













This product contains components with multiple storage conditions



Store the First Strand Primer Mix, First Strand Enzyme Mix, and SPIA Primer Mix at  $-80^{\circ}\text{C}$  and the Agencourt Beads at  $4^{\circ}\text{C}$

### 1 FSGPSNBODF4QD

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  $\mu\text{g}$  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.

### 2 VBM

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

### 4 UPSBHFBOE4UBC

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  $4^{\circ}\text{C}$ .

The vial labeled First Strand Enzyme Mix (blue: A3) should be removed from the kit upon delivery and stored separately at  $-80^{\circ}\text{C}$ .

All remaining components should be stored at  $-20^{\circ}\text{C}$  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.

### 4 BGFUZ

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

## 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](http://www.beckmangenomics.com)
- MP Biomedicals, [www.mpbio.com](http://www.mpbio.com)
- New England BioLabs, [www.neb.com](http://www.neb.com)
- QIAGEN Inc., [www.qiagen.com](http://www.qiagen.com)
- Sigma-Aldrich, Inc., [www.sigmaaldrich.com](http://www.sigmaaldrich.com)
- Sorenson Biosciences, Inc., [www.sorbio.com](http://www.sorbio.com)
- Thermo Fisher Scientific, [www.thermofisher.com](http://www.thermofisher.com)
- Zymo Research, [www.zymoresearch.com](http://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.

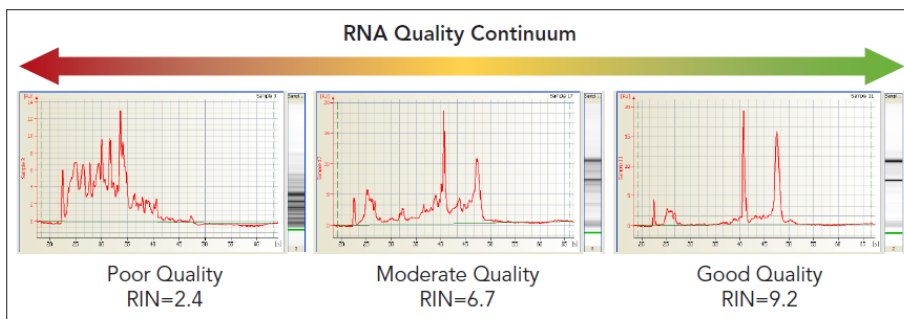
#### 3. RNA Integrity

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  $-80^{\circ}\text{C}$ . Avoid frequent freeze/thaw cycles or RNA degradation may result.

### D. SPIA<sup>®</sup> cDNA Storage

The SPIA cDNA produced by the Ovation PicoSL WTA System V2 may be stored at  $-20^{\circ}\text{C}$ .

## 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:

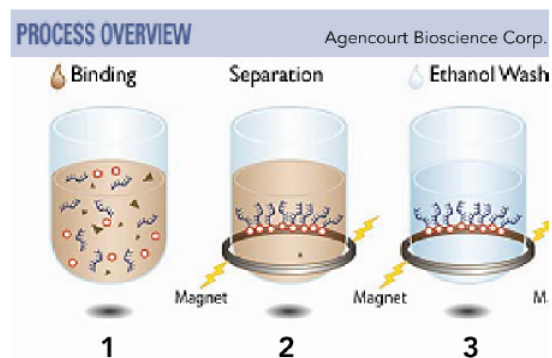
There are significant modifications to the Beckman Coulter Genomics' Agencourt SPRI Bead Kit standard procedure; therefore, you must follow the procedures outlined in this user guide for the use of these beads with the Ovation PicoSL WTA System V2.

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 µL) 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



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  $-20^{\circ}\text{C}$  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 $\mu\text{L}$	10 $\mu\text{L}$	10 $\mu\text{L}$

5. Add 40  $\mu\text{L}$  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](mailto: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  $-20^{\circ}\text{C}$  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 A<sub>320</sub> value from both A<sub>260</sub> and A<sub>280</sub> values. The adjusted  $(A_{260} - A_{320}) / (A_{280} - A_{320})$  ratio should be  $>1.8$ .
4. Yield: Assume 1 A<sub>260</sub> 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  $-20^{\circ}\text{C}$ .

## 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](mailto: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](mailto: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

### 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  $\mu\text{L}$  and 90  $\mu\text{L}$  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  $\mu\text{L}$  of the Hybridization Master Mix per sample.
3. Add 25  $\mu\text{L}$  of the biotin labeled SPIA cDNA (2.5  $\mu\text{g}$ ) to each Hybridization Master Mix Aliquot.
4. Add 40  $\mu\text{L}$  of the 25X WT Hyb Add 6 to each mixture to complete the Hybridization Cocktail. The total volume of each Hybridization Cocktail is 100  $\mu\text{L}$ .
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 $\mu\text{L}$	352 $\mu\text{L}$	528 $\mu\text{L}$	2112 $\mu\text{L}$	1X
2	Control Oligonucleotide B2 (3nM)	1.0 $\mu\text{L}$	17.6 $\mu\text{L}$	26.4 $\mu\text{L}$	105.6 $\mu\text{L}$	30 pM
3	20X Eukaryotic Hybridization Controls (bioB, bioC, bioD, cre)	5 $\mu\text{L}$	88 $\mu\text{L}$	132 $\mu\text{L}$	528 $\mu\text{L}$	1.5, 5, 25 and 100 pM
4	15X WT Hyb Add 4	6.7 $\mu\text{L}$	117.9 $\mu\text{L}$	176.9 $\mu\text{L}$	707.5 $\mu\text{L}$	1X
5	Nuclease-free Water	2.3 $\mu\text{L}$	40.5 $\mu\text{L}$	60.7 $\mu\text{L}$	242.9	
Total Volume		35 $\mu\text{L}$	616 $\mu\text{L}$	924 $\mu\text{L}$	3696 $\mu\text{L}$	
6	Labeled SPIA cDNA (2.5 $\mu\text{g}$ )	25 $\mu\text{L}$				25 ng/ $\mu\text{L}$
7	25X WT Hyb Add 6	40 $\mu\text{L}$				1X
Total Volume		100 $\mu\text{L}$				

\*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 µL	29.9 µL	44.9 µL	179.5 µL	30 pM
2	20X Eukaryotic Hybridization Controls (bioB, bioC, bioD, cre)	5 µL	88 µL	132 µL	528 µL	1.5, 5, 25 and 100 pM
3	1.3X Hybridization Solution A	26.9 µL	473.4 µL	710.2 µL	2841 µL	1X
4	1.3X Hybridization Solution B	50 µL	880 µL	1320 µL	5280 µL	
Total Volume		83.6 µL	1471 µL	2207 µL	8828 µL	
5	Labeled SPIA cDNA (1.6 µg)	16.4 µL				16 ng/µL
Total Volume		100 µL				

\*10% overage has been included in these volumes.

8. Place 90 µL 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

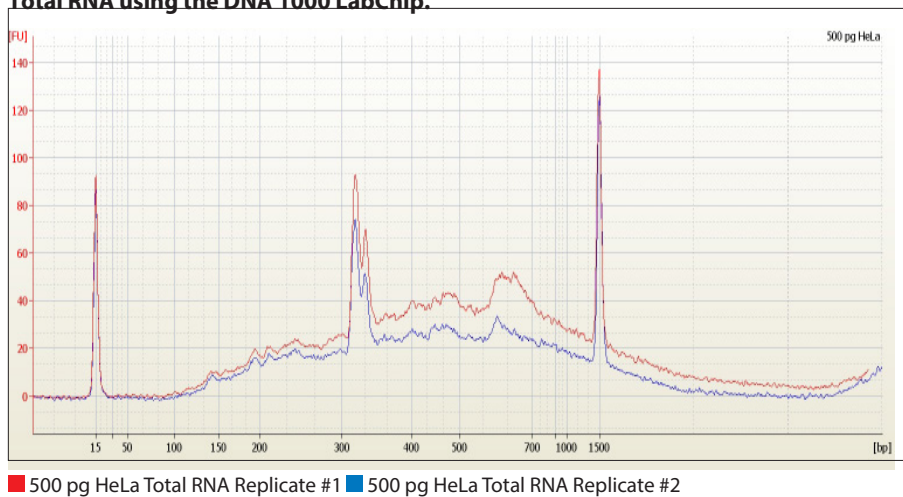
There are several methods which we can recommend for quality control analysis of the SPIA cDNA including spectrophotometric quantification of amplification yield (as described in Section IV, Protocol I on page 19) and fragment size distribution analysis using the Agilent Bioanalyzer.

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

- 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 Services 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.

## VI. Appendix

- 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.

## VI. Appendix

- 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.

##### 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](http://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](http://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.

