

cUMP is a cyclic nucleotide with a pyrimidine nucleobase.

*BLG-U001-250 (5 x 50 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

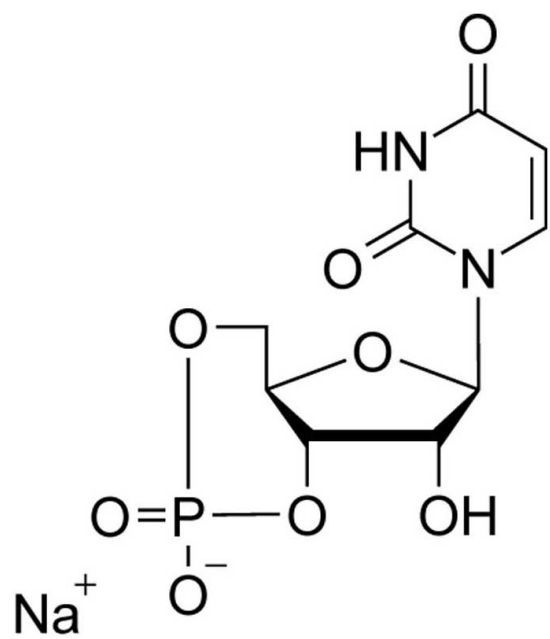
Ordering Information

[Order Online »](#)

BLG-U001-50	50 μ mol
BLG-U001-250	5x50 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 Uridine-3',5'-cyclic monophosphate . sodium salt

CAS 56632-58-7

Couple Type Modified nucleotides

Formula $C_9H_{10}N_2O_8P \cdot Na$

MW 328.2

Purity > 98% HPLC

Quantity 50µmol≈~16mg

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

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