

Adenosine 3',5'-cyclic Monophosphate . sodium salt

PKA activator

Key regulator of many cellular reactions. Naturally occurring activator of cAMP-dependent protein kinase (PKA).

Citations: 1

[View Online »](#)

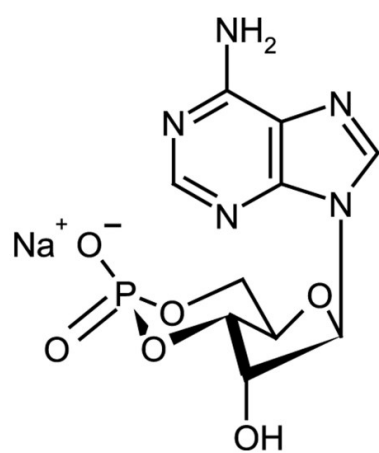
Ordering Information

[Order Online »](#)

ALX-480-011-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions should be stored in the refrigerator and should be lyophilized and frozen for longer storage periods.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Cyclic AMP . Na, cAMP . Na
Appearance	White to off-white powder.
CAS	37839-81-9
Couple Target	PKA
Couple Type	Activator
Formula	$C_{10}H_{11}N_5O_6P \cdot Na$
Identity	Identity determined by ESI-MS and UV/VIS
MI	14: 2708
MW	328.2 . 23.0
Purity	≥98% (HPLC)
Solubility	Soluble in water.

Last modified: May 29, 2024



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)