

8-(4-

Chlorophenylthio)guanosine

3',5'-cyclic

Monophosphothioate,

Rp-Isomer,

triethylammonium salt  
Potent selective inhibitor of cGMP-dependent protein kinase 1 $\alpha$ , 1 $\beta$  and type II. Resistant to mammalian cyclic nucleotide-dependent phosphodiesterases, 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).

(high purity)

Citations: 4

[View Online »](#)

PKG inhibitor

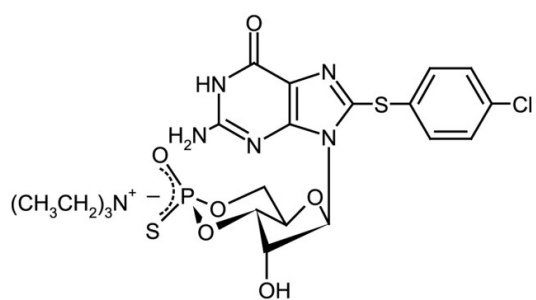
Ordering Information

[Order Online »](#)

|                  |             |
|------------------|-------------|
| ALX-480-032-M001 | 1mg         |
| ALX-480-032-C100 | 100 $\mu$ g |

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** Rp-8-pCPT-cGMPS . TEA, 8-pCPT-cGMPS . TEA, Rp-Isomer

**Appearance** White to off-white crystalline powder.

**CAS** 153660-04-9

**Couple Target** PKG

**Couple Type** Inhibitor

**Formula**  $C_{16}H_{14}ClN_5O_6PS_2 \cdot C_6H_{16}N$

**Identity** Determined by UV and ESI-MS.

**MW** 502.9 . 102.2

**Purity** ≥99% (HPLC)

**Solubility** Soluble in water and aqueous buffers.

**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)