

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 dibutyl-cAMP or 8-Br-cAMP. Concentration for half-maximal activation is 1.8 μ M. The product is not sterile. No metabolic side effects.

Citations: 11

[View Online »](#)

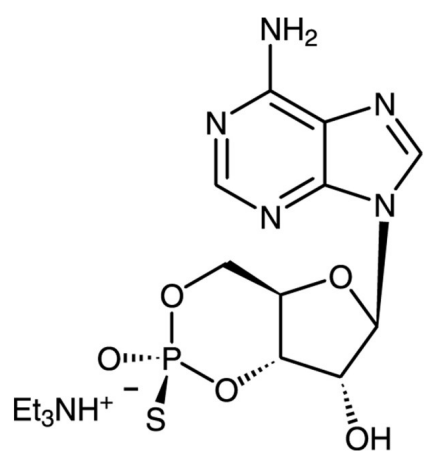
Ordering Information

[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



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)