GoScript[™] Reverse Transcription System

INSTRUCTIONS FOR USE OF PRODUCTS A5000 AND A5001.

First-Strand cDNA Synthesis

The following procedure can be used to convert up to 5µg of total RNA or up to 500ng of poly(A) RNA into first-strand cDNA.

1. Mix and briefly centrifuge each component before use. Combine the following:

Component Experimental RNA (up to 5ug/reaction)	Volume Xul
Primer [Oligo(dT) ₁₅ (0.5µg/reaction) and/or Random Primer (0.5µg/reaction) or	
gene-specific primer (10-20pmol/reaction)]	XμI
Nuclease-Free Water	XμI
Final volume	5µl

- 2. Heat in a 70°C heat block for 5 minutes. Immediately chill in ice water for at least 5 minutes. Centrifuge 10 seconds in a microcentrifuge. Store on ice until reverse transcription mix is added.
- 3. Prepare the reverse transcription reaction mix, 15µl for each cDNA reaction. Combine on ice, in the order listed.

Component	Volume
GoScript [™] 5X Reaction Buffer	4.0µI
MgCl ₂ (final concentration 1.5–5.0mM) ¹	1.2–6.4µI
PCR Nucleotide Mix (final concentration 0.5mM each dNTP) ²	1.0µI
Recombinant RNasin [®] Ribonuclease Inhibitor (optional)	20units
GoScript™ Reverse Transcriptase	1.0µI
Nuclease-Free Water (to a final volume of 15µl)	XμI
Final volume	15µl

¹Mg²⁺ concentration should be optimized to 1.5–5.0mM (MgCl₂ provided at 25mM). ²If isotopic or nonisotopic incorporation is desired for monitoring first-strand cDNA synthesis. $\alpha^{[32P]}$ -dCTP or other modified nucleotides may be supplemented into the PCR Nucleotide Mix. See Section 4.D, TM316, for analysis suggestions.

- 4. Combine 15µl of reverse transcription mix with 5µl of RNA and primer mix.
- 5. **Anneal** in a heat block at 25°C for 5 minutes.
- 6. Extend in a heat block at 42°C for up to one hour. Reactions can be stopped at this point for analysis of the cDNA or may be frozen for long-term storage.
- 7. Inactivate Reverse Transcriptase: Before proceeding with gPCR, inactivate the reverse transcriptase in a heat block at 70°C for 15 minutes.

(continued)

ORDERING/TECHNICAL INFORMATION:

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



2977MB

Place tube on ice. Centrifuge briefly. Assemble a 20µl reaction (Step 4). Incubate at \bigcirc 25°C (anneal) 25°C then at 42°C (extend). 42°C Store at 0-5°C. 0-5°C qPCR: Inactivate 70°C reverse transcriptase at 70°C. Proceed with aPCR (next page).



Incubate RNA and

primers at 70°C for

5 minutes.

 \bigcirc

70°C

GoScript[™] Reverse Transcription System

INSTRUCTIONS FOR USE OF PRODUCTS A5000 AND A5001.



cDNA Quantification Using qPCR

- 1. Quantify specific targets in samples of undiluted or diluted cDNA using GoTaq[®] qPCR Master Mix.
- 2. Alternatively, add a diluted sample of cDNA, as determined to be optimal (up to 20% of the reaction volume or 100ng of input RNA).

See additional protocol information in the *GoScript*[™] *Reverse Transcription System Technical Manual,* #TM316, available at: www.promega.com/tbs



