

mant-GDP, S pack

Ordering Information

[Order Online »](#)

JBS-NU-204S	150µl
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2'/3'-O-(N-Methyl-anthraniloyl)-guanosine-5'-diphosphate, Triethylammonium salt

Appearance colorless to slightly yellow solution

Application Notes Conformational dynamic: DnaB/C-proteins, Csk
Fluorescence stop-flow kinetics: release factor, eIF5B, EF-1B, EF-Tu
FRET: AC

CAS 148821-02-7

Couple Type Modified nucleotides

Formula $C_{18}H_{22}N_6O_{12}P_2$ (free acid)

Formulation solution in water

MW 576.35 (free acid)

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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)