

c[A(2',5')pU(3',5')p]

Hybrid purine-pyrimidine cyclic dinucleotide

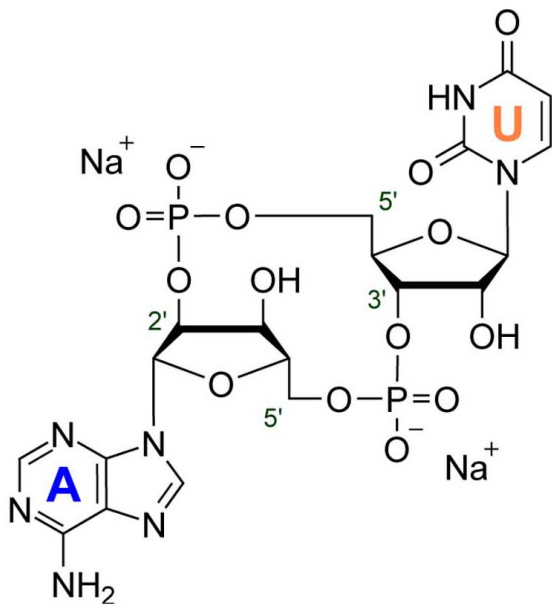
c[A(2',5')pU(3',5')p] is a hybrid purine-pyrimidine cyclic dinucleotide and stereoisomer of the second messengers c-(ApUp) (Cat. No. C 357) and c[U(2',5')pA(3',5')p] (Cat. No. C 399) which were found to be biosynthesized in bacteria and oyster species, respectively.

*BLG-C398-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

Ordering Information [Order Online »](#)

BLG-C398-01	1 µmol
BLG-C398-05	5x1 µmol

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 (adenosine-(2'→5')-monophosphate-uridine-(3'→5')-monophosphate) . sodium salt, 2'3'-cAUMP, 2',5'-3',5'-cAUMP

Couple Type Modified nucleotides

Formula $C_{19}H_{23}N_7O_{14}P_2$ (free acid)

MW 635.4 (free acid)

Purity > 95% HPLC

Quantity 0.5µmol is equivalent to ~318 µg

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-fr@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