

2'-O-MB-cGMP

Metabolically activatable PKG agonist

2'-O-MB-cGMP is a lipophilic precursor of cGMP. During metabolic activation by esterases cGMP and butyrate are released. Significantly more lipophilic and membrane permeant compared to cGMP.

Caution: This analogue releases butyrate due to intracellular and extracellular esterase action. Butyrate was shown to have distinct biological effects.

BLG-M006-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

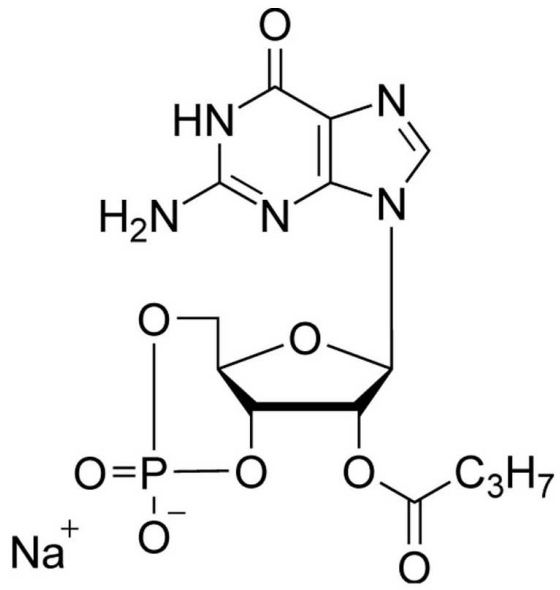
Ordering Information

[Order Online »](#)

BLG-M006-50	5x10 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-Monobutyrylguanosine-3',5'-cyclic monophosphate . sodium salt

CAS 58329-72-9

Couple Type Modified nucleotides

Formula $C_{14}H_{17}N_5O_8P \cdot Na$

MW 437.3

Purity > 97% HPLC

Quantity 10µmol≈4.4mg

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