

NTP Bundle, S pack

4x 100 mM (ATP, CTP, GTP, UTP)

Ordering Information

[Order Online »](#)

JBS-NU-1014S	4x200µl
--------------	---------

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 Adenosine-5'-triphosphate, Sodium salt; Cytidine-5'-triphosphate, Sodium salt; Guanosine-5'-triphosphate, Sodium salt; Uridine-5'-triphosphate, Sodium salt

Appearance clear solution

Couple Type Modified nucleotides

Formula
ATP: $C_{10}H_{16}N_5O_{13}P_3$ (free acid)
CTP: $C_9H_{16}N_3O_{14}P_3$ (free acid)
GTP: $C_{10}H_{16}N_5O_{14}P_3$ (free acid)
UTP: $C_9H_{15}N_2O_{15}P_3$ (free acid)

Formulation aqueous solution

MW ATP: 507.18 (free acid) CTP: 483.16 (free acid) GTP: 523.18 (free acid) UTP: 484.14 (free acid)

Purity ≥ 99 % (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-fr@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