

A multiplexed SNP panel using oligonucleotide ligation assays run on the N-PLEX™ platform for the allelic assignment of genetic risk factors of lung cancer development

Seth B. Harkins, Timothy J. Break, Yui Machida, Mariluz Rodriguez, Galina Nikolenko, John Kenten, Robert M. Umek, and Jacob N. Wohlstadter
Meso Scale Discovery, Rockville, Maryland, USA

1 Introduction

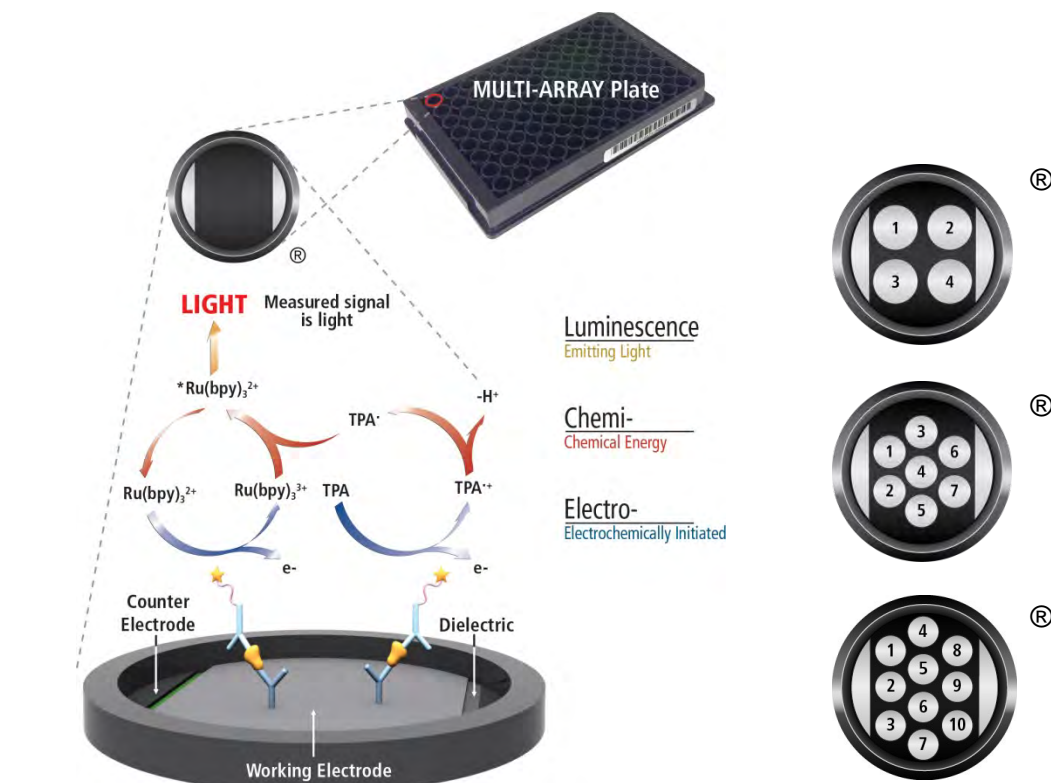
Lung cancer is the leading cause of cancer death worldwide, with an estimate of more than 1 million lung cancer deaths per year. Aggressive treatment strategies are almost always needed to completely eradicate the disease; however, even with aggressive treatment, many lung cancer patients have low 5-year survival expectancies. As a result, efforts are being made to be able to predict which individuals may be more at risk for development of lung cancer (and other cancers) on a genetic level. This information, along with lifestyle and environmental factors like smoking, pollution, etc., may enhance the ability of healthcare professionals to catch this deadly disease at an earlier stage, thus improving patient outcomes.

Genomic variations are common in the general population, with most of these changes being benign or silent. However, in some cases, even individual nucleotide changes can have a profound impact on homeostasis and disease. Consequently, nucleic acid assays for single nucleotide polymorphisms (SNPs) have become a mainstay in most disease-related research. Unfortunately, the routine analysis for genomic variants can be challenging and tedious, making the development of new technology to simplify this process and allow for high-throughput assessment of SNPs in large sample sets highly beneficial. MSD's N-PLEX platform takes advantage of the oligonucleotide ligation assay (OLA) to afford single-nucleotide resolution and pairs it with the robustness of MSD's electrochemiluminescence technology to allow for multiplexed measurements of SNPs in DNA samples. In this study, the N-PLEX platform was used to provide a simple and scalable work-flow for the establishment of a multiplexed panel of assays for nine SNPs that are genetic risk factors for lung cancer development.

2 Methods

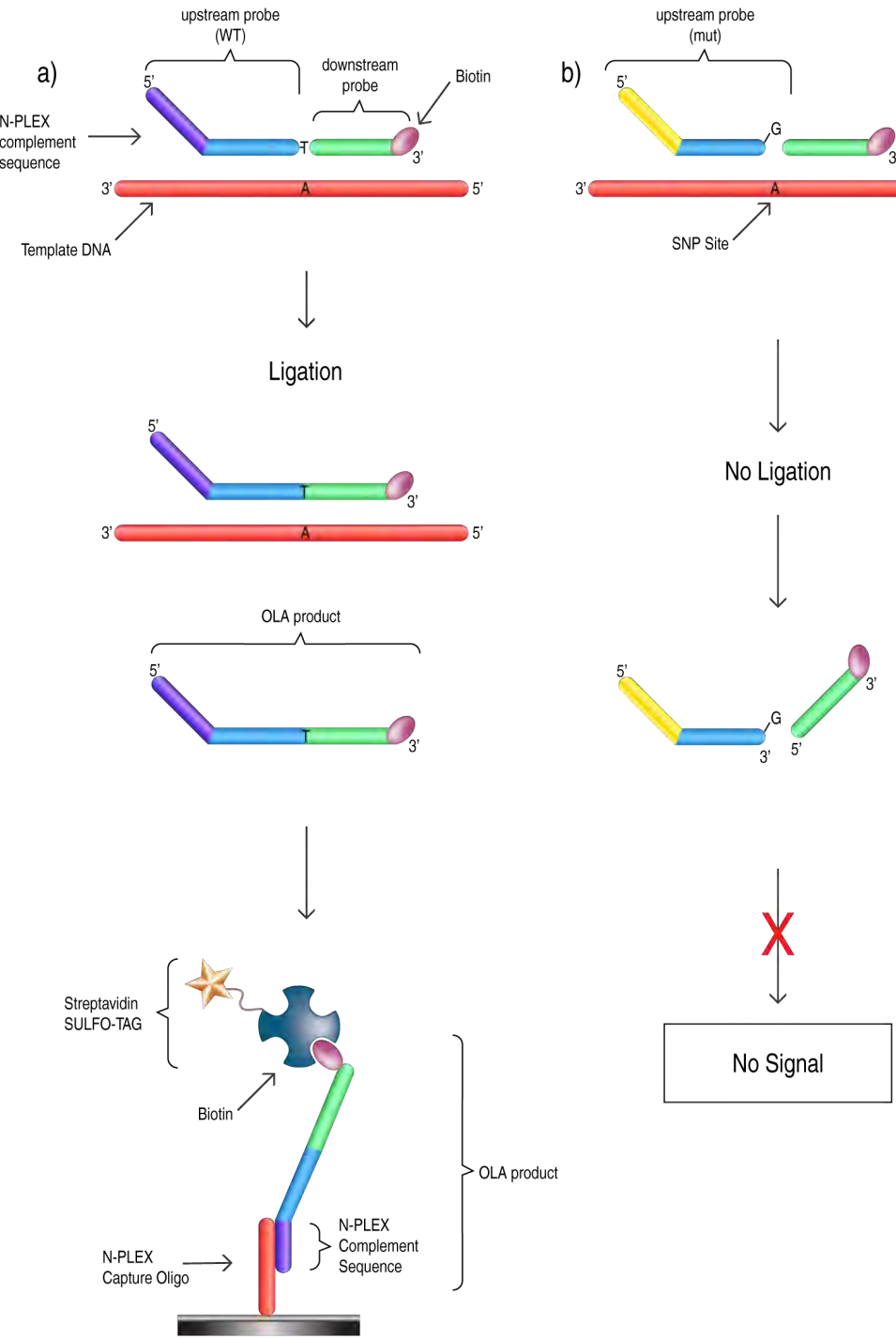
MSD Technology:

MSD's electrochemiluminescence detection technology uses SULFO-TAG™ labels that emit light upon electrochemical stimulation initiated at the electrode surfaces of MULTI-ARRAY® and MULTI-SPOT™ microplates.



SNP detection with OLA and N-PLEX:

The oligonucleotide ligation assay (OLA) used probes that were specific for DNA sequences upstream and downstream of the SNP of interest, with only exact matches at the SNP site allowing for the ligation of the two probes. The ligated probes were hybridized to spot-specific capture oligos on the N-PLEX plates to allow for detection.



Electrochemiluminescence Technology

- Minimal non-specific background and strong responses to analyte yield high signal-to-background ratios.
- The stimulation mechanism (electricity) is decoupled from the response (light signal), minimizing matrix interference.
- Only labels bound near the electrode surface are excited, enabling non-washed assays.
- Labels are stable, non-radioactive, and directly conjugated to biological molecules.
- Emission at ~620 nm eliminates problems with color quenching.
- Multiple rounds of label excitation and emission enhance light levels and improve sensitivity.
- Carbon electrode surface has 10X greater binding capacity than polystyrene wells.
- Surface coatings can be customized.

SNP Detection

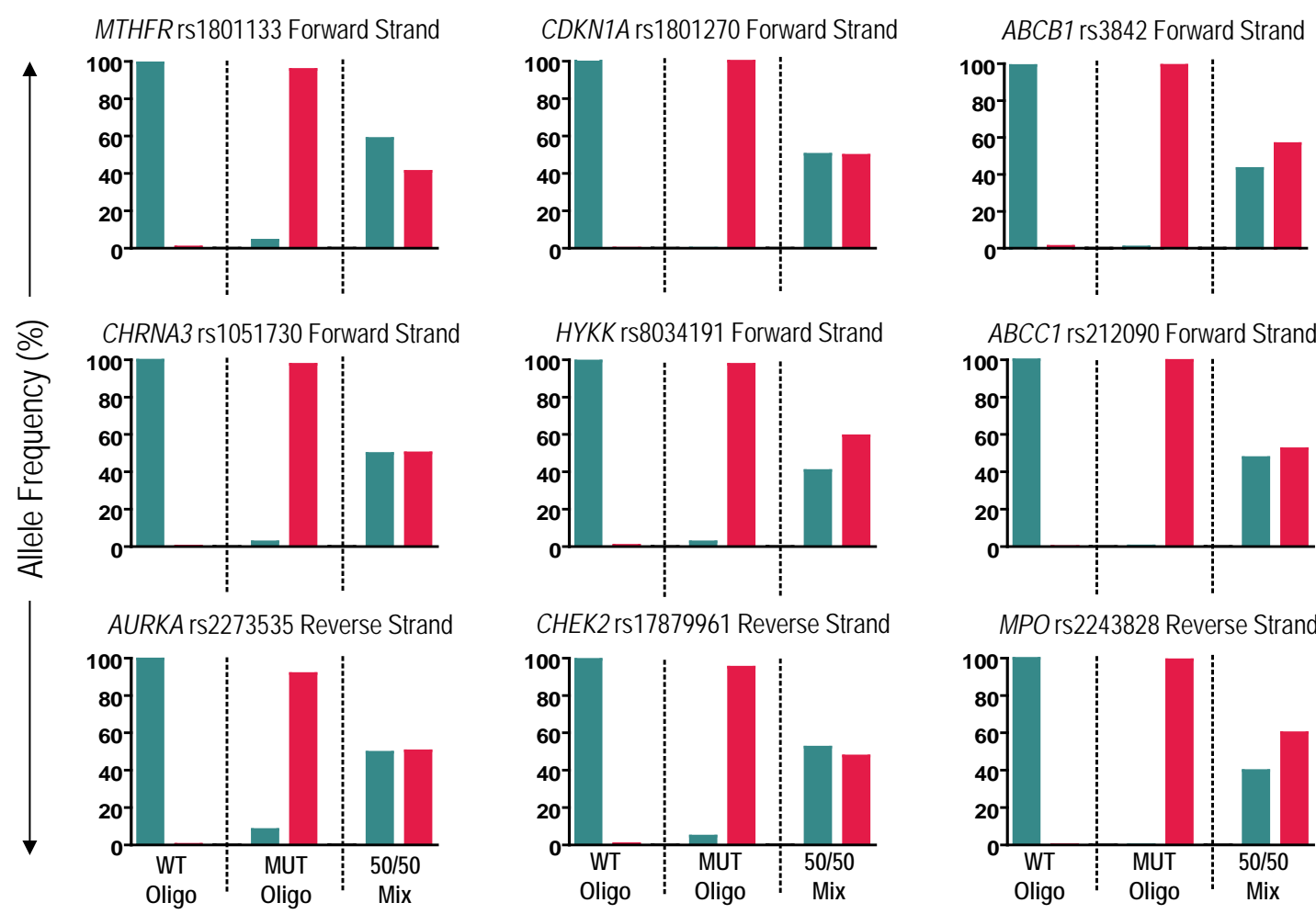
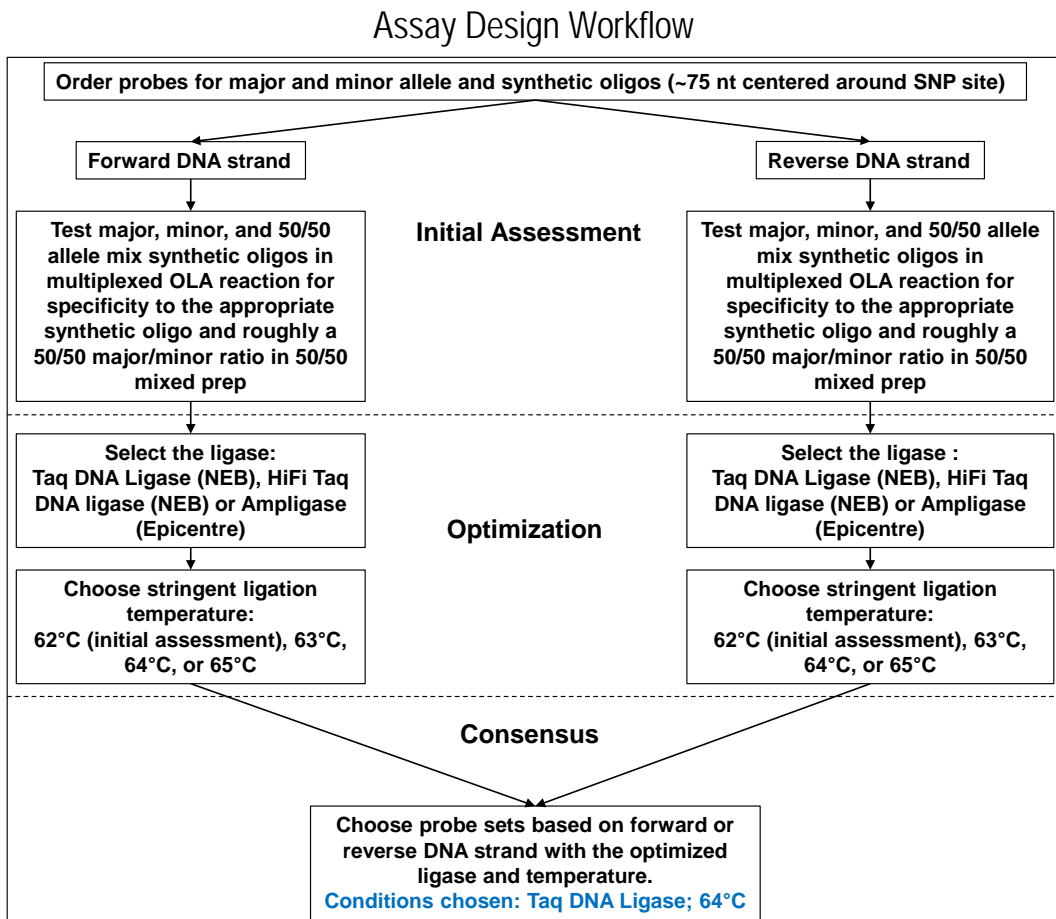
- Each well in an N-PLEX 96-well plate has an array of 10 unique capture oligos attached to the surface of the plate, allowing for the detection of up to five SNPs per well.
- Three probes were needed per target: a biotinylated (downstream) probe and two (upstream) probes that recognized either polymorphic base and contained a sequence complementary to a specific capture oligo.
- Probe characteristics: 17-33 nucleotides in length; optimal probe ligation temperatures between 62-67 °C for a given pair.
- Taq DNA ligase was used to join upstream and downstream probes that aligned correctly on a given DNA sample. Fragments of unmodified template complements were added to prevent bridging of unligated probes.
- OLA products (from synthetic oligos or PCR products from DNA samples) were hybridized to the appropriate capture oligo on the N-PLEX plate, bound by SULFO-TAG labeled streptavidin, and analyzed using an MSD® instrument.
- Synthetic oligos were obtained from IDT
- Blood samples were obtained from BioIVT and DNA was extracted using the QIAamp DNA Blood Mini Kit (Qiagen).
- DNA was extracted from the HL-60 cell line using the PureGene Cell and Tissue Kit (Qiagen).
- For concordance, SNP determination for three SNPs was compared with Taqman SNP Genotyping Assays (Thermo Fisher Scientific), using manufacturer's guidelines.

3 Selected SNPs

SNP	Gene	Chromosome	Global minor allele frequency (GMAF) – 1000 genomes project	Protective/ Deleterious Mutation?
rs1801133	MTHFR	1	0.2454	Deleterious
rs1801270	CDKN1A	6	0.2562	Deleterious
rs3842	ABCB1	7	0.1879	Deleterious
rs1051730	CHRNA3	15	0.1681	Deleterious
rs8034191	HYKK	15	0.1889	Deleterious
rs212090	ABCC1	16	0.2833	Deleterious
rs2273535	AURKA	20	0.3097	Protective
rs17879961	CHEK2	22	0.0010	Protective
rs2243828	MPO	17	0.2270	Protective

SNPs that have been shown to be genetic risk factors (either protective or deleterious) for lung cancer development were chosen that were 1) well-distributed throughout the human genome and 2) well-represented in the general population for the majority of the targets.

4 Results – Assay Design & Optimization



Synthetic oligo templates show high specificity for the appropriate allele using the optimized ligase and temperature conditions of Taq DNA ligase and 64 °C, respectively. WT = Wild-type (major allele), MUT = Mutant (minor allele). All nine assays fit well within our allele frequency guidelines for calling SNPs: Homozygous ≥ 80%, Heterozygous = 30-70%, and Not Present ≤ 20%.



5 Results – Sample Testing

Sample	Allele Frequency % (WT / MUT)									Wild-type	Heterozygous	Mutant	Minor Allele Frequency (MAF)	
	MTHFR	CDKN1A	ABCB1	CHRNA3	HYKK	ABCC1	AURKA	CHEK2	MPO				Number of Alleles	
													Experimental	Global
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11	0.25		
99.4/0.6	46.9/53.2	47.2/52.8	99.0/1.0	99.3/0.7	49.5/50.5	53.9/46.1	99.5/0.5	99.5/0.5	99.5/0.5	99.5/0.5	0.11			