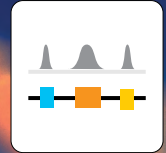


RNA-Seq Solutions.



PRODUCT SELECTION GUIDE

	Revelo™ High Sensitivity RNA-Seq	Universal Plus™ mRNA-Seq	Universal Plus™ Total RNA-Seq	Ovation® RNA-Seq V2	SoLo® RNA-Seq
Applications	<ul style="list-style-type: none"> Whole transcriptome profiling Differential gene expression Rare transcript detection Infectious disease 	<ul style="list-style-type: none"> mRNA Seq Splice variant analysis Differential gene expression Fusion detection 	<ul style="list-style-type: none"> Whole transcriptome profiling Differential gene expression Transcript discovery Splice variant and isoform detection 	<ul style="list-style-type: none"> cDNA preparation from total RNA 	<ul style="list-style-type: none"> Whole transcriptome profiling Differential gene expression Transcript discovery Splice variant and isoform detection
Sample Type	<ul style="list-style-type: none"> Total RNA FFPE Nasal swabs Microbial RNA 	<ul style="list-style-type: none"> Total RNA (polyadenylated) 	<ul style="list-style-type: none"> Total RNA (any organism) 	<ul style="list-style-type: none"> Total RNA FFPE Nasal swabs Microbial RNA 	<ul style="list-style-type: none"> Total RNA FFPE RNA cfRNA (liquid biopsies) Single cell RNA Exosome RNA Microbial RNA
Input Amount	250 pg - 10 ng	10 ng - 1 µg	10 ng - 500 ng	500 pg - 100 ng	10 pg - 10 ng
Integrated Fragmentation	Yes	Yes	Yes	—	Yes
Workflow Time	6.5 hrs	7.5-8.5 hrs	7.5 hrs	4.5 hrs	14.75 hrs
Index Adaptors (barcode length, number of barcodes)	Unique Dual Index (UDI) (8 nt, up to 96)	UDI (8 nt, up to 384)	UDI (8 nt, up to 384)	N/A	Single index (8 nt, up to 96)
UMI (Unique Molecular Identifiers)	—	Yes	Yes	N/A	Yes
Strand Specific	—	Yes	Yes	N/A	Yes
Automation Readiness	DreamPrep® NGS (Fluent® 780 workstation), and other platforms				—
Sequencing Platform	Illumina®			Any	Illumina®
Tecan Patented Technologies:					
NuQuant®	✓	✓	✓	N/A	—
DimerFree®	✓	✓	✓	N/A	✓
SPIA® (Single Primer Isothermal Amplification)	✓	—	—	✓	—
Transcript Depletion Technology	SPIABoost™	AnyDeplete®	AnyDeplete	N/A	AnyDeplete

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, Revelo, Universal Plus, SoLo, Ovation, DreamPrep, Fluent, NuQuant, DimerFree, SPIA, SPIABoost, and AnyDeplete are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA. Illumina® is a registered trademark of Illumina, Inc., San Diego, CA, US.

© 2023 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com

