

HighYield T7 RNAi Kit

dsRNA synthesis for RNAi applications via
T7 RNA Polymerase-mediated in vitro
transcription

HighYield T7 RNAi Kit is designed to produce large amounts of double-stranded RNA (dsRNA) > 200 bp via T7 RNA Polymerase-mediated in vitro transcription. The resulting dsRNA can subsequently be used for RNA interference (RNAi) applications. Preparation of small interfering dsRNA, known as siRNA (approximately 20 bp long) is principally feasible as well however, we recommend a chemical siRNA synthesis approach for short dsRNAs. The kit contains sufficient reagents for 50 reactions of 20 µl each (7.5 mM each NTP). A 20 µl reaction yields > 40 µg dsRNA (> 2 mg/ml) after 30 min incubation (1 µg each T7 control template I & II, 0.5 kbp RNA transcripts). Yields may however vary depending on the template (promotor design, sequence length, secondary structure formation).

Ordering Information

[Order Online »](#)

JBS-RNT-134	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)