

Table 1

Tube	Index ID	Index
1	iPCRtagT1	ATCACGTT
2	iPCRtagT2	CGATGTTT
3	iPCRtagT3	TTAGGCAT
4	iPCRtagT4	TGACCACT
5	iPCRtagT5	ACAGTGGT
6	iPCRtagT6	GCCAATGT
7	iPCRtagT7	CAGATCTG
8	iPCRtagT8	ACTTGATG
9	iPCRtagT9	GATCAGCG
10	iPCRtagT10	TAGCTTGT
11	iPCRtagT11	GGCTACAG
12	iPCRtagT12	CTTGTA CT