

Rp-2'-O-MB-cAMPS

Metabolically activatable Form of the PKA inhibitor Rp-cAMPS

Lipophilic precursor of Rp-cAMPS, a competitive inhibitor of protein kinase A type I and II (cyclic AMP antagonist). During metabolic activation by esterases the inhibitor and butyrate are released. Significantly more lipophilic and membrane permeant compared to Rp-cAMPS. Resistant against mammalian cyclic nucleotide-dependent phosphodiesterases.

BLG-M004-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

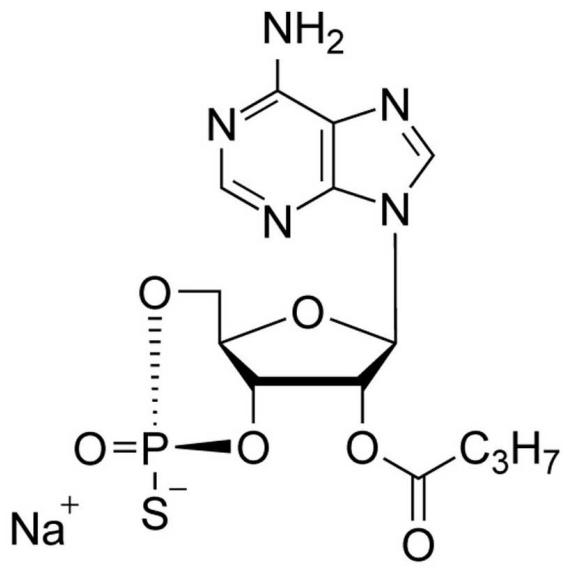
Ordering Information

[Order Online »](#)

BLG-M004-05	5 μ mol
BLG-M004-25	5x5 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-Monobutyryl-adenosine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

Appearance Lyophilized or crystallized sodium salt.

CAS 152218-23-0

Couple Type Modified nucleotides

Formula $C_{14}H_{17}N_5O_6PS \cdot Na$

MW 437.4

Purity > 97% HPLC

Quantity 5µmol≈2.2mg

Reconstitution Rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

Solubility Soluble in water or buffer.

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