

Sp-5,6-Dichloro-cBIMPS

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 »](#)

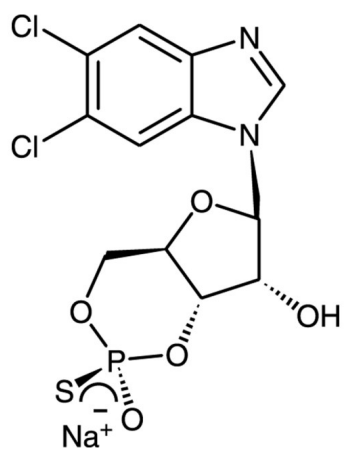
Ordering Information

[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	$C_{12}H_{10}N_2O_5PSCl_2 \cdot Na$
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 phosphate 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 Na ₂ HPO ₄ .
Source	Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)