

# 5-Methyl-UTP, L pack

## Ordering Information

[Order Online »](#)

|             |              |
|-------------|--------------|
| JBS-NU-880L | 5x10 $\mu$ l |
|-------------|--------------|

## 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** m5UTP, 5-Methyluridine-5'-O-triphosphate, Sodium salt, 5-Me-UTP

**Appearance** colorless to slightly yellow solution

**CAS** 23198-01-8 (free acid), 1801968-79-5 (Tetrasodium salt)

**Couple Type** Modified nucleotides

**Formula**  $C_{10}H_{17}N_2O_{15}P_3$  (free acid)

**Formulation** solution in water

**MW** 498.17 (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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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