

ddCTP, L pack

Ordering Information

[Order Online »](#)

JBS-NU-1016L	5x100µl
--------------	---------

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'-Dideoxycytidine-5'-triphosphate, Lithium salt

Appearance colorless to slightly yellow solution

Application Notes
SNP: multiplexing by minisequencing
Multiplexed detection of alternatively spliced transcripts
Effect on reverse transcriptase of HIV-1
Molecular dynamic calculations of DNA-polymerase beta-ddCTP complex
Crystal structure with DNA- polymerase-I

CAS 66004-77-1

Concentration 10 mM

Couple Type Modified nucleotides

Formula $C_9H_{16}N_3O_{12}P_3$ (free acid)

Formulation solution in water

MW 451.16 (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