

Methyl-Seq Solutions.



PRODUCT SELECTION GUIDE

	Ovation® RRBS Methyl-Seq with TrueMethyl® oxBS	Ultralow™ Methyl Seq with TrueMethyl® OXBS	TrueMethyl® oxBS Module
Applications	<ul style="list-style-type: none"> Reduced Representation Bisulfite Sequencing 	<ul style="list-style-type: none"> Methyl-Seq Bisulfite sequencing Oxidative bisulfite sequencing 	<ul style="list-style-type: none"> Bisulfite conversion Oxidative bisulfite conversion Methyl-Seq DNA methylation arrays
Sample Type	<ul style="list-style-type: none"> High quality gDNA 	<ul style="list-style-type: none"> Pre-fragmented or degraded DNA FFPE 	<ul style="list-style-type: none"> Pre-fragmented or degraded DNA FFPE
Input Amount	100 ng	10 ng - 300 ng	100 ng - 1 µg
Enzymatic Fragmentation	Yes	—	—
Workflow Time	9 hrs	9.25 hrs	2.25 hrs
Index Adaptor (barcode length, number of barcodes)	Single Index (8 nt, up to 96)	Single Index (8 nt, up to 96)	N/A
Diversity adaptors	Yes	—	N/A
UMI (Unique Molecular Identifiers)	Yes	—	N/A
Oxidative Bisulfite Reagents Included	Yes (5mC and 5hmC differentiation)	Yes (5mC and 5hmC differentiation)	Yes (5mC and 5hmC differentiation)
Automation Readiness	<ul style="list-style-type: none"> Beckman Coulter® Biomek i7 Agilent® Bravo® 		—
Sequencing platform	Illumina®		Any

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 Ultralow are registered trademarks and trademarks of Tecan Group Ltd., Männedorf, Switzerland or of Tecan Genomics, Inc., Redwood City, USA. TrueMethyl® is a registered trademark of Cambridge EpiGenetix, Ltd., London, GB. Illumina® is a registered trademark of Illumina, Inc., San Diego, CA, US.

© 2023 Tecan Genomics, Inc., all rights reserved. For disclaimer and trademarks please visit www.tecan.com

