

Dibutyryl-cGMP

PKG activator

A cell-permeable cGMP analog which preferentially activates cGMP-dependent protein kinase. Inhibits thrombin stimulated arachidonic acid release in human platelets. Cell permeable.

Citations: 7

[View Online »](#)

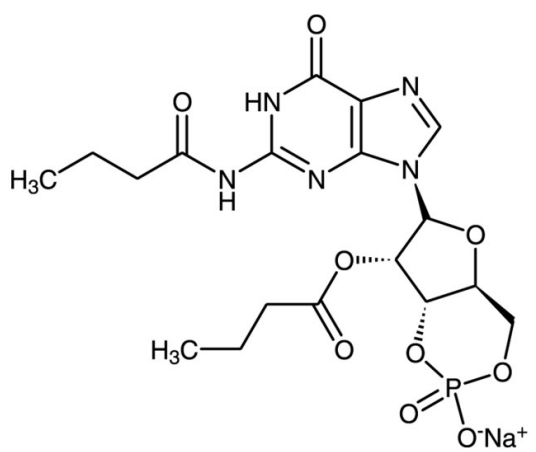
Ordering Information

[Order Online »](#)

BML-CN215-0010	10mg
BML-CN215-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Make solutions fresh daily. Avoid pH >8.5.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	51116-00-8
Couple Target	PKG
Couple Type	Activator
Formula	$C_{18}H_{23}N_5O_9P \cdot Na$
MW	484.4 . 23.0
Purity	≥96%
Solubility	Soluble in water (100mg/ml). Also soluble in 100% ethanol, DMSO, dimethyl formamide or methanol.

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-
be@enzolifesciences.com](mailto:info-be@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)