

# Microarray Solutions.



## PRODUCT SELECTION GUIDE

	Input Amount and Quality	
	500 pg - 50 ng High quality or degraded RNA	50 - 100 ng FFPE or degraded RNA
Platform	Recommended Kit	
<b>Affymetrix GeneChip 3' Arrays</b>	Ovation® Pico WTA System V2 with Encore™ Biotin Module	Ovation® FFPE WTA System with Encore™ Biotin Module
<b>Affymetrix GeneChip Exon/ Gene ST Array</b>	Ovation® Pico WTA System V2 with Encore™ Biotin Module	Ovation® FFPE WTA System with Encore™ Biotin Module
<b>Agilent® Dual-Mode Gene Expression Arrays</b>	Ovation® Pico WTA System V2	Ovation® FFPE WTA System*

\*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, and Encore are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA. Agilent is a registered trademark of Agilent Technologies Ltd., Santa Clara, CA, US.

© 2023 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit [www.tecan.com](http://www.tecan.com)

