

pppA(2',5')pA

In pppA(2',5')pA a 5'-ATP unit is connected with a 5'-AMP unit via a 2' → 5' linkage to form a linear dinucleotide. pppA(2',5')pA is considered to be a linear intermediate in the enzymatic production of 2' → 5'-linked oligoadenylates by oligoadenylate synthase 1 (OAS1).

*BLG-T073-05 (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-T073-01	1 μmol
BLG-T073-05	5x1 μmol

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5'-Triphosphoadenylyl-(2'→5')-adenosine

CAS 65954-94-1

Couple Type Modified nucleotides

Formula $C_{20}H_{28}N_{10}O_{19}P_4$ (free acid)

Formulation 10 mM aqueous solution of pH 7.6

MW 836.4 (free acid)

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.8 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