

8-Br-cGMP

PKG Agonist

Activator of cGMP-dependent protein kinases I α , I β & type II and cGMP-dependent ion channels (cyclic GMP agonist). Increased metabolic stability compared to cGMP, but not completely stable as often stated. 8-pCPT-cGMP (Prod. No. [BLG-C009](#)), a highly membrane permeant and metabolically resistant cGMP-agonists seems to be superior for most applications.

BLG-B004-250 (5 x 50 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

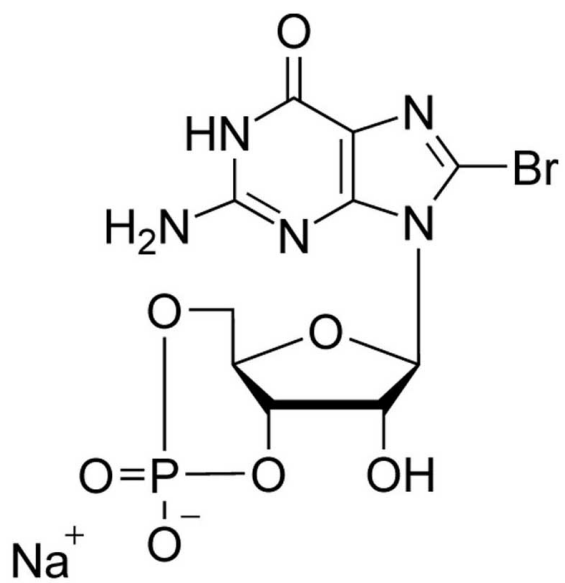
Ordering Information

[Order Online »](#)

BLG-B004-250	5x50 μ mol
--------------	----------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-Bromoguanosine-3',5'-cyclic monophosphate . sodium salt

CAS 51116-01-9

Couple Type Activator, Modified nucleotides

Formula $C_{10}H_{10}BrN_5O_7P \cdot Na$

MW 446.1

Purity > 99% HPLC

Quantity 50µmol≈22mg

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



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-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@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