

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-01	A01	HQP005890	NM_004119	FLT3
QG070-01	A02	HQP011547	NM_005957	MTHFR
QG070-01	A03	HQP008849	NM_002116	HLA-A
QG070-01	A04	HQP054047	BC008403	HLA-DRB1
QG070-01	A05	HQP018175	NM_000546	TP53
QG070-01	A06	HQP003772	NM_000499	CYP1A1
QG070-01	A07	HQP000615	NM_004364	CEBPA
QG070-01	A08	HQP004317	NM_000903	NQO1
QG070-01	A09	HQP009850	NM_004972	JAK2
QG070-01	A10	HQP013100	NM_000927	ABCB1
QG070-01	A11	HQP018562	NM_006297	XRCC1
QG070-01	A12	HQP018546	NM_000378	WT1
QG070-01	B01	HQP003841	NM_000777	CYP3A5
QG070-01	B02	HQP011914	NM_002524	NRAS
QG070-01	B03	HQP011135	NM_002392	MDM2
QG070-01	B04	HQP004976	NM_000400	ERCC2
QG070-01	B05	HQP009351	NM_005896	IDH1
QG070-01	B06	HQP010099	NM_000222	KIT
QG070-01	B07	HQP018213	NM_000367	TPMT
QG070-01	B08	HQP018342	NM_001071	TYMS
QG070-01	B09	HQP018141	NM_000594	TNF
QG070-01	B10	HQP011241	NM_005933	MLL
QG070-01	B11	HQP003817	NM_000773	CYP2E1
QG070-01	B12	HQP003814	NM_000106	CYP2D6
QG070-01	C01	HQP013696	NM_017628	KIAA1546
QG070-01	C02	HQP009353	NM_002168	IDH2
QG070-01	C03	HQP021347	NM_001754	RUNX1
QG070-01	C04	HQP018564	NM_005432	XRCC3
QG070-01	C05	HQP016212	NM_000657	BCL2
QG070-01	C06	HQP012100	NM_002562	P2RX7
QG070-01	C07	HQP010133	NM_004985	KRAS
QG070-01	C08	HQP009781	NM_002460	IRF4
QG070-01	C09	HQP009685	NM_000572	IL10
QG070-01	C10	HQP008757	NM_000410	HFE
QG070-01	C11	HQP011235	NM_000249	MLH1
QG070-01	C12	HQP001136	NM_000015	NAT2
QG070-01	D01	HQP017490	NM_003057	SLC22A1
QG070-01	D02	HQP011687	NM_002485	NBN
QG070-01	D03	HQP011309	NM_000250	MPO
QG070-01	D04	HQP010095	NM_013289	KIR3DL1
QG070-01	D05	HQP010083	NM_014218	KIR2DL1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-01	D06	HQP004985	NM_000123	ERCC5
QG070-01	D07	HQP018475	NM_001025366	VEGFA
QG070-01	D08	HQP016801	NM_022162	NOD2
QG070-01	D09	HQP016204	NM_053056	CCND1
QG070-01	D10	HQP011555	NM_002454	MTRR
QG070-01	D11	HQP010086	NM_015868	KIR2DL3
QG070-01	D12	HQP009851	NM_000215	JAK3
QG070-01	E01	HQP009640	NM_000575	IL1A
QG070-01	E02	HQP004975	NM_202001	ERCC1
QG070-01	E03	HQP004287	NM_015338	ASXL1
QG070-01	E04	HQP003811	NM_000771	CYP2C9
QG070-01	E05	HQP003809	NM_000769	CYP2C19
QG070-01	E06	HQP023467	NM_000662	NAT1
QG070-01	E07	HQP022337	NM_014207	CD5
QG070-01	E08	HQP020297	NM_004346	CASP3
QG070-01	E09	HQP020207	NM_001223	CASP1
QG070-01	E10	HQP018966	NM_001080124	CASP8
QG070-01	E11	HQP018563	NM_005431	XRCC2
QG070-01	E12	HQP018149	NM_001066	TNFRSF1B
QG070-01	F01	HQP010092	NM_012313	KIR2DS3
QG070-01	F02	HQP010089	NM_002255	KIR2DL4
QG070-01	F03	HQP009662	NM_000589	IL4
QG070-01	F04	HQP009641	NM_000576	IL1B
QG070-01	F05	HQP009101	NM_001168	BIRC5
QG070-01	F06	HQP009061	NM_001641	APEX1
QG070-01	F07	HQP005280	NM_021642	FCGR2A
QG070-01	F08	HQP001810	NM_153443	KIR3DL3
QG070-01	F09	HQP000397	NM_006060	IKZF1
QG070-01	F10	HQP010096	NM_006737	KIR3DL2
QG070-01	F11	HQP022870	NM_001775	CD38
QG070-01	F12	HQP021557	NM_003844	TNFRSF10A
QG070-01	G01	HQP020515	NM_032199	ARID5B
QG070-01	G02	HQP020481	NM_033338	CASP7
QG070-01	G03	HQP018565	NM_003401	XRCC4
QG070-01	G04	HQP018556	NM_004628	XPC
QG070-01	G05	HQP018474	NM_000376	VDR
QG070-01	G06	HQP018466	NM_080682	VCAM1
QG070-01	G07	HQP018340	NM_003331	TYK2
QG070-01	G08	HQP017753	NM_000059	BRCA2
QG070-01	G09	HQP017616	NM_000636	SOD2
QG070-01	G10	HQP016418	NM_004327	BCR

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-01	G11	HQP016260	NM_004050	BCL2L2
QG070-01	G12	HQP011492	NM_002439	MSH3
QG070-01	H01	HGDC		
QG070-01	H02	HGDC		
QG070-01	H03	HQP006940	NM_002046	GAPDH
QG070-01	H04	HQP016381	NM_001101	ACTB
QG070-01	H05	HQP015171	NM_004048	B2M
QG070-01	H06	HQP006171	NM_012423	RPL13A
QG070-01	H07	HQP009026	NM_000194	HPRT1
QG070-01	H08	HQP054253	NR_003286	RN18S1
QG070-01	H09	RT		
QG070-01	H10	RT		
QG070-01	H11	PCR		
QG070-01	H12	PCR		
QG070-02	A01	HQP011491	NM_000251	MSH2
QG070-02	A02	HQP009849	NM_002227	JAK1
QG070-02	A03	HQP009678	NM_000584	IL8
QG070-02	A04	HQP009671	NM_000639	FASLG
QG070-02	A05	HQP009645	NM_000577	IL1RN
QG070-02	A06	HQP008934	NM_001160	APAF1
QG070-02	A07	HQP008620	NM_013371	IL19
QG070-02	A08	HQP007767	NM_002075	GNB3
QG070-02	A09	HQP005951	NM_001460	FMO2
QG070-02	A10	HQP004986	NM_000124	ERCC6
QG070-02	A11	HQP004984	NM_005236	ERCC4
QG070-02	A12	HQP004948	NM_000120	EPHX1
QG070-02	B01	HQP003808	NM_000767	CYP2B6
QG070-02	B02	HQP001475	NM_001005176	SP140
QG070-02	B03	HQP000370	NM_058195	CDKN2A
QG070-02	B04	HQP054690	NM_001099436	ULK3
QG070-02	B05	HQP022699	NM_006139	CD28
QG070-02	B06	HQP022202	NM_004213	SLC28A1
QG070-02	B07	HQP021906	NM_001239	CCNH
QG070-02	B08	HQP021834	NM_005092	TNFSF18
QG070-02	B09	HQP021553	NM_003842	TNFRSF10B
QG070-02	B10	HQP021526	NM_003824	FADD
QG070-02	B11	HQP021430	NM_005188	CBL
QG070-02	B12	HQP021408	NM_022845	CBFB
QG070-02	C01	HQP020384	NM_004347	CASP5
QG070-02	C02	HQP019982	NM_001017388	TLR10
QG070-02	C03	HQP018802	NM_001008540	CXCR4

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-02	C04	HQP018568	NM_021141	XRCC5
QG070-02	C05	HQP018113	NM_003263	TLR1
QG070-02	C06	HQP018044	NM_000660	TGFB1
QG070-02	C07	HQP017767	NM_003150	STAT3
QG070-02	C08	HQP017764	NM_007315	STAT1
QG070-02	C09	HQP017663	NM_003120	SPI1
QG070-02	C10	HQP016747	NM_003006	SELPLG
QG070-02	C11	HQP016746	NM_003005	SELP
QG070-02	C12	HQP016670	NM_001033886	CXCL12
QG070-02	D01	HQP016621	NM_002982	CCL2
QG070-02	D02	HQP016125	NM_002890	RASA1
QG070-02	D03	HQP016076	NM_002874	RAD23B
QG070-02	D04	HQP013422	NM_002690	POLB
QG070-02	D05	HQP013152	NM_002648	PIM1
QG070-02	D06	HQP012889	NM_002609	PDGFRB
QG070-02	D07	HQP012441	NM_016123	IRAK4
QG070-02	D08	HQP012078	NM_000607	ORM1
QG070-02	D09	HQP011866	NM_000625	NOS2A
QG070-02	D10	HQP011807	NM_003998	NFKB1
QG070-02	D11	HQP011774	NM_000267	NF1
QG070-02	D12	HQP011554	NM_000254	MTR
QG070-02	E01	HQP011308	NM_005373	MPL
QG070-02	E02	HQP010613	NM_002312	LIG4
QG070-02	E03	HQP010609	NM_000234	LIG1
QG070-02	E04	HQP010070	NM_002253	KDR
QG070-02	E05	HQP009718	NM_001562	IL18
QG070-02	E06	HQP009697	NM_002188	IL13
QG070-02	E07	HQP009694	NM_005535	IL12RB1
QG070-02	E08	HQP009692	NM_000882	IL12A
QG070-02	E09	HQP009681	NM_001557	IL8RB
QG070-02	E10	HQP009675	NM_175767	IL6ST
QG070-02	E11	HQP009670	NM_000600	IL6
QG070-02	E12	HQP009667	NM_000564	IL5RA
QG070-02	F01	HQP009642	NM_000877	IL1R1
QG070-02	F02	HQP009556	NM_000041	APOE
QG070-02	F03	HQP009472	NM_005534	IFNGR2
QG070-02	F04	HQP009469	NM_000416	IFNGR1
QG070-02	F05	HQP009399	NM_004509	SP110
QG070-02	F06	HQP009085	NM_002155	HSPA6
QG070-02	F07	HQP008690	NM_004131	GZMB
QG070-02	F08	HQP008565	NM_001039877	STRN4

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-02	F09	HQP008483	NM_000849	GSTM3
QG070-02	F10	HQP008402	NM_001020825	NR3C1
QG070-02	F11	HQP008362	NM_020469	ABO
QG070-02	F12	HQP007755	NM_000516	GNAS
QG070-02	G01	HQP006202	NM_012276	LILRA4
QG070-02	G02	HQP005058	NM_000130	F5
QG070-02	G03	HQP005022	NM_001987	ETV6
QG070-02	G04	HQP004991	NM_001014431	AKT1
QG070-02	G05	HQP003908	NM_000500	CYP21A2
QG070-02	G06	HQP003775	NM_000104	CYP1B1
QG070-02	G07	HQP003499	NM_001037631	CTLA4
QG070-02	G08	HQP003173	NM_000759	CSF3
QG070-02	G09	HQP003159	NM_000758	CSF2
QG070-02	G10	HQP003120	NM_001618	PARP1
QG070-02	G11	HQP002222	NM_005201	CCR8
QG070-02	G12	HQP002217	NM_001838	CCR7
QG070-02	H01	HGDC		
QG070-02	H02	HGDC		
QG070-02	H03	HQP006940	NM_002046	GAPDH
QG070-02	H04	HQP016381	NM_001101	ACTB
QG070-02	H05	HQP015171	NM_004048	B2M
QG070-02	H06	HQP006171	NM_012423	RPL13A
QG070-02	H07	HQP009026	NM_000194	HPRT1
QG070-02	H08	HQP054253	NR_003286	RN18S1
QG070-02	H09	RT		
QG070-02	H10	RT		
QG070-02	H11	PCR		
QG070-02	H12	PCR		
QG070-03	A01	HQP001228	NM_007027	TOPBP1
QG070-03	A02	HQP000459	NM_006092	NOD1
QG070-03	A03	HQP000415	NM_006068	TLR6
QG070-03	A04	HQP000145	NM_005732	RAD50
QG070-03	A05	HQP021498	NM_003808	TNFSF13
QG070-03	A06	HQP000648	NM_001805	CEBPE
QG070-03	A07	HQP054701	NM_001035	RYR2
QG070-03	A08	HQP054688	NM_032453	ZNF527
QG070-03	A09	HQP054326	NM_001099287	ICHTHYIN
QG070-03	A10	HQP054058	NM_005041	PRF1
QG070-03	A11	HQP054030	NM_130398	EXO1
QG070-03	A12	HQP023464	NM_012115	CASP8AP2
QG070-03	B01	HQP023409	NM_001790	CDC25C

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-03	B02	HQP023354	NM_001254	CDC6
QG070-03	B03	HQP023203	NM_001785	CDA
QG070-03	B04	HQP023168	NM_004356	CD81
QG070-03	B05	HQP023155	NM_000626	CD79B
QG070-03	B06	HQP023145	NM_021601	CD79A
QG070-03	B07	HQP023095	NM_012291	ESPL1
QG070-03	B08	HQP023074	NM_153029	N4BP1
QG070-03	B09	HQP023038	NM_014641	MDC1
QG070-03	B10	HQP023012	NM_006536	CLCA2
QG070-03	B11	HQP022998	NM_004917	KLK4
QG070-03	B12	HQP022955	NM_001250	CD40
QG070-03	C01	HQP022884	NM_004881	TP53I3
QG070-03	C02	HQP022490	NM_000591	CD14
QG070-03	C03	HQP022361	NM_004747	DLG5
QG070-03	C04	HQP022323	NM_004217	AURKB
QG070-03	C05	HQP022316	NM_000616	CD4
QG070-03	C06	HQP022241	NM_003856	IL1RL1
QG070-03	C07	HQP022197	NM_004212	SLC28A2
QG070-03	C08	HQP022170	NM_004708	PDCD5
QG070-03	C09	HQP021889	NM_003955	SOCS3
QG070-03	C10	HQP021725	NM_003921	BCL10
QG070-03	C11	HQP021636	NM_032454	STK19
QG070-03	C12	HQP021604	NM_003879	CFLAR
QG070-03	D01	HQP021602	NM_003877	SOCS2
QG070-03	D02	HQP021569	NM_003854	IL1RL2
QG070-03	D03	HQP021550	NM_003839	TNFRSF11A
QG070-03	D04	HQP021524	NM_003821	RIPK2
QG070-03	D05	HQP021522	NM_003820	TNFRSF14
QG070-03	D06	HQP021502	NM_003810	TNFSF10
QG070-03	D07	HQP021492	NM_003804	RIPK1
QG070-03	D08	HQP021484	NM_033274	ADAM19
QG070-03	D09	HQP021462	NM_003789	TRADD
QG070-03	D10	HQP021459	NM_003786	ABCC3
QG070-03	D11	HQP021399	NM_003745	SOCS1
QG070-03	D12	HQP021385	NM_003739	AKR1C3
QG070-03	E01	HQP021290	NM_003682	MADD
QG070-03	E02	HQP021260	NM_033035	TSLP
QG070-03	E03	HQP021130	NM_003632	CNTNAP1
QG070-03	E04	HQP020808	NM_003593	FOXN1
QG070-03	E05	HQP020707	NM_003579	RAD54L
QG070-03	E06	HQP020499	NM_032169	ACAD11

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-03	E07	HQP020427	NM_001226	CASP6
QG070-03	E08	HQP020332	NM_001225	CASP4
QG070-03	E09	HQP020071	NM_000070	CAPN3
QG070-03	E10	HQP020070	NM_001748	CAPN2
QG070-03	E11	HQP019941	NM_030931	DEFB126
QG070-03	E12	HQP019768	NM_130445	COL18A1
QG070-03	F01	HQP019462	NM_024945	RMI1
QG070-03	F02	HQP018854	NM_024051	C7orf24
QG070-03	F03	HQP018768	NM_001954	DDR1
QG070-03	F04	HQP018739	NM_006300	ZNF230
QG070-03	F05	HQP018682	NM_001018011	ZBTB16
QG070-03	F06	HQP018552	NM_000379	XDH
QG070-03	F07	HQP018411	NM_003357	SCGB1A1
QG070-03	F08	HQP018391	NM_000606	C8G
QG070-03	F09	HQP018380	NM_000066	C8B
QG070-03	F10	HQP018353	NM_000587	C7
QG070-03	F11	HQP018343	NM_000372	TYR
QG070-03	F12	HQP018326	NM_001736	C5AR1
QG070-03	G01	HQP018276	NM_000716	C4BPB
QG070-03	G02	HQP018259	NM_000715	C4BPA
QG070-03	G03	HQP018237	NM_004620	TRAF6
QG070-03	G04	HQP018236	NM_001033910	TRAF5
QG070-03	G05	HQP018234	NM_003300	TRAF3
QG070-03	G06	HQP018233	NM_021138	TRAF2
QG070-03	G07	HQP018226	NM_000063	C2
QG070-03	G08	HQP018176	NM_005657	TP53BP1
QG070-03	G09	HQP018172	NM_001067	TOP2A
QG070-03	G10	HQP018170	NM_172369	C1QC
QG070-03	G11	HQP018115	NM_003265	TLR3
QG070-03	G12	HQP018106	NM_001212	C1QBP
QG070-03	H01	HGDC		
QG070-03	H02	HGDC		
QG070-03	H03	HQP006940	NM_002046	GAPDH
QG070-03	H04	HQP016381	NM_001101	ACTB
QG070-03	H05	HQP015171	NM_004048	B2M
QG070-03	H06	HQP006171	NM_012423	RPL13A
QG070-03	H07	HQP009026	NM_000194	HPRT1
QG070-03	H08	HQP054253	NR_003286	RN18S1
QG070-03	H09	RT		
QG070-03	H10	RT		
QG070-03	H11	PCR		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-03	H12	PCR		
QG070-04	A01	HQP018022	NM_001211	BUB1B
QG070-04	A02	HQP017964	NM_003203	C2orf3
QG070-04	A03	HQP017895	NM_003189	TAL1
QG070-04	A04	HQP017816	NM_000456	SUOX
QG070-04	A05	HQP017811	NM_177536	SULT1A1
QG070-04	A06	HQP017770	NM_003151	STAT4
QG070-04	A07	HQP017766	NM_005419	STAT2
QG070-04	A08	HQP017702	NM_001725	BPI
QG070-04	A09	HQP017625	NM_005633	SOS1
QG070-04	A10	HQP017615	NM_000454	SOD1
QG070-04	A11	HQP017549	NM_030807	SLC2A11
QG070-04	A12	HQP017482	NM_194255	SLC19A1
QG070-04	B01	HQP017022	NM_001079533	CPEB1
QG070-04	B02	HQP016982	NM_001033855	DCLRE1C
QG070-04	B03	HQP016844	NM_000057	BLM
QG070-04	B04	HQP016763	NM_022127	SLC28A3
QG070-04	B05	HQP016744	NM_000450	SELE
QG070-04	B06	HQP016738	NM_003004	SECTM1
QG070-04	B07	HQP016644	NM_002991	CCL24
QG070-04	B08	HQP016639	NM_004591	CCL20
QG070-04	B09	HQP016637	NM_002988	CCL18
QG070-04	B10	HQP016630	NM_005408	CCL13
QG070-04	B11	HQP016629	NM_002986	CCL11
QG070-04	B12	HQP016626	NM_002985	CCL5
QG070-04	C01	HQP016620	NM_002981	CCL1
QG070-04	C02	HQP016581	NM_006919	SERPINB3
QG070-04	C03	HQP016565	NM_001710	CFB
QG070-04	C04	HQP016500	NM_001032295	SERPING1
QG070-04	C05	HQP016391	NM_002947	RPA3
QG070-04	C06	HQP016389	NM_002945	RPA1
QG070-04	C07	HQP016344	NM_001024808	BCL7A
QG070-04	C08	HQP016223	NM_002912	REV3L
QG070-04	C09	HQP016214	NM_006509	RELB
QG070-04	C10	HQP016213	NM_021975	RELA
QG070-04	C11	HQP016208	NM_002908	REL
QG070-04	C12	HQP016172	NM_005611	RBL2
QG070-04	D01	HQP016107	NM_002885	RAP1GAP
QG070-04	D02	HQP016064	NM_018890	RAC1
QG070-04	D03	HQP015947	NM_002853	RAD1
QG070-04	D04	HQP015917	NM_001188	BAK1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-04	D05	HQP015625	NM_020706	SFRS15
QG070-04	D06	HQP015565	NM_020661	AICDA
QG070-04	D07	HQP015537	NM_000953	PTGDR
QG070-04	D08	HQP015530	NM_000264	PTCH1
QG070-04	D09	HQP015248	NM_001080452	GPR108
QG070-04	D10	HQP015240	NM_020162	DHX33
QG070-04	D11	HQP015102	NM_001012965	KLK6
QG070-04	D12	HQP015096	NM_005046	KLK7
QG070-04	E01	HQP015061	NM_019619	PARD3
QG070-04	E02	HQP015023	NM_054027	ANKH
QG070-04	E03	HQP014732	NM_018416	FOXJ2
QG070-04	E04	HQP014462	NM_017509	KLK15
QG070-04	E05	HQP014292	NM_018315	FBXW7
QG070-04	E06	HQP013991	NM_017944	USP47
QG070-04	E07	HQP013596	NM_001184	ATR
QG070-04	E08	HQP013496	NM_019009	TOLLIP
QG070-04	E09	HQP013447	NM_017545	HAO1
QG070-04	E10	HQP013425	NM_006231	POLE
QG070-04	E11	HQP013388	NM_017442	TLR9
QG070-04	E12	HQP013133	NM_005025	SERPINI1
QG070-04	F01	HQP013071	NM_006212	PFKFB2
QG070-04	F02	HQP013015	NM_000442	PECAM1
QG070-04	F03	HQP012755	NM_001037872	REV1
QG070-04	F04	HQP012651	NM_016639	TNFRSF12A
QG070-04	F05	HQP012584	NM_016546	C1RL
QG070-04	F06	HQP012393	NM_016009	SH3GLB1
QG070-04	F07	HQP012335	NM_015895	GMNN
QG070-04	F08	HQP012166	NM_022047	DEF6
QG070-04	F09	HQP012155	NM_002575	SERPINB2
QG070-04	F10	HQP012154	NM_000602	SERPINE1
QG070-04	F11	HQP012079	NM_000608	ORM2
QG070-04	F12	HQP012073	NM_002552	ORC4L
QG070-04	G01	HQP012062	NM_001008505	OPRM1
QG070-04	G02	HQP011938	NM_001007792	NTRK1
QG070-04	G03	HQP011868	NM_000603	NOS3
QG070-04	G04	HQP011858	NM_000904	NQO2
QG070-04	G05	HQP011813	NM_004556	NFKBIE
QG070-04	G06	HQP011810	NM_020529	NFKBIA
QG070-04	G07	HQP011808	NM_001077493	NFKB2
QG070-04	G08	HQP011741	NM_006154	NEDD4
QG070-04	G09	HQP011728	NM_004550	NDUFS2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-04	G10	HQP011694	NM_000631	NCF4
QG070-04	G11	HQP011693	NM_000433	NCF2
QG070-04	G12	HQP011663	NM_005967	NAB2
QG070-04	H01	HGDC		
QG070-04	H02	HGDC		
QG070-04	H03	HQP006940	NM_002046	GAPDH
QG070-04	H04	HQP016381	NM_001101	ACTB
QG070-04	H05	HQP015171	NM_004048	B2M
QG070-04	H06	HQP006171	NM_012423	RPL13A
QG070-04	H07	HQP009026	NM_000194	HPRT1
QG070-04	H08	HQP054253	NR_003286	RN18S1
QG070-04	H09	RT		
QG070-04	H10	RT		
QG070-04	H11	PCR		
QG070-04	H12	PCR		
QG070-05	A01	HQP011620	NM_000488	SERPINC1
QG070-05	A02	HQP011616	NM_001040113	MYH11
QG070-05	A03	HQP011603	NM_002468	MYD88
QG070-05	A04	HQP011597	NM_002467	MYC
QG070-05	A05	HQP011325	NM_019899	ABCC1
QG070-05	A06	HQP011320	NM_005590	MRE11A
QG070-05	A07	HQP011268	NM_004995	MMP14
QG070-05	A08	HQP011264	NM_002425	MMP10
QG070-05	A09	HQP011263	NM_004994	MMP9
QG070-05	A10	HQP011262	NM_002424	MMP8
QG070-05	A11	HQP011219	NM_002415	MIF
QG070-05	A12	HQP011113	NM_002389	CD46
QG070-05	B01	HQP011109	NM_006739	MCM5
QG070-05	B02	HQP011108	NM_005914	MCM4
QG070-05	B03	HQP011097	NM_005912	MC4R
QG070-05	B04	HQP011085	NM_001025081	MBP
QG070-05	B05	HQP011073	NM_015844	MBD1
QG070-05	B06	HQP011024	NM_016835	MAPT
QG070-05	B07	HQP011004	NM_002372	MAN2A1
QG070-05	B08	HQP010996	NM_022438	MAL
QG070-05	B09	HQP010966	NM_005904	SMAD7
QG070-05	B10	HQP010932	NM_005582	CD180
QG070-05	B11	HQP010660	NM_005570	LMAN1
QG070-05	B12	HQP010612	NM_013975	LIG3
QG070-05	C01	HQP010584	NM_002303	LEPR
QG070-05	C02	HQP010560	NM_004139	LBP

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-05	C03	HQP010497	NM_000426	LAMA2
QG070-05	C04	HQP010104	NM_000892	KLKB1
QG070-05	C05	HQP010100	NM_002257	KLK1
QG070-05	C06	HQP009819	NM_000213	ITGB4
QG070-05	C07	HQP009805	NM_002208	ITGAE
QG070-05	C08	HQP009782	NM_002200	IRF5
QG070-05	C09	HQP009776	NM_001570	IRAK2
QG070-05	C10	HQP009745	NM_005538	INHBC
QG070-05	C11	HQP009716	NM_001561	TNFRSF9
QG070-05	C12	HQP009708	NM_000585	IL15
QG070-05	D01	HQP009696	NM_001559	IL12RB2
QG070-05	D02	HQP009693	NM_002187	IL12B
QG070-05	D03	HQP009687	NM_000628	IL10RB
QG070-05	D04	HQP009679	NM_000634	IL8RA
QG070-05	D05	HQP009672	NM_000565	IL6R
QG070-05	D06	HQP009666	NM_000879	IL5
QG070-05	D07	HQP009664	NM_000418	IL4R
QG070-05	D08	HQP009649	NM_000586	IL2
QG070-05	D09	HQP009639	NM_001556	IKBKB
QG070-05	D10	HQP009460	NM_000874	IFNAR2
QG070-05	D11	HQP009458	NM_000629	IFNAR1
QG070-05	D12	HQP009251	NM_002163	IRF8
QG070-05	E01	HQP009201	NM_001039132	ICAM4
QG070-05	E02	HQP009177	NM_176822	NLRP14
QG070-05	E03	HQP009139	NM_001643	APOA2
QG070-05	E04	HQP009091	NM_001167	BIRC4
QG070-05	E05	HQP009011	NM_014621	HOXD4
QG070-05	E06	HQP008846	NM_001607	ACAA1
QG070-05	E07	HQP008808	NM_021155	CD209
QG070-05	E08	HQP008801	NM_001010931	HGF
QG070-05	E09	HQP008799	NM_014593	CXXC1
QG070-05	E10	HQP008700	NM_005323	HIST1H1T
QG070-05	E11	HQP008554	NM_012092	ICOS
QG070-05	E12	HQP008498	NM_005513	GTF2E1
QG070-05	F01	HQP008490	NM_001513	GSTZ1
QG070-05	F02	HQP008476	NM_000846	GSTA2
QG070-05	F03	HQP008285	NM_002085	GPX4
QG070-05	F04	HQP008282	NM_002084	GPX3
QG070-05	F05	HQP008281	NM_002083	GPX2
QG070-05	F06	HQP008198	NM_173681	ATG9B
QG070-05	F07	HQP007821	NM_019844	SLCO1B3

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-05	F08	HQP007722	NM_014905	GLS
QG070-05	F09	HQP007689	NM_013274	POLL
QG070-05	F10	HQP007152	NM_032638	GATA2
QG070-05	F11	HQP007100	NM_015670	SENP3
QG070-05	F12	HQP006547	NM_001469	XRCC6
QG070-05	G01	HQP006427	NM_001039130	ALOX15B
QG070-05	G02	HQP006425	NM_001140	ALOX15
QG070-05	G03	HQP006396	NM_153289	DEFB119
QG070-05	G04	HQP006375	NM_001629	ALOX5AP
QG070-05	G05	HQP006359	NM_000698	ALOX5
QG070-05	G06	HQP006356	NM_000697	ALOX12
QG070-05	G07	HQP006320	NM_080671	KCNE4
QG070-05	G08	HQP006296	NM_015364	LY96
QG070-05	G09	HQP006249	NM_012381	ORC3L
QG070-05	G10	HQP006233	NM_012114	CASP14
QG070-05	G11	HQP006131	NM_014294	TRAM1
QG070-05	G12	HQP005959	NM_001002294	FMO3
QG070-05	H01	HGDC		
QG070-05	H02	HGDC		
QG070-05	H03	HQP006940	NM_002046	GAPDH
QG070-05	H04	HQP016381	NM_001101	ACTB
QG070-05	H05	HQP015171	NM_004048	B2M
QG070-05	H06	HQP006171	NM_012423	RPL13A
QG070-05	H07	HQP009026	NM_000194	HPRT1
QG070-05	H08	HQP054253	NR_003286	RN18S1
QG070-05	H09	RT		
QG070-05	H10	RT		
QG070-05	H11	PCR		
QG070-05	H12	PCR		
QG070-06	A01	HQP005589	NM_012072	CD93
QG070-06	A02	HQP005511	NM_001033053	NLRP1
QG070-06	A03	HQP005405	NM_002007	FGF4
QG070-06	A04	HQP005403	NM_002006	FGF2
QG070-06	A05	HQP005395	NM_002005	FES
QG070-06	A06	HQP005242	NM_002002	FCER2
QG070-06	A07	HQP005240	NM_004106	FCER1G
QG070-06	A08	HQP005238	NM_002001	FCER1A
QG070-06	A09	HQP005052	NM_000506	F2
QG070-06	A10	HQP004998	NM_000125	ESR1
QG070-06	A11	HQP004983	NM_000122	ERCC3
QG070-06	A12	HQP004950	NM_001979	EPHX2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-06	B01	HQP004934	NM_145007	NLRP11
QG070-06	B02	HQP004864	NM_001975	ENO2
QG070-06	B03	HQP004605	NM_005228	EGFR
QG070-06	B04	HQP004599	NM_001963	EGF
QG070-06	B05	HQP004557	NM_001955	EDN1
QG070-06	B06	HQP004526	NM_004091	E2F2
QG070-06	B07	HQP004426	NM_001381	DOK1
QG070-06	B08	HQP004411	NM_022552	DNMT3A
QG070-06	B09	HQP004309	NM_000791	DHFR
QG070-06	B10	HQP004242	NM_194320	ZNF169
QG070-06	B11	HQP004186	NM_021010	DEFA5
QG070-06	B12	HQP004182	NM_001925	DEFA4
QG070-06	C01	HQP004115	NM_000790	DDC
QG070-06	C02	HQP004081	NM_000789	ACE
QG070-06	C03	HQP004066	NM_000788	DCK
QG070-06	C04	HQP003855	NM_000779	CYP4B1
QG070-06	C05	HQP003826	NM_001017420	ESCO2
QG070-06	C06	HQP003813	NM_000772	CYP2C18
QG070-06	C07	HQP003810	NM_000770	CYP2C8
QG070-06	C08	HQP003539	NM_001904	CTNNB1
QG070-06	C09	HQP003420	NM_182919	TICAM1
QG070-06	C10	HQP003408	NM_134444	NLRP4
QG070-06	C11	HQP003393	NM_144685	HIPK4
QG070-06	C12	HQP003151	NM_172211	CSF1
QG070-06	D01	HQP003069	NM_139074	DEFB127
QG070-06	D02	HQP002873	NM_000098	CPT2
QG070-06	D03	HQP002869	NM_004377	CPT1B
QG070-06	D04	HQP002353	NM_176811	NLRP8
QG070-06	D05	HQP002209	NM_005508	CCR4
QG070-06	D06	HQP002198	NM_001295	CCR1
QG070-06	D07	HQP002181	NM_080867	SOCS4
QG070-06	D08	HQP001954	NM_054112	DEFB118
QG070-06	D09	HQP001862	NM_000082	ERCC8
QG070-06	D10	HQP001708	NM_001278	CHUK
QG070-06	D11	HQP001608	NM_007284	TWF2
QG070-06	D12	HQP001457	NM_004778	GPR44
QG070-06	E01	HQP001408	NM_007199	IRAK3
QG070-06	E02	HQP001399	NM_007195	POLI
QG070-06	E03	HQP001396	NM_001005735	CHEK2
QG070-06	E04	HQP001384	NM_003465	CHIT1
QG070-06	E05	HQP001351	NM_001008396	WDHD1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-06	E06	HQP001341	NM_007083	NUDT6
QG070-06	E07	HQP001288	NM_007051	FAF1
QG070-06	E08	HQP001287	NM_013441	DSCR1L2
QG070-06	E09	HQP000973	NM_006664	CCL27
QG070-06	E10	HQP000888	NM_006615	CAPN9
QG070-06	E11	HQP000657	NM_006407	ARL6IP5
QG070-06	E12	HQP000586	NM_006363	SEC23B
QG070-06	F01	HQP000421	NM_006072	CCL26
QG070-06	F02	HQP000274	NM_001259	CDK6
QG070-06	F03	HQP000245	NM_000075	CDK4
QG070-06	F04	HQP000020	NM_133484	TANK
QG070-06	F05	HQP054687	NM_005961	MUC6
QG070-06	F06	HQP018996	NM_021160	BAT5
QG070-06	F07	HQP018180	NM_005427	TP73
QG070-06	F08	HQP018159	NM_000491	C1QB
QG070-06	F09	HQP017618	NM_003102	SOD3
QG070-06	F10	HQP016651	NM_002995	XCL1
QG070-06	F11	HQP015394	NM_020375	C12orf5
QG070-06	F12	HQP015091	NM_001031849	MASP1
QG070-06	G01	HQP012805	NM_016463	CXXC5
QG070-06	G02	HQP011865	NM_000620	NOS1
QG070-06	G03	HQP011571	NM_152291	MUC7
QG070-06	G04	HQP010851	NM_006151	LPO
QG070-06	G05	HQP010101	NM_001002231	KLK2
QG070-06	G06	HQP009192	NM_002162	ICAM3
QG070-06	G07	HQP005588	NM_207341	ZP1
QG070-06	G08	HQP004414	NM_006892	DNMT3B
QG070-06	G09	HQP004189	NM_001926	DEFA6
QG070-06	G10	HQP002354	NM_153447	NLRP5
QG070-06	G11	HQP002352	NM_176810	NLRP13
QG070-06	G12	HQP001979	NM_001285	CLCA1
QG070-06	H01	HGDC		
QG070-06	H02	HGDC		
QG070-06	H03	HQP006940	NM_002046	GAPDH
QG070-06	H04	HQP016381	NM_001101	ACTB
QG070-06	H05	HQP015171	NM_004048	B2M
QG070-06	H06	HQP006171	NM_012423	RPL13A
QG070-06	H07	HQP009026	NM_000194	HPRT1
QG070-06	H08	HQP054253	NR_003286	RN18S1
QG070-06	H09	RT		
QG070-06	H10	RT		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG070-06	H11	PCR		
QG070-06	H12	PCR		