

8-(4-

Chlorophenylthio)guanosine

3',5'-cyclic

Monophosphothioate,

Rp-Isomer,

triethylammonium salt
Potent selective inhibitor of cGMP-dependent protein kinase 1 α , 1 β 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

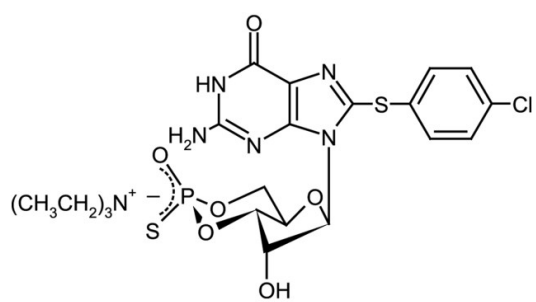
Ordering Information

[Order Online »](#)

ALX-480-032-M001	1mg
ALX-480-032-C100	100 μ 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)