

GTP – Solid – Purity

85 % (100 g)

Ordering Information

[Order Online »](#)

JBS-NU-1047-100G	100g
------------------	------

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 Guanosine 5'-triphosphate, Disodium salt trihydrate

Appearance colorless to slightly yellow solid

Application Notes Assembly of ribosomal units. Microdomain formation by small GTPases. Antiviral activity of large GTPases (dynamin superfamily). Regulation of exocytosis by Rho GTPases. Mechanism of hydrolysis by ADP-ribosylation factors.

CAS 56001-37-7

Couple Type Modified nucleotides

Formula $C_{10}H_{16}N_5O_{14}P_3$ (free acid)

MW 523.18 (free acid)

Purity ≥ 85 % (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)