



QUICK PROTOCOL

Ovation[®] Ultralow System V2

Part Nos. 0344, 0344NB

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:

Kit Lot No:

Number of Samples[†]:

THERMAL CYCLER PROGRAMS

End Repair

Program 1: End Repair

25 °C - 30 min, 70 °C - 10 min, hold at 4 °C

Ligation

Program 2: Ligation

25 °C - 30 min, 70 °C - 10 min, hold at 4 °C

Amplification

Program 3: Amplification

72 °C - 2 min, 95 °C - 3 min, 7-15* cycles (98 °C - 20 s, 65 °C - 30 s, 72 °C - 30 s) 72 °C - 1 min, hold 4 °C

***Important Note:** The number of cycles used for PCR amplification depends on the starting amount of genomic DNA. Alternatively, real-time PCR can be used to determine the appropriate number of PCR cycles. For more information, contact Tecan NGS Technical Support.

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



Remove reagents for each section from storage as needed. Return to storage promptly after use. Continue immediately to the next section unless otherwise directed.

END REPAIR

Thaw ER1 , ER2 , ER3 and D1 . Mix as needed and spin down. Place ER1 , ER2 and ER3 on ice, and D1 at room temperature.				
Prepare End Repair Master Mix (calculation allows for appropriate overfill). Per sample combine: 3.5 µL End Repair Buffer Mix ER1 + 0.5 µL End Repair Enzyme Mix ER2 + 1.0 µL End Repair Enhancer ER3 Mix thoroughly in a 0.5 mL capped tube.	No. of Samples	ER1	ER2	ER3
	1	3.5 µL	0.5 µL	1.0 µL
Add 5 µL of End Repair Master Mix to each tube containing 10 µL fragmented DNA sample (10 pg-100 ng), mix by pipetting, spin and place on ice.				
Place tube in a thermal cycler running Program 1: (25 °C - 30 min, 70 °C - 10 min, hold at 4 °C)				
Once thermal cycler reaches 4 °C, remove tubes, spin and place on ice.	Leave D1 at room temp for use in the next step.			

LIGATION

Thaw L1 , L2 , L3 and D1 . Mix as needed and spin down. Place L1 , L2 and L3 on ice, and D1 at room temperature.				
If using adaptors from tubes (O344NB-08), add 6 µL of appropriate Ligation Adaptor Mix L2 to each sample and mix thoroughly. If using an adaptor plate (O344-32 or O344NB-32), add the entire 15 µL End Repair mixture to the appropriate adaptor well, mix well, then transfer the reaction to a PCR tube.	No. of Samples	D1	L1	L3
	1	1.5 µL	6.0 µL	1.5 µL
Prepare Ligation Master Mix (calculation allows for appropriate overfill). Per sample combine: 1.5 µL Nuclease-free Water D1 + 6.0 µL Ligation Buffer Mix L1 + 1.5 µL Ligation Enzyme Mix L3 Mix thoroughly in a 0.5 mL capped tube.				
Add 9 µL of Ligation Master Mix to each tube, mix thoroughly by pipetting slowly and gently, spin and place on ice.				
Place tube in a thermal cycler running Program 2: (25 °C - 30 min, 70 °C - 10 min, hold at 4 °C)				
Once thermal cycler reaches 4 °C, remove tubes, spin and place on ice.	Leave D1 at room temp for use in the next step.			

LIGATION PURIFICATION

Ensure the Agencourt beads have reached room temperature.	
Add 70 µL of room temperature Nuclease-free Water (D1) to each ligation reaction.	
Mix the beads by inverting several times. At room temperature, add 80 µL of Agencourt XP 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 the binding buffer and discard it.	
Add 200 µL of freshly prepared 70% ethanol and let stand for 30 seconds. Remove the ethanol using a pipette.	
Repeat the ethanol wash one more time for a total of two washes.	
Remove all excess ethanol after the final wash and let beads air dry for 10 minutes.	
Ensure the tubes have completely dried and no residual ethanol is left.	
Add 40 µL 1xTE buffer (low EDTA), mix thoroughly by pipetting and let stand for 3 minutes.	
Transfer the tubes to the magnet and let stand for 3 minutes.	
Remove 35 µL of eluate and transfer to a fresh set of PCR tubes and place on ice.	



*Optional stopping point:
store samples at -20 °C.*

LIBRARY AMPLIFICATION

Thaw P1 , P2 and P3 . Mix as needed, spin down and place on ice.													
Prepare Amplification Master Mix (calculation allows for appropriate overfill). Per sample combine: 12.75 µL Amplification Buffer Mix P1 + 1.25 µL Amplification Primer Mix P2 + 1.0 µL Amplification Enzyme Mix P3 Mix thoroughly in a 0.5 mL capped tube.	<table border="1"> <thead> <tr> <th>No. of Samples</th> <th>P1</th> <th>P2</th> <th>P3</th> </tr> </thead> <tbody> <tr> <td>1</td> <td>12.75 µL</td> <td>1.25 µL</td> <td>1.0 µL</td> </tr> <tr> <td></td> <td></td> <td></td> <td></td> </tr> </tbody> </table>	No. of Samples	P1	P2	P3	1	12.75 µL	1.25 µL	1.0 µL				
	No. of Samples	P1	P2	P3									
	1	12.75 µL	1.25 µL	1.0 µL									
<table border="1"> <thead> <tr> <th>Starting Input</th> <th>PCR cycles</th> </tr> </thead> <tbody> <tr> <td>< 1 ng</td> <td>may require optimization</td> </tr> <tr> <td>1 - 10 ng</td> <td>13 - 15</td> </tr> <tr> <td>10 - 50 ng</td> <td>10 - 12</td> </tr> <tr> <td>50 - 100 ng</td> <td>7 - 9</td> </tr> </tbody> </table>	Starting Input	PCR cycles	< 1 ng	may require optimization	1 - 10 ng	13 - 15	10 - 50 ng	10 - 12	50 - 100 ng	7 - 9			
Starting Input	PCR cycles												
< 1 ng	may require optimization												
1 - 10 ng	13 - 15												
10 - 50 ng	10 - 12												
50 - 100 ng	7 - 9												
On ice, add 15 µL of Amplification Master Mix to each tube, mix by pipetting, spin and place on ice.													
Place the tubes in a thermal cycler programmed to run Program 3. Recommended PCR cycle numbers are given to the right. 72 °C - 2 min, 95 °C - 3 min, 7-15 cycles (98 °C - 20 s, 65 °C - 30 s, 72 °C - 30 s) 72 °C - 1 min, hold 4 °C													
Once thermal cycler reaches 4 °C, remove tubes, spin and place on ice.													

AMPLIFIED LIBRARY PURIFICATION

Retrieve Agencourt beads and 70% ethanol set aside previously at room temperature.

Mix the beads by inverting several times.

At room temperature, add 50 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 binding buffer and discard it.

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 10 minutes.

Ensure the tubes have completely dried and no residual ethanol is left.

Add 33 μL 1xTE buffer (low EDTA), mix thoroughly by pipetting.

Transfer the tubes to the magnet and let stand for 2 minutes.

Remove 30 μL of eluate and transfer to a fresh set of PCR tubes.

Proceed to Qualitative and Quantitative Assessment of the library.



*Optional stopping point:
store samples at $-20\text{ }^{\circ}\text{C}$.*

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

© 2020 Tecan Genomics, Inc., all rights reserved. For disclaimer and trademarks please visit www.tecan.com

www.tecan.com

