

pA(2',5')pA

Potential metabolite of 2' → 5'-linked oligoadenylates (2-5As)

pA(2',5')pA is a potential metabolite of 2' → 5'-linked oligoadenylates (2-5As).

BLG-P112-05 (5 x 1 μmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

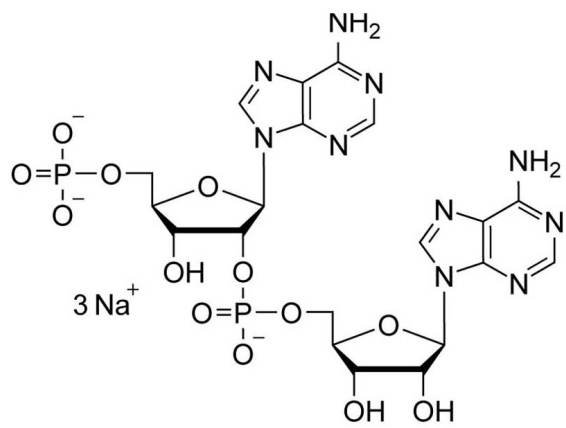
Ordering Information

[Order Online »](#)

BLG-P112-01	1 μmol
BLG-P112-05	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 5'-Phosphoadenylyl-(2'→5')-adenosine . sodium salt, 5'-p(rA)(2',5')p(rA)

Appearance Solid

CAS 20307-28-2

Couple Type Modified nucleotides

Formula $C_{20}H_{26}N_{10}O_{13}P_2$ (free acid)

MW 676.5 (free acid)

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.7 mg



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