Sp-5,6-DichlorocBIMPS

PKA activator

A cAMP analog which is a potent and specific activator of cAMP-dependent protein kinase. It is cell permeable and is not significantly hydrolyzed by type I, II or III phosphodiesterases. It may be used to distinguish the cAMP-versus the cGMP-mediated signal transduction pathways. Surpasses the widely used but problematic dibutyryl-cAMP (Prod. No. BML-CN125) or 8-CPT-cAMP (Prod. No. BML-CN130).

Citations: 5

View Online »

Ordering Information

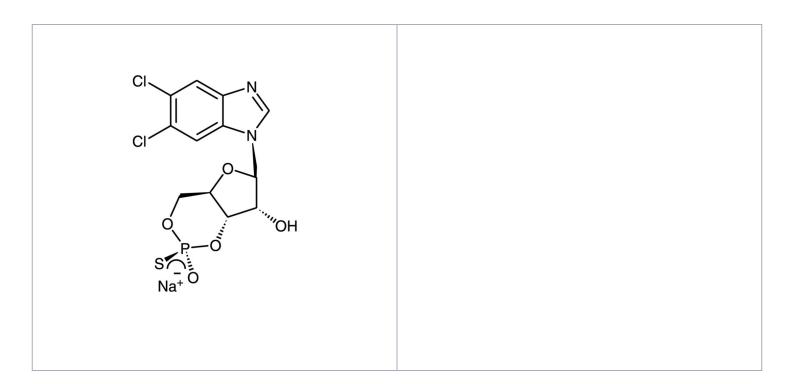
Order Online »

BML-CN120-0001

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 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 White solid.

CAS 120912-54-1

Couple Target PKA

Couple Type Activator

Formula $C_{12}H_{10}N_2O_5PSCI_2$. Na

MW 419.2

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (167 mM), dimethyl formamide (167

mM), 96 % ethanol (167 mM), methanol (25 mM), water (25 mM) and aqueous buffers such as phostphate buffered saline (pH 7.4; 1.4 mM), 25 mM TRIS (pH 7.4-; 25 mM), 25

mM HEPES (pH 7.2; 35.7 mM) or Na2HPO4.

Source Synthetic.

Last modified: May 29, 2024

