

Rp-8-Cl-cAMPS

Membrane-permeant analogue of the cyclic AMP antagonist Rp-cAMPS

Rp-8-Cl-cAMPS is a novel combination of the well accepted protein kinase inhibitor Rp-cAMPS with 8-chloro cyclic AMP, resulting in a membrane permeant cyclic AMP antagonist which is not metabolized by mammalian cyclic nucleotide phosphodiesterases.

Rp-8-Cl-cAMPS is about 1.3 times more lipophilic compared to 8-Cl-cAMP, and 1.7 times and 2.2 times more compared to Rp-cAMPS and cAMP respectively.

In contrast to common ATP-site inhibitors or to its parent compound Rp-cAMPS, Rp-8-Cl-cAMPS discriminates between both isozymes of protein kinase A, preferring type I¹³ and thus provides additional selectivity. By occupying cAMP binding sites Rp-8-Cl-cAMPS prevents the kinase holoenzyme from dissociation and thus from activation.

Due to this working principle preincubation of the inhibitor prior to the activation step is necessary for optimal results.

BLG-C003-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

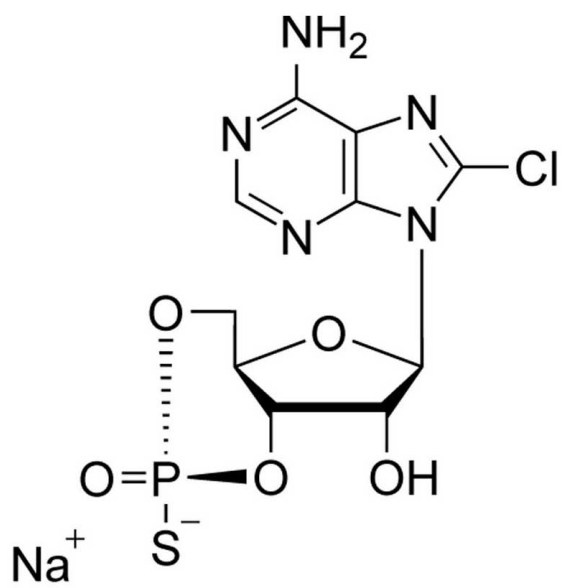
Ordering Information

[Order Online »](#)

BLG-C003-05	5µmol
BLG-C003-25	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-Chloroadenosine-3',5'-cyclic monophosphorothioate . sodium salt, Rp-isomer

Appearance Lyophilized or crystallized sodium salt.

CAS 142754-27-6

Couple Type Modified nucleotides

Formula $C_{10}H_{10}ClN_5O_5PS \cdot Na$

MW 401.7

Purity > 99% HPLC

Quantity 5µmol≈2mg

Reconstitution Please 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)