

Sp-8-pCPT-2'-OMe-cAMPS

PDE-resistant, membrane-permeant Epac agonist

Analog of the natural signal molecule cyclic AMP in which the hydrogen in position 8 of the heterocyclic nucleobase is replaced by the lipophilic 4-chlorophenylthio moiety, and the ribose 2'-hydroxyl group has been methylated. In addition, the axial one of the two exocyclic oxygen atoms in the cyclic phosphate moiety is modified by sulfur. The suffix "p" indicates that the R/S nomenclature refers to phosphorus atom.

Hydrolysis-resistant form of 8-pCPT-2'-O-Me-cAMP (Prod. no. BLG-C041), a specific activator of the exchange protein directly activated by cyclic AMP (Epac) or cAMP-GEF, respectively. The compound does not activate protein kinase A, has high lipophilicity and membrane permeability and is not metabolized by phosphodiesterases.

BLG-C052-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

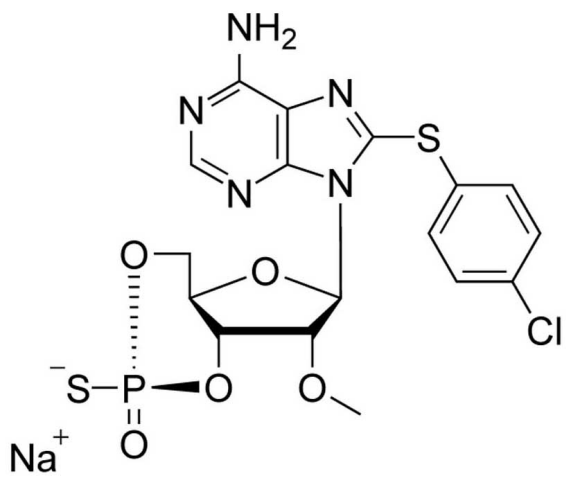
Ordering Information

[Order Online »](#)

BLG-C052-01	1µmol
BLG-C052-05	5x1µmol

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-(4-Chlorophenylthio)-2'-O-methyl-cAMPS, Sp isomer, sodium salt

CAS 634208-37-0

Couple Type Activator, Modified nucleotides

Formula $C_{17}H_{16}ClN_5O_5PS_2 \cdot Na$

MW 523.9

Purity > 98% HPLC

Quantity 1µmol≈0.5mg

Reconstitution 5-10mM stock solutions in water can be obtained. Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

Solubility Limited solubility in water.

Technical Info / Product For the Original Manufacturer's data sheet please [click here](#).

Notes



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