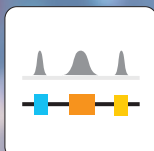


Ovation[®] SoLo[™] RNA-Seq.

PRODUCT SHEET



Whole transcriptome analysis at ultra-low input levels

Ovation SoLo RNA-Seq System is a complete end-to-end sample to library preparation solution for stranded, whole transcriptome RNA-Seq analysis of ultra-low inputs from exosomes, formalin-fixed, paraffin-embedded (FFPE) samples and cell-free RNA. The highly flexible and customizable **AnyDeplete[™]** technology offers depletion of rRNA and other highly abundant, unwanted transcripts. This depletion increases the dynamic range that can be studied, reduces sequencing costs and simplifies data analysis.

Why use the Ovation SoLo RNA-Seq System?

1. Fully integrated workflow for whole transcriptome analysis of as little as 10 pg of precious samples, providing access to complete biological information.
2. Customizable, targeted transcript depletion with AnyDeplete, allowing elimination of unwanted reads after library preparation, even from ultra low inputs.
3. Direct integration with cell lysis protocols, providing an easy-to-use workflow for ultra-low RNA-Seq.

Benefits

- Whole transcriptome analysis
- 10 pg to 10 ng of isolated RNA or cell lysate
- Efficient stranded library preparation without adaptor titration
- Pre-plated adaptors with unique barcodes included for every sample
- Integrated molecular tag for accurate detection of PCR duplicates
- Customizable transcript depletion post-library preparation

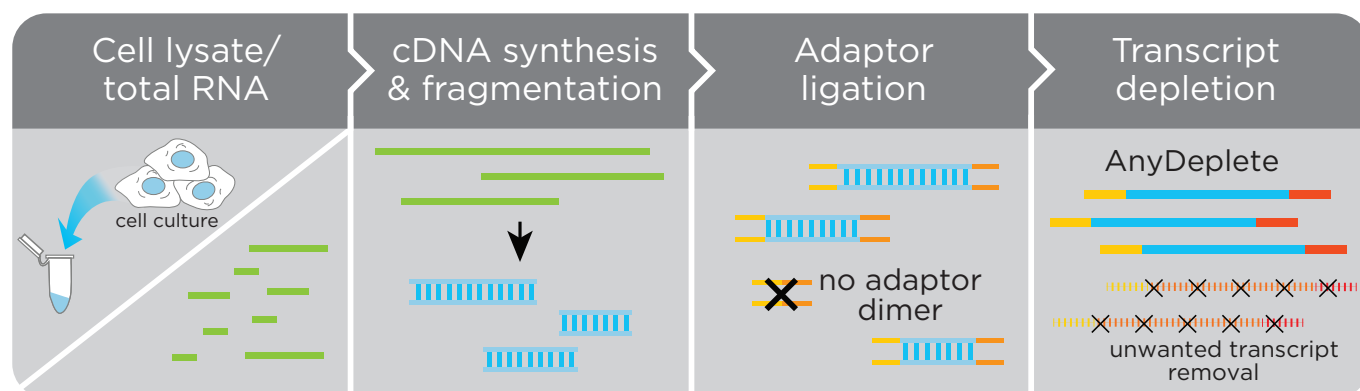


Figure 1: Ovation SoLo RNA-Seq is a complete workflow enabling whole transcriptome analysis from ultra-low inputs.



Technical details

- Input range: 10 pg to 10 ng total RNA or cell lysate
- Excellent correlation down to 10 pg sample input
- Excellent correlation and data from low quality samples

Applications

- Whole transcriptome profiling
- Gene expression
- RNA-Seq from a wide variety of sample sources: liquid biopsy, FFPE samples, exosomes, cell-free RNA, and tissues

Why use AnyDeplete?

- Remove unwanted transcripts after library preparation
- Add new probes to existing probe sets without re-optimization
- Customizable depletion of species- or experiment-specific transcripts

Highly correlated data from 1 ng to 10 pg

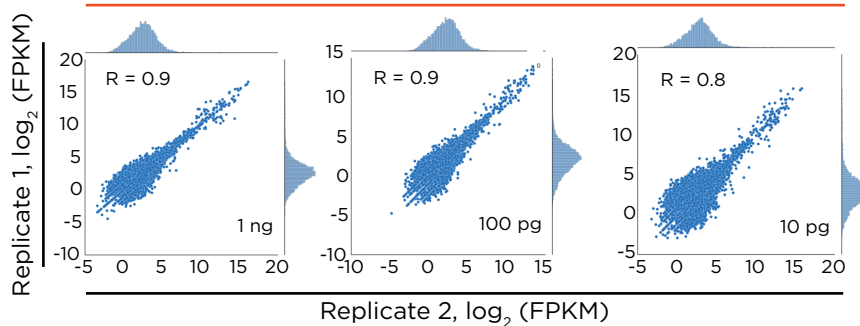


Figure 2: High correlation between technical replicates from inputs as low as 10 pg of total RNA, giving you confidence in your data.

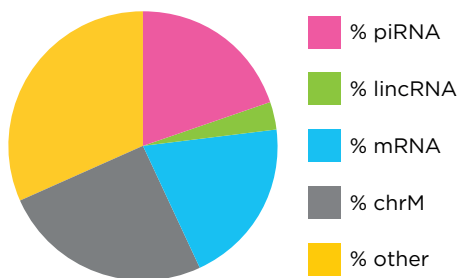


Figure 3: Libraries generated from 5 ng of cell-free plasma RNA with a DV200 score of 10-20 %. Analysis of the whole transcriptome provides complete information from coding, non-coding and regulatory transcripts.

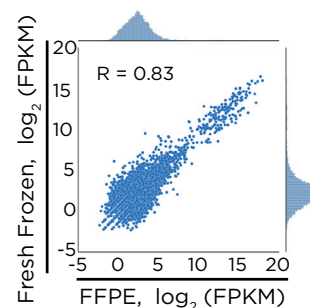


Figure 4: Ovation SoLo RNA Seq provides highly correlated data between fresh frozen and FFPE adenocarcinoma samples.

Ordering information

Ordering information	Part no.	No. of reactions
Ovation SoLo RNA-Seq, Human AnyDeplete	0500	32, 96
Ovation SoLo RNA-Seq, Mouse AnyDeplete	0501	32, 96
Ovation SoLo RNA-Seq, Drosophila AnyDeplete	0502	32, 96
Ovation SoLo RNA-Seq, Custom AnyDeplete	Contact your account executive	

AnyDeplete probe sets can be customized to any transcript from any organism. For custom probe sets, contact your account executive or request a quote via our website.

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

www.tecan.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, SoLo and AnyDeplete 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.

