

Introduction

Next-generation sequencing (NGS) target enrichment protocols are in high demand for preparation of Illumina sequencing-ready libraries from samples of varying quality and input. Automation of these complex protocols eliminates the need for highly trained specialists while considerably improving the NGS library prep turnaround time, quality, and reproducibility. A fully automated walk-away protocol on the Agilent Magnis NGS Prep system was developed and optimized for a wide range of DNA inputs (10 to 200 ng) and sample types (for example, high quality, formalin-fixed paraffin-embedded [FFPE], circulating tumor DNA samples). It includes different shearing methods within the automated protocol (mechanical or enzymatic fragmentation) using 192 unique dual indexes (UDI) to minimize the effects of possible index hopping and contamination. Further, incorporation of inline duplex molecular barcodes filters out PCR or sequencing errors by making consensus calls using barcode information. The automated workflow on Magnis, referred to as the SSEL-DNA-XTHS2-ILM protocol, is easily set up from pre-aliquoted, single-use reagents to minimize contamination risk. It delivers up to eight target-enriched NGS libraries per run, in nine to 10 hours, without the need of user intervention. To optimize library yields, catalog and custom small, medium, and large probes can be used and PCR cycles can easily be adjusted from the touch screen. The expanded Magnis SureSelect XT HS2 DNA protocol now includes dual indexing capabilities and a new method to process initial and poor-quality FFPE RNA. This secondary automated procedure, the SSEL-RNA-XTHS2-ILM protocol, can be set up and run separately using RNA that is first converted to cDNA and then further processed using similar steps as the SSEL-DNA-XTHS2-ILM protocol to create a final NGS library.

Methods

SureSelect XT HS2 DNA libraries were generated from over 80 runs (640 libraries) on six Magnis instruments using high-quality HapMap and FFPE DNA at 10 to 200ng inputs with small, medium, and large bait panels. Ultrasonication methods were used to manually shear the DNA prior to the runs, and short and long fragments were created by changing the shearing time. In addition, the enzymatic fragmentation option was used and is programmed in the protocol to shear the genomic DNA on-deck. All sequencing results shown in Figure 2 are from libraries generated from DNA using the large capture V8 probe set and were down sampled to 100X raw sequencing depth. Single nucleotide polymorphism (SNP) calling was completed using the Agilent SureCall Software (Figure 3). Inter- and intrarun contamination assays were performed using 50 ng of input DNA for all eight samples (15,000 copies). Every other sample was spiked with one million copies of Alien 6 DNA fragment (239 bp) which contains a specific DNA sequence absent from the human genome and can be explicitly detected using a sensitive qPCR assay. A Taqman qPCR assay specific to the Alien 6 DNA was used for the pre- and postcapture samples. It was used postrun to rule out intrarun contamination and in the following run to rule out inter-run contamination. Quantification of copies per μ L sample was completed using a standard curve (spiked samples were pre-diluted at 100X; data not shown). The sequencing data from each pool was aligned to all 192 indexes. When fewer than eight samples per run were sequenced, alternate samples were sequenced. Any sample-to-sample or index contamination would likely be higher over a shorter distance and thus more likely to contaminate neighboring wells. Therefore, sequencing alternate samples allowed for capture of contamination throughout the process, from mixing the index pairs and plating into the strip tubes, through the Magnis library prep and target enrichment steps, sample collection, and quantification (Figure 4).

