

USER GUIDE

Ovation® PicoSL WTA System V2

PART NO. 3312

Patents, Licensing and Trademarks

©2019 Tecan Genomics, Inc. All rights reserved. The Encore®, Ovation® and Applause® families of products and methods of their use are covered by several issued U.S. and International patents and pending applications (www.nugen.com). NuGEN, Allegro, Celero, NuQuant, Solo, Metaplex, DimerFree, AnyDeplete, Ovation, SPIA, Ribo-SPIA, Applause, Encore and Imagine More From Less are trademarks or registered trademarks of Tecan Genomics, Inc. Other marks appearing in these materials are marks of their respective owners.

The purchase of this product conveys to the buyer the limited, non-exclusive, non-transferable right (without the right to modify, reverse engineer, resell, repackage or further sublicense) under these patent applications and any patents issuing from these patent applications to use this product and methods, accompanying this user guide, for research and development purposes solely in accordance with the intended use described and the written instructions provided in this user guide. No license to make or sell products by use of this product is granted to the buyer whether expressly, by implication, by estoppels or otherwise. In particular, the purchase of this product does not include or carry any right or license to use, develop or otherwise exploit this product commercially and no rights are conveyed to the buyer to use the product or components of the product for purposes including commercial services or clinical diagnostics.

For information on purchasing a license to Tecan Genomics' patents for uses other than in conjunction with this product or to use this product for purposes other than research, please contact Tecan Genomics, Inc., 900 Chesapeake Drive, Redwood City, CA 94063. Phone 888-654-6544 or 650-590-3600; FAX 888-296-6544 or 650-590-3630.

Warranty

Tecan Genomics warrants that this product meets the performance standards described in the Company's product and technical literature for a period of six months from the date of purchase, provided that the product is handled and stored according to published instructions, and that the product is not altered or misused. If the product fails to meet these performance standards, Tecan Genomics will replace the product free of charge or issue a credit for the purchase price. Tecan Genomics' liability under this warranty shall not exceed the purchase price of the product. Tecan Genomics shall assume no liability for direct, indirect, consequential or incidental damages arising from the use, results of use or inability to use its products. Tecan Genomics reserves the right to change, alter or modify any product to enhance its performance and design.

Tecan Genomics' products are developed, designed and sold FOR RESEARCH USE ONLY. This product is not to be used for diagnostic or therapeutic purposes, nor is it to be administered to humans or animals.

Except as expressly set forth herein, no right to modify, reverse engineer, distribute, offer to sell or sell Tecan Genomics' product is conveyed or implied by buyer's purchase of this Tecan Genomics product. The buyer agrees to use Tecan Genomics products accompanying the product insert in accordance with the intended use and the written instructions provided.

Technical Support

For help with any of our products, please contact Tecan's NGS Technical Support at 650.590.3674 (direct) or 888.654.6544, option 2 (toll-free, U.S. only). You may also email Tecan's NGS Technical Support at techserv-gn@tecan.com.

In Europe contact Tecan NGS Technical Support at +31 13 5780215 (Phone) or +31 13 5780216 (Fax) or europe-gn@tecan.com.

In all other locations, contact your Tecan distributor for technical support.

Table of Contents

Contents

I.	Introduction	1
A.	Background.....	1
B.	Ribo-SPIA [®] Technology	2
C.	Performance Specifications.....	3
D.	Quality Control	3
E.	Storage and Stability	3
F.	Safety Data Sheet (SDS).....	3
II.	Kit Components	4
A.	Reagents and Supplies Provided.....	4
B.	Additional Equipment, Reagents and Labware	5
III.	Planning the Experiment	7
A.	Input RNA Requirements	7
B.	Using RNase-free Techniques	8
C.	RNA Storage.....	9
D.	SPIA [®] cDNA Storage.....	9
IV.	Protocol	10
A.	Overview.....	10
B.	Agencourt [®] Purification Beads.....	12
C.	Programming the Thermal Cycler.....	14
D.	First Strand cDNA Synthesis	15
E.	Second Strand cDNA Synthesis	16
F.	Purification of cDNA	17
G.	SPIA Amplification	18
H.	Purification of SPIA cDNA	19
I.	Measuring SPIA cDNA Yield and Purity.....	19
V.	Technical Support	21
VI.	Appendix	22
A.	Purification Protocols for SPIA cDNA	22
B.	Guidelines for Use in Microarray Experiments	26
C.	Use with Affymetrix [®] GeneTitan [®] Systems	27
D.	Performing Quantitative PCR on SPIA cDNA	31
E.	Quality Control of the SPIA cDNA Product	32
F.	DNase Treatment of RNA.....	34
G.	Preventing Non-specific Amplification	37
H.	Frequently Asked Questions (FAQs)	39
I.	Update History	43

I. Introduction

A. Background

The Ovation® PicoSL WTA System V2 provides a fast and simple method for preparing amplified cDNA for gene expression analysis. Amplification is initiated at the 3' end as well as randomly throughout the transcriptome in the sample, making the Ovation PicoSL WTA System V2 ideal for amplification of partially degraded and compromised RNA samples. The amplified product of the Ovation PicoSL WTA System V2 is optimized for gene expression analysis applications on Affymetrix® Array Plates, Agilent Gene Expression microarrays and Illumina Genome-Wide Expression BeadChips utilizing the appropriate Tecan Genomics® labeling modules and protocols. The amplified cDNA can also be used as input for qPCR reactions. For details please visit the website.

The Ovation PicoSL WTA System V2 is powered by Ribo-SPIA® technology, a rapid, simple and sensitive strand-displacement amplification process developed by Tecan Genomics. This high fidelity method uses DNA/RNA chimeric primers (SPIA primers), DNA polymerase and RNase H in an isothermal assay. The amplification process is highly reproducible, maintaining the stoichiometry of the input RNA sample and thereby faithfully preserving biological information. Starting with 500 pg to 50 ng of total RNA, microgram quantities of cDNA can be prepared in approximately 4.8 hours.

The Ovation PicoSL WTA System V2 (Part No. 3312) provides optimized reagent mixes and a protocol to process total RNA samples. Control RNA is not provided with the Ovation PicoSL WTA System V2 but we recommend the use of a control RNA when first using this product.

I. Introduction

B. Ribo-SPIA® Technology

Ribo-SPIA is a three-step process, based on Tecan Genomics' Single Primer Isothermal Amplification (SPIA) technology, that generates amplified cDNA from as little as 500 picograms of total RNA.

1. Generation of First Strand cDNA (1.2 hours)

First strand cDNA is prepared from total RNA using a unique first strand DNA/RNA chimeric primer mix and reverse transcriptase (RT). The primer mix contains a unique mixture of random and oligo dT primers such that priming occurs across the whole transcript. RT extends the 3' DNA end of each primer generating first strand cDNA.

The resulting cDNA/mRNA hybrid molecule contains a unique RNA tag sequence (SPIA tag) at the 5' end of the cDNA strand which will be used as a priming site for the SPIA process.

2. Generation of a DNA/RNA Heteroduplex Double Strand cDNA (2 hours)

Fragmentation of the mRNA within the cDNA/mRNA complex creates priming sites for DNA polymerase to synthesize a second cDNA strand, which includes DNA complementary to the 5' SPIA tag sequence from the first strand chimeric primers. The result is a double-stranded cDNA with a DNA/RNA heteroduplex corresponding to the SPIA tag at one end.

3. SPIA® Amplification (1.6 hours)

SPIA is a rapid, simple and sensitive strand-displacement amplification process developed by Tecan Genomics. It uses a DNA/RNA chimeric primer (SPIA primer), DNA polymerase and RNase H in an isothermal assay. RNase H removes the RNA portion of the heteroduplex SPIA tag sequence, revealing a site for binding the SPIA primer. DNA polymerase synthesizes cDNA starting at the 3' end of the primer, displacing the existing forward strand. Priming with the chimeric SPIA primer recapitulates the heteroduplex SPIA tag, creating a new substrate for RNase H and the initiation of the next round of cDNA synthesis. The process of SPIA DNA/RNA primer binding, DNA replication, strand displacement and RNA cleavage is repeated in a highly processive manner, resulting in rapid accumulation of micrograms of amplified cDNA from picograms of total RNA.

I. Introduction

C. Performance Specifications

The Ovation PicoSL WTA System V2 synthesizes microgram quantities of SPIA cDNA starting with total cellular RNA inputs of 500 pg to 50 ng. In approximately 4.8 hours the Ovation PicoSL WTA System V2 can produce 2–4 µg of cDNA ready for qPCR or other analytical tests. When used with intact input RNA, the size of the majority of the cDNA products produced by the Ribo-SPIA amplification process is between 50 bases and 1.5 Kb. When used with degraded input RNA the size of the SPIA cDNA products may be smaller, in proportion to the degree of input RNA degradation. With a whole transcriptome amplification approach, the size distribution of the product is less important compared to a 3' amplification strategy, since it results in densely overlapping cDNA fragments representing the entire transcriptome.

D. Quality Control

Each Ovation PicoSL WTA System V2 lot is tested to meet specifications of yield, qPCR and array performance.

E. Storage and Stability

The Ovation PicoSL WTA System V2 is shipped on dry ice and should be unpacked immediately upon receipt.

The vial labeled Agencourt® Beads (clear cap) should be removed from the top of the shipping carton upon delivery and stored at 4°C.

The vial labeled First Strand Enzyme Mix (blue: A3) should be removed from the kit upon delivery and stored separately at –80°C.

All remaining components should be stored at –20°C in a freezer without a defrost cycle.

This product has been tested to perform to specifications after as many as six freeze/thaw cycles. Kits handled and stored according to the above guidelines will perform to specifications for at least six months. Tecan Genomics has not yet established long-term storage conditions for the Ovation PicoSL WTA System V2.

F. Safety Data Sheet (SDS)

An SDS for this product is available on the website at <http://www.nugen.com/products/ovation-picosl-wta-system-v2>.

 This product contains components with multiple storage conditions

 Store the First Strand Primer Mix, First Strand Enzyme Mix, and SPIA Primer Mix at –80°C and the Agencourt Beads at 4°C

II. Kit Components

A. Reagents and Supplies Provided

Table 1. First Strand cDNA Reagents

COMPONENT	3312-24 PART NUMBER	3312-48 PART NUMBER	VIAL CAP	VIAL NUMBER
First Strand Primer Mix	S01493	S01403	Blue	A1 VER 8
First Strand Buffer Mix	S01494	S01191	Blue	A2 VER 3
First Strand Enzyme Mix	S02271	S02251	Blue	A3 VER 7

Table 2. Second Strand cDNA Reagents

COMPONENT	3312-24 PART NUMBER	3312-48 PART NUMBER	VIAL CAP	VIAL NUMBER
Second Strand Buffer Mix	S01496	S01192	Yellow	B1 VER 3
Second Strand Enzyme Mix	S01377	S01193	Yellow	B2 VER 2

Table 3. SPIA Reagents

COMPONENT	3312-24 PART NUMBER	3312-48 PART NUMBER	VIAL CAP	VIAL NUMBER
SPIA Primer Mix	S01497	S01400	Red	C1 VER 9
SPIA Buffer Mix	S01498	S01401	Red	C2 VER 10
SPIA Enzyme Mix	S01499	S01402	Red	C3 VER 7

II. Kit Components

Table 4. Additional Reagents

COMPONENT	3312-24 PART NUMBER	3312-48 PART NUMBER	VIAL CAP	VIAL NUMBER
Nuclease-free Water	S01001	S01001	Green	D1
Agencourt® Beads	S01307	S01307	Clear	—

Note: The reagents in the Ovation PicoSL WTA System V2 are similar to reagents in our other kits; however, unless the part numbers are identical, these reagents do not have exactly the same composition and, therefore, are not interchangeable. Do not exchange or replace one reagent named, for example, A1 with another A1, as it will adversely affect performance.

B. Additional Equipment, Reagents and Labware

Required Materials

- Equipment
 - Microcentrifuge for individual 1.5 mL and 0.5 mL tubes
 - 0.5 to 10 µL pipette, 2 to 20 µL pipette, 20 to 200 µL pipette, and 200 to 1000 µL pipette
 - Vortexer
 - Thermal cycler with 0.2 mL tube heat block, heated lid, and 100 µL reaction capacity
 - Appropriate spectrophotometer and cuvettes, or Nanodrop® UV-Vis Spectrophotometer
- Reagents
 - Ethanol (Sigma-Aldrich, Cat. #E7023), for purification steps
- Supplies and Labware
 - Nuclease-free pipette tips
 - 1.5 mL and 0.5 mL RNase-free microcentrifuge tubes
 - Low-retention microcentrifuge tubes (SafeSeal Low Binding 0.65 mL Microcentrifuge Tubes, Sorenson Biosciences, Inc., Cat #11300)
 - 0.2 mL individual thin wall PCR tubes, 8 X 0.2 mL strip PCR tubes or 0.2 mL thin wall PCR plates
 - SPRIPlate® 96R, Ring Magnet Plate (Beckman Coulter Genomics, Cat. #A29164) or SPRIPlate Ring Super Magnet Plate, (Beckman Coulter Genomics, Cat. #A32782). Other magnetic stands may be used as well, although their performance has not been validated by Tecan Genomics.

II. Kit Components

- Purification options for final SPIA cDNA purification (select one option):
 - Agencourt® Beads (Beckman Coulter Genomics, Cat. #A63880 or A63987)
 - MinElute® Reaction Cleanup Kit (QIAGEN®, Cat. #28204)
 - QIAquick® PCR Purification Kit (QIAGEN, Cat. #28104)
 - DNA Clean & Concentrator™-25 (Zymo Research, Cat. #D4005/D4006)
- Disposable gloves
- Kimwipes
- Ice bucket
- Cleaning solutions such as RNase Zap® (Thermo Fisher Scientific, Cat.#AM9780) and DNA-OFF™ (MP Biomedicals, Cat.#QD0500)
- Optional Equipment
 - Agilent 2100 Bioanalyzer or materials and equipment for electrophoretic analysis of RNA
 - Real-time PCR system

To Order:

- Beckman Coulter Genomics, www.beckmangenomics.com
- MP Biomedicals, www.mpbio.com
- QIAGEN Inc., www.qiagen.com
- Sigma-Aldrich, Inc., www.sigmaaldrich.com
- Sorenson Biosciences, Inc., www.sorbio.com
- Thermo Fisher Scientific, www.thermofisher.com
- Zymo Research, www.zymoresearch.com

III. Planning the Experiment

A. Input RNA Requirements

It is important to assess the quality of your RNA sample prior to planning your amplification. While the Ovation PicoSL WTA System V2 will allow the amplification of many RNA samples of variable quality, use of highly degraded RNA samples can lead to lower yields and shorter SPIA cDNA. It is impossible to guarantee success with all degraded RNA samples. To assess RNA quality prior to using the Ovation PicoSL WTA System V2, follow the guidelines below.

1. RNA Quantity

Total RNA input must be between 500 pg and 50 ng. Inputs above 50 ng per reaction may inhibit amplification, while lower amounts of input will potentially result in insufficient yields depending on the requirements of the analytical platform. We strongly recommend quantification of total RNA to ensure the minimum input requirement is met.

2. RNA Purity

RNA samples must be free of contaminating proteins and other cellular material, organic solvents (including phenol and ethanol) and salts used in many RNA isolation methods. Use of a commercially available system for preparing small amounts of RNA that does not require organic solvents is recommended. If a method such as Trizol is used, we recommend using a column purification after isolation in order to remove any residual organic solvents that may be present. One measure of RNA purity is the ratio of absorbance readings at 260 and 280 nm. The A260:A280 ratio for RNA samples should be in excess of 1.8. RNA samples with lower ratios may result in low amplification yield.

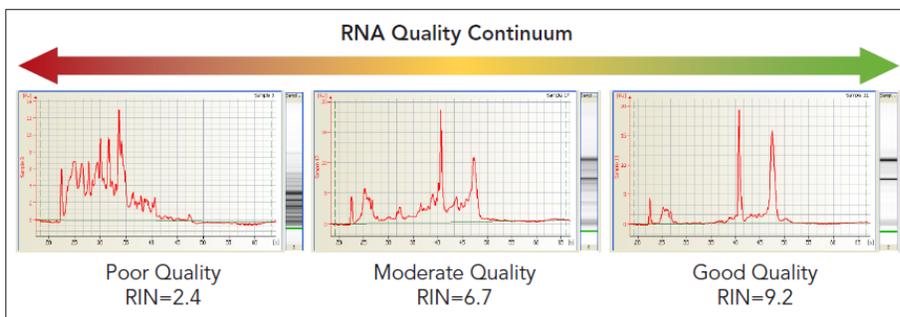
3. RNA Integrity

RNA samples of high molecular weight with little or no evidence of degradation will amplify very well with this product. Due to the whole transcriptome amplification approach, even lower quality, compromised RNA samples can be amplified successfully using the Ovation PicoSL WTA System V2.

RNA integrity can be determined using the Agilent 2100 Bioanalyzer, RNA 6000 Nano LabChip® or RNA 6000 Pico LabChip®, and the RNA Integrity Number (RIN) calculation available in the Bioanalyzer 2100 Expert Software. The instrument provides a sensitive and rapid way of confirming RNA integrity prior to amplification, both visually, with a detailed electrophoretic trace of the RNA, and computationally, by calculating a RIN score. While it is impossible to guarantee satisfactory results with all degraded samples, the Ovation PicoSL WTA System V2 can work with many samples that are moderately to severely degraded. In our tests using an RNA degradation model system, RNA samples that are severely degraded showing RIN scores of approximately 2 to 5 still amplify successfully and reproducibly. On occasions when the Bioanalyzer software fails to calculate a RIN score, we recommend viewing the electrophoretic trace to determine if the sample may still be of adequate integrity for use.

III. Planning the Experiment

Figure 1. This continuum of RNA quality shows Bioanalyzer traces of three different RNAs with varying levels of degradation, all of which have amplified successfully using the Ovation PicoSL WTA System V2 approach.



4. DNase Treatment

It is generally recommended to use DNase-treated RNA for amplification using the Ovation PicoSL WTA System V2. The presence of genomic DNA in the RNA sample may have adverse effects on downstream analytical platforms. Contaminating genomic DNA may be amplified along with the RNA. Additionally, if the total RNA sample contains a significant amount of contaminating genomic DNA, it will be difficult to accurately quantify the true RNA concentration. The RNA input quantity may, therefore, be over-estimated based on an absorbance measurement. Since it is important that RNA input be within the stated range of 500 pg to 50 ng, we recommend using a DNase treatment that will remove contaminating genomic DNA during RNA purification.

5. Carrier use for RNA isolation

We strongly recommend against the use of nucleic acid based carriers during RNA purification because many have been shown to produce cDNA product in first strand synthesis. We also advise against the use of glycogen in RNA isolation, as it inhibits reverse transcription. For the latest information regarding other carriers, contact the Tecan Technical Services Team.

B. Using RNase-free Techniques

RNase contamination through reagents and work environment will lead to experimental failure. Follow these guidelines to minimize contamination:

- Wear disposable gloves and change them frequently.
- Avoid touching surfaces or materials that could introduce RNases.
- Use reagents provided. Substitutions may introduce RNases.
- Clean and decontaminate work areas and instruments, including pipettes, with commercially available cleaning reagents, such as RNaseZap.
- Use only new RNase-free pipette tips and microcentrifuge tubes.
- Use a work area specifically designated for RNA work and do not use other high copy number materials in the same area.

III. Planning the Experiment

C. RNA Storage

RNA samples for use with the Ovation PicoSL WTA System V2 must be stored at -80°C . Avoid frequent freeze/thaw cycles or RNA degradation may result.

D. SPIA[®] cDNA Storage

The SPIA cDNA produced by the Ovation PicoSL WTA System V2 may be stored at -20°C .

IV. Protocol

A. Overview

The Ribo-SPIA amplification process used in the Ovation PicoSL WTA System V2 is performed in three stages:

1. First strand cDNA synthesis	1.2 hours
2. Second strand cDNA synthesis and purification	2 hours
3. SPIA amplification and purification	1.6 hours
<hr/>	
Total time to prepare SPIA cDNA	~4.8 hours

Ovation PicoSL WTA System V2 components are color coded, with each color linked to a specific stage of the process. Performing each stage requires making a master mix then adding it to the reaction, followed by incubation. Master mixes are prepared by mixing components provided for that stage.

The SPIA cDNA must be purified following amplification if you intend to use the cDNA for labeling using a Tecan Genomics Encore™ labeling module or other supported labeling protocol for applications such as microarray analysis.

The Ovation PicoSL WTA System V2 may also be used as a method of pre-amplification prior to qPCR. Although for qPCR applications it is not absolutely necessary, we recommend purifying the cDNA after SPIA. If quantification of the cDNA product is desired, it must be purified first. Spectrophotometric quantification of unpurified amplification products will result in artificially high readings due to amplification components present in the sample.

Protocol Notes

- We recommend the routine use of a positive control RNA. Especially the first time you set up an amplification reaction, the use of a positive control RNA will allow the establishment of a baseline of performance and provide the opportunity to become familiar with the bead purification step. This step may be unfamiliar to many users and can be especially prone to handling variability in using the magnet, so a practice run with the magnet is highly recommended.
- In working with very small, picogram amounts of RNA we strongly recommend the use of low retention tubes for storage and dilution of the samples in order to reduce the loss of RNA samples due to adhesion to polypropylene surfaces.
- Due to the high sensitivity inherent in this amplification system we strongly recommend taking measures to minimize the potential for the carryover of previously amplified SPIA cDNA into new amplification reactions. The two steps to accomplish this are: 1. Designating separate workspaces for “pre-amplification” and “post-amplification” steps and materials and 2. Implementing routine clean-up protocols for workspaces as standard operating procedure. A detailed set of these recommendations is given in the Appendix.
- Use the Nuclease-free Water provided with the kit (green: D1) or an alternate source of nuclease-free water. We do not recommend the use of DEPC-treated water with this protocol.

IV. Protocol

- Setting up a minimum of four reactions at a time ensures that you are not pipetting very small volumes (see the second strand synthesis section).
- The reagent volumes recovered greatly depend on the number of batches processed with each kit. Set up no fewer than four reactions at a time with a 24-reaction kit, and no fewer than 8 reactions with a 48-reaction kit. This ensures there will be sufficient reagent volumes to perform the full number of reactions specified for each kit size.
- Thaw components used in each step and immediately place them on ice. It is best to not thaw all reagents at once.
- Always keep thawed reagents and reaction tubes on ice unless otherwise instructed.
- After thawing and mixing buffer mixes, if any precipitate is observed, re-dissolve it completely prior to use. You may gently warm the buffer mix for two minutes at room temperature followed by brief vortexing. Do not warm any enzyme or primer mixes.
- When placing small amounts of reagents into the reaction mix, pipet up and down several times to ensure complete transfer.
- When instructed to pipet mix, gently aspirate and dispense a volume that is at least half of the total volume of the reaction mix.
- Always allow the thermal cycler to reach the initial incubation temperature prior to placing the tubes in the block.
- When preparing master mixes, use the minimal amount of extra material to ensure obtaining the full number of reactions in the kit. The Ovation PicoSL WTA System V2 Quick Protocol has been designed to automatically calculate an appropriate overfill volume, based on the desired number of reactions, which can be used as a guideline in setting up master mixes.
- Components and reagents from other Tecan Genomics products should not be interchanged with the components supplied with this product.
- Use only fresh ethanol stocks to make 70% ethanol used in the post-second strand bead purification, and ethanol for washes in the SPIA cDNA purification protocols. Make the ethanol mixes fresh as well, carefully measuring both the ethanol and water with pipettes. Lower concentrations of ethanol in wash solutions will result in loss of yield as the higher aqueous content will dissolve the cDNA and wash it off the beads or column.

IV. Protocol

B. Agencourt® Purification Beads

Tips and notes Relevant to the Second Strand cDNA Cleanup, Protocol G:

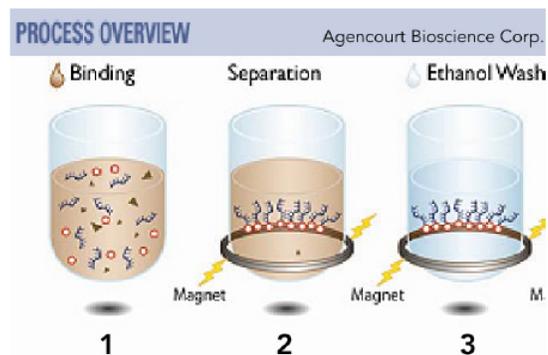
There are significant modifications to the Beckman Coulter Genomics' Agencourt SPRI Bead Kit standard procedure; therefore, you must follow the procedures outlined in this user guide for the use of these beads with the Ovation PicoSL WTA System V2.

The bead purification process used for cDNA purification before amplification consists of the following steps:

1. Binding of cDNA to beads
2. Magnetic separation of beads from supernatant
3. Ethanol wash of bound beads to remove contaminants

Elution takes place upon addition of the SPIA Master Mix. At this stage the beads are left in the reaction tube and removed only after amplification.

Figure 2. Bead purification process overview.



Additional Tips and Notes

- Remove beads from 4°C and leave at room temperature for at least 15 minutes before use. Ensure that they have completely reached room temperature. Cold beads will result in reduced recovery.
- Fully resuspend the beads by inverting and tapping before adding to the sample.
- Note that we recommend using 1.6 volumes (32 μ L) of beads. This is different from the standard Agencourt protocol.
- It is critical to let the beads separate on the magnet for a full five minutes. Removing the binding buffer before the beads have completely separated will impact cDNA yields.
- After the binding step has been completed, it is important to minimize bead loss when removing the binding buffer. With the samples placed on the magnet,

IV. Protocol

remove only 45 μ L of the binding buffer from each sample. Some liquid will remain at the bottom of the tube but this will minimize bead loss.

- Any significant loss of beads during the ethanol washes will impact cDNA yields, so make sure the beads are not lost with the wash.
- Ensure that the ethanol wash is freshly prepared from fresh ethanol stocks at the indicated concentration. Lower percent ethanol mixes will reduce recovery.
- During the ethanol washes, keep the samples on the magnet. The beads should not be allowed to disperse; the magnet will keep the beads on the walls of the sample tubes.
- It is critical that all residual ethanol be removed prior to continuing with the SPIA amplification. Therefore, when removing the final ethanol wash, first remove most of the ethanol, then allow the excess to collect at the bottom of the tube before removing the remaining ethanol. This reduces the required bead air drying time.
- After drying the beads for 15 to 20 minutes, inspect each tube carefully and make certain that all the ethanol has evaporated before proceeding with the amplification step.
- It is strongly recommended that strip tubes or partial plates are firmly placed when used with the magnetic plate. We do not advise the use of individual tubes as they are not very stable on the magnet.

IV. Protocol

C. Programming the Thermal Cycler

Use a thermal cycler with a heat block designed for 0.2 mL tubes, equipped with a heated lid, and with a capacity of 100 μ L reaction volume. Prepare the programs shown in Table 5, following the operating instructions provided by the manufacturer. For thermal cyclers with an adjustable heated lid, set the lid temperature at 100°C. For thermal cyclers with a fixed temperature heated lid (e.g., ABI GeneAmp[®] PCR 9600 and 9700 models) use the default settings (typically 100 to 105°C).

Table 5. Thermal Cycler Programming

FIRST STRAND cDNA SYNTHESIS	
Program 1 Primer Annealing	65°C – 2 min, hold at 4°C
Program 2 First Strand Synthesis	4°C – 2 min, 25°C – 30 min, 42°C – 15 min, 70°C – 15 min, hold at 4°C
SECOND STRAND cDNA SYNTHESIS	
Program 3 Second Strand Synthesis	4°C – 1 min, 25°C – 10 min, 50°C – 30 min, 80°C – 20 min, hold at 4°C
SPIA AMPLIFICATION	
Program 4 SPIA Amplification	4°C – 1 min, 47°C – 75 min, 95°C – 5 min, hold at 4°C

IV. Protocol

D. First Strand cDNA Synthesis

Important Note: Carry out Protocol D (First-Strand cDNA Synthesis) through Protocol G, step 7 (SPIA Amplification) in a pre-amplification workspace using dedicated pre-amplification consumables and equipment. Wipe all surfaces, equipment and instrumentation with a DNA removal solution such as DNA-OFF (MP Biomedicals, Cat. #Q0500) to avoid the potential introduction of previously amplified cDNA into new amplifications. For more information on our recommendations for workflow compartmentalization and routine lab cleanup please refer to Appendix G. If you have any questions on this important topic, please contact the Tecan's NGS Technical Support at 650.590.3674 (direct) or 888.654.6544, option 2 (toll-free, U.S. only). You may also email Tecan's NGS Technical Support at techserv-gn@tecan.com.



Do not vortex any enzyme mixes.

1. Obtain the First Strand Primer Mix (blue: A1 VER 8), First Strand Buffer Mix (blue: A2 VER 3) and the Nuclease-free Water (green: D1) from the -20°C storage and the First Strand Enzyme Mix (blue: A3 VER 7) from the -80°C storage.
2. Spin down the contents of A3 and place on ice.
3. Thaw the other reagents at room temperature, mix by vortexing, spin and place on ice. Leave the Nuclease-free Water at room temperature.
4. Add 2 μL of A1 to a 0.2 mL PCR tube.
5. Add 5 μL of total RNA sample (500 pg to 50 ng) to the primer.
6. Mix by pipetting 5 times, spin and place on ice.
7. Place the tubes in a pre-warmed thermal cycler programmed to run Program 1 (Primer Annealing; see Table 5):
65°C – 2 min, hold at 4°C
8. Remove the tubes from the thermal cycler and place on ice.
9. Once Primer Annealing (Step 7) is complete, prepare a master mix by combining A2 and A3 in a 0.5 mL capped tube, according to the volumes shown in Table 6.

Table 6. First Strand Master Mix (volumes listed are for a single reaction)

FIRST STRAND BUFFER MIX (BLUE: A2 <small>VER 3</small>)	FIRST STRAND ENZYME MIX (BLUE: A3 <small>VER 7</small>)
2.5 μL	0.5 μL



Mix by pipetting and spin down the master mix briefly. Immediately place on ice.

10. Add 3 μL of the First Strand Master Mix to each tube.

IV. Protocol



The second strand reagents may be thawed and put on ice 10 minutes before the completion of First Strand cDNA Synthesis



The purification beads should be removed from 4°C storage and left on the bench top to reach room temperature well before the start of purification.



In order to ensure accurate measurement of B2, do not make this mix for fewer than 4 reactions.



Mix by pipetting and spin down the master mix briefly. Immediately place on ice.

11. Mix by pipetting 5 times, spin and place on ice.
12. Place the tubes in a pre-cooled thermal cycler programmed to run Program 2 (First Strand cDNA Synthesis; see Table 5):
4°C – 2 min, 25°C – 30 min, 42°C – 15 min, 70°C – 15 min, hold at 4°C
13. Remove the tubes from the thermal cycler, spin to collect condensation and place on ice.
14. Continue immediately with the Second Strand cDNA Synthesis protocol.

E. Second Strand cDNA Synthesis

1. Remove the Agencourt Beads (supplied with the Ovation PicoSL WTA System V2) from 4°C storage and place on the bench top to reach room temperature for use in the next step.
2. Obtain the Second Strand Buffer Mix (yellow: B1 ver 3) and Second Strand Enzyme Mix (yellow: B2 ver 2) from –20°C storage.
3. Spin down the contents of B2 and place on ice.
4. Thaw reagent B1 at room temperature, mix by vortexing, spin and place on ice.
5. Make a master mix by combining B1 and B2 in a 0.5 mL capped tube, according to the volumes shown in Table 7.

Table 7. Second Strand Master Mix (volumes listed are for a single reaction)

SECOND STRAND BUFFER MIX (YELLOW: B1 ver 3)	SECOND STRAND ENZYME MIX (YELLOW: B2 ver 2)
9.7 µL	0.3 µL

6. Add 10 µL of the Second Strand Master Mix to each First Strand reaction tube.
7. Mix by pipetting 5 times, spin and place on ice.
8. Place the tubes in a pre-cooled thermal cycler programmed to run Program 3 (Second Strand cDNA Synthesis; see Table 5):
4°C – 1 min, 25°C – 10 min, 50°C – 30 min, 80°C – 20 min, hold at 4°C
9. Remove the tubes from the thermal cycler and spin to collect condensation. Place in a rack on the bench top.
10. Continue immediately with the Purification of cDNA protocol.

IV. Protocol

F. Purification of cDNA

1. Ensure the Agencourt Beads have completely reached room temperature before proceeding.
2. Prepare a 70% ethanol wash solution. It is critical that this solution be prepared fresh on the same day of the experiment from a recently opened stock container. Measure both the ethanol and the water components carefully prior to mixing. Failure to do so can result in a higher than anticipated aqueous content, which may reduce amplification yield.
3. Resuspend the beads by inverting and tapping the tube. Ensure that the beads are fully resuspended before adding to the sample. After resuspending do not spin the beads. A large excess of beads is provided; therefore, it is not necessary to recover any trapped in the cap.
4. At room temperature, add 32 μL (1.6 volumes) of the bead suspension to each reaction and mix by pipetting 10 times.
5. Incubate at room temperature for 10 minutes.
6. Transfer the tubes to the magnet and let stand 5 minutes to completely clear the solution of beads.
7. Keeping the tubes on the magnet, carefully remove only 45 μL of the binding buffer and discard it. Leaving some of the volume behind minimizes bead loss at this step.
8. With the tubes still on the magnet, add 200 μL of freshly prepared 70% ethanol and allow to stand for 30 seconds.
Note: The beads should not disperse; instead they will stay on the walls of the tubes. Significant loss of beads at this stage will impact cDNA yields, so ensure beads are not removed with the binding buffer or the washes.
9. Remove the 70% ethanol wash using a pipette.
10. Repeat the wash 2 more times.
Note: With the final wash, it is critical to remove as much of the ethanol as possible. Use at least 2 pipetting steps and allow excess ethanol to collect at the bottom of the tubes after removing most of the ethanol in the first pipetting step.
11. Air dry the beads on the magnet for 15 to 20 minutes. Inspect each tube carefully to ensure that all the ethanol has evaporated. It is critical that all residual ethanol be removed prior to continuing with SPIA Amplification.
12. Continue immediately with the SPIA Amplification protocol with the cDNA still bound to the dry beads.

! Minimize bead loss by leaving a residual volume of binding buffer after completion of the binding step.

! Best results are obtained by using fresh 70% EtOH in wash step.

IV. Protocol



Ensure the enzyme is well mixed without introducing bubbles.



Use SPIA Master Mix immediately after preparation.



Mix by pipetting and spin down the master mix briefly. Immediately place on ice.

G. SPIA Amplification

1. Obtain the SPIA Primer Mix (red: C1 ver 9), SPIA Buffer Mix (red: C2 ver 10) and SPIA Enzyme Mix (red: C3 ver 7) from -20°C storage.
2. Thaw C3 on ice and mix the contents by inverting gently 5 times, spin and place on ice. Ensure the enzyme is well mixed without introducing bubbles.
3. Thaw reagents C1 and C2 at room temperature, mix by vortexing, spin and place on ice.
4. Make a master mix by sequentially combining C2, C1 and C3 in an appropriately sized capped tube, according to the volumes shown in Table 8.

Note: Make sure the addition of C3 is at the last moment.

Table 8. SPIA Master Mix (volumes listed are for a single reaction)

SPIA BUFFER MIX (RED:C2 ver 10)	SPIA PRIMER MIX (RED:C1 ver 9)	SPIA ENZYME MIX (RED:C3 ver 7)
20 μL	10 μL	10 μL

5. Add 40 μL of the SPIA Master Mix to each tube containing the double-stranded cDNA bound to the dried beads. Mix thoroughly by pipetting at least 8 to 10 times. Attempt to get the majority of the beads in suspension and remove most of the beads from the tube walls.
Note: Beads may not form a perfectly uniform suspension, but this will not affect the reaction. The addition of SPIA master mix will elute the cDNA from the beads.
6. Place the tubes in a pre-cooled thermal cycler programmed to run Program 4 (SPIA Amplification, see Table 5):
4 $^{\circ}\text{C}$ – 1 min, 47 $^{\circ}\text{C}$ – 75 min, 95 $^{\circ}\text{C}$ – 5 min, hold at 4 $^{\circ}\text{C}$
7. Remove the tubes from the thermal cycler, spin to collect condensation and place on ice. Do not re-open the tubes in the pre-amplification workspace.

IV. Protocol

Important Note: At this point the tubes should be removed from the pre-amplification workspace. Carry out all remaining steps in a post-amplification workspace using dedicated post-amplification consumables and equipment. Take care to avoid the introduction of previously amplified cDNA into your pre-amplification workspace. For more information on our recommendations for workflow compartmentalization and routine lab cleanup, please refer to Appendix G of this user guide. If you have any questions on this important topic, please contact Tecan's NGS Technical Support at 650.590.3674 (direct) or 888.654.6544, option 2 (toll-free, U.S. only). You may also email Tecan's NGS Technical Support at techserv-gn@tecan.com.

Note: If using the Agencourt Beads for final SPIA cDNA cleanup, skip steps 8–10 below and go directly to the Agencourt bead protocol on 22. It is not necessary to remove the beads at this point.

8. Transfer the tubes to the magnet and let stand for 5 minutes to completely clear the solution of beads.
9. Carefully remove all of the cleared supernatant containing the eluted SPIA cDNA and transfer to a fresh tube. The beads may now be discarded.
10. Continue immediately with the Purification of SPIA cDNA protocol or store the reaction products at -20°C prior to continuing.

H. Purification of SPIA cDNA

The SPIA cDNA product can be purified using various methods listed in Appendix A. Purification is required if the SPIA cDNA is intended for use in an Encore labeling module or other supported labeling protocol. We recommend that the SPIA cDNA be purified prior to qPCR analysis.

I. Measuring SPIA cDNA Yield and Purity

1. Mix the purified SPIA cDNA sample by brief vortexing and spinning prior to checking the concentration.
2. Measure the absorbance of the SPIA cDNA at 260, 280 and 320 nm. You may need to make a 1:20 dilution of the cDNA in water prior to measuring the absorbance.
3. Purity: Subtract the A₃₂₀ value from both A₂₆₀ and A₂₈₀ values. The adjusted $(A_{260} - A_{320}) / (A_{280} - A_{320})$ ratio should be >1.8 .
4. Yield: Assume 1 A₂₆₀ unit = 33 $\mu\text{g}/\text{mL}$ for single-stranded cDNA.

IV. Protocol

To calculate: $(A_{260} - A_{320} \text{ of diluted sample}) \times (\text{dilution factor}) \times 33$ (concentration in $\mu\text{g}/\text{mL}$ of a 1 A260 unit solution) $\times 0.03$ (final volume in mL) = total yield in micrograms

5. Alternatively, you may measure the concentration and purity of the SPIA cDNA with a Nanodrop, using the ssDNA setting or using 1 A260 unit = 33 $\mu\text{g}/\text{mL}$ as the constant.
6. The purified SPIA cDNA may be stored at -20°C .

V. Technical Support

For support in the U.S., please contact Tecan's NGS Technical Support at 650.590.3674 (direct) or 888.654.6544, option 2 (toll-free, U.S. only). You may also email Tecan's NGS Technical Support at techserv-gn@tecan.com.

In Europe contact Tecan NGS Technical Support at +31 13 5780215 (Phone) or +31 13 5780216 (Fax) or europe-gn@tecan.com

In all other locations, contact your Tecan distributors Technical Support Team.

VI. Appendix

A. In all other locations, contact your Tecan distributors Technical Support Team. Purification Protocols for SPIA cDNA

There are two supported alternatives for carrying out the final purification of SPIA cDNA: 1) the Agencourt Beads and 2) the QIAGEN MinElute Reaction Cleanup Kit.

Selection of the optimum purification option can depend on many factors. Generally speaking, the column based option provides for superior recovery (higher yield) while the Agencourt bead method is more scalable for larger batch sizes and can be easily automated. We have found that the Agencourt bead method may result in somewhat lower amplification yields, so this option may not be appropriate for the most demanding applications where small and/or degraded RNA samples are anticipated. Please contact the Tecan Technical Support Team for assistance in selecting the appropriate purification option for your application.

The procedures given below are specifically adapted for use with Tecan Genomics products and may differ significantly from the protocols published by the manufacturers. Failure to follow the purification procedures as given below may negatively impact your results.

Agencourt Beads (instructions for a single reaction)

Important notes:

- Stop after step 7 on 18. It is not necessary to remove the beads from the SPIA reactions at this point.
 - Prepare a room temperature 80% ethanol wash solution. It is critical that this solution be prepared fresh on the same day of the experiment from a recently opened stock container. Measure both the ethanol and the water components carefully prior to mixing. Failure to do so can result in a higher than anticipated aqueous content which may reduce amplification yield.
 - The use of 96-well microplates and multi-channel pipettes is recommended for processing large batches with this procedure.
1. Obtain the AMPure XP or RNAClean XP bead bottle from 4°C storage. Allow the bead solution to reach room temperature.
 2. Invert the bead bottle several times to ensure the beads are fully in suspension. It may be necessary to remix the bead stock from time to time to ensure beads remain in suspension while in use.
 3. At room temperature, add 72 µL of resuspended beads (1.8 times the sample volume) to the 40 µL SPIA reaction.
 4. Mix the sample and beads thoroughly by pipetting 10 times.
 5. Incubate the sample/bead mixture at room temperature for 5 minutes.
 6. Place the sample/bead mixture on the magnet for 10 minutes to completely clear the solution of beads.
 7. Keeping the tube on the magnet, carefully remove and discard the supernatant from the first set of samples. Do not disturb the ring of beads.

VI. Appendix

Note: Leaving several microliters of supernatant behind at this step can help mini-mize bead loss.

8. Keeping the tube on the magnet, add 200 μL of freshly prepared 80% ethanol to each sample and incubate for 30 seconds or until the solution clears. Add slowly so as not to disturb the separated magnetic beads.
9. Carefully remove and discard the ethanol.
10. Repeat the 80% ethanol wash once more. Ensure as much ethanol as possible is removed from the tube.

Note: With the final wash, it is important to remove as much of the ethanol as possible. Use at least 2 pipetting steps and allow excess ethanol to collect at the bottom of the tubes after removing most of the ethanol in the first pipetting step.

11. Remove the tube from the magnet and air dry on the bench top for no more than 2 minutes. If the beads dry too long, they are difficult to resuspend.
12. Add 30 μL of room temperature, Nuclease-free Water (green: D1) to the tube.
13. Resuspend the beads by repeated pipetting. Alternatively, the beads may be resuspended by carefully vortexing the tube for 30 seconds or using a plate shaker set to medium speed. Ensure the beads are fully resuspended. Vortex longer if necessary.
14. Replace the tube on the magnet and allow the beads to separate for 5 minutes or until the solution clears.
15. Carefully remove the eluted sample and transfer to a fresh tube.
Note: Small amounts of bead carryover may interfere with sample quantification. Take care to minimize bead carryover.
16. Continue with the Measuring SPIA cDNA Yield and Purity protocol or store the purified SPIA cDNA at -20°C .



Use nuclease-free water at room temperature to elute sample.

VI. Appendix

QIAGEN MinElute Reaction Cleanup Kit (instructions for a single reaction)

Important notes:

- Buffer ERC is considered hazardous according to QIAGEN, and an MSDS may be consulted.
- Add the appropriate amount of 100% ethanol to Buffer PE before use (see bottle label for volume).
- All centrifuge steps are carried out at maximum speed in a conventional table-top microcentrifuge at room temperature.

1. Into a clean, labeled 1.5 mL microcentrifuge tube, add 300 μ L of Buffer ERC from the QIAGEN kit.
2. Add the entire volume (40 μ L) of the SPIA reaction to the tube.
3. Vortex for 5 seconds, then spin briefly.
4. Obtain and label a MinElute spin column and place it into a collection tube.
5. Load the sample/buffer mixture onto the column.
6. Centrifuge for 1 minute at maximum speed in a microcentrifuge.
7. Discard the flow-through and replace the column in the same collection tube.
8. Add 750 μ L of Buffer PE to the column.
9. Centrifuge for 1 minute at maximum speed.
10. Discard the flow-through and replace the column in the same collection tube.
11. Centrifuge the column for an additional 2 minutes at maximum speed to remove all residual Buffer PE.
Note: Residual ethanol from the wash buffer will not be completely removed unless the flow-through is discarded before this additional centrifugation.
12. Discard the flow-through with the collection tube. Blot the column onto clean, absorbant paper to remove any residual wash buffer from the tip of the column.
Note: Blotting the column tip prior to transferring it to a clean tube is necessary to prevent any wash buffer transferring to the eluted sample.
13. Place the column into a clean, labeled 1.5 mL microcentrifuge tube.
14. Add 20 μ L of room temperature, Nuclease-free Water (green: D1) from the Tecan Genomics® kit to the center of each column.
Note: Ensure that the water is dispensed directly onto the membrane for complete elution of the bound cDNA.
15. Let the column stand for 1 minute at room temperature.



100% ethanol must be added to the QIAGEN Buffer PE upon first use. Failure to do so will result in low amplification yields.



Use nuclease-free water at room temperature to elute sample

VI. Appendix

16. Centrifuge for 1 minute at maximum speed.
17. Discard the column and measure the volume recovered.
18. Mix the sample by vortexing, then spin briefly.
19. Continue with the Measuring SPIA cDNA Yield and Purity protocol or store purified SPIA cDNA at -20°C .

VI. Appendix

B. Guidelines for Use in Microarray Experiments

The SPIA cDNA produced by the Ovation PicoSL WTA System V2 may be labeled for use in microarray analysis. Tecan Genomics has developed specific labeling products and protocols for use with various microarray platforms. Workflows suitable for use with specific microarray platforms are described below.

Affymetrix GeneChip Arrays

The Ovation PicoSL WTA System V2 is suitable for use only with mini (169 format) GeneChip Arrays, GeneChip Array Strips and Array Plates due to restrictions on cDNA yield. The Encore Biotin Module (Part No. 4200) is designed to prepare the SPIA cDNA produced with the Ovation PicoSL WTA System V2 for hybridization to these arrays. To label the cDNA it will be necessary to carry out the Encore Biotin Module protocol using half the normal volumes given in the user guide and a 2.5 µg input of SPIA cDNA. The SPIA cDNA sample must have a concentration of at least 200 ng/µL for use in the Encore Biotin Module half reaction since there is only 12.5 µL for sample input. For this reason we strongly recommend using the QIAGEN MinElute method given in this user guide to avoid the necessity to concentrate the SPIA cDNA prior to labeling.

Refer to Appendix C for more information about the use of the Ovation PicoSL WTA System V2 with GeneChip Array Strips and Array Plates.

Illumina Genome-Wide Expression BeadChips

The Encore BiotinIL Module (Part No. 4210) is designed to prepare the SPIA cDNA produced with the Ovation PicoSL WTA System V2 for hybridization to Illumina Genome-Wide Expression BeadChips. The Encore BiotinIL Module user guide provides an optimized protocol for use with the Ovation PicoSL WTA System V2. 0.75-1.5 µg of SPIA cDNA is required for labeling with this method, depending on the specific array format.

Agilent Dual-mode Gene Expression Arrays

We have developed a protocol for the fluorescent labeling of SPIA cDNA for use on Agilent Dual-mode Gene Expression Arrays, which is described in the Agilent Solution technical report. This document may be found at www.nugen.com/products/ovation-picosl-wta-system-v2. 2 µg of SPIA cDNA is required for labeling with this method.

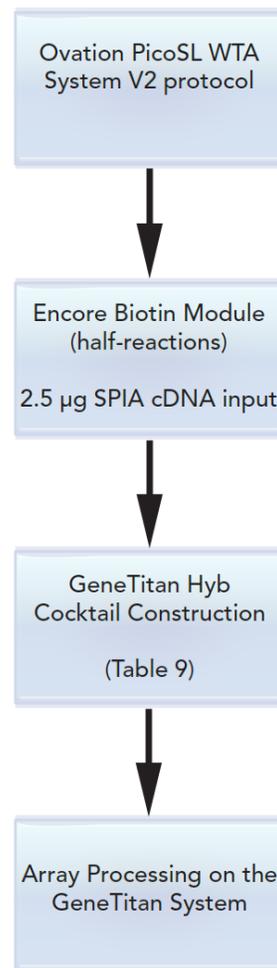
VI. Appendix

C. Use with Affymetrix® GeneTitan® Systems

The GeneTitan from Affymetrix is a family of instruments designed to enable the automated processing of GeneChip® Array Plates. The Ovation PicoSL WTA System V2 can be used with the Encore Biotin Module (Part No. 4200) as part of an integrated workflow with a GeneTitan System in gene expression applications.

Figure 3 shows the workflow recommended when using the Ovation PicoSL WTA System V2 on a GeneTitan System. The Ovation PicoSL WTA System V2 protocol is carried out as described in this user guide. After appropriate QC analysis, 2.5 µg of the SPIA cDNA is biotin labeled using a half-scale Encore Biotin Module protocol. (Follow the protocol given in the Encore Biotin Module user guide, but use only one-half the recommended volumes.)

Figure 3. Ovation PicoSL WTA System V2 workflow with GeneTitan Systems.



VI. Appendix

Processing GeneChip WT Array Plates on the GeneTitan

The Hybridization Cocktail construction for use with Affymetrix WT Array Plates is described below. The total volume of the cocktail is 100 μL and 90 μL of that volume is then used for each array.

1. Make a Hybridization Master Mix by combining the first 5 components in Table 9 below. Volumes for 16-, 24- and 96-Array Plates have been provided. These volumes include a 10% overage to allow for pipetting error and handling loss.
2. Aliquot 35 μL of the Hybridization Master Mix per sample.
3. Add 25 μL of the biotin labeled SPIA cDNA (2.5 μg) to each Hybridization Master Mix Aliquot.
4. Add 40 μL of the 25X WT Hyb Add 6 to each mixture to complete the Hybridization Cocktail. The total volume of each Hybridization Cocktail is 100 μL .
5. Mix the Hybridization Cocktail by vortexing for 5 seconds, then spin briefly.

Table 9. GeneTitan Hybridization Cocktail Formulation for WT Array Plates.

Add Order	Component	Volume per Array	16-Array Plate*	24-Array Plate*	96-Array Plate*	Final Conc.
1	5X WT Hyb Add 1	20 μL	352 μL	528 μL	2112 μL	1X
2	Control Oligonucleotide B2 (3nM)	1.0 μL	17.6 μL	26.4 μL	105.6 μL	30 pM
3	20X Eukaryotic Hybridization Controls (bioB, bioC, bioD, cre)	5 μL	88 μL	132 μL	528 μL	1.5, 5, 25 and 100 pM
4	15X WT Hyb Add 4	6.7 μL	117.9 μL	176.9 μL	707.5 μL	1X
5	Nuclease-free Water	2.3 μL	40.5 μL	60.7 μL	242.9	
Total Volume		35 μL	616 μL	924 μL	3696 μL	
6	Labeled SPIA cDNA (2.5 μg)	25 μL				25 ng/ μL
7	25X WT Hyb Add 6	40 μL				1X
Total Volume		100 μL				

*10% overage has been included in these volumes.

VI. Appendix

6. Denature the Hybridization Cocktail by heating to 95°C if using PCR plates or 99°C if using microcentrifuge tubes.
7. Incubate at 45°C for 5 minutes.
8. Spin for 1 minute at 5,000 RPM to remove any insoluble material from the Hybridization Cocktail.
9. Place 90 µL of the Hybridization Cocktail into the appropriate well of the hybridization tray.
10. Follow the manufacturer's recommendations to process the Array Plate using a GeneTitan System.

Processing GeneChip 3' Plate Arrays on the GeneTitan

The Hybridization Cocktail construction for use with Affymetrix 3' Array Plates is described below. The total volume of the cocktail is 100 µL and 90 µL of that volume is then used for each array.

1. Make a Hybridization Master Mix by combining the first 4 components in Table 10 below. Volumes for 16-, 24- and 96-Array Plates have been provided. These volumes include a 10% overage to allow for pipetting error and handling loss.
2. Aliquot 83.6 µL of the Hybridization Master Mix per sample.
3. Add 16.4 µL of the biotin labeled SPIA cDNA (1.6 µg) to each Hybridization Master Mix. The total volume of each Hybridization Cocktail is 100 µL.
4. Mix the Hybridization Cocktail by vortexing for 5 seconds, then spin briefly.
5. Denature the Hybridization Cocktail by heating to 95°C if using PCR plates or 99°C if using microcentrifuge tubes.
6. Incubate at 45°C for 5 minutes.
7. Spin for 1 minute at 5,000 RPM to remove any insoluble material from the Hybridization Cocktail.

VI. Appendix

Table 10. GeneTitan Hybridization Cocktail Formulation for 3' Array Plates.

Add Order	Component	Volume per Array	16-Array Plate*	24-Array Plate*	96-Array Plate*	Final Conc.
1	Control Oligonucleotide B2 (3nM)	1.7 μ L	29.9 μ L	44.9 μ L	179.5 μ L	30 pM
2	20X Eukaryotic Hybridization Controls (bioB, bioC, bioD, cre)	5 μ L	88 μ L	132 μ L	528 μ L	1.5, 5, 25 and 100 pM
3	1.3X Hybridization Solution A	26.9 μ L	473.4 μ L	710.2 μ L	2841 μ L	1X
4	1.3X Hybridization Solution B	50 μ L	880 μ L	1320 μ L	5280 μ L	
Total Volume		83.6 μ L	1471 μ L	2207 μ L	8828 μ L	
5	Labeled SPIA cDNA (1.6 μ g)	16.4 μ L				16 ng/ μ L
Total Volume		100 μ L				

*10% overage has been included in these volumes.

8. Place 90 μ L of the Hybridization Cocktail into the appropriate well of the hybridization tray.
9. Follow the manufacturer's recommendations to process the Array Plate using a GeneTitan System.

VI. Appendix

D. Performing Quantitative PCR on SPIA cDNA

It is recommended that the amplified cDNA (SPIA cDNA) generated from the Ovation PicoSL WTA System V2 be purified prior to use in quantitative PCR reactions (qPCR). Since different SPIA cDNA samples will vary in concentration, the purified products may be quantified and mass normalized to ensure the cDNA input into the qPCR reaction is equal for all samples. Purified SPIA cDNA produced with this kit has been successfully used as template for qPCR systems, including TaqMan® and SYBR® Green. Note that RT-PCR master mixes containing the enzyme Uracil N-Glycosylase (UNG) are not compatible with the Ovation PicoSL WTA System V2.

We can recommend the following reagents for qPCR:

- TaqMan: Absolute qPCR Mix plus ROX (Thermo Fisher Scientific, Cat. #AB-1136/B), Fast Universal PCR Master Mix 2x (Thermo Fisher Scientific, Cat. #4352042)
- SYBR: QuantiTect™ SYBR Green PCR Kit (QIAGEN, Cat. #204143), iQ SYBR Green Supermix (BioRad, Cat. #170-8880), FastStart SYBR Green Master (ROX) (Roche, Cat. #04 673 514 001)

Recommendations to Achieve Optimal Results

1. Dilute the SPIA cDNA

After purification and quantification of the SPIA cDNA, it can be diluted to an appropriate concentration for qPCR reaction. We recommend using 20 ng of SPIA cDNA in a 20 µL TaqMan reaction and 2 ng of cDNA for a 25 µL SYBR Green reaction. Depending on the abundance of the transcripts of interest you may wish to use more or less cDNA.

2. Primer Design

We recommend using primers and probes designed with amplicon sizes of less than 200 nt. Primers may be designed at any position along a transcript since the Ovation PicoSL WTA System V2 amplification covers the whole transcriptome.

VI. Appendix

E. Quality Control of the SPIA cDNA Product

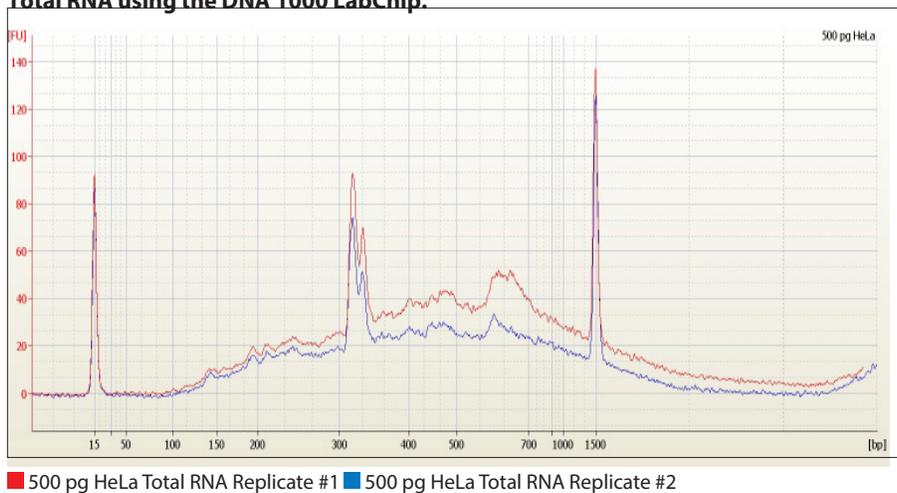
There are several methods which we can recommend for quality control analysis of the SPIA cDNA including spectrophotometric quantification of amplification yield (as described in Section IV, Protocol I on page 19) and fragment size distribution analysis using the Agilent Bioanalyzer.

When performing fragment size distribution analysis on the Bioanalyzer, we recommend using either the Agilent DNA 1000 LabChip (Agilent Cat. #5067-1504) or the RNA 6000 Nano LabChip (Agilent Cat. #5065-4476) following the manufacturer's instructions and the specific guidelines given below.

Guidelines for using the Agilent DNA 1000 LabChip

When using the DNA 1000 LabChip (Agilent Cat. # 5065-1504) for analysis of SPIA cDNA fragment distribution, use the DNA 1000 Series II program and follow the manufacturer's instructions. A typical size distribution trace may look like the one obtained from HeLa Total RNA (Figure 4). Note that the shape of this distribution trace is highly dependent on the input RNA integrity as well as RNA source, and can vary significantly.

Figure 4. Bioanalyzer trace of SPIA cDNA product obtained from 500 pg of HeLa Total RNA using the DNA 1000 LabChip.

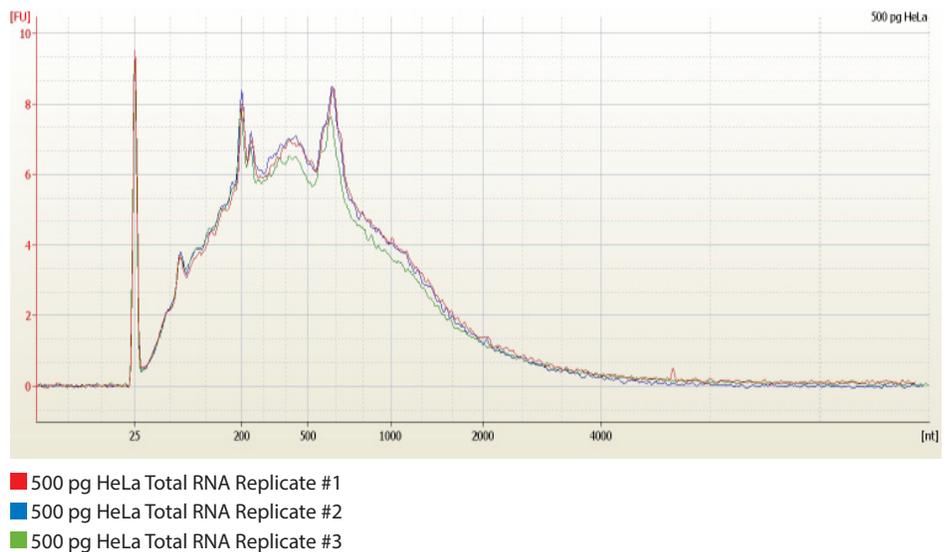


VI. Appendix

Guidelines for using the Agilent RNA 6000 Nano LabChip

When using the RNA 6000 Nano LabChip (Agilent Cat. #5065-4476) for analysis of SPIA cDNA fragment distribution, use the Eukaryotic Total RNA Nano program and follow the manufacturer's instructions. Denaturation of the SPIA cDNA sample (as described in the Agilent RNA 6000 Nano protocol) is required for optimum resolution and reproducibility using this option. A typical size distribution trace may look like the one obtained from HeLa Total RNA (Figure 5). Note that the shape of this distribution trace is highly dependent on the input RNA integrity as well as RNA source, and can vary significantly.

Figure 5. Bioanalyzer trace of SPIA cDNA product obtained from 500 pg of HeLa Total RNA using the RNA 6000 Nano LabChip.



VI. Appendix

F. DNase Treatment of RNA

DNase Treatment During Purification: Using the QIAGEN RNase-free DNase Set and the RNeasy Mini RNA Purification Kit.

1. Homogenize sample in Buffer RLT including β -mercaptoethanol according to the type of sample as described in the RNeasy Mini Kit protocol.
2. Add 1X volume of 70% ethanol to the homogenized lysate, pipet up and down to mix sample well. Do not centrifuge.
3. Place an RNeasy mini column in a 2 mL collection tube.
4. Apply the sample (up to 700 μ L), including any precipitate that may have formed, to the column.
5. Close the tube gently and centrifuge for 15 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
6. For volumes greater than 700 μ L, load aliquots onto the RNeasy column successively and centrifuge as before.
7. Add 350 μ L Buffer RW1 into the RNeasy mini column to wash, and centrifuge for 15 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
8. Add 10 μ L DNase I to 70 μ L Buffer RDD. Gently invert the tube to mix.
9. Pipet the DNase I incubation mix (80 μ L) directly onto the membrane inside the RNeasy mini column. Incubate on the bench top ($\sim 25^\circ\text{C}$) for 15 min.
10. Add 350 μ L Buffer RW1 into the RNeasy mini column, and centrifuge for 15 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm) to wash. Discard the flow-through.
11. Transfer the RNeasy column to a fresh 2 mL collection tube. Add 500 μ L Buffer RPE (with the added ethanol) to the RNeasy column.
12. Close the tube gently and centrifuge for 15 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
13. Add another 500 μ L Buffer RPE to the RNeasy column.
14. Close the tube gently and centrifuge for 2 minutes at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
15. Transfer the RNeasy column to a new 1.5 mL collection tube.
16. Pipet 30 to 50 μ L RNase-free water directly onto the RNeasy membrane.
17. Close the tube gently and centrifuge for 1 minute at $\geq 8000 \times g$ ($\geq 10,000$ rpm) to elute.
18. If yields of greater than 30 μ g are expected, repeat elution step and collect in the same collection tube.

VI. Appendix

DNase Treatment of RNA Post-purification: Using RNase-free DNase and either the RNA Clean & Concentrator™-5 Columns or the RNeasy MinElute Columns

Note: If you are unable to quantify your RNA because the sample is contaminated with DNA, we recommend DNase treatment followed by purification.

1. On ice, mix together 2.5 µL 10 X DNase I Reaction buffer (Roche Cat. #04716728001) with 1 µL rDNase (10 Units Roche Cat. #04716728001).
2. Add RNA sample (up to 500 ng) and add RNase-free water (D1, green cap) to bring the final volume to 25 µL.
3. Incubate at 25°C for 15 minutes followed by 37°C for 15 minutes and return to ice.
4. After the DNase treatment, the sample must be purified. We recommend either of the two purification procedures below:

Purification with RNA Clean & Concentrator-5 (Zymo Research, Cat. #R1015)

1. Add 4 volumes (100 µL) of RNA binding buffer to the sample.
2. Obtain one RNA Clean & Concentrator Kit-5 column and apply sample to column.
3. Spin column for 30 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
4. Add 200 µL wash buffer (with ethanol added as per vendor's specifications).
5. After closing the column spin for 30 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
6. Add 200 µL fresh 80% ethanol, close cap and spin for 30 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
7. Place the RNA Clean & Concentrator Kit-5 column in a fresh 1.5 mL collection tube.
8. Add 10 µL nuclease-free water (green: D1) directly to the center of the filter in the tube and close the cap.

Important: Do not use cold water!
9. Spin for 1 minute at $\geq 8000 \times g$ ($\geq 10,000$ rpm) to collect the purified RNA.

Purification with QIAGEN RNeasy MinElute Cleanup Columns (QIAGEN, Cat. #74204)

1. Add 80 µL ice-cold RNase-free water (D1, green cap) to the sample on ice.
2. Add 350 µL Buffer RLT and mix by pipetting.
3. Add 250 µL 96-100% ethanol and mix thoroughly by pipetting.



Use nuclease-free water at room temperature to elute sample.

VI. Appendix



Best results can be obtained by using fresh 80% ethanol in the wash step. Lower percent ethanol mixes will reduce recovery.



Use nuclease-free water at room temperature to elute sample.

4. Place an RNeasy MinElute Spin Column into a 2 mL collection tube (one column per sample) and apply the 700 μ L sample to the column.
5. After closing the column, spin for 15 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
6. Place the RNeasy MinElute Spin Column into a fresh 2 mL collection tube. Add 500 μ L Buffer RPE to the column and close the tube. Spin for 15 seconds at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through, keeping the same collection tube.
7. Add 500 μ L 80% ethanol to the RNeasy MinElute Spin Column and close the tube.
Note: Use fresh 80% ethanol. Lower percent ethanol mixes will reduce recovery.
8. Spin for 2 minutes at $\geq 8000 \times g$ ($\geq 10,000$ rpm). Discard the flow-through.
9. Place the RNeasy MinElute Spin Column in a fresh 2 mL collection tube and place in the microcentrifuge with the cap open. Spin for 5 minutes at $\geq 8000 \times g$ ($\geq 10,000$ rpm) and discard the flow-through.
10. Place the RNeasy MinElute Spin Column in a fresh 1.5 mL collection tube.
11. Add 14 μ L nuclease-free water (D1, green cap) directly to the center of the filter in the tube and close the cap. Do not use cold water!
12. Spin for 1 minute at $\geq 8000 \times g$ ($\geq 10,000$ rpm) to collect the purified RNA.

VI. Appendix

G. Preventing Non-specific Amplification

Due to the high sensitivity inherent in our amplification systems, we have developed a set of recommendations designed to minimize the potential for the generation of non-specific amplification products through the carry-over of previously amplified SPIA cDNA. We strongly recommend implementing these procedures, especially for the high-throughput and low-RNA input environments typical in today's gene expression laboratories. We have two general recommendations. First, designate separate workspaces for "pre-amplification" and "post-amplification" steps and materials. This provides the best work environment for processing RNA using our highly sensitive amplification protocols. Our second recommendation is to implement routine clean-up protocols for workspaces as standard operating procedure. This will prevent non-specific amplification products from spreading through the laboratory. Details regarding establishing and maintaining a suitable work environment are listed below:

1. Designate a pre-amplification workspace separate from the post-amplification workspace or general lab areas:
 - a. Pre-amplification includes all steps and materials related to RNA sample handling and dilution, Tecan Genomics' first strand reaction, second strand reaction, second strand cleanup and SPIA amplification reaction setup. After SPIA incubation the reactions are immediately removed from the pre-amplification workspace and opened only in the post-amplification area.
 - b. Post-amplification includes all steps and materials related to the handling of the final amplified cDNA product including bead removal, final purification, post-SPIA modification, array hybridization and any other analytical work.
 - c. Ideally the pre-amplification workspace will be in a separate room. If this is not possible, ensure the pre-amplification area is sufficiently isolated from post-amplification work.
 - d. PCR Workstation enclosures with UV illumination for use as pre-amplification workspaces can be an option in situations where conditions preclude physical separation of pre- and post-amplification activities.
2. Establish and maintain a clean work environment:
 - a. Initially clean the entire lab thoroughly with DNA-OFF. Follow this treatment with a thorough rinse with water to ensure no residual cleaning agents are left behind.
 - b. In the pre-amplification area, remove all small equipment, and then clean every surface that may have been exposed to amplified SPIA cDNA (surfaces, drawer handles, key pads, etc.). Before reintroducing any equipment, clean every piece of equipment thoroughly.

Clean thermal cycler blocks by heating to 99°C for 15 minutes, then wipe down exposed surfaces and keypad with cleaning solution.

Clean magnets by immersion in cleaning solution or use a cotton swab.

VI. Appendix

- c. Carry out a thorough external and internal cleaning of all pipettes with DNA-OFF. Carefully follow the manufacturer's instructions for this process to avoid damaging the pipettes. It is a good idea to keep a clean set of pipettes as a backup.
 - d. Always wear gloves and don fresh gloves upon entry into this controlled area. Frequently change gloves while working in the pre-amplification area, especially prior to handling stock reagents, reactions and RNA samples.
 - e. Stock this area with clean (preferably new) equipment (pipettes, racks, consumables) that has not been exposed to post-amplification workspace.
 - f. Make it a policy to carry out regular cleaning of all workspaces.
 - g. Capture waste generated in both pre- and post-amplification areas (tips, columns, wash solutions from beads and columns, tubes, everything) in sealable plastic bags and dispose of promptly after each experiment to avoid waste spillage.
 - h. Do not open amplified product reaction vessels in the pre-amplification workspace.
3. Avoid running negative controls (i.e., no RNA input reactions). Instead use low-template controls (inputs of 50 pg to 100 pg) in order to detect and monitor any non-specific amplification issues. The clearest indication that non-specific amplification is taking place is the appearance of higher than expected yields or irregular Bioanalyzer traces in a low template control (LTC) reaction.
 - a. Typical amplification performance:
 - i. LTC yields for Ovation PicoSL WTA System V2 amplifications should be significantly lower than yields for RNA inputs within the recommended input range of 500 pg to 50 ng.
 - ii. The Bioanalyzer trace of the LTC amplification product is consistent with that seen with higher input.
 - b. Atypical amplification performance:
 - i. LTC yields may be similar to those obtained using higher inputs of total RNA.
 - ii. The Bioanalyzer traces of amplification products may look significantly different than the typical Ovation PicoSL WTA reaction traces. The LTC reaction is designed to be an especially sensitive indicator of atypical amplification performance.
 - iii. Sensitivity on arrays may be lower than expected.
 - iv. Contact Tecan Technical Support when atypical performance is suspected.

H. Frequently Asked Questions (FAQs)

Q1. What materials are provided with the Ovation PicoSL WTA System V2?

The Ovation PicoSL WTA System V2 provides all necessary buffers, primers and enzymes for first strand synthesis, second strand synthesis and amplification. The kit also provides nuclease-free water and Agencourt Beads for double-stranded cDNA purification.

Q2. Does the Ovation PicoSL WTA System V2 provide any labeling reagents?

No. The Ovation PicoSL WTA System V2 is used to generate cDNA from small amounts of total RNA for qPCR analysis or cDNA storage; however, the cDNA output of this kit may be processed further using other Tecan Genomics products such as the Encore labeling modules (for labeling cDNA for analysis on Affymetrix GeneChip or Illumina BeadChip arrays) or other supported labeling protocols.

Q3. How much of the SPIA cDNA from the Ovation PicoSL WTA System V2 should I use for array labeling?

We recommend following the guidelines given for the original Ovation PicoSL WTA System (Part No. 3310) for use in an Encore labeling module or other supported labeling protocol. Refer to the Encore labeling module user guide or labeling protocol for specific guidance.

Q4. Will I need to use the WT-Ovation Exon Module when the cDNA will be labeled for use on an Affymetrix GeneChip Gene ST array?

No. The cDNA produced by the Ovation PicoSL WTA System V2 is suitable for labeling and use with both Affymetrix GeneChip Gene ST Arrays as well as certain GeneChip 3' Arrays without further processing.

Q5. What equipment is required or will be useful?

Required equipment includes a microcentrifuge, pipettes, vortexer, a thermal cycler, a spectrophotometer and a magnetic plate. An Agilent Bioanalyzer may also be useful for optional analytical tests.

Q6. What additional consumables does the user need?

For the SPIA cDNA purification step, purification columns or beads are required.

Q7. Can I do reactions in smaller batches than four?

We recommend a minimum batch size of four reactions. Smaller batch sizes may result in difficulty with pipetting small volumes, as well as obtaining fewer than 12 reactions in total.

Q8. Has Tecan Genomics performed reproducibility studies on the Ovation PicoSL WTA System V2?

Yes. Sample-to-sample, lot-to-lot and operator-to-operator reproducibility tests are conducted.

Q9. Is the Ovation PicoSL WTA System V2 3' biased?

In this system, oligo dT primers are mixed with random primers for the first strand synthesis of cDNA products. This allows the product to be analyzed on 3' expression arrays and whole transcriptome arrays when used with an appropriate Encore labeling module, or other supported labeling protocol. Additionally, the random primers allow the detection of the entire transcripts when used as a pre-qPCR amplification system.

Q10. How much total RNA do I need for amplification?

We recommend staying within the range of 500 pg to 50 ng total RNA as starting material. Input amounts outside this range may produce unsatisfactory and variable results.

Q11. Do I need to use high-quality total RNA?

RNA samples of high molecular weight with little or no evidence of degradation, as expected, will amplify very well with this product. However, due to the whole transcriptome amplification approach, lower quality RNA samples and transcripts with a compromised poly(A) tail can also be amplified successfully using the Ovation PicoSL WTA System V2. The RNA should have high purity, however, and be free of contaminants.

Q12. Can contaminating genomic DNA interfere with the amplification performance?

This system is designed to amplify RNA but large amounts of contaminating genomic DNA may amplify during the process. For this reason we recommend DNase treatment during RNA purification.

Q13. Do you recommend DNase treatment of my total RNA sample?

Yes. For an explanation of DNase requirements see section III.A.5. You may also find recommended procedures for DNase treatment in Appendix D.

Q14. Can I use the Ovation PicoSL WTA System V2 on bacterial RNA samples?

The Ovation PicoSL WTA System V2 amplification process has been shown to work with some bacterial RNAs. However, the kit has not been optimized for this purpose.

Q15. Are there any tissues that will not work with the Ovation PicoSL WTA System V2?

We have not encountered any specific RNA sources that will not work with the Ovation PicoSL WTA System V2. The RNA should have high purity and be free of contaminants.

Q16. How much cDNA can I expect from a single reaction?

You should expect 2–4 µg of cDNA from input of 500 pg to 50 ng total RNA.

Q17. Is the cDNA yield dependent upon the quantity of total RNA input?

Yes, higher RNA inputs will produce higher yields. However, at inputs of above 50 ng, the yields may become variable without increasing.

Q18. What size cDNA is generated by the Ovation PicoSL WTA System V2?

The SPIA cDNA size distribution is somewhat dependent on the input RNA integrity. A representative size distribution is given in Figure 4.

Q19. Does the Ovation PicoSL WTA System V2 generate product in the absence of RNA input?

In the complete absence of input RNA, approximately 0.5 µg or less of non-specific product is generated. However, in the presence of even a very small amount of RNA the amplified cDNA has been demonstrated to be specific.

Q20. How many rounds of amplification are performed with the Ovation PicoSL WTA System V2?

This System has a single round of amplification. It cannot be used for multiple rounds.

Q21. Can I use the Ovation PicoSL WTA System V2 for archiving cDNA?

SPIA cDNA may be stored at -20°C for at least six months.

Q22. Do I have to use your DNA/RNA primers, or can I order specific primers for the amplification?

The Ovation PicoSL WTA System V2 will not work properly with other primers. There is no need to order any primers, as DNA/RNA primers provided in the Ovation PicoSL WTA System V2 are universal.

Q23. What purification methods do you recommend?

- For the Second Strand cDNA purification step (pre-amplification) we require the use of the Agencourt Beads provided with the kit.
- Several purification options are available for the final SPIA cDNA cleanup step. These are described in Appendix A of this user guide. Selection of the optimum purification option can depend on many factors. Please contact the Tecan Technical Support team for assistance in selecting the appropriate option for your application. Refer to section II.B. for ordering information.

Q24. How do I measure my SPIA cDNA product yield?

You may use a standard spectrophotometer or a Nanodrop. Refer to Section IV, Protocol I for specific guidance.

Q25. Where can I safely stop in the protocol?

The SPIA cDNA can be stored at -20°C prior to performing the purification. We do not recommend stopping at any intermediate stage of the protocol.

Q26. How many qPCR reactions will I get from one Ovation PicoSL WTA System V2 amplification?

The number of qPCR reactions depends on the abundance level of the genes being interrogated. For medium- to high-copy genes, the cDNA may be diluted as much as 400-fold, enough for thousands of qPCR reactions. For very-low-copy genes you will need to use more cDNA per reaction. The user will need to determine how much cDNA to use per reaction depending on the abundance of the gene being interrogated. Note that we recommend purification of the SPIA cDNA prior to qPCR analysis.

VI. Appendix

Q27. Do you recommend purification of the cDNA prior to qPCR analysis?

Yes. Although this is not absolutely necessary, it is important to be able to quantify the SPIA cDNA. This allows assessment of amplification success based on the amplification yields. It also allows mass normalization of the cDNA into qPCR.

Q28. Where in my target sequence can I design my qPCR primers?

The Ovation PicoSL WTA System V2 does not have a 3' bias and, therefore, primers can be designed at any location within the mRNA. In order to avoid qPCR interference from possible genomic DNA contamination, we recommend treating your RNA with DNase and designing your amplicons to span an intron.

VI. Appendix

I. Update History

This document, the Ovation Pico WTA System V2 user guide (M01224 v6), is an update to address the following topics:

Description	Section	Page(s)
Update Technical support contact	Throughout	Throughout
NuGEN to Tecan Genomics	Throughout	Throughout



Tecan Genomics Inc.

Headquarters USA

900 Chesapeake Dr.
Redwood City, CA 94063 USA
Toll Free Tel: 888.654.6544
Toll Free Fax: 888.296.6544
techserv-gn@tecan.com

Europe

P.O. Box 109
9350 AC Leek
The Netherlands
Tel: +31-13-5780215
Fax: +31-13-5780216
europe-gn@tecan.com

For our international distributors contact
information, visit our website

www.nugen.com

©2019 Tecan Genomics, Inc. All rights reserved. The Encore®, Ovation® and Applause® families of products and methods of their use are covered by several issued U.S. and International patents and pending applications (www.nugen.com). NuGEN, Allegro, Celero, NuQuant, SoLo, Metaplex, DimerFree, AnyDeplete, Ovation, SPIA, Ribo-SPIA, Applause, Encore and Imagine More From Less are trademarks or registered trademarks of Tecan Genomics, Inc. Other marks appearing in these materials are marks of their respective owners.

