

# HighYield T7 P&L RNA Synthesis Kit

RNA synthesis via in vitro transcription with a modified T7 RNA Polymerase (P266L)

HighYield T7 P&L RNA Synthesis Kit is designed to produce large amounts of RNA via in vitro transcription with a modified T7 RNA Polymerase. Proline 266 has been replaced by leucine (P266L) which has been associated with decreased abortive transcription[1].increased 5' homogeneity of transcripts synthesized from A-initiating phi2.5 promoter[2]increased 5' incorporation efficiency of GTP analogs[3] The resulting RNA can subsequently be used for a variety of applications such as microinjection, transfection or in vitro translation experiments, structure/function studies or probes for in situ hybridization and Northern Blot experiments. One kit contains sufficient reagents for 50 reactions of 20 µl each (7.5 mM each NTP). A 20 µl reaction yields about 140-160 µg RNA after 30 min incubation (1 µg T7 control template, 1.4 kb RNA transcript). Yields may however vary depending on the template (promoter design, sequence length, secondary structure formation).

## Ordering Information

[Order Online »](#)

JBS-RNT-201	50Reactions
-------------	-------------

## 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)