High Throughput RNA Amplification and Labeling for Array Analysis

Part Number AM1821



#### (Part Number AM1821)

#### Protocol

I.	Introduction
	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 Invitrogen
II.	aRNA Amplification Procedure
	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.	Assessing aRNA Yield and Size
	A. aRNA Quantitation
	B. Analysis of aRNA Size with a Bioanalyzer
IV.	Dye Coupling and Cleanup
	A. Concentrate the Amino Allyl aRNA
	B. aRNA:Dye Coupling Reaction
	C. Dye Labeled aRNA Purification
	D. Spectrophotometric Analysis of Dye Incorporation
	E. Preparing Labeled aRNA for Hybridization
V.	Troubleshooting
	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
	E. Troubleshooting Dye Coupling
VI.	Appendix
	A. References
	B. Quality Control
	C. Safety Information

#### P/N 1821M Revision G

For research use only. Not intended for animal or human therapeutic or diagnostic. 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.

Revision Date: September 6, 2010

Information in this document is subject to change without notice. Ambion, a part of Life Technologies Corporation, assumes no responsibility for any errors that may appear in this document.

Ambion 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 Ambion 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, please refer to it as the Amino Allyl MessageAmp $^{TM}$  II-96 Kit.

**Warranty and Liability:** Ambion, a part of Life Technologies Corporation, 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 Ambion if you have any questions about our warranties or would like information about post-warranty support.

Patents and Licensing Notifications: The Amino Allyl 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 Life Technologies 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. DIS-**CLAIMER OF WARRANTIES. GE HEALTHCARE BIO-SCIENCES CORP. PROVIDES NO WARRANTIES** TO END USER (STATUTORY OR IMPLIED), INCLUDING WITHOUT LIMITATION, AS TO PROD-UCT QUALITY, CONDITION, DESCRIPTION, MERCHANTABILITY OR FITNESS FOR A PARTICU-LAR PURPOSE, AND ALL SUCH WARRANTIES ARE HEREBY EXPRESSLY DISCLAIMED. HEALTHCARE BIO-SCIENCES CORP. HEREBY EXPRESSLY DISCLAIMS ANY WARRANTY REGARD-ING RESULTS OBTAINED THROUGH THE USE OF THE PRODUCTS, INCLUDING WITHOUT LIMI-TATION 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:** Trademarks of Life Technologies Corporation and its affiliated companies: Applied Biosystems®, AB (Design)®, AlexaFluor®, Ambion®, ArrayControl™, ArrayScript™, MagMAX™, MEGAclear™, MEGAscript®, MessageAmp™, RiboGreen®, RNAqueous®, RNase*Zap*®, and SlideHyb™. Cy and CyDye are trademarks of GE Healthcare. TRI Reagent is a registered trademark of Molecular Research Center, Inc. All other trademarks are the sole property of their respective owners.

#### I. Introduction

#### A. Product Description and Background

The Amino Allyl MessageAmp™ II-96 Kit (patent pending) 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 2 on page 4; Van Gelder et al. 1990). It consists of reverse transcription of RNA using an oligo(dT) primer bearing a T7 promoter and ArrayScript<sup>™</sup> reverse transcriptase (RT), which is engineered to produce higher yields of first strand cDNA than wild type enzymes. ArrayScript RT 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 cleanup to become a template for in vitro transcription with T7 RNA Polymerase. To maximize aRNA yield, Ambion proprietary MEGAscript<sup>®</sup> in vitro transcription (IVT) technology is used in the kit to generate hundreds to thousands of antisense RNA copies of each mRNA in a sample. (In this protocol the antisense amplified RNA is referred to as aRNA, in scientific literature it is also commonly called cRNA.) The IVT is configured to incorporate the modified nucleotide, 5-(3-aminoallyl)-UTP (aaUTP) into the aRNA during in vitro transcription, aaUTP contains a reactive primary amino group on the C5 position of uracil that can be chemically coupled to N-hydroxysuccinimidyl ester-derivatized reactive dyes (NHS ester dyes) in a simple, efficient reaction (see Figure 1). Once purified, the labeled 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 RNA samples of very limited amounts 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).

Figure 1. Amino Allyl Labeling Reaction

The Amino Allyl MessageAmp II-96 advantage

#### Optimized protocols

All reagents and protocols have been optimized to produce high quality amino allyl-modified aRNA with maximum yield and reproducibility. The simple, rapid aRNA purification procedure prepares the aRNA for post-synthesis labeling reactions.

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

#### **Built-in flexibility**

The NTPs are provided in three separate tubes: a mixture of ATP, CTP and GTP, a UTP Solution, and an aaUTP Solution. Separating the UTP and the aaUTP provides the flexibility to make unlabeled aRNA or to incorporate aaUTP during the amplification. Unlike most modified nucleotides, the inclusion of aaUTP in an IVT reaction has only a minor effect on the reaction efficiency and yield. Additionally, since the incorporation of aaUTP by IVT will be virtually identical in different samples and since the dye coupling reaction is efficient and reproducible, labeled samples will not have the biases that can result from direct incorporation of modified nucleotides.

#### 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:

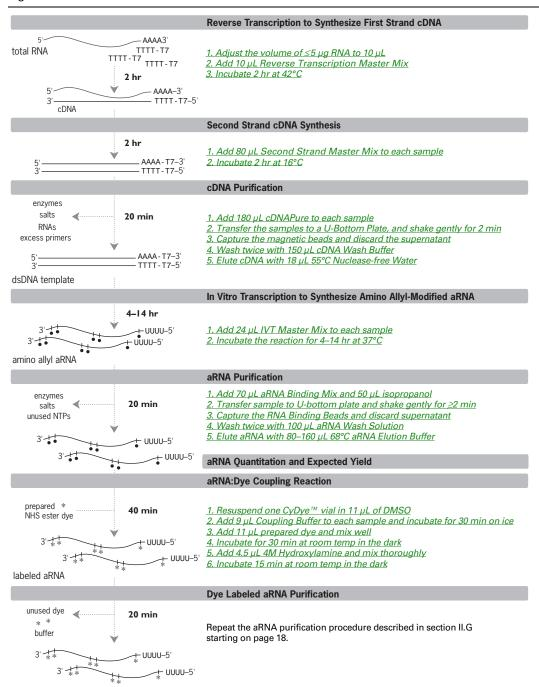
www4.appliedbiosystems.com/techlib/automation

#### 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 (P/N AM1751).

- 1. For the first round amplification, use the MessageAmp II Kit to make unmodified aRNA. (You will need more than one kit to process more than 20 samples.)
- For the second round of amplification, use the Amino Allyl MessageAmp II-96 Kit components plus the Second Round Primers from the MessageAmp II Kit, and follow instructions from both protocols as detailed below:
  - For second round reverse transcription and second strand cDNA synthesis, follow the instructions in the MessageAmp II protocol. (You will need Second Round Primers from more than one MessageAmp II Kit to process more than 20 samples.)
  - For second round cDNA purification, follow either the MessageAmp II Kit or the Amino Allyl MessageAmp II-96 protocol, depending on your preference and available kit components.
  - For the second round amplification IVT reaction, follow the instructions in the Amino Allyl MessageAmp II-96 Kit protocol.

Figure 2. Procedure Overview



#### B. Materials Provided with the Kit and Storage

The Amino Allyl MessageAmp II-96 Kit includes reagents for amplification and labeling 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™ RT
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
220 µL	T7 UTP Solution (75 mM)
220 µL	aaUTP Solution (75 mM) [5-(3-amino allyl)-UTP]
1.32 mL	ATP, CTP, GTP Mix

## cDNA and aRNA purification components

Amount	Component	Storage
50 mL	Nuclease-free Water	any temp*
20 mL	cDNAPure†	4°C ±
2.2 mL	RNA Binding Beads	4°C‡
40 mL	cDNA Wash Buffer Concentrate Add 32 mL 100% ethanol before use	room temp
50 mL	aRNA Wash Solution Concentrate Add 40 mL 100% ethanol before use	room temp
900 μL	Bead Resuspension Solution	room temp
22 mL	aRNA Elution Buffer	room temp
11 mL	aRNA Binding Buffer Concentrate	room temp
3	U-Bottom Plate	room temp
5	PCR Plate	room temp
1	Reservoir	room temp

<sup>\*</sup> Store Nuclease-free Water at room temperature, 4°C, or -20°C.

<sup>†</sup> cDNAPure is manufactured by Agencourt Bioscience Corporation.

<sup>‡</sup> Do not freeze.

#### Dye labeling reagents

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

Amount	Component
1 mL	4M Hydroxylamine
2.6 mL	DMSO
2 mL	Coupling Buffer

#### C. Materials Not Provided with the Kit

#### Lab equipment and supplies

- ACS grade (or higher quality) alcohol:
  - -100% ethanol
  - -100% isopropanol
- Additional 100% DMSO may be needed depending on how the dye is supplied.
- Thermal cycler with adjustable-temperature heated lid (recommended), hybridization oven, or fixed temperature air incubators set at 70°C, 42°C, 37°C, and 16°C. (See "Procedural notes" on page 10 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, P/N AM10050, for its high-strength magnets and quality design)
- Microcentrifuge
- Centrifugal vacuum concentrator with plate adaptor
- RNase-free pipettors and tips, positive-displacement type recommended to increase the accuracy and precision of reaction inputs
- 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)

## NHS esters of fluorescent dyes

This kit was developed using CyDye™ fluorescent dyes from GE Healthcare (Amersham Biosciences), but mono-reactive NHS esters of any label moiety should be capable of coupling to the amino allyl modified aRNA generated with this kit. Some suppliers of compatible dyes are listed below:

- Amersham Biosciences CyDye fluors
- Molecular Probes Alexa Fluor® Succinimidyl Esters
- Pierce Biotechnology NHS-Fluorescein and -Rhodamine
- Denovo Biolabels GmbH Oyster® Dyes

#### D. Related Products Available from Invitrogen

96-well Magnetic-Ring Stand P/N AM10050	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.
Magnetic Stand-96 P/N AM10027	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 (P/N AM10050).
MagMAX™-96 Total RNA Isolation Kit P/N AM1830	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.
MagMAX™-96 for Microarrays Total RNA Isolation Kit P/N AM1839	The MagMAX-96 for Microarrays Total RNA Isolation Kit (patent pending) 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.
RNAqueous®-96 Automated Kit P/N AM1812	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.

ArrayControl™ Spikes P/N AM1780	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.
RNA Fragmentation Reagents P/N 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 web or print catalog for P/Ns	There are 3 unique SlideHyb Glass Array Hybridization Buffers; they have identical salt and formamide compositions, but differ in hybridization kinetics and blocking reagents. The Glass Array Hybridization Cassette is also available, for incubation of glass microarray hybridization reactions.

#### II. aRNA Amplification Procedure

#### Α. 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 Protocol, please contact our high throughput experts by emailing:

automation@ambion.com

#### RNA quantity and IVT incubation time

The Amino Allyl 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 ug of total RNA should be used in an Amino Allyl 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 RNA samples of limited amounts (e.g. -250 ng or less), however, incubating the IVT reaction for 14 hr will maximize the amount of aRNA produced.



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

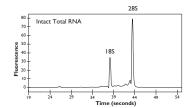
#### RNA purity

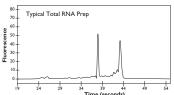
RNA samples should be free of contaminating proteins, DNA, and other cellular material as well as phenol, ethanol, and salts associated with RNA isolation procedures. Impurities can lower the efficiency of reverse transcription and subsequently reduce the level of amplification. An effective measure 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 1.7-2.1.

#### **RNA** integrity

The integrity of the RNA sample, or the proportion of full-length mRNA, can have a significant impact on 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 3).





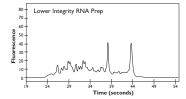


Figure 3. Bioanalyzer Images of Increasingly Degraded Total RNA

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**

#### Thermal cycler recommended

The Amino Allyl MessageAmp II-96 procedure is very sensitive to temperature; variable or inaccurate incubation temperatures can limit aRNA synthesis. 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, which can greatly reduce yield.

### A thermal cycler with a temperature adjustable heated lid is recommended.

A calibrated thermal cycler, with a temperature-adjustable heated lid, is recommended, for the greatest temperature control and stability during Amino Allyl MessageAmp II-96 reaction incubations. Allow the thermal cycler to equilibrate to the required temperature before placing the plates in the block for incubation. Follow the recommended settings for the lid temperatures. Too high a lid setting may inhibit the reaction; too low a setting may cause condensation.

#### aRNA Amplification Procedure



Even if you use a hybridization oven or incubator for most of the Amino Allyl MessageAmp II-96 reactions, a thermal cycler is strongly recommended for the 16°C second-strand synthesis reaction incubation (step II.D.2 on page 15). Turn off the heated lid if it cannot be adjusted to match the 16°C block temperature.

If your thermal cycler does not have a temperature-adjustable lid, or a thermal cycler is unavailable, calibrated hybridization ovens or incubators (at constant temperature) may also be used. Preheat incubators so that the correct temperature has stabilized before reactions are placed in the incubator.

 Heat blocks or water baths are not recommended for Amino Allyl MessageAmp II-96 reaction incubations.

#### 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 µL 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 <2 mL in volume, we recommend that you minimize the void volume by aliquoting the master mix into a PCR plate instead of using a reservoir.

A web-based interactive calculator to calculate master mix quantities is available at:

www4.appliedbiosystems.com/tools/aama296

We recommend using the web calculator to determine reagent volumes needed for all the master mixes in the procedure and printing the results to accompany this 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 40 mL 100% ethanol to the aRNA Wash Solution Concentrate

Add 40 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 temperature for one week; prepare only the amount needed for all samples in the experiment plus ~10% overage to cover pipetting error. A master mix calculator is available at:

www4.appliedbiosystems.com/tools/aama296

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

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

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

<sup>\*</sup> Mix the RNA Binding Beads by vortexing before dispensing.

#### b. Next, add aRNA Binding Buffer Concentrate

At room temperature, 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 μL	
RNA Binding Bead Mixture (from step 3.a)	20 μL	

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

#### C. Reverse Transcription to Synthesize First Strand cDNA

#### Incubator needed

Thermal cycler programmed as follows:

Temp	Time	Cycles
42°C (50°C lid)	2 hr	1
4°C	hold	

It is important to follow the lid temperature recommendation; higher lid temperatures may inhibit reverse transcription, while lower temperatures may cause condensation, resulting in changes to the reaction composition.

- 1. Adjust the volume of ≤5 µg RNA to 10 µL
- a. Place up to 5  $\mu g$  of total RNA (typically 100 ng–5  $\mu g$  ) into the wells of a PCR Plate.
- b. Add Nuclease-free Water to a final volume of 10  $\mu$ L.
- 2. Add 10 µL Reverse Transcription Master Mix
- a. Prepare *Reverse Transcription Master Mix* in a nuclease-free tube at room temperature. Assemble enough master mix for all samples in the experiment plus ~10% overage to allow for pipetting error. A master mix calculator to calculate reagent amounts is available at:

www4.appliedbiosystems.com/tools/aama296

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

Reverse Transcription Master Mix (for a single 20 $\mu L$ reaction)		
Amount	Component	
1 μL	Nuclease-free Water	
1 μL	T7 Oligo(dT) Primer	
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 10  $\mu$ L 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.
- d. Place the samples in the thermal cycler and start the run.

#### 3. 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 and immediately proceed to the second strand cDNA synthesis (next).

#### D. Second Strand cDNA Synthesis

#### Incubator needed

Thermal cycler programmed as follows:

Temp	Time	Cycles
16°C (heat-disabled lid, or no lid)	2 hr	1
4°C	hold	

Disable the heated lid of the thermocycler if the lid temperature cannot be set in the range 16°C to room temperature. If the lid is too hot, inhibition of the reaction may occur.

# 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. A master mix calculator to calculate reagent amounts is available at:

www4.appliedbiosystems.com/tools/aama296

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

Second Strand Master Mix (for a single 100 $\mu L$ 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.

#### aRNA Amplification Procedure

#### 2. Incubate 2 hr at 16°C

Place the plate in the 16°C thermal cycler, start the run, and incubate for 2 hr.

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. (Do not incubate in a heat block in a 4°C refrigerator because the temperature may fluctuate.)

3. Place reactions on ice briefly or freeze immediately After the 2 hr incubation at 16°C, place the reactions on ice and proceed to section <u>E. cDNA Purification</u> (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 temperature for 30 min before use.
- 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.
- Preheat a minimum of 20 µL per sample of Nuclease-free Water to 55°C.



#### **IMPORTANT**

Preheat the Nuclease-free Water to a maximum of 55°C; temperatures above 58°C can partially denature the cDNA, compromising final aRNA yield.

- 1. Add 180 µL cDNAPure to each sample
- Add 180  $\mu L$  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
- a. Transfer the samples from the PCR Plate to a U-Bottom Plate.



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

b. 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.

#### Capture the magnetic beads and discard the supernatant

- a. 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 (P/N AM10050).
- b. Carefully aspirate the supernatant without disturbing the RNA Binding Beads. Discard the supernatant.
- c. 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 μL cDNA Wash Buffer

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

a. Add  $150\,\mu\text{L}$  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 <u>a-c</u> to wash a second time with 150  $\mu L$  of cDNA Wash Buffer.
- 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 µL 55°C Nuclease-free Water

- a. Elute the cDNA from the magnetic beads by adding 18  $\mu$ L preheated (55°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  $\mu$ L of the supernatant, which contains the eluted cDNA, to a new PCR Plate.



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

#### F. In Vitro Transcription to Synthesize aRNA

#### Incubator needed

Thermal cycler programmed as follows:

Temp	Time	Cycles
37°C (default lid; 100–105°C)	4–14 hr; see step 2	1
4°C	hold	

It is important to use a thermal cycler with a heated lid (100–105°C), to maintain uniform temperature, and because it is extremely important that condensation does not form inside the tubes; this would change the reagent concentrations and reduce yield.

Because the amino allyl UTP (aaUTP) does not have a bulky modification group, it is possible to replace 100% of the UTP with aaUTP in the IVT reaction. We recommend, however, using a 1:1 ratio of UTP to aaUTP (i.e. 50%) for most samples. If desired, it is possible to adjust the level of aaUTP incorporation by adjusting the UTP:aaUTP ratio. Increasing the amount of aaUTP in the ratio will produce aRNA with more dye-coupling sites, which in some cases will lead to brighter signal on microarrays. This kit contains enough 75 mM aaUTP to set up 100 IVT reactions at the 1:1 ratio recommended. If more aaUTP is needed, it can be purchased from Ambion (P/N AM8437).

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

 a. Assemble the *IVT Master Mix* at room temperature in a nuclease-free tube. Prepare enough master mix for all samples in the experiment plus ~10% overage to allow for pipetting error. A master mix calculator to calculate reagent amounts is available at:

www4.appliedbiosystems.com/tools/aama296

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

IVT Master Mix (for a single 40 $\mu\text{L}$ reaction)		
Amount	Component	
2 μL	aaUTP Solution (75 mM)	
12 µL	ATP, CTP, GTP Mix	
2 μL	T7 UTP Solution (75 mM)	
4 μL	T7 10X Reaction Buffer	
4 μL	T7 Enzyme Mix	

- 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  $\mu L$  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

Place the plate in the thermal cycler and start the run.

The minimum recommended incubation time is 4 hr. The maximum incubation time is 14 hr.

#### G. aRNA Purification

#### Before beginning the aRNA purification:

Preheat the aRNA Elution Buffer to 68°C.

- Add 70 μL aRNA Binding Mix and 50 μL isopropanol
- a. Add 70  $\mu L$  aRNA Binding Mix (prepared as described in section II.B.3 on page 12) to each sample.
- b. Add 50 μL of 100% isopropanol to each sample.

2. Transfer sample to U-bottom plate and shake gently for ≥2 min Transfer the 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
- a. Move the plate to a magnetic stand and capture the magnetic beads, until the mixture becomes transparent.
- b. Carefully aspirate and discard the supernatant without disturbing the magnetic beads; then remove the plate from the magnetic stand.
- 4. Wash twice with 100 μL aRNA Wash Solution
- a. Add 100 µL aRNA Wash Solution (prepared as described in step II.B.2 on page 12) to each sample, and shake at moderate speed for 1 min.



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

- b. Move the plate to a magnetic stand and capture the RNA Binding Beads until the mixture becomes transparent.
- c. Carefully aspirate and discard the supernatant without disturbing the RNA Binding Beads and remove the plate from the magnetic stand.
- d. Repeat steps <u>a-c</u> to wash a second time with 100  $\mu L$  of aRNA Wash Solution.
- e. Move the plate to a shaker and shake the plate vigorously for 1 min to evaporate residual ethanol from the beads.
   Do not overdry the beads as this may make them difficult to resuspend in the next step.

#### aRNA Amplification Procedure

- 5. Elute aRNA with 80–160 µL 68°C aRNA Elution Buffer
- a. Elute the purified aRNA from the RNA Binding Beads by adding  $80{\text -}160~\mu\text{L}$  preheated (68°C) aRNA Elution Buffer to each sample.
- b. 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.
- c. 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.
- Store the aRNA at -20°C or determine the aRNA concentration and continue



#### STOPPING POINT

The purified aRNA can be stored at –20°C at this point.

The purified aRNA can be stored at -20°C if desired. Alternatively, determine its concentration using one of the methods described in section *III.A. aRNA Quantitation* on page 20. The aRNA concentration must be known in order to continue to the dye coupling reaction.

### III. Assessing aRNA Yield and Size

#### A. aRNA Quantitation

### 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 because no dilutions or cuvettes are needed. Follow the manufacturer's instructions.

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  $\mu$ g/mL using the equation shown below. (1  $A_{260}$  = 40  $\mu$ g RNA/mL)

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

### Assessing aRNA yield with the RiboGreen® assay

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 the RiboGreen assay.

#### 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 (P/N 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  $\mu$ g of high quality input total RNA, the expected aRNA profile for unmodified aRNA is a distribution of sizes from 250–5500 nt with a peak centered at 1000–1500 nt (see Figure 4 on page 26). Reaction products may be smaller when less input RNA or lower quality input RNA is used. When comparing unmodified aRNA with amino allyl modified aRNA, you will notice that amino allyl aRNA migrates at a slightly higher molecular weight than unmodified aRNA.

#### IV. Dye Coupling and Cleanup

This kit was developed using  $Cy^{\infty}3$  and  $Cy^{\infty}5$  dyes from Amersham Biosciences, but mono-reactive NHS esters of any label moiety should be capable of coupling to the amino allyl modified aRNA generated with this kit. For a partial list of suppliers of NHS ester dyes, see section *L.C. Materials Not Provided with the Kit* on page 6. The choice of dye will depend on your preference and on the type of microarray scanning equipment that will be used.

#### A. Concentrate the Amino Allyl aRNA

- 1. Determine the aRNA concentration
- (See section III.A on page 20 for instructions on determining the aRNA concentration.)
- 2. Vacuum dry 5–20 μg aRNA to completion
- Place 5–20 µg of amino allyl aRNA into separate wells of a PCR Plate and vacuum dry using the low heat setting until no liquid remains.
- Check the progress of drying every 5–10 min and remove the sample from the concentrator as soon as it is dry; do not overdry.

#### B. aRNA: Dye Coupling Reaction

Resuspend one CyDye™ vial in 11 µL of DMSO

Prepare dye immediately before starting the dye coupling procedure; add 11  $\mu$ L of DMSO to one tube of Amersham Biosciences Cy3 or Cy5 Post Labelling Reactive Dye, and vortex to mix thoroughly.

Keep the resuspended dye in the dark at room temperature for up to 1 hr until you are ready to use it.



#### **IMPORTANT**

The preparation and storage of solubilized dye is important for the efficient labeling of amino allyl modified aRNA. It is imperative that the dye compounds remain dry both before and after dissolving in DMSO because any water that is introduced will cause hydrolysis of the NHS esters, lowering the efficiency of coupling.

- 2. Add 9 µL Coupling Buffer to each sample and incubate for 30 min on ice
- a. To each well containing dried amino allyl aRNA, add 9  $\mu L$  Coupling Buffer.
- b. Resuspend samples by shaking vigorously for 1 min. Let the mixture equilibrate for 30 min on ice to ensure full resuspension.
- 3. Add 11 µL prepared dye and mix well

Add 11  $\mu L$  of NHS ester dye in DMSO to each well containing aRNA: Coupling Buffer mixture. Mix well.

4. Incubate for 30 min at room temp in the dark

This 30 min incubation at room temperature allows the dye coupling reaction to occur. To keep the samples in the dark, cover the plate with aluminum foil or incubate it in a closed drawer.

Add 4.5 μL
 4M Hydroxylamine and mix thoroughly

To quench the reaction, add 4.5  $\mu$ L 4M Hydroxylamine to each sample and mix thoroughly.

6. Incubate 15 min at room temp in the dark

Incubate the reaction at room temp in the dark for 15 minutes.

During this incubation, the large molar excess of hydroxylamine will quench the amine-reactive groups on the unreacted dye molecules.

#### C. Dye Labeled aRNA Purification

Repeat the aRNA purification procedure described in section <u>II.G</u> starting on page 18.

Use 80 µL preheated aRNA Elution Buffer in step II.G.5.

After the purification, if the sample will not be used the same day, store it at  $-20^{\circ}$ C in the dark.

#### D. Spectrophotometric Analysis of Dye Incorporation

The Nanodrop ND-1000A UV-Vis Spectrophotometer is recommended for UV-Vis spectrophotometry measurements because there is no need to dilute samples. Measure 1.5  $\mu$ L of undiluted sample.

1. Determine the appropriate dilution for spectrophometer readings

To use a conventional spectrophotometer, dilute labeled aRNA 1:20 by mixing 5  $\mu L$  labeled aRNA in 95  $\mu L$  TE (10 mM Tris-HCl, 1 mM EDTA). Mix well and measure the  $A_{260}$  in a UV-Vis spectrophotometer.

To obtain an accurate reading at both 260 nm and at the maximum absorbance wavelength for the dye, the aRNA dilution should result in absorbance readings (A $_{260}$  and A $_{\rm dye}$ )between 0.1 and 1.0. If absorbance readings are below 0.1, reduce the dilution factor and check the A $_{260}$  and A $_{\rm dye}$  again.

2. Measure the aRNA absorbance at 260 nm and at the absorbance max for the dye

Blank the instrument with the TE used for making dilutions. Measure the absorbance of each sample at 260 nm ( $A_{260}$ ) and also at the maximum absorbance wavelength for the dye used in the coupling reaction ( $A_{dye}$ ).

Dye type	Absorbance maximum	Extinction coefficient*
fluorescein	491 nm	66,000
Cy3	550 nm	150,000
Cy5	650 nm	250,000
Alexa Fluor 546	546 nm	104,000
Alexa Fluor 555	555 nm	150,000
Alexa Fluor 660	663 nm	132,000
Alexa Fluor 647	650 nm	239,000
Oyster 556	556 nm	155,000
Oyster 645	645 nm	250,000
Oyster 656	656 nm	220,000

<sup>\*</sup> Extinction coefficient ( $\epsilon$ ) at  $\lambda_{max}$  in cm<sup>-1</sup>M<sup>-1</sup>

- 3. Calculate the number of dye molecules incorporated per 1000 nt
- With the absorbance values ( $A_{\rm dye}$  and  $A_{\rm 260}$ ) in hand, use the convenient dye incorporation calculator on our website to calculate the number of dye molecules incorporated per 1000 nt:

www4.appliedbiosystems.com/techlib/misc/aama\_dye\_calc.html

 Alternatively, use this formula to estimate the number of dye molecules incorporated per 1000 nt of labeled aRNA.

$$\frac{\text{# dye molecules}}{1000 \text{ nt}} = \frac{A_{\text{dye}}}{A_{\text{260}}} \times \frac{X}{\text{dye extinction coefficient}} \times 1000$$

The expected incorporation rate is 20–50 dye molecules per 1000 nt.

4. (optional) Calculate the RNA concentration

Use the A<sub>260</sub> to calculate the RNA concentration if desired (see section *III.A. Assessing aRNA yield by UV absorbance* on page 20).

#### E. Preparing Labeled aRNA for Hybridization

1. Concentrate the labeled aRNA if necessary

Typically, microarrays are hybridized with 25–100  $\mu L$  hybridization solution. Since the labeled aRNA is eluted in about 40  $\mu L$ , it may need to be concentrated to use it for microarray hybridization. We recommend vacuum drying the labeled aRNA in the dark until the volume is reduced to 1–10  $\mu L$  (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  $\mu$ 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
- 3. Dilute the labeled aRNA into the hybridization solution

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 (P/N AM8740) for this procedure. Follow the procedure associated with the fragmentation method used.

#### 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 µg 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 (P/N AM8861– AM8863).

### V. Troubleshooting

#### A. Positive Control Reaction

### Control RNA amplification instructions

To establish if the reaction conditions are optimal, a tube of Control RNA is provided.

- 1. Use 1  $\mu$ L (1  $\mu$ g) of the Control RNA in an Amino Allyl MessageAmp II-96 reaction; follow the protocol starting at step II.C.1 on page 13.
- 2. At step <u>II.F.2</u> on page 18, use a *14 hr* incubation for the IVT reaction.
- Continue with the procedure for making amino allyl aRNA through section II.G.

# 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. on page 20; the positive control reaction should produce  $\geq 80$  µg of aRNA.

#### Qualitative analysis

Also, run 100-200 ng of the reaction products on a bioanalyzer as described in section <u>III.B</u> on page 20; the expected aRNA profile for unmodified aRNA is a distribution from 250-5500 nt with a peak centered at 1000-1500 nt (Figure 4).

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

#### Labeling the aRNA produced in the positive control amplification reaction

Use 15  $\mu$ g of aRNA (made from Control RNA) in a dye coupling reaction following the instructions in section  $\underline{IV.A}$  through  $\underline{IV.D}$  starting on page  $\underline{21}$ .

Evaluate the positive control reaction by determining the number of dye molecules incorporated per 1000 nt as described in section <u>IV.D</u> on page 22.

The expected incorporation rate is 30–60 dye molecules per 1000 nt.

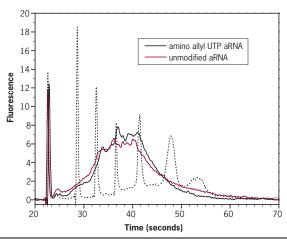


Figure 4. Bioanalyzer Electropherogram of aa-UTP Modified aRNA and Unmodified aRNA.

Double stranded cDNA produced with the Amino Allyl MessageAmp™ II-96 Kit was used as template for in vitro transcription reactions containing either 50% aaUTP or 100% UTP. Approximately 500 ng of aRNA was analyzed on an Agilent bioanalyzer with the RNA 6000 Ladder. The Ladder consists of six RNA transcripts with lengths of 0.2, 0.5, 1, 2, 4, and 6 kb. aRNA of similar size and yield are produced in reactions containing either 50% aaUTP or 100% UTP.

#### 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.

# 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  $1-2~\mu L$  of condensate in an IVT reaction changes the concentrations of the nucleotides and magnesium which can reduce yields.

*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 result in cleavage of 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 (P/N 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.

- 1. Be sure to resuspend RNA Binding Beads just before each step in preparing the aRNA Binding Mix (section <u>II.B.3</u> on page 12).
- 2. Make sure that magnetic bead mixtures are fully resuspended before adding them to cDNA or aRNA samples for purification procedures (steps <u>II.E.1</u> on page 15, and <u>II.G.1</u> on page 18).
- 3. For efficient elution of purified reaction products, make sure that magnetic beads are fully resuspended in the elution solution (steps <u>II.E.5</u> on page 16 and <u>II.G.5</u> on page 19).
- 4. Avoid overdrying magnetic beads before eluting the RNA (in steps II.E.4.e on page 16 and II.G.4.e on page 18) 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 <u>II.E.</u>, 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 MEGAclear™ Kit (P/N 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 RNA sample by determining the size of the 18S and 28S rRNA bands and the relative abundance of 28S to 18S rRNA (See section <u>II.A. RNA integrity</u> on page 9 for more information).

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

Different RNA samples contain different amounts of mRNA. In healthy cells, mRNA constitutes  $1{\text -}3\%$  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  $10{\text -}30$  ng per  $\mu g$  of total RNA (assuming good RNA quality). Most total RNA samples can be amplified up to 1000 fold producing  $10{\text -}30$   $\mu g$  of aRNA from 1  $\mu g$  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.

#### E. Troubleshooting Dye Coupling

NHS ester dye coupling is a simple and robust chemical reaction, but unreacted NHS ester dye solutions are susceptible to hydrolysis and photobleaching. Dye esters that have been improperly handled or stored may lose much of their reactivity. Refer to information provided by the NHS ester dye supplier for more details.



### VI. Appendix

#### A. 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.

Kacharmina 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 tanscription 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.

#### **B.** Quality Control

#### Functional testing

The Control RNA (2  $\mu g$ ) is used in an aRNA amplification experiment following the instructions in section  $\underline{V.A}$  on page 25. The amino allyl aRNA yield is assessed by measuring the  $A_{260}$  on the NanoDrop ND1000A spectrophotometer. Cy5 is used to label the 10  $\mu g$  of the amino allyl aRNA following the instructions in section IV starting on page 21; 20–50 dye molecules per 1000 nt are incorporated.

#### **Nuclease testing**

Relevant kit components are tested in the following nuclease assays:

#### RNase activity

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

#### Nonspecific endonuclease activity

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

#### **Exonuclease activity**

Meets or exceeds specification when a sample is incubated with labeled double-stranded DNA, followed by PAGE analysis.

#### Protease testing

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

#### C. Safety Information

#### Chemical safety guidelines

To minimize the hazards of chemicals:

- Read and understand the Material Safety Data Sheets (MSDS) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety goggles, gloves, or protective clothing). For additional safety guidelines, consult the MSDS.
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood). For additional safety guidelines, consult the MSDS.
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended on the MSDS.
- Comply with all local, state/provincial, or national laws and regulations related to chemical storage, handling, and disposal.

#### **About MSDSs**

Chemical manufacturers supply current Material Safety Data Sheets (MSDSs) with shipments of hazardous chemicals to new customers. They also provide MSDSs with the first shipment of a hazardous chemical to a customer after an MSDS has been updated. MSDSs provide the safety information you need to store, handle, transport, and dispose of the chemicals safely.

Each time you receive a new MSDS packaged with a hazardous chemical, be sure to replace the appropriate MSDS in your files.

#### **Obtaining the MSDS**

To obtain Material Safety Data Sheets (MSDSs) for any chemical product supplied by Applied Biosystems or Ambion:

 At www.appliedbiosystems.com, select Support, then MSDS. Search by chemical name, product name, product part number, or MSDS part number. Right-click to print or download the MSDS of interest.

- At www.ambion.com, go to the web catalog page for the product of interest. Click MSDS, then right-click to print or download.
- E-mail (MSDS\_Inquiry\_CCRM@appliedbiosystems.com), telephone (650-554-2756; USA), or fax (650-554-2252; USA) your request, specifying the catalog or part number(s) and the name of the product(s). The associated MSDSs will be e-mailed unless you request fax or postal delivery. Requests for postal delivery require 1 to 2 weeks for processing.

Note: For the MSDSs of chemicals not distributed by Applied Biosystems or Ambion, contact the chemical manufacturer.