

DNA-Seq Solutions.

PRODUCT SELECTION GUIDE

Input Amount	10ng - 500ng	10ng - 500ng	100ng - 500ng	200ng - 500ng	10pg - 100ng	10ng - 100ng
Input Type	High quality gDNA	High quality gDNA	High quality gDNA	High quality gDNA	High Quality gDNA, FFPE, cfDNA or degraded DNA	High quality gDNA
Enzymatic Fragmentation	NO	YES	NO	YES	NO	YES
PCR-Free	NO	NO	YES	YES	NO	NO
Unique Dual Index	YES	YES	YES	YES	YES	NO
High-throughput Multiplexing	YES	YES	NO	NO	NO	YES
Molecular Tags	NO	NO	NO	NO	YES	NO
Custom Target Enrichment Probes	NO	NO	NO	NO	NO	YES
NuQuant®	YES	YES	NO	NO	NO	NO
	Celero™ DNA-Seq	Celero EZ DNA-Seq	Ovation® Rapid DNA-Seq	Rapid EZ DNA-Seq™	Ovation Ultralow DNA-Seq V2	Allegro® Targeted Genotyping V2

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, Celero, NuQuant, Rapid EZ DNA-Seq and Allegro are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA.

© 2020 Tecan Trading AG, Switzerland, all rights reserved. For disclaimer and trademarks please visit www.tecan.com

