

8-CPT-cAMP

Lipophilic activator of both cAMP- and cGMP-dependent protein kinase and of Epac

Lipophilic activator of both cAMP- and cGMP- dependent protein kinase. Higher activation potential compared to cAMP; excellent cell membrane permeability, improved phosphodiesterase stability and high site selectivity preferring site B of cAK type II.

Caution: This analogue is not selective but activates cGMP-dependent protein kinase as well! In addition, it increases basal cGMP level by inhibition of PDE V.

For specific cAMP pathway activation the highly permeant Sp-5,6-DCI-cBIMPS (Prod. No. [BLG-D014](#)) is the better choice.

BLG-C010-500 (5 x 100 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

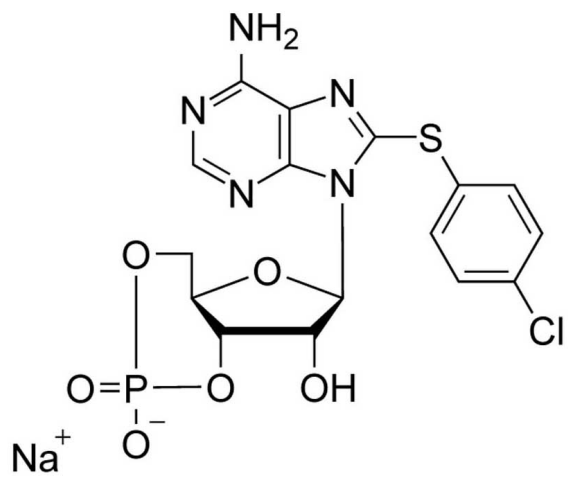
Ordering Information

[Order Online »](#)

BLG-C010-500	5x100µmol
--------------	-----------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-(4-Chlorophenylthio)adenosine-3',5'-cyclic monophosphate . sodium salt

CAS 93882-12-3

Couple Type Modified nucleotides

Formula $C_{16}H_{14}ClN_5O_6PS \cdot Na$

MW 493.8

Purity > 99% HPLC

Quantity 100µmol≈49mg

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