

8-Bromo-cAMP

PKA activator

A cell-permeable cAMP analog which is more resistant to phosphodiesterases than cAMP and which preferentially activates cAMP-dependent protein kinase (PKA). Has increased metabolic stability and membrane permeability compared with cAMP.

Citations: 5

[View Online »](#)

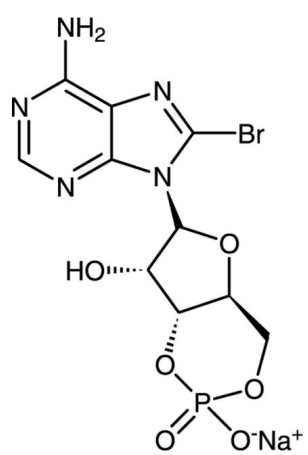
Ordering Information

[Order Online »](#)

BML-CN115-0010	10mg
BML-CN115-0050	50mg

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 to grey powder
CAS	76939-46-3
Couple Target	PKA
Couple Type	Activator
Formula	$C_{10}H_{10}BrN_5O_6P \cdot Na$
Identity	Determined by 1H -NMR, ^{31}P -NMR, ESI-MS and UV.
MW	430.1
Purity	≥98% (HPLC)
Solubility	Soluble in water (100mg/ml).

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)