

# Sp-2',3'-cGMPS

BLG-G071-25 (5 x 5  $\mu$ mol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

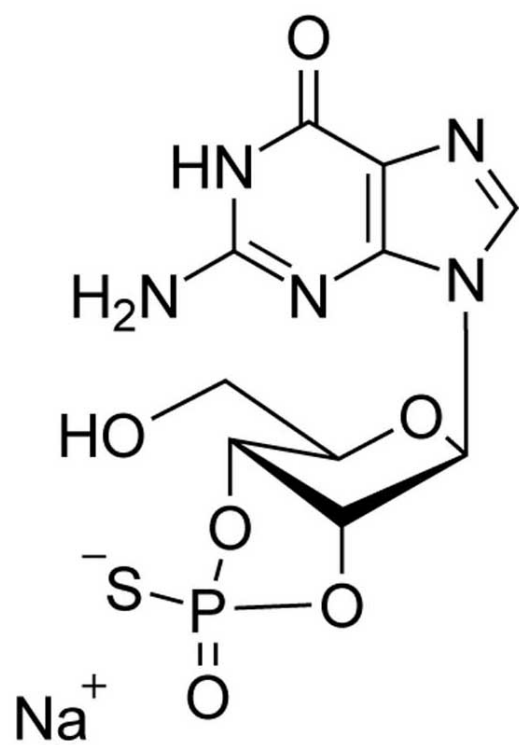
## Ordering Information

[Order Online »](#)

|             |               |
|-------------|---------------|
| BLG-G071-05 | 5 $\mu$ mol   |
| BLG-G071-25 | 5x5 $\mu$ 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** Guanosine- 2',3'-cyclic monophosphorothioate, endo / Sp- isomer . sodium salt

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{11}N_5O_6PS \cdot Na$

**MW** 383.3

**Purity** > 97% HPLC

**Quantity** 5µmol is equivalent to ~1.9 mg



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)