

pG-3'-p / pGp

Nucleoside bisphosphate suitable for the characterisation of ribonucleases and nucleotidases

pG-3'-p is a nucleoside bisphosphate that is suitable for characterising ribonucleases, nucleotidases and other nucleotide-binding enzymes. pG-3'-p has been successfully employed for the crystallization of ribonuclease T1.

BLG-G081-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

[Order Online »](#)

BLG-G081-50	5x10µmol
-------------	----------

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 Guanosine-3',5'-O-bisphosphate . sodium salt

Appearance Solid

CAS 110972-29-7

Couple Type Modified nucleotides

Formula $C_{10}H_{15}N_5O_{11}P_2$ (free acid)

MW 443.2 (free acid)

Purity > 95 % HPLC

Quantity 10 µmol are equivalent to ~4.4 mg

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