

# dNTP Mix / 25 mM, L pack

Equimolar Mix of 25 mM dATP, dCTP, dGTP and dTTP

## Ordering Information

[Order Online »](#)

JBS-NU-1023L	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'-Deoxythymidine 5'-triphosphate, Sodium salt

**Appearance** clear solution

**Application Notes** For standard PCR applications a final concentration of 200 µM each dNTP is recommended.

**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)  
dTTP:  $C_{10}H_{17}N_2O_{14}P_3$  (free acid)

**Formulation** aqueous solution

**MW** dATP: 491.18 (free acid) dCTP: 467.15 (free acid) dGTP: 507.18 (free acid) dTTP: 482.17 (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](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)