

pppA(2',5')pG

In pppA(2',5')pG a 5'-ATP unit is connected with a 5'-GMP unit via a 2' → 5' linkage to form a linear dinucleotide. pppA(2',5')pG is the linear intermediate in the two-step enzymatic production of 3',2'-cGAMP (Cat. No. C 238) by the *Drosophila melanogaster* enzyme cGLR1 in response to dsRNA.

BLG-T074-05 (5 x 1 μmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

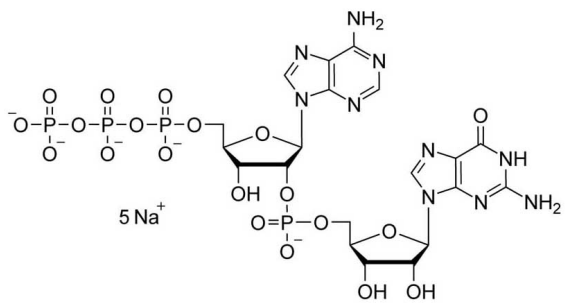
Ordering Information

[Order Online »](#)

BLG-T074-01	1 μmol
BLG-T074-05	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')-guanosine

Couple Type Modified nucleotides

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

Formulation 10 mM aqueous solution of pH 7.6

MW 852.4

Purity > 95% HPLC

Quantity 1µmol is equivalent to ~0.9 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