

NGS Constitutional Disease Research Workflow

1

Ordering a Catalog Bait Library or Creating a Custom Panel



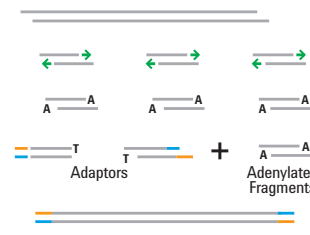
Order an NGS catalog panel or easily design a custom panel in minutes using Agilent's SureDesign web application

SureSelect Hybridization-Based Target Enrichment

Ligation-based

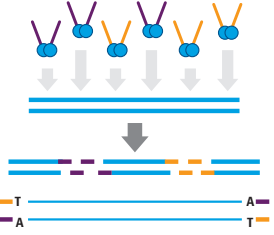
Transposase-based

Preparation of DNA Libraries



1. Shear DNA into smaller fragments
2. Repair fragment ends
3. Adenylate fragments
4. Ligate adaptors
5. Amplify pre-capture library

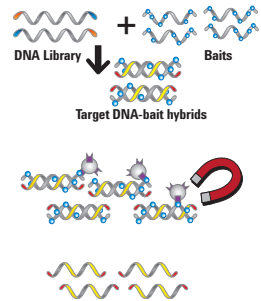
Preparation of DNA Libraries



Genomic DNA is enzymatically fragmented and tagged with the adaptors
Adaptor-tagged fragments are amplified to generate the pre-capture libraries

3

Hybridization & Capture

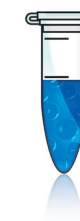


1. DNA library is allowed to hybridize with biotinylated SureSelect RNA baits (90 min for QXT, 16 hr for XT or XT2)
2. DNA-RNA hybrids are pulled down using streptavidin-coated magnetic beads
3. Stringent washes are applied to eliminate non-specific targets

HaloPlex Next-Generation PCR

2

Digest & Denature Sample DNA



Digest genomic DNA with 16 restriction enzymes and QC the reactions using a 4200 TapeStation or 2100 Bioanalyzer System

3

Hybridize Probes to DNA Targets



Mix the HaloPlex probes, indexing primer cassettes and digested DNA; hybridize for 3 hours (up to 500 kb, if >500 kb hyb time = 16 hours)

4

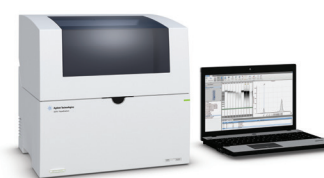
Amplify Targets



Amplify enriched and ligated DNA fragments using a single primer pair

5

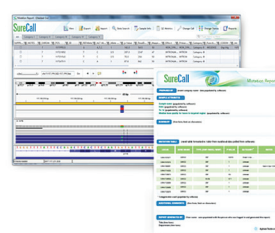
Final Quality Control



Quantify and QC prepared libraries using a 4200 TapeStation or 2100 Bioanalyzer System

6

Sequence & Analysis



Sequence libraries, analyze and obtain report of mutations using SureCall data analysis software



For Research Use Only.
Not for use in diagnostic procedures.

PR7000-0187
© Agilent Technologies, Inc. 2016
Printed in USA, May 19, 2016
5991-5909EN