

2'-O-Me-cAMP

Polar Epac agonist

Analogue of the natural signal molecule cyclic AMP in which the 2'-hydroxyl group of the ribose has been methylated.

Since a free 2'-ribose hydroxyl group in cyclic AMP is essential for stimulation of cAMP-dependent protein kinase (PKA), the methylated analog 2'-OMe-cAMP does not activate PKA, whereas it is a selective stimulator of the exchange factors directly activated by cAMP (Epac or cAMP-GEF).

2'-O-Me-cAMP is also suitable for testing the 2'-position of the cAMP skeleton in receptor mapping studies. In comparison with potent Epac-activating analogues such as 8-pCPT-2'-O-Me-cAMP (Prod. No. BLG-C041) or 8-pMeOPT-2'-OMe-cAMP (Prod. No. BLG-M034) 2'-O-Me-cAMP shows considerably reduced potency as well as lipophilicity which makes it much less membrane-permeant. Furthermore, it is metabolized rapidly by cyclic nucleotide-dependent phosphodiesterases. For these reasons, 2'-O-Me-cAMP is not recommended for investigating the role of Epac in cell culture experiments.

BLG-M050-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

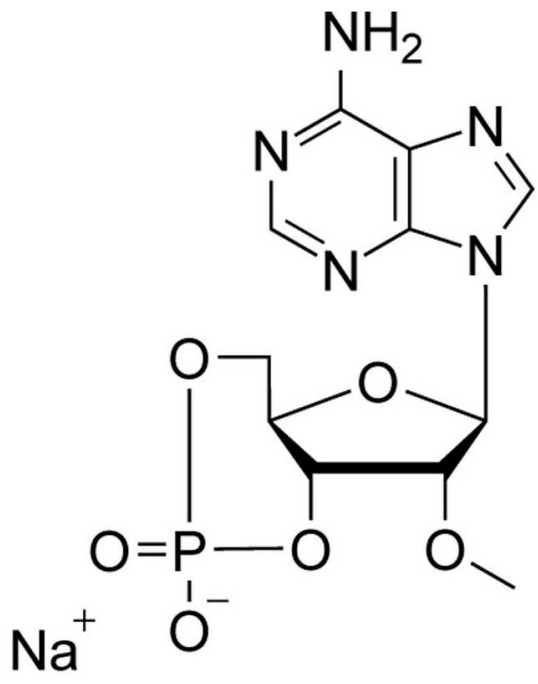
Ordering Information

[Order Online »](#)

BLG-M050-25	5x5µmol
-------------	---------

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 2'-O-Methyladenosine-3',5'-cyclic monophosphate . sodium salt

Appearance Lyophilized or crystallized sodium salt.

CAS 40269-29-2

Couple Type Modified nucleotides

Formula $C_{11}H_{13}N_5O_6P \cdot Na$

MW 365.2

Purity > 98% HPLC

Quantity 5µmol≈1.8mg

Reconstitution Soluble in water or in buffers. Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

Solubility Soluble in water or in buffers.

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)