

# Access RT-PCR System

INSTRUCTIONS FOR USE OF PRODUCTS A1250, A1260 AND A1280.

# Quick PROTOCOL

## Assemble Reactions

1. Combine the reagents below in a thin-walled 0.5ml reaction tube on ice.

Reagents	Volume	Final Concentration
Nuclease-Free Water (to a final volume of 50 $\mu$ l)	X $\mu$ l	
AMV/T7/5X Reaction Buffer	10 $\mu$ l	1X
dNTP Mix (10mM each dNTP)	1 $\mu$ l	0.2mM
Downstream primer	50pmol	1 $\mu$ M
Upstream primer	50pmol	1 $\mu$ M
25mM MgSO <sub>4</sub>	2 $\mu$ l	1mM

2. Mix by pipetting. Add the remaining components.

AMV Reverse Transcriptase (5u/ $\mu$ l)	1 $\mu$ l	0.1u/ $\mu$ l
T7/ DNA Polymerase (5u/ $\mu$ l)	1 $\mu$ l	0.1u/ $\mu$ l

3. Gently vortex. Initiate the reaction by adding:

RNA template	<u>Y</u> $\mu$ l	10 <sup>3</sup> –10 <sup>6</sup> copies
	50 $\mu$ l	

4. Overlay the reactions with 1 or 2 drops of mineral oil.

## First Strand cDNA Synthesis

These PCR cycling profiles are only a guideline. Cycling conditions should be optimized for each RNA template.

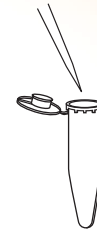
1 cycle	45°C for 45 minutes	reverse transcription
1 cycle	94°C for 2 minutes	AMV RT inactivation and RNA/cDNA/primer denaturation

## Second Strand Synthesis and PCR Amplification

40 cycles	94°C for 30 seconds	denaturation
	60°C for 1 minute	annealing
	68°C for 2 minutes	extension
1 cycle (optional)	68°C for 7 minutes	final extension
1 cycle	4°C	soak

Analyze 2.5 $\mu$ l of the reaction products by agarose gel electrophoresis. Store the remainder of the reaction at –20°C.

Additional protocol information is available in Technical Bulletin #TB220, available online at: [www.promega.com](http://www.promega.com)



Assemble reactions.



First strand cDNA synthesis.



Second strand cDNA synthesis and PCR amplification.



Analyze products by agarose gel electrophoresis.

2946MA05\_0A

### ORDERING/TECHNICAL INFORMATION:

[www.promega.com](http://www.promega.com) • Phone 608-274-4330 or 800-356-9526 • Fax 608-277-2601



**Promega**

Printed in USA, Revised 3/09  
Part# 9FB034

© 2000–2009 Promega Corporation. All Rights Reserved.