

| Plate | Position | Catalog No. of Primer | Accession No. of Gene | Symbol |
|----------|----------|--------------------------|--------------------------|----------|
| QG005-01 | A01 | HQP009653 | NM_152872 | FAS |
| QG005-01 | A02 | HQP022974 | NM_001001390 | CD44 |
| QG005-01 | A03 | HQP009026 | NM_000194 | HPRT1 |
| QG005-01 | A04 | HQP009802 | NM_001011645 | AR |
| QG005-01 | A05 | HQP018481 | NM_003376 | VEGFA |
| QG005-01 | A06 | HQP018175 | NM_000546 | TP53 |
| QG005-01 | A07 | HQP018172 | NM_001067 | TOP2A |
| QG005-01 | A08 | HQP018144 | NM_006291 | TNFAIP2 |
| QG005-01 | A09 | HQP018091 | NM_005424 | TIE1 |
| QG005-01 | A10 | HQP018069 | NM_003247 | THBS2 |
| QG005-01 | A11 | HQP018068 | NM_003246 | THBS1 |
| QG005-01 | A12 | HQP018043 | NM_003236 | TGFA |
| QG005-01 | B01 | HQP018035 | NM_003225 | TFF1 |
| QG005-01 | B02 | HQP021341 | NM_003714 | STC2 |
| QG005-01 | B03 | HQP017678 | NM_003125 | SPRR1B |
| QG005-01 | B04 | HQP012154 | NM_000602 | SERPINE1 |
| QG005-01 | B05 | HQP013125 | NM_002639 | SERPINB5 |
| QG005-01 | B06 | HQP002590 | NM_001085 | SERPINA3 |
| QG005-01 | B07 | HQP011193 | NM_002407 | SCGB2A1 |
| QG005-01 | B08 | HQP016429 | NM_000988 | RPL27 |
| QG005-01 | B09 | HQP016067 | NM_002872 | RAC2 |
| QG005-01 | B10 | HQP015598 | NM_000963 | PTGS2 |
| QG005-01 | B11 | HQP015535 | NM_000314 | PTEN |
| QG005-01 | B12 | HQP013204 | NM_002658 | PLAU |
| QG005-01 | C01 | HQP013099 | NM_000926 | PGR |
| QG005-01 | C02 | HQP012189 | NM_002581 | PAPPA |
| QG005-01 | C03 | HQP011853 | NM_000269 | NME1 |
| QG005-01 | C04 | HQP011828 | NM_002507 | NGFR |
| QG005-01 | C05 | HQP011827 | NM_002506 | NGFB |
| QG005-01 | C06 | HQP011825 | NM_006166 | NFYB |
| QG005-01 | C07 | HQP011559 | NM_001018016 | MUC1 |
| QG005-01 | C08 | HQP011539 | NM_005954 | MT3 |
| QG005-01 | C09 | HQP011232 | NM_002417 | MKI67 |
| QG005-01 | C10 | HQP014926 | NM_145185 | MAP2K7 |
| QG005-01 | C11 | HQP010233 | NM_002276 | KRT19 |
| QG005-01 | C12 | HQP010197 | NM_000224 | KRT18 |
| QG005-01 | D01 | HQP054014 | NM_012427 | KLK5 |
| QG005-01 | D02 | HQP017898 | NM_001730 | KLF5 |
| QG005-01 | D03 | HQP010099 | NM_000222 | KIT |
| QG005-01 | D04 | HQP009853 | NM_002228 | JUN |
| QG005-01 | D05 | HQP009819 | NM_000213 | ITGB4 |

| | | | | |
|----------|-----|-----------|-----------|---------|
| QG005-01 | D06 | HQP009774 | NM_000210 | ITGA6 |
| QG005-01 | D07 | HQP009674 | NM_002184 | IL6ST |
| QG005-01 | D08 | HQP009672 | NM_000565 | IL6R |
| QG005-01 | D09 | HQP009670 | NM_000600 | IL6 |
| QG005-01 | D10 | HQP009650 | NM_000417 | IL2RA |
| QG005-01 | D11 | HQP009541 | NM_000597 | IGFBP2 |
| QG005-01 | D12 | HQP009273 | NM_002166 | ID2 |
| QG005-01 | E01 | HQP009089 | NM_001540 | HSPB1 |
| QG005-01 | E02 | HQP008883 | NM_002128 | HMGB1 |
| QG005-01 | E03 | HQP007166 | NM_002051 | GATA3 |
| QG005-01 | E04 | HQP006660 | NM_014211 | GABRP |
| QG005-01 | E05 | HQP019708 | NM_005438 | FOSL1 |
| QG005-01 | E06 | HQP005400 | NM_000800 | FGF1 |
| QG005-01 | E07 | HQP009671 | NM_000639 | FASLG |
| QG005-01 | E08 | HQP009651 | NM_000043 | FAS |
| QG005-01 | E09 | HQP005002 | NM_001437 | ESR2 |
| QG005-01 | E10 | HQP004998 | NM_000125 | ESR1 |
| QG005-01 | E11 | HQP004969 | NM_004448 | ERBB2 |
| QG005-01 | E12 | HQP004605 | NM_005228 | EGFR |
| QG005-01 | F01 | HQP000462 | NM_006094 | DLC1 |
| QG005-01 | F02 | HQP003904 | NM_000103 | CYP19A1 |
| QG005-01 | F03 | HQP003599 | NM_001909 | CTSD |
| QG005-01 | F04 | HQP003596 | NM_001908 | CTSB |
| QG005-01 | F05 | HQP003539 | NM_001904 | CTNNB1 |
| QG005-01 | F06 | HQP002549 | NM_001848 | COL6A1 |
| QG005-01 | F07 | HQP002037 | NM_001831 | CLU |
| QG005-01 | F08 | HQP000369 | NM_000077 | CDKN2A |
| QG005-01 | F09 | HQP000342 | NM_004064 | CDKN1B |
| QG005-01 | F10 | HQP000331 | NM_000389 | CDKN1A |
| QG005-01 | F11 | HQP022972 | NM_000610 | CD44 |
| QG005-01 | F12 | HQP021819 | NM_001238 | CCNE1 |
| QG005-01 | G01 | HQP016204 | NM_053056 | CCND1 |
| QG005-01 | G02 | HQP021701 | NM_001237 | CCNA2 |
| QG005-01 | G03 | HQP021692 | NM_003914 | CCNA1 |
| QG005-01 | G04 | HQP018238 | NM_000064 | C3 |
| QG005-01 | G05 | HQP016260 | NM_004050 | BCL2L2 |
| QG005-01 | G06 | HQP016211 | NM_000633 | BCL2 |
| QG005-01 | G07 | HQP015574 | NM_004323 | BAG1 |
| QG005-01 | G08 | HQP015538 | NM_004322 | BAD |
| QG005-01 | G09 | HQP009801 | NM_000044 | AR |
| QG005-01 | G10 | HQP023466 | NM_004360 | CDH1 |
| QG005-01 | G11 | HQP002852 | NM_001307 | CLDN7 |
| QG005-01 | G12 | HQP006343 | NM_013280 | FLRT1 |

| | | | | |
|----------|-----|-----------|-----------|--------|
| QG005-01 | H01 | HGDC | | |
| QG005-01 | H02 | HGDC | | |
| QG005-01 | H03 | HQP006940 | NM_002046 | GAPDH |
| QG005-01 | H04 | HQP016381 | NM_001101 | ACTB |
| QG005-01 | H05 | HQP015171 | NM_004048 | B2M |
| QG005-01 | H06 | HQP006171 | NM_012423 | RPL13A |
| QG005-01 | H07 | HQP009026 | NM_000194 | HPRT1 |
| QG005-01 | H08 | HQP054253 | NR_003286 | RN18S1 |
| QG005-01 | H09 | RT | | |
| QG005-01 | H10 | RT | | |
| QG005-01 | H11 | PCR | | |
| QG005-01 | H12 | PCR | | |