



SureSelect

Your Complete Target
Enrichment Solution

Christina Chiu

Global Product Manager

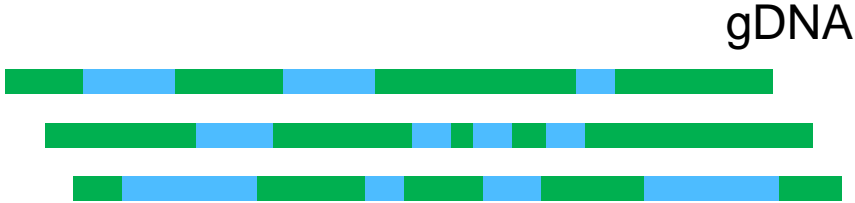
Focus on what matters

Target Enrichment



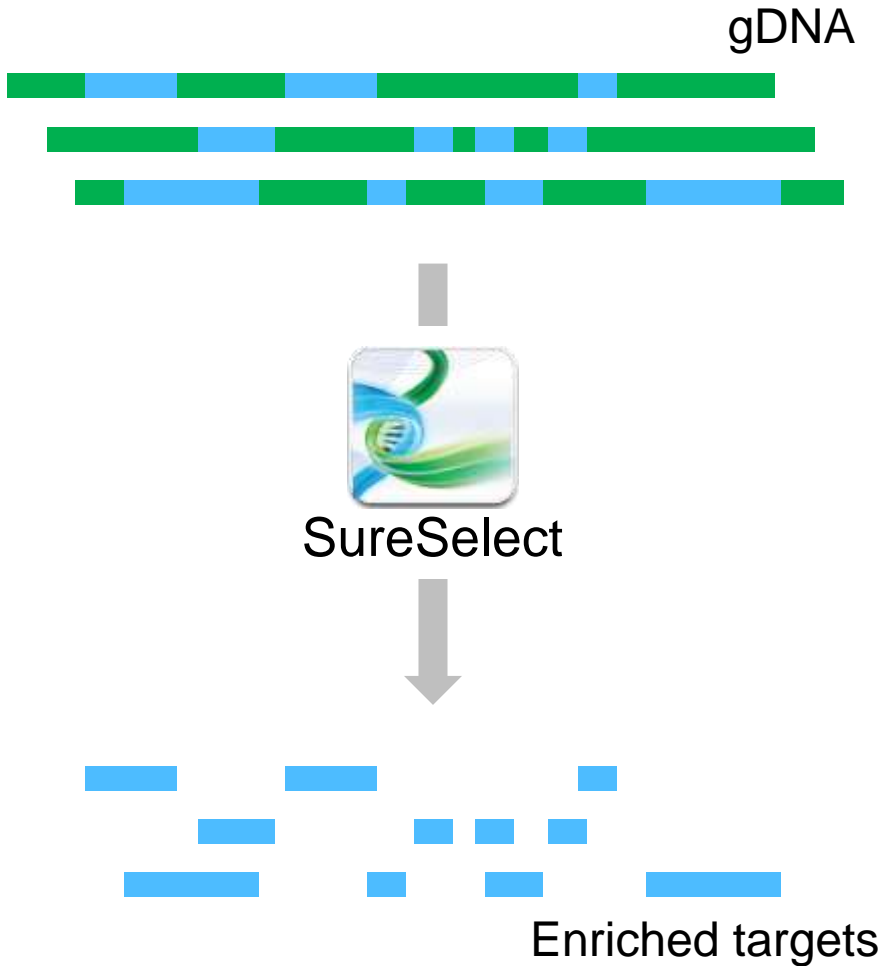
Focus on what matters

Target Enrichment



Focus on what matters

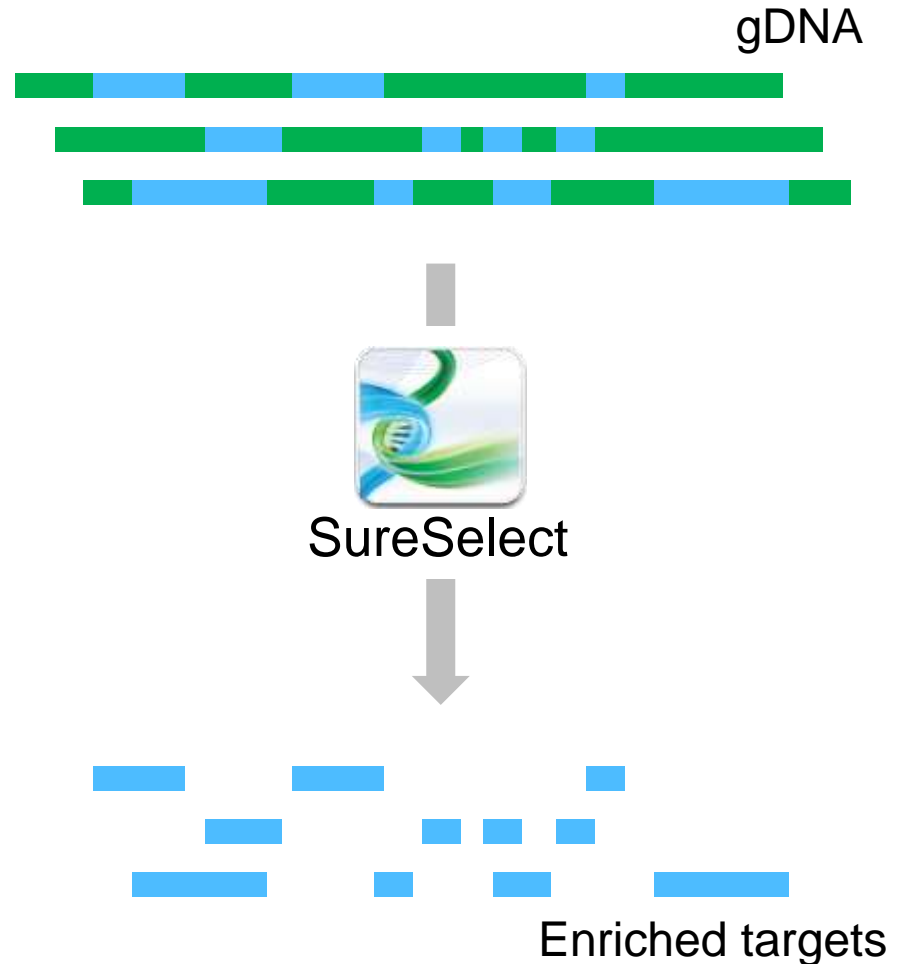
Target Enrichment



Focus on what matters

Target Enrichment

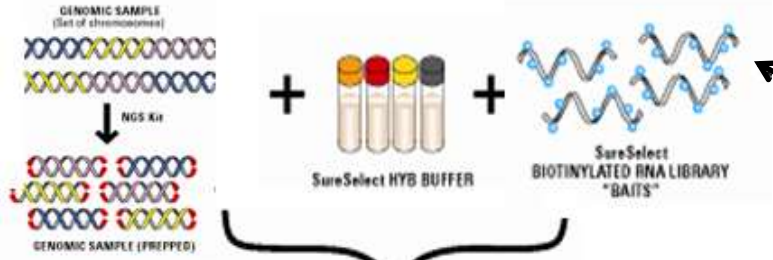
- Focus on a subset of the genome/transcriptome
- Achieve greater depth in targeted regions – discover more mutations, SNPs and Indels
- Comprehensive detection of genome-wide CNVs and LOH
- Lower sequencing costs and obtain faster time to result



SureSelect Technology

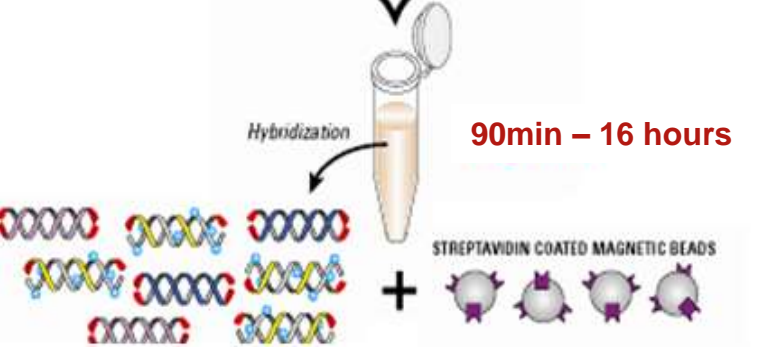
Simple Target Enrichment Workflow

Library prep
Hyb set-up

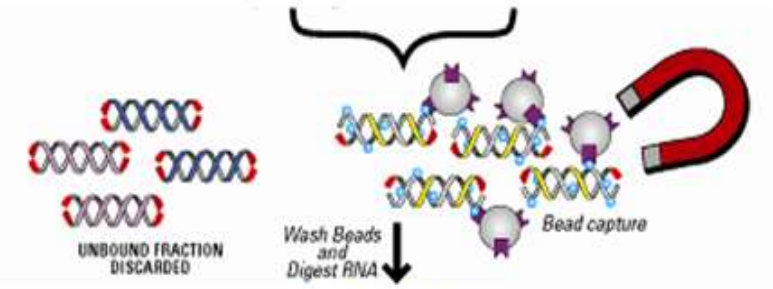


- Baits:**
- cRNA probes
 - Long (120bp)
 - Biotin labeled
 - User-defined

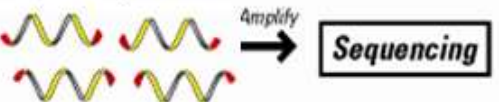
Hybridization



Capture



Wash / Elution / Amp



SureSelect

The Leader In Target Enrichment

COMPLETE
Solution

Best
PERFORMANCE

FLEXIBLE
Designs

SureSelect – Complete Solution

The Leader In Target Enrichment

Library
Prep Kits



TE Kits



Automation

DNA

XT

XT2

QXT

Exomes (Clinical Research Exome, Human, Non-human)

Panels (Inherited Disease, DNA Kinome)

Human Methyl-Seq

Whole Genome

Custom Solutions

RNA

Stranded RNA RNA-Seq

Stranded RNA RNA capture

Panels (RNA Kinome)



OR



SureSelect – Complete Solution

Complete workflow solution from sample to data

1. Create a design



SureDesign – human or nonhuman organisms

2. Qualify samples



Tapestation DIN or NGS FFPE QC

3. Enrich libraries



SureSelect Workflows

4. Quantify and qualify libraries



2100 BioAnalyzer or 4200 TapeStation

5. Analyze and Interpret



SureCall – single sample, trio or pair analysis and Cartagenia Bench

SureSelect

The Leader In Target Enrichment

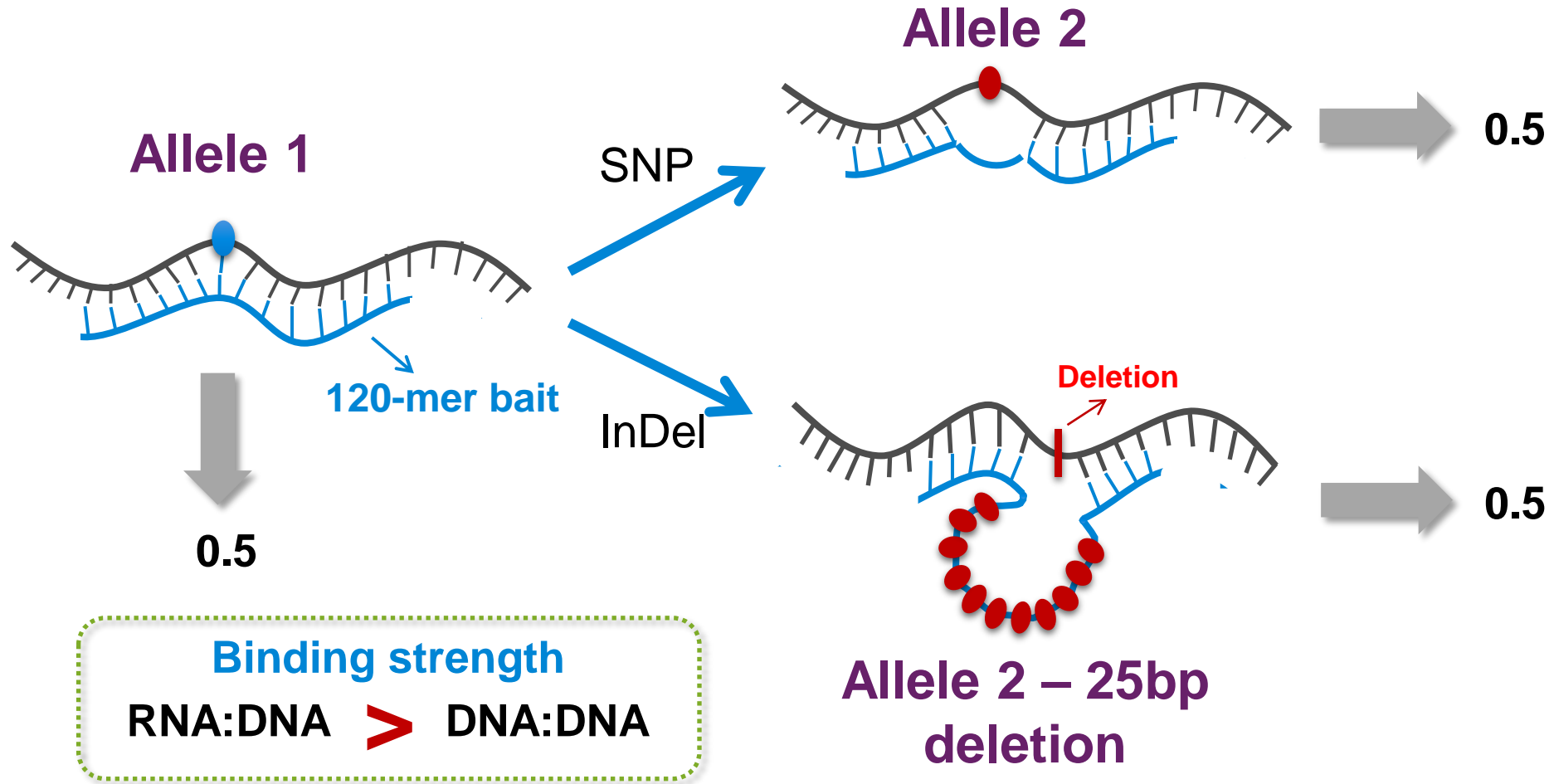
COMPLETE
Solution

Best
PERFORMANCE

FLEXIBLE
Designs

SureSelect – Best Performance

Longer baits = Better sensitivity



SureSelect – Best Performance

The most published target enrichment platform

- Over 2000 publications describe research enabled by Agilent NGS technology

SureSelect

The Leader In Target Enrichment

COMPLETE
Solution

Best
PERFORMANCE

FLEXIBLE
Designs

SureSelect – Flexible Designs

Customize your target enrichment with Custom SureSelect

**Capture only
what you want**

Custom content from <200Kb up to
24Mb+

Cost effective

Minimize enrichment and sequencing
costs

**Faster time from
sample to results**

Greater sample throughput, easier data
analysis

SureSelect – Flexible Designs

Customize your target enrichment with Custom SureSelect



Create custom designs specific to your research needs

- Choose from pre-defined content or create your own
- Review Agilent catalog and custom designs for purchase
- Collaborate within your workgroup or other researchers



Assistance with complex designs

- Access to Field Application Scientists
- Bioinformatics, commercial and industry experience