APPLICATION NOTE

Optimized protocol for human whole-exome sequencing

Introduction

Since the first human genome was sequenced in 2003, new sequencing technologies have been evolving rapidly. Even though genome sequencing costs have decreased drastically over the years due to the emergence of next-generation sequencing (NGS), it has yet to become affordable and efficient enough for routine use in diagnostics research. In contrast to whole-genome sequencing (WGS), whole-exome sequencing (WES) and targeted sequencing provide a balance between cost and benefit. WES targets approximately 2% of the whole genome, corresponding to protein-coding genes [1]. It is a proven and cost-effective method to detect disease-causing variants and discover gene targets. This application note describes a workflow solution for NGS library preparation of human whole-exome samples using the Invitrogen™ Collibri™ ES DNA Library Prep Kit for Illumina[™] Systems with unique dual indexes (UDIs), from library preparation to sequencing (Figure 1).

Methods

Samples

Well-characterized human genomic DNA (gDNA) samples obtained from the GM12878 human cell line (Ceph/Utah pedigree) were purchased from the Coriell Institute for Medical Research (Cat. No. NA12878) for use in whole-exome library preparation.

Library generation

The gDNA samples were converted into sequencing-ready libraries via enzymatic shearing (ES) for fragmentation, using the Collibri ES DNA Library Prep Kit for Illumina Systems with UDIs. The standard protocol of the Collibri ES DNA Library Prep Kit for Illumina Systems was used with the following modifications: 31 μ L of EDTA-free double-stranded DNA (recommended input of 50 ng). A buffer/DNA mixture (36 μ L) was prepared by adding 5 μ L of 10X Fragmentation and dA-Tailing Buffer to the DNA mixture. The 5X Fragmentation and dA-Tailing Enzyme Mix (14 μ L) was then added to the buffer/DNA mixture for a 50 μ L total volume.

Fragmentation was performed for 20 min at 37°C. The adaptor-ligated library was purified according to a customized cleanup protocol (**Pub. No. MAN0025489**) for the Collibri ES DNA Library Prep Kit for Illumina Systems. The library was then amplified for 8 PCR cycles before the enrichment step and subsequent cleanup (**Pub. No. MAN0025490**). The complete manual for the human whole-exome library preparation protocol described in this application note is also available (**Pub. No. MAN0025534**).



Figure 1. NGS workflow for human whole-exome sequencing.



Table 1. Composition of hybridization kits (Long and Fast hybridization protocols).

F	ast hybridization protoc	col
Name	Cat. No.	Kit composition
	101175	Twist fast hybridization reagents:
		Fast hybridization mix
Todas Forest I belondations and		Hybridization enhancer
Twist Fast Hybridization and Wash Kit (2 boxes),		Amplification primers
96 reactions		2. Twist fast wash buffers:
		Fast binding buffer
		Fast wash buffer 1
		Wash buffer 2
Twist Comprehensive Exome, 96 reactions	102033 (1.1 rev)/ 101953 (1 rev)	Comprehensive panel
Twist Comprehensive Exome, 12 reactions	102032 (1.1 rev)/ 101952 (1 rev)	Comprehensive panel
Twist Comprehensive Exome, 2 reactions	102031(1.1 rev)/ 101951 (1 rev)	Comprehensive panel
Twist Binding and		Streptavidin binding beads
Purification Beads (96 reactions/12 reactions)	100984/100983	DNA purification beads
Twist Universal Blockers	400707/400570	Universal blockers
(96 reactions/12 reactions)	100767/100578	Blocker solution

Long hybridization protocol			
Name	Cat. No.	Kit composition	
	101026	Twist hybridization reagents:	
		Hybridization mix	
T. dat III de delle etter en el		Hybridization enhancer	
Twist Hybridization and Wash Kit (2 boxes),		Amplification primers	
96 reactions		2. Twist wash buffers:	
		Binding buffer	
		Wash buffer 1	
		Wash buffer 2	
Twist Comprehensive Exome, 96 reactions	102033 (1.1 rev)/ 101953 (1 rev)	Comprehensive panel	
Twist Comprehensive Exome, 12 reactions	102032 (1.1 rev)/ 101952 (1 rev)	Comprehensive panel	
Twist Comprehensive Exome, 2 reactions	102031 (1.1 rev)/ 101951 (1 rev)	Comprehensive panel	
Twist Binding and Purification Beads (96 reactions/12 reactions)		Streptavidin binding beads	
	100984/100983	DNA purification beads	
Twist Universal Blockers	100707/100570	Universal blockers	
(96 reactions/12 reactions)	100767/100578	Blocker solution	

Enrichment

The WGS libraries were quantified using the Invitrogen™ Qubit™ Fluorometer and pooled. The exome targets were then enriched using a synthetic probe panel (Twist Human Core Exome Panel, from Twist Bioscience). Two separate hybridization reactions were performed:

- "Fast" hybridization—using the Twist Fast Hybridization and Wash Kit; 2 hr target enrichment
- "Long" hybridization—using the Twist Standard Hybridization and Wash Kit; overnight (~16 hr) target enrichment

The composition of each hybridization kit is noted in Table 1. The enriched libraries were amplified for 8 PCR cycles using the 2X Library Amplification Master Mix from the Collibri ES DNA Library Prep Kit for Illumina Systems instead of the recommended KAPA™ HiFi HotStart ReadyMix from Roche (Table 2). The DNA purification beads from Twist hybridization kits can be replaced with Invitrogen™ Dynabeads™ MyOne™ Streptavidin T1 magnetic beads (Cat. No. 65601), if required.

Library quantification and sequencing

The final concentrations of the sequencing libraries were determined using the Invitrogen™ Collibri™ Library Quantification Kit. Libraries were sequenced on the Illumina™ NovaSeq™ 6000 Sequencing System by paired-end sequencing of 2 x 150 bp.

NGS analysis

The Illumina™ BaseSpace™ Enrichment App v3.1.0 was used to align a subsample of 30M passing filter (PF) clusters (60M reads) from each sample to the hg19 reference genome. The enrichment parameters were calculated using the Picard CollectHsMetrics tool. Variant calling in the Variant Calling Assessment Tool v4.0.2 was limited by the target regions. Gold-standard data for NA12878 from the Genome in a Bottle (GIAB) consortium hosted by NIST were used to determine the precision and recall parameters.

Results

Some of the main parameters that are usually calculated in WES include enrichment efficiency, coverage of enriched targets, and single nucleotide variant (SNV) analysis. In this study, we were able to compare the parameters of the Long and Fast enrichment protocols so that customers may identify and choose the performance they desire for their experiments.

Table 2. Recommended protocol optimization for library generation using Collibri ES DNA Library Prep Kits.

Step	Standard recommendation		Recommended changes for WES samples		
Removal of EDTA from DNA	(If needed)		If the DNA is ≤1 ng, no need to remove EDTA from DNA samples		
samples	Input: 1-500 ng		Input: 50 ng		
	On ice or a cooling rack, assemble the fragmentation and dA-tailing reaction for each DNA sample in a sterile 0.2 mL thin-wall PCR tube. Add the reagents in the order given.		On ice or a cooling rack, assemble the fragmentation and dA-tailing reaction for each DNA sample in a sterile 0.2 mL thin-wall PCR tube. Add the reagents in the order given.		
	Component	Volume	Component	Volume	
	Double-stranded DNA (1-500 ng)	XμL	Double-stranded DNA	ΧμL	
	10 mM Tris-HCl, pH 7.5-8.5	to 40 µL	10 mM Tris-HCI, pH 7.5-8.5	to 36 µL	
	10X Fragmentation and dA-Tailing Buffer (blue)	5 μL	10X Fragmentation and dA-Tailing Buffer (blue)	5 μL	
	Total volume (light blue mixture)	40 μL	Total volume (light blue mixture)	36 µL	
	Add 5X Fragmentation and dA-Tailing Enzyme Mix to the sample.		Add 5X Fragmentation and dA-Tailing Enzyme Mix to the sample.		
2. DNA fragmentation and addition of dA tails	Component	Volume	Component	Volume	
addition of da tails	Buffer-DNA mixture from above (light blue mixture)	40 μL	Buffer-DNA mixture from above (light blue mixture)	36 µL	
	5X Fragmentation and dA-Tailing Enzyme Mix (colorless)	10 μL	5X Fragmentation and dA-Tailing Enzyme Mix (colorless)	14 µL	
	Total volume (light blue mixture)	50 μL	Total volume (light blue mixture)	50 μL	
	Recommended fragmentation time		Recommended fragmentation time		
	(at 37°C)		(at 37°C)		
	Desired fragment size	Fragmentation time (optimization range)	Fragment for 20 minutes		
	150–300 bp	20 min (20–30 min)			
	300-500 bp	10 min (10–20 min)			
	500–700 bp	5 min (5–10 min)			
3. Adaptor ligation	Described in the "Dual-Indexed Adaptor ligation" section of the workflow provided in the Collibri ES DNA Library Prep Kit for Illumina User Guide		Described in the "Dual-Indexed Adaptor ligation" section in the Collibri ES DNA Library Prep Kit for Illumina User Guide		
4. Post-ligation double-sided size selection	Double-sided size selection to match target insert size		Check product page for the protocol (Pub. No. MAN0025489)		
5. PCR amplification of the DNA library	The number of PCR cycles depends on the starting amount of input DNA		Amplify the enriched library for 8 PCR cycles		
6. Amplified DNA library purification	Post-amplification cleanup of the DNA library		Check product page for the protocol (Pub. No. MAN0025490)		
7. Human exome enrichment	Not applicable		Target enrichment protocol		
8. Post-capture library amplification with Library Amplification Master Mix			Component	Volume	
	Not applicable		Enriched library from the Target Enrichment and Amplification Beads slurry (brown)	20 μL	
			2X Collibri Library Amplification Master Mix (blue)	25 μL	
			Collibri Primer Mix	5 μL	
			Total volume (brown mixture)	50 μL	
			Amplify the enriched DNA library for 8 PCR cycles		
9. Purification of the amplified bost-capture libraries	Not applicable		Purification of the amplified post-capture libraries		

Single-nucleotide variants

Accurate SNV determination is one of the key parameters in WES. Using the modified Collibri ES DNA library prep protocol described in Table 2 in combination with the Twist Bioscience enrichment kits (for Long and Fast hybridization protocols, Table 1), we were able to reach an SNV precision greater than 99%. An SNV recall of 96.8% was achieved using the Fast hybridization protocol, whereas the Long enrichment protocol yielded an SNV recall of 98.1% (Figure 2).

Coverage and enrichment

Efficient target enrichment and high coverage are key to successful WES. In addition, high coverage uniformity is required for efficient sequencing so that the targeted regions are covered as uniformly as possible. In this study, the enrichment and coverage parameters were calculated using the 150 bp padded target. High coverage uniformity was achieved using a modified Collibri ES DNA library prep protocol in combination with the Twist Bioscience enrichment module. The results showed 96.7% coverage uniformity when using the Fast hybridization protocol and 97.8% when target enrichment was performed using the Long hybridization protocol (Figure 3).

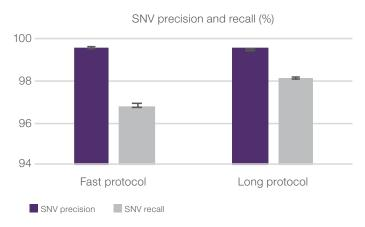
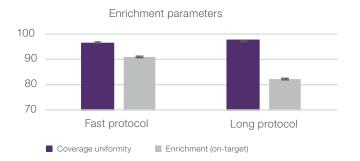


Figure 2. SNV precision and recall for the Fast and Long enrichment protocols.



The overall target enrichment of 91.4% was achieved when using the Fast enrichment protocol, compared to the 82.2% target enrichment achieved with the Long protocol. We also observed that the percentage of targets with coverage greater than 20x is 98.2% when using the Long enrichment module, but only 96.2% when the Fast hybridization module is used (Table 4).

Table 3. Comparison of whole-exome sequencing performance parameters for Fast and Long enrichment protocols. All of the parameters were calculated using 16 repetitions of the same sample preparation.

Parameter	Fast	Long	
Panel	Twist Comprehensive Exome		
Panel size	36.8 Mb		
Probe size	120 bp		
Padded read enrichment (on-target)	91.4 ± 0.2%	82.5 ± 0.4%	
Coverage at 20x	97.1 ± 0.1%	98.1 ± 0.1%	
Uniformity of coverage	96.7 ± 0.1%	97.8 ± 0.1%	
Read depth per sample	30M PF clusters		
SNV precision	99.6 ± 0.0%	99.6 ± 0.0%	
SNV recall	96.8 ± 0.1%	98.1 ± 0.1%	

Summary

Strong target coverage and sensitive variant detection for the human whole exome were accomplished using two different enrichment methods (Long and Fast) and a modified protocol for the Collibri ES DNA Library Prep Kit for Illumina Systems. Over 97% of exome bases were covered with only 20x coverage. Enrichment quality corresponded to the percentage of on-target reads, and all of the samples were sufficiently enriched to detect single-nucleotide polymorphisms (SNPs). The optimized protocol for the Collibri ES DNA Library Prep Kit for Illumina Systems, when used in combination with the Twist Bioscience hybridization kit, was extremely precise (>99%) and resulted in outstanding recall rates (>96%) for detecting SNPs.

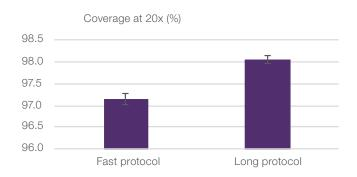


Figure 3. Coverage and enrichment parameters for the Fast and Long enrichment modules.

invitrogen

Conclusions

- The optimized method described in this application note for the Collibri ES DNA Library Prep Kit enables highly sensitive detection of variants for human whole-exome samples.
- Uniformity and high on-target rates show the efficiency of targeted sequencing using the optimized Collibri ES DNA Library Prep Kit.
- The Collibri Library Amplification Master Mix containing Invitrogen™ Platinum™ SuperFi™ DNA Polymerase provides uniform exome amplification and minimal error rates similar to those of PCR-free methods; this is crucial for detecting SNPs accurately.

Reference

 Suwinski P, Ong C, Ling MHT et al. (2019) Advancing personalized medicine through the application of whole exome sequencing and big data analytics. Front Genet 10:49. doi: 10.3389/fgene.2019.00049

Ordering information

Product	Quantity	DNA index type	Cat. No.
	24 preps	CDI*	A38605024
Collibri ES DNA Library Prep Kit for Illumina Systems, with indexes	96 preps	CDI	A38607096
	24 preps	UDI* set A (1-24)	A38606024
		UDI set B (25-48)	A43605024
		UDI set C (49-72)	A43606024
		UDI set D (73-96)	A43607024
	96 preps	UDI set A-D (1-96)	A38607196
Collibri ES DNA Library Prep Core Kit for Illumina Systems	96 preps	_	A38607096W
Collibri Library Quantification Kit	100 reactions	_	A38524100
Collibri Library Quantification Kit	500 reactions	_	A38524500
Collibri Library Amplification Master Mix with Primer Mix**	50 reactions	_	A38540050
Collibri Library Amplification Master Mix with Primer Mix**	250 reactions	_	A38540250

^{*} CDI = combinatorial dual indexes, UDI = unique dual indexes.



^{**} Collibri Library Amplification Master Mix with Primer Mix can be purchased separately for use with the PCR-free version of the Collibri DNA Library Prep Kit.