

# DNA-Seq Solutions.



## PRODUCT SELECTION GUIDE

	Ovation® Ultralow System V2	Celero™ DNA-Seq	Celero™ EZ DNA-Seq	Allegro® Targeted Genotyping V2
<b>Applications</b>	<ul style="list-style-type: none"> <li>WGS</li> <li>WES</li> <li>Targeted sequencing</li> <li>Microbial sequencing</li> <li>Amplicon sequencing</li> </ul>	<ul style="list-style-type: none"> <li>WGS</li> <li>WES</li> <li>Targeted sequencing</li> <li>Microbial sequencing</li> <li>Amplicon sequencing</li> </ul>	<ul style="list-style-type: none"> <li>WGS</li> <li>WES</li> <li>Targeted sequencing</li> <li>Microbial sequencing</li> <li>Amplicon sequencing</li> </ul>	<ul style="list-style-type: none"> <li>Targeted Genotyping</li> </ul>
<b>Sample Type</b>	<ul style="list-style-type: none"> <li>cfDNA</li> <li>FFPE</li> <li>Pre-fragmented gDNA</li> <li>ds cDNA</li> </ul>	<ul style="list-style-type: none"> <li>FFPE</li> <li>Pre-fragmented gDNA</li> <li>ds cDNA</li> <li>Amplicons</li> </ul>	<ul style="list-style-type: none"> <li>High quality gDNA</li> <li>Large DNA fragments</li> </ul>	<ul style="list-style-type: none"> <li>High quality gDNA</li> </ul>
<b>Input Amount</b>	10 pg - 100 ng	10 ng - 500 ng	10 ng - 500 ng	10 ng - 100 ng
<b>Enzymatic Fragmentation</b>	—	—	Yes	Yes
<b>Workflow Time</b>	4 hrs	3.5 hrs	3.5 hrs	6 hrs over 1.5 days
<b>Index Adaptors</b> (barcode length, number of barcodes)	<ul style="list-style-type: none"> <li>Single Index (8 nt, up to 96)</li> <li>Unique Dual Index (UDI), (8 nt, up to 96)</li> </ul>	<ul style="list-style-type: none"> <li>Combinatorial Dual Index (8 nt, up to 384)</li> <li>UDI (10 nt, up to 384)</li> </ul>	<ul style="list-style-type: none"> <li>Combinatorial Dual Index (8 nt, up to 384)</li> <li>UDI (10 nt, up to 384)</li> </ul>	<ul style="list-style-type: none"> <li>Single Index (8 nt, up to 192)</li> <li>Combinatorial Dual Index (when used with Metaplex module, 8 nt, up to 3072)</li> </ul>
<b>UMI</b> (Unique Molecular Identifiers)	Yes, with UDI adaptors	—	—	—
<b>Automation Readiness</b>	DreamPrep® NGS (Fluent® 780 workstation), and other platforms			
<b>Sequencing platform</b>	Illumina®			
<b>Tecan Patented Technologies:</b>				
<b>NuQuant®</b>		✓	✓	
<b>DimerFree®</b>	✓	✓	✓	✓
<b>SPET™</b> (Single Primer Enrichment Technology)				✓

**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, DreamPrep, Fluent, Allegro, NuQuant, DimerFree, and SPET 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](http://www.tecan.com)



**TECAN.**