

# 8-NBD-cGMP

## Fluorescent analogue of cyclic GMP

Potent membrane permeant fluorescent activator of both isozymes I alpha and I  $\beta$  of cGMP-dependent protein kinase. While 8-NBD-cGMP is almost nonfluorescent in aqueous solutions, it fluoresces strongly in hydrophobic environments such as in hydrophobic protein binding site. Excitation 475nm, Emission 525nm.

*BLG-N001-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

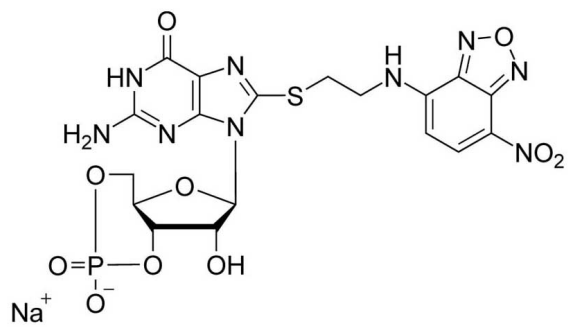
### Ordering Information

[Order Online »](#)

BLG-N001-01	1 $\mu$ mol
BLG-N001-05	5x1 $\mu$ mol

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	8-(2-[7-Nitro-4-benzofurazanyl]aminoethylthio)guanosine-3',5'-cyclic monophosphate . sodium salt
CAS	115993-88-9
Couple Type	Modified nucleotides
Formula	$C_{18}H_{17}N_9O_{10}PS \cdot Na$
MW	605.4
Purity	> 95% HPLC
Quantity	1µmol≈0.7mg
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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)