

T7 P&L RNA Polymerase HC

**Highly concentrated and modified T7 RNA
Polymerase for set-up of high yield in vitro
transcription reactions**

Bacteriophage T7 RNA Polymerase is a DNA-dependent RNA polymerase (99 kDa) that catalyzes in vitro RNA synthesis from a DNA sequence containing a T7 phage promoter. The highly concentrated version and modified T7 RNA Polymerase HC is ideally suited for set-up and optimization of high yield in vitro transcription reactions. Proline 266 has been replaced by leucine (P266L) which has been associated with decreased abortive transcription[1], increased 5' homogeneity of transcripts synthesized from A-initiating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 provided T7 IVT Set-up Buffer, MgCl₂ and DTT solutions enable testing of different reaction conditions such as nucleotide/ MgCl₂ concentrations. Recommendation for assay set-up with minimal optimization: HighYield T7 P&L RNA Synthesis Kit (#RNT-201). Activity: 1 µl T7 P&L RNA Polymerase HC generates \geq 30 µg RNA in 10 min at 37°. This is comparable to 1000 U of NEB's T7 RNA Polymerase (High Concentration), #M0460T (activity assay conditions: 1x T7 IVT Set-up Buffer supplemented with 10 mM DTT, 12 mM MgCl₂, 3 mM NTPs, 1 µl T7 RNA Polymerase HC and 1 µg DNA template (1.4 kpb RNA transcript) in 20 µl. RNA yield was determined with a fluorescence microplate assay.)

Ordering Information

[Order Online »](#)

JBS-RNT-018

50µl

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance liquid

Purity ≥ 95 % (SDS-PAGE)

Quantity 2 µg/µl

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-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

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

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