

# HighYield T7 mRNA Synthesis Kit (m1ATP), S pack

## Synthesis of N1-methyladenosine-modified (m)RNA

HighYield T7 mRNA Synthesis Kit (N1-Methyl-ATP) is designed to produce large amounts of N1-methyladenosine-modified (m)RNA via in vitro transcription with T7 RNA polymerase. The internally modified (m)RNA can subsequently be used for microinjection, transfection or in vitro translation experiments. N1-methyladenosine modifications have been shown to increase (m)RNA stability and to reduce immunogenicity. A 5'-cap structure further increases translation efficiency. 5'-capping can be achieved by a) co-transcriptional capping with cap analogs (e.g. ARCA) or b) post-transcriptional capping with capping enzymes. The kit contains sufficient reagents for 15 reactions (S-Pack) or 50 reactions (L-Pack) à 20 µl (7.5 mM GTP, 7.5 mM UTP, 7.5 mM CTP, 3.75 mM ATP, 3.75 mM N1-Methyl-ATP). An individual optimization of N1-Methyl-ATP concentration can easily be achieved with the single nucleotide format. A 20 µl reaction yields about 80-100 µg RNA after 30 min incubation (1 µg T7 control template, 1.4 kb RNA transcript). Yields may however vary depending on the template (promotor design, sequence length, secondary structure formation)).

### Ordering Information

[Order Online »](#)

JBS-RNT-113-S	15Reactions
---------------	-------------

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping                      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

**Technical Info / Product**      For the Original Manufacturer's data sheet please [click here](#).

**Notes**



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)