

ddNTP bundle, S pack

4x 10 mM (ddATP, ddCTP, ddGTP, ddTTP)

Ordering Information

[Order Online »](#)

JBS-NU-1019S	4x100µl
--------------	---------

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'-Dideoxyadenosine-5'-triphosphate, Lithium salt; 2',3'-Dideoxycytidine-5'-triphosphate, Lithium salt; 2',3'-Dideoxyguanosine-5'-triphosphate, Lithium salt; 2',3'-Dideoxythymidine-5'-triphosphate, Lithium salt

Appearance colorless to slightly yellow solution

Couple Type Modified nucleotides

Formula
ddATP: $C_{10}H_{16}N_5O_{11}P_3$ (free acid)
ddCTP: $C_9H_{16}N_3O_{12}P_3$ (free acid)
ddGTP: $C_{10}H_{16}N_5O_{12}P_3$ (free acid)
ddTTP: $C_{10}H_{17}N_2O_{13}P_3$ (free acid)

Formulation solution in water

MW ddATP: 475.18 (free acid)ddCTP: 451.16 (free acid)ddGTP: 491.18 (free acid)ddTTP: 466.17 (free acid)

Purity ≥ 98 % (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-be@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