## 8-(4-Chlorophenylthio)guanosin 3',5'-cyclic Monophosphate .

SGE LILLION c Statement protein kinase (PKG) Iα, Iβ, and type II as well as of cGMP-gated ion channels. Higher activation potential as compared to cGMP; excellent cell membrane permeability and prosphetically preferred alternative to unsuitable analogs such as dibutyryl-cGMP.

Citations: 2

View Online »

**Ordering Information** 

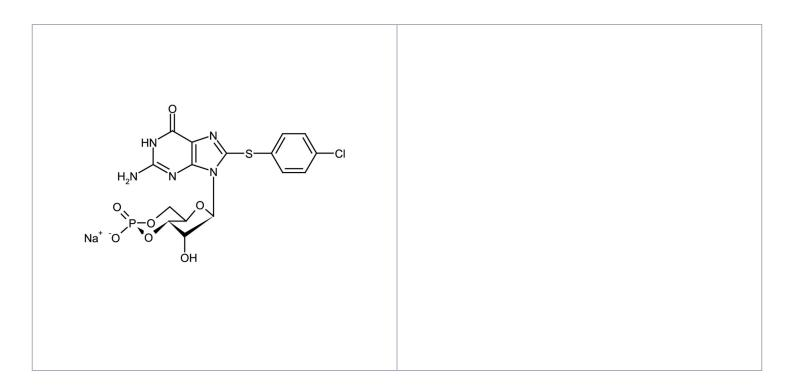
Order Online »

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  ${
m C_{16}H_{14}CIN_5O_7PS}$  . Na

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

**MW** 486.8 . 23.0

Purity ≥99% (HPLC)

**Solubility** Soluble in water.

eu@enzolifesciences.com

Last modified: May 29, 2024