Sp-8-Br-cGMPS

cGMP agonist

Cell-permeable, phosphodiesterase-resistant cGMP agonist. This analog activates PKA and PKG with similar potency. Much more lipophilic and membrane permeant as compared to cyclic GMP (Prod. No. ALX-480-059) or Sp-cGMPS (Prod. No. ALX-480-073)

Citations: 1

View Online »

Ordering Information

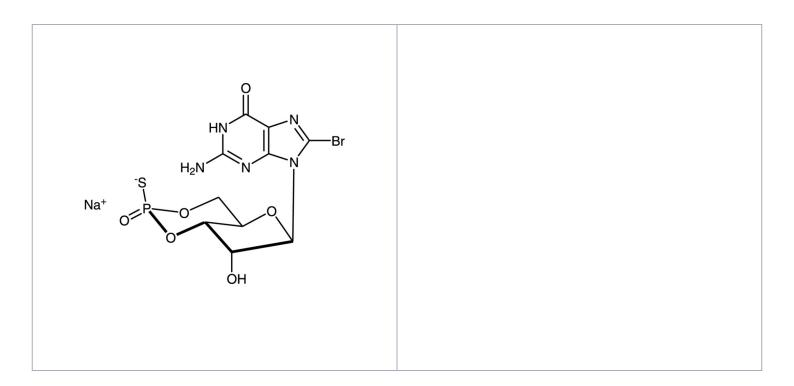
Order Online »

BML-CN217-0500

500µg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Lyophilized solid.

CAS 153660-03-8

Couple Target PKA, PKG

Couple Type Inhibitor

 $\mathbf{Formula} \qquad \qquad \mathbf{C_{10}H_{10}BrN_5O_6PS} \text{ . Na}$

MW 462.2

Purity ≥98% (HPLC)

Solubility Soluble in water (25mg/ml).

Last modified: May 29, 2024

