

2'-O-MB-cAMP

Metabolically activatable prodrug form of cyclic AMP

2'-O-MB-cAMP is a lipophilic precursor of cAMP. During metabolic activation by esterases cAMP and butyrate are released.

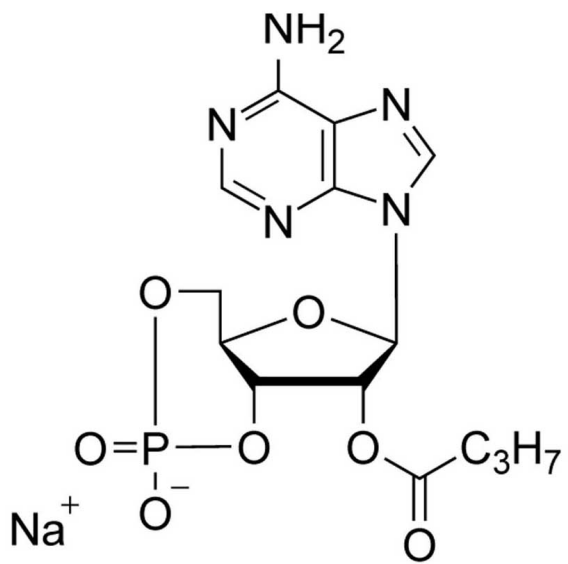
Caution: Butyrate was shown to have distinct biological effects! Significantly more lipophilic and membrane permeant compared to cAMP.

BLG-M007-250 (5 x 50 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-M007-250	5x50 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 2'-O-Monobutyryladenine-3',5'-cyclic monophosphate . sodium salt

CAS 55443-13-5

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