

Rp-8-CPT-cAMPS

Membrane-permeant and PDE-resistant PKA inhibitor

Rp-8-CPT-cAMPS is a membrane permeant, isozyme-selective inhibitor of protein kinase A with special preference for cAK II. Its metabolic stability avoids potential side effects through active metabolites. In contrast to common ATP-site inhibitors, Rp-8-CPT-cAMPS occupies cAMP binding sites at the regulatory subunit of PKA and 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-C011-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

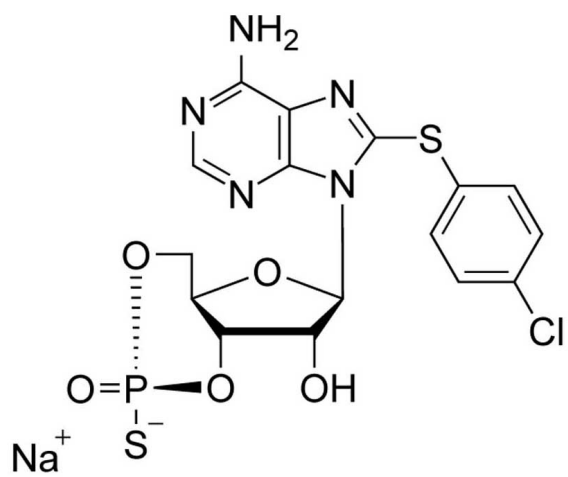
Ordering Information

[Order Online »](#)

BLG-C011-05	5µmol
BLG-C011-25	5x5µmol

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 monophosphorothioate . sodium salt, Rp-isomer
Appearance	Lyophilized or crystallized sodium salt.
CAS	129735-01-9
Couple Type	Inhibitor, Modified nucleotides
Formula	$C_{16}H_{14}ClN_5O_5PS_2 \cdot Na$
MW	509.8
Purity	> 99% HPLC
Quantity	5µmol≈2.5mg
Reconstitution	Rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.
Solubility	Limited soluble in water.
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