

Rp-8-pCPT-cGMPS . sodium salt

PKG inhibitor

A cell-permeable cGMP antagonist which is resistant to hydrolysis by cyclic nucleotide phosphodiesterases ($K_i=0.5 \mu\text{M}$). No metabolic side effects. Significantly more lipophilic and membrane permeant as compared to Rp-cGMPS (Prod. No. ALX-480-072) or Rp-8-Br-cGMPS (Prod. No. BML-CN216)

Citations: 8

[View Online »](#)

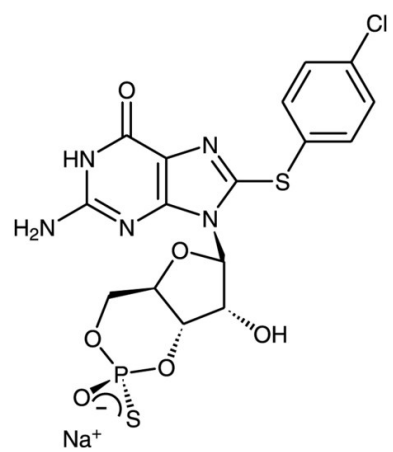
Ordering Information

[Order Online »](#)

BML-CN206-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	153660-04-9
Couple Target	PKG
Couple Type	Inhibitor
Formula	$C_{16}H_{14}N_5O_6PS_2Cl \cdot Na$
MW	525.9
Purity	≥98% (HPLC)
Solubility	Soluble in water (25 mg/ml).

Last modified: May 29, 2024



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)