

# Rp-cAMPS

## PKA inhibitor

Rp-cAMPS is a potent and specific competitive inhibitor of the activation of cAMP-dependent protein kinases by cAMP ( $IC_{50}=4.9\text{ }\mu\text{M}$ ). Its cell permeability and complete resistance to cyclic nucleotide phosphodiesterases make it a unique tool for studying cAMP-dependent signaling. Significantly more lipophilic as compared to cAMP and thus membrane permeable in many biological systems. This product is not sterile.

Citations: 14

[View Online »](#)

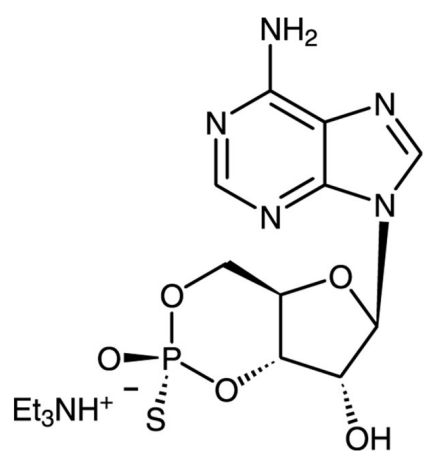
## Ordering Information

[Order Online »](#)

BML-CN135-0001	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	Clear film.
CAS	151837-09-1
Couple Target	PKA
Couple Type	Inhibitor
Formula	$C_{10}H_{11}N_5O_5PS \cdot C_6H_{16}N$
MW	446.5
Purity	≥98% (HPLC)
Solubility	Soluble in water (25mg/ml), aqueous buffers, methanol, 100% ethanol or DMSO.
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-  
fr@enzolifesciences.com](mailto:info-fr@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)