

GpA can be useful in kinetic studies on RNA polymerases or ribonucleases. GpA is a metabolite of the linear RNA dinucleotide pGpA (Cat. No. P 120) and GpAp (Cat. No. G 056).

BLG-G040-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

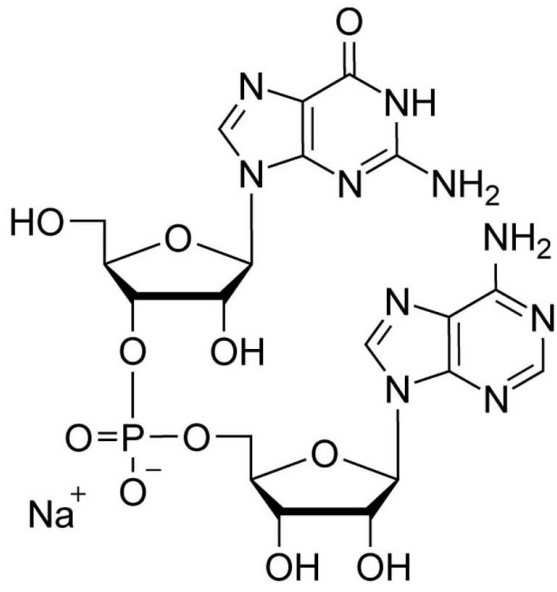
Ordering Information

[Order Online »](#)

| | |
|--------------------|---------------|
| BLG-G040-01 | 1 μ mol |
| BLG-G040-05 | 5x1 μ mol |

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 Guanylyl-(3'→5')-adenosine . sodium salt

CAS 103192-56-9

Couple Type Modified nucleotide

Formula $C_{20}H_{25}N_{10}O_{11}P$ (free acid)

MW 612.5

Purity > 95% HPLC

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