

8-(4-

Chlorophenylthio)guanosine

3',5'-cyclic

Monophosphate .

sodium salt

Selective activator of cAMP-dependent protein kinase (PKG) I α , I β , and type II as well as of cGMP-gated ion channels. Higher activation potential as compared to cGMP; excellent cell membrane permeability and phosphodiesterase stability. Preferred alternative to unsuitable analogs such as dibutyl-cGMP.

Activator of PKG and CNG channels

Citations: 2

[View Online »](#)

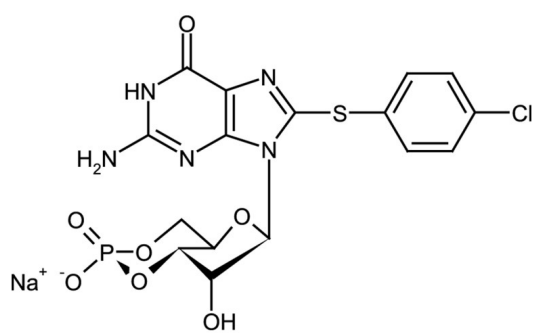
Ordering Information

[Order Online »](#)

ALX-480-088-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-pCPT-cGMP . Na

Appearance White to off-white powder.

CAS 51239-26-0

Couple Target CNG channel, PKG

Couple Type Activator

Formula $C_{16}H_{14}ClN_5O_7PS \cdot Na$

Identity Determined by ESI-MS and UV/VIS.

MW 486.8 . 23.0

Purity ≥99% (HPLC)

Solubility Soluble in water.



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-
fr@enzolifesciences.com](mailto:info-fr@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)