

Rp-8-Br-MB-cAMPS

Metabolically activatable form of the PKA inhibitor Rp-8-Br-cAMPS

Rp-8-Br-MB-cAMPS is a lipophilic precursor of the cyclic AMP antagonist Rp-8-Br-cAMPS (Prod. No. [BLG-B001](#)). The butyryl group masks the polar 2' hydroxyl group and facilitate membrane permeability. During metabolic activation by intracellular esterases the inhibitor and butyrate are released. As observed with dibutyryl cAMP, release of butyrate can already start in the medium if it contains serum esterases. Please note that butyrate can have its own biochemical effects, therefore a control experiment with sodium butyrate is necessary.

Significantly more lipophilic and membrane permeant compared to Rp-8-Br-cAMPS.

BLG-B010-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

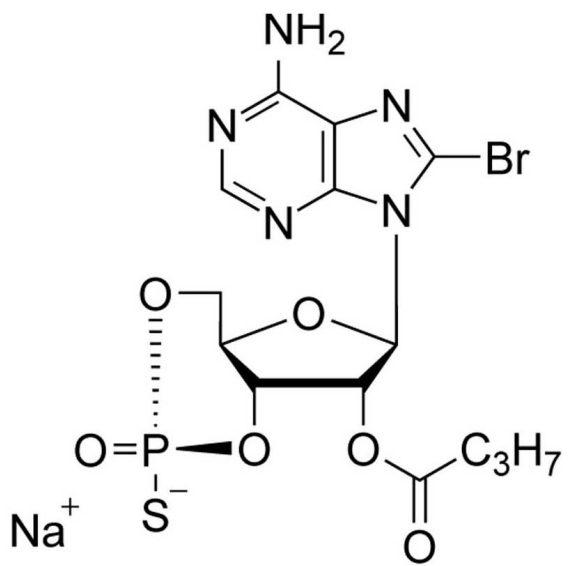
Ordering Information

[Order Online »](#)

BLG-B010-05	5 μ mol
BLG-B010-25	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 8-Bromo-2'-O-monobutryladenine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

Appearance Lyophilized or crystallized sodium salt.

CAS 788807-32-9

Couple Type Modified nucleotides

Formula $C_{14}H_{16}BrN_5O_6PS \cdot Na$

MW 516.2

Purity > 97% HPLC

Quantity 5µmol≈2.6mg

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-fr@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