

Sp-8-pCPT-cGMPS

. Na

Membrane-permeant and PDE-resistant PKG agonist

Lipophilic activator of both protein kinase G (I α , I β & type II) and protein kinase A type II. Excellent cell membrane permeability and phosphodiesterase stability.

BLG-C014-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

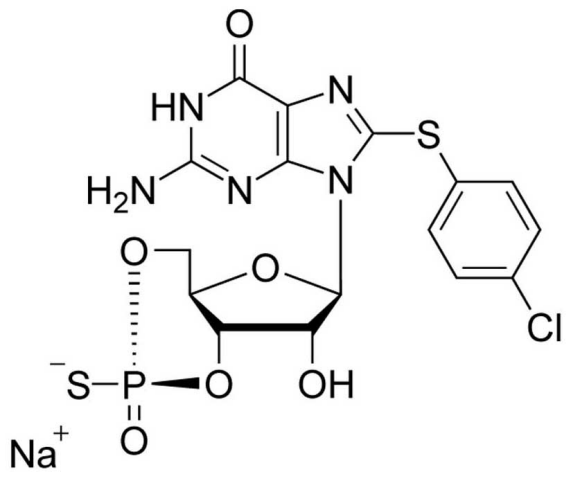
Ordering Information

[Order Online »](#)

BLG-C014-05	5 μ mol
BLG-C014-25	5x5 μ mol

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)guanosine-3',5'-cyclic monophosphorothioate . sodium salt, Sp-isomer
CAS	160385-87-5
Couple Type	Activator, Modified nucleotides
Formula	$C_{16}H_{14}ClN_5O_6PS_2 \cdot Na$
MW	525.9
Purity	> 99% HPLC
Quantity	5µmol≈2.6mg
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)