

c-(CpUp) / cCMP-UMP

Pyrimidine cyclic dinucleotide

c-(CpUp) is a pyrimidine cyclic dinucleotide that was found to be biosynthesized by an alternative cyclic dinucleotide synthase identified in *Legionella pneumophila*, LpCdnE02. c-(CpUp) is a minor product of LpCdnE02 while c-diUMP (Cat. No. BLG-C256) is the major product of this cyclic dinucleotide synthase.

BLG-C375-025 (5 x 0.5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

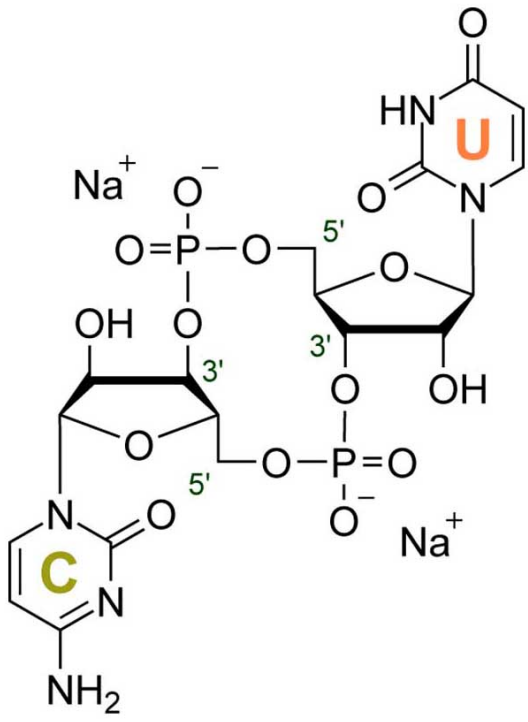
Ordering Information

[Order Online »](#)

BLG-C375-025	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 (cytidine monophosphate-uridine monophosphate), c[C(3',5')pU(3',5')]

Couple Type Modified nucleotides

Formula $C_{18}H_{23}N_5O_{15}P_2$ (free acid)

MW 611.4 (free acid)

Purity > 95% HPLC

Quantity 0.5 μ mol \approx 306 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