

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG037-01	A01	HQP004202	NM_213566	DFFA
QG037-01	A02	HQP020250	NM_032983	CASP2
QG037-01	A03	HQP009026	NM_000194	HPRT1
QG037-01	A04	HQP022754	NM_152942	TNFRSF8
QG037-01	A05	HQP018234	NM_003300	TRAF3
QG037-01	A06	HQP018233	NM_021138	TRAF2
QG037-01	A07	HQP018232	NM_005658	TRAF1
QG037-01	A08	HQP022769	NM_001244	TNFSF8
QG037-01	A09	HQP023108	NM_001252	CD70
QG037-01	A10	HQP015306	NM_020232	TNFSF5IP1
QG037-01	A11	HQP018329	NM_003326	TNFSF4
QG037-01	A12	HQP021834	NM_005092	TNFSF18
QG037-01	B01	HQP021496	NM_003807	TNFSF14
QG037-01	B02	HQP000821	NM_006573	TNFSF13B
QG037-01	B03	HQP021501	NM_003809	TNFSF12
QG037-01	B04	HQP021502	NM_003810	TNFSF10
QG037-01	B05	HQP009716	NM_001561	TNFRSF9
QG037-01	B06	HQP022753	NM_001243	TNFRSF8
QG037-01	B07	HQP022667	NM_001242	CD27
QG037-01	B08	HQP007613	NM_014452	TNFRSF21
QG037-01	B09	HQP018149	NM_001066	TNFRSF1B
QG037-01	B10	HQP018148	NM_001065	TNFRSF1A
QG037-01	B11	HQP021069	NM_032871	TNFRSF19L
QG037-01	B12	HQP014424	NM_018647	TNFRSF19
QG037-01	C01	HQP021536	NM_004195	TNFRSF18
QG037-01	C02	HQP016367	NM_001192	TNFRSF17
QG037-01	C03	HQP021522	NM_003820	TNFRSF14
QG037-01	C04	HQP006148	NM_012452	TNFRSF13B
QG037-01	C05	HQP012651	NM_016639	TNFRSF12A
QG037-01	C06	HQP012049	NM_002546	TNFRSF11B
QG037-01	C07	HQP021550	NM_003839	TNFRSF11A
QG037-01	C08	HQP021551	NM_003840	TNFRSF10D
QG037-01	C09	HQP021552	NM_003841	TNFRSF10C
QG037-01	C10	HQP021553	NM_003842	TNFRSF10B
QG037-01	C11	HQP021557	NM_003844	TNFRSF10A
QG037-01	C12	HQP018145	NM_006290	TNFAIP3
QG037-01	D01	HQP018141	NM_000594	TNF
QG037-01	D02	HQP017688	NM_003127	SPTAN1
QG037-01	D03	HQP016131	NM_000321	RB1
QG037-01	D04	HQP003120	NM_001618	PARP1
QG037-01	D05	HQP012170	NM_002577	PAK2

QG037-01	D06	HQP012156	NM_002576	PAK1
QG037-01	D07	HQP011828	NM_002507	NGFR
QG037-01	D08	HQP011810	NM_020529	NFKBIA
QG037-01	D09	HQP011807	NM_003998	NFKB1
QG037-01	D10	HQP014886	NM_002750	MAPK8
QG037-01	D11	HQP017891	NM_003188	MAP3K7
QG037-01	D12	HQP053966	NM_005921	MAP3K1
QG037-01	E01	HQP016830	NM_003010	MAP2K4
QG037-01	E02	HQP021290	NM_003682	MADD
QG037-01	E03	HQP010915	NM_002342	LTBR
QG037-01	E04	HQP010907	NM_000595	LTA
QG037-01	E05	HQP020961	NM_032737	LMNB2
QG037-01	E06	HQP010680	NM_005573	LMNB1
QG037-01	E07	HQP009853	NM_002228	JUN
QG037-01	E08	HQP021140	NM_003639	IKBKG
QG037-01	E09	HQP009639	NM_001556	IKBKB
QG037-01	E10	HQP009042	NM_004504	HRB
QG037-01	E11	HQP009671	NM_000639	FASLG
QG037-01	E12	HQP009651	NM_000043	FAS
QG037-01	F01	HQP021526	NM_003824	FADD
QG037-01	F02	HQP016300	NM_021783	EDA2R
QG037-01	F03	HQP004498	NM_004417	DUSP1
QG037-01	F04	HQP004201	NM_004401	DFFA
QG037-01	F05	HQP021493	NM_003805	CRADD
QG037-01	F06	HQP001708	NM_001278	CHUK
QG037-01	F07	HQP022962	NM_000074	CD40LG
QG037-01	F08	HQP022955	NM_001250	CD40
QG037-01	F09	HQP020297	NM_004346	CASP3
QG037-01	F10	HQP020249	NM_032982	CASP2
QG037-01	F11	HQP018928	NM_004341	CAD
QG037-01	F12	HQP022871	NM_004874	BAG4
QG037-01	G01	HQP010610	NM_001175	ARHGDIB
QG037-01	G02	HQP020548	NM_001228	CASP8
QG037-01	G03	HQP021525	NM_003823	TNFRSF6B
QG037-01	G04	HQP023417	NM_005118	TNFSF15
QG037-01	G05	HQP021469	NM_148970	TNFRSF25
QG037-01	G06	HQP021141	NM_003640	IKBKAP
QG037-01	G07	HQP010673	NM_005572	LMNA
QG037-01	G08	HQP010908	NM_002341	LTB
QG037-01	G09	HQP022956	NM_152854	CD40
QG037-01	G10	HQP021321	NM_003701	TNFSF11
QG037-01	G11	HQP021293	NM_130472	MADD
QG037-01	G12	HQP021465	NM_003790	TNFRSF25

QG037-01	H01	HGDC		
QG037-01	H02	HGDC		
QG037-01	H03	HQP006940	NM_002046	GAPDH
QG037-01	H04	HQP016381	NM_001101	ACTB
QG037-01	H05	HQP015171	NM_004048	B2M
QG037-01	H06	HQP006171	NM_012423	RPL13A
QG037-01	H07	HQP009026	NM_000194	HPRT1
QG037-01	H08	HQP054253	NR_003286	RN18S1
QG037-01	H09	RT		
QG037-01	H10	RT		
QG037-01	H11	PCR		
QG037-01	H12	PCR		