# Sp-cAMPS

#### **PKA** activator

Sp-cAMPS is a potent and specific activator of cAMP-dependent protein kinases. Its cell permeability and complete resistance to cyclic nucleotide phosphodiesterases make it a unique tool for studying cAMP-dependent signaling. Has a considerably higher resistance against cyclic nucleotide-dependent phosphodiesterases as compared to dibutyryl-cAMP or 8-Br-cAMP. Concentration for half-maximal activation is 1.8µM. The product is not sterile. No metabolic side effects.

Citations: 11

**View Online »** 

**Ordering Information** 

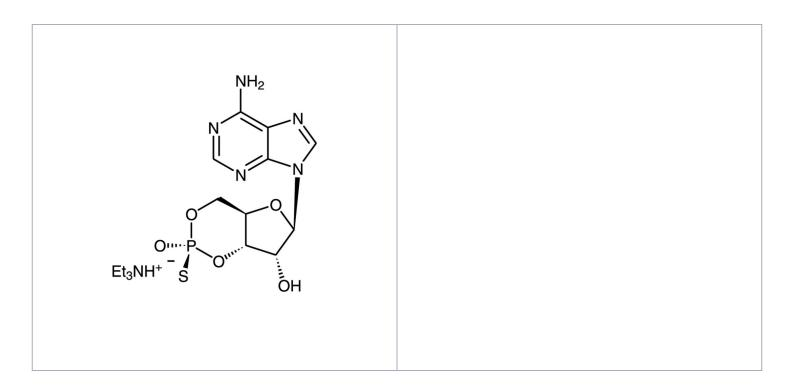
Order Online »

BML-CN136-0001

1mg

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 one 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** 93602-66-5

Couple Target PKA

Couple Type Activator

Formula  $C_{10}H_{11}N_5O_5PS \cdot C_6H_{16}N$ 

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

**MW** 446.5

Purity ≥98% (HPLC)

**Solubility** Soluble in water (25mg/ml), DMSO, dimethyl formamide or

methanol

Last modified: May 29, 2024

