

Clontech Laboratories, Inc.

# SMARTer® Ultra® Low RNA Kit for Illumina® Sequencing User Manual

Cat. No. 634936  
PT5163-1  
(010616)

---

**Clontech Laboratories, Inc.**

A Takara Bio Company  
1290 Terra Bella Avenue, Mountain View, CA 94043, USA  
U.S. Technical Support: [tech@clontech.com](mailto:tech@clontech.com)

United States/Canada 800.662.2566	Asia Pacific +1.650.919.7300	Europe +33.(0)1.3904.6880	Japan +81.(0)77.543.6116
---	---------------------------------	------------------------------	-----------------------------

Page 1 of 21

Table of Contents

I. List of Components ..... 4

II. Additional Materials Required ..... 5

III. Introduction ..... 6

IV. SMARTer cDNA Synthesis ..... 9

    A. Requirements for Preventing Contamination ..... 9

    B. General Requirements ..... 9

    C. Sample Recommendations ..... 10

    D. Sample Requirements ..... 10

    E. Protocol: First-Strand cDNA Synthesis ..... 11

    F. Protocol: Purification of First-Strand cDNA using SPRI Ampure Beads ..... 13

    G. Protocol: ds cDNA Amplification by LD PCR ..... 14

V. Amplified cDNA Purification & Validation ..... 15

    A. Protocol: Purification of ds cDNA using SPRI Ampure Beads ..... 15

    B. Validation Using the Agilent 2100 BioAnalyzer ..... 16

VI. Covaris Shearing ..... 17

    A. Protocol: Covaris Shearing of Full-length cDNA ..... 17

VII. Appendix A: PCR Clean Work Station Guidelines ..... 17

    A. Equipment Needed in the PCR Clean Work Station ..... 17

    B. PCR Clean Work Station Operation Instructions ..... 18

VIII. Appendix B: Working with Cells ..... 19

    A. Protocol: Screen for Cell Culture Media ..... 19

    B. Cell Preparation for First-Strand cDNA Synthesis ..... 19

IX. Appendix C: PCR Optimization ..... 19

X. Appendix D: Troubleshooting Guide ..... 20

XI. References ..... 21

**Table of Figures**

Figure 1. Protocol Overview ..... 6  
Figure 2. Flowchart of SMARTer cDNA synthesis..... 7  
Figure 3. SMART Adapter in Primer 2 Read ..... 8  
Figure 4. Optional setup to ensure proper and steady positioning of tubes containing first-strand cDNA..... 13  
Figure 5. Electropherogram example results from Agilent 2100 Bioanalyzer..... 16

**Table of Tables**

Table 1. Sample Preparation Guidelines ..... 11  
Table 2. Cycling Guidelines Based on Amount of Starting Material ..... 14  
Table 3. Process Configuration Panel Set Up ..... 17  
Table 4. Validated Media..... 19  
Table 5. Troubleshooting Guide for SMARTer cDNA Synthesis & Amplification..... 20

## I. List of Components

The following components have been specifically designed to work together and are optimized for this particular protocol. Please do not make any substitutions. The substitution of reagents in the kit and/or a modification of the protocol may lead to unexpected results.

The SMARTer Ultra Low RNA Kit for Illumina Sequencing (Cat. No. 634936) consists of:

- **The Advantage® 2 PCR Kit (Cat. No. 639206)**, which has been specially formulated to provide automatic hot-start PCR (Kellogg *et al.*, 1994)—and can efficiently amplify full-length cDNAs with a significantly lower error rate than that of conventional PCR (Barnes, 1994).
- **The SMARTer Ultra Low RNA Kit for Illumina Sequencing Components (Cat. No. 634948)**, which includes:

### Box 1:

- 10 x 20 µl SMARTer II A Oligonucleotide (12 µM)
- 5 µl Control Total RNA (1 µg/µl)

### Box 2:

- 200 µl 3' SMART CDS Primer II A (12 µM)
- 250 µl IS PCR Primer (12 µM)
- 400 µl 5X First-Strand Buffer (RNase-Free)
  - 250 mM Tris-HCl (pH 8.3)
  - 375 mM KCl
  - 30 mM MgCl<sub>2</sub>
- 400 µl dNTP Mix (dATP, dCTP, dGTP, and dTTP, each at 10 mM)
- 100 µl Dithiothreitol (DTT; 100 mM)
- 200 µl SMARTScribe™ Reverse Transcriptase (100 U/µl)
- 10 ml Nuclease-Free Water
- 550 µl RNase Inhibitor (40 U/µl)
- 10 ml Dilution Buffer
- 10 ml Purification Buffer (10 mM Tris-Cl, pH 8.5)

## Storage Conditions

- Store Control Total RNA and SMARTer IIA Oligonucleotide at  $-70^{\circ}\text{C}$ .
- Store Dilution Buffer at  $-20^{\circ}\text{C}$ . Once thawed, the buffer can be stored at  $4^{\circ}\text{C}$ .
- Store Purification Buffer at  $-20^{\circ}\text{C}$ . Once thawed, the buffer can be stored at Room Temperature.
- Store all other reagents at  $-20^{\circ}\text{C}$ .

## II. Additional Materials Required

The following reagents are required but not supplied. These materials have been validated to work with this protocol. Please do not make any substitutions because you may not obtain the expected results:

- Single channel pipette: 10 µl, 20 µl and 200 µl, one each
- Eight channel pipette: 20 µl and 200 µl, one each
- Filter pipette tips: 10 µl, 20 µl and 200 µl, one box each
- One QuickSpin minicentrifuge for 1.5-ml tubes
- One QuickSpin minicentrifuge for 0.2-ml tubes
- One marker pen
- Small 1.5-ml tube rack
- One bag of 8-strip nuclease-free 0.2-ml thin wall PCR tubes with caps
- DNA-OFF™ Solution (Takara® Cat. No. 9036)

### **For PCR Amplification & Validation:**

- One dedicated PCR thermal cycler used only for first-strand synthesis
- High Sensitivity DNA Kit (Agilent Cat No. 5067-4626)
- IsoFreeze Flipper Rack (MIDSCI Cat. No. TFL-20)
- IsoFreeze PCR Rack (MIDSCI Cat. No. 5640-T4)
- Nuclease-free thin-wall PCR tubes (0.2 ml; USA Scientific Cat. No.1402-4700)
- Nuclease-free nonsticky 1.5-ml tubes (USA Scientific Cat. No. 1415-2600)

### **For SPRI Bead Purification:**

- Agencourt AMPure PCR Purification Kit  
(5 ml Beckman Coulter Part No. A63880; 60 ml Beckman Coulter Part No. A63881)  
Use this kit for the SPRI Purifications (Sections IV.F & V.A).
- MagnaBot II Magnetic Separation Device (Promega Part No. V8351)  
Use this stand for the first purification (Section IV.F).
- Magnetic Stand-96 (Ambion Part No. AM10027)  
Use this stand for the second purification (Section V.A).
- 96-well V-bottom Plate (500 µl; VWR Cat. No. 47743-996)
- MicroAmp Clean Adhesive Seal (AB Part No. 4306311)
- 80% Ethanol

### **For Sequencing Library Generation:**

- Low Input Library Prep Kit (Cat. No. 634947)
- Covaris Instrument and Related Materials for DNA Shearing

### III. Introduction

#### SMARTer cDNA Synthesis for the Illumina Sequencing Platform

The SMARTer Ultra Low RNA Kit allows high-quality cDNA synthesis starting from as little as 10 pg of total RNA or cells. The kit has been designed and validated to prepare cDNA samples for sequencing and quantitation with the Illumina HiSeq® and Genome Analyzer™ sequencing instruments. The entire library construction protocol can be completed within two days (see Figure 1 for an overview of the protocol). SMART® technology offers unparalleled sensitivity and unbiased amplification of cDNA transcripts, enabling a direct start from your sample. Most importantly, SMART technology enriches for full-length transcripts and maintains the true representation of the original mRNA transcripts; these factors are critical for transcriptome sequencing and gene expression analysis.

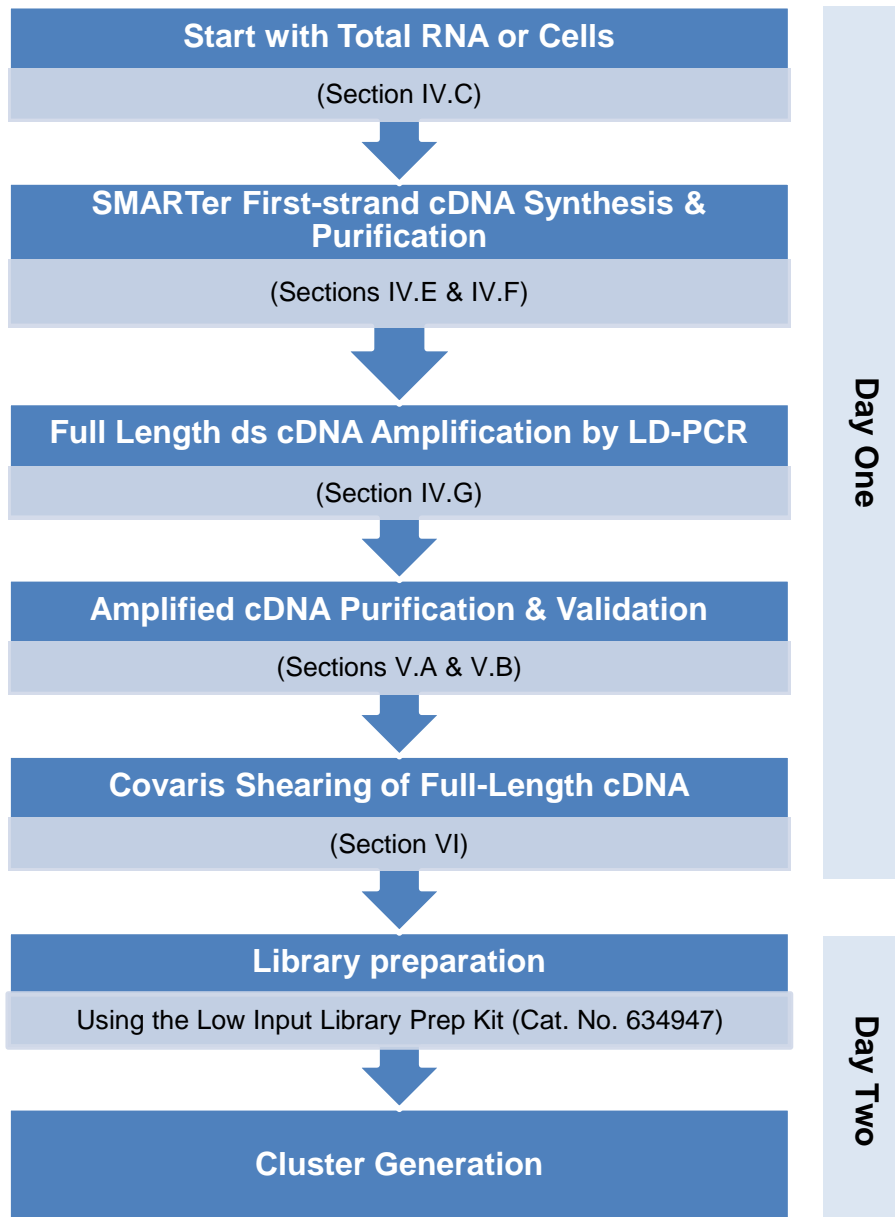
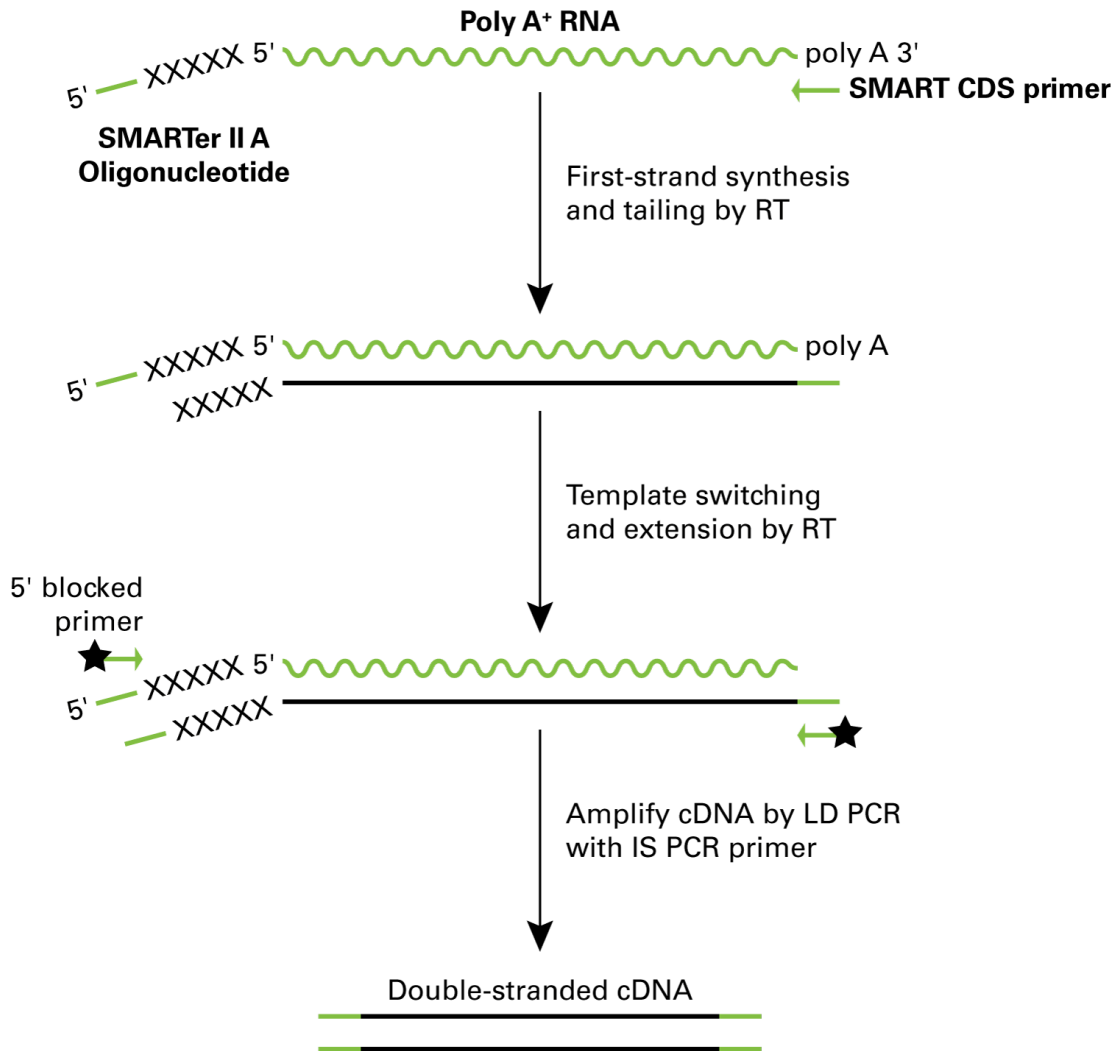


Figure 1. Protocol Overview.

SMARTer cDNA synthesis starts with picogram amounts of total RNA. A modified oligo(dT) primer (the SMART CDS Primer) primes the first-strand synthesis reaction (Figure 2). When SMARTScribe Reverse Transcriptase reaches the 5' end of the mRNA, the enzyme's terminal transferase activity adds a few additional nucleotides to the 3' end of the cDNA. The carefully-designed SMARTer Oligonucleotide base-pairs with the non-template nucleotide stretch, creating an extended template to enable SMARTScribe RT continue replicating to the end of the oligonucleotide (Chenchik *et al.*, 1998). The resulting full-length, single-stranded (ss) cDNA contains the complete 5' end of the mRNA, as well as sequences that are complementary to the SMARTer Oligonucleotide. In cases where the RT pauses before the end of the template, the addition of non-template nucleotides is much less efficient than with full-length cDNA/RNA hybrids, thus the overhang needed for base-pairing with the SMARTer Oligonucleotide is absent. The SMARTer anchor sequence and the poly A sequence serve as universal priming sites for end-to-end cDNA amplification. In contrast, cDNA without these sequences, such as prematurely terminated cDNAs, contaminating genomic DNA, or cDNA transcribed from poly A-RNA, will not be exponentially amplified. However, truncated RNAs with poly A tails that are present in poor quality RNA starting material will be amplified yielding shorter cDNA fragments.



**Figure 2. Flowchart of SMARTer cDNA synthesis.** The SMARTer II A Oligonucleotide, 3' SMART CDS Primer II A, and IS PCR Primer all contain a stretch of identical sequence (see Section I for sequence information).

## SMART Adapter in Primer 2 Read

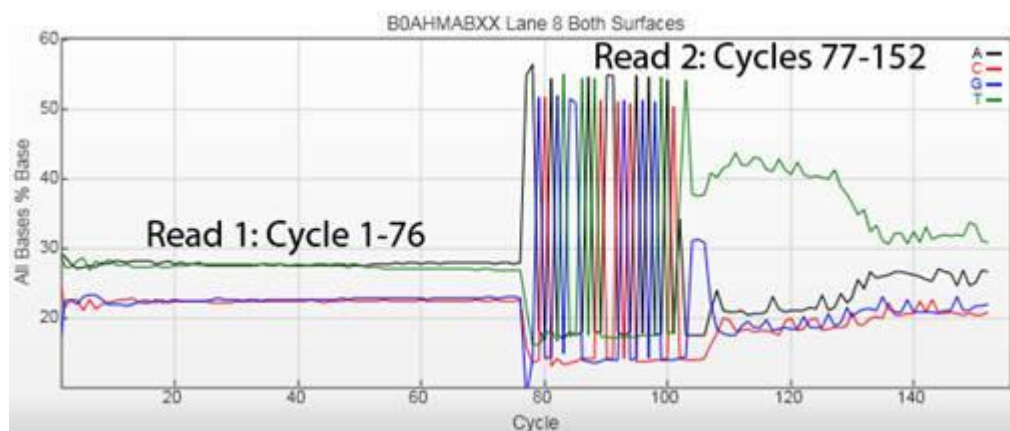
The IS primer used for amplification of the double-stranded cDNA is blocked, which stops Illumina sequencing adapter ligation at the 5' ends of the ds cDNA fragments containing the SMART sequence. In many library preparation methods, the double-stranded Illumina adapters are added to the cDNA fragments through ligation. Unfortunately, in these reactions, ligation takes place between the bottom strand of the cDNA fragment and the Illumina adapter containing Read Primer 2, at a low and somewhat variable rate. If ligation is also successful on the other, unblocked side of the same cDNA fragment, this bottom strand can be amplified by the subsequent PCR and can ultimately form clusters for sequencing on Illumina machines.

Upon sequencing these clusters, the SMART adapter will be present in the first 30 cycles in Read 2 of an Illumina sequencing run. In addition, the dT30 sequence from the CDS primer will also be present after the adapter in a subset of these clusters. The presence of the SMART adapter in Read 2 commonly occurs at a high enough rate to be observed in the base distribution by cycle graph generated by the run analysis (Figure 3, cycles 77-106), as does the dT30 sequence (Figure 3, cycles 107-136).

If you are interested in avoiding sequencing the SMART adapter, there are three options:

1. Sequence only from Read Primer 1.
2. Use the Nextera® or Nextera XT library preparation method.
3. Use the Low Input Library Prep Kit (Cat. No. 634947). This unique adapter addition method does not allow for erroneous ligation.

If you have already sequenced with Read Primer 2, the SMART adapter sequence can be trimmed from reads prior to mapping to your transcriptome.



**Figure 3. SMART Adapter in Primer 2 Read.** The presence of the SMART adapter in Read 2 commonly occurs at a high enough rate to be observed in the base distribution by cycle graph generated by the run analysis (cycles 77-106), as does the dT30 sequence (cycles 107-136).



## IV. SMARTer cDNA Synthesis

**NOTE:** Please read the entire protocol before starting. This protocol is optimized for the generation of cDNA starting from ultra-low amounts of total RNA using Clontech's SMART technology. The protocol also works starting from cells. Due to the sensitivity of the protocol, the input material (total RNA or cells) needs to be collected and purified under clean room conditions to avoid contamination. The whole process of SMARTer cDNA Synthesis should be carried out in a PCR Clean Work Station under clean room conditions (see Appendix A: PCR Clean Work Station Guidelines).

### A. Requirements for Preventing Contamination

- **Before you set up the experiment, make sure you have two physically separated work stations. The first should be the PCR Clean Work Station (see Appendix A) for all Pre-PCR experiments that require clean room conditions such as first-strand cDNA synthesis (Section IV.E.) and purification of first-strand cDNA (Section IV.F). The second work station can be in the general laboratory where you will perform PCR, measure cDNA concentration, and work on any other experiments involving the PCR-amplified cDNA.**
- **It is absolutely required to have a PCR work station in a clean room with positive air flow, as contamination may occur very easily. Once contamination occurs it can be difficult to remove.**
- You need to strictly follow the clean room operation rules, i.e., you can only move materials/ supplies from the clean room to the general lab, NOT the other way around. Don't share any equipment/reagents between the clean room and the general lab. You need a separate PCR machine inside the PCR workstation for cDNA synthesis.
- Wear gloves and sleeve covers throughout the procedure to protect your RNA samples from degradation by contaminants and nucleases. Be sure to change gloves and sleeve covers between each section of the protocol.

### B. General Requirements

- **The success of your experiment depends on the quality of your starting sample of RNA. Prior to cDNA synthesis, please make sure that your RNA is intact and free of contaminants.**
- The assay is very sensitive to variations in pipette volume, etc. Please make sure all your pipettes are calibrated for reliable delivery, and make sure nothing is attached to the outside of the tips.
- All lab supplies related to SMARTer cDNA synthesis need to be stored in a DNA-free, closed cabinet. Reagents for SMARTer cDNA synthesis need to be stored in a freezer/refrigerator that has not previously been used to store PCR amplicons.
- Add enzymes to reaction mixtures last, and thoroughly incorporate them by gently pipetting the reaction mixture up and down.
- Do not increase (or decrease) the amount of enzyme added or the concentration of DNA in the reactions. The amounts and concentrations have been carefully optimized for the SMARTer amplification reagents and protocol.
- If you are using this protocol for the first time we strongly recommend that you perform negative and positive control reactions to verify that kit components are working properly.

### C. Sample Recommendations

The sequence complexity and the average length of SMARTer cDNA is noticeably dependent on the quality of starting RNA material. Due to the limiting sample size, most traditional RNA isolation methods may not be applicable. There are several commercially available products that enable purification of total RNA preparations from extremely small samples [e.g. Clontech offers the NucleoSpin RNA XS Kit (Cat. No. 740902.10) for purification of RNA from  $10^2$  cells]. When choosing a purification method (kit) ensure that it is appropriate for your sample amount.

Though SMART Technology is sensitive enough to generate cDNA from as little as 10 pg of total RNA, the use of a higher amount of starting material (100 pg to 10 ng) is recommended for reproducible amplification of low abundance mRNA transcripts.

After RNA extraction, if your sample size is not limiting, we recommend evaluating total RNA quality using the Agilent RNA 6000 Pico Kit (Cat. No. 5067-1513).

### D. Sample Requirements

#### Total RNA

This protocol has been optimized for cDNA synthesis starting from 10 pg of total RNA. However, if your RNA sample is not limiting, we recommend that you start with more total RNA (up to 10 ng). Purified total RNA should be in nuclease-free water. If the total RNA volume is larger than 1  $\mu$ l, perform ethanol precipitation and resuspend the pellet in 1  $\mu$ l of nuclease-free water.

#### Cells

Although this protocol was optimized for cDNA synthesis starting from total RNA, the protocol has also been validated to work starting from cells. See **Appendix B: Working with Cells** for instructions on selecting the appropriate media and preparing your cells for first-strand cDNA synthesis.

**E. Protocol: First-Strand cDNA Synthesis  
(Perform in PCR Clean Work Station)**

**IMPORTANT:** To avoid introducing contaminants to your RNA sample, the first part of the cDNA synthesis protocol (Sections E–G) requires use of a PCR work station in a clean room. Standard clean room procedure should be followed. If no clean room is available, you may work with just a PCR Clean Work Station on a temporary basis. We strongly recommend putting the PCR work station in a clean room to avoid contamination. It is critical to have an air blower in the PCR work station turned “on” during the whole process. Refer to Appendix A for PCR Clean Work Station Guidelines.

1. Prepare a stock solution of Reaction Buffer by mixing the Dilution Buffer with the RNase Inhibitor as indicated below, scale-up as needed:

19 µl	Dilution Buffer
1 µl	RNase Inhibitor
<hr/>	
20 µl	Total Volume

2. See Table 1 for guidelines on setting up your control & test samples. Transfer each whole volume of 3.5 µl to individual 0.2 ml RNase-free PCR tubes in an 8-well strip.

**Table 1. Sample Preparation Guidelines**

Components	Negative Control	Positive Control	Test Sample
Reaction Buffer	2.5 µl	2.5 µl	2.5 µl
Nuclease-free water	1 µl	–	–
Diluted Control RNA*	–	1 µl	–
Sample	–	–	1 µl
<b>Total Volume</b>	<b>3.5 µl</b>	<b>3.5 µl</b>	<b>3.5 µl</b>

\*The Control RNA is supplied at a concentration of 1 µg/µl. The Control RNA should be diluted in nuclease-free water to match the concentration of your test sample. Perform serial dilutions on the Control RNA until you obtain the appropriate concentration.

3. Place the samples on a –20°C prechilled IsoFreeze PCR rack in a PCR clean station, and add 1 µl of 3’ SMART CDS Primer II A (12 µM). Mix the contents and spin the tube(s) briefly in a microcentrifuge:

3.5 µl	Cell/Total RNA in Reaction Buffer (from Table 1)
1 µl	3’ SMART CDS Primer II A (12 µM)
<hr/>	
4.5 µl	Total Volume

4. Incubate the tube(s) at 72°C in a hot-lid thermal cycler for 3 min, and then put the samples on the IsoFreeze PCR rack.

**NOTE:** The initial reaction steps (Step 5-7) are critical for first-strand synthesis and should not be delayed after Step 4. You can prepare your master mix (for Step 5) while your tubes are incubating (Step 4) in order to jump start the cDNA synthesis.

5. Meanwhile, prepare a Master Mix for all reactions plus one by combining the following reagents in the order shown at room temperature:

2 $\mu$ l	5X First-Strand Buffer
0.25 $\mu$ l	DTT (100 mM)
1 $\mu$ l	dNTP Mix (10 mM)
1 $\mu$ l	SMARTer IIA Oligonucleotide (12 $\mu$ M)
0.25 $\mu$ l	RNase Inhibitor
1 $\mu$ l	SMARTScribe Reverse Transcriptase (100 U)*
<hr/>	
5.5 $\mu$ l	Total Volume added per reaction

\* Add the reverse transcriptase to the master mix just prior to use. Mix well by gently vortexing and spin the tube(s) briefly in a microcentrifuge.

6. Add 5.5  $\mu$ l of the Master Mix to each reaction tube from Step 4. Mix the contents of the tube(s) by gently pipetting, and spin the tube(s) briefly to collect the contents at the bottom.
7. Incubate the tubes at 42°C for 90 min.
8. Terminate the reaction by heating the tube(s) at 70°C for 10 min.

## F. Protocol: Purification of First-Strand cDNA using SPRI Ampure Beads (Perform in PCR Clean Work Station)

The first-strand cDNA selectively binds to SPRI beads leaving contaminants in solution which is removed by a magnetic separation. The beads are then directly used for PCR amplification.

### NOTES:

- Before use, beads should be brought to room temperature and mixed well to disperse.
- In order to ensure proper and steady positioning of the tubes containing first-strand cDNA (from Protocol E), you may place the tubes in the top part of an inverted P20 or P200 Rainin Tip Holder which is taped to the MagnaBot II Magnetic Separator.

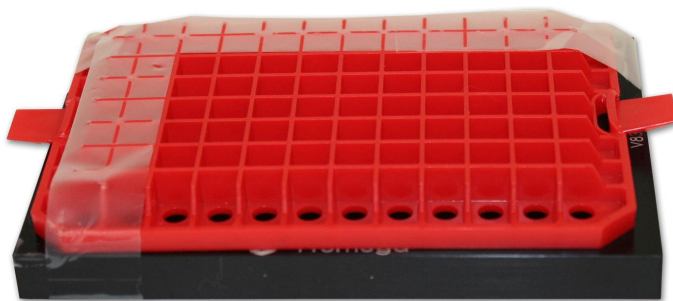


Figure 4. Optional setup to ensure proper and steady positioning of tubes containing first-strand cDNA.

To purify the SMART cDNA from unincorporated nucleotides and small (<0.1 kb) cDNA fragments, follow this procedure for each reaction tube:

1. Add 25  $\mu$ l of SPRI Ampure XP beads to each sample using a 200  $\mu$ l pipetter. Adjust the pipetter to 35  $\mu$ l, and pipette the entire volume up and down 10 times to mix thoroughly. The beads are viscous; suck the entire volume up, and push it out slowly. Incubate at room temperature for 8 minutes to let DNA bind to the beads.
2. Briefly spin the sample tubes to collect the liquid from the side of the wall. Place the sample tubes on the Promega MagnaBot II Magnetic Separation Device for 5 min or longer, until the solution is completely clear.
3. While samples are still on the Magnetic Separation Device pipette out the solution and discard. Briefly spin the tubes to collect the liquid at the bottom.
4. Place the tubes back in the Promega MagnaBot II Magnetic Separation Device for 2 min or longer to let beads separate from the liquid completely. Pipette out the residual liquid from the beads using a 10  $\mu$ l pipetter and discard. Make sure that there is no supernatant remaining in the tube. **Be careful not to take out any beads with the supernatant.**

**G. Protocol: ds cDNA Amplification by LD PCR  
(Perform steps 1 & 2 in PCR Clean Work Station)**

We strongly recommend use of the **Advantage 2 PCR Kit** (included in this kit; also sold separately as Cat. Nos. 639206 & 639207) for PCR amplification. The Advantage 2 Polymerase Mix has been specially formulated for efficient and accurate amplification of cDNA templates by long-distance PCR (LD PCR; Barnes, 1994). Table 2 provides guidelines for optimizing your PCR, depending on the amount of total RNA used in the first-strand synthesis.

**IMPORTANT:** Optimal parameters may vary with different templates and thermal cyclers. To determine the optimal number of cycles for your sample and conditions, we strongly recommend that you perform a range of cycles. Do not exceed the recommended cycle numbers in the table below for different starting amounts of material. Refer to Appendix C for PCR optimization guidelines.

**Table 2. Cycling Guidelines Based on Amount of Starting Material**

Input Amount, Total RNA	Input Amount, Cells	Typical No. of PCR Cycles
10 ng	1,000 cells	12
1 ng	100 cells	12
500 pg	50 cells	13
100 pg	10 cells	15
10 pg	1 cell	18

1. Prepare a PCR Master Mix for all reactions, plus one additional reaction. Combine the following reagents in the order shown, then mix well by vortexing and spin the tube briefly in a microcentrifuge:

5 µl	10X Advantage 2 PCR Buffer
2 µl	dNTP Mix (10 mM)
2 µl	IS PCR Primer (12 µM)
2 µl	50X Advantage 2 Polymerase Mix
39 µl	Nuclease-Free Water
<hr/>	
50 µl	Total Volume per reaction

2. Add 50 µl of PCR Master Mix to each tube containing DNA bound to the beads from Section IV.F., Step 4. Mix well and briefly spin down.

**Important:** Transfer the samples from the PCR Clean Work Station to the general lab. All downstream processes will be performed in the general lab.

3. Place the tube in a preheated thermal cycler with a heated lid. Commence thermal cycling using the following program:

95°C	1 min	
X <sup>a</sup> cycles:		
95°C	15 sec	]
65°C	30 sec	
68°C	6 min	
72°C	10 min	
4°C	forever	

<sup>a</sup>Consult Table 2 for guidelines. Determine the optimal number of PCR cycles using the protocol in **Appendix C: PCR Optimization**.

## V. Amplified cDNA Purification & Validation

### A. Protocol: Purification of ds cDNA using SPRI Ampure Beads

PCR-amplified cDNA is purified by immobilizing it onto SPRI beads. The beads are then washed with 80% Ethanol and eluted in Purification Buffer.

1. Take out a 96-well Axygen V-bottom plate and cover all the wells with a MicroAmp Clean Adhesive Seal. Uncover only the wells that you want to use. Vortex SPRI beads till even, and then add 90  $\mu$ l of SPRI Ampure XP Beads to the wells of the 96-well plate.
2. Transfer the entire PCR product including the SPRI beads (from Section IV.G, Step 3) to the wells of the plate containing the SPRI beads (from Step 1 above). Pipette the entire volume up and down 10 times to mix thoroughly. Incubate at room temperature for 8 min to let the DNA bind to the beads.

**NOTE:** The beads are viscous; suck the entire volume up, and push it out slowly.

3. Place the 96-well plate on the Ambion Magnetic Stand-96 for ~5 min or longer, until the liquid appears completely clear, and there are no beads left in the supernatant.
4. While the plate is sitting on the magnetic stand, pipette out the supernatant.
5. Keep the plate on the magnetic stand. Add 200  $\mu$ l of freshly made 80% Ethanol to each sample without disturbing the beads to wash away contaminants. Wait for 30 seconds and carefully pipette out the supernatant. DNA will remain bound to the beads during the washing process.
6. Repeat Step 5 one more time.
7. Seal the sample wells on the plate and briefly spin down for 10 seconds at 1,000 rpm to collect the liquid at the bottom of the well.
8. Place the 96-well plate on the magnetic stand for 30 seconds, then remove all the remaining Ethanol.
9. Place the plate at room temperature for ~3-5 min until the pellet appears dry. You may see a tiny crack in the pellet.

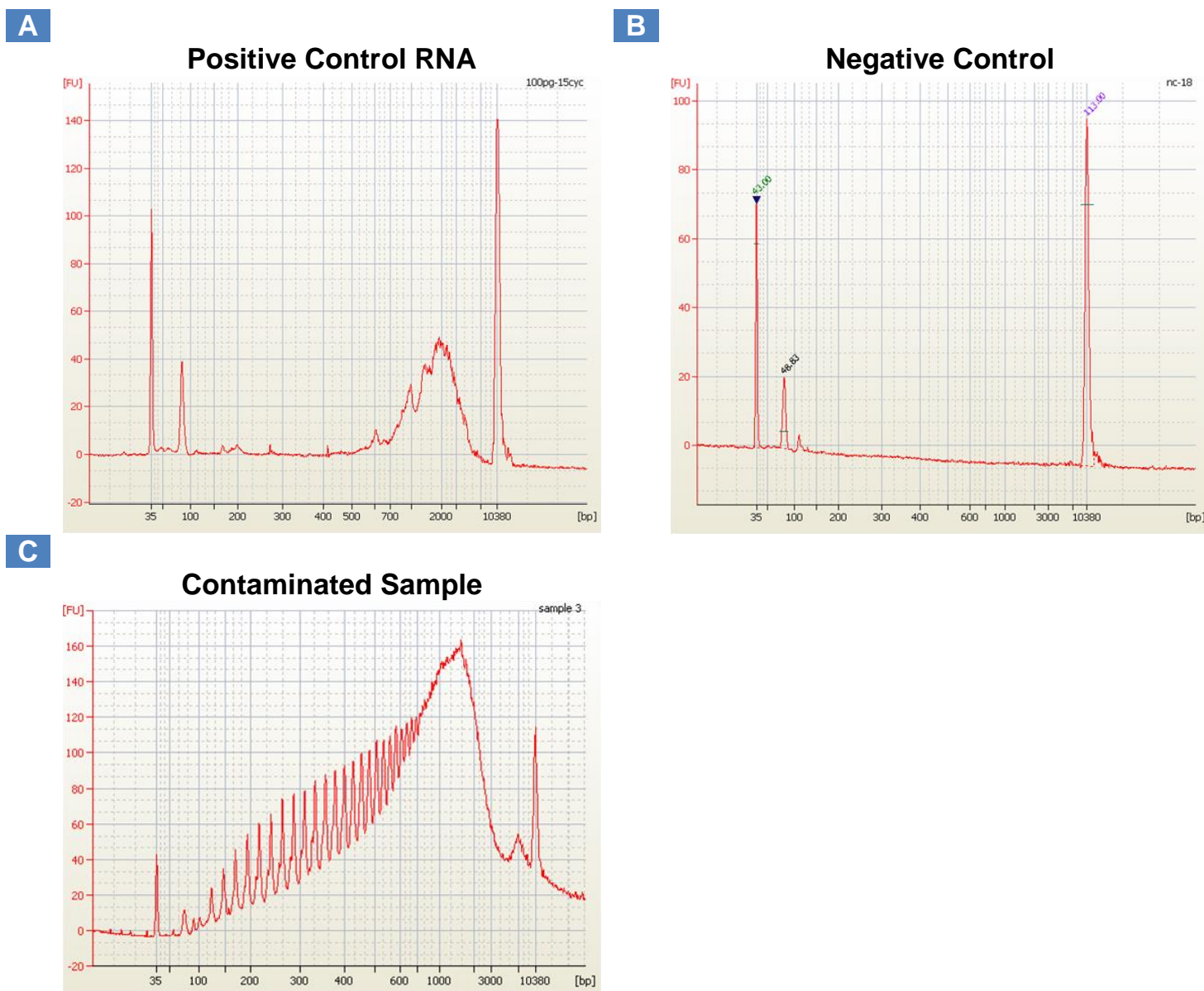
**NOTE:** If you over-dried the beads, you will see many cracks in the pellet. If it is under-dried, the DNA recovery rate will be lower because of the remaining Ethanol.

10. Once the beads are dried, add 12  $\mu$ l of Purification Buffer to cover the beads. Remove the plate from the magnetic stand and incubate at room temperature for 2 min to rehydrate.
11. Mix the pellet by pipetting up and down 10 times to elute DNA from the beads, then put the plate back on the magnetic stand for 1 minute or longer until the solution is completely clear.
12. Transfer clear supernatant containing purified cDNA from each well to a nuclease-free nonsticky tube. Label each tube with sample information and store at  $-20^{\circ}\text{C}$ .



## B. Validation Using the Agilent 2100 BioAnalyzer

1. Aliquot 1  $\mu$ l of the amplified cDNA for validation using the Agilent 2100 BioAnalyzer and the High Sensitivity DNA Chip from Agilent's High Sensitivity DNA Kit (Cat. No. 5067-4626). See the user manual for the Agilent High Sensitivity DNA Kit for instructions.
2. Compare the results for your samples & controls (see Figure 5) to verify whether the sample is suitable for further processing. Successful cDNA synthesis and amplification should yield no product in the negative control (Figure 5, Panel B), and a distinct peak spanning 400 bp to 9000 bp, peaked at  $\sim$ 2000 bp for the control RNA sample (Figure 5, Panel A), yielding approximately 2–7 ng of cDNA (depending on the input). Contaminated samples will have a broader peak, and abnormally high yield (Figure 5, Panel C).
3. Proceed to **Section VI: Covaris Shearing**.



**Figure 5. Electropherogram example results from Agilent 2100 Bioanalyzer.** All samples were subjected to SMARTer cDNA synthesis and amplification as described in the protocol. FU = fluorescence absorption units. **Panel A (top left).** Clean SMARTer Amplification Product (15 PCR cycles) **Panel B (top right).** Clean SMART Negative Control (18 PCR cycles). **Panel C (bottom).** Example of Contaminated SMARTer Amplification Product.



## VI. Covaris Shearing

### A. Protocol: Covaris Shearing of Full-length cDNA

Prior to generating the final library for Illumina sequencing, the Covaris AFA system is used for controlled DNA shearing. The resulting DNA will be in the 200–500 bp size range.

**NOTE:** The full-length cDNA output of the SMARTer Ultra Low Input RNA Kit for Illumina Sequencing can be processed using either the following protocol or [our modified protocol for the Nextera DNA Sample Preparation kits from Illumina](#).

1. Turn power ON for the Covaris system and the main cooler. Add about 1.9 L of distilled or deionized water to the water bath. The water level in the cooler should be within +/- 3 mm of the “FULL” waterline when the transducer is submerged. If needed, add distilled or deionized water to the water bath until the “FULL” line is reached.

**Important:** Never run a process without the water bath. This will permanently damage the transducer.

2. Close the door and open the Sonolab software. Click “ON” for the degassed button, and degas the water bath for ½ hour (30 minutes).
3. Add 65 µl of Purification Buffer to the DNA from Section V.A., Step 12. Transfer 75 µl of the Purification Buffer + DNA mixture into the 130-µl Covaris tube. Put the sample tubes into the appropriate location on the Sample holder.
4. Set up the process configuration panel based on the following table:

Table 3. Process Configuration Panel Set Up

Duty %	Intensity	Burst Cycle	Time (min)	Mode
10	5	200	5 min	Frequency Sweeping

5. Save the file and click return to go back to the main page.
6. Open the door. Place the tube holder with sample tubes on the transducer positioning system.
7. Close the door.
8. Click “START” on the main page to run the process.
9. After shearing is complete, transfer 75 µl of sheared DNA to 1.5 ml tubes.
10. Proceed to generate an Illumina Sequencing Library with the Low Input Library Prep Kit (Cat. No. 634947). Dispose all tubes and pipettes that have been exposed to amplicons in a sealed trash bag.

## VII. Appendix A: PCR Clean Work Station Guidelines

### A. Equipment Needed in the PCR Clean Work Station

**IMPORTANT:** The following equipment must be kept in the PCR Clean Work Station only. Don’t share any equipment/reagents between the clean room and the general lab. You can only move materials/supplies from the clean room to the general lab, NOT the other way around. Do not use equipment from the general lab for the portions of the protocol that require a PCR Clean Work Station.

- Single channel pipette: 10 µl, 20 µl and 200 µl, one each
- Eight channel pipette: 20 µl and 200 µl, one each
- Filter pipette tips: 10 µl, 20 µl and 200 µl, one box each
- One dedicated PCR thermal cycler used only for first-strand synthesis

- One QuickSpin minicentrifuge for 1.5-ml tube & one QuickSpin minicentrifuge for 0.2-ml tube
- MagnaBot II Magnetic Separation Device • (Promega Part No. V8351)
- One marker pen
- Small 1.5-ml tube rack
- Small bottle of nuclease-free water
- One bag of 1.5-ml nuclease-free tubes
- One bag of 8-strip nuclease-free 0.2-ml thin wall PCR tubes with caps
- DNA-OFF Solution (Takara Cat. No. 9036)

### B. PCR Clean Work Station Operation Instructions

**IMPORTANT:** At the beginning of each section of the protocol listed below, be sure to put on a clean pair of gloves and sleeve covers, then turn on the light and blower on the PCR clean station and lift the cover window. Bring a clean trash bag into the PCR clean station as well as two pieces of kimwipes to cover the work area. When done using the PCR Clean Work Station (after Section IV.G.), wipe the working area dry with Kimwipes, and clean with 70% EtOH. Once a week, clean the working area with DNA-OFF Solution.

#### Clean Work Station Instructions for First-Strand cDNA Synthesis (Section IV.E):

1. Bring the IsoFreeze PCR Rack and the Flipper Rack containing the SMARTer cDNA synthesis reagents from the  $-20^{\circ}\text{C}$  freezer to the PCR clean work station, and put them on top of the kimwipes.
2. Take the cell or RNA samples from  $-80^{\circ}\text{C}$  and put them on the IsoFreeze PCR Rack in the PCR clean work station.
3. Set up the cDNA synthesis reaction(s), and put the sample(s) in the PCR thermal cycler for cDNA synthesis.
4. Place all the used tips and tubes in a trash bag, remove from the PCR clean work station, and dispose.
5. Put the IsoFreeze PCR Rack and the Flipper Rack containing the SMARTer cDNA synthesis reagents back in the  $-20^{\circ}\text{C}$  freezer.
6. Close the cover window of the PCR clean work station, leave the blower on, and dispose of gloves and sleeve covers.

#### Clean Work Station Instructions for Purification of fs cDNA & LD-PCR (Sections IV.F & IV.G):

1. Bring the Ampure XP beads from  $4^{\circ}\text{C}$  to the PCR clean work station. Aliquot the beads into 2 tubes: one for the first-strand cDNA purification (Section IV.F) and one for the PCR product purification (Section V.A) prior to the experiment to avoid a cross-contamination.
2. Bring the IsoFreeze Flipper Rack containing the SMARTer PCR reagents from the  $-20^{\circ}\text{C}$  freezer to the PCR clean work station, and put it on top of the kimwipes.
3. Set up the PCR reactions.
4. Take the PCR reaction tubes out of the PCR clean work station, and perform PCR in a thermal cycler in the general lab.
5. Place used tips and tubes in a trash bag, remove from the PCR clean work station, and dispose.
6. Put the IsoFreeze Flipper Rack containing the SMARTer PCR reagents back in the  $-20^{\circ}\text{C}$  freezer and put the Ampure XP beads back in the  $4^{\circ}\text{C}$  refrigerator.
7. Close the cover window of the PCR clean work station. Turn off the light and blower, and turn on the UV light.

## VIII. Appendix B: Working with Cells

### A. Protocol: Screen for Cell Culture Media

**Important:** When working with cultured cells you need to ensure that the cell culture medium does not inhibit first-strand cDNA synthesis. Therefore it is important to select a medium with the least inhibitory effect.

We recommend testing your cells' compatibility with the media listed in Table 4. These media have been tested with this protocol.

Table 4. Validated Media

Superblock (Pierce Cat. No. 37515)
0.1 ml of [DMEM F12 + Glutamax (Invitrogen Cat. No. 10565)] + 3.6 µl of 25% BSA (Invitrogen Cat. No. A10008-01)
For 1 L of PBS Buffer (0.2 micron filtered): 0.2 g of KCl 0.24 g of KH <sub>2</sub> PO <sub>4</sub> (anhydrous) 8.00 g of NaCl 1.44 g of Na <sub>2</sub> HPO <sub>4</sub> (anhydrous) add dH <sub>2</sub> O to 1 L

### B. Cell Preparation for First-Strand cDNA Synthesis

**Important:** The cDNA synthesis protocol has been tested with suspension cells without internal labeling. Any cell that has gone through fixation won't work (the RNA won't release efficiently).

1. Pick cell(s) in 1 µl of validated media (see Section A. Screen for Cell Culture Media), and then transfer to a 0.2 mL RNase-free PCR tube containing 2.5 µl of Reaction Buffer. Put the tube(s) on dry ice immediately.

#### NOTES:

- The total volume of suspended cells should not exceed 1–2 µl. If necessary, you can spin down your cells and resuspend them in Reaction Buffer.
  - If you choose to transfer your cell(s) in 2 µl of media, you need to use double the amount of Reaction Buffer (5 µl). You must also double the volumes for all of the components in the first-strand synthesis reaction (Section IV.E, Steps 1–6), as well as double the volume of beads for the first SPRI purification (Section IV.F, Step 1).
2. Proceed with **Section IV.E: Protocol: First-Strand cDNA Synthesis**. If you will not be performing cDNA synthesis immediately, store the tube on dry ice or at –80°C until use.

## IX. Appendix C: PCR Optimization

If you have a sufficient amount of starting material (>1 ng total RNA), we recommend optimizing the PCR cycling parameters for your experiment. If you have a very limited amount of material or your sample is unique, use a similar source of RNA or cells to perform PCR cycle optimization prior to using the actual sample. Choosing the optimal number of PCR cycles ensures that the ds cDNA will remain in the exponential phase of amplification. When the yield of PCR products stops increasing with more cycles, the reaction has reached its plateau. Overcycled cDNA can result in a less representative sample. Undercycling, on the other hand, results in a lower yield of cDNA. The optimal number of cycles for your experiment is one cycle fewer than is needed to reach the plateau. Be conservative: when in doubt, it is better to use fewer cycles than too many.

To perform PCR cycle optimization, prepare several tubes containing an amount of RNA equal to your sample amount. Subject each tube to a different range of cycles. For example if you have 1 ng of RNA, subject one tube to the recommended a number of cycles. Subject the other two tubes to 2–3 cycles fewer or more than the first tube, i.e. 15, 12 (recommended for 1 ng), and 10 cycles.

1. Use the following program for thermal cycling:

95°C	1 min	
X cycles:		
95°C	15 sec	}
65°C	30 sec	
68°C	6 min	
72°C	10 min	
4°C	forever	

2. Perform **Purification of ds cDNA using SPRI Ampure Beads** (Section V.A).
3. Run samples on an Agilent High Sensitivity DNA Chip using the Agilent 2100 Bioanalyzer to evaluate DNA output. See the user manual for the Agilent High Sensitivity DNA Kit for instructions.
4. Determine the optimal number of cycles required for each experimental and control sample. We recommend using the lowest PCR cycle number that generates enough material for Illumina library construction.
5. Apply the optimal number of PCR cycles to your sample material.

## X. Appendix D: Troubleshooting Guide

**Table 5. Troubleshooting Guide for SMARTer cDNA Synthesis & Amplification**

Problem	Cause	Solution
No cDNA yield with the control RNA	Control RNA is degraded	The control RNA is vigorously tested for RNase contaminations and provided in high concentrations (1 µg/µl) to reserve its integrity. It is essential that any item that could contact the control RNA is RNase-free.
	Control RNA is too diluted	Make sure to use recently calibrated pipettes to accurately perform serial dilutions of the control RNA. Prepare fresh dilutions of the Control RNA and try again. The diluted RNA is less stable. Therefore, avoid using previously diluted low concentration RNA samples.
	cDNA synthesis reaction failed	Carefully check the protocol and make sure all components are added in a right order; Failure of any components of the reaction will result in a failed cDNA synthesis. Make sure to prepare all the necessary items prior cDNA synthesis and to spin down tubes prior opening them. Do not delay cDNA synthesis reaction after RNA is denatured to jump start cDNA synthesis.
	PCR failed	Make sure to use the recommended PCR enzyme. Before using the PCR enzyme check it with previously tested primers and template.
	cDNA was not efficiently bound to the SPRI beads during purification steps	Make sure beads are mixed correctly after adding them to DNA samples. Vortexing the beads once they are added to samples can shear the DNA or break it free from the beads. Make sure the SPRI beads are collected correctly on the magnetic stand and given enough time to completely separate the liquid phase. Make sure not to disturb the beads while removing the supernatant. Supernatant that is not completely removed from SPRI beads may inhibit downstream reactions.

## SMARTer Ultra Low RNA Kit for Illumina Sequencing User Manual

Problem	Cause	Solution
No cDNA yield with the experimental RNA, but control is OK	RNA is low quality, degraded or has impurities that inhibit the cDNA synthesis reaction	If possible, check RNA quality prior to cDNA synthesis. Make sure RNA is good quality. Certain tissues, such as plants, have a high level of polysaccharides that interfere with first-strand synthesis. Make sure to use appropriate RNA isolation kits for each given tissue/species.
Low cDNA yield with the control RNA	Control RNA is too diluted	Make sure to use recently calibrated pipettes to accurately perform serial dilutions of the control RNA. Prepare fresh dilutions of the Control RNA and try again. Diluted RNA is less stable therefore avoid using previously diluted low concentration RNA samples.
	cDNA was not efficiently bound to the SPRI beads during purification steps	Make sure the SPRI beads are collected correctly on the magnetic stand and given enough time to completely separate the liquid phase. Make sure not to disturb beads while removing the supernatant. Repeat cDNA synthesis with the freshly diluted control RNA.
	SPRI beads were overdried during purification of amplified cDNA	Don't allow the post-wash bead pellet to be overdried. It's important to allow any residual ethanol to dry, but overdrying (too long, or drying at a temperature higher than room temperature) will make it difficult to elute the DNA from the beads.
Low cDNA yield with the experimental RNA, but control is OK	Insufficient amount of starting material or RNA is degraded or has impurities that inhibit the cDNA synthesis reaction	If more RNA sample is available, quantitate RNA prior to cDNA synthesis. Follow PCR cycle guidelines appropriate for the starting amount of sample you are using.

## XI. References

- Barnes, W. M. (1994) PCR amplification of up to 35-kb DNA with high fidelity and high yield from  $\lambda$  bacteriophage templates. *Proc. Natl. Acad. Sci. USA* **91**:2216–2220.
- Chenchik, A., Zhu, Y., Diatchenko, L., Li, R., Hill, J. & Siebert, P. (1998) Generation and use of high-quality cDNA from small amounts of total RNA by SMART PCR. In *RT-PCR Methods for Gene Cloning and Analysis*. Eds. Siebert, P. & Larrick, J. (BioTechniques Books, MA), pp. 305–319.
- Kellogg, D. E., Rybalkin, I., Chen, S., Mukhamedova, N., Vlasik, T., Siebert, P. & Chenchik, A. (1994) TaqStart Antibody: Hot start PCR facilitated by a neutralizing monoclonal antibody directed against *Taq* DNA polymerase. *BioTechniques* **16**:1134–1137.

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: 800.424.1350 (toll-free)
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.

Illumina®, Genome Analyzer, HiSeq®, and Nextera® are trademarks of Illumina, Inc. Takara®, the Takara logo, and DNA-OFF are trademarks of Takara Holdings, Inc. Clontech®, the Clontech logo, Advantage®, SMART®, SMARTer®, SMARTScribe, and Ultra® are trademarks of Clontech Laboratories, Inc. All other trademarks are the property of their respective owners. Certain trademarks may not be registered in all jurisdictions. Clontech Laboratories, Inc. is a Takara Bio Company. ©2016 Clontech Laboratories, Inc.

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