

5'-O-MB-2',3'-cAMP

Analogue of 2',3'-cAMP

5'-O-MB-2',3'-cAMP is a membrane permeable analogue and prodrug / precursor of 2',3'-cAMP (BIOLOG Cat. No. A 307) and was used to identify pathways regulated by 2',3'-cNMPs, including flagellar motility and biofilm formation in *E. coli*. After entering a target cell by passive diffusion, the butyrate group is split off by endogenous esterases to release the parent compound 2',3'-cAMP. Caution: Also extracellular esterases activities were reported in literature and premature release of 2',3'-cAMP strongly reduces membrane permeability. 2',3'-cAMP is a positional isomer of the natural second messenger 3',5'-cAMP (BIOLOG Cat. No. A 001) and was found to be produced by organ systems such as rat and mouse kidney and mouse brain from RNA degradation. It was also recently reported to promote the formation of stress granules in *Arabidopsis thaliana*.

BLG-M176-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

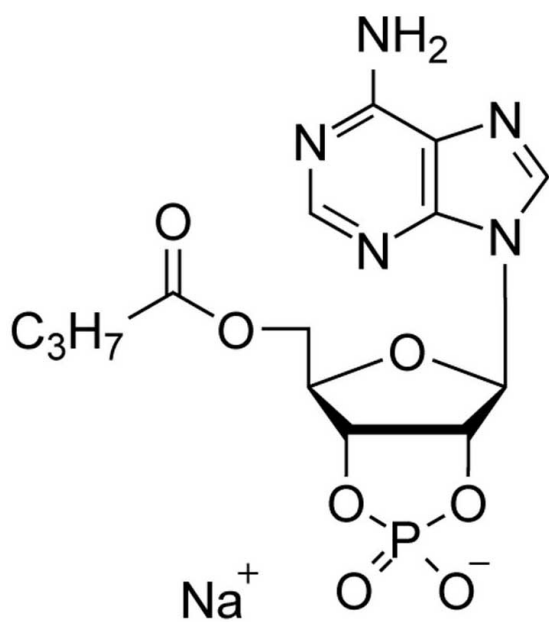
Ordering Information

[Order Online »](#)

BLG-M176-05	5µmol
BLG-M176-25	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 5'-O-Monobutyryl-adenosine-2',3'-cyclic monophosphate . sodium salt

Formula $C_{14}H_{17}N_5O_7 \cdot Na$

MW 421.3 (free acid)

Purity > 97% HPLC



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