

c-di-2'-d-2''-O-Me-GMP

Analogue of the bacterial second messenger
c-diGMP

c-di-2'-d-2''-O-Me-GMP is an analogue of the bacterial second messenger c-diGMP (Cat. No. BLG-C057) with modified ribose 2'-hydroxy positions.

BLG-D142-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

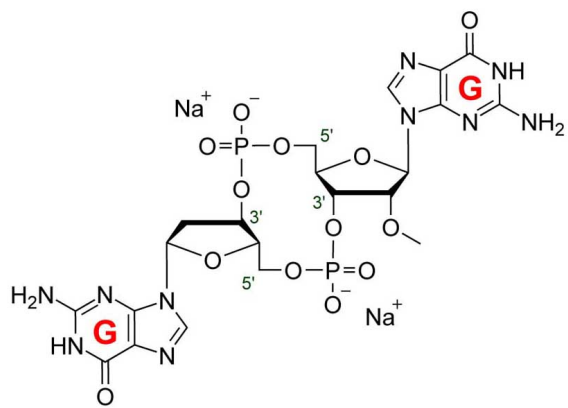
Ordering Information

[Order Online »](#)

BLG-D142-001	0.1µmol
BLG-D142-005	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 c-(2'-dGp-2'-O-Me-Gp), 2'-Deoxy-2''-O-methyl-cyclic diguanosine monophosphate . sodium salt

Appearance Solid

Formula $C_{21}H_{26}N_{10}O_{13}P_2$ (free acid)

MW 688.4 (free acid)

Purity > 95% HPLC



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