



QUICK PROTOCOL

Ovation[®] Whole Blood Solution

Part No. 1300

Enter the following information to automatically calculate the volumes needed to prepare each reaction. The calculated volumes include an appropriate overfill in excess of the nominal volume requirements to allow for volume 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:

Kit Lot No:

Number of Samples*:

THERMAL CYCLER PROGRAMS

First Strand cDNA Synthesis

Program 1: First Strand Primer Annealing

65 °C - 5 min, hold at 4 °C

Program 2: First Strand Synthesis

48 °C - 60 min, 70 °C - 15 min, hold at 4 °C

Second Strand cDNA Synthesis

Program 3: Second Strand Synthesis

37 °C - 30 min, 75 °C - 15 min, hold at 4 °C

SPIA[®] Amplification

Program 4: SPIA Amplification 1

48 °C - 30 min, hold at 4 °C

Program 5: SPIA Amplification 2

48 °C - 30 min, 95 °C - 5 min, hold at 4 °C

ENCORE BIOTIN MODULE

Refer to the Encore Biotin Module (Part No. 4200) documentation for guidance with this portion of the protocol.

*Number of samples field ties into embedded logic to calculate master mix volumes, number of reactions.



FIRST STRAND cDNA SYNTHESIS

Thaw the First Strand Reagents (blue) . Mix each reagent, spin and place on ice.			
For each assay place 5 µL of total RNA into 0.2 mL PCR tube and place on ice.			
Add 2 µL of First Strand Primer Mix A1 VER 1 , flick tubes to mix and spin.			
Place the tubes in a thermal cycler running Program 1 (65 °C - 5 min, hold at 4 °C).			
After 5 minutes at 65 °C, immediately snap cool tubes on ice.			
Prepare First Strand Master Mix (calculation allows for appropriate overfill). Please be sure to pipet A3 VER 1 enzyme slowly and rinse out tip at least five times into buffer. Per sample combine: 12 µL Buffer Mix A2 VER 4 + 1 µL Enzyme Mix A3 VER 1 . Mix well.	No. of Samples	A2	A3
	1	12 µL	1 µL
Add 13 µL of the First Strand Master Mix to each tube, mix and spin.			
Place the tubes in a thermal cycler running Program 2 (48 °C - 60 min, 70 °C - 15 min, hold at 4 °C).			

SECOND STRAND cDNA SYNTHESIS

Thaw the Second Strand Reagents (yellow) . Mix each reagent, spin and place on ice.			
Once the thermal cycler reaches 4 °C, remove tubes, spin and place on ice.			
Prepare Second Strand Master Mix (calculation allows for appropriate overfill). Per sample combine: 18 µL Buffer Mix B1 VER 4 + 2 µL Enzyme Mix B2 VER 1 . Mix well.	No. of Samples	B1	B2
	1	18 µL	2 µL
Add 20 µL of Second Strand Master Mix to each first strand reaction tube, mix and spin.			
Place the tubes in a thermal cycler running Program 3 (37 °C - 30 min, 75 °C - 15 min, hold at 4 °C).			
Once the thermal cycler reaches 4 °C, spin and place tubes on ice.			

SPIA AMPLIFICATION

Thaw the **SPIA Amplification Reagents (red)**. Vortex **C1 VER 1**, **C2 VER 6** and **C4 VER 1**. Invert **C3 VER 5** 5 times. Spin all, place on ice.

Prepare **SPIA Master Mix** (calculation allows for appropriate overfill). Per sample combine:
72 µL **C2 VER 6** + 2 µL **C1 VER 1** + 4 µL **D1** + 40 µL **C3 VER 5**. **Mix well.**

No. of Samples	C2	C1	D1	C3
1	72 µL	2 µL	4 µL	40 µL

On ice, add 118 µL of **SPIA Master Mix** to each second strand reaction tube, mix and spin.

Place half of the reaction volume (79 µL) into a separate 0.2 mL PCR tube, cap tightly and spin.

Place tubes in a thermal cycler running Program 4 (48 °C - 30 min, hold at 4 °C).

Add 3 µL of **C4 VER 1** to each reaction tube, mix and spin.

Return tubes to the thermal cycler running Program 5 (48 °C - 30 min, 95 °C - 5 min, hold at 4 °C).

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

Recombine the split reactions (total volume of 164 µL).

Proceed immediately to purification step or store SPIA cDNA at -20 °C.

PURIFICATION OF AMPLIFIED SPIA cDNA

Refer to the user guide and follow the method of choice for purification:

Purification Kit Part No.

Purification Kit Lot No.

Add Binding Buffer in volume of:

Spin at speed:

For a duration of:

Add Wash Buffer in volume of:

Spin at speed:

For a duration of:

Number of washes:

To elute sample use **Nuclease-free Water D1** provided with the Ovation RNA Amplification System V2 kit.

Add **Nuclease-free Water D1** 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

