

Table 1

Well	Index ID	Index
A1	iPCRtagT1	ATCACGTT
A2	iPCRtagT2	CGATGTTT
A3	iPCRtagT3	TTAGGCAT
A4	iPCRtagT4	TGACCACT
A5	iPCRtagT5	ACAGTGGT
A6	iPCRtagT6	GCCAATGT
A7	iPCRtagT7	CAGATCTG
A8	iPCRtagT8	ACTTGATG
A9	iPCRtagT9	GATCAGCG
A10	iPCRtagT10	TAGCTTGT
A11	iPCRtagT11	GGCTACAG
A12	iPCRtagT12	CTTGTA CT
B1	iPCRtagT13	TGGTTGTT
B2	iPCRtagT14	TCTCGGTT
B3	iPCRtagT15	TAAGCGTT
B4	iPCRtagT16	TCCGTCTT
B5	iPCRtagT17	TGTACCTT
B6	iPCRtagT18	TTCTGTGT
B7	iPCRtagT19	TCTGCTGT
B8	iPCRtagT20	TTGGAGGT
B9	iPCRtagT21	TCGAGCGT
B10	iPCRtagT22	TGATACGT
B11	iPCRtagT99	GTGCTACC
B12	iPCRtagT101	GGTTGGAC
C1	iPCRtagT25	TGCGATCT
C2	iPCRtagT26	TTCTGCT
C3	iPCRtagT27	TAGTGA CT
C4	iPCRtagT28	TACAGGAT
C5	iPCRtagT29	TCCTCAAT

C6	iPCRtagT30	TGTGGTTG
C7	iPCRtagT31	TAGTCTTG
C8	iPCRtagT32	TTCCATTG
C9	iPCRtagT33	TCGAAGTG
C10	iPCRtagT34	TAACGCTG
C11	iPCRtagT35	TTGGTATG
C12	iPCRtagT36	TGAACTGG
D1	iPCRtagT102	GGCACAAC
D2	iPCRtagT38	TCTCACGG
D3	iPCRtagT39	TCAGGAGG
D4	iPCRtagT40	TAAGTTCG
D5	iPCRtagT41	TCCAGTCG
D6	iPCRtagT42	TGTATGCG
D7	iPCRtagT43	TCATTGAG
D8	iPCRtagT44	TGGCTCAG
D9	iPCRtagT45	TATGCCAG
D10	iPCRtagT46	TCAGATTC
D11	iPCRtagT47	TACTAGTC
D12	iPCRtagT48	TTCAGCTC