

ApA(2',3'-cp)

Degradation product of the bacterial second messenger c-tetraAMP

ApA(2',3'-cp) is considered to be the product of the degradation of the bacterial second messenger c-tetraAMP (Biolog Cat. No. C 335) by ring nucleases via the intermediate ApApApA(2',3'-cp) (Biolog Cat. No. A 373).

BLG-A399-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

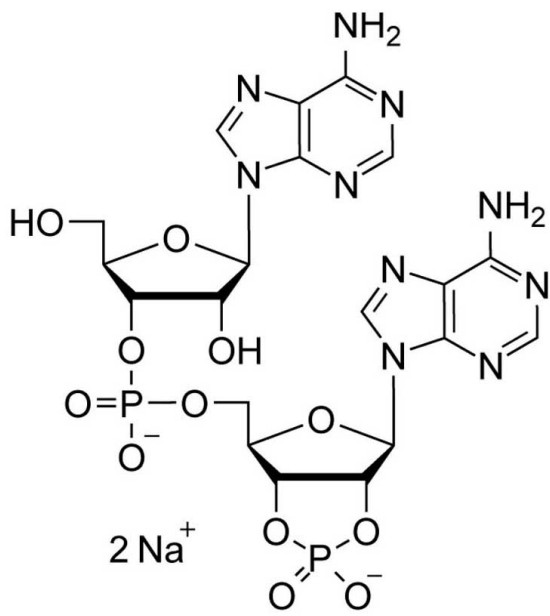
Ordering Information

[Order Online »](#)

BLG-A399-005	0.5µmol
BLG-A399-025	5x0.5µmol

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')-adenosine-2',3'-cyclic monophosphate . sodium salt, A2>P, ApA>p

Appearance Crystallized or lyophilized solid

CAS 4527-93-9

Couple Type Modified nucleotides

Formula $C_{20}H_{24}N_{10}O_{12}P_2$ (free acid)

MW 658.4

Purity > 95% HPLC

Quantity 0.5µmol are equivalent to ~0.3 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