

8-CPT-cAMP

PKA activator

A lipophilic, cell-permeable cAMP analog which activates cAMP- and cGMP-dependent protein kinase. More potent than cAMP and more resistant to phosphodiesterases, with good selectivity for site B of PKA. For specific activation of the cAMP pathway the highly membrane permeable Sp-5,6-DCI-cBIMPS (Prod. No. BML-CN135).

Citations: 13

[View Online »](#)

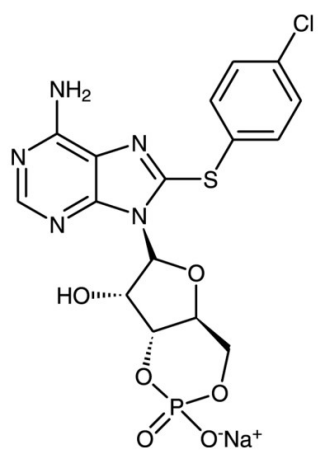
Ordering Information

[Order Online »](#)

BML-CN130-0020	20mg
BML-CN130-0100	100mg

Manuals, SDS & CoFA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store as supplied (solid) at -20°C for 1 year. Once dissolved in water, it can be stored frozen at -20°C up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	93882-12-3
Couple Target	PKA
Couple Type	Activator
Formula	$C_{16}H_{14}ClN_5O_6PS \cdot Na$
MW	493.8
Purity	≥96% (HPLC)
Solubility	Soluble in water (25mg/ml).
Source	Synthetic.

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)