

6-Bnz-8-PIP-cAMP

Membrane-permeant activator of protein kinase A which does not activate Epac. Increased metabolic stability. Selects the A site of cAK I and thus can be used for synergistic activation when combined with the site B I-selective analogues 2-Cl-8-HA-cAMP (Cat. No. C 071) or 2-Cl-8-MA-cAMP (Cat. No. C 080) or 8-HA-cAMP (Cat. No. H 006).

BLG-B055-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

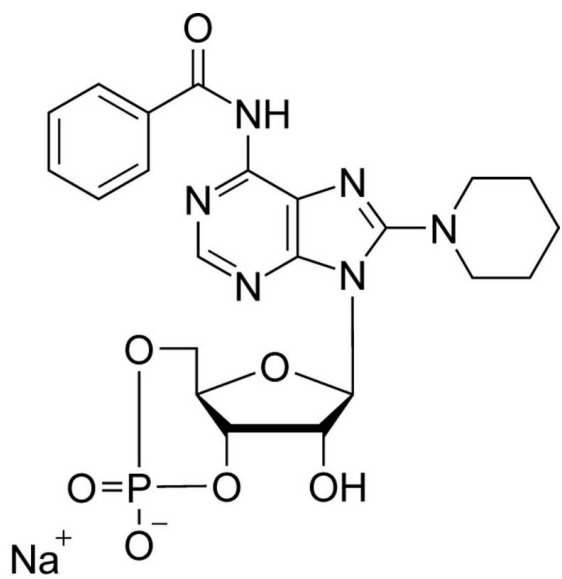
Ordering Information

[Order Online »](#)

BLG-B055-05	5 μ mol
BLG-B055-25	5x5 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 N6-Benzoyl-8-piperidinoadenosine-3',5'-cyclic monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{22}H_{26}N_6O_7P \cdot Na$

MW 540.4

Purity > 98% HPLC

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