

Plate	Position	Catalog No.of Primer	Accession No.of Gene	Symbol
QG099-01	A01	HQP004991	NM_001014431	AKT1
QG099-01	A02	HQP008934	NM_001160	APAF1
QG099-01	A03	HQP021206	NM_006595	API5
QG099-01	A04	HQP015389	NM_020371	AVEN
QG099-01	A05	HQP015538	NM_004322	BAD
QG099-01	A06	HQP015574	NM_004323	BAG1
QG099-01	A07	HQP022872	NM_004281	BAG3
QG099-01	A08	HQP022871	NM_004874	BAG4
QG099-01	A09	HQP015917	NM_001188	BAK1
QG099-01	A10	HQP015964	NM_004324	BAX
QG099-01	A11	HQP007522	NM_014417	BBC3
QG099-01	A12	HQP021725	NM_003921	BCL10
QG099-01	B01	HQP016211	NM_000633	BCL2
QG099-01	B02	HQP016222	NM_004049	BCL2A1
QG099-01	B03	HQP016237	NM_001191	BCL2L1
QG099-01	B04	HQP000028	NM_006538	BCL2L11
QG099-01	B05	HQP019021	NM_030766	BCL2L14
QG099-01	B06	HQP016260	NM_004050	BCL2L2
QG099-01	B07	HQP016328	NM_001706	BCL6
QG099-01	B08	HQP012590	NM_016561	BFAR
QG099-01	B09	HQP016683	NM_001197	BIK
QG099-01	B10	HQP009072	NM_001166	BIRC2
QG099-01	B11	HQP009084	NM_001165	BIRC3
QG099-01	B12	HQP009091	NM_001167	BIRC4
QG099-01	C01	HQP009099	NM_001012270	BIRC5
QG099-01	C02	HQP017593	NM_001205	BNIP1
QG099-01	C03	HQP017605	NM_004330	BNIP2
QG099-01	C04	HQP017619	NM_004052	BNIP3
QG099-01	C05	HQP017631	NM_004331	BNIP3L
QG099-01	C06	HQP017642	NM_032515	BOK
QG099-01	C07	HQP008522	NM_014550	CARD10
QG099-01	C08	HQP020249	NM_032982	CASP2
QG099-01	C09	HQP020297	NM_004346	CASP3

QG099-01	C10	HQP020480	NM_001227	CASP7
QG099-01	C11	HQP018966	NM_001080124	CASP8
QG099-01	C12	HQP020648	NM_001229	CASP9
QG099-01	D01	HQP022955	NM_001250	CD40
QG099-01	D02	HQP002037	NM_001831	CLU
QG099-01	D03	HQP003943	NM_001344	DAD1
QG099-01	D04	HQP003975	NM_004394	DAP
QG099-01	D05	HQP003979	NM_004938	DAPK1
QG099-01	D06	HQP006265	NM_014326	DAPK2
QG099-01	D07	HQP004127	NM_004083	DDIT3
QG099-01	D08	HQP004201	NM_004401	DFFA
QG099-01	D09	HQP015116	NM_019887	DIABLO
QG099-01	D10	HQP016218	NM_006268	DPF2
QG099-01	D11	HQP004524	NM_005225	E2F1
QG099-01	D12	HQP004526	NM_004091	E2F2
QG099-01	E01	HQP021526	NM_003824	FADD
QG099-01	E02	HQP009651	NM_000043	FAS
QG099-01	E03	HQP009671	NM_000639	FASLG
QG099-01	E04	HQP005434	NM_000142	FGFR3
QG099-01	E05	HQP005759	NM_001455	FOXO3A
QG099-01	E06	HQP001042	NM_006705	GADD45G
QG099-01	E07	HQP008279	NM_000581	GPX1
QG099-01	E08	HQP008469	NM_002093	GSK3B
QG099-01	E09	HQP008831	NM_001530	HIF1A
QG099-01	E10	HQP008358	NM_022740	HIPK2
QG099-01	E11	HQP021494	NM_003806	HRK
QG099-01	E12	HQP009077	NM_005345	HSPA1A
QG099-01	F01	HQP007710	NM_013247	HTRA2
QG099-01	F02	HQP009518	NM_000618	IGF1
QG099-01	F03	HQP010907	NM_000595	LTA
QG099-01	F04	HQP010915	NM_002342	LTBR
QG099-01	F05	HQP011104	NM_021960	MCL1
QG099-01	F06	HQP011670	NM_004536	NAIP
QG099-01	F07	HQP011807	NM_003998	NFKB1
QG099-01	F08	HQP012197	NM_002583	PAWR

QG099-01	F09	HQP016753	NM_022121	PERP
QG099-01	F10	HQP018000	NM_005809	PRDX2
QG099-01	F11	HQP015535	NM_000314	PTEN
QG099-01	F12	HQP016074	NM_006265	RAD21
QG099-01	G01	HQP021492	NM_003804	RIPK1
QG099-01	G02	HQP006080	NM_012238	SIRT1
QG099-01	G03	HQP017615	NM_000454	SOD1
QG099-01	G04	HQP017771	NM_003152	STAT5A
QG099-01	G05	HQP017774	NM_012448	STAT5B
QG099-01	G06	HQP018044	NM_000660	TGFB1
QG099-01	G07	HQP018141	NM_000594	TNF
QG099-01	G08	HQP018145	NM_006290	TNFAIP3
QG099-01	G09	HQP021557	NM_003844	TNFRSF10A
QG099-01	G10	HQP012049	NM_002546	TNFRSF11B
QG099-01	G11	HQP018148	NM_001065	TNFRSF1A
QG099-01	G12	HQP018149	NM_001066	TNFRSF1B
QG099-01	H01	HGDC		
QG099-01	H02	HGDC		
QG099-01	H03	HQP006940	NM_002046	GAPDH
QG099-01	H04	HQP016381	NM_001101	ACTB
QG099-01	H05	HQP015171	NM_004048	B2M
QG099-01	H06	HQP006171	NM_012423	RPL13A
QG099-01	H07	HQP009026	NM_000194	HPRT1
QG099-01	H08	HQP054253	NR_003286	RN18S1
QG099-01	H09	RT		
QG099-01	H10	RT		
QG099-01	H11	PCR		
QG099-01	H12	PCR		