

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG051-01	A01	HQP002912	NM_001880	ATF2
QG051-01	A02	HQP015964	NM_004324	BAX
QG051-01	A03	HQP016211	NM_000633	BCL2
QG051-01	A04	HQP016222	NM_004049	BCL2A1
QG051-01	A05	HQP016238	NM_138578	BCL2L1
QG051-01	A06	HQP011670	NM_004536	NAIP
QG051-01	A07	HQP009072	NM_001166	BIRC2
QG051-01	A08	HQP009084	NM_001165	BIRC3
QG051-01	A09	HQP017333	NM_001200	BMP2
QG051-01	A10	HQP053910	NM_130851	BMP4
QG051-01	A11	HQP017713	NM_007294	BRCA1
QG051-01	A12	HQP016621	NM_002982	CCL2
QG051-01	B01	HQP016639	NM_004591	CCL20
QG051-01	B02	HQP016204	NM_053056	CCND1
QG051-01	B03	HQP022337	NM_014207	CD5
QG051-01	B04	HQP000225	NM_001798	CDK2
QG051-01	B05	HQP000331	NM_000389	CDKN1A
QG051-01	B06	HQP000342	NM_004064	CDKN1B
QG051-01	B07	HQP000369	NM_000077	CDKN2A
QG051-01	B08	HQP003159	NM_000758	CSF2
QG051-01	B09	HQP011220	NM_002416	CXCL9
QG051-01	B10	HQP003904	NM_000103	CYP19A1
QG051-01	B11	HQP004612	NM_001964	EGR1
QG051-01	B12	HQP009651	NM_000043	FAS
QG051-01	C01	HQP009671	NM_000639	FASLG
QG051-01	C02	HQP005134	NM_004104	FASN
QG051-01	C03	HQP006022	NM_002026	FN1
QG051-01	C04	HQP006188	NM_005252	FOS
QG051-01	C05	HQP008906	NM_021784	FOXA2
QG051-01	C06	HQP004125	NM_001924	GADD45A
QG051-01	C07	HQP023078	NM_014668	GREB1
QG051-01	C08	HQP008653	NM_002103	GYS1
QG051-01	C09	HQP008843	NM_000189	HK2
QG051-01	C10	HQP008966	NM_005522	HOXA1
QG051-01	C11	HQP009068	NM_005526	HSF1
QG051-01	C12	HQP009089	NM_001540	HSPB1
QG051-01	D01	HQP009184	NM_000201	ICAM1
QG051-01	D02	HQP009544	NM_000598	IGFBP3
QG051-01	D03	HQP009639	NM_001556	IKBKB
QG051-01	D04	HQP009640	NM_000575	IL1A
QG051-01	D05	HQP009649	NM_000586	IL2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG051-01	D06	HQP009662	NM_000589	IL4
QG051-01	D07	HQP009664	NM_000418	IL4R
QG051-01	D08	HQP009678	NM_000584	IL8
QG051-01	D09	HQP009778	NM_002198	IRF1
QG051-01	D10	HQP009853	NM_002228	JUN
QG051-01	D11	HQP012480	NM_016269	LEF1
QG051-01	D12	HQP010581	NM_000230	LEP
QG051-01	E01	HQP010907	NM_000595	LTA
QG051-01	E02	HQP011135	NM_002392	MDM2
QG051-01	E03	HQP011264	NM_002425	MMP10
QG051-01	E04	HQP011258	NM_002423	MMP7
QG051-01	E05	HQP011597	NM_002467	MYC
QG051-01	E06	HQP011807	NM_003998	NFKB1
QG051-01	E07	HQP011866	NM_000625	NOS2A
QG051-01	E08	HQP020042	NM_003489	NRIP1
QG051-01	E09	HQP012013	NM_002539	ODC1
QG051-01	E10	HQP013015	NM_000442	PECAM1
QG051-01	E11	HQP013634	NM_015869	PPARG
QG051-01	E12	HQP014706	NM_002737	PRKCA
QG051-01	F01	HQP015530	NM_000264	PTCH1
QG051-01	F02	HQP015598	NM_000963	PTGS2
QG051-01	F03	HQP016190	NM_002899	RBP1
QG051-01	F04	HQP016744	NM_000450	SELE
QG051-01	F05	HQP016747	NM_003006	SELPLG
QG051-01	F06	HQP000019	NM_004180	TANK
QG051-01	F07	HQP017958	NM_003202	TCF7
QG051-01	F08	HQP018041	NM_003234	TFRC
QG051-01	F09	HQP015258	NM_020182	TMEPAI
QG051-01	F10	HQP018141	NM_000594	TNF
QG051-01	F11	HQP018175	NM_000546	TP53
QG051-01	F12	HQP022884	NM_004881	TP53I3
QG051-01	G01	HQP053975	NM_001078	VCAM1
QG051-01	G02	HQP018481	NM_003376	VEGFA
QG051-01	G03	HQP021610	NM_003882	WISP1
QG051-01	G04	HQP018528	NM_005430	WNT1
QG051-01	G05	HQP018529	NM_003391	WNT2
QG051-01	G06	HQP009026	NM_000194	HPRT1
QG051-01	G07	HQP000020	NM_133484	TANK
QG051-01	G08	HQP017959	NM_201632	TCF7
QG051-01	G09	HQP011136	NM_006878	MDM2
QG051-01	G10	HQP000623	NM_005194	CEBPB

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG051-01	G11	HQP004846	NM_001426	EN1
QG051-01	G12	HQP014742	NM_005400	PRKCE
QG051-01	H01	HGDC		
QG051-01	H02	HGDC		
QG051-01	H03	HQP006940	NM_002046	GAPDH
QG051-01	H04	HQP016381	NM_001101	ACTB
QG051-01	H05	HQP015171	NM_004048	B2M
QG051-01	H06	HQP006171	NM_012423	RPL13A
QG051-01	H07	HQP009026	NM_000194	HPRT1
QG051-01	H08	HQP054253	NR_003286	RN18S1
QG051-01	H09	RT		
QG051-01	H10	RT		
QG051-01	H11	PCR		
QG051-01	H12	PCR		