



# Genotyping Polymorphic MicroRNA Target Sites in Individuals with Schizophrenia

Michael Moreau  
Marco Azaro  
Anthony Marcketta



# Overview of Presentation

- Schizophrenia background
- Experimental design:
  - Samples, selecting candidate genes
  - Identifying polymorphic microRNA targets sites
  - An improved bead-based SNP genotyping method
- Results of family-based genetic association analysis



# Schizophrenia

- Schizophrenia is a type of psychosis characterized by disturbed thought, perception, emotion, and behavior
  - Prevalence is about 0.5-1% in the general population
- The Psychotic Symptoms: Active manifestation of abnormal behavior, distortions of normal behavior
  - Delusions, hallucinations
- The Negative Symptoms: Absence or lack of normal behavior
  - Apathy, alogia, anhedonia, affective flattening
- The Cognitive Symptoms: Severe disruptions in speech, attention and concentration, psychomotor function, learning and memory

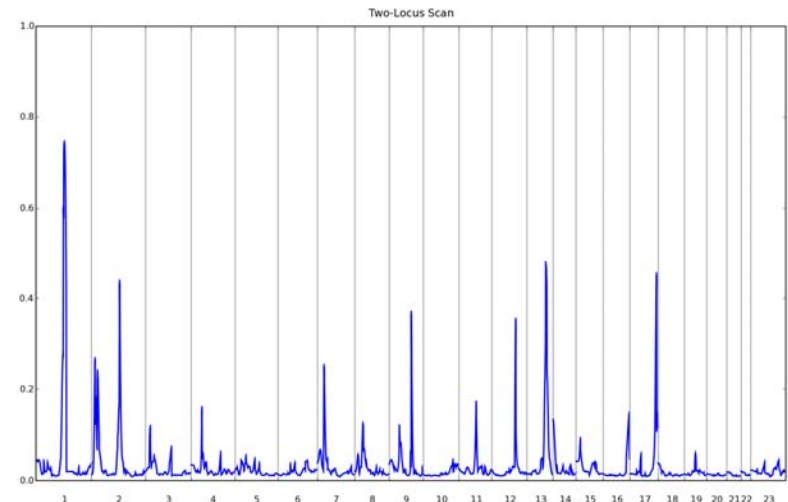
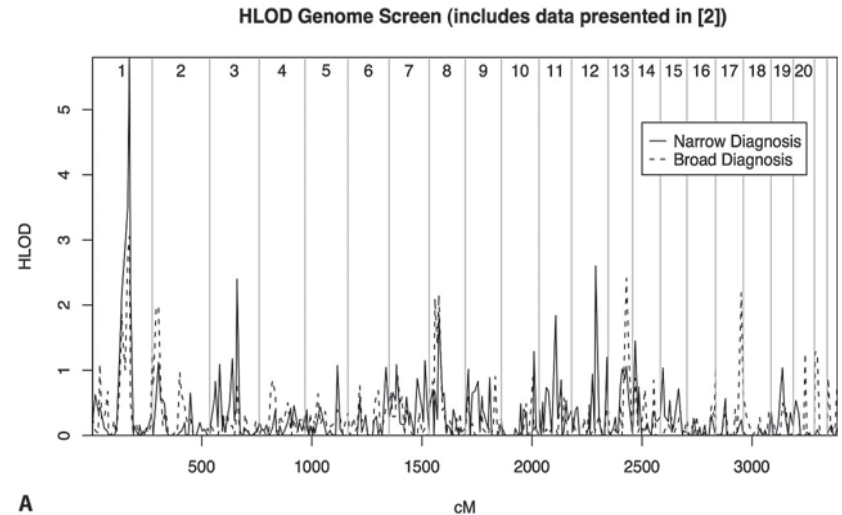


# Study Design - Samples

- Family-based genetic association study
  - 392 subjects from 25 extended pedigrees of Canadian Celtic ancestry
  - On average, 13.8 individuals per family and 3.6 affected with schizophrenia or schizoaffective disorder
  - Illness appeared to segregate in a unilineal, autosomal dominant manner

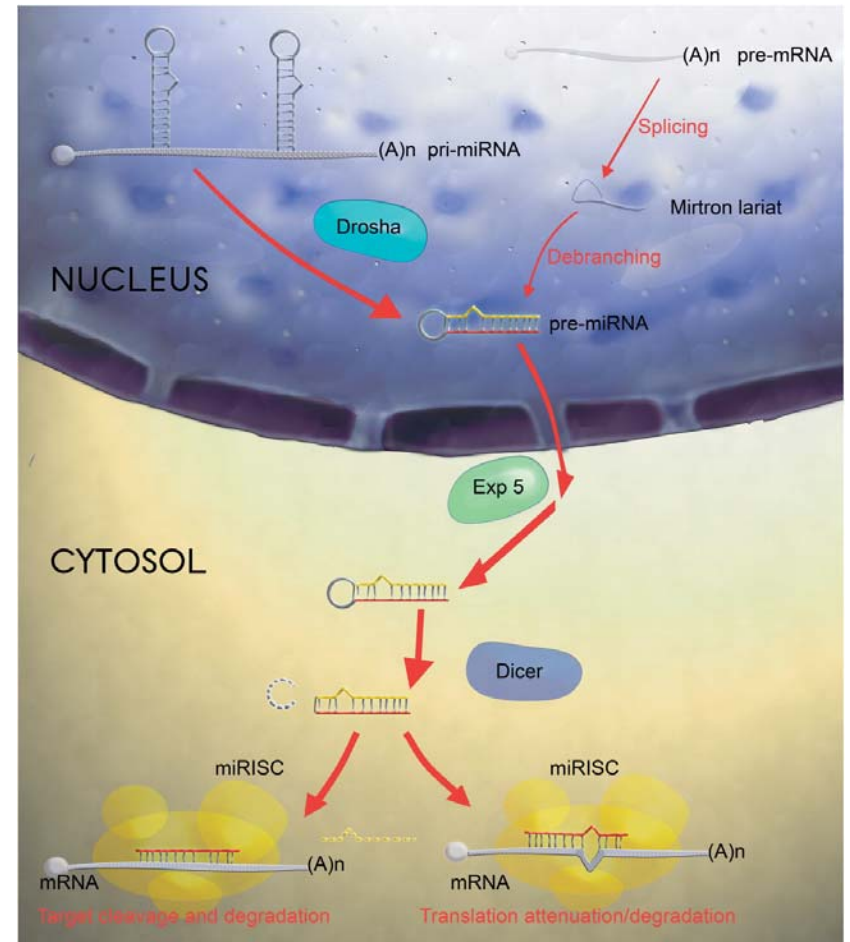
# Study Design – Candidate Gene Selection

- In 2000, a genome-wide linkage scan revealed strong linkage to 1q21-q22
- Reanalyzed using the PPL in 2006
- Fine mapping and mutation analysis identified putative functional variant in NOS1AP
- NOS1AP risk allele used to define liability classes, conditioned linkage scan performed to find factors that interact epistatically
- Peaks from conditional scan used to define genomic boundaries, mined for polymiRTs



# microRNAs

- microRNAs are ~21 nt non-coding RNA molecules that regulate gene expression post-transcriptionally
- miRNAs biosynthesis consists of two enzymatic processing steps
- Bind to 3' end of mRNA target
- Sequence complementarity affects efficacy of gene repression and mechanism of action



# Variation in miRNA genes

- Mutations in miRNA genes and targets have very different implications
  - Mutations in miRNA genes have broad implications
  - One miRNA can target hundreds of transcripts
  - Evolutionary selection against disrupting entire regulatory networks

Redundancy may ameliorate impact:

## A – Redundant genes encode miR-30c

Homo sapiens miR-30c-1 stem-loop (MI0000736)

```

a      cu      ugu u      u      aca      ---g a
ccaug  guag  g guaaaca ccu  cucucagcu  ug g
|||||  ||||  | ||||| |||  ||||| |||  ||
gguac  cguc  c cauuugu ggg  gagggucgg  ac c
a      --      uuc u      u      --a      ugga u
  
```

Homo sapiens miR-30c-2 stem-loop (MI0000254)

```

      uacu      u      aca      guggaa
aga      guaaaca ccu  cucucagcu  a
|||      ||||| |||  ||||| |||  g
ucu      cauuugu gga  gagggucga  g
uucu      c      --a      aagaau
  
```

## B – Multiple paralogues of miR-30

← Seed →

```

>hsa-miR-30a  UGUAAACAUCCUCGACUGGAAG
>hsa-miR-30b  UGUAAACAUCCUACACUCAGCU
>hsa-miR-30d  UGUAAACAUCCCCGACUGGAAG
>hsa-miR-30e  UGUAAACAUCCUUGACUGGAAG
>hsa-miR-30c  UGUAAACAUCCUACACUCUCAGC
  
```

↓  
A\*

\*C>A SNP in mir-30c from Iwai & Naraba, (2005)

# Variation in miRNA target sites

- Variants in miRNA target sites are more likely
  - Biological effects can be subtle or pronounced
  - SNPs may create, strengthen, weaken, or destroy miRNA target sites



Clop, A., F. Marcq, et al. (2006). "A mutation creating a potential illegitimate microRNA target site in the myostatin gene affects muscularity in sheep." *Nat Genet* **38**(7): 813-8.

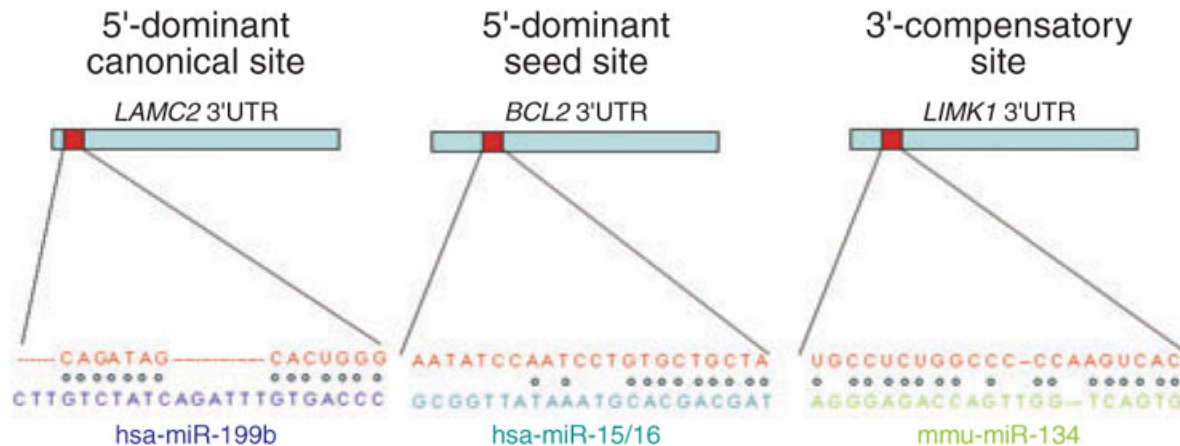


## Prediction of miRNA targets that overlap validated SNPs

- Given a list of genes, what do we need to identify polymorphic target sites?
  - in this particular case, we are interested in a set of 6467 genes that fall beneath linkage peaks
- We need:
  - 1) a source of accurately annotated mRNAs (AceView)
  - 2) a database of SNPs with accurate positional and validation status
  - 3) a high-throughput method to predict miRNA targets that does not rely on interspecies conservation
    - PROBLEM: this requirement invalidates the majority of existing target prediction programs!

# Why is miRNA target prediction so challenging?

miRNA-target heteroduplexes fall into 3 general classes:



- although perfect complementarity in the seed region (positions 2-7) is characteristic of a large class of alignments, the number and position of remaining matches, mismatches and bulges is highly variable
- standard signature-based alignment algorithms prioritize perfect seed matching
  - Many chance alignments, improved signal/noise by factoring in context determinants, evolutionary conservation
- may be better to look for conserved patterns that are not necessarily contiguous



# Identifying Polymorphic miRNA Target Sites

- *rna22* is a pattern-based approach
  - Training set from (Rfam) rel 7.1
  - Teiresias algorithm used to discover variable-length motifs
    - e.g. [AT][CG].TTTTT[CG]G..[AT]
  - Map reverse complements and define “target islands”
  - Pair each target with a candidate miRNA
- Experimentally demonstrated repressions of 30% or more for 168 of 226 tested targets



## Our procedure is a mutant offspring of Rna22

- with a few differences:
  - 1) We have newer and greatly expanded sets of mature miRNAs
    - Release 7.1 (10/06) had 319 human miRNAs in a total of 3102
    - Release 12.0 (9/08) has 692 human miRNAs in a total of 9169
      - we are considering patterns derived from human miRNAs only!
  - 2) We consider all alternative splice variants curated by AceView
    - NOTE: most target prediction programs ignore tissue-specific splice variants and alt. poly-A termini
  - 3) Our background probability computation uses a random sample of human genomic fragments that reflect size distribution of miRNA set
  - **4) We substitute 3'UTRs with all relevant valid SNPs**

# The Use of A Novel Genotyping Method

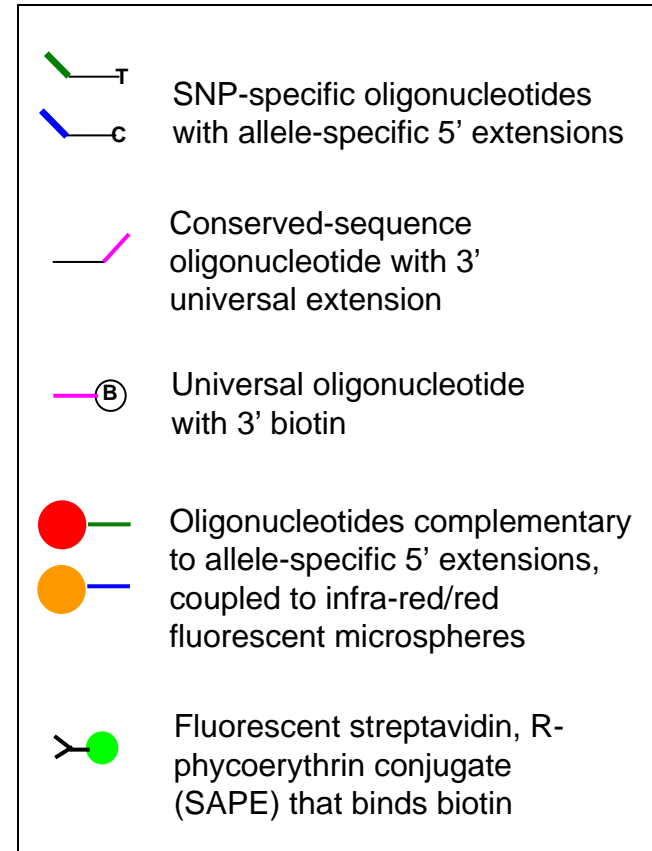
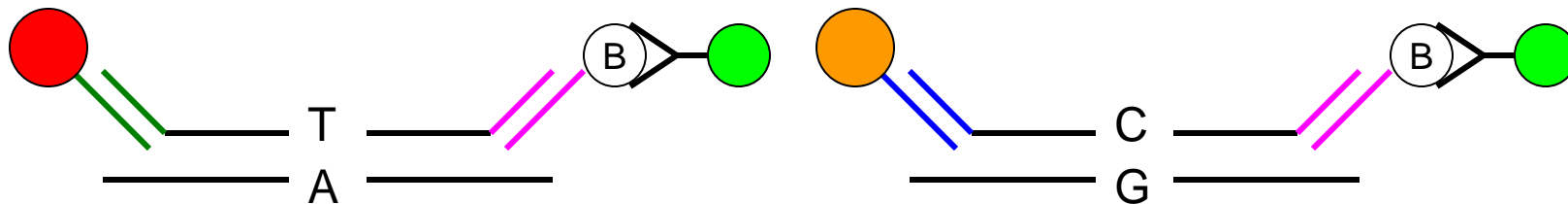
## PCR Product



**Step #1** – Single tube oligonucleotide ligation assay (OLA) to interrogate SNP



**Step #2** – Denature and hybridize to microspheres and universal oligos; incubate with SAPE



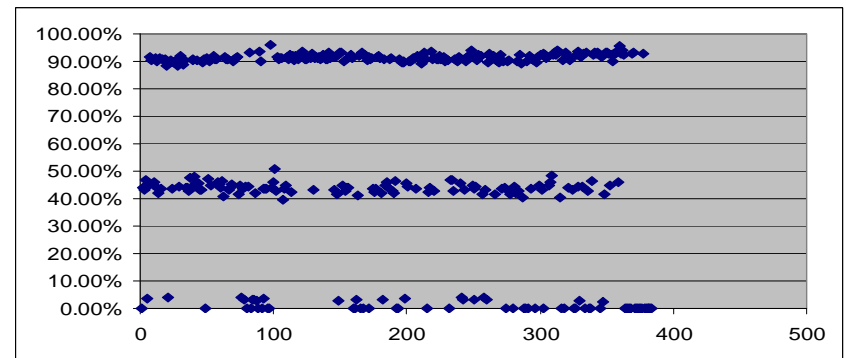
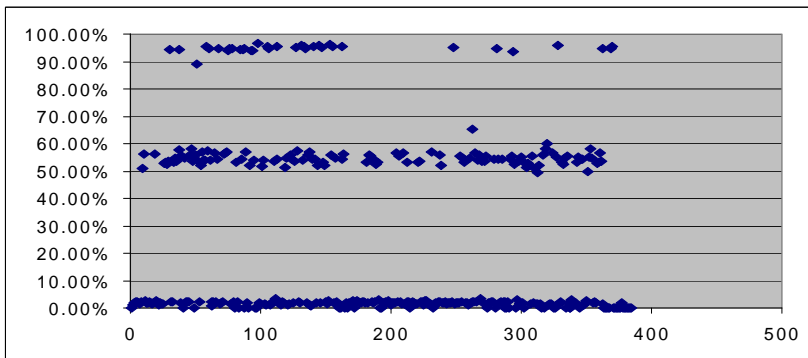
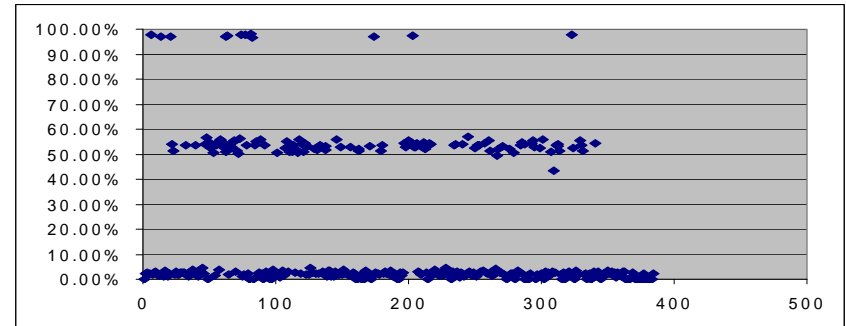
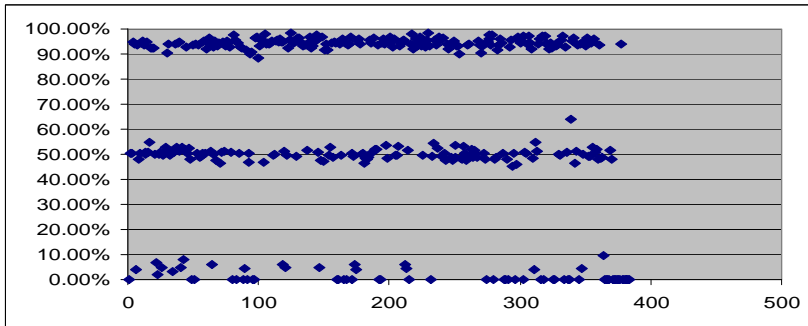
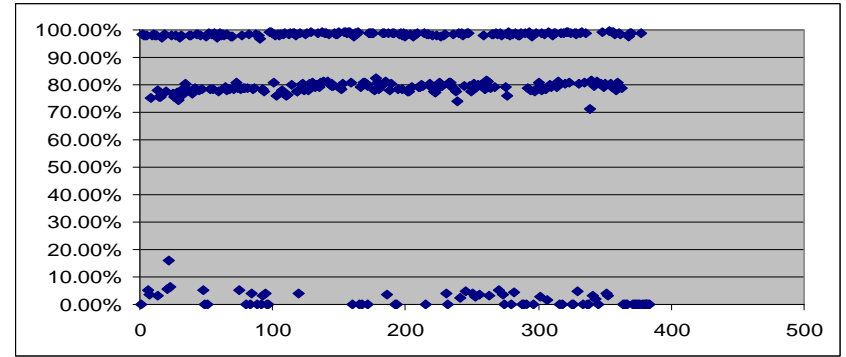
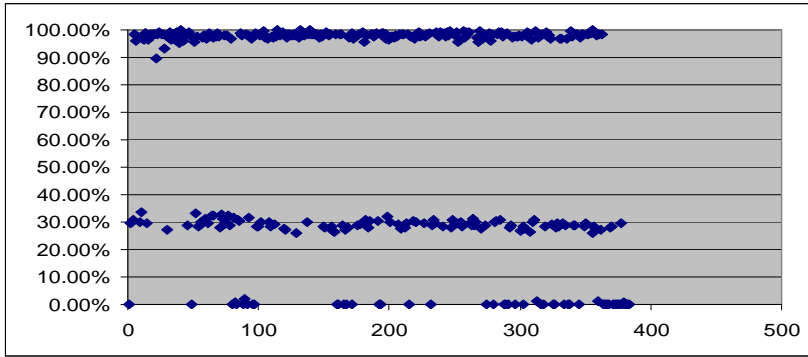
# The Use of A Novel Genotyping Method

- **Step #3** Detect using two color flow cytometric analysis (Luminex 100)



|   | rs75797_C | rs75797_G | rs79974_C | rs79974_T | rs80104_C | rs80104_G | rs98566_C | rs98566_T | rs98932_C | rs98932_G |
|---|-----------|-----------|-----------|-----------|-----------|-----------|-----------|-----------|-----------|-----------|
| 5 | 417       | 6202      | 188       | 91.5      | 69        | 44        | 52        | 34        | 53        | 82.5      |
| 3 | 119       | 4758      | 1707      | 1670      | 4668      | 6168.5    | 8399.5    | 10675.5   | 4883      | 504       |
| 5 | 207       | 4903      | 1878      | 1845      | 4548.5    | 6147      | 8849      | 11570     | 4965      | 333       |
| 5 | 233.5     | 4871.5    | 3370      | 192.5     | 4271      | 6266      | 9006      | 10236     | 4487.5    | 530       |
| 5 | 81        | 5132      | 3552      | 226       | 6777.5    | 240       | 599.5     | 15745     | 4371.5    | 347.5     |
| 3 | 304       | 4495      | 143       | 3463      | 7870.5    | 325.5     | 8815      | 10833     | 4218.5    | 405       |
| 5 | 230.5     | 5223      | 3675      | 257       | 5035      | 6047      | 13317     | 1217.5    | 5319      | 352.5     |
| 3 | 187       | 4734.5    | 1911      | 2084      | 4381      | 5751      | 12484.5   | 1352      | 4733      | 376       |
| 3 | 248       | 5213.5    | 1803      | 1782      | 4778      | 6661      | 9092.5    | 11030     | 5367      | 297       |
| 5 | 218.5     | 5115      | 3075      | 190       | 4486      | 6518      | 9744.5    | 11392     | 5074      | 454.5     |
| 0 | 170       | 4486      | 3367      | 172       | 7660      | 342       | 13771     | 1351      | 4980      | 476       |
| 5 | 248       | 5346      | 3525      | 247.5     | 8015      | 304.5     | 12279.5   | 1337      | 5556.5    | 532       |
| 5 | 365       | 4641      | 1841.5    | 1790      | 7793      | 151       | 8356.5    | 11476     | 5217.5    | 407       |
| 3 | 300       | 5306      | 1873      | 1871      | 4166.5    | 5967      | 8544      | 11016     | 4435      | 325.5     |
| 3 | 264       | 4867      | 3560      | 195.5     | 329.5     | 9530      | 12861     | 1248      | 4291      | 344       |
| 7 | 118       | 5231      | 2204      | 2126.5    | 4627      | 6374      | 9197      | 11821     | 4806      | 430       |
| 3 | 167       | 4936.5    | 2398      | 1985.5    | 8078      | 271.5     | 14065     | 1428      | 4998      | 393.5     |

# Genotype Calling





## Improvements to bead-based oligonucleotide ligation SNP genotyping assays

- Refinement and validation of a robust method for multiplexed SNP genotyping on the Luminex platform:
  - Oligonucleotide ligation assay (OLA)
  - Use of fewer beads
  - Incorporation of biotin using a universal oligo
  - Multiplexed PCR
  - No wash steps



# Improvements to bead-based oligonucleotide ligation SNP genotyping assays

**Table 1. Inter-well CV and MFI Range Using Low and High Number of Beads in Control Experiment**

| <b>200 input/20 count</b>   | <b>Positive control</b> | <b>Negative control</b> |
|---|-------------------------|-------------------------|
| Avg. inter-well CV  | 2%                      | 26%                     |
| Avg. MFI  | 21,968                  | 192                     |
| MFI range   | 21,003–22,842           | 62–299                  |
| 5000 input/300 count  | Positive control        | Negative control        |
| Avg. inter-well CV  | 2%                      | 20%                     |
| Avg. MFI  | 22,619                  | 162                     |
| MFI range   | 21,496–23,022           | 115–205                 |
| Median fluorescence intensity (MFI) was assessed using the high PMT setting of the Luminex 100 flow cytometer. Inter-well CVs were calculated manually and are based on the median. |                         |                         |



# Concordance between Luminex OLA and other genotyping methods

- Comparison of 8 SNPs in 14 human genomic DNA samples using direct sequencing; 100% concordance
- Comparison of 6 SNPs in 84 samples using a gel-based version of the OLA; 100% concordance
- Comparison using pyrosequencing:

| SNP               | concordant calls | % concordant  |
|-------------------|------------------|---------------|
| <b>ID345549</b>   | 231/232          | 99.60%        |
| <b>rs1123005</b>  | 287/289          | 99.30%        |
| <b>rs11806859</b> | 168/170          | 98.80%        |
| <b>rs12122048</b> | 58/58            | 100.00%       |
| <b>rs4657179</b>  | 290/290          | 100.00%       |
| <b>rs4657187</b>  | 167/168          | 99.40%        |
| <b>rs905720</b>   | 263/264          | 99.62%        |
| Total             | <b>1464/1471</b> | <b>99.52%</b> |



# Cost breakdown

|                       | Cost/rxn     | Cost/<br>genotype |
|-----------------------|--------------|-------------------|
| PCR reagents          | 0.156        |                   |
| OLA reagents          | 0.085        |                   |
| Luminex<br>reagents   | 0.053        |                   |
| <b>Fixed Cost</b>     | <b>0.294</b> |                   |
| Primers               | 0.30         | 0.02              |
| Beads<br>(200/allele) | 0.6          | 0.04              |
| <b>Total</b>          | <b>1.194</b> | <b>0.080</b>      |



# Results Summary

- 48 SNPs genotyped in 392 samples
  - 3 multiplexed assays: 15-plex, 16-plex, 17-plex, total of 12 luminex reaction plates
  - All genotyping was completed within 2 days
  - 45/48 (95.6%) converted to working assays, standard conditions with no optimization required
  - 2 markers were non-polymorphic in the Canadian pedigrees
  - 43 markers analyzed for genetic association to schizophrenia or schizoaffective disorder



# Association Analysis

- Posterior probability of linkage disequilibrium (PPLD)
  - A Bayesian method which estimates the probability of linkage disequilibrium between a marker and a causal mutation
  - Integrates out nuisance parameters related to mode of inheritance (MOI) of disease, allowing for a “model-free” analysis
  - The prior probability is set to 2%; values above 2% represent evidence for LD, below 2% represent evidence against LD
  - Simulations indicate that 95% of PPLD values are below 4%; >99% are below 10%

| Marker  | PPLD L narrow | PPLD L broad | Marker  | PPLD L narrow | PPLD L broad |
|---------|---------------|--------------|---------|---------------|--------------|
| rs01271 | 0.0190        | 0.0178       | rs64149 | 0.0132        | 0.0131       |
| rs07539 | 0.0181        | 0.0190       | rs75427 | 0.0300        | 0.0185       |
| rs11768 | 0.0153        | 0.0166       | rs77494 | 0.0175        | 0.0181       |
| rs20360 | 0.0234        | 0.0226       | rs82136 | 0.0248        | 0.0151       |
| rs24834 | 0.0148        | 0.0152       | rs88303 | 0.0175        | 0.0165       |
| rs35575 | 0.0144        | 0.0165       | rs89885 | 0.0147        | 0.0131       |
| rs36147 | 0.0400        | 0.0221       | rs11895 | 0.0168        | 0.0181       |
| rs53550 | 0.0159        | 0.0180       | rs12311 | 0.0158        | 0.0150       |
| rs69391 | 0.0167        | 0.0177       | rs15793 | 0.0154        | 0.0197       |
| rs75797 | 0.0177        | 0.0177       | rs17507 | 0.0161        | 0.0149       |
| rs79974 | 0.0185        | 0.0165       | rs18910 | 0.0102        | 0.0135       |
| rs80104 | 0.0145        | 0.0121       | rs23444 | 0.0600        | 0.0197       |
| rs98566 | 0.0154        | 0.0197       | rs24663 | 0.0188        | 0.0179       |
| rs98932 | 0.0151        | 0.0114       | rs28687 | 0.0115        | 0.0041       |
| rs02294 | 0.0152        | 0.0155       | rs32305 | 0.0181        | 0.0154       |
| rs11218 | 0.0179        | 0.0171       | rs37530 | 0.0180        | 0.0187       |
| rs11988 | 0.0149        | 0.0132       | rs50892 | 0.0165        | 0.0158       |
| rs25003 | 0.0196        | 0.0182       | rs60120 | 0.0500        | 0.1700       |
| rs34228 | 0.0219        | 0.0213       | rs72916 | 0.0195        | 0.0201       |
| rs43293 | 0.0126        | 0.0144       | rs84430 | 0.0157        | 0.0151       |
| rs57007 | 0.0159        | 0.0160       | rs98931 | 0.0125        | 0.0097       |
| rs63196 | 0.0172        | 0.0165       |         |               |              |



# Results Cont.

| <b>Marker</b>     | <b>PPLD Narrow</b> | <b>PPLD Broad</b> | <b>Gene</b>  |
|-------------------|--------------------|-------------------|--------------|
| <b>rs3736147</b>  | <b>0.0400</b>      | <b>0.0221</b>     | <b>BIN3</b>  |
| <b>rs10923444</b> | <b>0.0600</b>      | <b>0.0197</b>     | <b>GDAP2</b> |
| <b>rs1060120</b>  | <b>0.0500</b>      | <b>0.1700</b>     | <b>H3F3B</b> |



# Conclusions

- We have developed an efficient, robust, and economical medium throughput genotyping assay for the Luminex platform
- Ideally suited for:
  - Candidate gene studies - tagSNP approach
  - Follow-up, high density SNP genotyping after WGAS
  - Custom panels of putative functional variants
    - Functional elements such as PolymiRTS
    - Polymorphisms in drug metabolizing enzymes



# Acknowledgements

- Dr. Linda Brzustowicz
- Dr. Marco Azaro
- Anthony Marcketta
- Dan Lee
- Academic Collaborators
  - Dr. Anne Bassett
  - Dr. Veronica Vieland
- Luminex Corp.
  - Tim Betit
- Perkin Elmer
  - Brian Polizzi
  - Nicole Flaherty