

# 2',2''-Di-O-Me-c-diGMP

2',2''-Di-O-Me-c-diGMP is an analogue of the bacterial second messenger c-diGMP (Cat. No. BLG-C057) which selectively binds the c-diGMP class II riboswitch over the class I riboswitch. It may be useful for discriminating between these two classes of c-diGMP-binding RNAs and for studying RNA-based c-diGMP signalling.

*BLG-D136-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

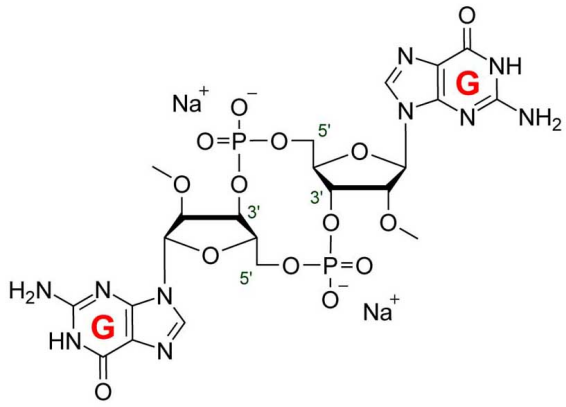
## Ordering Information

[Order Online »](#)

BLG-D136-001	0.1µmol
BLG-D136-005	5x0.1µmol

## 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** 2', 2"-O-(Dimethyl)-cyclic diguanosine monophosphate . sodium salt

**CAS** 849214-02-4

**Couple Type** Modified nucleotides

**Formula**  $C_{22}H_{28}N_{10}O_{14}P_2$  (free acid)

**MW** 718.5 (free acid)

**Purity** >95% HPLC

**Quantity** 0.1µmol≈72µg

**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)