

C8-Alkyne-dCTP (L pack)

Ordering Information

[Order Online »](#)

JBS-CLK-T06-L	5x5µl
---------------	-------

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 5-(Octa-1,7-diynyl)-2'-deoxycytidine 5'-triphosphate, Sodium salt

Appearance colorless to slightly yellow solution

Application Notes Incorporation into DNA by PCR with family B polymerases, Incorporation into DNA by Primer Extension with family A polymerase Taq and family B polymerases , Incorporation into DNA by 3'-End Labeling with Terminal deoxynucleotidyl Transferase(TdT)

Couple Type Modified nucleotides

Formula $C_{17}H_{24}N_3O_{13}P_3$ (free acid)

Formulation solution in water

MW 571.31 (free acid)

Purity > 95 %

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