

ApApA(2',3'-cp)

Degradation product of the bacterial second messenger c-hexaAMP

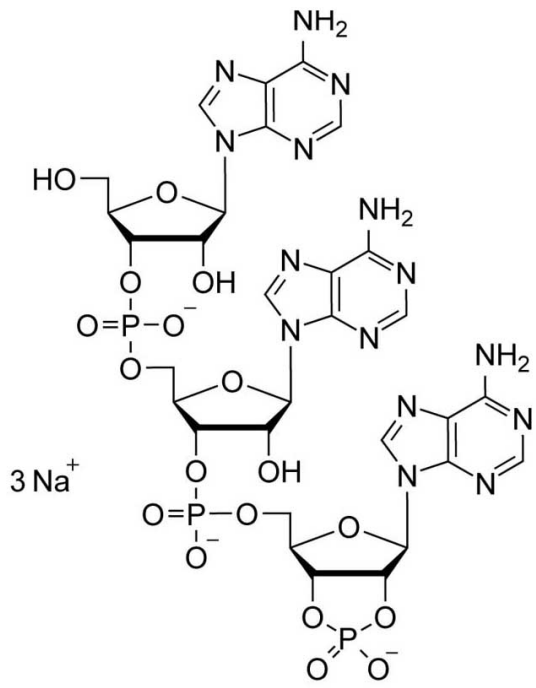
ApApA(2',3'-cp) is considered to be the product of the degradation of the bacterial second messenger c-hexaAMP (Biolog Cat. No. C 332) by certain Csm6 enzymes that degrade their own activator, thereby providing a self-limiting regulation for the c-hexaAMP signalling.

BLG-A425-005 (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-A425-001	0.1µmol
BLG-A425-005	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')-adenosine-2',3'-cyclic monophosphate . sodium salt, A3>P, ApApA>p

Appearance Crystallized or lyophilized solid

CAS 4527-94-0

Couple Type Modified nucleotides

Formula $C_{30}H_{36}N_{15}O_{18}P_3$ (free acid)

MW 987.6

Purity > 95% HPLC

Quantity 0.1 μmol are equivalent to ~99 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