

dNTP Mix incl. dUTP, L pack

Premix of 10 mM dATP, dCTP, dGTP and 20 mM dUTP

Ordering Information

[Order Online »](#)

JBS-NU-1020L	5x1ml
--------------	-------

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'-Deoxyadenosine 5'-triphosphate, Sodium salt 2'-Deoxycytidine 5'-triphosphate, Sodium salt 2'-Deoxyguanosine 5'-triphosphate, Sodium salt 2'-Deoxyuridine 5'-triphosphate, Sodium salt

Appearance clear solution

Couple Type Modified nucleotides

Formula
dATP: $C_{10}H_{16}N_5O_{12}P_3$ (free acid)
dCTP: $C_9H_{16}N_3O_{13}P_3$ (free acid)
dGTP: $C_{10}H_{16}N_5O_{13}P_3$ (free acid)
dUTP: $C_9H_{15}N_2O_{14}P_3$ (free acid)

Formulation aqueous solution

MW dATP: 491.18 (free acid) dCTP: 467.15 (free acid) dGTP: 507.18 (free acid) dUTP: 468.14 (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