

6-MB-cAMP

Site-selective PKA agonist

Membrane permeant, site-selective activator of protein kinase A (cyclic AMP agonist). Together with e.g. 8-HA-cAMP (Prod. No. [BLG-H006](#)), 8-ABA-cAMP (Prod. No. [BLG-A017](#)) or 8-AHA-cAMP (Prod. No. [BLG-A011](#)) selective stimulation of type I of cAK can be achieved.

*BLG-M003-250 (5 x 50 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

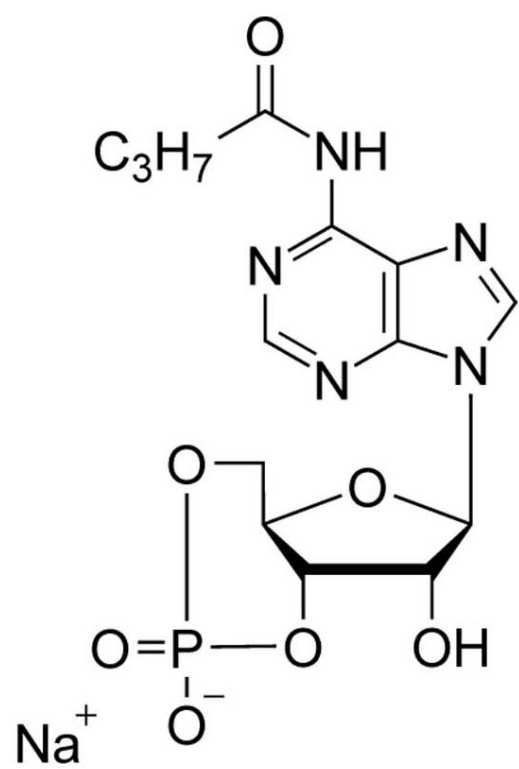
Ordering Information

[Order Online »](#)

BLG-M003-250	5x50µmol
---------------------	----------

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-Monobutyryladenosine-3',5'-cyclic monophosphate . sodium salt

CAS 70253-67-7

Couple Type Modified nucleotides

Formula $C_{14}H_{17}N_5O_7P \cdot Na$

MW 421.3

Purity > 97% HPLC

Quantity 50µmol≈21mg

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