

8-Cl-cAMP

Site-selective PKA agonist

8-Cl-cAMP is a potent stimulator of cAMP-dependent protein kinases with high affinity for sites A and B of PKA type I. In addition, it shows high site selectivity for site B of PKA type II. The analogue is degraded by cyclic nucleotide phosphodiesterases more slowly compared to natural cyclic AMP.

The increased lipophilicity which resembles 8-bromo cyclic AMP and Sp-cAMPS enables membrane permeability in several biosystems.

Due to its ability to stop growth of several cancer cell lines at very low doses, its therapeutic use is under investigation. However, recent reports on active 8-Cl- metabolites showed that also degradation products could be responsible for cytostatic properties.

BLG-C007-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

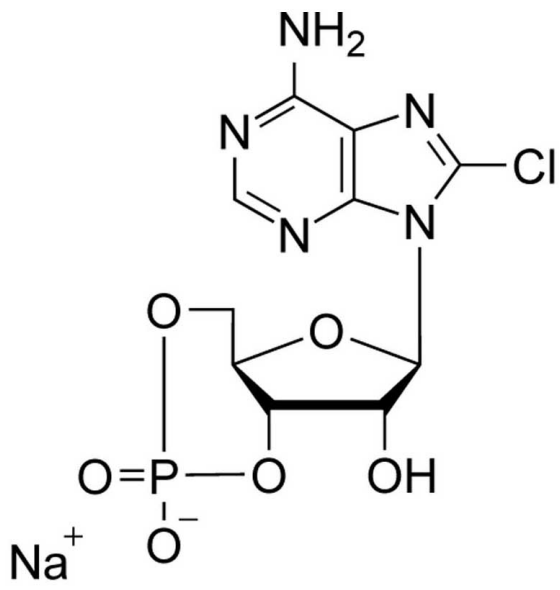
Ordering Information

[Order Online »](#)

BLG-C007-50	5x10µmol
-------------	----------

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 monophosphate . sodium salt

Appearance Lyophilized or crystallized sodium salt.

CAS 124705-03-9

Couple Type Activator, Modified nucleotides

Formula $C_{10}H_{10}ClN_5O_6P \cdot Na$

MW 385.6

Purity > 99% HPLC

Quantity 10µmol≈3.9mg

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