

| Plate | Position | Catalog No. of Primer | Accession No. of Gene | Symbol |
|----------|----------|--------------------------|--------------------------|--------|
| QG057-01 | A01 | HQP003654 | NM_181780 | BTLA |
| QG057-01 | A02 | HQP021438 | NM_170662 | CBLB |
| QG057-01 | A03 | HQP053971 | NM_021006 | CCL3L1 |
| QG057-01 | A04 | HQP002209 | NM_005508 | CCR4 |
| QG057-01 | A05 | HQP022667 | NM_001242 | CD27 |
| QG057-01 | A06 | HQP022699 | NM_006139 | CD28 |
| QG057-01 | A07 | HQP022955 | NM_001250 | CD40 |
| QG057-01 | A08 | HQP022962 | NM_000074 | CD40LG |
| QG057-01 | A09 | HQP023108 | NM_001252 | CD70 |
| QG057-01 | A10 | HQP000225 | NM_001798 | CDK2 |
| QG057-01 | A11 | HQP000245 | NM_000075 | CDK4 |
| QG057-01 | A12 | HQP002142 | NM_001836 | CMA1 |
| QG057-01 | B01 | HQP003149 | NM_000757 | CSF1 |
| QG057-01 | B02 | HQP003159 | NM_000758 | CSF2 |
| QG057-01 | B03 | HQP003500 | NM_005214 | CTLA4 |
| QG057-01 | B04 | HQP004613 | NM_000399 | EGR2 |
| QG057-01 | B05 | HQP004616 | NM_004430 | EGR3 |
| QG057-01 | B06 | HQP020132 | NM_005442 | EOMES |
| QG057-01 | B07 | HQP009651 | NM_000043 | FAS |
| QG057-01 | B08 | HQP009671 | NM_000639 | FASLG |
| QG057-01 | B09 | HQP006188 | NM_005252 | FOS |
| QG057-01 | B10 | HQP022662 | NM_014491 | FOXP2 |
| QG057-01 | B11 | HQP012269 | NM_014009 | FOXP3 |
| QG057-01 | B12 | HQP007166 | NM_002051 | GATA3 |
| QG057-01 | C01 | HQP008690 | NM_004131 | GZMB |
| QG057-01 | C02 | HQP009184 | NM_000201 | ICAM1 |
| QG057-01 | C03 | HQP008554 | NM_012092 | ICOS |
| QG057-01 | C04 | HQP009467 | NM_000619 | IFNG |
| QG057-01 | C05 | HQP009685 | NM_000572 | IL10 |
| QG057-01 | C06 | HQP009686 | NM_001558 | IL10RA |
| QG057-01 | C07 | HQP009697 | NM_002188 | IL13 |
| QG057-01 | C08 | HQP009708 | NM_000585 | IL15 |
| QG057-01 | C09 | HQP009717 | NM_002190 | IL17A |
| QG057-01 | C10 | HQP009640 | NM_000575 | IL1A |
| QG057-01 | C11 | HQP009649 | NM_000586 | IL2 |
| QG057-01 | C12 | HQP009650 | NM_000417 | IL2RA |
| QG057-01 | D01 | HQP009658 | NM_000878 | IL2RB |
| QG057-01 | D02 | HQP010154 | NM_001014336 | IL31 |
| QG057-01 | D03 | HQP009662 | NM_000589 | IL4 |
| QG057-01 | D04 | HQP009666 | NM_000879 | IL5 |
| QG057-01 | D05 | HQP009670 | NM_000600 | IL6 |

| | | | | |
|----------|-----|-----------|-----------|-----------|
| QG057-01 | D06 | HQP009677 | NM_002185 | IL7R |
| QG057-01 | D07 | HQP012449 | NM_016162 | ING4 |
| QG057-01 | D08 | HQP009781 | NM_002460 | IRF4 |
| QG057-01 | D09 | HQP020315 | NM_031483 | ITCH |
| QG057-01 | D10 | HQP009793 | NM_181501 | ITGA1 |
| QG057-01 | D11 | HQP009849 | NM_002227 | JAK1 |
| QG057-01 | D12 | HQP009851 | NM_000215 | JAK3 |
| QG057-01 | E01 | HQP009853 | NM_002228 | JUN |
| QG057-01 | E02 | HQP007463 | NM_014387 | LAT |
| QG057-01 | E03 | HQP010581 | NM_000230 | LEP |
| QG057-01 | E04 | HQP010590 | NM_002306 | LGALS3 |
| QG057-01 | E05 | HQP010907 | NM_000595 | LTA |
| QG057-01 | E06 | HQP011149 | NM_005587 | MEF2A |
| QG057-01 | E07 | HQP011788 | NM_172390 | NFATC1 |
| QG057-01 | E08 | HQP011789 | NM_012340 | NFATC2 |
| QG057-01 | E09 | HQP011792 | NM_004555 | NFATC3 |
| QG057-01 | E10 | HQP011807 | NM_003998 | NFKB1 |
| QG057-01 | E11 | HQP011873 | NM_017617 | NOTCH1 |
| QG057-01 | E12 | HQP012662 | NM_005018 | PDCD1 |
| QG057-01 | F01 | HQP054058 | NM_005041 | PRF1 |
| QG057-01 | F02 | HQP014751 | NM_002739 | PRKCG |
| QG057-01 | F03 | HQP015544 | NM_000956 | PTGER2 |
| QG057-01 | F04 | HQP015598 | NM_000963 | PTGS2 |
| QG057-01 | F05 | HQP016745 | NM_000655 | SELL |
| QG057-01 | F06 | HQP017767 | NM_003150 | STAT3 |
| QG057-01 | F07 | HQP017775 | NM_003153 | STAT6 |
| QG057-01 | F08 | HQP008682 | NM_013351 | TBX21 |
| QG057-01 | F09 | HQP018044 | NM_000660 | TGFB1 |
| QG057-01 | F10 | HQP021557 | NM_003844 | TNFRSF10A |
| QG057-01 | F11 | HQP021522 | NM_003820 | TNFRSF14 |
| QG057-01 | F12 | HQP021536 | NM_004195 | TNFRSF18 |
| QG057-01 | G01 | HQP018332 | NM_003327 | TNFRSF4 |
| QG057-01 | G02 | HQP022753 | NM_001243 | TNFRSF8 |
| QG057-01 | G03 | HQP009716 | NM_001561 | TNFRSF9 |
| QG057-01 | G04 | HQP021502 | NM_003810 | TNFSF10 |
| QG057-01 | G05 | HQP021496 | NM_003807 | TNFSF14 |
| QG057-01 | G06 | HQP022769 | NM_001244 | TNFSF8 |
| QG057-01 | G07 | HQP016381 | NM_001101 | ACTB |
| QG057-01 | G08 | HQP015171 | NM_004048 | B2M |
| QG057-01 | G09 | HQP006940 | NM_002046 | GAPDH |
| QG057-01 | G10 | HQP009026 | NM_000194 | HPRT1 |
| QG057-01 | G11 | HQP054097 | NM_001002 | RPLP0 |
| QG057-01 | G12 | HQP021151 | NM_003646 | DGKZ |

| | | | | |
|----------|-----|-----------|-----------|--------|
| QG057-01 | H01 | HGDC | | |
| QG057-01 | H02 | HGDC | | |
| QG057-01 | H03 | HQP006940 | NM_002046 | GAPDH |
| QG057-01 | H04 | HQP016381 | NM_001101 | ACTB |
| QG057-01 | H05 | HQP015171 | NM_004048 | B2M |
| QG057-01 | H06 | HQP006171 | NM_012423 | RPL13A |
| QG057-01 | H07 | HQP009026 | NM_000194 | HPRT1 |
| QG057-01 | H08 | HQP054253 | NR_003286 | RN18S1 |
| QG057-01 | H09 | RT | | |
| QG057-01 | H10 | RT | | |
| QG057-01 | H11 | PCR | | |
| QG057-01 | H12 | PCR | | |