyl ester

**Couple Type** Modified nucleotides

**Formula**  $C_{20}H_{20}N_5O_9P$

**MW** 505.4

**Purity** > 97% HPLC for mixture of isomers

**Quantity** 1µmol≈0.5 mg

**Solubility** Due to its rather high lipophilicity, the solubility of 6-Bnz-cAMP-AM in water or buffers is limited. It is suggested to use a small amount of anhydrous organic solvent such as anhydrous DMSO or DMF for dissolution at 1-100mM, and to dilute with water or buffer down to the concentrations required. In some cases, especially at high concentrations (~1mM), Pluronic® F-127 (Molecular Probes) can be useful to facilitate solubilization in physiological media. Please keep in mind that due to the high potency of 6-Bnz-cAMP-AM relatively low concentrations (0.005-0.1mM) should be sufficient, and be sure to check for DMSO/DMF tolerance in your system.

**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)