



Automated NGS Library Construction Methods for Biomek i-Series Workstations

DEMONSTRATED METHOD		BIOMEK i7 HYBRID	BIOMEK i7/i5 MULTICHANNEL	BIOMEK i5 SPAN-8
CANCER PANELS	Illumina AmpliSeq Cancer HotSpot Panel v2	●		●
	Illumina AmpliSeq Comprehensive Panel v3	●		●
	Illumina AmpliSeq Focus Panel	●		●
	Illumina TruSight Oncology 500 HT Panel	★		
	Illumina TruSight Tumor 170 Panel	●		●
	NEB NEBNext Direct Cancer HotSpot Panel	■		
	Illumina TruSight RNA Pan Cancer Panel	■		
DNA SEQUENCING	ATAC-Seq	■		
	IDT xGEN cfDNA and FFPE DNA Lib Prep Kit	■		
	Illumina AmpliSeq for Illumina NGS Kit	●		●
	Illumina DNA PCR-Free Prep kit	■		■
	Illumina Nextera DNA Flex Library Prep Kit	●		
	Illumina Nextera XT Library Prep Kit	●		
	Illumina TruSeq DNA PCR-Free	■		■
	Illumina TruSeq Nano DNA kit	■		
	NEB NEBNext Ultra II DNA Library Prep Kit for Illumina NGS	■		
	NEB NEBNext Ultra II FS DNA Library Prep Kit for Illumina NGS	■		
	NEB NEBNext Enzymatic Methyl Seq kit	■		
	Roche KAPA EvoPlus Library Prep kit		▲	
	Roche KAPA HyperPrep Kit for Illumina NGS Library Prep	■▲		
	Roche KAPA HyperPlus Kit for Illumina NGS Library Prep	■▲		
	Swift Biosciences Accel - NGS 2S Plus DNA Library Kit for Illumina NGS	■		
Twist Library Preparation Kit (Mechanical & enzymatic fragmentation)	■		■	
HLA TYPING	GenDx NGSgo Kit	■		
	Immucor MIA FORA Flex HLA Kit	▲	▲	
RNA SEQUENCING	Illumina AmpliSeq for Illumina NGS kit	●		●
	Illumina RNA Prep with Enrichment Kit	■		
	Illumina TruSeq Stranded mRNA Library Prep Kit	■		
	Illumina TruSeq Stranded Total RNA Library Prep Kit	■		
	Illumina Stranded mRNA Library Prep kit	■		
	Illumina Stranded Total RNA Library Prep kit	■		
	Lexogen QuantSeq 3' mRNA -Seq Library Prep Kit	■		
	NEBNext ARTIC SARS CoV-2 FS Library Prep Kit	■		
	NEB NEBNext Ultra II Directional RNA Library Prep Kit for Illumina NGS	■		
	NuGEN Universal Plus mRNA-Seq Library Preparation Kit	■		
	Oxford Nanopore Midnight Covid Seq Kit	■		
	Qiagen QIAseq DIRECT SARS-Cov-2 Kit	■		■
	Qiagen QIAseq miRNA UDI Library kit	■		
	Roche KAPA mRNA HyperPrep Kit	■		
	Roche KAPA RNA HyperPrep Kit with RiboErase (HMR)	■		
SINGLE-CELL SEQ	10X Genomics Single Cell 3' kit- Library Preparation	■		
	NEB NEBNext Single Cell/ Low Input RNA Library Prep Kit	■		
	Takara SMART-Seq HT Kit	■		
TARGET/EXOME CAPTURE	Agilent SureSelect XT Kit	■		
	IDT xGen Hybridization Capture Kit	■		
	Illumina Nextera Flex Enrichment Kit	●		
	Illumina TruSeq Exome Kit	■		
	NEBNext Direct Genotyping Solution kit for Illumina	■		
	Twist Fast Hybridization and Wash Kit	■		
Twist Standard Hybridization and Wash Kit	■			



Automated Nucleic Acid Sample Preparation Methods for Biomek i-Series Workstations

DEMONSTRATED METHOD		BIOMEK 17 HYBRID	BIOMEK 17/15 MULTICHANNEL	BIOMEK 15 SPAN-8
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 kit	■	■	
	Beckman CleanSEQ - Dye Terminator Removal kit	■	■	
	Beckman SPRIselect - Size Selection kit	■		■
	COVARIS DNA Shearing Plate Setup for E220 Focused Ultrasonicators	■		■
	Promega Wizard MagneSil Tfx Plasmid Purification System		■	
	Quantitation/Normalization procedures	■		■
	Roche Kapa Biosystems Library Quantification Kit - Illumina (96 and 384)	■		■
NUCLEIC ACID ISOLATION	Apostle Minimax™ High Efficiency cfDNA Isolation kit - Isolation of cfDNA from Plasma	■	■	
	Beckman DNAdvance kit - DNA Isolation from Tissue System	■	■	
	Beckman FormaPure XL DNA Kit - DNA Isolation from FFPE Tissue	■	■	■
	Beckman FormaPure XL RNA Kit - RNA Isolation from FFPE Tissue	■	■	
	Beckman FormaPure XL Total Kit -DNA and RNA Isolation from a Single FFPE Tissue Sample	■	■	
	Beckman GenFind v2 Kit- DNA Isolation from Blood	■	■	
	Beckman GenFind v3 Kit - DNA Isolation from Blood, Cells & Serum	■	■	■
	Beckman RNAdvance Blood Kit- RNA Isolation from Blood	■	■	
	Beckman RNAdvance Cell Kit- RNA Isolation from Cells		■	
	Beckman RNAdvance Viral Kit- RNA Isolation from Viral samples		■	
	Beckman RNAdvance Tissue Kit- RNA Isolation from Tissue	■	■	
	Cytiva Viral Pathogen RNA 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.

© 2022 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.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://www.beckman.com)

2022-GBL-EN-100049-v1