

## Biomek-Automated NGS Library Construction Methods

	DEMONSTRATED METHOD	BIOMEK i7 HYBRID	BIOMEK i7/i5 MULTICHANNEL	BIOMEK i5 SPAN-8	BIOMEK FXP HYBRID	BIOMEK FXP/NXP MULTICHANNEL	BIOMEK NXP SPAN-8	BIOMEK 4000
	Illumina AmpliSeq Cancer HotSpot Panel v2 Illumina AmpliSeq Comprehensive Panel v3			•				
	Illumina AmpliSeq Focus Panel			•				
	Illumina TruSight Oncology 500 DNA/RNA					•		
	Illumina TruSight Oncology 500 HT	*						
CANCER PANELS	Illumina TruSight Cancer							•
	Illumina TruSight Tumor 15 Illumina TruSight Tumor 170			•				•
	Illumina TruSight Tumor 500						•	
	NEB NEBNext Direct Cancer HotSpot Panel	-			•			
	TruSight RNA Pan Cancer							
	ATAC-Seq							
	Illumina AmpliSeq for Illumina NGS Illumina DNA PCR-Free Prep			•				
	Illumina ForenSeq			_				
	Illumina Nextera DNA Flex Library Prep Kit	•			•			
	Illumina Nextera XT	•			•			
	Illumina TruSeq Custom Amplicon Low Input Kit				•			
	Illumina TruSeq DNA PCR-Free	-		•			_	-
DNA	Illumina TruSeq Nano DNA  NEB NEBNext Ultra DNA for Illumina NGS (ChIP-seq and HLA)	•			•		•	_
SEQUENCING	NEB NEBNext Ultra II DNA Kit for Illumina NGS				•			
	NEB NEBNext Ultra II FS DNA							
	NuGEN Ovation Ultralow Library System v2				<b>A</b>			
	NuGEN Ovation Rapid Library System (PCR-Free)				<b>A</b>			
	Roche KAPA HyperPrep Library Prep Kit for Illumina NGS  Roche KAPA HyperPlus Library Prep Kit for Illumina NGS				<b>A</b>			
	Swift Biosciences Accel - NGS 2S Plus DNA Library Kit for Illumina NGS							
	Takara SMARTer ThruPLEX DNA-seq				•			
	Twist Library Preparation Kit (Mechanical & enzymatic fragmentation)	-		•				
	GenDx NGSgo							•
HLA TYPING	Illumina TruSight HLA Sequencing Panel v2 Immucor MIA FORA Flex HLA	<b>A</b>			<b>A</b>		•	
	Omixon Holotype HLA X2							
	Illumina AmpliSeq for Illumina NGS			•				
	Illumina RNA Prep with Enrichment	-						
	Illumina TruSeq RNA v2				•			
	Illumina TruSeq RNA Exome							
	Illumina TruSeq Stranded mRNA Illumina TruSeq Stranded Total RNA				•			
	Lexogen QuantSeq 3' mRNA -Seq Library Prep Kit							
RNA SEQUENCING	NEBNext ARTIC SARS-CoV-2 FS Express	-						
	NEB NEBNext Small RNA Kit for Illumina NGS							-
	NEB NEBNext Ultra II Directional RNA Library Kit for Illumina NGS				•			
	NuGEN Ovation Universal RNA-Seq  NuGEN Trio RNA-Seq				<b>A</b>			
	NuGEN Universal Plus mRNA-Seq							
	Roche KAPA mRNA HyperPrep Kit				<b>A</b>			
	Roche KAPA RNA HyperPrep Kit with RiboErase (HMR)	-			<b>A</b>			
SINGLE-CELL SEQ	10X Genomics Single Cell 3' kit- Library Preparation							
	NEB NEBNext Single Cell/ Low Input RNA Library Prep Kit	•						
	Takara SmartSeq HT  Agilent HaloPlexTarget Enrichment - Ion Torrent	-			•			_
	Agilent SureSelect XT				-		_	_
	IDT xGen Hybridization and capture of DNA Libraries	-						
	Illumina Nextera DNA Exome Kit				•			
	Illumina Nextera Flex Enrichment	•			•			
TARGET/EXOME CAPTURE	Illumina Nextera Rapid Capture Illumina TruSeq Exome				•			•
CAPTURE	NEBNext Direct Genotyping Solution				•			
	NuGEN Allegro Targeted Genotyping				<b>A</b>			
	NuGEN Allegro Targeted Genotyping  Roche Nimblegen SeqCap EZ for Illumina NGS				•			



## Biomek-Automated Nucleic Acid Sample Preparation Methods

	DEMONSTRATED METHOD	BIOMEK i7 HYBRID	BIOMEK i7/i5 MULTICHANNEL	BIOMEK i5 SPAN-8	BIOMEK FXP HYBRID	BIOMEK FXP/NXP MULTICHANNEL	BIOMEK NXP SPAN-8	BIOMEK 4000
DNA AND RNA PURIFICATION AND CLEAN UP/ SIZE SELECTION/ QUANTITATION/ NORMALIZATION/ POOLING	Agilent TapeStation 2200 Plate Setup for analysis of DNA Fragments			-				
	Beckman AMPure XP - PCR Purification		•		•	•	•	•
	Beckman CleanSEQ - Dye Terminator Removal	•	•		•	•		•
	Beckman RNAClean XP - Post cDNA Purification and Post-IVT cRNA Purification				•	•		•
	Beckman SPRIselect - Size Selection	•		•				
	COVARIS DNA Shearing Plate Setup for E220 Focused Ultrasonicators			•				
	Promega Wizard MagneSil Tfx Plasmid Purification System		•					
	qPCR/PCR Setup 384				-		•	•
	qPCR/PCR Setup 96				•		•	
	Quantitation/Normalization			•	•			•
	Roche Kapa Biosystems Library Quantification Kit - Illumina (96 and 384)			•	•		•	
	Sample Pooling for Multiplexing				•			
NUCLEIC ACID ISOLATION	Apostle Minimax™ High Efficiency cfDNA Isolation kit - Isolation of cFDA from Plasma	•	•					
	Beckman CosMCPrep - High/Low Copy Plasmid Purification				•	•		
	Beckman DNAdvance - DNA Isolation from Tissue	•	•		•	•		•
	Beckman Formapure XL DNA - DNA Isolation from FFPE Tissue		•	•	•			•
	Beckman Formapure XL RNA - RNA Isolation from FFPE Tissue	•	•					
	Beckman Formapure XL Total -DNA and RNA Isolation from a Single FFPE Tissue Sample	•	•					
	Beckman GenFind v2 - DNA Isolation from Blood	•	•			•	•	
	Beckman GenFind v3 - DNA Isolation from Blood, Cells & Serum	•		•				
	Beckman RNAdvance Blood - RNA Isolation from Blood	•			•		•	
	Beckman RNAdvance Cell - RNA Isolation from Cells						•	
	Beckman RNAdvance Viral - RNA Isolation from Viral samples							
	Beckman RNAdvance Tissue - RNA Isolation from Tissue	•			•			
	Mo Bio PowerMag Soil DNA Isolation Kit							•

## Method Key

- Demonstrated Method: This indicates that the method was developed for a sample preparation kit following a vendor's published manual protocol. Each one is tested with scientifically relevant samples and has yielded results that meet the kit vendor's specifications either in a customer lab (customer demonstrated) or in a Beckman Coulter Life Sciences Lab (Beckman demonstrated). Beckman Coulter makes no claims or warranties regarding the use or performance of these methods.
- Illumina Qualified Method: This indicates that Illumina's analysis of libraries prepared with the Biomek-automated method has shown the libraries to perform comparably to those prepared manually. This method is not an Illumina product, and Illumina does not directly support this product. Illumina makes no representations or warranties with respect to this product.
- ▲ Method created and provided by kit vendor
- ★TruSight Qualified method: These methods are co-developed by Illumina and the automation partner, undergo verification testing at Illumina and are alpha tested at a user site.

  The methods are installed and supported by the partner, but Illumina FAS will offer method training and will provide front line support. Technical support will also offer front line support in order to determine if the customer issue is the method/automation or the kit.

Biomek Automated Workstation Stations and methods are not intended or validated for use in the diagnosis of disease or other conditions. In some cases, method data was generated on pre-production automated workstations.

The automated methods identified are demonstrated methods only and are not validated by Beckman Coulter. Beckman Coulter makes no warranties of any kind whatsoever express or implied, with respect to this protocol, including but not limited to warranties of fitness for a particular purpose or merchantability or that the protocol is non-infringing. All warranties are expressly disclaimed. Your use of the method is solely at your own risk, without recourse to Beckman Coulter. Not intended or validated for use in the diagnosis of disease or other conditions. The demonstrated methods are available through Beckman Coulter automation field applications team.



© 2021 Beckman Coulter Life Sciences. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All additional trademarks used throughout this document are the property of their respective owners.