

5F-UTP, L pack

Ordering Information

[Order Online »](#)

JBS-NU-1115L	5x10µl
--------------	--------

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 5F-UTP, 5-Fluoro-uridine-5'-triphosphate, Sodium salt

Appearance colorless to slightly yellow solution

Application Notes Inhibition of viral application
Mutagenic activity
Conformational studies (NMR-NOE) of HIV-2 TAR-RNA
Use as NMR-label (F)[2] X-ray of viral RNA polymerase

CAS 3828-96-4

Couple Type Modified nucleotides

Formula $C_9H_{14}N_2O_{15}P_3F$ (free acid)

Formulation solution in water

MW 502.13 (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