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

treitethayrbammenbehmsesaltd 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 attendable to unsuitable analogs such as dibutyryl-cGMP.

Citations: 2

View Online »

Ordering Information

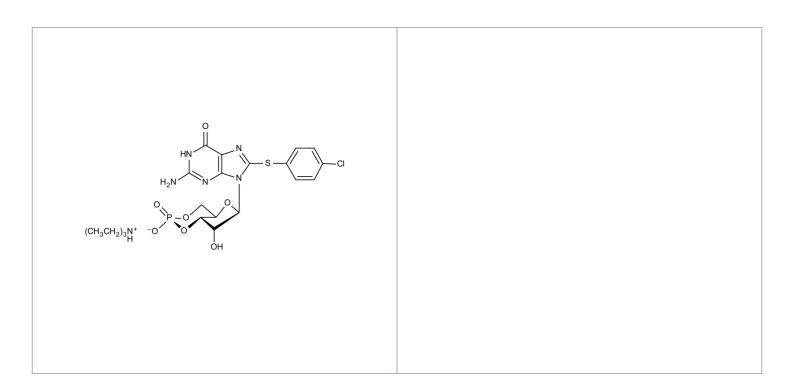
Order Online »

ALX-480-031-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. TEA

Appearance White to off-white powder.

CAS 54364-02-2

CNG channel, PKG **Couple Target**

Couple Type Activator

 $C_{16}H_{14}CIN_5O_7PS \cdot C_6H_{16}N$ **Formula**

Identity Determined by MS and UV.

MW 486.8 . 102.2

Purity ≥99% (HPLC)

Solubility Soluble in water.

Last modified: May 29, 2024

info-