

# 8-Fluo-cGMP

## Fluorescent cyclic GMP analogue

Fluorescent cGMP analogue. Excitation 494nm, Emission 520 nm.  
Potent activator of cGMP-dependent protein kinase and cGMP-gated ion channels.

*BLG-F001-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

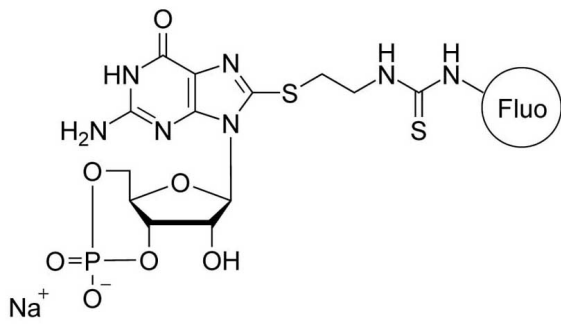
### Ordering Information

[Order Online »](#)

BLG-F001-01	1 $\mu$ mol
BLG-F001-05	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-[(Fluoresceinylthioureido)amino]ethyl]thio]guanosine-3',5'-cyclic monophosphate . sodium salt
Appearance	Crystallized or lyophilized sodium salt.
Couple Type	Modified nucleotides
Formula	$C_{33}H_{27}N_7O_{12}PS_2 \cdot Na$
MW	831.7
Purity	> 95% HPLC
Quantity	1µmol≈0.8mg
Solubility	Soluble in DMSO and ethanol.

**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](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)