

| Plate | Position | Catalog No. of Primer | Accession No. of Gene | Symbol |
|----------|----------|--------------------------|--------------------------|---------|
| QG094-01 | A01 | HQP014510 | NM_017677 | MTMR8 |
| QG094-01 | A02 | HQP014378 | NM_018498 | PPP4R1L |
| QG094-01 | A03 | HQP013910 | NM_017857 | SSH3 |
| QG094-01 | A04 | HQP013874 | NM_017823 | DUSP23 |
| QG094-01 | A05 | HQP013834 | NM_017762 | MTMR10 |
| QG094-01 | A06 | HQP013561 | NM_001040446 | MTMR12 |
| QG094-01 | A07 | HQP013463 | NM_018984 | SSH1 |
| QG094-01 | A08 | HQP012967 | NM_016226 | VPS29 |
| QG094-01 | A09 | HQP012940 | NM_016086 | STYXL1 |
| QG094-01 | A10 | HQP012781 | NM_016396 | CTDSPL2 |
| QG094-01 | A11 | HQP012516 | NM_016361 | ACP6 |
| QG094-01 | A12 | HQP008416 | NM_014188 | SSU72 |
| QG094-01 | B01 | HQP008395 | NM_014172 | PHPT1 |
| QG094-01 | B02 | HQP007815 | NM_013239 | PPP2R3B |
| QG094-01 | B03 | HQP007528 | NM_001002837 | PIB5PA |
| QG094-01 | B04 | HQP007478 | NM_014395 | DAPP1 |
| QG094-01 | B05 | HQP007244 | NM_014369 | PTPN18 |
| QG094-01 | B06 | HQP007113 | NM_012411 | PTPN22 |
| QG094-01 | B07 | HQP006898 | NM_015466 | PTPN23 |
| QG094-01 | B08 | HQP006041 | NM_015319 | TENC1 |
| QG094-01 | B09 | HQP005901 | NM_194449 | PHLPP |
| QG094-01 | B10 | HQP005691 | NM_015020 | PHLPPL |
| QG094-01 | B11 | HQP005535 | NM_014937 | INPP5F |
| QG094-01 | B12 | HQP005496 | NM_014906 | PPM1E |
| QG094-01 | C01 | HQP001479 | NM_007240 | DUSP12 |
| QG094-01 | C02 | HQP001419 | NM_007207 | DUSP10 |
| QG094-01 | C03 | HQP001332 | NM_007079 | PTP4A3 |
| QG094-01 | C04 | HQP001285 | NM_007050 | PTPRT |
| QG094-01 | C05 | HQP001261 | NM_007039 | PTPN21 |
| QG094-01 | C06 | HQP001227 | NM_007026 | DUSP14 |
| QG094-01 | C07 | HQP023452 | NM_001042388 | PPP4R1 |
| QG094-01 | C08 | HQP023249 | NM_014787 | DNAJC6 |
| QG094-01 | C09 | HQP023200 | NM_014750 | DLG7 |
| QG094-01 | C10 | HQP023028 | NM_014634 | PPM1F |
| QG094-01 | C11 | HQP022913 | NM_004897 | MINPP1 |
| QG094-01 | C12 | HQP022301 | NM_014241 | PTPLA |
| QG094-01 | D01 | HQP022191 | NM_004715 | CTDP1 |
| QG094-01 | D02 | HQP022092 | NM_004687 | MTMR4 |
| QG094-01 | D03 | HQP022089 | NM_004686 | MTMR7 |
| QG094-01 | D04 | HQP022088 | NM_004685 | MTMR6 |
| QG094-01 | D05 | HQP021686 | NM_016156 | MTMR2 |

| | | | | |
|----------|-----|-----------|--------------|--------|
| QG094-01 | D06 | HQP021683 | NM_021090 | MTMR3 |
| QG094-01 | D07 | HQP021548 | NM_003837 | FBP2 |
| QG094-01 | D08 | HQP021531 | NM_003828 | MTMR1 |
| QG094-01 | D09 | HQP021487 | NM_003800 | RNGTT |
| QG094-01 | D10 | HQP021340 | NM_003713 | PPAP2B |
| QG094-01 | D11 | HQP021335 | NM_003711 | PPAP2A |
| QG094-01 | D12 | HQP021276 | NM_003672 | CDC14A |
| QG094-01 | E01 | HQP021112 | NM_003626 | PPFIA1 |
| QG094-01 | E02 | HQP021090 | NM_015053 | PPFIA4 |
| QG094-01 | E03 | HQP021053 | NM_003620 | PPM1D |
| QG094-01 | E04 | HQP020754 | NM_003584 | DUSP11 |
| QG094-01 | E05 | HQP019733 | NM_080391 | PTP4A2 |
| QG094-01 | E06 | HQP019066 | NM_001018041 | EPM2A |
| QG094-01 | E07 | HQP018760 | NM_003463 | PTP4A1 |
| QG094-01 | E08 | HQP018223 | NM_199259 | TPTE |
| QG094-01 | E09 | HQP018167 | NM_022648 | TNS1 |
| QG094-01 | E10 | HQP017809 | NM_145251 | STYX |
| QG094-01 | E11 | HQP016574 | NM_002972 | SBF1 |
| QG094-01 | E12 | HQP015943 | NM_002851 | PTPRZ1 |
| QG094-01 | F01 | HQP015939 | NM_002850 | PTPRS |
| QG094-01 | F02 | HQP015937 | NM_002849 | PTPRR |
| QG094-01 | F03 | HQP015933 | NM_002848 | PTPRO |
| QG094-01 | F04 | HQP015929 | NM_002847 | PTPRN2 |
| QG094-01 | F05 | HQP015927 | NM_002845 | PTPRM |
| QG094-01 | F06 | HQP015926 | NM_002844 | PTPRK |
| QG094-01 | F07 | HQP015925 | NM_002843 | PTPRJ |
| QG094-01 | F08 | HQP015924 | NM_002842 | PTPRH |
| QG094-01 | F09 | HQP015921 | NM_002840 | PTPRF |
| QG094-01 | F10 | HQP015919 | NM_006504 | PTPRE |
| QG094-01 | F11 | HQP015912 | NM_001040712 | PTPRD |
| QG094-01 | F12 | HQP015908 | NM_002838 | PTPRC |
| QG094-01 | G01 | HQP015907 | NM_002837 | PTPRB |
| QG094-01 | G02 | HQP015897 | NM_006264 | PTPN13 |
| QG094-01 | G03 | HQP015890 | NM_002835 | PTPN12 |
| QG094-01 | G04 | HQP015878 | NM_002834 | PTPN11 |
| QG094-01 | G05 | HQP015873 | NM_002833 | PTPN9 |
| QG094-01 | G06 | HQP015863 | NM_002832 | PTPN7 |
| QG094-01 | G07 | HQP015856 | NM_002830 | PTPN4 |
| QG094-01 | G08 | HQP015854 | NM_002829 | PTPN3 |
| QG094-01 | G09 | HQP015839 | NM_002828 | PTPN2 |
| QG094-01 | G10 | HQP015828 | NM_002827 | PTPN1 |
| QG094-01 | G11 | HQP015535 | NM_000314 | PTEN |
| QG094-01 | G12 | HQP014907 | NM_002755 | MAP2K1 |

| | | | | |
|----------|-----|-----------|--------------|----------|
| QG094-01 | H01 | HGDC | | |
| QG094-01 | H02 | HGDC | | |
| QG094-01 | H03 | HQP006940 | NM_002046 | GAPDH |
| QG094-01 | H04 | HQP016381 | NM_001101 | ACTB |
| QG094-01 | H05 | HQP015171 | NM_004048 | B2M |
| QG094-01 | H06 | HQP006171 | NM_012423 | RPL13A |
| QG094-01 | H07 | HQP009026 | NM_000194 | HPRT1 |
| QG094-01 | H08 | HQP054253 | NR_003286 | RN18S1 |
| QG094-01 | H09 | RT | | |
| QG094-01 | H10 | RT | | |
| QG094-01 | H11 | PCR | | |
| QG094-01 | H12 | PCR | | |
| QG094-02 | A01 | HQP014381 | NM_002721 | PPP6C |
| QG094-02 | A02 | HQP014377 | NM_006247 | PPP5C |
| QG094-02 | A03 | HQP014356 | NM_000945 | PPP3R1 |
| QG094-02 | A04 | HQP014346 | NM_005605 | PPP3CC |
| QG094-02 | A05 | HQP014334 | NM_021132 | PPP3CB |
| QG094-02 | A06 | HQP014321 | NM_002720 | PPP4C |
| QG094-02 | A07 | HQP014309 | NM_000944 | PPP3CA |
| QG094-02 | A08 | HQP014301 | NM_006246 | PPP2R5E |
| QG094-02 | A09 | HQP014282 | NM_006245 | PPP2R5D |
| QG094-02 | A10 | HQP014262 | NM_002719 | PPP2R5C |
| QG094-02 | A11 | HQP014251 | NM_006244 | PPP2R5B |
| QG094-02 | A12 | HQP014244 | NM_006243 | PPP2R5A |
| QG094-02 | B01 | HQP014212 | NM_002718 | PPP2R3A |
| QG094-02 | B02 | HQP014202 | NM_020416 | PPP2R2C |
| QG094-02 | B03 | HQP014174 | NM_002717 | PPP2R2A |
| QG094-02 | B04 | HQP014161 | NM_002716 | PPP2R1B |
| QG094-02 | B05 | HQP014115 | NM_002715 | PPP2CA |
| QG094-02 | B06 | HQP014033 | NM_005398 | PPP1R3C |
| QG094-02 | B07 | HQP013980 | NM_002710 | PPP1CC |
| QG094-02 | B08 | HQP013970 | NM_002709 | PPP1CB |
| QG094-02 | B09 | HQP013953 | NM_001008709 | PPP1CA |
| QG094-02 | B10 | HQP013893 | NM_021003 | PPM1A |
| QG094-02 | B11 | HQP013674 | NM_006240 | PPEF1 |
| QG094-02 | B12 | HQP013643 | NM_006239 | PPEF2 |
| QG094-02 | C01 | HQP013613 | NM_021129 | PPA1 |
| QG094-02 | C02 | HQP013073 | NM_004567 | PFKFB4 |
| QG094-02 | C03 | HQP013072 | NM_004566 | PFKFB3 |
| QG094-02 | C04 | HQP012011 | NM_000276 | OCRL |
| QG094-02 | C05 | HQP011657 | NM_002481 | PPP1R12B |
| QG094-02 | C06 | HQP011625 | NM_001002841 | MYL4 |
| QG094-02 | C07 | HQP011549 | NM_000252 | MTM1 |

| | | | | |
|----------|-----|-----------|--------------|--------|
| QG094-02 | C08 | HQP010565 | NM_001042771 | LCK |
| QG094-02 | C09 | HQP009837 | NM_033453 | ITPA |
| QG094-02 | C10 | HQP009756 | NM_001567 | INPPL1 |
| QG094-02 | C11 | HQP009752 | NM_005539 | INPP5A |
| QG094-02 | C12 | HQP009750 | NM_001566 | INPP4A |
| QG094-02 | D01 | HQP009747 | NM_002194 | INPP1 |
| QG094-02 | D02 | HQP009727 | NM_014214 | IMPA2 |
| QG094-02 | D03 | HQP009726 | NM_005536 | IMPA1 |
| QG094-02 | D04 | HQP005224 | NM_000507 | FBP1 |
| QG094-02 | D05 | HQP005048 | NM_001990 | EYA3 |
| QG094-02 | D06 | HQP005044 | NM_005244 | EYA2 |
| QG094-02 | D07 | HQP005041 | NM_000503 | EYA1 |
| QG094-02 | D08 | HQP004507 | NM_004420 | DUSP8 |
| QG094-02 | D09 | HQP004506 | NM_001947 | DUSP7 |
| QG094-02 | D10 | HQP004504 | NM_001946 | DUSP6 |
| QG094-02 | D11 | HQP004503 | NM_004419 | DUSP5 |
| QG094-02 | D12 | HQP004501 | NM_001394 | DUSP4 |
| QG094-02 | E01 | HQP004500 | NM_004090 | DUSP3 |
| QG094-02 | E02 | HQP004498 | NM_004417 | DUSP1 |
| QG094-02 | E03 | HQP004337 | NM_004087 | DLG1 |
| QG094-02 | E04 | HQP000418 | NM_005192 | CDKN3 |
| QG094-02 | E05 | HQP023409 | NM_001790 | CDC25C |
| QG094-02 | E06 | HQP023399 | NM_004358 | CDC25B |
| QG094-02 | E07 | HQP023385 | NM_001789 | CDC25A |
| QG094-02 | E08 | HQP006448 | NM_031313 | ALPPL2 |
| QG094-02 | E09 | HQP006440 | NM_000478 | ALPL |
| QG094-02 | E10 | HQP006433 | NM_001631 | ALPI |
| QG094-02 | E11 | HQP004724 | NM_001623 | AIF1 |
| QG094-02 | E12 | HQP023341 | NM_138448 | ACYP2 |
| QG094-02 | F01 | HQP023213 | NM_001107 | ACYP1 |
| QG094-02 | F02 | HQP014890 | NM_001099 | ACPP |
| QG094-02 | F03 | HQP013957 | NM_001611 | ACP5 |
| QG094-02 | F04 | HQP013357 | NM_001610 | ACP2 |
| QG094-02 | F05 | HQP021106 | NM_003625 | PPFIA2 |
| QG094-02 | F06 | HQP013165 | NM_001040649 | ACP1 |
| QG094-02 | F07 | HQP021272 | NM_001077181 | CDC14B |
| QG094-02 | F08 | HQP020019 | NM_001222 | CAMK2G |
| QG094-02 | F09 | HQP004499 | NM_004418 | DUSP2 |
| QG094-02 | F10 | HQP004508 | NM_001395 | DUSP9 |
| QG094-02 | F11 | HQP004980 | NM_004100 | EYA4 |
| QG094-02 | F12 | HQP013069 | NM_002625 | PFKFB1 |
| QG094-02 | G01 | HQP013070 | NM_001018053 | PFKFB2 |
| QG094-02 | G02 | HQP013904 | NM_001033556 | PPM1B |

| | | | | |
|----------|-----|-----------|--------------|---------|
| QG094-02 | G03 | HQP013921 | NM_002707 | PPM1G |
| QG094-02 | G04 | HQP014048 | NM_006242 | PPP1R3D |
| QG094-02 | G05 | HQP014125 | NM_001009552 | PPP2CB |
| QG094-02 | G06 | HQP014187 | NM_004576 | PPP2R2B |
| QG094-02 | G07 | HQP014229 | NM_021131 | PPP2R4 |
| QG094-02 | G08 | HQP015523 | NM_004577 | PSPH |
| QG094-02 | G09 | HQP015859 | NM_002831 | PTPN6 |
| QG094-02 | G10 | HQP015901 | NM_005401 | PTPN14 |
| QG094-02 | G11 | HQP021337 | NM_003712 | PPAP2C |
| QG094-02 | G12 | HQP015923 | NM_002841 | PTPRG |
| QG094-02 | H01 | HGDC | | |
| QG094-02 | H02 | HGDC | | |
| QG094-02 | H03 | HQP006940 | NM_002046 | GAPDH |
| QG094-02 | H04 | HQP016381 | NM_001101 | ACTB |
| QG094-02 | H05 | HQP015171 | NM_004048 | B2M |
| QG094-02 | H06 | HQP006171 | NM_012423 | RPL13A |
| QG094-02 | H07 | HQP009026 | NM_000194 | HPRT1 |
| QG094-02 | H08 | HQP054253 | NR_003286 | RN18S1 |
| QG094-02 | H09 | RT | | |
| QG094-02 | H10 | RT | | |
| QG094-02 | H11 | PCR | | |
| QG094-02 | H12 | PCR | | |