

DB-cGMP

Metabolically activatable PKG agonist

We do not recommend to use this cGMP analogue due to its lacking potency and selectivity. The metabolically released N²-monobuteryl cGMP is only a poor activator of cGMP-dependent protein kinase, while butyrate can disturb seriously with own effects. 8-pCPT-cGMP (Prod. No. [BLG-C009](#)) or other cGMP analogues are much better and selective agonists.

BLG-D010-250 (5 x 50 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

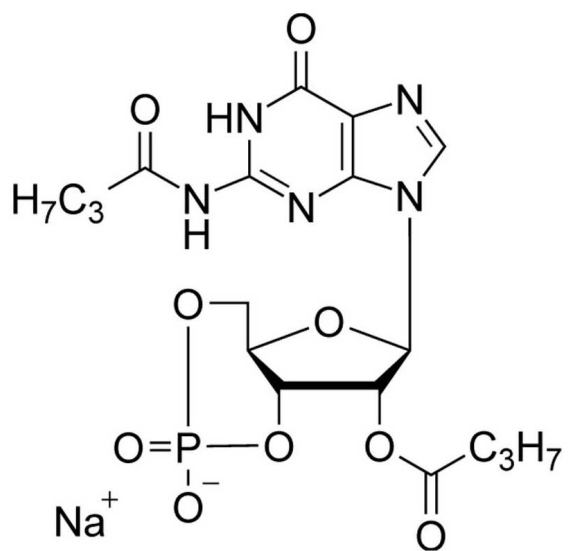
Ordering Information

[Order Online »](#)

BLG-D010-50	50µmol
BLG-D010-250	5x50µmol

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 N2,2'-O-Dibutyrylguanosine-3',5'-cyclic monophosphate . sodium salt

CAS 51116-00-8

Couple Type Modified nucleotides

Formula $C_{18}H_{23}N_5O_9P \cdot Na$

MW 507.4

Purity > 97% HPLC

Quantity 50µmol≈25mg

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