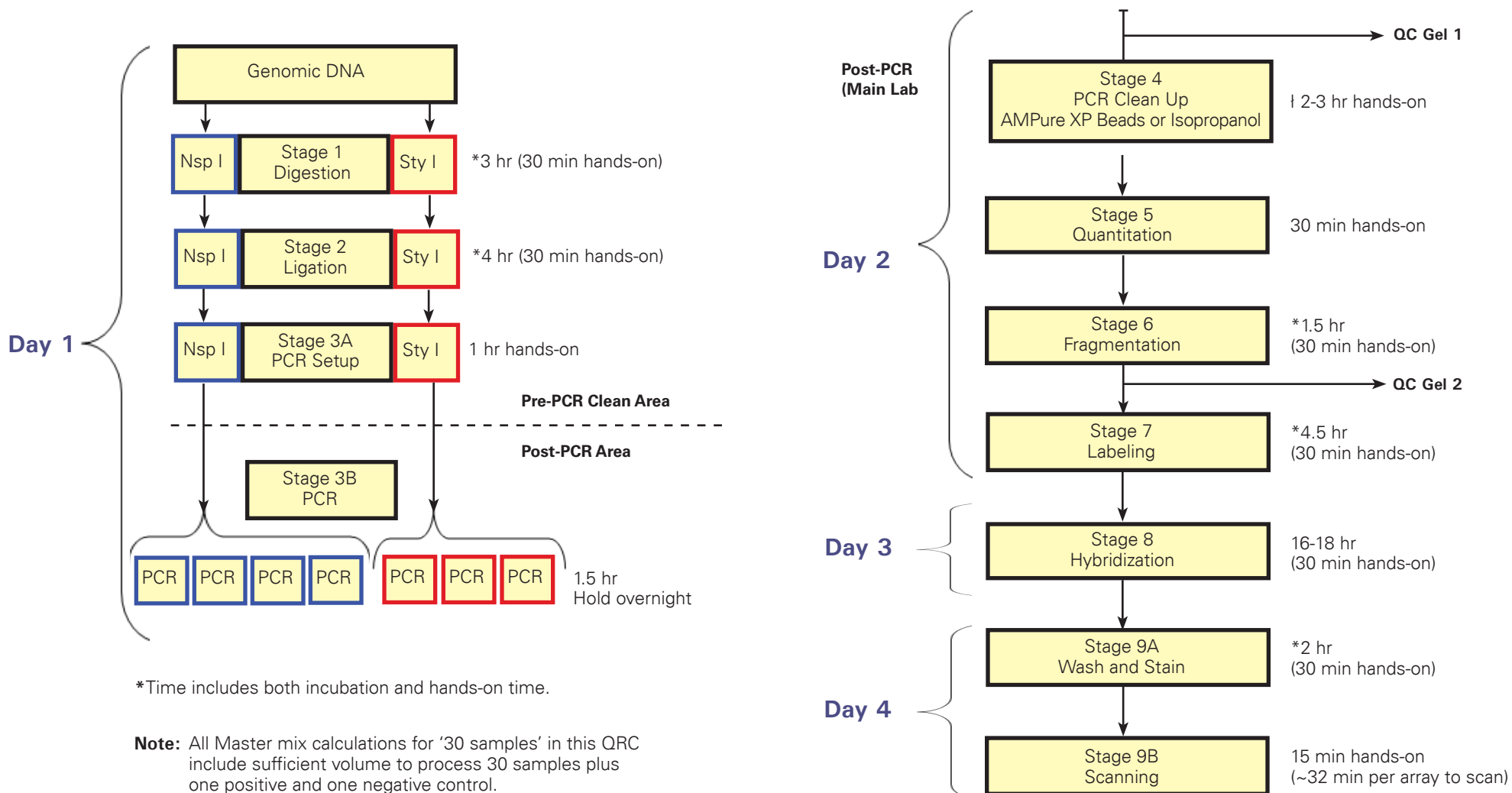


# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Workflow Overview

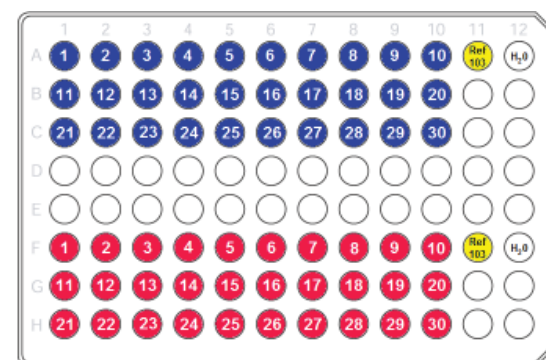


# Quick Reference Card

SNP 6.0 Cytogenetics Copy Number Assay  
Core Reagent 100 reaction Kit, 3x30 Sample  
Stage 1 – Digestion

DIGESTION MASTER MIX						
Reagent	Per Sample	30 Samples Nsp	✓	30 Samples Sty	✓	Lot Number
Water, molecular biology grade	11.55 µL	423.5 µL		423.5 µL		
Nsp Buffer	2.00 µL	73.3 µL		—		
Sty Buffer	2.00 µL	—		73.3 µL		
BSA (100X; 10 mg/mL)	0.20 µL	7.3 µL		7.3 µL		
Nsp I (10 U/ µL)	1.00 µL	36.7 µL		—		
Sty I (10 U/ µL)	1.00 µL	—		36.7 µL		
<b>Total Volume</b>	<b>14.75 µL</b>	<b>540.8 µL</b>	<b>—</b>	<b>540.8 µL</b>	<b>—</b>	<b>—</b>

## Suggested Digestion and Ligation Plate Map



Nsp Samples	Volume	Sty Samples	Volume
gDNA (50 ng/ µL)	5.00 µL (250 ng)	gDNA (50 ng/ µL)	5.00 µL (250 ng)
Nsp Master Mix	14.75 µL	Sty Master Mix	14.75 µL
<b>Total Volume</b>	<b>19.75 µL</b>	<b>Total Volume</b>	<b>19.75 µL</b>

Cyto Digest	
Temp	Time
37 °C	2 hr
65 °C	20 min
4 °C	Hold

1. Seal plate with adhesive film.
2. Vortex plate at high speed for 3 sec.
3. Spin down at 2000 rpm for 30 sec.
4. Ensure lid of thermal cycler is preheated.
5. Load plate onto thermal cycler and run the *Cyto Digest* program.

Proceed to Ligation

### Important Points

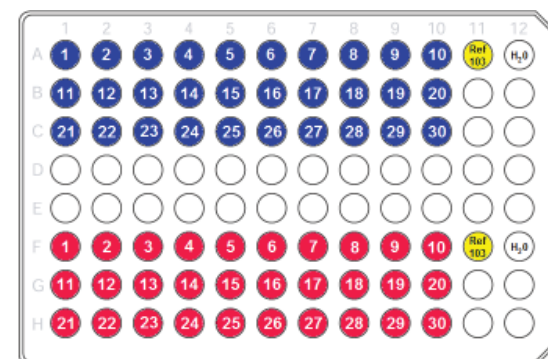
- Aliquot genomic DNA (gDNA) to opposite ends of the plate to lessen the chance of pipetting errors.
- Add gDNA to all sample wells.
- Two digestion master mixes are prepared (Nsp and Sty).
  - Be sure to use the correct enzyme for each master mix (Nsp or Sty)
  - Leave Nsp and Sty enzymes at -20 °C until ready to use.
- Add 5 µL Ref103 DNA as positive control to wells marked “Ref 103”.
- Add 5 µL water (molecular biology grade) as negative control to wells marked “H<sub>2</sub>O”.

# Quick Reference Card

SNP 6.0 Cytogenetics Copy Number Assay  
Core Reagent 100 reaction Kit, 3x30 Sample  
Stage 2 – Ligation

LIGATION MASTER MIX						
Reagent	Per Sample	30 Samples Nsp	✓	30 Samples Sty	✓	Lot Number
10XT4 DNA Ligase Buffer	2.50 µL	116.0 µL		116.0 µL		
Nsp Adaptor	0.75 µL	34.8 µL		—		
Sty Adaptor	0.75 µL	—		34.8 µL		
T4 DNA Ligase (400 U/µL)	2.00 µL	92.8 µL		92.8 µL		
<b>Total Volume</b>	<b>5.25 µL</b>	<b>243.6 µL</b>	—	<b>243.6 µL</b>	—	—

## Suggested Digestion and Ligation Plate Map



Nsp Samples	Volume	Sty Samples	Volume
Nsp Digested Sample	19.75 µL	Sty Digested Sample	19.75 µL
Nsp Master Mix	5.25 µL	Sty Master Mix	5.25 µL
<b>Total Volume</b>	<b>25.00 µL</b>	<b>Total Volume</b>	<b>25.00 µL</b>

Cyto Ligate	
Temp	Time
16 °C	3 hr
70 °C	20 min
4 °C	Hold

1. Seal plate with adhesive film.
2. Vortex plate at high speed for 3 sec.
3. Spin down at 2000 rpm for 30 sec.
4. Ensure lid of thermal cycler is preheated.
5. Load plate onto thermal cycler and run the *Cyto Ligate* program.

Proceed to PCR Setup

### Important Points

- Sample plate used for digestion and ligation.
- Two ligation master mixes are prepared (Nsp and Sty).
  - Be sure to use the correct adaptor for each master mix (Nsp or Sty).
  - Leave T4 DNA Ligase at -20 °C until ready to use.
  - Thaw T4 DNA Ligase Buffer on ice. Vortex to ensure any precipitate is resuspended and buffer is clear.

# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Stage 3a – PCR

1. Dilute ligated samples.

Nsp Samples	Volume Per Sample	Sty Samples	Volume Per Sample
Nsp Ligated Sample	25.0 µL	Sty Ligated Sample	25.0 µL
Water, molecular biology grade	75.0 µL	Water, molecular biology grade	75.0 µL
<b>Total Volume</b>	<b>100 µL</b>	<b>Total Volume</b>	<b>100 µL</b>

2. Seal plate with adhesive film.

3. Vortex at high speed for 3 sec; spin down at 2000 rpm for 30 sec.

4. Transfer four 10 µL aliquots of each Nsp sample to the PCR plate.

5. Transfer three 10 µL aliquots of each Sty sample to the PCR plate.



6. Prepare the PCR Master Mix.

7. Add PCR Master Mix to samples.

Sample	Volume
Nsp or Sty Sample	10 µL
PCR Master Mix	90 µL
<b>Total Volume</b>	<b>100 µL</b>

8. Seal PCR plate, vortex at high speed for 3 sec, spin down at 2000 rpm for 30 sec.

9. Keep plate on ice and move to Post-PCR Room/Area.

10. Ensure thermal cycler lid is preheated.

11. Load plate onto thermal cycler and run the *Cyto PCR* program.

12. Hold overnight.

PCR MASTER MIX				
Reagent	Per Sample	30 Samples	✓	Lot Number
Water, molecular biology grade	39.5 µL	9,730 µL		
TITANIUM™ Taq PCR Buffer (10X)	10.0 µL	2,468 µL		
GC-Melt/Beatine (5M)	20.0 µL	4,918 µL		
dNTPs (2.5 mM each)	14.0 µL	3,448 µL		
PCR Primer 002 (100 µM)	4.5 µL	1,103 µL		
TITANIUM™ Taq Polymerase	2.0 µL	490 µL		
<b>Total Volume</b>	<b>90.0 µL</b>	<b>22,157 µL</b>		

Pre-PCR Area

Post-PCR Area

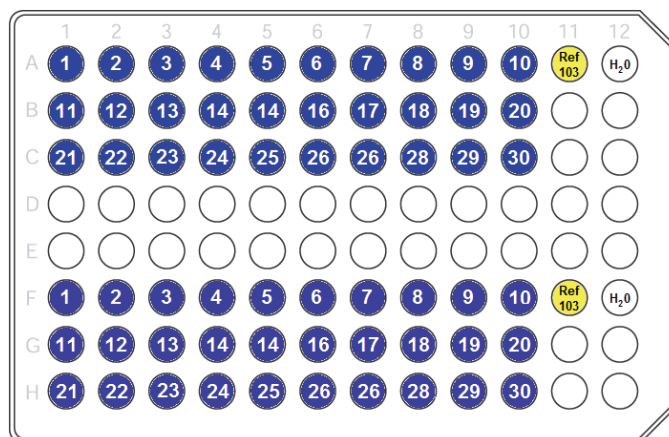
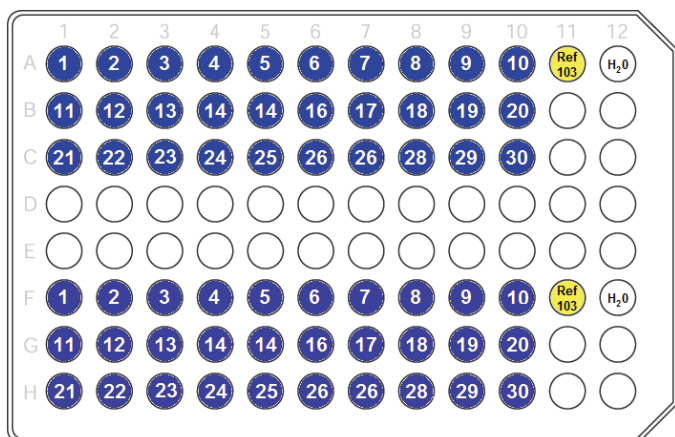
# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay

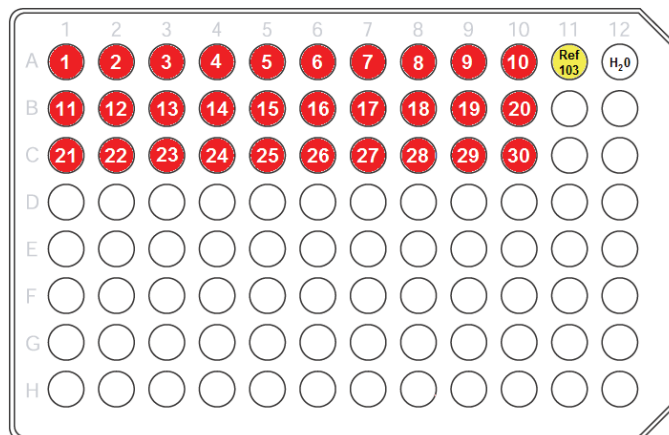
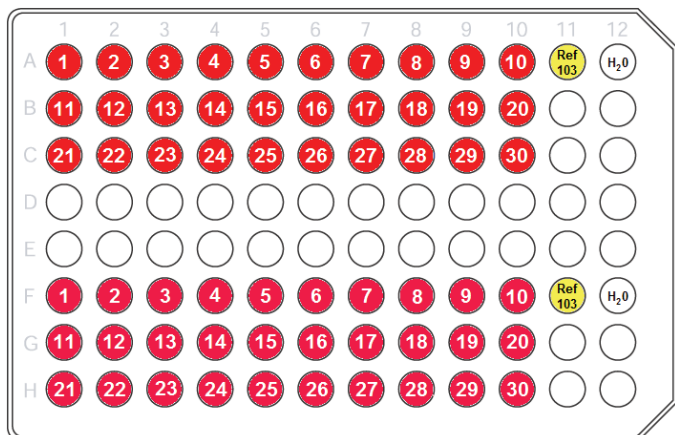
Core Reagent 100 reaction Kit, 3x30 Sample

Stage 3a – PCR Plate Maps and Thermocycler Programs

### Suggested Nsp PCR Plate Map



### Suggested Sty PCR Plate Map



### Thermocycling Programs

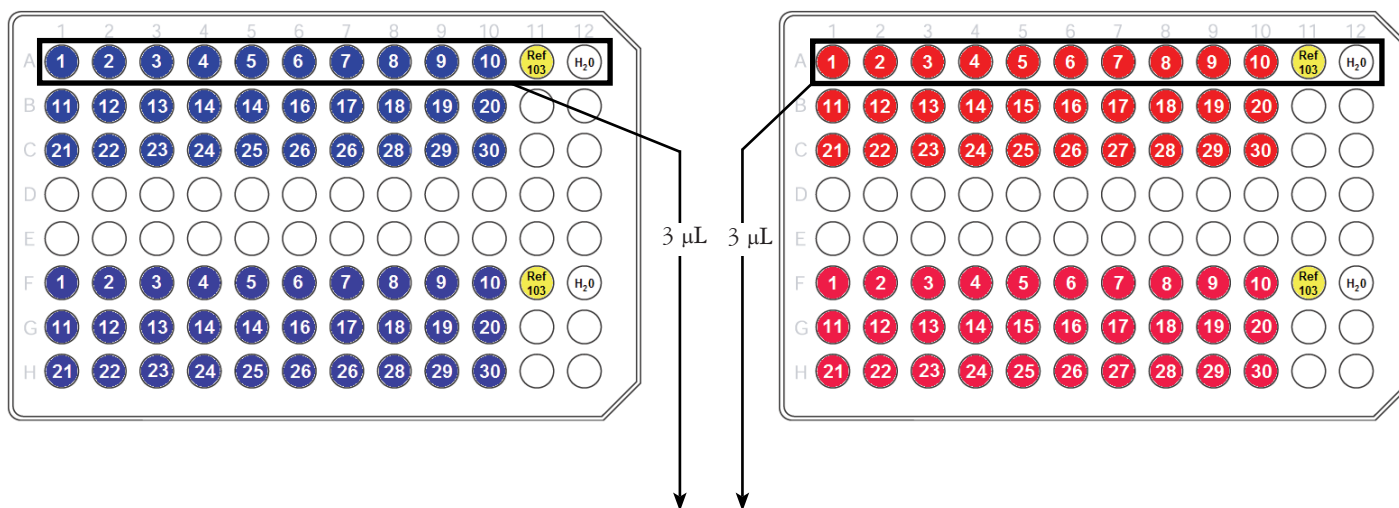
Cyto PCR – ABI 9700		
Temp	Time	Cycles
94 °C	3 min	—
94 °C	30 sec	30
60 °C	45 sec	
68 °C	15 sec	
68 °C	7 min	—
4 °C	Hold	—

Cyto PCR – MJ Tetrad PTC-225		
Temp	Time	Cycles
94 °C	3 min	—
94 °C	30 sec	30
60 °C	30 sec	
68 °C	15 sec	
68 °C	7 min	—
4 °C	Hold	—

# Quick Reference Card

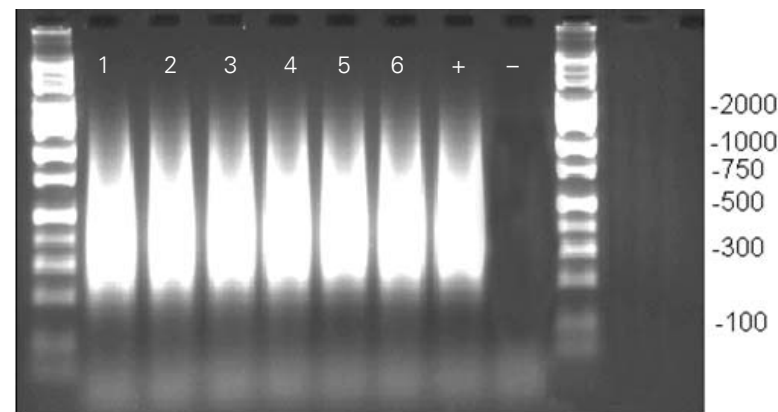
SNP 6.0 Cytogenetics Copy Number Assay  
Core Reagent 100 reaction Kit, 3x30 Sample  
Stage 3b – QC Gel #1, PCR Product Quality Check

## Suggested PCR Plate Maps



Transfer to corresponding wells of 96-well plate.  
These wells should already contain 3  $\mu$ L Gel Loading Dye.

1. Aliquot 3  $\mu$ L of 2X Gel Loading Dye to a new 96-well plate labeled "PCR gel". Mark rows A-D as Nsp and rows F-H as Sty.
2. Transfer 3  $\mu$ L of each representative Nsp PCR product to the PCR gel plate (rows A-D).
3. Transfer 3  $\mu$ L of each representative Sty PCR product to the PCR gel plate (rows F-H). Be sure to cover reactions across thermal cyclers. Be sure to include the positive and negative control.
4. Seal the gel plate.
5. Vortex on high speed for 3 sec; spin down at 2000 rpm for 30 sec.
6. Load the total volume of each well from the PCR gel plate onto a 2% TBE gel with ladder.
7. Run the gel at 120V for one hour.
8. Confirm that the product distribution is between ~200 to 1100bp.
9. Save the image.
10. Continue to stage 4 – PCR purification.



Example of PCR products run on a 2% TBE gel at 120 volts for 1 hour.  
Average product size is between 200 and 1100 bp.

# Quick Reference Card

SNP 6.0 Cytogenetics Copy Number Assay  
Core Reagent 100 reaction Kit, 3x30 Sample  
Stage 4 – Post-PCR Cleanup

**IMPORTANT:** There are two methods that can be used to clean up the PCR product. See the tables below for method comparison.

## Bead-Based Protocol

Stage 4A (page 8-10)

Stage 4A – DNA Clean Up Using AMPure XP Beads Duration ~ 180 min		
Steps in Process	Step Duration (min)	Hands on Time (min)
Sample/magnetic bead incubation	10	5
Initial vacuum step	40 - 60	5
First ethanol vacuum step (wash step)	10 - 20	5
Final ethanol vacuum step (dry step)	15	5
Elute DNA in Elution Buffer (EB)	30	< 2
Resuspend beads in Elution Buffer (EB)	30	5
Elution on vacuum manifold	5 - 15	< 2
<b>Total Time</b>	<b>140 - 180</b>	<b>&lt; 45</b>

## Isopropanol Method

Stage 4B (page 11)

Stage 4B – DNA Clean Up Using Isopropanol Duration ~ 130 min		
Step	Step Duration (min)	Hands on Time (min)
EDTA incubation	10 - 15	5
Isopropanol precipitation	30 - 40	< 10
Centrifugation	30	< 2
Pour off isopropanol	5	< 2
Ethanol wash	10	5
Pour off ethanol and dry	5	5
Resuspend pellet in Elution Buffer (EB)	30 - 50	5
<b>Total Time</b>	<b>90 - 155</b>	<b>&lt; 35</b>



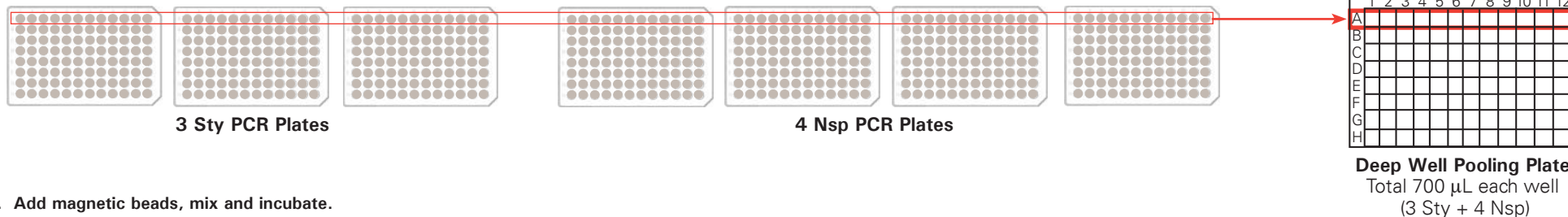
# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay

Core Reagent 100 reaction Kit, 3x30 Sample

Stage 4a – Post PCR Clean-up Using AMPure XP Beads

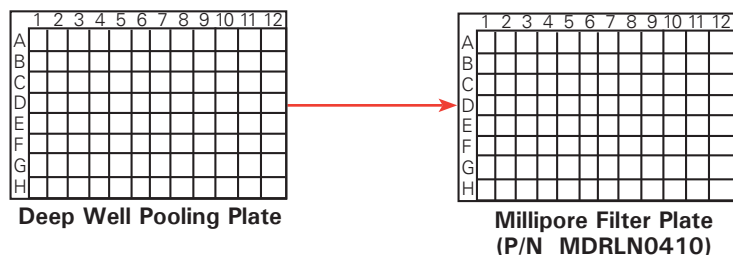
### 1. Pool Sty and Nsp PCR products into deep well plate using 12-channel P200 set to 110 $\mu$ L.



### 2. Add magnetic beads, mix and incubate.

- Shake bottle of magnetic beads vigorously until all beads are resuspended.
- Add 1.0 mL magnetic beads to each pooled sample.
- Mix well 5X by pipetting up and down using 12-channel P1200.
- Visually confirm that each well is mixed.
- Cover plate and incubate for 10 min at room temperature.

### 3. Place Millipore filter plate on vacuum manifold; then transfer incubated samples to filter plate (2 x 900 $\mu$ L). Seal the empty wells.



#### Note:

- Take care not to mix up samples while pooling. Work uninterrupted.
- Do not touch filters with pipet tips.
- Preheat hybridization oven for 30 min at 50  $^{\circ}$ C



## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Stage 4a – Post PCR Clean-up Using AMPure XP Beads

### 4. Apply vacuum until all of the liquid has been pulled through the filter (~40 to 60 min).

Visually check that all wells are dry, they should appear dull (matte).

#### Ensure all liquid is removed from the filter plate.

- With the vacuum on, tap the top of the plate twice with the palm of your hand.
- Turn the vacuum off and remove the plate.
- Firmly blot the plate on lint-free tissue until no wet spots are observed.
- Place the plate back on the manifold.
- Turn on the vacuum for 3 min.
- With the vacuum on, tap the top of the plate twice with the palm of your hand. (Repeat step 4B - 4D, 1x.)

### 5. Wash each sample with 1.8 ml 75% EtOH (room temperature) using a P1200 as follows:

- With the vacuum on, add 900  $\mu$ l to each sample.
- After ~2 min under vacuum, add another 900  $\mu$ l to each sample.
- Continue to apply vacuum until all of the liquid has been pulled through the filter (~10-15 min).  
*Visually check that all wells are dry. Maximum time 20 min.*

### 6. Ensure all EtOH is removed from the filter plate

- With the vacuum on, tap the top of the plate twice with the palm of your hand.
- Turn the vacuum off and remove the plate.
- Firmly blot the plate on lint-free tissue until no wet spots are observed.
- Return the plate to the manifold and apply vacuum for 2 - 5 min.
- Turn the vacuum off and remove the plate.
- Firmly blot the plate on lint-free tissue until no wet spots are observed.
- Let sit at room temperature for 10 min.

### 7. Elute DNA from beads at 50C for 30 min.

- Attach elution catch plate using lab tape.  
Seal on two sides only – do not seal completely.
- Add 60  $\mu$ l buffer EB to each sample. Buffer EB at 4 °C.  
*Tap plate stack to move all buffer EB onto the filters at the bottom of each well.*
- Loosely cover with plate lid or clear lid from pipet tip box.
- Place the covered plate stack in a 50 °C hyb oven for 30 min  
*If using an Affymetrix hybridization oven, place plate on the right front side (away from the air vent).*
- Remove the plate cover and seal all wells tightly with adhesive film.
- Place on Jitterbug for 30 min at setting 7.  
*Inspect each well to ensure all beads are resuspended.*
- Remove the seal and place plate stack INSIDE the vacuum manifold.  
*Ensure empty wells are sealed.*
- Apply vacuum until all liquid has been pulled through (~5 to 15 min).  
*Visually check that all wells are dry. If any wells are still wet or shiny, continue to apply vacuum for 15 min.*
- Turn off vacuum, seal entire plate and centrifuge the plate stack for 5 min at 1400 RCF.  
*Centrifuge at room temperature.*
- Remove the catch plate containing the eluate and seal.  
There should be ~ 50  $\mu$ l in each well.

#### NOTE:

- Do not extend vacuum steps beyond the maximum time listed.  
**UNDER DRYING CAN RESULT IN INCOMPLETE ELUTION.**
- Once samples are applied to the filter plate, keep at room temperature. If stored at 4 °C, yields will be reduced and array results compromised.

# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay

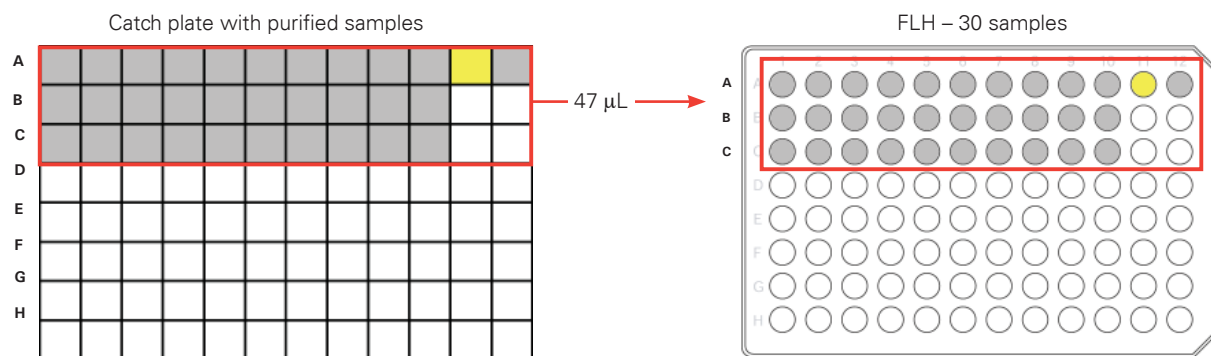
Core Reagent 100 reaction Kit, 3x30 Sample

Stage 4a – Post PCR Clean-up Using AMPure XP Beads

### 8. Prepare fragmentation/label/hyb plate (FLH).

- A. Mix eluted DNA before transfer for fragmentation and quantification on the Jitterbug for 5 min at setting 7.
- B. Spin briefly, then remove seal.
- C. Transfer 47  $\mu$ L to a new PCR plate using a 12-channel P200 (this is the fragmentation plate).

**NOTE:** To avoid mixing up your samples, re-label top 4 rows of plate FLH (Frag, Label, Hyb plate).

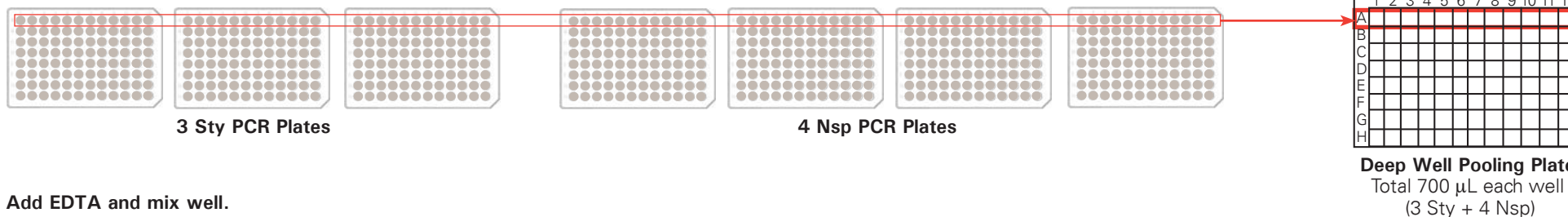


**OD is measured in Stage 5. Proceed to page 12 for instructions.**

# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Stage 4b – Post PCR Clean up Using Isopropanol

1. Pool 700  $\mu$ L of Sty and Nsp PCR products into deep well plate, row by row (see diagram).



2. Add EDTA and mix well.

- Pipette 1 mL of 0.5 M EDTA into a reservoir.
- From this reservoir, pipette 12  $\mu$ L of 0.5 M EDTA into each well using a multi-channel pipette.
- Mix the sample thoroughly by aspirating up and down 5 - 10 times using a P200 multi-channel pipette set at 200  $\mu$ L. Change tips between rows.

3. Cover the plate incubate for 10 min at room temperature.  
Inspect the plate and make sure the samples have turned clear.

4. Prepare Master Mix: 200  $\mu$ L of NH<sub>4</sub>OAC (7.5M) and 700  $\mu$ L of isopropanol for each sample.  
Transfer 900  $\mu$ L of the Master Mix to each pooled PCR product.  
Change tips and move to the next row until all the samples are done.

5. Cover the plate with a loose fitting lid and leave at room temperature for 30 min.  
**Note:** Turn on the centrifuge and cool to 4 °C.

6. Seal the plate with clear adhesive seal. Centrifuge the plate at 2,250 RCF for 30 minutes at 4 °C.  
**Note:** RCF is NOT the same as RPM. Use RCF (relative centrifugal force)

7. Remove the plate seal and **carefully** decant the supernatant by slowly inverting the plate upside down without disturbing the pellets. Discard supernatant in a waste container.

8. Place the plate upside down on lab tissue for 2 min on a bed of Kimwipes® tissues. Do not tap the plate!

9. Wash the pellets by adding 1.6 mL of 75% ethanol. The ethanol **MUST** be at room temperature.  
Pipet directly onto the pellets to help dislodge them from the bottom of the plate.  
Once the pellets are dislodged, leave the plate at room temperature on the bench top for 2 min.

10. Seal the plate with clear adhesive seal.  
Spin the samples again at 2,250 RCF for 5 min at 4 °C.

11. Remove the plate seal and carefully pour off the ethanol wash by slowly inverting the plate upside down without disturbing the pellets.

12. Place the plate upside down on Kimwipes® tissue for 2 min. Do NOT immediately tap the plate on the Kimwipes tissues! Change the tissues if too wet. After the 2 min time is up, tap the plate gently on the Kimwipes until no wet spots are observed on the tissues. Then flip the plate upright and tap on the bench to bring the pellets to the bottom of each well. Let dry another 2 min on the bench.

13. Add 55  $\mu$ L EB buffer to each sample. Check that pellets are immersed.

14. Seal the plate. Shake gently for 30 min (Jitterbug setting 5).  
Vortex until pellets are dissolved.

15. Using a multi-channel pipette, transfer 47  $\mu$ L to a new PCR plate.

**Proceed to Stage 5 - Quantitation on page 12.**

# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Stage 5 – QC #2, Quantitation

### Plate Spectrophotometer

1. Dilute purified PCR 1:100.
  - Pipette 2  $\mu$ L from the Fragmentation plate into 198  $\mu$ L water.
  - Thoroughly mix the samples and water using one of these methods:
    - Seal the plate, vortex, and spin down OR pipet up and down 5 times.
2. Measure the OD of each PCR product at 260, 280 and 320 nm.
3. Determine the OD<sub>260</sub> measurement for the water blank and average.
4. Calculate one OD reading for every sample:  $OD = (\text{sample OD}) - (\text{average water blank OD})$
5. Calculate the undiluted concentration for each sample in  $\mu\text{g}/\mu\text{L}$ :  $OD \times 0.05 \mu\text{g}/\mu\text{L} \times 100$ .

### NanoDrop

1. Dilute purified PCR 1:10.
  - Pipette 2  $\mu$ L from the Fragmentation plate into 18  $\mu$ L water.
  - Thoroughly mix the samples and water using one of these methods:
    - Seal the plate, vortex, and spin down OR pipet up and down 5 times.
2. Blank the NanoDrop using water.
3. Take 2  $\mu$ L of 1:10 diluted sample and measure the OD of each PCR product at 260, 280 and 320 nm.
4. Calculate the undiluted concentration of each sample in  $\mu\text{g}/\mu\text{L}$ :  $OD \text{ reading} \times 0.05 \mu\text{g}/\mu\text{L} \times 10$ .

### Assess OD Readings

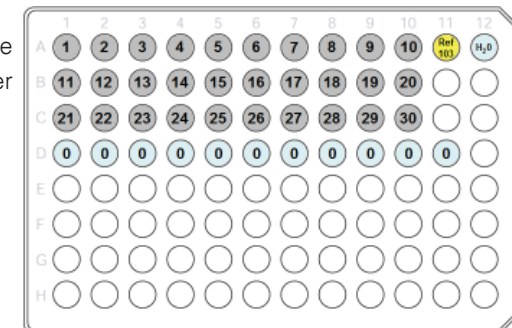
- An acceptable OD should fall within the range of 0.9 to 1.4.
- DNA yield equivalent = 4.5 to 7.0  $\mu\text{g}/\mu\text{L}$ .
- The OD<sub>260</sub>/OD<sub>280</sub> ratio should be between 1.8 and 2.0.
- The OD<sub>320</sub> measurement should be very close to zero ( $< 0.1$ ).
- If metrics fall outside of these ranges, refer to the *Affymetrix® Cytogenetics Copy Number Assay User Guide* for more information.

**NOTE:** Optimal PCR yield for downstream steps is above 5.5  $\mu\text{g}/\mu\text{L}$ .

### Example showing one row of samples

#### Plate Spectrophotometer

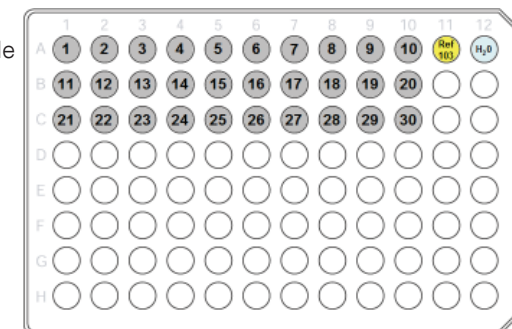
**Row A:** 198  $\mu$ L water + 2  $\mu$ L sample  
**Row B:** 200  $\mu$ L water



	1	2	3	4	5	6	7	8	9	10	11	12
A	1	2	3	4	5	6	7	8	9	10	100	H <sub>2</sub> O
B	11	12	13	14	15	16	17	18	19	20		
C	21	22	23	24	25	26	27	28	29	30		
D	0	0	0	0	0	0	0	0	0	0		
E												
F												
G												
H												

#### NanoDrop

**Row A:** 18  $\mu$ L water + 2  $\mu$ L sample



	1	2	3	4	5	6	7	8	9	10	11	12
A	1	2	3	4	5	6	7	8	9	10	100	H <sub>2</sub> O
B	11	12	13	14	15	16	17	18	19	20		
C	21	22	23	24	25	26	27	28	29	30		
D												
E												
F												
G												
H												

# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Stage 6 – Fragmentation

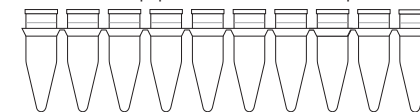
FRAGMENTATION MASTER MIX					
Reagent	2 U/ $\mu$ L	2.5 U/ $\mu$ L	3.0 U/ $\mu$ L	✓	Lot Number
Water, molecular biology grade	170 $\mu$ L	215 $\mu$ L	260 $\mu$ L		
10X Fragmentation Buffer	20 $\mu$ L	25 $\mu$ L	30 $\mu$ L		
Fragmentation Reagent	10 $\mu$ L	10 $\mu$ L	10 $\mu$ L		
<b>Total Volume</b>	<b>200 <math>\mu</math>L</b>	<b>250 <math>\mu</math>L</b>	<b>300 <math>\mu</math>L</b>		

**NOTE:** Please refer to manual if the U/ $\mu$ L of this reagent is different from the table above or if processing less than 30 samples.

1. Add 5  $\mu$ L of 10X Fragmentation Buffer to each sample. ....

Purified Sample	45 $\mu$ L
10X Fragmentation Buffer	5 $\mu$ L
<b>Total Volume</b>	<b>50 <math>\mu</math>L</b>

Aliquot Fragmentation Master Mix equally to strip tubes. Use a multi-channel pipet to add to samples.



- Prepare the Fragmentation Master Mix from table above.
- Aliquot the master mix equally to one set of strip tubes.
- Using a multi-channel pipet, add 5  $\mu$ L of Fragmentation Master Mix to each sample.
- Seal the plate with adhesive film.
- Vortex at high speed for 3 sec.
- Spin down at 2000 rpm for 30 sec.
- Start the Cyto Frag program and pause when it reaches temperature.
- Load plate onto thermal cycler and un-pause to run program.

Diluted sample	50 $\mu$ L
Fragmentation Master Mix	5 $\mu$ L
<b>Total Volume</b>	<b>55 <math>\mu</math>L</b>

Cyto Fragment	
Temp	Time
37 °C	35 min
95 °C	15 min
4 °C	Hold

Proceed immediately to Labeling.

### Important Points – Fragmentation Master Mix Preparation

- Check concentration of Fragmentation Reagent (enzyme; varies between 2 and 3 U/ $\mu$ L).
- Leave Fragmentation Reagent (enzyme) at  $-20$  °C until ready to use.
- Thaw 10X Fragmentation Buffer on ice.
- Keep all reagents, including water, on ice. Perform all additions on ice.
- Preheat thermal cycler block to 37 °C.

# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay

Core Reagent 100 reaction Kit, 3x30 Sample

Stage 7 – Labeling and QC #3, Fragmentation QC Gel

### Labeling

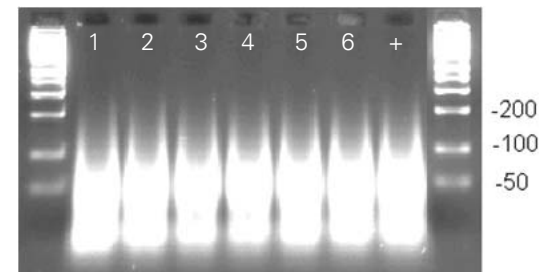
LABELING MASTER MIX				
Reagent	Per Sample 2 U/ $\mu$ L	30 Samples	✓	Lot Number
5xTdT Buffer	14.0 $\mu$ L	537 $\mu$ L		
30 mM DNA Labeling Reagent	2.0 $\mu$ L	77 $\mu$ L		
30U/ $\mu$ L TdT Enzyme (30 U/ $\mu$ L)	3.5 $\mu$ L	134 $\mu$ L		
<b>Total Volume</b>	<b>19.5 <math>\mu</math>L</b>	<b>748 <math>\mu</math>L</b>		

1. Transfer 2  $\mu$ L of each fragmented sample to the corresponding well of a fresh 96-well plate (the Fragmentation QC Gel Plate).
2. Prepare the Labeling Master Mix.
3. Add 19.5  $\mu$ L of master mix to each sample.
4. Tightly seal the plate, and vortex at high speed for 3 sec.
5. Spin down at 2000 rpm for 30 sec.
6. Load plate onto thermal cycler and run the *Cyto Label* program.

Cyto Label	
Temp	Time
37 °C	4 hr
95 °C	15 min
4 °C	Hold

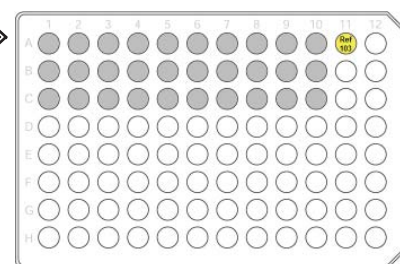
If possible, store the Label plate overnight at  $-20$  °C. Otherwise, OK to hold at 4 °C overnight.

### Fragmentation QC Gel



Example of fragmented samples run on a 4% TBE gel at 120 volts for 1 hr. Average fragment size is < 180 bp.

2  $\mu$ L



Fragmentation QC Gel Plate

1. While the *Cyto Label* program is running, finish preparing the gel plate by adding 4  $\mu$ L of Gel Loading Dye to each sample.
2. Seal the plate, vortex, and spin down.
3. Onto a 4% TBE gel, load 10  $\mu$ L of BioNexus All Purpose Hi-Lo Ladder to the first and last lanes.
4. Load the samples and run the gel.
5. Inspect the gel and compare against the figure shown here.

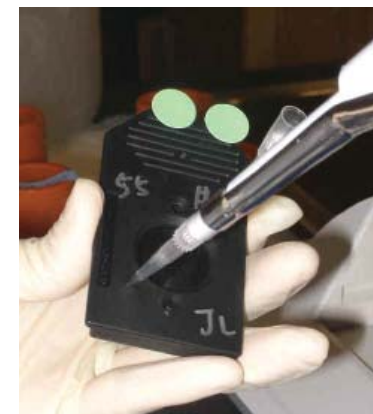
#### Important Points

- Leave the TdT enzyme at  $-20$  °C until ready to use.
- Thaw the 5X TdT Buffer and DNA Labeling Reagent on ice.
- Ensure the plate is tightly sealed to avoid evaporation while on the thermal cycler.

# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Stage 8 – Hybridization

HYBRIDIZATION MASTER MIX				
Reagent	Per Sample	30 Samples	✓	Lot Number
Hyb Buffer Part 1	165 µL	6,325 µL		
Hyb Buffer Part 2	15 µL	575 µL		
Hyb Buffer Part 3	7 µL	268.3 µL		
Hyb Buffer Part 4	1 µL	38.3 µL		
OCR	2 µL	76.7		
<b>Total Volume</b>	<b>190 µL</b>	<b>7,283.3 µL</b>		



Use a P200 pipet to load arrays.

1. Unpackage the arrays and allow to equilibrate to room temperature prior to use.
2. Preheat the hybridization ovens for at least 1 hr at 50 °C with the rotation turned on.
3. Prepare the Hybridization Master Mix.
4. Add 190 µL of master mix to each sample.
5. Tightly seal the plate, vortex at high speed for 30 sec, and spin down at 2000 rpm for 30 sec.
6. Load plate onto thermal cycler and run the *Cyto Hyb* program.

7. Leaving the samples on the thermal cycler, load 200 µL of sample onto each array using a single-channel P200 pipet.
8. Clean any excess fluid from around the septa.
9. Apply Tough-Spots to the septa and press firmly.
10. Load arrays into the hybridization oven four at a time.

Hyb arrays  
16 to 18 hr at 50 °C

Cyto Hyb	
Temp	Time
95 °C	10 min
49 °C	Hold

### Important Points

- Samples must remain on the thermal cycler while loading the arrays.
- To avoid damaging the septa, use a single-channel P200 pipet to load the arrays.
- Shake arrays a few times to ensure bubbles are not visible through the window.
- When 4 arrays are loaded, immediately place them into the hybridization oven.



# Quick Reference Card

## SNP 6.0 Cytogenetics Copy Number Assay Core Reagent 100 reaction Kit, 3x30 Sample Stage 9 – Washing and Staining

All buffers are prepared for you. You will need to aliquot the following from the bottles:



**Stain Cocktail 1**  
600  $\mu$ L/sample



**Stain Cocktail 2**  
600  $\mu$ L/sample



**Array Holding Buffer**  
1000  $\mu$ L/sample

### Washing and Staining Arrays

1. Remove the hybridization solution from each array.
2. Fill the arrays with 270  $\mu$ L **1X Array Holding Buffer**.
3. Load arrays onto the Fluidics Station.
4. Using AGCC, run the *GenomeWideSNP6\_450* protocol.

### Before Scanning

1. Ensure no bubbles are visible through the window.
2. Cover the septa with Tough-Spots; then load onto the scanner.

### Important Points

- The hybridization solution removed from the arrays can be stored long term at  $-80^{\circ}\text{C}$ .
- Stain Cocktail 1 and the Array Holding Buffer are light sensitive and must be stored at  $4^{\circ}\text{C}$ .
- If necessary, the array can be stored in Array Holding Buffer at  $4^{\circ}\text{C}$  for up to 3 hr before washing and staining.