

3TCTP, L pack

Ordering Information

[Order Online »](#)

JBS-NU-1606L	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 Lamivudine triphosphate, Sodium Salt (L-Isomer), 2', 3'-Dideoxy-3'-thia-cytidine-5'-triphosphate, Sodium salt (L isomer)

Appearance colorless to slightly yellow solution

Application Notes Monitoring in mononuclear cells by ion-pair LC and MS
Influence on incorporation efficiency
Inhibition of DNA/RNA-dependent DNA-polymerase

CAS 143188-53-8

Couple Type Modified nucleotides

Formula $C_8H_{14}N_3O_{12}P_3S$ (free acid)

Formulation solution in water

MW 469.19 (free acid)

Purity ≥ 95 % (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-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com