

ApApApA(2',3'-cp)

Intermediate in the degradation of the bacterial second messenger c-tetraAMP

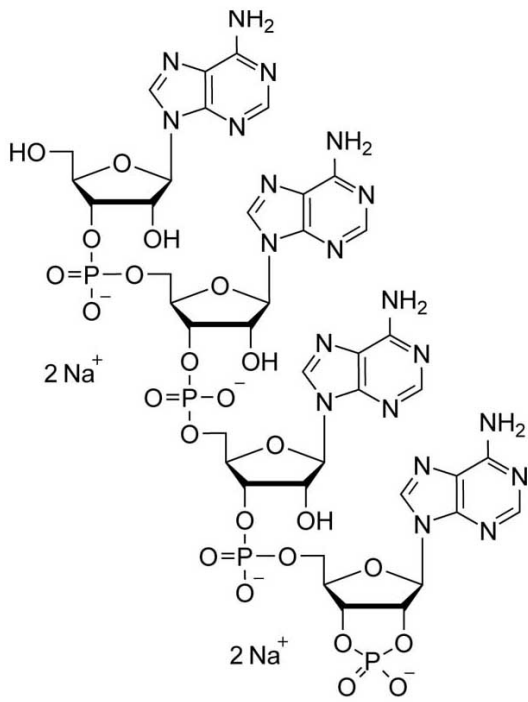
ApApApA(2',3'-cp) is considered to be an intermediate in the degradation of the bacterial second messenger c-tetraAMP (Biolog Cat. No. C 335) by ring nucleases. ApApApA(2',3'-cp) was also found to have activating effects on Csm6/Csx1-family ribonucleases.

BLG-A373-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-A373-001	0.1µmol
BLG-A373-005	5x0.1µmol

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 Adenylyl-(3'→5')-adenylyl-(3'→5')-adenylyl-(3'→5')-adenosine-2',3'-cyclic monophosphate . sodium salt, A4>P, ApApApA>p

Appearance Crystallized or lyophilized solid

CAS 62222-55-3

Couple Type Modified nucleotides

Formula $C_{40}H_{48}N_{20}O_{24}P_4$ (free acid)

MW 1316.8

Purity > 95% HPLC

Quantity 0.1 μmol are equivalent to ~0.1 mg

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