

Dibutyryl-cAMP

PKA activator

Dibutyryl-cAMP (DB-cAMP) is a cell-permeable cAMP analog that preferentially activates cAMP-dependent protein kinase (e.g. protein kinase A) and inhibits phosphodiesterases. Used together with other compounds, DB-cAMP increases survival rate and promote differentiation of neural stem/progenitor cells into neurons *in vivo*. DB-cAMP also has anti-inflammatory properties and can be used to facilitate wound healing. Caution: Due to extra- and intra-cellular esterases, DB-cAMP releases butyrate, which may have distinct biological effects.

Citations: 20

[View Online »](#)

Ordering Information

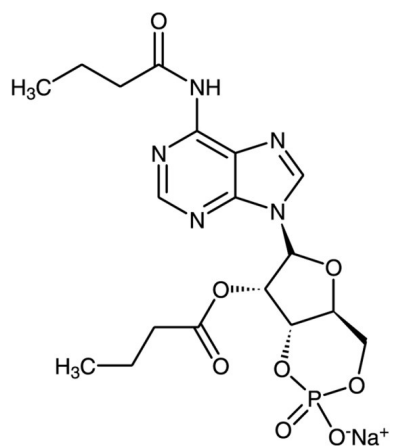
[Order Online »](#)

BML-CN125-0030	30mg
BML-CN125-0100	100mg

Manuals, SDS & CofA

[View Online »](#)

- Cell-permeable PKA activator
- Compound used to study neurogenesis



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solution must 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	Dibutyl cyclic AMP sodium salt, N6-2'-O-dibutyryladenosine-3',5'-cyclic monophosphate, DB-cAMP
Appearance	White crystalline solid.
CAS	16980-89-5
Couple Target	PKA
Couple Type	Activator
Formula	$C_{18}H_{23}N_5O_8P \cdot Na$
Identity	Identity determined by MS, ¹ H-NMR, ³¹ P-NMR, and UV spectrum.
MW	491.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (200mM), 96% ethanol (25mM), or water (100mg/ml).
Source	Synthetic.



Last modified: May 29, 2024

ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com