Reverse Transcription System

INSTRUCTIONS FOR USE OF PRODUCT A3500.

Reverse Transcription Protocol

Reverse Transcription Reaction (First-Strand cDNA Synthesis)

- Place 1µg (2µl) of 1.2kb Kanamycin Positive Control RNA, poly(A)+ mRNA or total RNA in a microcentrifuge tube, and incubate at 70°C for 10 minutes. Centrifuge briefly in a microcentrifuge, then place on ice.
- Prepare a 20µl reaction by adding the following reagents in the order listed (this reaction can be scaled up or down, depending on the amount of RNA):

Component	Amount
MgCl ₂ , 25mM	4µl
Reverse Transcription 10X Buffer	2µl
dNTP Mixture, 10mM	2µl
Recombinant RNasin [®] Ribonuclease Inhibitor	0.5µl
AMV Reverse Transcriptase (High Conc.)	15u
Oligo(dT) ₁₅ Primer OR Random Primers	0.5µg
1.2kb Kanamycin Positive Control RNA (2µI)	
OR poly(A)+ mRNA OR total RNA	<u> 1µg</u>
Nuclease-Free Water to a final volume of	20µl

- When using Oligo(dT)₁₅ Primer, incubate the reaction at 42°C for 15 minutes. When using Random Primers, incubate the reaction at room temperature for 10 minutes, then incubate at 42°C for 15 minutes.
- 4. Heat the sample at 95°C for 5 minutes, then incubate at 0–5°C for 5 minutes.

Dilution of the Reaction for Amplification

- 1. Dilute the first-strand cDNA synthesis reaction to 100µl with TE buffer or Nuclease-Free Water.
- Prepare a 100µl PCR amplification mix by combining the following reagents. Template-specific upstream and downstream primers should be used for this reaction.

Component	Amount
first-strand cDNA reaction	10–20µl
dNTP Mixture, 10mM	1.8µI
MgCl ₂ , 25mM	7.5µl
Reverse Transcription 10X Buffer	9.8µI
upstream primer	50pmol
downstream primer	50pmol
<i>Taq</i> DNA polymerase	<u>2.5 units</u>
Nuclease-Free Water to a final volume of	100µl

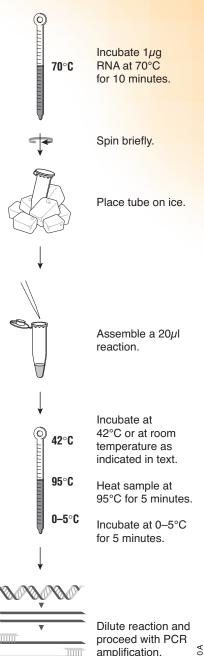
3. Proceed to thermal cycling according to your own specific experiment.

See additional protocol information in Technical Bulletin #TB099, available online at: www.promega.com

ORDERING/TECHNICAL INFORMATION:

www.promega.com • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601







reaction and d with PCR cation.



