

## HiScribe™ T7 Quick High Yield RNA Synthesis Kit NEB #E2050S/L

50/250 reactions  
Version 7.0\_12/25

### Table of Contents

Introduction.....	<a href="#">2</a>
DNA Template Preparation .....	<a href="#">3</a>
Plasmid Templates .....	<a href="#">3</a>
PCR Templates .....	<a href="#">3</a>
Synthetic DNA Oligonucleotides .....	<a href="#">3</a>
RNA Synthesis Protocols .....	<a href="#">4</a>
Standard RNA Synthesis .....	<a href="#">5</a>
RNA Synthesis with Modified Nucleotides.....	<a href="#">6</a>
Capped RNA Synthesis.....	<a href="#">7</a>
Purification of Synthesized RNA .....	<a href="#">9</a>
LiCl Precipitation.....	<a href="#">9</a>
Phenol:Chloroform Extraction and Ethanol Precipitation .....	<a href="#">9</a>
Spin Column Chromatography .....	<a href="#">10</a>
Gel Purification .....	<a href="#">10</a>
Evaluation of Reaction Products .....	<a href="#">10</a>
Quantification by UV Light Absorbance .....	<a href="#">10</a>
Analysis of Transcription Products by Gel Electrophoresis .....	<a href="#">10</a>
Troubleshooting .....	<a href="#">11</a>
Control Reaction .....	<a href="#">11</a>
Low Yield of Full-length RNA .....	<a href="#">11</a>
Addition of DTT .....	<a href="#">11</a>
Low Yield of Short Transcript .....	<a href="#">11</a>
RNA Transcript Smearing on Denaturing Gel .....	<a href="#">11</a>
RNA Transcript of Larger Size than Expected .....	<a href="#">11</a>
RNA Transcript of Smaller Size than Expected .....	<a href="#">11</a>
Ordering Information .....	<a href="#">12</a>
Revision History .....	<a href="#">13</a>

## The HiScribe T7 Quick High Yield RNA Synthesis Kit Includes:

All kit components should be stored at  $-20^{\circ}\text{C}$ . Each kit contains sufficient reagents for 50 x 20  $\mu\text{l}$  reactions (S size) or 250 x 20  $\mu\text{l}$  reactions (L size). Each standard reaction yields up to 180  $\mu\text{g}$  of unmodified RNA from 1  $\mu\text{g}$  control template.

NTP Buffer Mix (20 mM each NTP)

T7 RNA Polymerase Mix

FLuc Control Template (0.5  $\mu\text{g}/\mu\text{l}$ )

DNase I (2 units/ $\mu\text{l}$ )

LiCl Solution (7.5 M LiCl, 10 mM EDTA)

Dithiothreitol (DTT) (0.1 M)

## Required Materials Not Included:

DNA Template: The DNA template must be linear and contain the T7 RNA Polymerase promoter with correct orientation in relation to target sequence to be transcribed.

Cap Analogs: Dinucleotide Cap Analogs: NEB #S1411, #S1405, #S1406 and #S1407

Modified-NTP: N1-Methyl-Pseudouridine-5'-Triphosphate (NEB #N0431)  
5-Methyl-Cytidine-5'-Triphosphate (NEB #N0432)  
Pseudouridine-5'-Triphosphate (NEB #N0433)  
5-Methoxy-Uridine-5'-Triphosphate (NEB# N0434)  
Biotin-, Fluorescein-, Digoxigenin-, or Aminoallyl-NTP

General:  $37^{\circ}\text{C}$  incubator or thermocycler, nuclease-free water

DNase: DNase I (RNase-free) (NEB #M0303) or DNase I-XT (NEB #M0570)

Purification: Buffer- or water-saturated phenol/chloroform, ethanol and 3M sodium acetate, pH 5.2, or Monarch<sup>®</sup> RNA Spin Cleanup Kit (50, or 500  $\mu\text{g}$  capacity) (#T2040, or #T2050)

Gel Analysis: Gels and running buffers, gel apparatus, power supply

## Introduction

The HiScribe T7 Quick High Yield RNA Synthesis Kit is designed for quick set-up and production of large amounts of RNA *in vitro*. The reaction can be set up conveniently by combining the NTP buffer mix, T7 RNA Polymerase mix, and a suitable DNA template. The kit also allows for capped RNA or partially modified RNA synthesis by incorporation of cap analog (ARCA, NEB #S1411) or base-modified nucleotides. RNA synthesized with the kit can be used for RNA structure and function studies, ribozyme biochemistry, as probes for RNase protection assays and hybridization-based blots, anti-sense RNA and RNAi experiments, microarray analysis and microinjection, as well as *in vitro* translation and RNA vaccines.

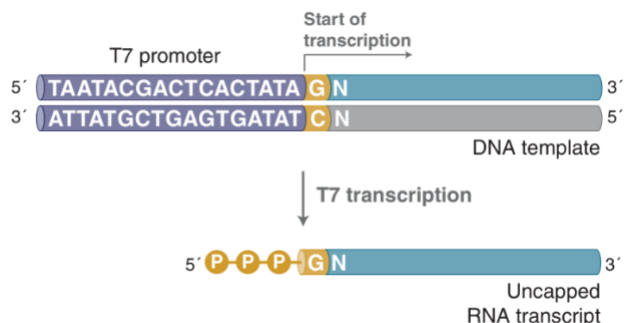
To synthesize high specific activity radioactive RNA probes or RNA with 100% substitution of one or more modified nucleotides we recommend using the T7 High Yield RNA Synthesis Kit (NEB #E2040), in which the four nucleotides are supplied separately.

The kit contains sufficient reagents for 50 or 250 reactions of 20  $\mu\text{l}$  each. Each standard reaction yields up to 180  $\mu\text{g}$  of RNA from 1  $\mu\text{g}$  control template. Each kit can yield up to 9 mg RNA (S size) or 45 mg RNA (L size). Reactions may be scaled linearly to increase yield proportional to reaction volume.

## DNA Template Preparation

Linearized plasmid DNA, PCR products or synthetic DNA oligonucleotides can be used as templates for *in vitro* transcription with the HiScribe T7 Quick High Yield RNA Synthesis Kit, provided that they contain a double-stranded T7 promoter region upstream of the sequence to be transcribed. Figure 1 illustrates the minimal T7 promoter sequence and the start of transcription as well as a run-off transcript after T7 transcription. The standard protocol results in 5' triphosphorylated RNA.

**Figure 1. Transcription by T7 RNA Polymerase**



### Plasmid Templates

Completely linearized plasmid template of highest purity is critical for successful use of the HiScribe T7 High Yield RNA Synthesis Kit. Quality of the template DNA affects transcription yield and the integrity of RNA synthesized. The highest transcription yield is achieved with the highest purity template. Plasmid purified by many laboratory methods can be successfully used, provided it contains mostly supercoiled form, and is free from contaminating RNase, protein, RNA, and salts.

To produce an RNA transcript of defined length, plasmid DNA must be completely linearized with a restriction enzyme, downstream of the insert to be transcribed. In contrast, circular plasmid templates will generate long heterogeneous RNA transcripts in higher quantities because of the high processivity of T7 RNA Polymerase. NEB has a large selection of restriction enzymes for this purpose; we recommend selecting restriction enzymes that generate blunt ends or 5'-overhangs.

After linearization, we recommend purifying the template DNA by phenol:chloroform extraction:

1. Extract DNA with an equal volume of 1:1 phenol:chloroform mixture, and repeat, if necessary.
2. Extract twice with an equal volume of chloroform to remove residual phenol.
3. Precipitate the DNA by adding 1/10th volume of 3 M sodium acetate, pH 5.2, and two volumes of ethanol. Incubate at  $-20^{\circ}\text{C}$  for at least 30 minutes.
4. Pellet the DNA in a microcentrifuge for 15 minutes at top speed. Carefully remove the supernatant.
5. Rinse the pellet by adding 500  $\mu\text{l}$  of 70% ethanol and centrifuging for 15 minutes at top speed. Carefully remove the supernatant.
6. Air dry the pellet and resuspend it in nuclease-free water at a concentration of 0.5–1  $\mu\text{g}/\mu\text{l}$ .

### PCR Templates

PCR products containing T7 RNA Polymerase promoter in the correct orientation can be transcribed. We recommend using Q5<sup>®</sup> Hot Start High-Fidelity DNA Polymerase (NEB #M0493/M0494). Though PCR mixture can be used directly, better yields will be obtained with purified PCR products. PCR products can be purified according to the protocol for plasmid restriction digests above, or by using commercially available spin columns (we recommend Monarch Spin PCR & DNA Cleanup Kit, NEB #T1130). PCR products should be examined on an agarose gel to estimate concentration and to confirm amplicon size prior to its use as a template in the HiScribe T7 Quick High Yield RNA Synthesis Kit. Depending on the PCR products, 0.1–0.5  $\mu\text{g}$  of PCR fragments can be used in a 20  $\mu\text{l}$  *in vitro* transcription reaction.

### Synthetic DNA Oligonucleotides

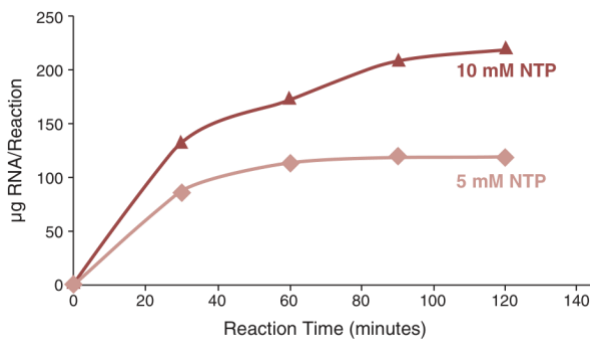
Synthetic DNA Oligonucleotides, which are either entirely double-stranded or mostly single-stranded, with a double-stranded T7 promoter sequence can be used in the T7 Quick transcription reaction. In general, the yields are relatively low and also variable depending upon the sequence, purity and preparation of the synthetic oligonucleotides. We do not recommend using synthetic oligonucleotides directly as template for IVT RNA greater than 100 nucleotides.

## RNA Synthesis Protocols

We strongly recommend wearing gloves and using nuclease-free tubes and reagents to avoid RNase contamination. Reactions are typically 20  $\mu$ l but can be scaled up linearly as needed. Reactions should be assembled in nuclease-free microfuge tubes or PCR strip tubes.

Figure 2 shows how the final yield is proportional to the amount of input nucleotides, so incorporation efficiency remains the same when different amounts of NTP are used; the amount of NTP Buffer Mix in a standard 20  $\mu$ l reaction can vary from 2 to 10  $\mu$ l. Figure 2 shows the time course of standard RNA synthesis from 1  $\mu$ g control DNA template coding for a 1.8 kb RNA transcript with the T7 Quick Kit using 10  $\mu$ l and 5  $\mu$ l NTP Buffer Mix in a 20  $\mu$ l reaction.

**Figure 2. RNA synthesis with different amounts of NTP**



*Time course of standard RNA synthesis from 1  $\mu$ g control DNA template coding for a 1.8 kb RNA transcript with the HiScribe T7 Quick Kit using 10  $\mu$ l and 5  $\mu$ l NTP Buffer Mix in a 20  $\mu$ l reaction. Reactions were incubated at 37°C in a thermocycler. Transcripts were purified by spin columns and quantified on a NanoDrop® Spectrophotometer.*

## Standard RNA Synthesis

1. Thaw the necessary components at room temperature. Keep the T7 RNA Polymerase Mix on ice.
2. Mix and pulse-spin in a microfuge to collect the solutions to the bottom of the tubes.
3. Set up the reaction at room temperature in the order listed in the table below:

COMPONENTS	20 $\mu$ l REACTION	FINAL AMOUNT
Nuclease-free Water	X $\mu$ l	
NTP Buffer Mix	10 $\mu$ l	10 mM each NTP
Template DNA	X $\mu$ l	1 $\mu$ g
DTT (0.1M)	1 $\mu$ l	5 mM
T7 RNA Polymerase Mix	2 $\mu$ l	

\* For reactions that synthesize short transcripts (< 0.3 kb), follow the general reaction above, except add an additional 10  $\mu$ l of nuclease-free water to bring the final reaction volume to 30  $\mu$ l. Increase template DNA up to 2  $\mu$ g.

4. Mix thoroughly by pipetting and pulse-spin in a microfuge. Incubate at 37°C for 2 hours in a dry air incubator or thermocycler to prevent evaporation. For reaction times of 60 minutes or less, a water bath or heating block may be used. The yield will not be compromised if the incubation temperature is within the range of 35–40°C.

Reactions for short RNA transcripts (< 0.3 kb) should be incubated for 4 hours or longer. It is safe to incubate the reaction for 16 hours (overnight).

*Optional:* Standard reactions normally generate large amounts of RNA at concentrations up to 10 mg/ml. As a result, the reaction mixture is quite viscous. It is easier to perform DNase treatment after the reaction mixture is diluted. To remove template DNA, add 30  $\mu$ l nuclease-free water and 2  $\mu$ l of DNase I (RNase-free) (NEB #M0303), mix, and incubate for 15 minutes at 37°C. Alternatively, 2  $\mu$ l of DNase I-XT (NEB# M0570) can be added directly to the IVT product and incubated for 15 minutes at 37°C.

5. Proceed with [purification of synthesized RNA](#) and/or [evaluation of transcription products](#) yield and/or length.

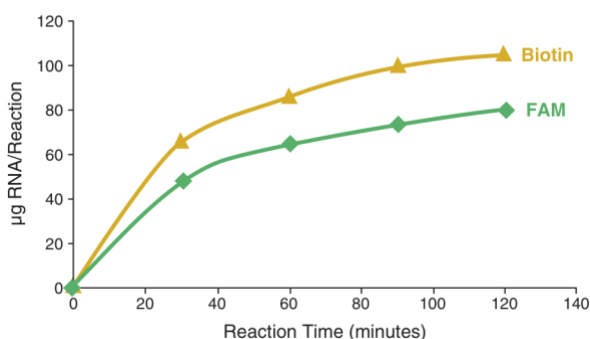
*Optional:* If a poly(A) tail is desired and is not encoded in the template plasmid, you can add one post-transcriptionally with *E. coli* Poly(A) Polymerase (NEB# M0276).

## RNA Synthesis with Modified Nucleotides

This protocol can be used to synthesize 5' triphosphorylated RNA containing partial substitutions of canonical nucleotides with base-modified nucleotides (not provided) from a DNA template containing the T7 RNA Polymerase promoter sequence immediately followed by guanosine. For complete modified nucleotide substitution, we recommended using the T7 High Yield RNA Synthesis Kit (NEB #E2040), in which the four nucleotides are supplied separately.

The ratio of modified nucleotide to standard nucleotide can be adjusted by varying the amount of the NTP Buffer Mix and modified nucleotide. The recommended molar ratio of user-supplied base-modified NTP (e.g. N1-Methyl-Pseudo-UTP, Biotin-, Fluorescein-, Digoxigenin-, or Aminoallyl-NTP) to standard NTP is 1/3. Modified ribonucleotides reduce transcription efficiency; therefore, lower transcription yields may be expected as compared to transcription using unmodified NTPs. In general, Biotin-NTP and Aminoallyl-NTP have an insignificant effect on yields, while lower yields can be expected for transcription reactions containing Fluorescein-NTP or Cy-NTP (Figure 3). In addition, transcripts containing modified ribonucleotides may have reduced electrophoretic mobility due to higher molecular weight.

**Figure 3. Labeled RNA synthesis with modified NTPs**



*Time course of labeled RNA synthesis using 1 µg control template with Biotin-16-UTP and Fluorescein-12-UTP according to the protocol below. Reactions were incubated at 37°C in a thermocycler. Transcripts were purified by spin columns and quantified on a NanoDrop Spectrophotometer.*

### Protocol

1. Thaw the necessary components at room temperature. Keep the T7 RNA Polymerase Mix on ice.
2. Mix and pulse-spin in a microfuge to collect the solutions to the bottom of the tubes.
3. Set up the reaction at room temperature in the order listed in the table below:

COMPONENTS	20 µl REACTION	FINAL AMOUNT
Nuclease-free water	X µl	
NTP Buffer Mix	5 µl	5 mM each NTP
10 mM Modified NTP	5 µl	2.5 mM
Template DNA	X µl	1 µg
DTT (0.1M)	1 µl	5 mM
T7 RNA Polymerase Mix	2 µl	

4. Mix thoroughly by pipetting and pulse-spin in a microfuge. Incubate at 37°C for 2 hours in a dry air incubator or thermocycler to prevent evaporation. Reactions for short RNA transcripts (< 0.3 kb) should be incubated for 4 hours or longer. It is safe to incubate the reaction for 16 hours (overnight).

*Optional:* To remove template DNA, add 30 µl nuclease-free water and 2 µl of DNase I (RNase-free) (NEB #M0303), mix, and incubate at 37°C for 15 minutes. Alternatively, 2 µl of DNase I-XT (NEB# M0570) can be added directly to the IVT product and incubated at 37°C for 15 minutes.

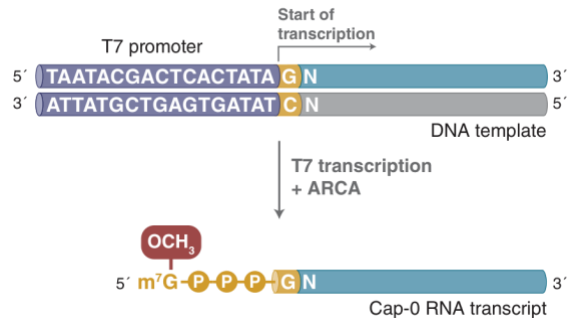
5. Proceed with [purification of synthesized RNA](#) and/or [evaluation of transcription products](#) yield and/or length.

*Optional:* If a poly(A) tail is desired and is not encoded in the template plasmid, you can add one post-transcriptionally with *E. coli* Poly(A) Polymerase (NEB# M0276).

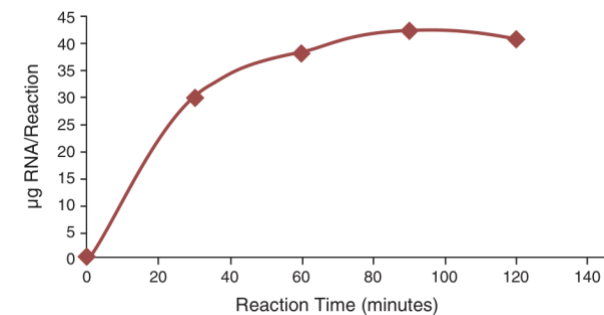
### Capped RNA Synthesis Using a Dinucleotide Cap Analog (i.e., ARCA)

This protocol can be used to synthesize Cap-0 RNA utilizing a dinucleotide cap analog, such as ARCA (NEB #S1411; Figure 3); cap analogs are sold separately. The recommended ratio of dinucleotide cap analog to GTP is 4:1, resulting in approximately 80% capped RNA transcripts. Increasing the ratio of cap analog to GTP will increase the proportion of capped RNA transcripts; however, it also significantly decreases the yield of the reaction. Figure 4 shows that most reactions will be complete in one hour and the yield per reaction is 30–40 µg RNA.

**Figure 3: T7 RNA Polymerase co-transcriptionally caps RNA with anti-reverse cap analog (ARCA)**



**Figure 4. Capped RNA Synthesis with ARCA**



*Time course of capped RNA synthesis from 1 µg control template. Reactions were incubated at 37°C in a thermocycler. Transcripts were purified by spin columns and quantified on a NanoDrop Spectrophotometer.*

#### Protocol

1. Thaw the necessary components at room temperature. Keep the T7 RNA Polymerase Mix on ice.
2. Mix and pulse-spin in a microfuge to collect the solutions to the bottom of the tubes.
3. Prepare a 40 mM solution of dinucleotide cap analog.
4. Set up the reaction at room temperature in the order listed in the table below:

COMPONENTS	20 µl REACTION	FINAL AMOUNT
Nuclease-free Water	X µl	
NTP Buffer Mix	2 µl	2 mM each NTP
40 mM Dinucleotide Cap Analog (NEB #S1411, S1405, S1406 or S1407)	4 µl	8 mM
Template DNA	X µl	1 µg
DTT (0.1M)	1 µl	5 mM
T7 RNA Polymerase Mix	2 µl	

5. Mix thoroughly by pipetting and pulse-spin in a microfuge. Incubate at 37°C for 2 hours in a dry air incubator or thermocycler to prevent evaporation.

*Optional:* To remove template DNA, add 30 µl nuclease-free water and 2 µl of DNase I (RNase-free) (NEB #M0303), mix, and incubate at 37°C for 15 minutes. Alternatively, 2 µl of DNase I-XT (NEB# M0570) can be added directly to the IVT product and incubated at 37°C for 15 minutes.

6. Proceed with [purification of synthesized RNA](#) and/or [evaluation of transcription products](#) yield and/or length.

*Optional:* If a poly(A) tail is desired and is not encoded in the template plasmid, you can add one post-transcriptionally with *E. coli* Poly(A) Polymerase (NEB# M0276).



## Purification of Synthesized RNA

In general, RNA synthesized by *in vitro* transcription can be purified by LiCl precipitation, phenol-chloroform extraction followed by ethanol precipitation, or by using a spin column-based method. If absolute full-length RNA is required, we recommend gel purification. For capped RNA, non-radioactively labeled RNA, or high specific activity radiolabeled RNA probes, spin column chromatography is the preferred method.

### LiCl Precipitation

LiCl precipitation can be used for quick recovery of the synthesized RNA and is an effective method for removing the majority of unincorporated NTPs and enzymes. However, RNAs shorter than 300 bases or at concentrations lower than 0.1 mg/ml do not precipitate well. In such cases, other purification methods may be used.

#### Protocol

1. Adjust the reaction volume to 50  $\mu$ l by adding nuclease-free water.
2. To the 50  $\mu$ l reaction, add 25  $\mu$ l LiCl solution and mix well.
3. Incubate at  $-20^{\circ}\text{C}$  for 30 minutes
4. Centrifuge at  $4^{\circ}\text{C}$  for 15 minutes at top speed to pellet the RNA.
5. Remove the supernatant carefully.
6. Rinse the pellet by adding 500  $\mu$ l of cold 70% ethanol and centrifuge at  $4^{\circ}\text{C}$  for 10 minutes.
7. Remove the ethanol carefully. Spin the tube briefly to bring down any liquid on the wall.
8. Remove residual liquid carefully using a sharp tip (e.g., loading tip).
9. Air dry the pellet and resuspend the mRNA in 50-200  $\mu$ l of 0.1 mM EDTA or a suitable RNA storage solution. Mix well.
10. Store the RNA at  $-20^{\circ}\text{C}$  or below.

### Phenol-chloroform Extraction and Ethanol Precipitation

For removal of proteins and most of the free nucleotides, phenol-chloroform extraction and ethanol precipitation of RNA transcripts is the preferred method.

#### Protocol

1. Adjust the reaction volume to 180  $\mu$ l by adding nuclease-free water. Add 20  $\mu$ l of 3 M sodium acetate (pH 5.2) or 20  $\mu$ l of 5 M ammonium acetate and mix thoroughly.
2. Extract with an equal volume of 1:1 phenol:chloroform mixture, followed by two extractions with chloroform. Collect the aqueous phase and transfer it to a new tube.
3. Precipitate the RNA by adding 2 volumes of ethanol. Incubate at  $-20^{\circ}\text{C}$  for at least 30 minutes and collect the pellet by centrifugation.
4. Remove the supernatant carefully.
5. Rinse the pellet by adding 500  $\mu$ l of cold 70% ethanol and centrifuge at  $4^{\circ}\text{C}$  for 10 minutes.
6. Remove the ethanol carefully. Spin the tube briefly to bring down any liquid on the wall.
7. Remove residual liquid carefully using a sharp tip (e.g., loading tip).
8. Air dry the pellet and resuspend the mRNA in 50-200  $\mu$ l of 0.1 mM EDTA or a suitable RNA storage solution.
9. Store the mRNA at  $-20^{\circ}\text{C}$  or below.

## Spin Column Chromatography

Spin columns will remove unincorporated nucleotides, proteins, and salts. We recommend using a Monarch Spin RNA Cleanup Kit (10 µg capacity NEB #T2030, 50 µg capacity #T2040, or 500 µg capacity #T2050) with the Monarch® Spin RNA Cleanup Kit Protocol.

Adjust the volume of the reaction mixture to 100 µl by adding nuclease-free water to the IVT product and mix well. Purify the RNA by following the manufacturer's instructions. Be aware that some reactions could produce 180 µg of RNA, which may exceed individual column capacity and require additional columns.

## Gel Purification

When high purity RNA transcript is desired (such as making RNA probe for RNase protection assays), we recommend gel purification of the transcription product. Gel purification of mRNA tailed by Poly(A) Polymerase is not recommended because the tailed mRNA is heterogeneous in length and will result in a smear on the gel. If the poly(A) tail is encoded in the DNA template, the RNA should resolve into a single band for gel purification.

The Monarch Spin RNA Cleanup Kits (NEB #T2030, #T2040 or #T2050) can be used for extraction of RNA from gels, although this is not their primary application, so recoveries may range from 40-70% (see protocol included in NEB #T2030, #T2040 or #T2050 product manual).

## Evaluation of Reaction Products

### Quantification by UV Light Absorbance

RNA concentration can be determined by measuring the ultraviolet light absorbance at 260 nm. However, any unincorporated nucleotides and template DNA in the mixture will affect the reading, so must be removed before the RNA concentration can be quantified.

A Nanodrop spectrophotometer can directly read RNA concentrations from 10 ng/µl to 3000 ng/µl; it may be necessary to dilute your RNA prior to measurement. For single-stranded RNA, 1 A<sub>260</sub> is equivalent to an RNA concentration of 40 µg/ml. The RNA concentration can be calculated as follows: A<sub>260</sub> x dilution factor x 40 = \_\_\_ µg/ml RNA

### Analysis of Transcription Products by Gel Electrophoresis

To evaluate transcript length, integrity and quantity, an aliquot of the transcription reaction should be run on an appropriate denaturing agarose or polyacrylamide gel. Transcripts larger than 0.3 kb can be run on agarose gels, whereas denaturing polyacrylamide gels (5–15%) are necessary for smaller transcripts. The gels should be run under denaturing conditions to minimize formation of secondary structures by the transcript.

#### Sample preparation

1. Prepare denatured samples by mixing 100–200 ng RNA sample with 10-20 µl of RNA Loading Dye, 2X (NEB #B0363).
2. Denature the RNA sample and an aliquot of RNA marker by heating at 70°C for 10 minutes.
3. Pulse-spin prior to loading onto the gel.

#### Preparation of denatured samples and gels

##### 4a. Denaturing agarose gel:

Load 100-200ng denatured RNA sample.

It is common practice to electrophorese RNA on a fully denaturing agarose gel, such as one containing formaldehyde. However, in many cases it is possible to run RNA on a native agarose gel and obtain suitable results. In fact, it has been demonstrated that treatment of RNA samples in a denaturing buffer (outlined below) maintains the RNA molecules in a denatured state for at least 3 hours during electrophoresis. The use of native agarose gels eliminates problems associated with toxic chemicals and the difficulties encountered when staining and blotting formaldehyde gels.

##### 4b. Denaturing PAGE/Urea Gel:

Load 50-100ng denatured RNA sample.

We recommend using commercially available premade gels and standard TBE gel running buffer (10X TBE buffer: 0.9 M Tris Base, 0.9 M Boric Acid, 20 mM EDTA).

#### Visualize RNA

5. Stain the gel post-electrophoresis with SYBR® Gold (preferred) or ethidium bromide.

## Troubleshooting

### Control Reaction

The FLuc control template DNA is a linearized plasmid containing the firefly luciferase gene under the transcriptional control of T7 promoter. The size of the runoff transcript is 1.76 kb. The control reaction should yield  $\geq 150$   $\mu$ g RNA transcript in 2 hours.

If the control reaction is not working, there may be technical problems during reaction set up. Repeat the reaction by following the protocol carefully and supplement with DTT to a final concentration of 5 mM; take any precaution to avoid RNase contamination. Contact NEB for technical assistance.

The control plasmid sequence can be found within the [DNA Sequences and Maps Tool](#) under the name “FLuc Control Plasmid”. The FLuc control template is generated by linearizing the plasmid with StuI.

### Low Yield of Full-length RNA

If the transcription reaction with your template generates full-length RNA, but the yield is significantly lower than expected, it is possible that contaminants in the DNA template are inhibiting the RNA polymerase, or the DNA concentration may be incorrect. Alternatively, additional purification of DNA template may be required. Phenol-chloroform extraction is recommended (see template DNA preparation section).

### Addition of DTT

The RNA polymerase in the kit is sensitive to oxidation and could result in lower RNA yield over time due to repeated handling etc. Adding DTT to the reaction may help restore the kit performance in such cases. Adding DTT will not compromise the reaction in any situation.

### Low Yield of Short Transcript

High yields of short transcripts (< 0.3 kb) are achieved by extending incubation time and increasing the amount of template. Incubation of reactions up to 16 hours (overnight) or using up to 2  $\mu$ g of template will help to achieve maximum yield.

### RNA Transcript Smearing on Denaturing Gel

If the RNA appears degraded (e.g. smeared) on a denaturing agarose or polyacrylamide gel, DNA template is contaminated with RNase. DNA templates contaminated with RNase can affect the length and yield of RNA synthesized (a smear below the expected transcript length). If the plasmid DNA template is contaminated with RNase, perform phenol/chloroform extraction, then ethanol precipitate and dissolve the DNA in nuclease-free water (see template DNA preparation section).

### RNA Transcript of Larger Size than Expected

If the RNA transcript appears larger than expected on a denaturing gel, template plasmid DNA may be incompletely digested. Even small amounts of undigested circular DNA can produce large amounts of long transcripts. Check template for complete digestion, if undigested plasmid is confirmed, repeat restriction enzyme digestion.

Larger size bands may also be observed when the RNA transcript is not completely denatured due to the presence of strong secondary structures.

### RNA Transcript of Smaller Size than Expected

If denaturing gel analysis shows the presence of smaller bands than the expected size, it is most likely due to premature termination by the polymerase. Some sequences which resemble T7 RNA Polymerase termination signals will cause premature termination. Incubating the transcription reaction at lower temperatures, for example at 30°C, may increase the proportion of full-length transcript, however the yield will be decreased. For GC rich templates, or templates with secondary structures, incubation at 42°C may improve yield of full-length transcript.

## Ordering Information

NEB #	PRODUCT	SIZE
E2050S/L	HiScribe T7 Quick High Yield RNA Synthesis Kit	50/250 reactions

## COMPANION PRODUCTS

NEB #	PRODUCT	SIZE
T2050S/L	Monarch RNA Cleanup Kit (500 µg)	10/100 preps
T2040S/L	Monarch RNA Cleanup Kit (50 µg)	10/100 preps
T2030S/L	Monarch RNA Cleanup Kit (10 µg)	10/100 preps
B0363S	RNA Loading Dye (2X)	4 x 1 ml
M0303S/L	DNase I (RNase-Free)	1,000/5,000 units
M0570S/L	DNase-XT	1,000/5,000 units
M0493S/L	Q5 Hot Start High-Fidelity DNA Polymerase	100/500 units
M0494S/L	Q5 Hot Start High-Fidelity 2X Master Mix	100/500 units
N0362S	ssRNA Ladder	25 gel lanes
N0364S	Low Range ssRNA Ladder	25 gel lanes
S1411S/L	3'-O-Me-m7G(5')ppp(5')G RNA Cap Structure Analog	1/5 µmol
S1404S/L	m7G(5')ppp(5')G RNA Cap Structure Analog	1/5 µmol
S1405S/L	m7G(5')ppp(5')A RNA Cap Structure Analog	1/5 µmol
S1406S/L	G(5')ppp(5')A RNA Cap Structure Analog	1/5 µmol
S1407S/L	G(5')ppp(5')G RNA Cap Structure Analog	1/5 µmol
M2080S	Vaccinia Capping System	400 units
M2081S/L	Faustovirus Capping Enzyme	500/2,500 units
M0366S	mRNA Cap 2'-O-Methyltransferase	2,000 units
M0276S/L	<i>E. coli</i> Poly(A) Polymerase	100/500 units
N0466S/L	Ribonucleotide Solution Mix	10/50 µmol of each
N0450S/L	Ribonucleotide Solution Set	10/50 µmol of each
N0431S	N1-Methyl-Pseudouridine-5'-Triphosphate	0.1 ml
N0432S	5-Methyl-Cytidine-5'-Triphosphate	0.1 ml
N0433S	Pseudouridine-5'-Triphosphate	0.1 ml
N0434S	5-Methoxy-Uridine-5'-Triphosphate	0.1 ml

## Revision History

REVISION #	DESCRIPTION	DATE
1.0	N/A	2/13
1.1		5/13
2.0		4/14
2.1		1/17
3.0		1/19
4.0	Applied new manual format.	4/20
5.0	Updated to include addition of DTT. Also updated location of control plasmid and LiCl Solution concentration. Also updated table formatting and legal footer.	7/23
6.0	Updated to include new L size.	1/24
6.1	Updated final concentration of NTP Buffer Mix in table on page 5 to "5 mM each NTP final".	9/24
7.0	Reorganized and updated protocols for consistency. Removed "optional" note about DTT.	12/25

This product is intended for research purposes only. This product is not intended to be used for therapeutic or diagnostic purposes in humans or animals.

Products and content are covered by one or more patents, trademarks and/or copyrights owned or controlled by New England Biolabs, Inc (NEB). The use of trademark symbols does not necessarily indicate that the name is trademarked in the country where it is being read; it indicates where the content was originally developed. See [www.neb.com/trademarks](http://www.neb.com/trademarks). The use of these products may require you to obtain additional third-party intellectual property rights for certain applications. For more information, please email [busdev@neb.com](mailto:busdev@neb.com).

B CORPORATION® is a registered trademark of B Lab IP, LLC, Inc.

NANODROP® is a registered trademark of NanoDrop Technologies, Inc.

SYBR® is a registered trademark of Life Technologies, Inc.

© Copyright 2025, New England Biolabs, Inc.; all rights reserved

