

## I. List of Components

Store all components at  $-20^{\circ}\text{C}$ .

- SMART MMLV Reverse Transcriptase
- 5X First-Strand Buffer
- 100 mM DTT

## II. Additional Materials Required

- RNase-free  $\text{H}_2\text{O}$
- dNTP Mix (10 mM each); we recommend Advantage® UltraPure PCR Deoxynucleotide Mix (Cat. No. 639125)
- Oligo (dT)<sub>12-18</sub>, random hexamer or gene-specific primers may be used.
- [Optional] 60 mM EDTA

## III. For Routine First-Strand cDNA Synthesis Reactions

This 20- $\mu\text{l}$  reaction is suitable for synthesizing first-strand cDNA from 5 ng–5  $\mu\text{g}$  of total RNA or 10 ng–1  $\mu\text{g}$  of poly A<sup>+</sup> RNA.

1. Add 2.5  $\mu\text{l}$  of 20  $\mu\text{M}$  primer stock (final concentration 2.5  $\mu\text{M}$ ) to your RNA sample. Add RNase-free  $\text{H}_2\text{O}$  to a final volume of 11.5  $\mu\text{l}$ .
2. Heat the mixture at  $70^{\circ}\text{C}$  for 3 min, and immediately cool on ice.
3. Centrifuge briefly, then add the following:
  - 4  $\mu\text{l}$  5X First-Strand Buffer
  - 2  $\mu\text{l}$  dNTP Mix
  - 2  $\mu\text{l}$  100 mM DTT
4. Mix the contents of the tube by gently pipetting up and down. Add 0.5  $\mu\text{l}$  SMART MMLV RT and mix again.
5. Incubate at  $42^{\circ}\text{C}$  for 60 min.

**NOTE:** Samples can be incubated for 50–90 min if necessary.

6. Terminate the reaction by heating at  $70^{\circ}\text{C}$  for 15 min, or by adding 4  $\mu\text{l}$  of 60 mM EDTA.

## IV. References

Ausubel, F. M., *et al.* (1995) In *Current Protocols in Molecular Biology* (Greene Publishing Associates, Inc. and John Wiley & Sons, Inc.), Supplement 29, Section 5.5.

Contact Us For Assistance	
Customer Service/Ordering	Technical Support
Telephone: 800.662.2566 (toll-free)	Telephone: 800.662.2566 (toll-free)
Fax: 800.424.1350 (toll-free)	Fax: 650.424.1064
Web: <a href="http://www.clontech.com">www.clontech.com</a>	Web: <a href="http://www.clontech.com">www.clontech.com</a>
E-mail: <a href="mailto:orders@clontech.com">orders@clontech.com</a>	E-mail: <a href="mailto:tech@clontech.com">tech@clontech.com</a>

## Notice to Purchaser

Our products are to be used for research purposes only. They may not be used for any other purpose, including, but not limited to, use in drugs, *in vitro* diagnostic purposes, therapeutics, or in humans. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without prior written approval of Clontech Laboratories, Inc.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product's web page at <http://www.clontech.com>. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Clontech®, the Clontech logo, Advantage, and SMART are trademarks of Clontech Laboratories, Inc. All other marks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. Clontech is a Takara Bio Company. ©2015 Clontech Laboratories, Inc.

This document has been reviewed and approved by the Clontech Quality Assurance Department.