

ApA

Adenylyl-(3→5')-adenosine (ApA) is a dinucleoside in which the 3'-hydroxy group of the adenosine moiety is connected to the 5'-hydroxy group of the other adenosine moiety by a phosphodiester bond.

*BLG-A083-25 (5 x 5 μmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

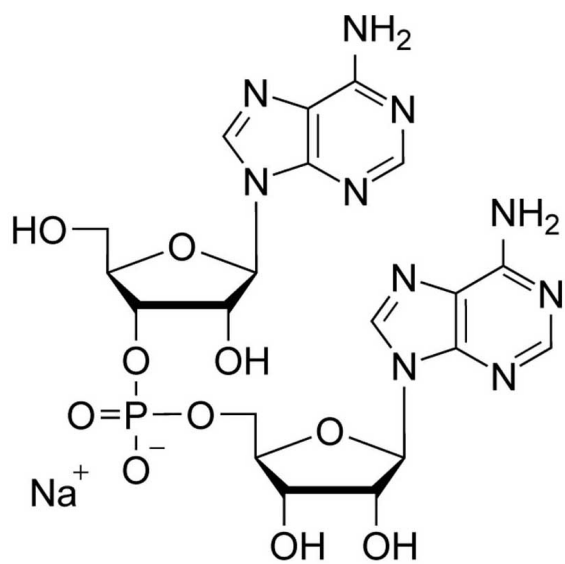
Ordering Information

[Order Online »](#)

BLG-A083-05	5μmol
BLG-A083-25	5x5μmol

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 . sodium salt

CAS 2391-46-0

Couple Type Modified nucleotides

Formula $C_{20}H_{25}N_{10}O_{10}P$ (free acid)

MW 596.5

Purity > 95% HPLC

Reconstitution Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

Solubility Very well soluble in water or aqueous buffers.

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