

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG050-01	A01	HQP011136	NM_006878	MDM2
QG050-01	A02	HQP020550	NM_033358	CASP8
QG050-01	A03	HQP020210	NM_033294	CASP1
QG050-01	A04	HQP009026	NM_000194	HPRT1
QG050-01	A05	HQP018563	NM_005431	XRCC2
QG050-01	A06	HQP018562	NM_006297	XRCC1
QG050-01	A07	HQP013614	NM_007120	UGT1A4
QG050-01	A08	HQP018175	NM_000546	TP53
QG050-01	A09	HQP021502	NM_003810	TNFSF10
QG050-01	A10	HQP018148	NM_001065	TNFRSF1A
QG050-01	A11	HQP018141	NM_000594	TNF
QG050-01	A12	HQP017616	NM_000636	SOD2
QG050-01	B01	HQP017615	NM_000454	SOD1
QG050-01	B02	HQP012154	NM_000602	SERPINE1
QG050-01	B03	HQP000145	NM_005732	RAD50
QG050-01	B04	HQP016075	NM_005053	RAD23A
QG050-01	B05	HQP015596	NM_000962	PTGS1
QG050-01	B06	HQP018000	NM_005809	PRDX2
QG050-01	B07	HQP012152	NM_002574	PRDX1
QG050-01	B08	HQP013504	NM_000941	POR
QG050-01	B09	HQP054038	NM_182649	PCNA
QG050-01	B10	HQP011866	NM_000625	NOS2A
QG050-01	B11	HQP011810	NM_020529	NFKBIA
QG050-01	B12	HQP011807	NM_003998	NFKB1
QG050-01	C01	HQP011538	NM_005953	MT2A
QG050-01	C02	HQP011219	NM_002415	MIF
QG050-01	C03	HQP011135	NM_002392	MDM2
QG050-01	C04	HQP010907	NM_000595	LTA
QG050-01	C05	HQP009670	NM_000600	IL6
QG050-01	C06	HQP009641	NM_000576	IL1B
QG050-01	C07	HQP009640	NM_000575	IL1A
QG050-01	C08	HQP009718	NM_001562	IL18
QG050-01	C09	HQP009555	NM_002178	IGFBP6
QG050-01	C10	HQP000946	NM_006644	HSPH1
QG050-01	C11	HQP009102	NM_002157	HSPE1
QG050-01	C12	HQP009098	NM_002156	HSPD1
QG050-01	D01	HQP009097	NM_007355	HSP90AB1
QG050-01	D02	HQP009089	NM_001540	HSPB1
QG050-01	D03	HQP009086	NM_006597	HSPA8
QG050-01	D04	HQP009085	NM_002155	HSPA6
QG050-01	D05	HQP009083	NM_005347	HSPA5

QG050-01	D06	HQP009081	NM_002154	HSPA4
QG050-01	D07	HQP009080	NM_021979	HSPA2
QG050-01	D08	HQP009079	NM_005527	HSPA1L
QG050-01	D09	HQP009077	NM_005345	HSPA1A
QG050-01	D10	HQP009068	NM_005526	HSF1
QG050-01	D11	HQP008898	NM_002133	HMOX1
QG050-01	D12	HQP008483	NM_000849	GSTM3
QG050-01	E01	HQP008473	NM_000637	GSR
QG050-01	E02	HQP008279	NM_000581	GPX1
QG050-01	E03	HQP022853	NM_004864	GDF15
QG050-01	E04	HQP004125	NM_001924	GADD45A
QG050-01	E05	HQP005979	NM_001461	FMO5
QG050-01	E06	HQP005939	NM_002021	FMO1
QG050-01	E07	HQP009671	NM_000639	FASLG
QG050-01	E08	HQP004983	NM_000122	ERCC3
QG050-01	E09	HQP004974	NM_001983	ERCC1
QG050-01	E10	HQP004950	NM_001979	EPHX2
QG050-01	E11	HQP004612	NM_001964	EGR1
QG050-01	E12	HQP004524	NM_005225	E2F1
QG050-01	F01	HQP001241	NM_007034	DNAJB4
QG050-01	F02	HQP009076	NM_001539	DNAJA1
QG050-01	F03	HQP004127	NM_004083	DDIT3
QG050-01	F04	HQP004111	NM_001923	DDB1
QG050-01	F05	HQP003859	NM_000780	CYP7A1
QG050-01	F06	HQP003817	NM_000773	CYP2E1
QG050-01	F07	HQP003772	NM_000499	CYP1A1
QG050-01	F08	HQP009746	NM_001565	CXCL10
QG050-01	F09	HQP003159	NM_000758	CSF2
QG050-01	F10	HQP000331	NM_000389	CDKN1A
QG050-01	F11	HQP021857	NM_004060	CCNG1
QG050-01	F12	HQP016204	NM_053056	CCND1
QG050-01	G01	HQP016625	NM_002984	CCL4
QG050-01	G02	HQP016622	NM_002983	CCL3
QG050-01	G03	HQP016640	NM_002989	CCL21
QG050-01	G04	HQP020709	NM_001230	CASP10
QG050-01	G05	HQP020208	NM_033292	CASP1
QG050-01	G06	HQP016238	NM_138578	BCL2L1
QG050-01	G07	HQP015964	NM_004324	BAX
QG050-01	G08	HQP011736	NM_000051	ATM
QG050-01	G09	HQP008829	NM_001154	ANXA5
QG050-01	G10	HQP020548	NM_001228	CASP8
QG050-01	G11	HQP020946	NM_001752	CAT
QG050-01	G12	HQP021735	NM_005190	CCNC

QG050-01	H01	HGDC		
QG050-01	H02	HGDC		
QG050-01	H03	HQP006940	NM_002046	GAPDH
QG050-01	H04	HQP016381	NM_001101	ACTB
QG050-01	H05	HQP015171	NM_004048	B2M
QG050-01	H06	HQP006171	NM_012423	RPL13A
QG050-01	H07	HQP009026	NM_000194	HPRT1
QG050-01	H08	HQP054253	NR_003286	RN18S1
QG050-01	H09	RT		
QG050-01	H10	RT		
QG050-01	H11	PCR		
QG050-01	H12	PCR		