



QUICK PROTOCOL

Ovation[®] RNA-Seq System V2

Part No. 7102

Enter the number of reactions you are running in the provided field to automatically calculate the volumes needed to prepare each master mix. The calculated volume includes an appropriate overfill in excess of the nominal volume requirements (typically 10%) to allow for loss due to handling. Simply print this document to create a working guide for your experiment, which can be kept as a record.

Operator's Name:

Date:

Kit Part No: 7102-

Kit Lot No:

Number of Samples*:

THERMAL CYCLER PROGRAMS

First Strand cDNA Synthesis

Program 1: First Strand Primer Annealing

(For RNA inputs ≤ 1 ng) 65 °C - 2 min, hold at 4 °C
(For RNA inputs > 1 ng) 65 °C - 5 min, hold at 4 °C

Program 2: First Strand Synthesis

4 °C - 1 min, 25 °C - 10 min, 42 °C - 10 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 - 60 min, 80 °C - 20 min, hold at 4 °C

*Number of samples field ties into embedded logic to calculate suggested master mix volumes.

Tecan recommends processing a minimum of 4 samples at a time.



FIRST STRAND cDNA SYNTHESIS

Thaw the **First Strand cDNA Synthesis reagents (blue)** and **Nuclease-free Water (green)**.

Spin **A3 VER 7** briefly and place on ice. Vortex **A1 VER 4** and **A2 VER 3**, spin and place on ice.

Leave **Nuclease-free Water** at room temperature.

On ice, mix 2 µL of **A1** and 5 µL of total RNA sample (500 pg to 100 ng) in a 0.2 mL PCR tube.

Place the tubes in a thermal cycler running Program 1 (65 °C - 2 min, hold at 4 °C or 65 °C - 5 min, hold at 4 °C).

Once the thermal cycler reaches 4 °C, remove tubes and place on ice.

Prepare **First Strand Master Mix** (calculated volumes allow for appropriate overfill). Be sure to pipet **A3** enzyme slowly and rinse out tip at least five times into buffer.

Per sample combine:

2.5 µL Buffer Mix **A2** + 0.5 µL Enzyme Mix **A3**.

Mix well.

No. of Samples	A2	A3
1	2.5 µL	0.5 µL

Add 3 µL of **First Strand Master Mix** to each tube, mix by pipetting, spin and place on ice.

Place the tubes in a thermal cycler running Program 2 (4 °C - 1 min, 25 °C - 10 min, 42 °C - 10 min, 70 °C - 15 min, hold at 4 °C).

Once the thermal cycler reaches 4 °C, remove tubes, spin and place on ice.

Continue immediately with Second Strand cDNA Synthesis.

SECOND STRAND cDNA SYNTHESIS

Resuspend the Agencourt® beads provided with the Ovation RNA-Seq System V2 and leave at room temperature for use in the next step.

Thaw the **Second Strand cDNA Synthesis reagents (yellow)**.

Spin **B2 VER 2** briefly and place on ice. Vortex **B1 VER 3**, spin and place on ice.

Prepare **Second Strand Master Mix**. Be sure to pipet **B2** enzyme slowly.

Per sample combine:

9.7 µL Buffer Mix **B1** + 0.3 µL Enzyme Mix **B2**.

Mix well.

No. of Samples	B1	B2
1	9.7 µL	0.3 µL

Add 10 µL of **Second Strand Master Mix** to each reaction tube, mix by pipetting, spin and place on ice.

Place the tubes in a thermal cycler running Program 3 (4 °C - 1 min, 25 °C - 10 min, 50 °C - 30 min, 80 °C - 20 min, hold at 4 °C).

Once the thermal cycler reaches 4 °C, remove tubes, spin and place on bench top.

Continue immediately with Purification of cDNA.

PURIFICATION OF cDNA

Ensure the Agencourt beads have reached room temperature.	
Mix the beads by inverting several times.	
At room temperature, add 32 µL of Agencourt beads to each reaction tube and mix by pipetting 10 times.	
Incubate at room temperature for 10 minutes.	
Transfer the tubes to the magnet and let stand for an additional 5 minutes	
Remove only 45 µL of the binding buffer.	
Add 200 µL of freshly prepared 70% ethanol and let stand for 30 seconds. Remove the ethanol using a pipette.	
Repeat the ethanol wash 2 more times.	
Remove all excess ethanol after the final wash and let beads air dry for 15 to 20 minutes.	
Ensure the tubes have completely dried and no residual ethanol is left.	
Continue immediately with SPIA Amplification, with the cDNA bound to the dry beads.	

SPIA AMPLIFICATION

Thaw the SPIA Amplification Reagents (red) .													
Invert C3 VER 7 5 times to mix, spin and place on ice. Vortex C1 VER 9 and C2 VER 11 , spin and place on ice.													
Prepare SPIA Master Mix . Per sample combine: 20 µL Buffer Mix C2 + 10 µL Primer Mix C1 + 10 µL Enzyme Mix C3 . Mix well.	<table border="1"> <thead> <tr> <th>No. of Samples</th> <th>C2</th> <th>C1</th> <th>C3</th> </tr> </thead> <tbody> <tr> <td>1</td> <td>20 µL</td> <td>10 µL</td> <td>10 µL</td> </tr> <tr> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>	No. of Samples	C2	C1	C3	1	20 µL	10 µL	10 µL				
No. of Samples	C2	C1	C3										
1	20 µL	10 µL	10 µL										
Add 40 µL of SPIA Master Mix to each reaction tube and resuspend beads thoroughly by pipetting. Place on ice.													
Place the tubes in a thermal cycler running Program 4 (4 °C - 1 min, 47 °C - 60 min, 80 °C - 20 min, hold at 4 °C).													
Once the thermal cycler reaches 4 °C, remove tubes, spin and place on ice.													
Transfer tubes to the magnet and let stand 5 minutes to completely clear the solution of beads.													
Carefully transfer 40 µL of the cleared supernatant to a fresh tube.													
Continue immediately with Purification of SPIA cDNA or store SPIA cDNA at -20 °C.													

PURIFICATION OF SPIA cDNA

We recommend using the QIAGEN® MinElute® Reaction Cleanup Kit for best results.	Purification Kit Part No.	Purification Kit Lot No.
Add Buffer ERC in volume of:	Spin at speed:	For a duration of:
Add Buffer PE in volume of:	Spin at speed:	For a duration of:
To elute sample use 1X TE or Buffer EB.		
Add 1X TE or Buffer EB in volume of:	Spin at speed:	For a duration of:

For Research Use Only. Not for use in diagnostic procedures.

Tecan Genomics, Inc.



USA

900 Chesapeake Drive
Redwood City, CA 94063 USA

Customer Service and

Technical Support:

Toll Free Tel: 888.654.6544
Toll Free Fax: 888.296.6544
cservice-gn@tecan.com
techserv-gn@tecan.com

Europe

P.O. Box 109,
9350 AC Leek
The Netherlands
**Customer Service and
Technical Support:**
Tel: +31-13-5780215
Fax: +31-13-5780216
europe-gn@tecan.com

Worldwide

For our international
distributors contact
information, visit
our website
www.nugen.com

.....
Australia +61 3 9647 4100 **Austria** +43 62 46 89 330 **Belgium** +32 15 42 13 19 **China** +86 21 220 63 206 **France** +33 4 72 76 04 80 **Germany** +49 79 51 94 170
Italy +39 02 92 44 790 **Japan** +81 44 556 73 11 **Netherlands** +31 18 34 48 17 4 **Nordic** +46 8 750 39 40 **Singapore** +65 644 41 886 **Spain** +34 93 595 25 31
Switzerland +41 44 922 89 22 **UK** +44 118 9300 300 **USA** +1 919 361 5200 **Other countries** +41 44 922 81 11
.....

Tecan Group Ltd. makes every effort to include accurate and up-to-date information within this publication; however, it is possible that omissions or errors might have occurred. Tecan Group Ltd. cannot, therefore, make any representations or warranties, expressed or implied, as to the accuracy or completeness of the information provided in this publication. Changes in this publication can be made at any time without notice. For technical details and detailed procedures of the specifications provided in this document please contact your Tecan representative. This brochure may contain reference to applications and products which are not available in all markets. Please check with your local sales representative.

Tecan, Ovation and SPIA are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA.

© 2019 Tecan Genomics, Inc., all rights reserved. For disclaimer and trademarks please visit www.tecan.com

www.tecan.com

