



# Biomek-Automated NGS Library Construction Methods

DEMONSTRATED METHOD		BIOMEK FX <sup>P</sup> HYBRID	BIOMEK FX <sup>P</sup> MULTI-CHANNEL	BIOMEK FX <sup>P</sup> SPAN 8	BIOMEK NX <sup>P</sup> SPAN 8	BIOMEK NX <sup>P</sup> MULTI-CHANNEL	BIOMEK 4000
CANCER PANELS	Illumina TruSight <sup>®</sup> Cancer						■
	Illumina TruSight <sup>®</sup> Tumor 15						■
DNA SEQUENCING	Illumina Nextera <sup>®</sup> XT	●					
	Illumina TruSeq <sup>®</sup> DNA PCR-Free	■					■
	Illumina TruSeq <sup>®</sup> Nano DNA	■			■		■
	Illumina TruSeq <sup>®</sup> Custom Amplicon Low Input Kit	●					
	KAPA Hyper Prep Library Prep Kit for Illumina NGS	▲					
	KAPA HyperPlus Library Prep Kit for Illumina NGS	▲					
	NEBNext <sup>®</sup> Ultra DNA for Illumina NGS (ChIP-seq and HLA)	■					
	NEBNext <sup>®</sup> Ultra DNA II Kit for Illumina NGS	■					
	Rubicon Genomics ThruPLEX <sup>®</sup> Plasma-seq Kit for Illumina NGS	■					
	Swift Biosciences Accel - NGS <sup>®</sup> 2S Plus DNA Library Kit for Illumina NGS	■					
HLA TYPING	Illumina TruSight <sup>®</sup> HLA Sequencing Panel v2				●		
	Immucor Mia Fora NGS	▲			▲		▲
	Omixon Holotype HLA X2						■
RNA SEQUENCING	Illumina TruSeq <sup>®</sup> RNA v2	■					
	Illumina TruSeq <sup>®</sup> RNA Access	■					
	Illumina TruSeq <sup>®</sup> Stranded mRNA	●					
	Illumina TruSeq <sup>®</sup> Stranded Total RNA	●					
	NEBNext <sup>®</sup> Small RNA Kit for Illumina NGS						■
	NEBNext <sup>®</sup> Ultra Directional RNA Library Kit for Illumina NGS	■					
TARGET/EXOME CAPTURE	Agilent HaloPlex <sup>™</sup> Target Enrichment - Ion Torrent	■			■		■
	Agilent SureSelect XT <sup>®</sup>	■					
	Epicentre ScriptSeq <sup>®</sup> Complete Gold Low Input	●					
	Illumina Nextera <sup>®</sup> Rapid Capture	●					
	Illumina TruSeq <sup>®</sup> Exome	■					
	Illumina TruSeq <sup>®</sup> Rapid Exome	■					
	Nimblegen SeqCap EZ <sup>®</sup> for Illumina NGS	■					



# Biomek-Automated Nucleic Acid Sample Preparation Methods

DEMONSTRATED METHOD		BIOMEK FX <sup>P</sup> HYBRID	BIOMEK FX <sup>P</sup> MULTI-CHANNEL	BIOMEK FX <sup>P</sup> SPAN 8	BIOMEK NX <sup>P</sup> SPAN 8	BIOMEK NX <sup>P</sup> MULTI-CHANNEL	BIOMEK 4000
DNA AND RNA PURIFICATION AND CLEAN UP/ SIZE SELECTION/ QUANTITATION/ NORMALIZATION/ POOLING	AMPure XP - PCR Purification	■	■		■	■	■
	CleanSEQ - Dye Terminator Removal	■	■			■	■
	KAPA Biosystems Library Quantification Kit - Illumina	■			■		
	qPCR setup 384	■			■		■
	qPCR setup 96	■			■		■
	Quantitation/Normalization	■			■		■
	RNAClean XP - Post cDNA Purification and Post-IVT cRNA Purification	■	■			■	■
	SPRIselect for DNA Size Selection	■			■		■
Sample Pooling for Multiplexing	■						
NUCLEIC ACID ISOLATION	CosMCPrep - High/Low Copy Plasmid Extraction	■	■			■	
	DNAdvance - DNA Isolation from Mammalian Tissue	■	■			■	■
	FormaPure DNA - DNA Isolation from FFPE Tissue						■
	FormaPure - RNA Isolation from FFPE Tissue	■		■	■		
	Mo Bio PowerMag Soil DNA Isolation Kit						■
	RNAdvance Tissue	■		■	■		
	RNAdvance Cell v2 - RNA Isolation from Tissues	■	■	■	■	■	

## Method Key

- Demonstrated Method developed for a sample preparation kit following a vendor's published manual protocol. Each one is tested with 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 regarding the use or performance of these methods.
- Methods currently in the data generation phase of development.
- Illumina Qualified Method, which 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 support this product. Illumina makes no representations or warranties with respect to this product.
- ▲ Method created and provided by kit vendor.



Methods are intended for molecular biology research applications. They are not intended, verified or validated, for use in the diagnosis of disease or other conditions.

\* All 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](http://beckman.com).

B2014-14915 DS-18966B beckman.com © 2015 Beckman Coulter, Inc. PRINTED IN U.S.A.