

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-01	A01	HQP004605	NM_005228	EGFR
QG071-01	A02	HQP018175	NM_000546	TP53
QG071-01	A03	HQP003772	NM_000499	CYP1A1
QG071-01	A04	HQP010133	NM_004985	KRAS
QG071-01	A05	HQP018562	NM_006297	XRCC1
QG071-01	A06	HQP004976	NM_000400	ERCC2
QG071-01	A07	HQP011309	NM_000250	MPO
QG071-01	A08	HQP004975	NM_202001	ERCC1
QG071-01	A09	HQP012021	NM_002542	OGG1
QG071-01	A10	HQP004948	NM_000120	EPHX1
QG071-01	A11	HQP004317	NM_000903	NQO1
QG071-01	A12	HQP001136	NM_000015	NAT2
QG071-01	B01	HQP011547	NM_005957	MTHFR
QG071-01	B02	HQP003817	NM_000773	CYP2E1
QG071-01	B03	HQP001643	NM_000745	CHRNA5
QG071-01	B04	HQP018556	NM_004628	XPC
QG071-01	B05	HQP011135	NM_002392	MDM2
QG071-01	B06	HQP001632	NM_000743	CHRNA3
QG071-01	B07	HQP018141	NM_000594	TNF
QG071-01	B08	HQP009061	NM_001641	APEX1
QG071-01	B09	HQP009641	NM_000576	IL1B
QG071-01	B10	HQP011255	NM_002421	MMP1
QG071-01	B11	HQP018564	NM_005432	XRCC3
QG071-01	B12	HQP015598	NM_000963	PTGS2
QG071-01	C01	HQP013100	NM_000927	ABCB1
QG071-01	C02	HQP009670	NM_000600	IL6
QG071-01	C03	HQP003775	NM_000104	CYP1B1
QG071-01	C04	HQP017616	NM_000636	SOD2
QG071-01	C05	HQP004985	NM_000123	ERCC5
QG071-01	C06	HQP008483	NM_000849	GSTM3
QG071-01	C07	HQP018475	NM_001025366	VEGFA
QG071-01	C08	HQP009678	NM_000584	IL8
QG071-01	C09	HQP018044	NM_000660	TGFB1
QG071-01	C10	HQP009685	NM_000572	IL10
QG071-01	C11	HQP017794	NM_000455	STK11
QG071-01	C12	HQP016204	NM_053056	CCND1
QG071-01	D01	HQP009671	NM_000639	FASLG
QG071-01	D02	HQP003841	NM_000777	CYP3A5
QG071-01	D03	HQP019829	NM_030782	CLPTM1L
QG071-01	D04	HQP011257	NM_002422	MMP3
QG071-01	D05	HQP004984	NM_005236	ERCC4

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-01	D06	HQP003814	NM_000106	CYP2D6
QG071-01	D07	HQP003809	NM_000769	CYP2C19
QG071-01	D08	HQP018180	NM_005427	TP73
QG071-01	D09	HQP018563	NM_005431	XRCC2
QG071-01	D10	HQP011868	NM_000603	NOS3
QG071-01	D11	HQP011687	NM_002485	NBN
QG071-01	D12	HQP011235	NM_000249	MLH1
QG071-01	E01	HQP004986	NM_000124	ERCC6
QG071-01	E02	HQP004446	NM_000795	DRD2
QG071-01	E03	HQP018342	NM_001071	TYMS
QG071-01	E04	HQP017811	NM_177536	SULT1A1
QG071-01	E05	HQP017753	NM_000059	BRCA2
QG071-01	E06	HQP013633	NM_005037	PPARG
QG071-01	E07	HQP011263	NM_004994	MMP9
QG071-01	E08	HQP011256	NM_004530	MMP2
QG071-01	E09	HQP010609	NM_000234	LIG1
QG071-01	E10	HQP009697	NM_002188	IL13
QG071-01	E11	HQP009645	NM_000577	IL1RN
QG071-01	E12	HQP009640	NM_000575	IL1A
QG071-01	F01	HQP003120	NM_001618	PARP1
QG071-01	F02	HQP002671	NM_000754	COMT
QG071-01	F03	HQP054030	NM_130398	EXO1
QG071-01	F04	HQP023467	NM_000662	NAT1
QG071-01	F05	HQP018983	NM_004639	BAT3
QG071-01	F06	HQP018966	NM_001080124	CASP8
QG071-01	F07	HQP018565	NM_003401	XRCC4
QG071-01	F08	HQP017733	NM_004333	BRAF
QG071-01	F09	HQP017482	NM_194255	SLC19A1
QG071-01	F10	HQP015535	NM_000314	PTEN
QG071-01	F11	HQP015192	NM_002789	PSMA4
QG071-01	F12	HQP011555	NM_002454	MTRR
QG071-01	G01	HQP011491	NM_000251	MSH2
QG071-01	G02	HQP011266	NM_002426	MMP12
QG071-01	G03	HQP011181	NM_000245	MET
QG071-01	G04	HQP009662	NM_000589	IL4
QG071-01	G05	HQP009544	NM_000598	IGFBP3
QG071-01	G06	HQP008849	NM_002116	HLA-A
QG071-01	G07	HQP005087	NM_000690	ALDH2
QG071-01	G08	HQP004658	NM_001621	AHR
QG071-01	G09	HQP003904	NM_000103	CYP19A1
QG071-01	G10	HQP003888	NM_000102	CYP17A1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-01	G11	HQP001667	NM_000750	CHRNA4
QG071-01	G12	HQP001396	NM_001005735	CHEK2
QG071-01	H01	HGDC		
QG071-01	H02	HGDC		
QG071-01	H03	HQP006940	NM_002046	GAPDH
QG071-01	H04	HQP016381	NM_001101	ACTB
QG071-01	H05	HQP015171	NM_004048	B2M
QG071-01	H06	HQP006171	NM_012423	RPL13A
QG071-01	H07	HQP009026	NM_000194	HPRT1
QG071-01	H08	HQP054253	NR_003286	RN18S1
QG071-01	H09	RT		
QG071-01	H10	RT		
QG071-01	H11	PCR		
QG071-01	H12	PCR		
QG071-02	A01	HQP000370	NM_058195	CDKN2A
QG071-02	A02	HQP054057	BC071181	TGFB1
QG071-02	A03	HQP021906	NM_001239	CCNH
QG071-02	A04	HQP021736	NM_003925	MBD4
QG071-02	A05	HQP021603	NM_003878	GGH
QG071-02	A06	HQP021557	NM_003844	TNFRSF10A
QG071-02	A07	HQP021518	NM_000071	CBS
QG071-02	A08	HQP021385	NM_003739	AKR1C3
QG071-02	A09	HQP020297	NM_004346	CASP3
QG071-02	A10	HQP018474	NM_000376	VDR
QG071-02	A11	HQP016076	NM_002874	RAD23B
QG071-02	A12	HQP013122	NM_000295	SERPINA1
QG071-02	B01	HQP011554	NM_000254	MTR
QG071-02	B02	HQP010612	NM_013975	LIG3
QG071-02	B03	HQP009693	NM_002187	IL12B
QG071-02	B04	HQP009692	NM_000882	IL12A
QG071-02	B05	HQP009679	NM_000634	IL8RA
QG071-02	B06	HQP009672	NM_000565	IL6R
QG071-02	B07	HQP009664	NM_000418	IL4R
QG071-02	B08	HQP009649	NM_000586	IL2
QG071-02	B09	HQP009460	NM_000874	IFNAR2
QG071-02	B10	HQP008479	NM_001512	GSTA4
QG071-02	B11	HQP004991	NM_001014431	AKT1
QG071-02	B12	HQP004081	NM_000789	ACE
QG071-02	C01	HQP003811	NM_000771	CYP2C9
QG071-02	C02	HQP002215	NM_001013619	LOC123688
QG071-02	C03	HQP011601	NM_005376	MYCL1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-02	C04	HQP054684	NM_001033	RRM1
QG071-02	C05	HQP054047	BC008403	HLA-DRB1
QG071-02	C06	HQP023203	NM_001785	CDA
QG071-02	C07	HQP018568	NM_021141	XRCC5
QG071-02	C08	HQP016418	NM_004327	BCR
QG071-02	C09	HQP013422	NM_002690	POLB
QG071-02	C10	HQP012420	NM_002592	PCNA
QG071-02	C11	HQP010613	NM_002312	LIG4
QG071-02	C12	HQP009472	NM_005534	IFNGR2
QG071-02	D01	HQP009064	NM_000413	HSD17B1
QG071-02	D02	HQP007304	NM_012419	RGS17
QG071-02	D03	HQP005440	NM_022963	FGFR4
QG071-02	D04	HQP004998	NM_000125	ESR1
QG071-02	D05	HQP004450	NM_000797	DRD4
QG071-02	D06	HQP003804	NM_000766	CYP2A13
QG071-02	D07	HQP003159	NM_000758	CSF2
QG071-02	D08	HQP001374	NM_007182	RASSF1
QG071-02	D09	HQP017618	NM_003102	SOD3
QG071-02	D10	HQP054045	BC004257	RET
QG071-02	D11	HQP023466	NM_004360	CDH1
QG071-02	D12	HQP023409	NM_001790	CDC25C
QG071-02	E01	HQP023003	NM_004296	RGS6
QG071-02	E02	HQP022884	NM_004881	TP53I3
QG071-02	E03	HQP021405	NM_003747	TNKS
QG071-02	E04	HQP020481	NM_033338	CASP7
QG071-02	E05	HQP018176	NM_005657	TP53BP1
QG071-02	E06	HQP018016	NM_005652	TERF2
QG071-02	E07	HQP018014	NM_003218	TERF1
QG071-02	E08	HQP016958	NM_000542	SFTPFB
QG071-02	E09	HQP016801	NM_022162	NOD2
QG071-02	E10	HQP016634	NM_001713	BHMT
QG071-02	E11	HQP016621	NM_002982	CCL2
QG071-02	E12	HQP016087	NM_134424	RAD52
QG071-02	F01	HQP015946	NM_000465	BARD1
QG071-02	F02	HQP013596	NM_001184	ATR
QG071-02	F03	HQP013152	NM_002648	PIM1
QG071-02	F04	HQP013150	NM_006218	PIK3CA
QG071-02	F05	HQP012755	NM_001037872	REV1
QG071-02	F06	HQP011866	NM_000625	NOS2A
QG071-02	F07	HQP011597	NM_002467	MYC
QG071-02	F08	HQP011325	NM_019899	ABCC1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-02	F09	HQP010868	NM_006152	LRMP
QG071-02	F10	HQP010581	NM_000230	LEP
QG071-02	F11	HQP009710	NM_002189	IL15RA
QG071-02	F12	HQP009708	NM_000585	IL15
QG071-02	G01	HQP009677	NM_002185	IL7R
QG071-02	G02	HQP009523	NM_000875	IGF1R
QG071-02	G03	HQP009518	NM_000618	IGF1
QG071-02	G04	HQP009059	NM_000862	HSD3B1
QG071-02	G05	HQP009024	NM_000038	APC
QG071-02	G06	HQP008493	NM_000179	MSH6
QG071-02	G07	HQP008490	NM_001513	GSTZ1
QG071-02	G08	HQP008476	NM_000846	GSTA2
QG071-02	G09	HQP008282	NM_002084	GPX3
QG071-02	G10	HQP008281	NM_002083	GPX2
QG071-02	G11	HQP006855	NM_001042594	POT1
QG071-02	G12	HQP005447	NM_002012	FHIT
QG071-02	H01	HGDC		
QG071-02	H02	HGDC		
QG071-02	H03	HQP006940	NM_002046	GAPDH
QG071-02	H04	HQP016381	NM_001101	ACTB
QG071-02	H05	HQP015171	NM_004048	B2M
QG071-02	H06	HQP006171	NM_012423	RPL13A
QG071-02	H07	HQP009026	NM_000194	HPRT1
QG071-02	H08	HQP054253	NR_003286	RN18S1
QG071-02	H09	RT		
QG071-02	H10	RT		
QG071-02	H11	PCR		
QG071-02	H12	PCR		
QG071-03	A01	HQP004983	NM_000122	ERCC3
QG071-03	A02	HQP004689	NM_001972	ELA2
QG071-03	A03	HQP004599	NM_001963	EGF
QG071-03	A04	HQP004309	NM_000791	DHFR
QG071-03	A05	HQP003871	NM_000781	CYP11A1
QG071-03	A06	HQP003499	NM_001037631	CTLA4
QG071-03	A07	HQP003173	NM_000759	CSF3
QG071-03	A08	HQP002590	NM_001085	SERPINA3
QG071-03	A09	HQP002413	NM_000669	ADH1C
QG071-03	A10	HQP001399	NM_007195	POLI
QG071-03	A11	HQP001282	NM_001274	CHEK1
QG071-03	A12	HQP000963	NM_012190	ALDH1L1
QG071-03	B01	HQP000261	NM_004935	CDK5

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-03	B02	HQP004972	NM_001042599	ERBB4
QG071-03	B03	HQP004414	NM_006892	DNMT3B
QG071-03	B04	HQP000320	NM_005845	ABCC4
QG071-03	B05	HQP053960	NM_001042	SLC2A4
QG071-03	B06	HQP053959	NM_001040	SHBG
QG071-03	B07	HQP023437	NM_005123	NR1H4
QG071-03	B08	HQP023414	NM_005847	SLC23A1
QG071-03	B09	HQP023399	NM_004358	CDC25B
QG071-03	B10	HQP023385	NM_001789	CDC25A
QG071-03	B11	HQP023240	NM_014781	RB1CC1
QG071-03	B12	HQP023195	NM_018834	MATR3
QG071-03	C01	HQP023168	NM_004356	CD81
QG071-03	C02	HQP022955	NM_001250	CD40
QG071-03	C03	HQP022863	NM_004870	MPDU1
QG071-03	C04	HQP022853	NM_004864	GDF15
QG071-03	C05	HQP022851	NM_004862	LITAF
QG071-03	C06	HQP022833	NM_005505	SCARB1
QG071-03	C07	HQP022747	NM_175862	CD86
QG071-03	C08	HQP022723	NM_004820	CYP7B1
QG071-03	C09	HQP022722	NM_005191	CD80
QG071-03	C10	HQP022580	NM_005454	CER1
QG071-03	C11	HQP022316	NM_000616	CD4
QG071-03	C12	HQP022202	NM_004213	SLC28A1
QG071-03	D01	HQP022170	NM_004708	PDCD5
QG071-03	D02	HQP021737	NM_003927	MBD2
QG071-03	D03	HQP021508	NM_001236	CBR3
QG071-03	D04	HQP021495	NM_001757	CBR1
QG071-03	D05	HQP021459	NM_003786	ABCC3
QG071-03	D06	HQP021388	NM_003742	ABCB11
QG071-03	D07	HQP021313	NM_001753	CAV1
QG071-03	D08	HQP021023	NM_003617	RGS5
QG071-03	D09	HQP020899	NM_000388	CASR
QG071-03	D10	HQP020707	NM_003579	RAD54L
QG071-03	D11	HQP020418	NM_001004426	PLA2G6
QG071-03	D12	HQP020384	NM_004347	CASP5
QG071-03	E01	HQP020136	NM_003507	FZD7
QG071-03	E02	HQP020116	NM_004655	AXIN2
QG071-03	E03	HQP020040	NM_006534	NCOA3
QG071-03	E04	HQP019768	NM_130445	COL18A1
QG071-03	E05	HQP019452	NM_001742	CALCR
QG071-03	E06	HQP018739	NM_006300	ZNF230

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-03	E07	HQP018550	NM_001079539	XBP1
QG071-03	E08	HQP018545	NM_000553	WRN
QG071-03	E09	HQP018488	NM_003379	VIL2
QG071-03	E10	HQP018466	NM_080682	VCAM1
QG071-03	E11	HQP018404	NM_003356	UCP3
QG071-03	E12	HQP018343	NM_000372	TYR
QG071-03	F01	HQP018269	NM_006292	TSG101
QG071-03	F02	HQP018149	NM_001066	TNFRSF1B
QG071-03	F03	HQP018060	NM_000359	TGM1
QG071-03	F04	HQP018035	NM_003225	TFF1
QG071-03	F05	HQP018013	NM_007110	TEP1
QG071-03	F06	HQP017764	NM_007315	STAT1
QG071-03	F07	HQP017745	NM_001051	SSTR3
QG071-03	F08	HQP017615	NM_000454	SOD1
QG071-03	F09	HQP017367	NM_003040	SLC4A2
QG071-03	F10	HQP017303	NM_003037	SLAMF1
QG071-03	F11	HQP016980	NM_003019	SFTPD
QG071-03	F12	HQP016876	NM_022437	ABCG8
QG071-03	G01	HQP016844	NM_000057	BLM
QG071-03	G02	HQP016819	NM_005410	SEPP1
QG071-03	G03	HQP016744	NM_000450	SELE
QG071-03	G04	HQP016626	NM_002985	CCL5
QG071-03	G05	HQP016527	NM_021976	RXRB
QG071-03	G06	HQP016526	NM_002957	RXRA
QG071-03	G07	HQP016380	NM_002944	ROS1
QG071-03	G08	HQP016303	NM_021133	RNASEL
QG071-03	G09	HQP016212	NM_000657	BCL2
QG071-03	G10	HQP016170	NM_021632	ZNF350
QG071-03	G11	HQP016064	NM_018890	RAC1
QG071-03	G12	HQP016003	NM_021213	PCTP
QG071-03	H01	HGDC		
QG071-03	H02	HGDC		
QG071-03	H03	HQP006940	NM_002046	GAPDH
QG071-03	H04	HQP016381	NM_001101	ACTB
QG071-03	H05	HQP015171	NM_004048	B2M
QG071-03	H06	HQP006171	NM_012423	RPL13A
QG071-03	H07	HQP009026	NM_000194	HPRT1
QG071-03	H08	HQP054253	NR_003286	RN18S1
QG071-03	H09	RT		
QG071-03	H10	RT		
QG071-03	H11	PCR		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-03	H12	PCR		
QG071-04	A01	HQP015917	NM_001188	BAK1
QG071-04	A02	HQP015596	NM_000962	PTGS1
QG071-04	A03	HQP015190	NM_020165	RAD18
QG071-04	A04	HQP014808	NM_018657	MYNN
QG071-04	A05	HQP014564	NM_017902	HIF1AN
QG071-04	A06	HQP014292	NM_018315	FBXW7
QG071-04	A07	HQP014243	NM_018272	CASC1
QG071-04	A08	HQP014089	NM_018081	WDR79
QG071-04	A09	HQP013614	NM_007120	UGT1A4
QG071-04	A10	HQP013581	NM_019077	UGT1A7
QG071-04	A11	HQP013423	NM_002691	POLD1
QG071-04	A12	HQP013388	NM_017442	TLR9
QG071-04	B01	HQP013352	NM_000535	PMS2
QG071-04	B02	HQP013312	NM_000534	PMS1
QG071-04	B03	HQP013262	NM_005030	PLK1
QG071-04	B04	HQP013205	NM_001005376	PLAUR
QG071-04	B05	HQP013204	NM_002658	PLAU
QG071-04	B06	HQP013194	NM_000300	PLA2G2A
QG071-04	B07	HQP013128	NM_006216	SERPINE2
QG071-04	B08	HQP013104	NM_002634	PHB
QG071-04	B09	HQP013099	NM_000926	PGR
QG071-04	B10	HQP012100	NM_002562	P2RX7
QG071-04	B11	HQP012062	NM_001008505	OPRM1
QG071-04	B12	HQP012008	NM_000275	OCA2
QG071-04	C01	HQP011852	NM_001678	ATP1B2
QG071-04	C02	HQP011836	NM_004148	NINJ1
QG071-04	C03	HQP011813	NM_004556	NFKBIE
QG071-04	C04	HQP011810	NM_020529	NFKBIA
QG071-04	C05	HQP011807	NM_003998	NFKB1
QG071-04	C06	HQP011693	NM_000433	NCF2
QG071-04	C07	HQP011685	NM_002484	NUBP1
QG071-04	C08	HQP011582	NM_002462	MX1
QG071-04	C09	HQP011492	NM_002439	MSH3
QG071-04	C10	HQP011219	NM_002415	MIF
QG071-04	C11	HQP011179	NM_002402	MEST
QG071-04	C12	HQP011073	NM_015844	MBD1
QG071-04	D01	HQP011007	NM_000240	MAOA
QG071-04	D02	HQP010924	NM_001668	ARNT
QG071-04	D03	HQP010878	NM_002335	LRP5
QG071-04	D04	HQP010877	NM_002336	LRP6

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-04	D05	HQP010847	NM_000237	LPL
QG071-04	D06	HQP010626	NM_000236	LIPC
QG071-04	D07	HQP010610	NM_001175	ARHGDI3
QG071-04	D08	HQP010584	NM_002303	LEPR
QG071-04	D09	HQP010577	NM_000527	LDLR
QG071-04	D10	HQP010564	NM_000229	LCAT
QG071-04	D11	HQP010070	NM_002253	KDR
QG071-04	D12	HQP010008	NM_198686	RAB15
QG071-04	E01	HQP009851	NM_000215	JAK3
QG071-04	E02	HQP009818	NM_000212	ITGB3
QG071-04	E03	HQP009788	NM_005544	IRS1
QG071-04	E04	HQP009780	NM_001571	IRF3
QG071-04	E05	HQP009778	NM_002198	IRF1
QG071-04	E06	HQP009777	NM_004136	IREB2
QG071-04	E07	HQP009764	NM_000208	INSR
QG071-04	E08	HQP009718	NM_001562	IL18
QG071-04	E09	HQP009694	NM_005535	IL12RB1
QG071-04	E10	HQP009687	NM_000628	IL10RB
QG071-04	E11	HQP009658	NM_000878	IL2RB
QG071-04	E12	HQP009642	NM_000877	IL1R1
QG071-04	F01	HQP009556	NM_000041	APOE
QG071-04	F02	HQP009555	NM_002178	IGFBP6
QG071-04	F03	HQP009551	NM_000599	IGFBP5
QG071-04	F04	HQP009541	NM_000597	IGFBP2
QG071-04	F05	HQP009539	NM_000596	IGFBP1
QG071-04	F06	HQP009529	NM_000612	IGF2
QG071-04	F07	HQP009469	NM_000416	IFNGR1
QG071-04	F08	HQP009458	NM_000629	IFNAR1
QG071-04	F09	HQP009171	NM_000482	APOA4
QG071-04	F10	HQP009139	NM_001643	APOA2
QG071-04	F11	HQP009138	NM_004507	HUS1
QG071-04	F12	HQP009117	NM_000864	HTR1D
QG071-04	G01	HQP009116	NM_000863	HTR1B
QG071-04	G02	HQP009101	NM_001168	BIRC5
QG071-04	G03	HQP009072	NM_001166	BIRC2
QG071-04	G04	HQP009067	NM_000414	HSD17B4
QG071-04	G05	HQP009066	NM_002153	HSD17B2
QG071-04	G06	HQP009060	NM_000198	HSD3B2
QG071-04	G07	HQP008934	NM_001160	APAF1
QG071-04	G08	HQP008891	NM_000859	HMGCR
QG071-04	G09	HQP008832	NM_181054	HIF1A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-04	G10	HQP008757	NM_000410	HFE
QG071-04	G11	HQP008718	NM_000182	HADHA
QG071-04	G12	HQP008659	NM_014579	SLC39A2
QG071-04	H01	HGDC		
QG071-04	H02	HGDC		
QG071-04	H03	HQP006940	NM_002046	GAPDH
QG071-04	H04	HQP016381	NM_001101	ACTB
QG071-04	H05	HQP015171	NM_004048	B2M
QG071-04	H06	HQP006171	NM_012423	RPL13A
QG071-04	H07	HQP009026	NM_000194	HPRT1
QG071-04	H08	HQP054253	NR_003286	RN18S1
QG071-04	H09	RT		
QG071-04	H10	RT		
QG071-04	H11	PCR		
QG071-04	H12	PCR		
QG071-05	A01	HQP008605	NM_013347	RPA4
QG071-05	A02	HQP008469	NM_002093	GSK3B
QG071-05	A03	HQP008464	NM_005314	GRPR
QG071-05	A04	HQP008285	NM_002085	GPX4
QG071-05	A05	HQP007737	NM_000481	AMT
QG071-05	A06	HQP007643	NM_014475	DHDH
QG071-05	A07	HQP007230	NM_000583	GC
QG071-05	A08	HQP007224	NM_014365	HSPB8
QG071-05	A09	HQP006945	NM_015515	KRT23
QG071-05	A10	HQP006733	NM_012134	LMOD1
QG071-05	A11	HQP006547	NM_001469	XRCC6
QG071-05	A12	HQP006425	NM_001140	ALOX15
QG071-05	B01	HQP006359	NM_000698	ALOX5
QG071-05	B02	HQP006356	NM_000697	ALOX12
QG071-05	B03	HQP006261	NM_014324	AMACR
QG071-05	B04	HQP006228	NM_001018078	FPGS
QG071-05	B05	HQP006199	NM_021115	SEZ6L
QG071-05	B06	HQP006189	NM_012429	SEC14L2
QG071-05	B07	HQP006188	NM_005252	FOS
QG071-05	B08	HQP006120	NM_018672	ABCA5
QG071-05	B09	HQP006119	NM_080284	ABCA6
QG071-05	B10	HQP005629	NM_001453	FOXC1
QG071-05	B11	HQP005459	NM_001004439	ITGA11
QG071-05	B12	HQP005436	NM_004466	GPC5
QG071-05	C01	HQP005081	NM_000135	FANCA
QG071-05	C02	HQP005002	NM_001437	ESR2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-05	C03	HQP004995	NM_001626	AKT2
QG071-05	C04	HQP004950	NM_001979	EPHX2
QG071-05	C05	HQP004849	NM_152686	DNAJC18
QG071-05	C06	HQP004727	NM_005502	ABCA1
QG071-05	C07	HQP004598	NM_001406	EFNB3
QG071-05	C08	HQP004557	NM_001955	EDN1
QG071-05	C09	HQP004445	NM_000794	DRD1
QG071-05	C10	HQP003923	NM_000785	CYP27B1
QG071-05	C11	HQP003917	NM_000782	CYP24A1
QG071-05	C12	HQP003818	NM_000774	CYP2F1
QG071-05	D01	HQP003741	NM_000397	CYBB
QG071-05	D02	HQP003691	NM_001337	CX3CR1
QG071-05	D03	HQP003620	NM_004390	CTSH
QG071-05	D04	HQP003539	NM_001904	CTNNB1
QG071-05	D05	HQP003493	NM_001902	CTH
QG071-05	D06	HQP003395	NM_001033521	CSTF1
QG071-05	D07	HQP003171	NM_006437	PARP4
QG071-05	D08	HQP003158	NM_005211	CSF1R
QG071-05	D09	HQP002210	NM_000579	CCR5
QG071-05	D10	HQP001449	NM_133491	SAT2
QG071-05	D11	HQP001262	NM_001818	AKR1C4
QG071-05	D12	HQP000956	NM_000735	CGA
QG071-05	E01	HQP000932	NM_006636	MTHFD2
QG071-05	E02	HQP000853	NM_000078	CETP
QG071-05	E03	HQP000720	NM_006446	SLCO1B1
QG071-05	E04	HQP000405	NM_006066	AKR1A1
QG071-05	E05	HQP000356	NM_000076	CDKN1C
QG071-05	E06	HQP000285	NM_001799	CDK7
QG071-05	E07	HQP000245	NM_000075	CDK4
QG071-05	E08	HQP000179	NM_147185	AKAP9
QG071-05	E09	HQP000134	NM_012225	NUBP2
QG071-05	E10	HQP015595	NM_000315	PTH
QG071-05	E11	HQP015245	NM_020168	PAK6
QG071-05	E12	HQP015091	NM_001031849	MASP1
QG071-05	F01	HQP012054	NM_000911	OPRD1
QG071-05	F02	HQP010701	NM_005574	LMO2
QG071-05	F03	HQP009537	NM_004970	IGFALS
QG071-05	F04	HQP005011	NM_000031	ALAD
QG071-05	F05	HQP003625	NM_144713	FAM82A
QG071-05	F06	HQP000394	NM_006058	TNIP1
QG071-05	F07	HQP054686	NM_001123396	LOC729230

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-05	F08	HQP053990	NM_003352	SUMO1
QG071-05	F09	HQP053978	NM_001300	KLF6
QG071-05	F10	HQP023261	NM_001786	CDC2
QG071-05	F11	HQP023233	NM_012289	KEAP1
QG071-05	F12	HQP023134	NM_014707	HDAC9
QG071-05	G01	HQP022764	NM_004832	GSTO1
QG071-05	G02	HQP022686	NM_004261	SEP15
QG071-05	G03	HQP022490	NM_000591	CD14
QG071-05	G04	HQP021670	NM_006773	DDX18
QG071-05	G05	HQP021630	NM_003889	NR1I2
QG071-05	G06	HQP021604	NM_003879	CFLAR
QG071-05	G07	HQP021417	NM_003756	EIF3S3
QG071-05	G08	HQP020994	NM_032793	MFSD2
QG071-05	G09	HQP020984	NM_032782	HAVCR2
QG071-05	G10	HQP020427	NM_001226	CASP6
QG071-05	G11	HQP020331	NM_001013836	MAD1L1
QG071-05	G12	HQP019278	NM_006528	TFPI2
QG071-05	H01	HGDC		
QG071-05	H02	HGDC		
QG071-05	H03	HQP006940	NM_002046	GAPDH
QG071-05	H04	HQP016381	NM_001101	ACTB
QG071-05	H05	HQP015171	NM_004048	B2M
QG071-05	H06	HQP006171	NM_012423	RPL13A
QG071-05	H07	HQP009026	NM_000194	HPRT1
QG071-05	H08	HQP054253	NR_003286	RN18S1
QG071-05	H09	RT		
QG071-05	H10	RT		
QG071-05	H11	PCR		
QG071-05	H12	PCR		
QG071-06	A01	HQP019201	NM_024670	SUV39H2
QG071-06	A02	HQP019141	NM_024608	NEIL1
QG071-06	A03	HQP019015	NM_030762	BHLHB3
QG071-06	A04	HQP018895	NM_001080432	FTO
QG071-06	A05	HQP018813	NM_001005914	SEMA3B
QG071-06	A06	HQP018755	NM_001007257	PRDM2
QG071-06	A07	HQP018747	NM_013309	SLC30A4
QG071-06	A08	HQP018744	NM_001004434	SLC30A2
QG071-06	A09	HQP018411	NM_003357	SCGB1A1
QG071-06	A10	HQP018379	NM_003345	UBE2I
QG071-06	A11	HQP018328	NM_000474	TWIST1
QG071-06	A12	HQP018326	NM_001736	C5AR1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-06	B01	HQP018022	NM_001211	BUB1B
QG071-06	B02	HQP017996	NM_003211	TDG
QG071-06	B03	HQP017767	NM_003150	STAT3
QG071-06	B04	HQP017492	NM_021977	SLC22A3
QG071-06	B05	HQP017246	NM_022902	SLC30A5
QG071-06	B06	HQP017181	NM_022820	CYP3A43
QG071-06	B07	HQP016830	NM_003010	MAP2K4
QG071-06	B08	HQP016670	NM_001033886	CXCL12
QG071-06	B09	HQP016572	NM_002970	SAT1
QG071-06	B10	HQP016538	NM_005978	S100A2
QG071-06	B11	HQP016252	NM_002923	RGS2
QG071-06	B12	HQP016224	NM_002913	RFC1
QG071-06	C01	HQP016131	NM_000321	RB1
QG071-06	C02	HQP016125	NM_002890	RASA1
QG071-06	C03	HQP016084	NM_014226	RAGE
QG071-06	C04	HQP015978	NM_002859	PXN
QG071-06	C05	HQP015925	NM_002843	PTPRJ
QG071-06	C06	HQP015897	NM_006264	PTPN13
QG071-06	C07	HQP015892	NM_021187	CYP4F11
QG071-06	C08	HQP015667	NM_020750	XPO5
QG071-06	C09	HQP015586	NM_000961	PTGIS
QG071-06	C10	HQP015544	NM_000956	PTGER2
QG071-06	C11	HQP014907	NM_002755	MAP2K1
QG071-06	C12	HQP014846	NM_019101	APOM
QG071-06	D01	HQP014362	NM_018407	LAPTM4B
QG071-06	D02	HQP014161	NM_002716	PPP2R1B
QG071-06	D03	HQP013708	NM_017646	TRIT1
QG071-06	D04	HQP013627	NM_006238	PPARD
QG071-06	D05	HQP013617	NM_001001928	PPARA
QG071-06	D06	HQP013473	NM_000446	PON1
QG071-06	D07	HQP013450	NM_018975	TERF2IP
QG071-06	D08	HQP013425	NM_006231	POLE
QG071-06	D09	HQP012484	NM_016299	HSPA14
QG071-06	D10	HQP012393	NM_016009	SH3GLB1
QG071-06	D11	HQP012169	NM_015721	GEMIN4
QG071-06	D12	HQP012155	NM_002575	SERPINB2
QG071-06	E01	HQP012154	NM_000602	SERPINE1
QG071-06	E02	HQP011920	NM_005173	ATP2A3
QG071-06	E03	HQP011914	NM_002524	NRAS
QG071-06	E04	HQP011853	NM_000269	NME1
QG071-06	E05	HQP011800	NM_006164	NFE2L2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-06	E06	HQP011656	NM_002480	PPP1R12A
QG071-06	E07	HQP011559	NM_001018016	MUC1
QG071-06	E08	HQP011523	NM_012331	MSRA
QG071-06	E09	HQP011304	NM_001015052	MPG
QG071-06	E10	HQP011278	NM_002431	MNAT1
QG071-06	E11	HQP011262	NM_002424	MMP8
QG071-06	E12	HQP011245	NM_004529	MLLT3
QG071-06	F01	HQP010138	NM_001172	ARG2
QG071-06	F02	HQP010127	NM_000045	ARG1
QG071-06	F03	HQP010099	NM_000222	KIT
QG071-06	F04	HQP009828	NM_002214	ITGB8
QG071-06	F05	HQP009801	NM_000044	AR
QG071-06	F06	HQP009681	NM_001557	IL8RB
QG071-06	F07	HQP009667	NM_000564	IL5RA
QG071-06	F08	HQP009666	NM_000879	IL5
QG071-06	F09	HQP009643	NM_002182	IL1RAP
QG071-06	F10	HQP009353	NM_002168	IDH2
QG071-06	F11	HQP009351	NM_005896	IDH1
QG071-06	F12	HQP009218	NM_000384	APOB
QG071-06	G01	HQP009091	NM_001167	BIRC4
QG071-06	G02	HQP009089	NM_001540	HSPB1
QG071-06	G03	HQP009086	NM_006597	HSPA8
QG071-06	G04	HQP009036	NM_005343	HRAS
QG071-06	G05	HQP009026	NM_000194	HPRT1
QG071-06	G06	HQP009022	NM_000860	HPGD
QG071-06	G07	HQP008522	NM_014550	CARD10
QG071-06	G08	HQP008502	NM_005316	GTF2H1
QG071-06	G09	HQP008482	NM_000848	GSTM2
QG071-06	G10	HQP008478	NM_000847	GSTA3
QG071-06	G11	HQP008474	NM_000178	GSS
QG071-06	G12	HQP008473	NM_000637	GSR
QG071-06	H01	HGDC		
QG071-06	H02	HGDC		
QG071-06	H03	HQP006940	NM_002046	GAPDH
QG071-06	H04	HQP016381	NM_001101	ACTB
QG071-06	H05	HQP015171	NM_004048	B2M
QG071-06	H06	HQP006171	NM_012423	RPL13A
QG071-06	H07	HQP009026	NM_000194	HPRT1
QG071-06	H08	HQP054253	NR_003286	RN18S1
QG071-06	H09	RT		
QG071-06	H10	RT		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG071-06	H11	PCR		
QG071-06	H12	PCR		