

# Guanosine 3',5'-cyclic Monophosphothioate, Rp-Isomer . sodium salt

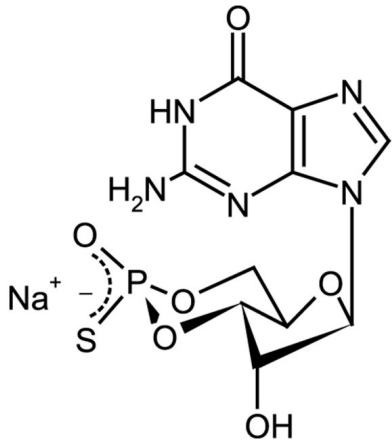
**PKG inhibitor**  
Inhibitor of cGMP-dependent protein kinase (PKG) 1 $\alpha$  (cGMP antagonist).  
Resistant against mammalian cyclic nucleotid-dependent  
phosphodiesterases. No metabolic side effects. Low membrane  
permeability.

Citations: 3

[View Online »](#)

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** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Rp-cGMPS . Na, cGMPS . Na, Rp-Isomer

**Appearance** White to off-white solid.

**CAS** 86562-09-6

**Couple Target** PKG

**Couple Type** Inhibitor

**Formula**  $C_{10}H_{11}N_5O_6SP \cdot Na$

**MW** 360.3 . 23.0

**Purity** ≥98% (HPLC)

**Solubility** Soluble in water (125mM), DMSO (2mM), DMF (1mM), methanol (7mM) or PBS, pH 7.4 (125mM).

**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)