

c-(GpUp) / cGMP-UMP

Bacterial hybrid purine-pyrimidine cyclic dinucleotide

c-(GpUp) is a hybrid purine-pyrimidine cyclic dinucleotide that was found to be biosynthesized as a minor product of alternative cyclic dinucleotide synthases identified in diverse bacteria.

BLG-C371-025 (5 x 0.5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

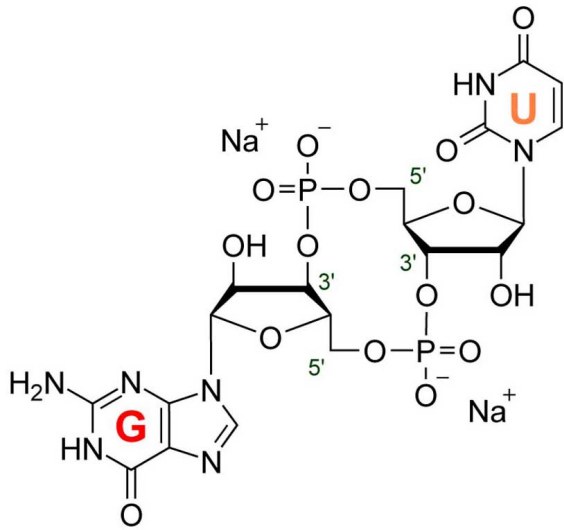
Ordering Information

[Order Online »](#)

BLG-C371-025	5x0.5 μ 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 Cyclic (guanosine monophosphate-uridine monophosphate), c[G(3',5')pU(3',5')]

CAS 232933-52-7

Couple Type Modified nucleotides

Formula $C_{19}H_{23}N_7O_{15}P_2$ (free acid)

MW 651.1 (free acid)

Purity > 95% HPLC

Quantity 0.5 μ mol \approx 326 mg



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