

c[C(3',5')pG(3',5')p]

Mixed purine/pyrimidine c-diNMP analogue

Analogue of the bacterial second messenger c-diGMP (Prod. No. BLG-C057) with mixed purine/pyrimidine nucleobases.

BLG-C199-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

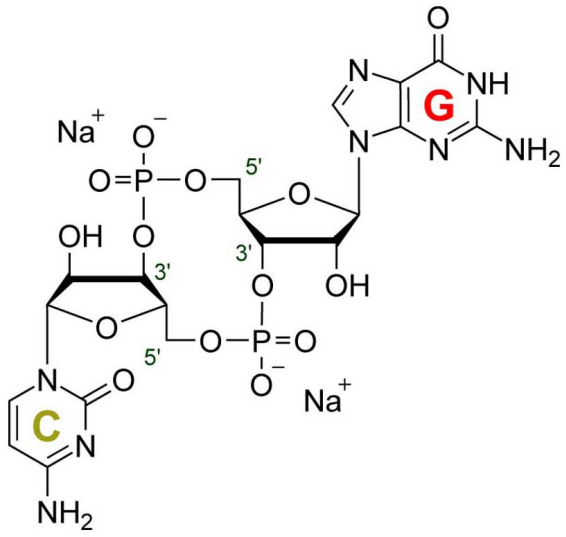
Ordering Information

[Order Online »](#)

BLG-C199-025	5x0.5µmol
--------------	-----------

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(CpGp), Cyclic (cytidine monophosphate-guanosine monophosphate) . sodium salt

Appearance Solid

CAS 132182-21-9

Formula $C_{19}H_{24}N_8O_{14}P_2$ (free acid)

MW 650.4 (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