

Microarray and qPCR Solutions.

PRODUCT SELECTION GUIDE

RNA Sample Preparation

Platform	Starting materials and sample sizes				
	Good Quality >50 ng	Good Quality 5-50 ng	Good Quality & Degraded 500 pg-50 ng	Whole Blood 5-50 ng	FFPE 50 ng
Affymetrix GeneChip® 3' Arrays	Ovation® RNA Amplification System V2 Encore Biotin Module	Ovation RNA Amp System V2 Encore Biotin Module	Ovation Pico WTA System V2 Encore Biotin Module	Ovation RNA Amp System V2 Ovation WB Reagent Encore Biotin Module	Ovation FFPE WTA System Encore Biotin Module
Affymetrix GeneChip Exon/Gene ST Arrays		Ovation Pico WTA System V2 Ovation PicoSL WTA System V2 Encore Biotin Module	Ovation Pico WTA System V2 Ovation PicoSL WTA System V2 Encore Biotin Module	Ovation Pico WTA System V2 Ovation PicoSL WTA System V2 Encore Biotin Module	Ovation FFPE WTA System Encore Biotin Module
Agilent Dual-Mode Gene Expression Arrays	Ovation RNA Amplification System V2	Ovation RNA Amp System V2	Ovation PicoSL WTA System V2	Ovation RNA Amp System V2	Ovation FFPE WTA System*
Quantitative PCR		Crescendo cDNA Synthesis for qPCR	Crescendo cDNA Synthesis for qPCR	Crescendo cDNA Synthesis for qPCR	Crescendo cDNA Synthesis for qPCR

*May require optimization of fragment sizes

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, Ovation, Encore, Applause, SPIA, Ribo-SPIA and Crescendo cDNA are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA.

© 2021 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com

