

Rp-8-PIP-cAMPS

Lipophilic, site-selective analogue of the protein kinase A inhibitor Rp-cAMPS

Lipophilic site selective analog of the protein kinase inhibitor Rp-cAMPS with extremely high preference for site B of cAK type II. Significantly more lipophilic and membrane permeant compared to Rp-cAMPS. Resistant against mammalian cyclic nucleotide-dependent phosphodiesterases.

BLG-P004-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

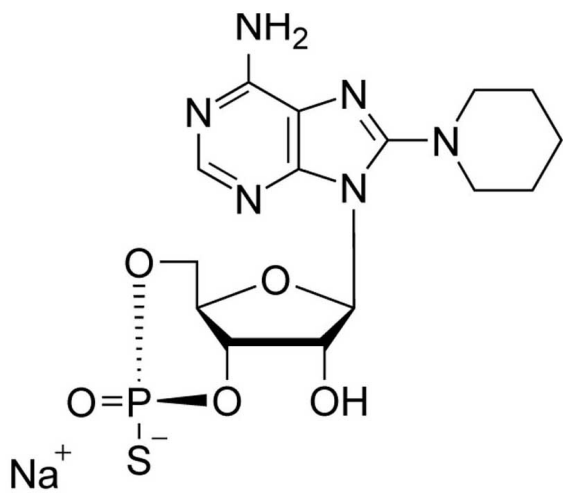
Ordering Information

[Order Online »](#)

BLG-P004-05	5µmol
BLG-P004-25	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 8-Piperidinoadenosine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

Appearance Lyophilized or crystallized sodium salt.

CAS 156816-36-3

Couple Type Modified nucleotides

Formula $C_{15}H_{20}N_6O_5PS \cdot Na$

MW 450.4

Purity > 99% HPLC

Quantity 5µmol≈2.3mg

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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)