

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-01	A01	HQP018175	NM_000546	TP53
QG068-01	A02	HQP003772	NM_000499	CYP1A1
QG068-01	A03	HQP018562	NM_006297	XRCC1
QG068-01	A04	HQP005087	NM_000690	ALDH2
QG068-01	A05	HQP003817	NM_000773	CYP2E1
QG068-01	A06	HQP004976	NM_000400	ERCC2
QG068-01	A07	HQP016204	NM_053056	CCND1
QG068-01	A08	HQP011547	NM_005957	MTHFR
QG068-01	A09	HQP002331	NM_000668	ADH1B
QG068-01	A10	HQP002413	NM_000669	ADH1C
QG068-01	A11	HQP001136	NM_000015	NAT2
QG068-01	A12	HQP018141	NM_000594	TNF
QG068-01	B01	HQP018564	NM_005432	XRCC3
QG068-01	B02	HQP015598	NM_000963	PTGS2
QG068-01	B03	HQP012021	NM_002542	OGG1
QG068-01	B04	HQP018475	NM_001025366	VEGFA
QG068-01	B05	HQP008849	NM_002116	HLA-A
QG068-01	B06	HQP004975	NM_202001	ERCC1
QG068-01	B07	HQP011255	NM_002421	MMP1
QG068-01	B08	HQP009641	NM_000576	IL1B
QG068-01	B09	HQP004948	NM_000120	EPHX1
QG068-01	B10	HQP004605	NM_005228	EGFR
QG068-01	B11	HQP000370	NM_058195	CDKN2A
QG068-01	B12	HQP018556	NM_004628	XPC
QG068-01	C01	HQP009685	NM_000572	IL10
QG068-01	C02	HQP004317	NM_000903	NQO1
QG068-01	C03	HQP018342	NM_001071	TYMS
QG068-01	C04	HQP011257	NM_002422	MMP3
QG068-01	C05	HQP011135	NM_002392	MDM2
QG068-01	C06	HQP008483	NM_000849	GSTM3
QG068-01	C07	HQP054047	BC008403	HLA-DRB1
QG068-01	C08	HQP018044	NM_000660	TGFB1
QG068-01	C09	HQP009670	NM_000600	IL6
QG068-01	C10	HQP010133	NM_004985	KRAS
QG068-01	C11	HQP009678	NM_000584	IL8
QG068-01	C12	HQP004599	NM_001963	EGF
QG068-01	D01	HQP023467	NM_000662	NAT1
QG068-01	D02	HQP017753	NM_000059	BRCA2
QG068-01	D03	HQP013150	NM_006218	PIK3CA
QG068-01	D04	HQP013100	NM_000927	ABCB1
QG068-01	D05	HQP011554	NM_000254	MTR

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-01	D06	HQP011263	NM_004994	MMP9
QG068-01	D07	HQP011256	NM_004530	MMP2
QG068-01	D08	HQP009662	NM_000589	IL4
QG068-01	D09	HQP003841	NM_000777	CYP3A5
QG068-01	D10	HQP003814	NM_000106	CYP2D6
QG068-01	D11	HQP018180	NM_005427	TP73
QG068-01	D12	HQP054045	BC004257	RET
QG068-01	E01	HQP018568	NM_021141	XRCC5
QG068-01	E02	HQP018563	NM_005431	XRCC2
QG068-01	E03	HQP017811	NM_177536	SULT1A1
QG068-01	E04	HQP011309	NM_000250	MPO
QG068-01	E05	HQP009718	NM_001562	IL18
QG068-01	E06	HQP009671	NM_000639	FASLG
QG068-01	E07	HQP003804	NM_000766	CYP2A13
QG068-01	E08	HQP003775	NM_000104	CYP1B1
QG068-01	E09	HQP004985	NM_000123	ERCC5
QG068-01	E10	HQP004414	NM_006892	DNMT3B
QG068-01	E11	HQP023466	NM_004360	CDH1
QG068-01	E12	HQP018565	NM_003401	XRCC4
QG068-01	F01	HQP018485	NM_000551	VHL
QG068-01	F02	HQP018474	NM_000376	VDR
QG068-01	F03	HQP017616	NM_000636	SOD2
QG068-01	F04	HQP016212	NM_000657	BCL2
QG068-01	F05	HQP009649	NM_000586	IL2
QG068-01	F06	HQP009645	NM_000577	IL1RN
QG068-01	F07	HQP004984	NM_005236	ERCC4
QG068-01	F08	HQP018176	NM_005657	TP53BP1
QG068-01	F09	HQP017733	NM_004333	BRAF
QG068-01	F10	HQP016689	NM_003000	SDHB
QG068-01	F11	HQP016670	NM_001033886	CXCL12
QG068-01	F12	HQP011807	NM_003998	NFKB1
QG068-01	G01	HQP011555	NM_002454	MTRR
QG068-01	G02	HQP009801	NM_000044	AR
QG068-01	G03	HQP009693	NM_002187	IL12B
QG068-01	G04	HQP009672	NM_000565	IL6R
QG068-01	G05	HQP009523	NM_000875	IGF1R
QG068-01	G06	HQP009061	NM_001641	APEX1
QG068-01	G07	HQP008832	NM_181054	HIF1A
QG068-01	G08	HQP005440	NM_022963	FGFR4
QG068-01	G09	HQP004986	NM_000124	ERCC6
QG068-01	G10	HQP004081	NM_000789	ACE

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-01	G11	HQP001396	NM_001005735	CHEK2
QG068-01	G12	HQP016706	NM_003002	SDHD
QG068-01	H01	HGDC		
QG068-01	H02	HGDC		
QG068-01	H03	HQP006940	NM_002046	GAPDH
QG068-01	H04	HQP016381	NM_001101	ACTB
QG068-01	H05	HQP015171	NM_004048	B2M
QG068-01	H06	HQP006171	NM_012423	RPL13A
QG068-01	H07	HQP009026	NM_000194	HPRT1
QG068-01	H08	HQP054253	NR_003286	RN18S1
QG068-01	H09	RT		
QG068-01	H10	RT		
QG068-01	H11	PCR		
QG068-01	H12	PCR		
QG068-02	A01	HQP018966	NM_001080124	CASP8
QG068-02	A02	HQP017673	NM_000582	SPP1
QG068-02	A03	HQP016131	NM_000321	RB1
QG068-02	A04	HQP016076	NM_002874	RAD23B
QG068-02	A05	HQP015897	NM_006264	PTPN13
QG068-02	A06	HQP015535	NM_000314	PTEN
QG068-02	A07	HQP013633	NM_005037	PPARG
QG068-02	A08	HQP011687	NM_002485	NBN
QG068-02	A09	HQP010613	NM_002312	LIG4
QG068-02	A10	HQP010612	NM_013975	LIG3
QG068-02	A11	HQP010609	NM_000234	LIG1
QG068-02	A12	HQP010581	NM_000230	LEP
QG068-02	B01	HQP009544	NM_000598	IGFBP3
QG068-02	B02	HQP009529	NM_000612	IGF2
QG068-02	B03	HQP009518	NM_000618	IGF1
QG068-02	B04	HQP009101	NM_001168	BIRC5
QG068-02	B05	HQP009036	NM_005343	HRAS
QG068-02	B06	HQP009024	NM_000038	APC
QG068-02	B07	HQP008476	NM_000846	GSTA2
QG068-02	B08	HQP006547	NM_001469	XRCC6
QG068-02	B09	HQP005447	NM_002012	FHIT
QG068-02	B10	HQP003811	NM_000771	CYP2C9
QG068-02	B11	HQP003810	NM_000770	CYP2C8
QG068-02	B12	HQP002671	NM_000754	COMT
QG068-02	C01	HQP000274	NM_001259	CDK6
QG068-02	C02	HQP000245	NM_000075	CDK4
QG068-02	C03	HQP011601	NM_005376	MYCL1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-02	C04	HQP054684	NM_001033	RRM1
QG068-02	C05	HQP022625	NM_004797	ADIPOQ
QG068-02	C06	HQP021736	NM_003925	MBD4
QG068-02	C07	HQP021542	NM_003834	RGS11
QG068-02	C08	HQP021023	NM_003617	RGS5
QG068-02	C09	HQP020481	NM_033338	CASP7
QG068-02	C10	HQP020297	NM_004346	CASP3
QG068-02	C11	HQP020116	NM_004655	AXIN2
QG068-02	C12	HQP019982	NM_001017388	TLR10
QG068-02	D01	HQP018983	NM_004639	BAT3
QG068-02	D02	HQP018263	NM_000368	TSC1
QG068-02	D03	HQP018149	NM_001066	TNFRSF1B
QG068-02	D04	HQP017900	NM_000544	TAP2
QG068-02	D05	HQP017899	NM_000593	TAP1
QG068-02	D06	HQP017812	NM_001056	SULT1C2
QG068-02	D07	HQP017744	NM_001050	SSTR2
QG068-02	D08	HQP016694	NM_001035511	SDHC
QG068-02	D09	HQP016389	NM_002945	RPA1
QG068-02	D10	HQP016252	NM_002923	RGS2
QG068-02	D11	HQP016087	NM_134424	RAD52
QG068-02	D12	HQP015667	NM_020750	XPO5
QG068-02	E01	HQP013581	NM_019077	UGT1A7
QG068-02	E02	HQP013473	NM_000446	PON1
QG068-02	E03	HQP013147	NM_002647	PIK3C3
QG068-02	E04	HQP013144	NM_002645	PIK3C2A
QG068-02	E05	HQP013143	NM_002644	PIGR
QG068-02	E06	HQP012996	NM_016362	GHRL
QG068-02	E07	HQP012501	NM_016341	PLCE1
QG068-02	E08	HQP012420	NM_002592	PCNA
QG068-02	E09	HQP012169	NM_015721	GEMIN4
QG068-02	E10	HQP011866	NM_000625	NOS2A
QG068-02	E11	HQP011597	NM_002467	MYC
QG068-02	E12	HQP011492	NM_002439	MSH3
QG068-02	F01	HQP011491	NM_000251	MSH2
QG068-02	F02	HQP011266	NM_002426	MMP12
QG068-02	F03	HQP011235	NM_000249	MLH1
QG068-02	F04	HQP010584	NM_002303	LEPR
QG068-02	F05	HQP009692	NM_000882	IL12A
QG068-02	F06	HQP009681	NM_001557	IL8RB
QG068-02	F07	HQP009679	NM_000634	IL8RA
QG068-02	F08	HQP009642	NM_000877	IL1R1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-02	F09	HQP009640	NM_000575	IL1A
QG068-02	F10	HQP009019	NM_005143	HP
QG068-02	F11	HQP007755	NM_000516	GNAS
QG068-02	F12	HQP007385	NM_000515	GH1
QG068-02	G01	HQP007287	NM_012199	EIF2C1
QG068-02	G02	HQP006460	NM_145043	NEIL2
QG068-02	G03	HQP006426	NM_004958	FRAP1
QG068-02	G04	HQP006356	NM_000697	ALOX12
QG068-02	G05	HQP005736	NM_015061	JMJD2C
QG068-02	G06	HQP004991	NM_001014431	AKT1
QG068-02	G07	HQP004658	NM_001621	AHR
QG068-02	G08	HQP003888	NM_000102	CYP17A1
QG068-02	G09	HQP003809	NM_000769	CYP2C19
QG068-02	G10	HQP003499	NM_001037631	CTLA4
QG068-02	G11	HQP003120	NM_001618	PARP1
QG068-02	G12	HQP002701	NM_000673	ADH7
QG068-02	H01	HGDC		
QG068-02	H02	HGDC		
QG068-02	H03	HQP006940	NM_002046	GAPDH
QG068-02	H04	HQP016381	NM_001101	ACTB
QG068-02	H05	HQP015171	NM_004048	B2M
QG068-02	H06	HQP006171	NM_012423	RPL13A
QG068-02	H07	HQP009026	NM_000194	HPRT1
QG068-02	H08	HQP054253	NR_003286	RN18S1
QG068-02	H09	RT		
QG068-02	H10	RT		
QG068-02	H11	PCR		
QG068-02	H12	PCR		
QG068-03	A01	HQP002659	NM_000672	ADH6
QG068-03	A02	HQP002469	NM_000670	ADH4
QG068-03	A03	HQP002210	NM_000579	CCR5
QG068-03	A04	HQP001416	NM_007204	DDX20
QG068-03	A05	HQP017743	NM_001049	SSTR1
QG068-03	A06	HQP011544	NM_002452	NUDT1
QG068-03	A07	HQP007397	NM_000823	GHRHR
QG068-03	A08	HQP054060	NM_021923	FGFRL1
QG068-03	A09	HQP054057	BC071181	TGFBR1
QG068-03	A10	HQP054030	NM_130398	EXO1
QG068-03	A11	HQP053999	NM_004955	SLC29A1
QG068-03	A12	HQP053998	NM_004873	BAG5
QG068-03	B01	HQP053978	NM_001300	KLF6

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-03	B02	HQP053963	NM_001048	SST
QG068-03	B03	HQP023414	NM_005847	SLC23A1
QG068-03	B04	HQP023409	NM_001790	CDC25C
QG068-03	B05	HQP023387	NM_007335	DLEC1
QG068-03	B06	HQP023385	NM_001789	CDC25A
QG068-03	B07	HQP023346	NM_006039	MRC2
QG068-03	B08	HQP023341	NM_138448	ACYP2
QG068-03	B09	HQP023240	NM_014781	RB1CC1
QG068-03	B10	HQP023203	NM_001785	CDA
QG068-03	B11	HQP023143	NM_014712	SETD1A
QG068-03	B12	HQP023005	NM_001778	CD48
QG068-03	C01	HQP023003	NM_004296	RGS6
QG068-03	C02	HQP022909	NM_004284	CHD1L
QG068-03	C03	HQP022884	NM_004881	TP53I3
QG068-03	C04	HQP022872	NM_004281	BAG3
QG068-03	C05	HQP022676	NM_004260	RECQL4
QG068-03	C06	HQP022490	NM_000591	CD14
QG068-03	C07	HQP022316	NM_000616	CD4
QG068-03	C08	HQP022170	NM_004708	PDCD5
QG068-03	C09	HQP022099	NM_004690	LATS1
QG068-03	C10	HQP022013	NM_004673	ANGPTL1
QG068-03	C11	HQP021906	NM_001239	CCNH
QG068-03	C12	HQP021829	NM_014240	LIMD1
QG068-03	D01	HQP021819	NM_001238	CCNE1
QG068-03	D02	HQP021737	NM_003927	MBD2
QG068-03	D03	HQP021727	NM_031966	CCNB1
QG068-03	D04	HQP021672	NM_006024	TAX1BP1
QG068-03	D05	HQP021630	NM_003889	NR1I2
QG068-03	D06	HQP021613	NM_006017	PROM1
QG068-03	D07	HQP021604	NM_003879	CFLAR
QG068-03	D08	HQP021603	NM_003878	GGH
QG068-03	D09	HQP021597	NM_003873	NRP1
QG068-03	D10	HQP021557	NM_003844	TNFRSF10A
QG068-03	D11	HQP021553	NM_003842	TNFRSF10B
QG068-03	D12	HQP021518	NM_000071	CBS
QG068-03	E01	HQP021502	NM_003810	TNFSF10
QG068-03	E02	HQP021459	NM_003786	ABCC3
QG068-03	E03	HQP021388	NM_003742	ABCB11
QG068-03	E04	HQP021385	NM_003739	AKR1C3
QG068-03	E05	HQP021313	NM_001753	CAV1
QG068-03	E06	HQP021205	NM_033345	RGS8

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-03	E07	HQP021125	NM_001005476	PKP4
QG068-03	E08	HQP020758	NM_003587	DHX16
QG068-03	E09	HQP020332	NM_001225	CASP4
QG068-03	E10	HQP020134	NM_012193	FZD4
QG068-03	E11	HQP020118	NM_006768	BRAP
QG068-03	E12	HQP020089	NM_000502	EPX
QG068-03	F01	HQP020070	NM_001748	CAPN2
QG068-03	F02	HQP020042	NM_003489	NRIP1
QG068-03	F03	HQP019829	NM_030782	CLPTM1L
QG068-03	F04	HQP019768	NM_130445	COL18A1
QG068-03	F05	HQP019568	NM_025113	C13orf18
QG068-03	F06	HQP019312	NM_024782	NHEJ1
QG068-03	F07	HQP019310	NM_024780	TMC5
QG068-03	F08	HQP019284	NM_024757	EHMT1
QG068-03	F09	HQP019267	NM_024743	UGT2A3
QG068-03	F10	HQP019238	NM_017412	FZD3
QG068-03	F11	HQP019189	NM_001006636	GTDC1
QG068-03	F12	HQP019141	NM_024608	NEIL1
QG068-03	G01	HQP018991	NM_033177	BAT4
QG068-03	G02	HQP018895	NM_001080432	FTO
QG068-03	G03	HQP018893	NM_001077198	ATG9A
QG068-03	G04	HQP018802	NM_001008540	CXCR4
QG068-03	G05	HQP018668	NM_001216	CA9
QG068-03	G06	HQP018622	NM_006965	ZNF24
QG068-03	G07	HQP018552	NM_000379	XDH
QG068-03	G08	HQP018545	NM_000553	WRN
QG068-03	G09	HQP018540	NM_004185	WNT2B
QG068-03	G10	HQP018539	NM_004626	WNT11
QG068-03	G11	HQP018425	NM_000373	UMPS
QG068-03	G12	HQP018372	NM_152653	UBE2E2
QG068-03	H01	HGDC		
QG068-03	H02	HGDC		
QG068-03	H03	HQP006940	NM_002046	GAPDH
QG068-03	H04	HQP016381	NM_001101	ACTB
QG068-03	H05	HQP015171	NM_004048	B2M
QG068-03	H06	HQP006171	NM_012423	RPL13A
QG068-03	H07	HQP009026	NM_000194	HPRT1
QG068-03	H08	HQP054253	NR_003286	RN18S1
QG068-03	H09	RT		
QG068-03	H10	RT		
QG068-03	H11	PCR		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-03	H12	PCR		
QG068-04	A01	HQP018265	NM_000548	TSC2
QG068-04	A02	HQP018144	NM_006291	TNFAIP2
QG068-04	A03	HQP018115	NM_003265	TLR3
QG068-04	A04	HQP018113	NM_003263	TLR1
QG068-04	A05	HQP018067	NM_000361	THBD
QG068-04	A06	HQP018022	NM_001211	BUB1B
QG068-04	A07	HQP017906	NM_004178	TARBP2
QG068-04	A08	HQP017897	NM_006755	TALDO1
QG068-04	A09	HQP017640	NM_138473	SP1
QG068-04	A10	HQP017631	NM_004331	BNIP3L
QG068-04	A11	HQP017563	NM_005631	SMO
QG068-04	A12	HQP017482	NM_194255	SLC19A1
QG068-04	B01	HQP017462	NM_000452	SLC10A2
QG068-04	B02	HQP017124	NM_022743	SMYD3
QG068-04	B03	HQP017001	NM_000337	SGCD
QG068-04	B04	HQP016980	NM_003019	SFTPD
QG068-04	B05	HQP016867	NM_022349	MS4A6A
QG068-04	B06	HQP016801	NM_022162	NOD2
QG068-04	B07	HQP016634	NM_001713	BHMT
QG068-04	B08	HQP016626	NM_002985	CCL5
QG068-04	B09	HQP016621	NM_002982	CCL2
QG068-04	B10	HQP016526	NM_002957	RXRA
QG068-04	B11	HQP016471	NM_003161	RPS6KB1
QG068-04	B12	HQP016470	NM_004586	RPS6KA3
QG068-04	C01	HQP016391	NM_002947	RPA3
QG068-04	C02	HQP016295	NM_002935	RNASE3
QG068-04	C03	HQP016281	NM_002931	RING1
QG068-04	C04	HQP016269	NM_002928	RGS16
QG068-04	C05	HQP016268	NM_002927	RGS13
QG068-04	C06	HQP016263	NM_001005339	RGS10
QG068-04	C07	HQP016262	NM_002924	RGS7
QG068-04	C08	HQP016260	NM_004050	BCL2L2
QG068-04	C09	HQP016259	NM_005613	RGS4
QG068-04	C10	HQP016253	NM_017790	RGS3
QG068-04	C11	HQP016223	NM_002912	REV3L
QG068-04	C12	HQP016160	NM_002895	RBL1
QG068-04	D01	HQP016125	NM_002890	RASA1
QG068-04	D02	HQP015983	NM_001017423	ALDH18A1
QG068-04	D03	HQP015946	NM_000465	BARD1
QG068-04	D04	HQP015925	NM_002843	PTPRJ



Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-04	D05	HQP015917	NM_001188	BAK1
QG068-04	D06	HQP015845	NM_020964	KIAA1632
QG068-04	D07	HQP015699	NM_020787	ZNF624
QG068-04	D08	HQP015669	NM_020752	GPR158
QG068-04	D09	HQP015596	NM_000962	PTGS1
QG068-04	D10	HQP015586	NM_000961	PTGIS
QG068-04	D11	HQP015190	NM_020165	RAD18
QG068-04	D12	HQP014905	NM_016205	PDGFC
QG068-04	E01	HQP014706	NM_002737	PRKCA
QG068-04	E02	HQP014564	NM_017902	HIF1AN
QG068-04	E03	HQP014454	NM_018725	IL17RB
QG068-04	E04	HQP014362	NM_018407	LAPTM4B
QG068-04	E05	HQP014089	NM_018081	WDR79
QG068-04	E06	HQP013874	NM_017823	DUSP23
QG068-04	E07	HQP013627	NM_006238	PPARD
QG068-04	E08	HQP013617	NM_001001928	PPARA
QG068-04	E09	HQP013614	NM_007120	UGT1A4
QG068-04	E10	HQP013596	NM_001184	ATR
QG068-04	E11	HQP013583	NM_205862	UGT1A6
QG068-04	E12	HQP013467	NM_000939	POMC
QG068-04	F01	HQP013431	NM_000938	POLR2B
QG068-04	F02	HQP013422	NM_002690	POLB
QG068-04	F03	HQP013388	NM_017442	TLR9
QG068-04	F04	HQP013352	NM_000535	PMS2
QG068-04	F05	HQP013312	NM_000534	PMS1
QG068-04	F06	HQP013262	NM_005030	PLK1
QG068-04	F07	HQP013204	NM_002658	PLAU
QG068-04	F08	HQP013154	NM_002649	PIK3CG
QG068-04	F09	HQP013151	NM_006219	PIK3CB
QG068-04	F10	HQP013149	NM_001695	ATP6V1C1
QG068-04	F11	HQP013135	NM_006217	SERPINI2
QG068-04	F12	HQP013101	NM_000443	ABCB4
QG068-04	G01	HQP013099	NM_000926	PGR
QG068-04	G02	HQP012958	NM_016169	SUFU
QG068-04	G03	HQP012889	NM_002609	PDGFRB
QG068-04	G04	HQP012857	NM_033016	PDGFB
QG068-04	G05	HQP012796	NM_016448	DTL
QG068-04	G06	HQP012755	NM_001037872	REV1
QG068-04	G07	HQP012725	NM_016218	POLK
QG068-04	G08	HQP012613	NM_016593	CYP39A1
QG068-04	G09	HQP012605	NM_016583	PLUNC

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-04	G10	HQP012444	NM_016133	INSIG2
QG068-04	G11	HQP012156	NM_002576	PAK1
QG068-04	G12	HQP012154	NM_000602	SERPINE1
QG068-04	H01	HGDC		
QG068-04	H02	HGDC		
QG068-04	H03	HQP006940	NM_002046	GAPDH
QG068-04	H04	HQP016381	NM_001101	ACTB
QG068-04	H05	HQP015171	NM_004048	B2M
QG068-04	H06	HQP006171	NM_012423	RPL13A
QG068-04	H07	HQP009026	NM_000194	HPRT1
QG068-04	H08	HQP054253	NR_003286	RN18S1
QG068-04	H09	RT		
QG068-04	H10	RT		
QG068-04	H11	PCR		
QG068-04	H12	PCR		
QG068-05	A01	HQP012131	NM_015713	RRM2B
QG068-05	A02	HQP011874	NM_000905	NPY
QG068-05	A03	HQP011868	NM_000603	NOS3
QG068-05	A04	HQP011841	NM_001677	ATP1B1
QG068-05	A05	HQP011812	NM_002503	NFKBIB
QG068-05	A06	HQP011810	NM_020529	NFKBIA
QG068-05	A07	HQP011706	NM_002490	NDUFA6
QG068-05	A08	HQP011609	NM_002470	MYH3
QG068-05	A09	HQP011325	NM_019899	ABCC1
QG068-05	A10	HQP011278	NM_002431	MNAT1
QG068-05	A11	HQP011268	NM_004995	MMP14
QG068-05	A12	HQP011264	NM_002425	MMP10
QG068-05	B01	HQP011241	NM_005933	MLL
QG068-05	B02	HQP011181	NM_000245	MET
QG068-05	B03	HQP011171	NM_000244	MEN1
QG068-05	B04	HQP011140	NM_002393	MDM4
QG068-05	B05	HQP011073	NM_015844	MBD1
QG068-05	B06	HQP010966	NM_005904	SMAD7
QG068-05	B07	HQP010960	NM_005902	SMAD3
QG068-05	B08	HQP010959	NM_001003652	SMAD2
QG068-05	B09	HQP010958	NM_001003688	SMAD1
QG068-05	B10	HQP010924	NM_001668	ARNT
QG068-05	B11	HQP010912	NM_206943	LTBP1
QG068-05	B12	HQP010868	NM_006152	LRMP
QG068-05	C01	HQP010070	NM_002253	KDR
QG068-05	C02	HQP009818	NM_000212	ITGB3

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-05	C03	HQP009803	NM_002207	ITGA9
QG068-05	C04	HQP009800	NM_002206	ITGA7
QG068-05	C05	HQP009788	NM_005544	IRS1
QG068-05	C06	HQP009780	NM_001571	IRF3
QG068-05	C07	HQP009756	NM_001567	INPPL1
QG068-05	C08	HQP009710	NM_002189	IL15RA
QG068-05	C09	HQP009708	NM_000585	IL15
QG068-05	C10	HQP009697	NM_002188	IL13
QG068-05	C11	HQP009694	NM_005535	IL12RB1
QG068-05	C12	HQP009666	NM_000879	IL5
QG068-05	D01	HQP009664	NM_000418	IL4R
QG068-05	D02	HQP009558	NM_001553	IGFBP7
QG068-05	D03	HQP009556	NM_000041	APOE
QG068-05	D04	HQP009551	NM_000599	IGFBP5
QG068-05	D05	HQP009541	NM_000597	IGFBP2
QG068-05	D06	HQP009539	NM_000596	IGFBP1
QG068-05	D07	HQP009472	NM_005534	IFNGR2
QG068-05	D08	HQP009469	NM_000416	IFNGR1
QG068-05	D09	HQP009460	NM_000874	IFNAR2
QG068-05	D10	HQP009140	NM_002160	TNC
QG068-05	D11	HQP009138	NM_004507	HUS1
QG068-05	D12	HQP009091	NM_001167	BIRC4
QG068-05	E01	HQP009067	NM_000414	HSD17B4
QG068-05	E02	HQP009066	NM_002153	HSD17B2
QG068-05	E03	HQP009065	NM_000197	HSD17B3
QG068-05	E04	HQP009064	NM_000413	HSD17B1
QG068-05	E05	HQP009060	NM_000198	HSD3B2
QG068-05	E06	HQP009059	NM_000862	HSD3B1
QG068-05	E07	HQP008934	NM_001160	APAF1
QG068-05	E08	HQP008859	NM_033554	HLA-DPA1
QG068-05	E09	HQP008808	NM_021155	CD209
QG068-05	E10	HQP008757	NM_000410	HFE
QG068-05	E11	HQP008683	NM_000180	GUCY2D
QG068-05	E12	HQP008669	NM_020804	PACSIN1
QG068-05	F01	HQP008554	NM_012092	ICOS
QG068-05	F02	HQP008493	NM_000179	MSH6
QG068-05	F03	HQP008490	NM_001513	GSTZ1
QG068-05	F04	HQP008469	NM_002093	GSK3B
QG068-05	F05	HQP008340	NM_000830	GRIK1
QG068-05	F06	HQP008152	NM_173827	COX18
QG068-05	F07	HQP008118	NM_173653	SLC9A9

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-05	F08	HQP007803	NM_000173	GP1BA
QG068-05	F09	HQP007767	NM_002075	GNB3
QG068-05	F10	HQP007702	NM_005270	GLI2
QG068-05	F11	HQP007613	NM_014452	TNFRSF21
QG068-05	F12	HQP007558	NM_012154	EIF2C2
QG068-05	G01	HQP007531	NM_148674	SMC1B
QG068-05	G02	HQP007506	NM_005268	GJB5
QG068-05	G03	HQP007476	NM_014394	GHITM
QG068-05	G04	HQP007396	NM_021081	GHRH
QG068-05	G05	HQP007391	NM_022556	GH2
QG068-05	G06	HQP007304	NM_012419	RGS17
QG068-05	G07	HQP007224	NM_014365	HSPB8
QG068-05	G08	HQP007098	NM_015668	RGS22
QG068-05	G09	HQP006767	NM_014351	SULT4A1
QG068-05	G10	HQP006764	NM_012473	TXN2
QG068-05	G11	HQP006583	NM_178510	ANKK1
QG068-05	G12	HQP006569	NM_001470	GABBR1
QG068-05	H01	HGDC		
QG068-05	H02	HGDC		
QG068-05	H03	HQP006940	NM_002046	GAPDH
QG068-05	H04	HQP016381	NM_001101	ACTB
QG068-05	H05	HQP015171	NM_004048	B2M
QG068-05	H06	HQP006171	NM_012423	RPL13A
QG068-05	H07	HQP009026	NM_000194	HPRT1
QG068-05	H08	HQP054253	NR_003286	RN18S1
QG068-05	H09	RT		
QG068-05	H10	RT		
QG068-05	H11	PCR		
QG068-05	H12	PCR		
QG068-06	A01	HQP006478	NM_152756	RICTOR
QG068-06	A02	HQP006441	NM_000350	ABCA4
QG068-06	A03	HQP006424	NM_145659	IL27
QG068-06	A04	HQP006359	NM_000698	ALOX5
QG068-06	A05	HQP006228	NM_001018078	FPGS
QG068-06	A06	HQP006078	NM_001017524	SIRT3
QG068-06	A07	HQP006010	NM_033071	SYNE1
QG068-06	A08	HQP005925	NM_015210	KIAA0802
QG068-06	A09	HQP005899	NM_182734	PLCB1
QG068-06	A10	HQP005879	NM_002019	FLT1
QG068-06	A11	HQP005747	NM_002015	FOXO1A
QG068-06	A12	HQP005629	NM_001453	FOXC1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-06	B01	HQP005416	NM_002010	FGF9
QG068-06	B02	HQP005408	NM_004464	FGF5
QG068-06	B03	HQP005403	NM_002006	FGF2
QG068-06	B04	HQP005391	NM_004111	FEN1
QG068-06	B05	HQP005280	NM_021642	FCGR2A
QG068-06	B06	HQP005084	NM_001018115	FANCD2
QG068-06	B07	HQP005081	NM_000135	FANCA
QG068-06	B08	HQP005069	NM_000129	F13A1
QG068-06	B09	HQP005058	NM_000130	F5
QG068-06	B10	HQP005052	NM_000506	F2
QG068-06	B11	HQP005049	NM_001991	EZH1
QG068-06	B12	HQP004998	NM_000125	ESR1
QG068-06	C01	HQP004995	NM_001626	AKT2
QG068-06	C02	HQP004992	NM_001433	ERN1
QG068-06	C03	HQP004983	NM_000122	ERCC3
QG068-06	C04	HQP004950	NM_001979	EPHX2
QG068-06	C05	HQP004526	NM_004091	E2F2
QG068-06	C06	HQP004494	NM_000029	AGT
QG068-06	C07	HQP004446	NM_000795	DRD2
QG068-06	C08	HQP004439	NM_000110	DPYD
QG068-06	C09	HQP004187	NM_174914	UGT3A2
QG068-06	C10	HQP004114	NM_000107	DDB2
QG068-06	C11	HQP003923	NM_000785	CYP27B1
QG068-06	C12	HQP003917	NM_000782	CYP24A1
QG068-06	D01	HQP003904	NM_000103	CYP19A1
QG068-06	D02	HQP003818	NM_000774	CYP2F1
QG068-06	D03	HQP003813	NM_000772	CYP2C18
QG068-06	D04	HQP003691	NM_001337	CX3CR1
QG068-06	D05	HQP003539	NM_001904	CTNNB1
QG068-06	D06	HQP003488	NM_001901	CTGF
QG068-06	D07	HQP003218	NM_004383	CSK
QG068-06	D08	HQP003209	NM_203394	E2F7
QG068-06	D09	HQP003171	NM_006437	PARP4
QG068-06	D10	HQP003013	NM_001884	HAPLN1
QG068-06	D11	HQP002612	NM_004370	COL12A1
QG068-06	D12	HQP002542	NM_000671	ADH5
QG068-06	E01	HQP002408	NM_016083	CNR1
QG068-06	E02	HQP001889	NM_052963	TOP1MT
QG068-06	E03	HQP001887	NM_133445	GRIN3A
QG068-06	E04	HQP001778	NM_178270	ATG4A
QG068-06	E05	HQP001712	NM_052917	GALNT13

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-06	E06	HQP001682	NM_052890	PGLYRP2
QG068-06	E07	HQP001429	NM_007215	POLG2
QG068-06	E08	HQP001414	NM_007202	AKAP10
QG068-06	E09	HQP001399	NM_007195	POLI
QG068-06	E10	HQP001374	NM_007182	RASSF1
QG068-06	E11	HQP001362	NM_021020	LZTS1
QG068-06	E12	HQP001282	NM_001274	CHEK1
QG068-06	F01	HQP001244	NM_022105	DIDO1
QG068-06	F02	HQP001111	NM_006828	ASCC3
QG068-06	F03	HQP001072	NM_015029	POP1
QG068-06	F04	HQP001051	NM_025256	EHMT2
QG068-06	F05	HQP001016	NM_013261	PPARGC1A
QG068-06	F06	HQP000963	NM_012190	ALDH1L1
QG068-06	F07	HQP000932	NM_006636	MTHFD2
QG068-06	F08	HQP000888	NM_006615	CAPN9
QG068-06	F09	HQP000771	NM_006480	RGS14
QG068-06	F10	HQP000765	NM_006476	ATP5L
QG068-06	F11	HQP000657	NM_006407	ARL6IP5
QG068-06	F12	HQP000508	NM_006320	PGRMC2
QG068-06	G01	HQP000462	NM_006094	DLC1
QG068-06	G02	HQP000459	NM_006092	NOD1
QG068-06	G03	HQP000415	NM_006068	TLR6
QG068-06	G04	HQP000406	NM_006067	COX4NB
QG068-06	G05	HQP000405	NM_006066	AKR1A1
QG068-06	G06	HQP000396	NM_001262	CDKN2C
QG068-06	G07	HQP000289	NM_013404	MSLN
QG068-06	G08	HQP000285	NM_001799	CDK7
QG068-06	G09	HQP000261	NM_004935	CDK5
QG068-06	G10	HQP000179	NM_147185	AKAP9
QG068-06	G11	HQP000018	NM_001792	CDH2
QG068-06	G12	HQP020119	NM_003503	CDC7
QG068-06	H01	HGDC		
QG068-06	H02	HGDC		
QG068-06	H03	HQP006940	NM_002046	GAPDH
QG068-06	H04	HQP016381	NM_001101	ACTB
QG068-06	H05	HQP015171	NM_004048	B2M
QG068-06	H06	HQP006171	NM_012423	RPL13A
QG068-06	H07	HQP009026	NM_000194	HPRT1
QG068-06	H08	HQP054253	NR_003286	RN18S1
QG068-06	H09	RT		
QG068-06	H10	RT		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG068-06	H11	PCR		
QG068-06	H12	PCR		