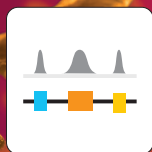


# Universal Plus Total RNA-Seq

with NuQuant<sup>®</sup>

## PRODUCT SHEET



### From total RNA to quantified libraries - a complete solution for whole transcriptome RNA-Seq

Universal Plus Total RNA-Seq with NuQuant provides a simple solution for the generation of stranded whole transcriptome RNA-Seq libraries ready for Illumina sequencing, and is compatible with varying quality input material from a range of sample types.

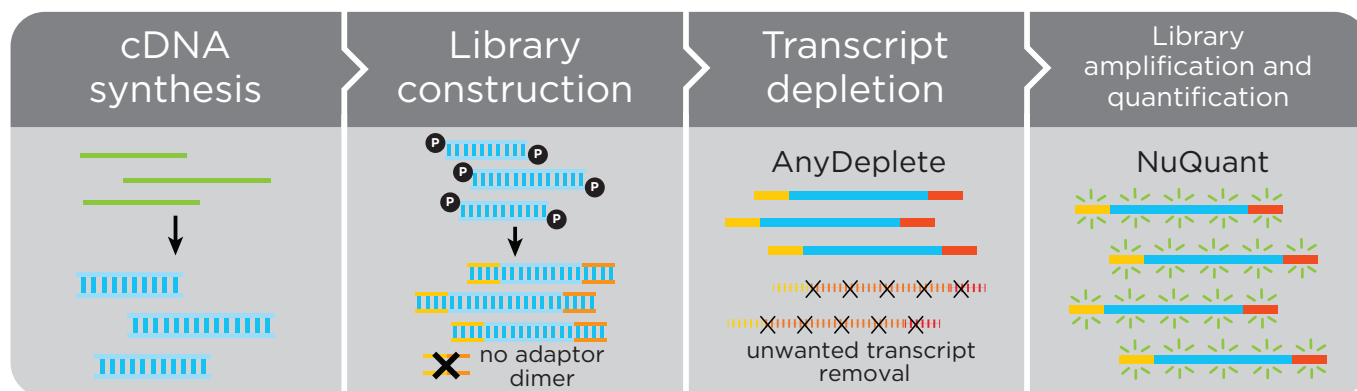


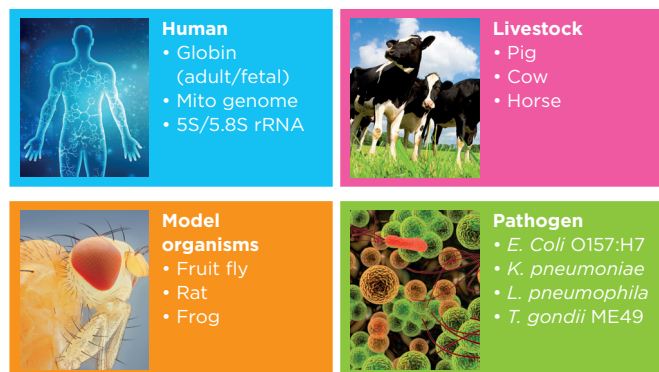
Figure 1: Universal Plus Total RNA-Seq library preparation kits provide end-to-end solutions for strand-specific libraries.



## Why use Universal Plus Total RNA-Seq

1. **Reduced hands-on time** – with fewer bead purifications and a simple workflow
2. **384 Unique Dual Index (UDI) adaptors** – for flexible multiplexing and sequencing
3. **AnyDeplete®** integrated targeted rRNA depletion after library construction:
  - Highly specific depletion maximizes informative reads
  - Lower input requirements than other depletion methods
  - Customizable for any transcript from any organism
4. **NuQuant®** library quantification – providing fast and easy determination of library molarity in minutes

### Available custom AnyDeplete probes



Contact Tecan NGS Technical Support for detailed target information for each design.

### Features

- Compatible with high quality fresh or FFPE/degraded total RNA samples
- Up to 384 unique, preplated UDI adaptors with DimerFree® technology included with the kit for every reaction
- Efficient target depletion of rRNA with AnyDeplete available for model organisms, or customizable to your desired targets
- Saves time and costs for library QC with NuQuant
- Automation-friendly



Figure 2: **Comparison of library quantification methods.** NuQuant is integrated into the kit and provides simple fluorescence-based determination of molar library concentration in minutes, without additional cost.

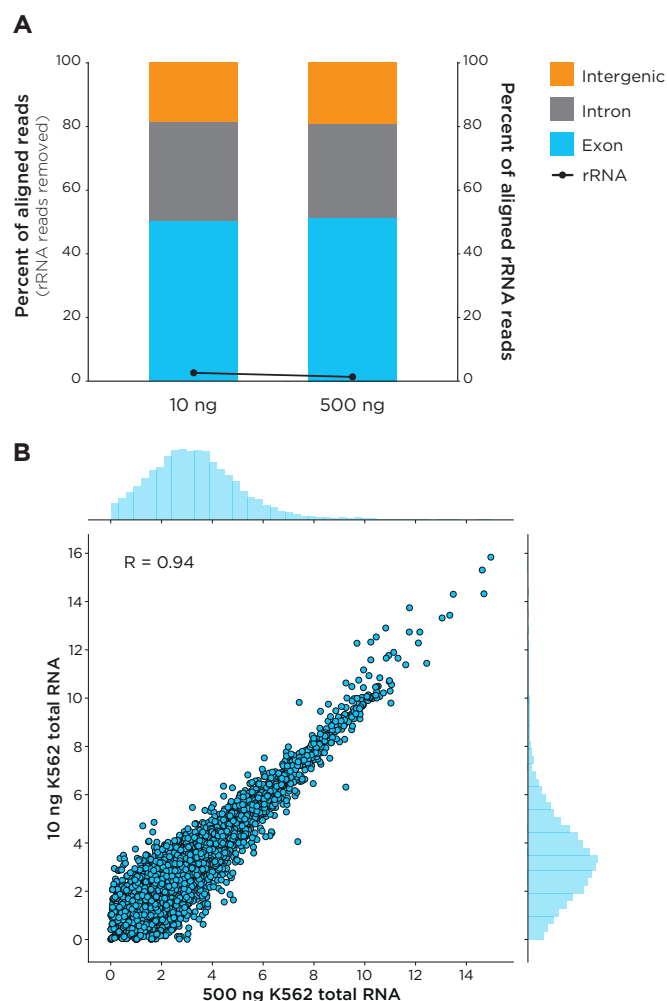


Figure 3: **Consistent data across a broad input range.** A) Total RNA-Seq libraries were generated with 10 ng and 500 ng of K562 total RNA (N=3). Metrics from the different libraries show similar results, demonstrating effective performance across the entire recommended input range. B) FPKM correlation between 10 ng and 500 ng input libraries shows an R value of 0.94, indicating similar data regardless of the 50-fold difference in input amounts.

Sample	DV200 (%)	Input (ng)	% aligned	% rRNA	% exon
Liver FFPE	64	100	95.3	13.8	31.8
Kidney FFPE	41	150	78.6	32.5	32.3

Table 1: **The Universal Plus Total RNA-Seq with NuQuant kit is compatible with FFPE and degraded RNA.** FFPE samples with a DV200 value (percentage of RNA fragments greater than 200 bases) of 64 % and 41 % were used to generate libraries (N=3). Even highly degraded FFPE samples generated high quality libraries with good alignment and a high percentage of exons, providing a single solution for whole transcriptome RNA-Seq for any sample type.

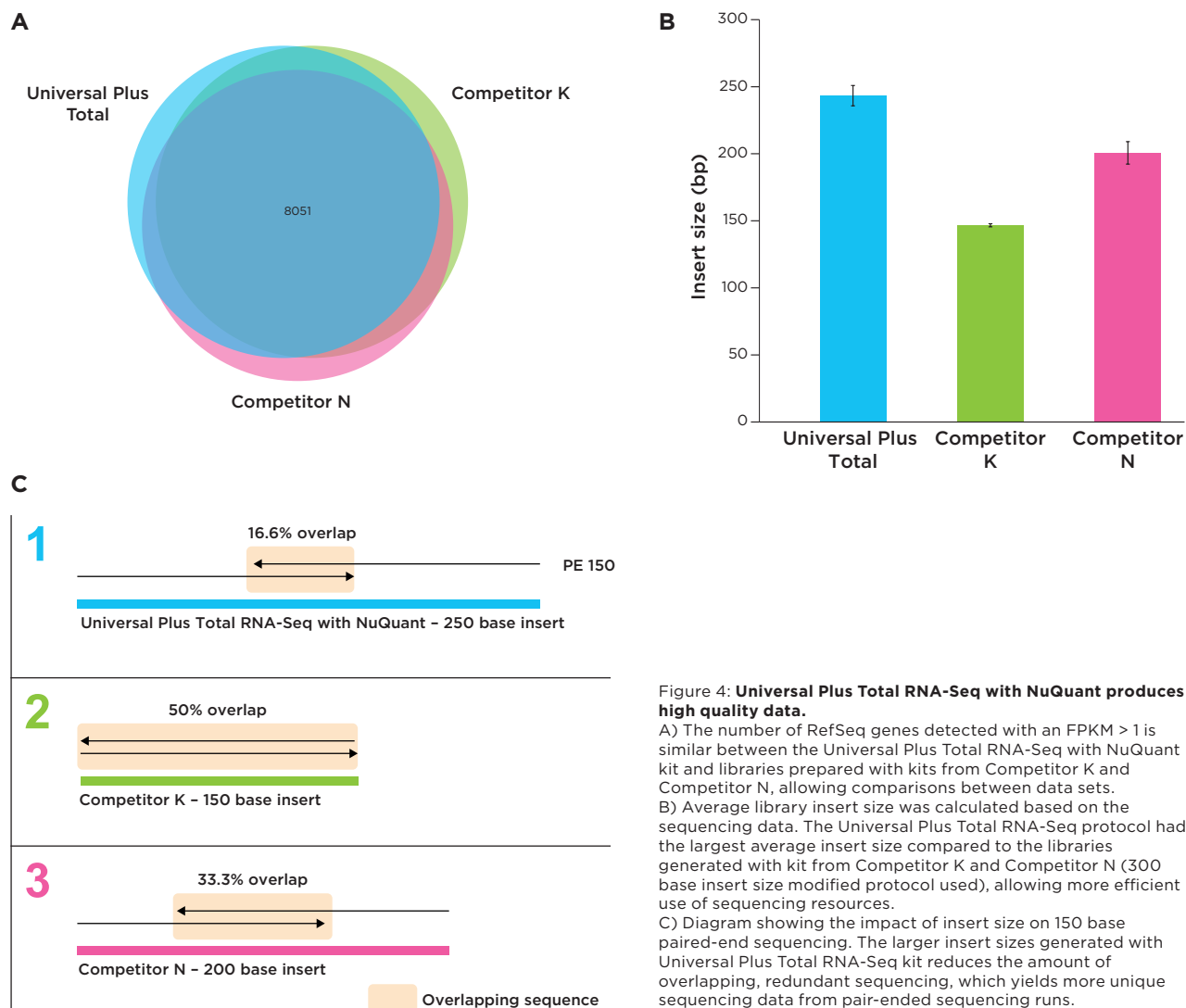


Figure 4: **Universal Plus Total RNA-Seq with NuQuant produces high quality data.**

A) The number of RefSeq genes detected with an FPKM > 1 is similar between the Universal Plus Total RNA-Seq with NuQuant kit and libraries prepared with kits from Competitor K and Competitor N, allowing comparisons between data sets.

B) Average library insert size was calculated based on the sequencing data. The Universal Plus Total RNA-Seq protocol had the largest average insert size compared to the libraries generated with kit from Competitor K and Competitor N (300 base insert size modified protocol used), allowing more efficient use of sequencing resources.

C) Diagram showing the impact of insert size on 150 base paired-end sequencing. The larger insert sizes generated with Universal Plus Total RNA-Seq kit reduces the amount of overlapping, redundant sequencing, which yields more unique sequencing data from pair-ended sequencing runs.

### Technical details

- 10-500 ng total RNA input range
- Robust chemical fragmentation
- Stranded libraries
- Consistent data across a wide input range
- Compatible with FFPE and degraded RNA samples
- Optimized protocol for larger insert size libraries

### Applications

- Whole transcriptome profiling
- Gene expression analysis
- Transcript discovery
- Splice variant detection

## Ordering information

Product name	AnyDeplete probe pool included	Part no.	Barcodes	No. of reactions
Universal Plus Total RNA-Seq with NuQuant, Human AnyDeplete	Human rRNA	9156-24	1-24	24
		9156-A01	1-96	96
		9156B-A01	97-192	96
		9156C-A01	193-288	96
		9156D-A01	289-384	96
Universal Plus Total RNA-Seq with NuQuant, Mouse AnyDeplete	Mouse rRNA, globin	9157-24	1-24	24
		9157-A01	1-96	96
		9157B-A0	97-19	96
		9157C-A01	193-28	96
		9157D-A01	289-384	96
Universal Plus Total RNA-Seq with NuQuant, Custom AnyDeplete*	Custom rRNA	9158-24	1-24	24
		9158-A01	1-96	96
		9158B-A01	97-192	96

\*Contact your local Tecan sales representative for more information

**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, NuQuant, DimerFree and AnyDeplete 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](http://www.tecan.com)

