

# 8-Bromo-cGMP

## PKG activator

A cell-permeable cGMP analog which is more resistant to phosphodiesterases than cGMP and which preferentially activates cGMP-dependent protein kinase. It inhibits thrombin stimulated arachidonic acid release in human platelets.

Citations: 11

[View Online »](#)

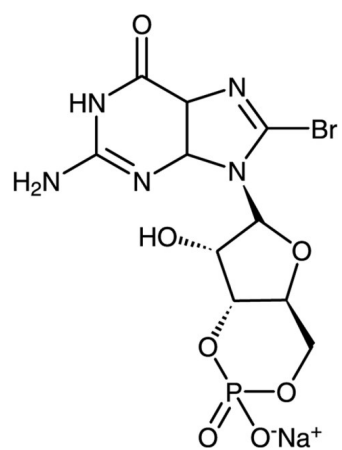
## Ordering Information

[Order Online »](#)

BML-CN205-0010	10mg
BML-CN205-0050	50mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Appearance</b>	White to off-white powder
<b>CAS</b>	51116-01-9
<b>Couple Target</b>	PKG
<b>Couple Type</b>	Activator
<b>Formula</b>	$C_{10}H_{10}BrN_5O_7PNa$
<b>Identity</b>	Determined by ESI-MS and UV/VIS.
<b>MW</b>	446.1
<b>Purity</b>	≥96% (HPLC)
<b>Solubility</b>	Soluble in water (25mg/ml) or aqueous buffers. Sparingly soluble in methanol (4mM). Insoluble in DMSO, dimethyl formamide and 96% ethanol.



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)