

# 8-BT-cAMP

Site selective activator of cAMP-dependent protein kinase. Increased hydrolytical stability against PDE and considerably higher membrane permeability compared to cyclic AMP. Activates membrane excitability in peptidergic neurons.

*BLG-B015-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

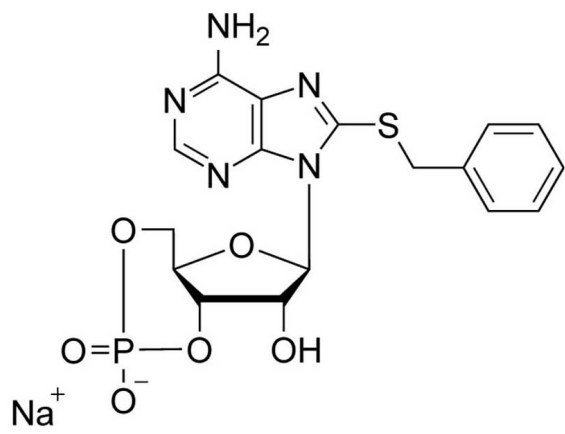
## Ordering Information

[Order Online »](#)

|             |          |
|-------------|----------|
| BLG-B015-50 | 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-Benzylthioadenosine-3',5'-cyclic monophosphate . sodium salt

**CAS** 50655-17-9

**Couple Type** Modified nucleotides

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

**MW** 473.4

**Purity** > 98% HPLC

**Quantity** 10µmol≈4.7mg

**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)