

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG021-01	A01	HQP021627	NM_003888	ALDH1A2
QG021-01	A02	HQP015964	NM_004324	BAX
QG021-01	A03	HQP009099	NM_001012270	BIRC5
QG021-01	A04	HQP017983	NM_001731	BTG1
QG021-01	A05	HQP018966	NM_001080124	CASP8
QG021-01	A06	HQP022955	NM_001250	CD40
QG021-01	A07	HQP003218	NM_004383	CSK
QG021-01	A08	HQP003488	NM_001901	CTGF
QG021-01	A09	HQP003539	NM_001904	CTNNB1
QG021-01	A10	HQP004524	NM_005225	E2F1
QG021-01	A11	HQP004528	NM_001950	E2F4
QG021-01	A12	HQP004557	NM_001955	EDN1
QG021-01	B01	HQP004559	NM_000114	EDN3
QG021-01	B02	HQP004564	NM_000115	EDNRB
QG021-01	B03	HQP004599	NM_001963	EGF
QG021-01	B04	HQP004605	NM_005228	EGFR
QG021-01	B05	HQP004968	NM_001005862	ERBB2
QG021-01	B06	HQP005001	NM_001040275	ESR2
QG021-01	B07	HQP005400	NM_000800	FGF1
QG021-01	B08	HQP005417	NM_004465	FGF10
QG021-01	B09	HQP005403	NM_002006	FGF2
QG021-01	B10	HQP005408	NM_004464	FGF5
QG021-01	B11	HQP005411	NM_002009	FGF7
QG021-01	B12	HQP005412	NM_006119	FGF8
QG021-01	C01	HQP005416	NM_002010	FGF9
QG021-01	C02	HQP005437	NM_000141	FGFR2
QG021-01	C03	HQP005434	NM_000142	FGFR3
QG021-01	C04	HQP005439	NM_002011	FGFR4
QG021-01	C05	HQP006188	NM_005252	FOS
QG021-01	C06	HQP005712	NM_021953	FOXM1
QG021-01	C07	HQP006930	NM_002045	GAP43
QG021-01	C08	HQP004493	NM_001945	HBEGF
QG021-01	C09	HQP008800	NM_000601	HGF
QG021-01	C10	HQP009026	NM_000194	HPRT1
QG021-01	C11	HQP000630	NM_006388	HTATIP
QG021-01	C12	HQP009518	NM_000618	IGF1
QG021-01	D01	HQP009523	NM_000875	IGF1R
QG021-01	D02	HQP009529	NM_000612	IGF2
QG021-01	D03	HQP009539	NM_000596	IGFBP1
QG021-01	D04	HQP009541	NM_000597	IGFBP2
QG021-01	D05	HQP009544	NM_000598	IGFBP3

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG021-01	D06	HQP009548	NM_001552	IGFBP4
QG021-01	D07	HQP009551	NM_000599	IGFBP5
QG021-01	D08	HQP009555	NM_002178	IGFBP6
QG021-01	D09	HQP009558	NM_001553	IGFBP7
QG021-01	D10	HQP010907	NM_000595	LTA
QG021-01	D11	HQP014848	NM_002745	MAPK1
QG021-01	D12	HQP011071	NM_005368	MB
QG021-01	E01	HQP011219	NM_002415	MIF
QG021-01	E02	HQP011774	NM_000267	NF1
QG021-01	E03	HQP011827	NM_002506	NGFB
QG021-01	E04	HQP011828	NM_002507	NGFR
QG021-01	E05	HQP012197	NM_002583	PAWR
QG021-01	E06	HQP012847	NM_002607	PDGFA
QG021-01	E07	HQP012889	NM_002609	PDGFRB
QG021-01	E08	HQP013099	NM_000926	PGR
QG021-01	E09	HQP013193	NM_000928	PLA2G1B
QG021-01	E10	HQP013194	NM_000300	PLA2G2A
QG021-01	E11	HQP013238	NM_002660	PLCG1
QG021-01	E12	HQP013627	NM_006238	PPARD
QG021-01	F01	HQP015530	NM_000264	PTCH1
QG021-01	F02	HQP015535	NM_000314	PTEN
QG021-01	F03	HQP016063	NM_006908	RAC1
QG021-01	F04	HQP016089	NM_000448	RAG1
QG021-01	F05	HQP016819	NM_005410	SEPP1
QG021-01	F06	HQP017098	NM_000193	SHH
QG021-01	F07	HQP006080	NM_012238	SIRT1
QG021-01	F08	HQP006284	NM_012444	SPO11
QG021-01	F09	HQP017698	NM_000348	SRD5A2
QG021-01	F10	HQP018043	NM_003236	TGFA
QG021-01	F11	HQP018044	NM_000660	TGFB1
QG021-01	F12	HQP018045	NM_001042454	TGFB111
QG021-01	G01	HQP018047	NM_003238	TGFB2
QG021-01	G02	HQP018048	NM_003239	TGFB3
QG021-01	G03	HQP018051	NM_004612	TGFB1
QG021-01	G04	HQP018060	NM_000359	TGM1
QG021-01	G05	HQP018141	NM_000594	TNF
QG021-01	G06	HQP018149	NM_001066	TNFRSF1B
QG021-01	G07	HQP022985	NM_004295	TRAF4
QG021-01	G08	HQP018336	NM_003330	TXNRD1
QG021-01	G09	HQP018474	NM_000376	VDR
QG021-01	G10	HQP018475	NM_001025366	VEGFA

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG021-01	G11	HQP005427	NM_015850	FGFR1
QG021-01	G12	HQP012204	NM_000438	PAX3
QG021-01	H01	HGDC		
QG021-01	H02	HGDC		
QG021-01	H03	HQP006940	NM_002046	GAPDH
QG021-01	H04	HQP016381	NM_001101	ACTB
QG021-01	H05	HQP015171	NM_004048	B2M
QG021-01	H06	HQP006171	NM_012423	RPL13A
QG021-01	H07	HQP009026	NM_000194	HPRT1
QG021-01	H08	HQP054253	NR_003286	RN18S1
QG021-01	H09	RT		
QG021-01	H10	RT		
QG021-01	H11	PCR		
QG021-01	H12	PCR		