

MessageAmp™ II-96 Kit

(Cat #AM1819)

Instruction Manual

I.	Introduction	1
	A. Product Description and Background	
	B. Materials Provided with the Kit and Storage	
	C. Materials Not Provided with the Kit	
	D. Related Products Available from Ambion	
II.	aRNA Amplification Protocol	8
	A. Planning the Experiment	
	B. Reagent Preparation	
	C. Reverse Transcription to Synthesize First Strand cDNA	
	D. Second Strand cDNA Synthesis	
	E. cDNA Purification	
	F. In Vitro Transcription to Synthesize aRNA	
	G. aRNA Purification	
III.	aRNA Evaluation and Preparation for Array Hybridization	19
	A. aRNA Quantitation	
	B. Analysis of aRNA Size with a Bioanalyzer	
	C. Preparing Labeled aRNA for Hybridization	
IV.	Troubleshooting	21
	A. Positive Control Reaction	
	B. Factors that Affect Both the Positive Control and Experimental Samples	
	C. Troubleshooting Low Yield and Small Average aRNA Size	
	D. Bead Carryover in Eluted Sample	
V.	Appendix	25
	A. IVT Reaction Using 10 mM Biotin-labeled CTP and UTP	
	B. References	
	C. MessageAmp II-96 Kit Contents	
	D. Quality Control	

Manual 1819M Revision C

Revision Date: December 7, 2010

For research use only. Not for use in diagnostic procedures. By use of this product, you accept the terms and conditions of all applicable Limited Label Licenses. For statement(s) and/or disclaimer(s) applicable to this product, see below.

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document.

Applied Biosystems disclaims all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. In no event shall Applied Biosystems be liable, whether in contract, tort, warranty, or under any statute or on any other basis for special, incidental, indirect, punitive, multiple or consequential damages in connection with or arising from this document, including but not limited to the use thereof.

Literature Citation: When describing a procedure for publication using this product, we would appreciate that you refer to it as the MessageAmp™ II-96 Kit.

If a paper that cites one of Ambion's products is published in a research journal, the author(s) may receive a free Ambion T-shirt by sending in the completed form at the back of this instruction manual, along with a copy of the paper.

Warranty and Liability: Applied Biosystems is committed to delivering superior product quality and performance, supported by industry-leading global service and technical support teams. Warranty information for the accompanying consumable product is available at www.ambion.com/info/warranty in "Limited Warranty for Consumables," which is subject to the exclusions, conditions, exceptions, and limitations set forth under the caption "EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS" in the full warranty statement. Please contact Applied Biosystems if you have any questions about our warranties or would like information about post-warranty support.

Patents and Licensing Notifications: The MessageAmp™ II-96 Kit is covered by US patents.

This product or portions thereof is manufactured and sold under license from GE Healthcare under U.S. Patent Nos. 5,523,231 and 5,681,946 and other foreign patents.

End User Terms And Conditions Acceptance. These terms and conditions shall govern the purchase, use, transfer and acceptance of the products described in the purchase order, quotation or invoice, which products are sold and distributed by Applied Biosystems Corporation to the buyer/transferee of such products (the "End User"). The transfer/sale of products to the End User is expressly conditional upon End User's acceptance of these terms and conditions. **Restrictions on Use.** End Users are specifically not authorized to and are forbidden from reselling, transferring or distributing any products either as a stand alone product or as a component of another product. The right to use the products does not, in and of itself, include or carry any right of the End User to any GE Healthcare Bio-Sciences Corp.'s technology or intellectual property other than expressly provided herein. End Users may not attempt to reverse engineer parameters of any of GE Healthcare Bio-Sciences Corp. proprietary products or services. **DISCLAIMER OF WARRANTIES.** GE HEALTHCARE BIO-SCIENCES CORP. PROVIDES NO WARRANTIES TO END USER (STATUTORY OR IMPLIED), INCLUDING WITHOUT LIMITATION, AS TO PRODUCT QUALITY, CONDITION, DESCRIPTION, MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, AND ALL SUCH WARRANTIES ARE HEREBY EXPRESSLY DISCLAIMED. GE HEALTHCARE BIO-SCIENCES CORP. HEREBY EXPRESSLY DISCLAIMS ANY WARRANTY REGARDING RESULTS OBTAINED THROUGH THE USE OF THE PRODUCTS, INCLUDING WITHOUT LIMITATION ANY CLAIM OF INACCURATE, INVALID OR INCOMPLETE RESULTS. **Exclusion of Liability.** GE Healthcare Bio-Sciences Corp. and its affiliates shall have no liability to an End User, including, without limitation, for any loss of use or profits, business interruption or any consequential, incidental, special or other indirect damages of any kind, regardless of how caused and regardless of whether an action in contract, tort, strict product liability or otherwise.

Trademarks: Applied Biosystems, AB (Design), Ambion, MEGAscript, RNAqueous, and RNaseZap are registered trademarks, and ArrayControl, ArrayScript, MagMAX, MEGAclean, MessageAmp, and SlideHyb are trademarks of Applied Biosystems Corporation or its subsidiaries in the US and/or certain other countries.

Cy and CyDye are trademarks of GE Healthcare. All other trademarks are the sole property of their respective owners.

All other trademarks are the sole property of their respective owners.

© 2007, 2010 Ambion, Inc. All Rights Reserved.

I. Introduction

A. Product Description and Background

The MessageAmp™ II-96 Kit is designed for amplification of RNA in a 96-well format. The procedure is based on the RNA amplification protocol developed in the laboratory of James Eberwine (Figure 1 on page 3) (Van Gelder et al. 1990). consists of reverse transcription with an oligo(dT) primer bearing a T7 promoter using Array-Script™, a reverse transcriptase (RT) engineered to produce higher yields of first strand cDNA than wild type enzymes. ArrayScript catalyzes the synthesis of virtually full-length cDNA, which is the best way to ensure production of reproducible microarray samples. The cDNA then undergoes second strand synthesis and clean-up to become a template for in vitro transcription with T7 RNA Polymerase. To maximize aRNA yield, Ambion proprietary MEGAscript® in vitro transcription (IVT) technology is used in the kit to generate hundreds to thousands of anti-sense RNA copies of each mRNA in a sample. (In this Instruction Manual the antisense amplified RNA is referred to as aRNA, in scientific literature it is also commonly called cRNA.) The IVT can be configured to synthesize either biotin labeled aRNA, or unlabeled aRNA that can subsequently be labeled by reverse transcription (for example with fluorescently labeled dNTPs). The resulting aRNA is suitable for use on most commercially available microarray gene expression systems.

Benefits of RNA amplification

RNA amplification was originally developed as a method to expand very small RNA samples to produce enough material for array hybridization (Yue et al. 2001). Several groups have conducted studies to determine whether amplification of RNA introduces bias, and they report that any bias is minimal (Li et al. 2004, Feldman et al. 2002, and Polacek et al. 2003). Additionally, among the benefits of amplification is a more reproducible expression profile from a wide range of RNA inputs. Some researchers conclude that amplification actually improves the reliability of array results regardless of whether it is needed for sample expansion (Feldman et al. 2002 and Polacek et al. 2003). RNA amplification has become the standard method for preparing RNA samples for array analysis (Kacharmina et al. 1999, Pabon et al. 2001).

The MessageAmp II-96 advantage

Optimized protocols

All reagents and protocols have been optimized to produce high quality aRNA with maximum yield and reproducibility. The simple, rapid aRNA purification procedure prepares the aRNA for downstream applications such as reverse transcription or post-synthesis labeling reactions.

The reagents used for enzymatic reactions in the MessageAmp II-96 Kit are the same as those used in the single prep version of the kit (Cat #AM1751). However, magnetic beads are used for cDNA and aRNA cleanup steps in the high throughput kits to streamline the amplification process.

Built-in flexibility

The NTPs for IVT are provided in separate tubes so that modified nucleotides, such as biotinylated UTP or CTP, or cyanine 3/cyanine 5 CTP and UTP, can be used in the reaction to synthesize labeled aRNA.

Ready for automation

This protocol was designed with both manual and robotic liquid handling systems in mind. It can easily be adapted for use on a robotic platform; downloadable protocols for the Biomec 3000 and NX workstations are available on our website at:

www.ambion.com/techlib/automation

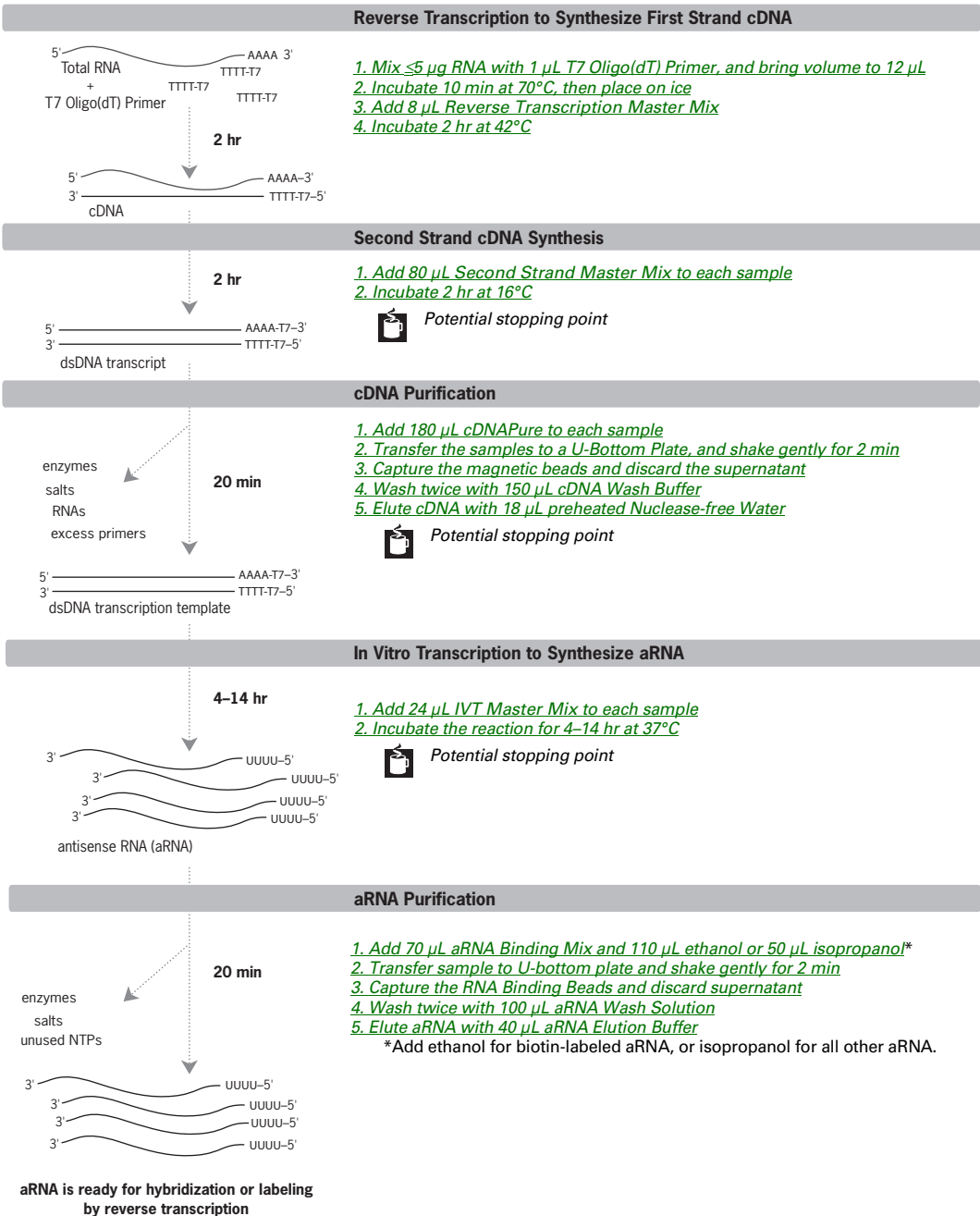
Compatible with second round amplification

This kit is designed for a single round of aRNA amplification, which provides sufficient aRNA for most uses. If more aRNA is needed, RNA can be subjected to two sequential rounds of amplification. If your experiments require two rounds of amplification, you will need to purchase the Ambion MessageAmp II aRNA Amplification Kit (Cat #AM1751). For the first round amplification, use the MessageAmp II Kit to make unmodified aRNA.

For the second round of amplification, use the MessageAmp II-96 Kit components plus the Second Round Primers from the MessageAmp II Kit, and follow instructions from both manuals as detailed below:

- For second round reverse transcription and second strand cDNA synthesis, follow the instructions in the MessageAmp II manual.
- For the second round amplification IVT reaction, follow the instructions in the MessageAmp II-96 Kit manual.

Figure 1. Procedure Overview



B. Materials Provided with the Kit and Storage

The MessageAmp II-96 Kit includes reagents for amplification of 96 samples using a single round of amplification in a 96-well plate. Reagents for 4 control reactions are also included.

cDNA synthesis and in vitro transcription reagents

Store at –20°C in a non-frost-free freezer.

Amount	Component
110 µL	T7 Oligo(dT) Primer*
110 µL	ArrayScript™
110 µL	RNase Inhibitor
220 µL	10X First Strand Buffer
880 µL	dNTP Mix
220 µL	DNA Polymerase
110 µL	RNase H
10 µL	Control RNA (1 µg/µL HeLa total RNA)
1.1 mL	10X Second Strand Buffer
440 µL	T7 Enzyme Mix
440 µL	T7 10X Reaction Buffer
440 µL	T7 ATP Solution (75 mM)
440 µL	T7 CTP Solution (75 mM)
440 µL	T7 GTP Solution (75 mM)
440 µL	T7 UTP Solution (75 mM)

* T7 Oligo(dT) Primer is also available separately (Cat #AM5710).

cDNA and aRNA purification components

Amount	Component	Storage
20 mL	Nuclease-free Water	any temp*
20 mL	cDNAPure†	4°C‡
1.1 mL	RNA Binding Beads	4°C‡
40 mL	cDNA Wash Buffer Concentrate Add 32 mL 100% ethanol before use	room temp
25 mL	aRNA Wash Solution Concentrate Add 20 mL 100% ethanol before use	room temp
450 µL	Bead Resuspension Solution	room temp
10 mL	aRNA Elution Buffer	room temp
5.5 mL	aRNA Binding Buffer Concentrate	room temp
2	U-Bottom Plate	room temp
4	PCR Plate	room temp
1	Reservoir	room temp

* Store Nuclease-free Water at room temp, 4°C, or –20°C.

† cDNAPure is manufactured by Agencourt Bioscience Corporation.

‡ Do not freeze.

C. Materials Not Provided with the Kit

Lab equipment and supplies

- ACS grade (or higher quality) alcohol:
 - 100% ethanol
 - 100% isopropanol
- Thermal cycler (recommended) or fixed temperature air incubators set at 70°C, 42°C, 37°C, and 16°C. (See Procedural notes on page [9](#) for more information.)
- Vortex mixer
- Orbital shaker for 96-well plates such as the Barnstead/Lab-Line Titer Plate Shaker (available from Fisher or VWR)
- Magnetic stand for 96-well plates (we recommend the Ambion 96-well Magnetic-Ring Stand, Cat #AM10050, for its high strength magnets and quality design)
- Microcentrifuge
- Aluminum sealing foil or tape for sealing reactions in 96-well plates during incubations

(optional) Materials and equipment for RNA analysis

- Spectrophotometer—such as the NanoDrop® ND-8000 UV-Vis Spectrophotometer
- RiboGreen® RNA quantitation reagent (Molecular Probes, Inc.)
- Bioanalyzer (Agilent) and RNA LabChip® Kits (Caliper Technologies)

(optional) Biotin labeled UTP

Biotin labeled UTP, for making biotinylated aRNA; Ambion recommends Biotin-11-UTP (Ambion Cat #AM8451, 75 mM) because it gives good incorporation, has minimal effect on aRNA recovery during purification, and results in high signal on most commercial microarrays. Biotin-16-UTP (Ambion Cat #AM8453, 75 mM) can also be used if it suits your experimental needs.

D. Related Products Available from Ambion

<p>96-well Magnetic-Ring Stand Cat #AM10050</p>	<p>The Ambion 96-well Magnetic-Ring Stand features 96 powerful ring-shaped magnets arranged to cradle each well of a 96-well plate for quick, thorough bead capture. Captured magnetic beads form evenly distributed donut-shaped pellets with a large hole in the center. This capture pattern facilitates both supernatant removal and subsequent bead resuspension. The stand is suitable for high throughput applications conducted with multichannel pipettors or with robotic liquid handlers. However, because the pellets will be evenly distributed around the edge of the wells, it may require practice for efficient manual removal of supernatants.</p>
<p>Magnetic Stand-96 Cat #AM10027</p>	<p>The Ambion Magnetic Stand-96 has powerful magnets positioned to capture beads to one side of the well. This capture pattern makes it very easy to remove supernatants manually without disturbing the beads, and therefore may be preferred by beginning users. In some applications, however, pellets formed with the Magnetic Stand-96 may be difficult to resuspend. If this occurs, we recommend the 96-well Magnetic-Ring Stand (Cat #AM10050).</p>
<p>MagMAX™ -96 Total RNA Isolation Kit Cat #AM1830</p>	<p>The MagMAX™-96 Total RNA Isolation Kit is a magnetic bead based total RNA purification system designed for rapid high throughput processing of cells in 96-well plates. High yield and high quality total RNA can be obtained from 100 to 500,000 cultured eukaryotic cells. The kit can also be used for total RNA isolation from small tissue samples.</p>
<p>MagMAX™ -96 for Microarrays Total RNA Isolation Kit Cat #AM1839</p>	<p>The MagMAX-96 for Microarrays Total RNA Isolation Kit is designed for rapid, high throughput isolation of total RNA from mammalian cells and tissues in 96-well plates. The procedure employs the robust and reliable lysis/denaturant, TRI Reagent® and Ambion MagMAX magnetic bead-based RNA purification technology to provide high yields of pure, intact RNA that can be used directly for quantitative reverse transcriptase PCR (qRT-PCR) and microarray analysis. The kit can also be used to process fewer than 96 samples at one time.</p>
<p>RNAqueous®-96 Automated Kit Cat #AM1812</p>	<p>The RNAqueous-96 Automated Kit employs a simple and rapid procedure to purify total RNA from source material without using organic solvents (such as phenol). It is formatted specifically for high throughput RNA isolation on robotic platforms.</p>
<p>ArrayControl™ Cat #AM1780</p>	<p>The ArrayControl Spikes are a set of eight control RNA transcripts designed for the normalization and validation of glass microarray experiments. The Spikes range in size from 750 to 2000 bases and each transcript has a 30-base 3' poly(A) tail. The precisely quantitated RNA Spikes are designed to be added to your RNA sample before labeling, to serve as internal controls for sample labeling and hybridization efficiency.</p>
<p>Amino Allyl cDNA Labeling Kit Cat #AM1705</p>	<p>The Amino Allyl cDNA Labeling Kit generates cDNA for secondary fluorescent dye labeling to be used for glass array analysis. It includes all the reagents, except the amine-reactive labeling moiety (e.g. cyanine dyes) for 2-step labeling of cDNA. The reaction produces more labeled cDNA, more efficiently than direct dye incorporation.</p>

5-(3-aminoallyl)-dUTP Cat #AM8439	This 50 μ M solution of amino allyl modified dUTP can be used with the RETROscript [®] Kit (Cat #AM1710) to synthesize amine-reactive cDNA from aRNA. The amine-reactive cDNA can then be postlabeled with any amine-reactive label moiety.
Biotin-11-UTP and Biotin-16-UTP Cat #AM8450, AM8451, AM8452, AM8453	Ambion biotinylated UTPs are ideal for use as substrates in vitro transcription reactions, and can be utilized by a variety of RNA polymerases, including T7, T3, and SP6 RNA polymerases. Biotinylated RNA can be used in place of radioactively labeled RNA in many applications with detection via one of a variety of streptavidin-based methods.
RNA Fragmentation Reagents Cat #AM8740	Amplified RNA is commonly fragmented prior to hybridization on oligonucleotide microarrays to improve the hybridization kinetics and signal produced on oligonucleotide microarrays. Ambion RNA Fragmentation Reagents include a 10X Fragmentation Reagent and a Stop Solution.
SlideHyb[™] Glass Array Hybridization Buffers and Glass Array Hybridization Cassette see our web or print catalog	There are 3 unique SlideHyb Glass Array Hybridization Buffers; they have identical salt and formamide compositions, but differ in hybridization kinetics and blocking reagents. Ambion also offers the Glass Array Hybridization Cassette for incubation of glass microarray hybridization reactions.

II. aRNA Amplification Protocol

A. Planning the Experiment

We understand that the needs of our high throughput customers are very diverse; if you have questions that are not covered in this Instruction Manual, please contact our high throughput experts by emailing: automation@ambion.com

RNA quantity and IVT incubation time

The MessageAmp II-96 procedure can accommodate a wide range of input RNA amounts, but for reproducible and comparable results, use a fixed amount of input RNA for all experiments. Typically **100 ng–5 µg of total RNA** should be used in a MessageAmp II-96 amplification reaction. *The recommended amount of RNA per reaction is 1 µg of total RNA.*

The amount of input RNA needed will depend on how much aRNA is required for your experiments; whether total RNA or poly(A) RNA is used; and how long the IVT amplification is incubated.

Using 1 µg or more total RNA and a 4 hr IVT incubation typically produces enough aRNA for several microarray hybridizations. When amplifying small RNA samples (e.g. ~250 ng or less), however, incubating the IVT reaction for 14 hr will maximize the amount of aRNA produced.



NOTE

Note that aRNA yield may vary considerably depending on the purity and integrity of the RNA used in the reaction and on the other parameters discussed in these instructions.

Other input RNA requirements

RNA purity

RNA samples should be free of contaminating proteins, DNA, and other cellular material as well as phenol, ethanol, and salts associated with RNA isolation procedures. Impurities can lower the efficiency of reverse transcription and subsequently reduce the level of amplification. An effective measure of RNA purity is the ratio of absorbance readings at 260 and 280 nm. The ratio of A_{260} to A_{280} values should fall in the range of 1.7–2.1.

RNA integrity

The integrity of the RNA sample, or the proportion of full-length mRNA, can significantly impact the amplification reaction. Reverse transcribing partially degraded mRNAs will typically generate relatively short cDNAs that lack the sequence upstream of the break in the RNA molecule.

The most common method for evaluating the relative integrity of mRNA is to measure the ratio of the 28S and 18S ribosomal RNAs (rRNA) using an RNA LabChip Kit with an Agilent 2100 bioanalyzer. The RNA LabChip fractionates RNA molecules according to size, and the relative amounts of 18S and 28S rRNA can be determined. The ratio of the 28S to 18S rRNA peaks will approach 2:1 in RNA samples comprising primarily full-length RNA (Figure 2).

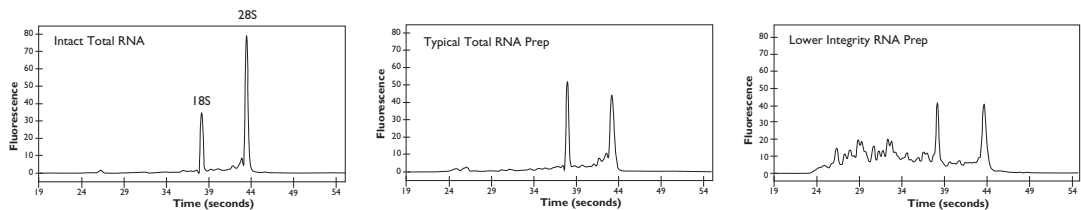


Figure 2. Bioanalyzer Images of Increasingly Degraded Total RNA

These electropherograms from the Agilent 2100 bioanalyzer show progressive RNA degradation. Notice that the ribosomal RNA (rRNA) peaks are at a ratio of about 2:1 (28S:18S) in an intact total RNA sample (left). Total RNA that is somewhat degraded shows nearly equal amounts of 28S and 18S rRNA (middle). Further degradation decreases the rRNA peaks, and degradation products become apparent in the lower molecular weight range (right).

Procedural notes

We recommend a thermal cycler for all enzymatic reaction incubations

Ideally, reactions should be incubated in a calibrated thermal cycler with a lid temperature that adjusts to the block temperature (typically 5°C above the block temperature). If the lid temperature is static (~100°C), turn off the lid heat or do not close the heated lid onto your samples. Otherwise, heat from the lid will raise the temperature of the solution in the tube, compromising the reaction.

If a thermal cycler is not available, you may use a calibrated hybridization oven or other constant temperature air incubator for second strand cDNA synthesis and IVT reactions (see next section). We do *not* recommend using heat blocks or water baths for any MessageAmp II-96 reaction incubations.

Use calibrated, preheated incubators for enzymatic reaction incubations (Alternative to thermal cycler)

The MessageAmp II-96 procedure is very sensitive to temperature; therefore use incubators that have been professionally calibrated according to the manufacturer's recommended schedule. Variable or inaccurate incubation temperatures can limit aRNA synthesis.

Preheat incubators if necessary so that the correct temperature has stabilized before reactions are placed in the incubator. It is also very important that condensation does not form in the plate wells during any of the incubations. Condensation changes the composition of reaction mixtures and can greatly reduce yield.

Orbital shaker settings

We recommend using an orbital shaker with a 2–3 mm orbit. Using too large an orbit will cause the liquid to spill out of the wells, while too small of an orbit may not mix the sample effectively.

Determine the maximum speed that can be used in this procedure by filling the wells of a 96-well plate with 290 μ L of water. Turn the shaker on and increase the speed slowly while observing the movement of the water in the wells. Note the maximum speed at which the plate can be shaken without causing the water to splash or spill out of the wells. When shaking experimental samples, always increase the orbital shaker speed slowly to avoid splashing. Always choose the lowest speed that is strong enough for complete mixing.

Master mixes

We strongly recommend preparing master mixes, instead of pipetting components individually, to improve pipetting consistency. When using a multichannel pipettor to dispense a master mix that consists of <2 mL, we recommend that you minimize the void volume by aliquoting the master mix into a PCR plate instead of using a reservoir.

Ambion provides a web-based interactive calculator to calculate master mix quantities at the following address:www.ambion.com/tools/ma296

We recommend using the web calculator to identify reagent volumes needed for all the master mixes in the procedure and printing the results to accompany this manual before starting the protocol.

B. Reagent Preparation

1. Add 32 mL 100% ethanol to the cDNA Wash Buffer Concentrate

Add 32 mL 100% ethanol (ACS grade or better) to the bottle labeled cDNA Wash Buffer Concentrate. Mix well and mark the label to indicate that the ethanol was added. This solution will be referred to as *cDNA Wash Buffer* in these instructions.

Store at room temperature.

2. Add 20 mL 100% ethanol to the aRNA Wash Solution Concentrate

Add 20 mL 100% ethanol (ACS grade or better) to the bottle labeled aRNA Wash Solution Concentrate. Mix well and mark the label to indicate that the ethanol was added. This solution will be referred to as *aRNA Wash Solution* in these instructions.

Store at room temperature.

3. Prepare aRNA Binding Mix for the experiment



IMPORTANT

Prepared aRNA Binding Mix can be stored at room temp for one week; prepare only the amount needed for all samples in the experiment plus ~10% overage to cover pipetting error. We provide a master mix calculator on our website to calculate reagent amounts:

www.ambion.com/tools/ma296

a. First, mix RNA Binding Beads with Bead Resuspension Solution and ethanol

Mixing beads with the Bead Resuspension Solution will facilitate bead dispersion in Binding Mix. At room temp, combine Bead Resuspension Solution with RNA Binding Beads in a nuclease-free tube and mix briefly. Then add 100% ethanol and mix thoroughly by vortexing.

a. Combine the following (for a single reaction):	
RNA Binding Beads*	10 μ L
Bead Resuspension Solution	4 μ L
b. Mix briefly, then add:	
100% ethanol	6 μ L
c. Mix well by vortexing.	

* Mix the RNA Binding Beads by vortexing before dispensing.

b. Next, add aRNA Binding Buffer Concentrate

At room temp, add aRNA Binding Buffer Concentrate to the bead mixture from the previous step in a nuclease-free container. Mix by vortexing:

aRNA Binding Mix (for a single reaction)	
Component	Amount
aRNA Binding Buffer Concentrate	50 μ L
RNA Binding Bead Mixture (from step 3.a)	20 μ L

This mixture is referred to as aRNA Binding Mix in these instructions.

C. Reverse Transcription to Synthesize First Strand cDNA

Incubators needed:

- 70°C: thermal cycler recommended
- 42°C: thermal cycler recommended

1. Mix $\leq 5 \mu$ g RNA with 1 μ L T7 Oligo(dT) Primer, and bring volume to 12 μ L

- Place up to 5 μ g of total RNA (typically 100 ng–5 μ g) into the wells of a PCR Plate.
- Add 1 μ L T7 Oligo(dT) Primer.

- c. Add Nuclease-free Water to a final volume of 12 µL and cover with aluminum sealing foil.



NOTE

If all samples use the same RNA volume, premix the Nuclease-free Water and T7 Oligo(dT) Primer to simplify the procedure.

2. Incubate 10 min at 70°C, then place on ice

- a. Incubate 10 min at 70°C.
- b. Place the samples on ice. If there is condensation on the foil, centrifuge briefly (~5 sec) to collect samples at the bottom of the wells.

3. Add 8 µL Reverse Transcription Master Mix

- a. Prepare **Reverse Transcription Master Mix** in a nuclease-free tube at room temp. Assemble enough master mix for all samples in the experiment plus ~10% overage to allow for pipetting error. We provide a Master Mix Calculator on our website to calculate reagent amounts: www.ambion.com/tools/ma296

At room temp, assemble the Reverse Transcription Master Mix in the order shown:

Reverse Transcription Master Mix (for a single 20 µL reaction)	
Amount	Component
2 µL	10X First Strand Buffer
4 µL	dNTP Mix
1 µL	RNase Inhibitor
1 µL	ArrayScript

- b. Mix well by gently vortexing. Centrifuge ~5 sec to collect the master mix at the bottom of the tube and place it on ice.
- c. Transfer 8 µL of the Reverse Transcription Master Mix to each RNA sample, mix by gently pipetting up and down 3–4 times. Cover the reactions with aluminum sealing foil.

To facilitate the transfer process, you can distribute the master mix into 8 wells of a PCR plate and use an 8-channel pipette to transfer the master mix.

4. Incubate 2 hr at 42°C

Incubate the reaction for 2 hr at 42°C. If condensation is apparent after the incubation, centrifuge the plate briefly (~5 sec) to collect the samples at the bottom of the wells.

Place the plate on ice, or change the thermal cycler temperature to 4°C, and immediately proceed to the second strand cDNA synthesis.

D. Second Strand cDNA Synthesis

Incubator needed:

16°C: thermal cycler recommended

1. Add 80 μL *Second Strand Master Mix* to each sample

- a. On ice, prepare a *Second Strand Master Mix* by adding the following reagents to a reservoir in the order listed below. Assemble enough master mix for all samples in the experiment plus ~10% overage to allow for pipetting error. We provide a master mix calculator on our website to calculate reagent amounts:

www.ambion.com/tools/ma296

Assemble the Second Strand Master Mix on ice in the order shown:

Second Strand Master Mix (for a single 100 μL reaction)	
Amount	Component
63 μL	Nuclease-free Water
10 μL	10X Second Strand Buffer
4 μL	dNTP Mix
2 μL	DNA Polymerase
1 μL	RNase H

- b. Mix gently by rocking the reservoir back and forth.
- c. Transfer 80 μL of Second Strand Master Mix to each sample in the PCR Plate. Mix by gently pipetting up and down 3–4 times. Cover the reactions with aluminum sealing foil.

2. Incubate 2 hr at 16°C

Incubate in a 16°C thermal cycler (do not incubate in a heat block in a 4°C refrigerator because the temperature may fluctuate). It is important to cool the thermal cycler to 16°C before placing the PCR Plate inside; subjecting the reactions to temperatures >16°C could compromise aRNA yield.

After the 2 hr incubation at 16°C, place the reactions on ice and proceed to section [E. cDNA Purification](#) (below), or immediately freeze reactions at –20°C. Do not leave the reactions on ice for more than 1 hr.



STOPPING POINT

This is a potential overnight stopping point (at –20°C), but it is better to complete the cDNA purification (next section) before stopping.

E. cDNA Purification

Before beginning the cDNA purification:

- Remove the cDNAPure from the refrigerator and allow it to equilibrate to room temp for 30 min before use.
- Preheat the bottle of Nuclease-free Water to 50–60°C for at least 10 min.
- Shake the container of cDNAPure to fully resuspend the magnetic beads before use. cDNAPure is a slurry of DNA binding magnetic beads in binding buffer. The magnetic beads will settle over time.

1. Add 180 µL cDNAPure to each sample

Add 180 µL of resuspended cDNAPure to each second strand cDNA synthesis reaction.

2. Transfer the samples to a U-Bottom Plate, and shake gently for 2 min

- Transfer the samples from the PCR Plate to a U-Bottom Plate.



NOTE

This transfer step facilitates good mixing of the cDNAPure with the cDNA.

- Gently shake for at least 2 min to thoroughly mix the sample with the cDNAPure. The cDNA in the samples will bind to the magnetic beads in the cDNAPure during this incubation.

3. Capture the magnetic beads and discard the supernatant

- Move the U-Bottom Plate to a magnetic stand to capture the magnetic beads. Leave the plate on the magnetic stand until the mixture becomes transparent, indicating that capture is complete. This may take 5–8 min with the Ambion 96-well Magnetic-Ring Stand (Cat #AM10050).
- Carefully aspirate the supernatant without disturbing the RNA Binding Beads. Discard the supernatant.
- Remove the plate from the magnetic stand.
For effective washing, it is critical to remove the plate from the magnetic stand before adding cDNA Wash Buffer (next step).

4. Wash twice with 150 μ L cDNA Wash Buffer

Make sure that ethanol has been added to the bottle of cDNA Wash Buffer Concentrate before using it.

- a. Add 150 μ L cDNA Wash Buffer to each sample and shake at moderate speed for 1 min.

**NOTE**

The beads will not resuspend at this point because of the low surface tension of the cDNA Wash Buffer.

- b. Move the plate to a magnetic stand, and capture the magnetic beads until the mixture becomes transparent. This takes 1–2 min with an Ambion magnetic stand.
- c. Carefully aspirate and discard the supernatant without disturbing the magnetic beads, and remove the plate from the magnetic stand.
- d. Repeat steps [a–c](#) to wash a second time with 150 μ L of cDNA Wash Buffer.
- e. Move the plate to a shaker and shake the plate vigorously for 2 min to evaporate residual ethanol from the plate.
Do not overdry the beads as this may make them difficult to resuspend in the next step.

5. Elute cDNA with 18 μ L preheated Nuclease-free Water

- a. Elute the cDNA from the magnetic beads by adding 18 μ L preheated (50–60°C) Nuclease-free Water to each sample.
- b. Vigorously shake the plate for 3 min, then check to make sure the magnetic beads are fully dispersed. If they are not, continue shaking until the beads are dispersed.
- c. Move the plate to the magnetic stand, and capture the magnetic beads until the mixture becomes transparent.
- d. Carefully transfer 16 μ L of the eluted cDNA into a new PCR Plate.

**NOTE**

The purified cDNA can be stored overnight at –20°C, if desired.

F. In Vitro Transcription to Synthesize aRNA

1. Add 24 μL *IVT Master Mix* to each sample

- a. Assemble the *IVT Master Mix* at room temp in a nuclease-free tube. Use the reaction components for the type of aRNA that you want to synthesize. Prepare enough master mix for all samples in the experiment plus ~10% overage to allow for pipetting error. We provide a master mix calculator on our website to calculate reagent amounts: www.ambion.com/tools/ma296

Assemble the IVT Master Mix at room temp in the order shown:

IVT Master Mix (for a single 40 μL reaction)*		
Biotin labeled	Unmodified	Component
4 μL	4 μL	T7 ATP Soln
4 μL	4 μL	T7 CTP Soln
4 μL	4 μL	T7 GTP Soln
2.6 μL	4 μL	T7 UTP Soln
1.4 μL	--	Biotin-11-UTP†, 75 mM
4 μL	4 μL	10X T7 Reaction Buffer
4 μL	4 μL	T7 Enzyme Mix

* These are our current recommendations for preparing biotin labeled aRNA, however, for continuity with previous experiments, our former instructions are provided in section [V.A](#) starting on page 25.

† Biotin-16-UTP can be used instead of Biotin-11-UTP, if needed.

- b. Mix well by gently vortexing. Centrifuge ~5 sec to collect the master mix at the bottom of the tube and place it on ice.
- c. Transfer 24 μL of IVT Master Mix to each sample; mix by pipetting up and down twice. Cover the reactions with aluminum sealing foil.

2. Incubate the reaction for 4–14 hr at 37°C

The minimum recommended incubation time is 4 hr. The maximum incubation time is 14 hr. It is important to maintain a constant 37°C incubation temperature. We recommend incubating in a thermal cycler with a heated lid (lid temp set at 42°C). An alternative is to use a mechanical convection oven, such as a hybridization oven, with the PCR Plate in an aluminum heat block (preferred), or on the oven shelf (rather than on the oven floor). It is extremely important that condensation does not form inside the tubes; this would change the concentration of the reaction components and reduce yield.

G. aRNA Purification

Before beginning the aRNA purification:

- Preheat the aRNA Elution Buffer to 50–60°C for at least 10 min.

1. Add 70 μ L aRNA Binding Mix and 110 μ L ethanol or 50 μ L isopropanol

Biotin-labeled aRNA requires slightly different conditions for optimal interaction with the RNA Binding Beads than unmodified aRNA or aRNA synthesized with fluorescent dye- or amino allyl-nucleotides. Follow the instructions below appropriate for the type of aRNA to be synthesized.

- Add 70 μ L aRNA Binding Mix (prepared as described in section [II.B.3](#) on page 11) to each sample.
- Biotin-labeled aRNA*: Add 110 μ L of 100% ethanol to each sample. *Unmodified aRNA, amino allyl aRNA, and CyDye-modified aRNA*: Add 50 μ L of 100% isopropanol to each sample.

2. Transfer sample to U-bottom plate and shake gently for 2 min

Transfer sample into a U-Bottom Plate. Gently shake for ≥ 2 min to thoroughly mix. The aRNA in the samples will bind to the RNA Binding Beads during this incubation.

3. Capture the RNA Binding Beads and discard supernatant

- Move the plate to a magnetic stand and capture the magnetic beads, until the mixture becomes transparent.
- Carefully aspirate and discard the supernatant without disturbing the magnetic beads; then remove the plate from the magnetic stand.

4. Wash twice with 100 μ L aRNA Wash Solution

- Add 100 μ L aRNA Wash Solution (prepared as described in step [II.B.2](#) on page 10) to each sample, and shake at moderate speed for 1 min.



NOTE

The RNA Binding Beads may not fully disperse during this step; this is expected and will not affect RNA purity or yield.

- Move the plate to a magnetic stand and capture the RNA Binding Beads until the mixture becomes transparent.
- Carefully aspirate and discard the supernatant without disturbing the RNA Binding Beads and remove the plate from the magnetic stand.
- Repeat steps [a–c](#) to wash a second time with 100 μ L of aRNA Wash Solution.
- Move the plate to a shaker and shake the plate vigorously for 1 min to evaporate residual ethanol from the beads.

5. Elute aRNA with 40 μ L aRNA Elution Buffer

- Elute the purified aRNA from the RNA Binding Beads by adding 40 μ L preheated (50–60°C) aRNA Elution Buffer to each sample.
- Vigorously shake the plate for 3 min, then check to make sure the RNA Binding Beads are fully dispersed. If they are not, continue shaking until the beads are dispersed.
- Move the plate to a magnetic stand, and capture the RNA Binding Beads until the mixture becomes transparent.

- d. Transfer the supernatant, which contains the eluted aRNA, to a nuclease-free PCR plate for storage.
- e. Store purified aRNA at -20°C overnight or at -80°C for long-term storage.

For long-term storage, keep the aRNA at -80°C and minimize repeated freeze-thawing. Splitting samples into 5–20 μg aliquots is a good way to prevent multiple freeze-thaws.

III. aRNA Evaluation and Preparation for Array Hybridization

A. aRNA Quantitation

1. Assessing aRNA yield by UV absorbance

The concentration of an aRNA solution can be determined by measuring its absorbance at 260 nm. We recommend using NanoDrop Spectrophotometers for convenience. No dilutions or cuvettes are needed; just measure 2 μL of the aRNA sample directly.

Alternatively, the aRNA concentration can be determined by diluting an aliquot of the preparation in TE (10 mM Tris-HCl pH 8, 1 mM EDTA) and reading the absorbance in a traditional spectrophotometer at 260 nm. Find the concentration in $\mu\text{g}/\text{mL}$ using the equation shown below. ($1 A_{260} = 40 \mu\text{g RNA}/\text{mL}$)

$$A_{260} \times \text{dilution factor} \times 40 = \mu\text{g RNA}/\text{mL}$$

2. Assessing aRNA yield with RiboGreen

If a fluorometer or a fluorescence microplate reader is available, the RiboGreen fluorescence-based assay for RNA quantitation (Invitrogen) is a convenient and sensitive way to measure RNA concentration. Follow the manufacturer's instructions for using RiboGreen.

B. Analysis of aRNA Size with a Bioanalyzer

The Agilent 2100 bioanalyzer with Caliper's RNA LabChip Kits provides better qualitative data than conventional gel analysis for characterizing aRNA. When used with the Ambion RNA 6000 Ladder (Cat #AM7152), this system can provide a fast and accurate size distribution profile of aRNA samples. Follow the manufacturer's instructions for performing the assay.

Expected results

Using 1 μg of high quality input total RNA, the expected aRNA profile is a distribution of sizes from 250–5500 nt with a peak centered at 1000–1500 nt (see Figure 3 on page 21). Reaction products may be smaller when less input RNA or lower quality input RNA is used.

C. Preparing Labeled aRNA for Hybridization

1. Concentrate the labeled aRNA if necessary

Typically, microarrays are hybridized with 25–100 μL hybridization solution. Since the labeled aRNA is eluted in about 40 μL , it may need to be concentrated to use it for microarray hybridization. We recommend vacuum drying the labeled aRNA in the dark (if appropriate) until the volume is reduced to 1–10 μL (to keep out ambient light cover the lid of the vacuum drier with aluminum foil, if necessary).

Bring the volume of the aRNA to 10 μ L with either hybridization buffer or Nuclease-free Water and calculate the aRNA concentration based on the amount of RNA coupled to dye or the A_{260} reading.

2. Fragment labeled RNA for hybridization to oligonucleotide microarrays

For microarrays printed with oligonucleotides, the labeled aRNA must typically be fragmented for proper hybridization. Any RNA hydrolysis method that produces RNA fragments in the 60–200 nucleotide size range can be used. We recommend Ambion RNA Fragmentation Reagents (Cat #AM8740) for this procedure. Follow the procedure associated with fragmentation method used.

3. Dilute the labeled aRNA into the hybridization solution

aRNA amount

The amount of aRNA to use for hybridization will depend on your microarray type and will have to be optimized for maximum sensitivity and minimal background.

Most arrays require 1–10 μ g of labeled aRNA.

Choice of hybridization solution

The choice of hybridization solution should be based on the type of microarray being hybridized. For glass arrays, we recommend Ambion SlideHyb™ Glass Array Hybridization Buffers (Cat #AM8861–AM8863).

IV. Troubleshooting

A. Positive Control Reaction

Control RNA amplification instructions

To establish if the reaction conditions are optimal, a tube of Control RNA is provided. *Use 1 μ L of the Control RNA* in a MessageAmp II-96 reaction; follow the protocol starting at section [II.C](#) step [1](#) on page 11. At step [II.F.2](#) on page 16, *use a 14 hr incubation for the IVT reaction*. Continue with the procedure for making unmodified aRNA through section [II.G](#) on page 16.

Analysis of the positive control amplification reaction

Quantitative analysis

After completing the aRNA purification, measure the A_{260} of the reaction product as described in section [III.A.1](#) on page 19; *the positive control reaction should produce $\geq 90 \mu$ g of aRNA*.

Qualitative analysis

Also, run 100–200 ng of the reaction products on a bioanalyzer as described in section [III.B](#) on page 19; *the expected aRNA profile is a distribution from 250–5500 nt with a peak centered at 1000–1500 nt* (Figure [3](#)).

Alternatively, if a bioanalyzer is not available, run 1 μ g of aRNA on a denaturing agarose gel.

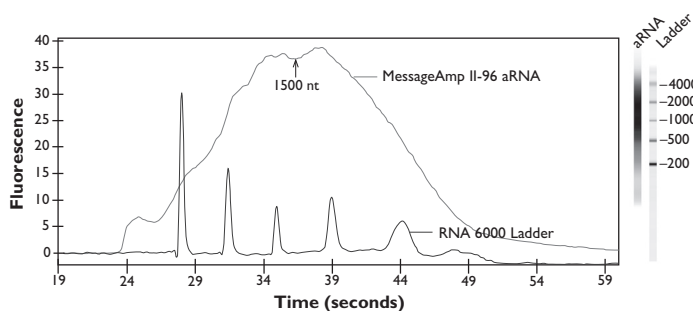


Figure 3. Bioanalyzer Electropherogram of MessageAmp II-96 aRNA.

A sample of the aRNA produced from the MessageAmp II-96 positive control reaction was analyzed on an Agilent 2100 Bioanalyzer.

B. Factors that Affect Both the Positive Control and Experimental Samples

If the positive control reaction yield or amplification product size does not meet expectations, consider the following possible causes and troubleshooting suggestions. These suggestions also apply to problems with amplification of experimental RNA.

Incubation temperature(s) were incorrect

The incubation temperatures are critical for effective RNA amplification.

- Check the temperatures of all incubators used in the procedure with a calibrated thermometer.
- If a thermal cycler is used for incubation, check the accuracy of the adjustable temperature lid. If the lid temperature cannot be adjusted to match the desired reaction temperature, use the lid with the heat turned off, or do not use it to cover the reaction vessel(s).

Condensation formed in the tube during the reaction incubation(s)

Condensation occurs when the cap of the reaction vessel is cooler (e.g. room temperature) than the bottom of the tube. As little as 1–2 μL of condensate in an IVT reaction changes the concentrations of the nucleotides and magnesium which can reduce yields.

The best way to avoid condensation is to use a thermal cycler for all incubations and to set the lid temp 5°C above the reaction temp.

If you see condensation, spin the plate briefly and mix the reaction gently. Move the plate to an incubator where condensation does not occur or is minimized.

Nuclease-contaminated tubes, tips, or equipment

Using pipettes, tubes, or equipment that are contaminated with nucleases can cleave the RNA or DNA being generated at each step in the procedure. This will reduce the size of the aRNA products and decrease aRNA yield. Both RNases and DNases can be removed from surfaces using Ambion RNaseZap® RNase Decontamination Solution (Cat #AM9780).

Absorbance readings were inaccurate

Confirm that your spectrophotometer is accurate by measuring the absorbance of an RNA or DNA sample of known concentration. Alternatively, assess the aRNA concentration by fractionating on an agarose gel adjacent to an RNA sample whose concentration is known. Comparing the ethidium bromide staining of the aRNA and control samples can approximate the concentration of the aRNA.

Incomplete resuspension or dispersion of magnetic beads

In general, the cDNAPure and the RNA Binding Beads will disperse more easily when the temperature of the mixture is warmer than 20°C.

- a. Be sure to resuspend RNA Binding Beads just before each step in preparing the aRNA Binding Mix (section [II.B.3](#) on page 11).
- b. Make sure that magnetic bead mixtures are fully resuspended before adding them to cDNA or aRNA samples for purification procedures (steps [II.E.1](#) on page 14, and [II.G.2](#) on page 17).
- c. For efficient elution of purified reaction products, make sure that magnetic beads are fully resuspended in the elution solution (steps [II.E.5](#) on page 15 and [II.G.5](#) on page 17).

- d. Avoid overdrying magnetic beads before eluting the RNA (in steps [II.E.4.e](#) on page 15 and [II.G.4.e](#) on page 17) because this may make the beads more difficult to resuspend. If the beads are inadvertently overdried, increase the mixing time (to 10 min) during the elution step to allow the beads to rehydrate.

Loss of the cDNAPure and/or the RNA Binding Beads

Since the reaction products are purified by immobilization on the cDNAPure and the RNA Binding Beads, any magnetic bead loss during the cDNA or aRNA purifications will result in lower final aRNA yield. Magnetic beads can be lost by incomplete magnetic bead capture and/or by inadvertently aspirating beads when removing supernatant from captured beads.

To determine whether cDNAPure or RNA Binding Beads have been aspirated along with supernatant, it may be helpful to collect all supernatants (except the final cDNA and aRNA-containing supernatant) in a single container. Observe the color of the collected supernatant, if magnetic beads are present, the solution will be light brown.

To prevent aspiration of magnetic beads in subsequent experiments, reduce the aspiration speed and make sure that the tip openings are not touching the bottom of the well when removing supernatant from captured cDNAPure and RNA Binding Beads.

To avoid incomplete magnetic bead capture, make sure that the solution becomes transparent before removing supernatant. This is especially important during the cDNA Purification in section [II.E](#), because the reaction mixture may be somewhat viscous.

C. Troubleshooting Low Yield and Small Average aRNA Size

Consider the following troubleshooting suggestions if the positive control reaction produced the expected results, but amplification of your experimental samples results in less aRNA than expected or average aRNA size below approximately 500 nt.

Impure RNA samples

RNA samples with significant amounts of contaminating DNA, protein, phenol, ethanol, or salts are reverse transcribed poorly and subsequently generate less aRNA than pure RNA samples. Phenol extract and ethanol precipitate your RNA, or use the Ambion MEGAclean™ Kit (Cat #AM1908) to further purify the samples before reverse transcription.

Lower than expected input RNA concentration

Take another A_{260} reading of your RNA sample or try using more RNA in the aRNA amplification procedure.

RNA sample integrity is poor

RNA that is partially degraded generates cDNA that is relatively short. This will reduce the average size of the aRNA population and subsequently reduce the yield of aRNA. You can assess the integrity of an

The mRNA content of your total RNA sample is lower than expected

RNA sample by determining the size of the 18S and 28S rRNA bands and the relative abundance of 28S to 18S rRNA (See section [II.A. RNA integrity](#) on page 8 for more information).

Different RNA samples contain different amounts of mRNA. In healthy cells, mRNA constitutes 1–3% of total cellular RNA. The actual amount of mRNA depends on the cell type and the physiological state of the sample. When calculating the amount of amplification, the starting mass of mRNA in a total RNA prep should always be considered a range from 10–30 ng per µg of total RNA (assuming good RNA quality). Most total RNA samples can be amplified up to 1000 fold producing 10–30 µg of aRNA from 1 µg of total RNA.

D. Bead Carryover in Eluted Sample

If cDNAPure or RNA Binding Beads are carried over into the eluate containing the RNA, they will cause the solution to be light brown in color.

To avoid bead carryover, allow sufficient time for bead capture and aspirate slowly and carefully when transferring eluent from the beads. Also avoid touching the bottom of the well with the pipette tip when removing solutions from captured beads.

V. Appendix

A. IVT Reaction Using 10 mM Biotin-labeled CTP and UTP

This is our former protocol for preparing biotin-labeled aRNA using the MessageAmp II-96 aRNA Amplification Kit. Our current recommendations are to use the procedure outlined in the main protocol (section [II.F. In Vitro Transcription to Synthesize aRNA](#) starting on page 16).

We include this protocol for the convenience of customers who choose to continue using our former recommendations for consistency with previous experiments.

Following is a list of biotin-labeled 10 mM CTP and UTP products recommended for this protocol.

- Biotin-11-CTP (10 mM): PerkinElmer Life Sciences is the preferred supplier of biotin-11-CTP (Cat #NEL542001EA), other biotin-labeled NTPs that may be used with this kit are available upon request. Alternatively, biotin-11-CTP is available from Enzo Biochemicals (Cat #42818).
- Biotin-16-UTP (10 mM) is available from Ambion Inc. (Cat #8452).

1. Protocol modifications

Follow the protocol detailed in section [II. aRNA Amplification Protocol](#) starting on page 8, but substitute these instructions for step [1](#) in section [II.F](#) starting on page 16.

1a. Biotin-labeled rxns: Mix biotin NTPs with the cDNA and concentrate to 16 μ L

For each sample, add the following to the eluted cDNA (16 μ L):

Amount	Component
7.5 μ L	10 mM biotin-11-CTP
7.5 μ L	10 mM biotin-16-UTP

Concentrate the mixture in a vacuum centrifuge concentrator until the volume is reduced to 16 μ L. This should take only a few minutes. Do not dry the mixture to completion. If volume is lower than 16 μ L, bring the mixture to 16 μ L with Nuclease-free Water.

1b. Unlabeled rxns: Proceed directly to step [2](#)

For synthesis of unmodified aRNA, proceed directly to step [2](#).

2. Add 24 μ L IVT Master Mix to each sample

- Assemble the IVT Master Mix at room temp in a nuclease-free tube. Use the reaction components for the type of aRNA that you want to synthesize. Prepare enough master mix for all samples in the experiment plus ~10% overage to allow for pipetting error.

Assemble the IVT Master Mix at room temp in the order shown:

IVT Master Mix (for a single 40 μ L reaction)		
Biotin labeled	Unmodified	Component
4 μ L	4 μ L	T7 ATP Soln
3 μ L	4 μ L	T7 CTP Soln
4 μ L	4 μ L	T7 GTP Soln
3 μ L	4 μ L	T7 UTP Soln
4 μ L	4 μ L	10X T7 Reaction Buffer
4 μ L	4 μ L	T7 Enzyme Mix
2 μ L	–	Nuclease-free Water

- b. Mix well by gently vortexing. Centrifuge \sim 5 sec to collect the master mix at the bottom of the tube and place it on ice.
- c. Transfer 24 μ L of *IVT Master Mix* to each sample; mix by pipetting up and down twice. Cover the reactions with aluminum sealing foil.
- d. Proceed to step [II.F.2](#) on page 16.

B. References

- Ausubel, FM, et al (editors) *Current Protocols in Molecular Biology*, John Wiley and Sons.
- Baugh LR, Hill EL, Hunter CP (2001) Quantitative analysis of mRNA amplification by in vitro transcription. *Nucl Acids Res* **29**(5): E29.
- Feldman AJ, Costouros NG, Wang E, Qian M, Marincola FM, Alexander HR, and Libutti SK (2002) Advantages of mRNA amplification for microarray analysis. *Biotechniques* **33**(4): 906–914.
- Li Y, Li T, Liu S, Qiu M, Han Z, Jiang Z, Li R, Ying K, Xie Y, Mao Y (2004) Systematic comparison of the fidelity of aRNA, mRNA and T-RNA on gene expression profiling using cDNA microarray. *J Biotechnology* **107**(1):19–28.
- Kacharina JE, Crino PB, Eberwine J (1999) Preparation of cDNA from single cells and subcellular regions. *Methods Enzymol*, **303**: 3–18.
- Pabon C, Modrusan Z, Ruvolo MV, Coleman IM, Daniel S, Yue H, Arnold LJ Jr. (2001) Optimized T7 amplification system for microarray analysis. *Biotechniques* **31**(4): 874–879.
- Polacek DC, Passerini AG, Shi C, Francesco NM, Manduchi E, Grant GR, Powell S, Bischof H, Winkler H, Stoeckert CJ Jr, Davies PF (2003) Fidelity of enhanced sensitivity of differential transcription profiles following linear amplification of nanogram amounts of endothelial mRNA. *Physiol Genomics* **13**: 147–156.
- Poirier G M-C, Erlander MG (1998) Postdifferential Display: Parallel Processing of Candidates Using Small Amounts of RNA. *Methods: A Companion to Method in Enzymology* **16**: 444–452.
- Van Gelder RN, von Xastrow ME, Yool A, Dement DC, Barchas JD, Eberwine JH (1990) Amplified RNA synthesized from limited quantities of heterogeneous cDNA. *Proc Natl Acad Sci USA*, **87**: 1663–1667.

Yue H, Eastman PS, Wang BB, Minor J, Doctolero MH, Nuttall RL, Stack R, Becker JW, Montgomery JR, Vainer M, Johnston R (2001) An evaluation of the performance of cDNA microarrays for detecting changes in global mRNA expression. *Nucleic Acids Res* 29(8):E41–1.

C. MessageAmp II-96 Kit Contents

cDNA synthesis and in vitro transcription reagents

Store at -20°C in a non-frost-free freezer.

Amount	Component
110 μL	T7 Oligo(dT) Primer*
110 μL	ArrayScript™
110 μL	RNase Inhibitor
220 μL	10X First Strand Buffer
880 μL	dNTP Mix
220 μL	DNA Polymerase
110 μL	RNase H
10 μL	Control RNA (1 $\mu\text{g}/\mu\text{L}$ HeLa total RNA)
1.1 mL	10X Second Strand Buffer
440 μL	T7 Enzyme Mix
440 μL	T7 10X Reaction Buffer
440 μL	T7 ATP Solution (75 mM)
440 μL	T7 CTP Solution (75 mM)
440 μL	T7 GTP Solution (75 mM)
440 μL	T7 UTP Solution (75 mM)

* T7 Oligo(dT) Primer is also available separately from Ambion (Cat #AM5710).

cDNA and aRNA purification components

Amount	Component	Storage
20 mL	Nuclease-free Water	any temp*
1.1 mL	RNA Binding Beads	$4^{\circ}\text{C}\dagger$
20 mL	cDNAPure‡	$4^{\circ}\text{C}\ddagger$
40 mL	cDNA Wash Buffer Concentrate Add 32 mL 100% ethanol before use	room temp
25 mL	aRNA Wash Solution Concentrate Add 20 mL 100% ethanol before use	room temp
450 μL	Bead Resuspension Solution	room temp
10 mL	aRNA Elution Buffer	room temp
5.5 mL	aRNA Binding Buffer Concentrate	room temp
2	U-Bottom Plate	room temp
4	PCR Plate	room temp
1	Reservoir	room temp

* Store Nuclease-free Water at room temp, 4°C , or -20°C .

† Do not freeze.

‡ cDNAPure is manufactured by Agencourt Bioscience Corporation.

To obtain Material Safety Data Sheets

- Material Safety Data Sheets (MSDSs) can be printed or downloaded from product-specific links on our website at the following address: www.ambion.com/techlib/msds
- Alternatively, e-mail your request to MSDS_Inquiry_CCRM@appliedbiosystems.com. Specify the catalog or part number(s) of the product(s), and we will e-mail the associated MSDSs unless you specify a preference for fax delivery.
- For customers without access to the internet or fax, our technical service department can fulfill MSDS requests placed by telephone or postal mail. (Requests for postal delivery require 1–2 weeks for processing.)

D. Quality Control**Functional testing**

The Control RNA is used in a MessageAmp II reaction following the instructions in section [IV.A](#) on page 21. The aRNA yield is assessed by measuring the A_{260} on the Nanodrop ND1000A spectrophotometer. The median size of the aRNA is assessed using the mRNA smear assay on the Agilent 2100 bioanalyzer.

Nuclease testing

Relevant kit components are tested in the following nuclease assays:

RNase activity

Meets or exceeds specification when a sample is incubated with 25 ng labeled RNA and analyzed by PAGE.

Nonspecific endonuclease activity

Meets or exceeds specification when a sample is incubated with 300 ng supercoiled plasmid DNA and analyzed by agarose gel electrophoresis.

Exonuclease activity

Meets or exceeds specification when a sample is incubated with 40 ng labeled *Sau3A* fragments of pUC19 and analyzed by PAGE.

Protease testing

Meets or exceeds specification when a sample is incubated with 1 μ g protease substrate and analyzed by fluorescence.