

Bn-cAMP

Membrane-permeant precursor of the second messenger cyclic AMP

Membrane-permeable precursor of cAMP. Hydrolysis slowly releases the polar cAMP which is trapped inside the cell and subsequently metabolized, resulting in an elevated cAMP level for some time.

*BLG-A065-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

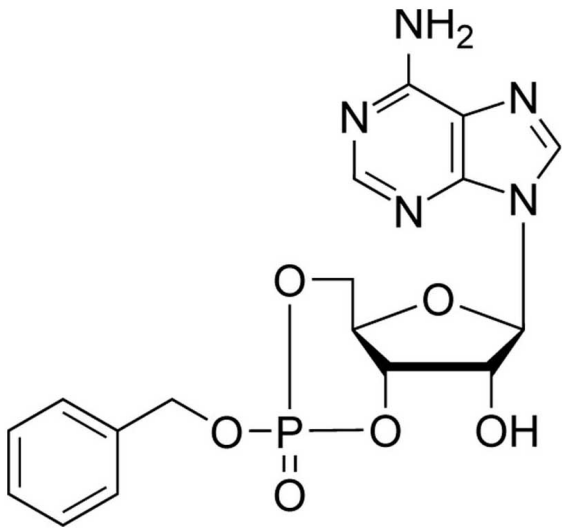
Ordering Information

[Order Online »](#)

BLG-A065-01	1 μ mol
BLG-A065-05	5x1 μ 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 Adenosine-3',5'-cyclic monophosphate . benzyl ester

CAS 62742-71-6

Couple Type Modified nucleotides

Formula $C_{17}H_{18}N_5O_6P$

MW 419.3

Purity > 95% HPLC for mixture of isomers

Quantity 1µmol≈0.4mg

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com