

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-01	A01	HQP011547	NM_005957	MTHFR
QG073-01	A02	HQP009685	NM_000572	IL10
QG073-01	A03	HQP018141	NM_000594	TNF
QG073-01	A04	HQP003772	NM_000499	CYP1A1
QG073-01	A05	HQP004317	NM_000903	NQO1
QG073-01	A06	HQP054047	BC008403	HLA-DRB1
QG073-01	A07	HQP011554	NM_000254	MTR
QG073-01	A08	HQP003817	NM_000773	CYP2E1
QG073-01	A09	HQP018562	NM_006297	XRCC1
QG073-01	A10	HQP018175	NM_000546	TP53
QG073-01	A11	HQP018342	NM_001071	TYMS
QG073-01	A12	HQP008849	NM_002116	HLA-A
QG073-01	B01	HQP004976	NM_000400	ERCC2
QG073-01	B02	HQP018213	NM_000367	TPMT
QG073-01	B03	HQP005280	NM_021642	FCGR2A
QG073-01	B04	HQP001136	NM_000015	NAT2
QG073-01	B05	HQP011309	NM_000250	MPO
QG073-01	B06	HQP009662	NM_000589	IL4
QG073-01	B07	HQP023467	NM_000662	NAT1
QG073-01	B08	HQP013100	NM_000927	ABCB1
QG073-01	B09	HQP011555	NM_002454	MTRR
QG073-01	B10	HQP016801	NM_022162	NOD2
QG073-01	B11	HQP009670	NM_000600	IL6
QG073-01	B12	HQP004948	NM_000120	EPHX1
QG073-01	C01	HQP003499	NM_001037631	CTLA4
QG073-01	C02	HQP018474	NM_000376	VDR
QG073-01	C03	HQP017616	NM_000636	SOD2
QG073-01	C04	HQP017482	NM_194255	SLC19A1
QG073-01	C05	HQP016224	NM_002913	RFC1
QG073-01	C06	HQP016212	NM_000657	BCL2
QG073-01	C07	HQP011687	NM_002485	NBN
QG073-01	C08	HQP011597	NM_002467	MYC
QG073-01	C09	HQP011135	NM_002392	MDM2
QG073-01	C10	HQP009664	NM_000418	IL4R
QG073-01	C11	HQP009645	NM_000577	IL1RN
QG073-01	C12	HQP009641	NM_000576	IL1B
QG073-01	D01	HQP003814	NM_000106	CYP2D6
QG073-01	D02	HQP003811	NM_000771	CYP2C9
QG073-01	D03	HQP021518	NM_000071	CBS
QG073-01	D04	HQP017753	NM_000059	BRCA2
QG073-01	D05	HQP014292	NM_018315	FBXW7

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-01	D06	HQP011868	NM_000603	NOS3
QG073-01	D07	HQP008483	NM_000849	GSTM3
QG073-01	D08	HQP003775	NM_000104	CYP1B1
QG073-01	D09	HQP000370	NM_058195	CDKN2A
QG073-01	D10	HQP012021	NM_002542	OGG1
QG073-01	D11	HQP004985	NM_000123	ERCC5
QG073-01	D12	HQP022955	NM_001250	CD40
QG073-01	E01	HQP018564	NM_005432	XRCC3
QG073-01	E02	HQP018475	NM_001025366	VEGFA
QG073-01	E03	HQP018044	NM_000660	TGFB1
QG073-01	E04	HQP016626	NM_002985	CCL5
QG073-01	E05	HQP016204	NM_053056	CCND1
QG073-01	E06	HQP011866	NM_000625	NOS2A
QG073-01	E07	HQP011807	NM_003998	NFKB1
QG073-01	E08	HQP010581	NM_000230	LEP
QG073-01	E09	HQP009693	NM_002187	IL12B
QG073-01	E10	HQP009692	NM_000882	IL12A
QG073-01	E11	HQP009678	NM_000584	IL8
QG073-01	E12	HQP009649	NM_000586	IL2
QG073-01	F01	HQP009640	NM_000575	IL1A
QG073-01	F02	HQP022490	NM_000591	CD14
QG073-01	F03	HQP021603	NM_003878	GGH
QG073-01	F04	HQP021385	NM_003739	AKR1C3
QG073-01	F05	HQP020515	NM_032199	ARID5B
QG073-01	F06	HQP020297	NM_004346	CASP3
QG073-01	F07	HQP018966	NM_001080124	CASP8
QG073-01	F08	HQP018563	NM_005431	XRCC2
QG073-01	F09	HQP018545	NM_000553	WRN
QG073-01	F10	HQP018466	NM_080682	VCAM1
QG073-01	F11	HQP018113	NM_003263	TLR1
QG073-01	F12	HQP017767	NM_003150	STAT3
QG073-01	G01	HQP017615	NM_000454	SOD1
QG073-01	G02	HQP016670	NM_001033886	CXCL12
QG073-01	G03	HQP015598	NM_000963	PTGS2
QG073-01	G04	HQP015535	NM_000314	PTEN
QG073-01	G05	HQP013388	NM_017442	TLR9
QG073-01	G06	HQP011491	NM_000251	MSH2
QG073-01	G07	HQP011320	NM_005590	MRE11A
QG073-01	G08	HQP010613	NM_002312	LIG4
QG073-01	G09	HQP010584	NM_002303	LEPR
QG073-01	G10	HQP009850	NM_004972	JAK2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-01	G11	HQP009781	NM_002460	IRF4
QG073-01	G12	HQP009697	NM_002188	IL13
QG073-01	H01	HGDC		
QG073-01	H02	HGDC		
QG073-01	H03	HQP006940	NM_002046	GAPDH
QG073-01	H04	HQP016381	NM_001101	ACTB
QG073-01	H05	HQP015171	NM_004048	B2M
QG073-01	H06	HQP006171	NM_012423	RPL13A
QG073-01	H07	HQP009026	NM_000194	HPRT1
QG073-01	H08	HQP054253	NR_003286	RN18S1
QG073-01	H09	RT		
QG073-01	H10	RT		
QG073-01	H11	PCR		
QG073-01	H12	PCR		
QG073-02	A01	HQP009666	NM_000879	IL5
QG073-02	A02	HQP009556	NM_000041	APOE
QG073-02	A03	HQP006228	NM_001018078	FPGS
QG073-02	A04	HQP005890	NM_004119	FLT3
QG073-02	A05	HQP005058	NM_000130	F5
QG073-02	A06	HQP004658	NM_001621	AHR
QG073-02	A07	HQP004309	NM_000791	DHFR
QG073-02	A08	HQP003908	NM_000500	CYP21A2
QG073-02	A09	HQP003888	NM_000102	CYP17A1
QG073-02	A10	HQP003841	NM_000777	CYP3A5
QG073-02	A11	HQP003691	NM_001337	CX3CR1
QG073-02	A12	HQP002210	NM_000579	CCR5
QG073-02	B01	HQP000963	NM_012190	ALDH1L1
QG073-02	B02	HQP000709	NM_006441	MTHFS
QG073-02	B03	HQP000405	NM_006066	AKR1A1
QG073-02	B04	HQP000145	NM_005732	RAD50
QG073-02	B05	HQP054686	NM_001123396	LOC729230
QG073-02	B06	HQP054058	NM_005041	PRF1
QG073-02	B07	HQP022870	NM_001775	CD38
QG073-02	B08	HQP022699	NM_006139	CD28
QG073-02	B09	HQP022527	NM_021950	MS4A1
QG073-02	B10	HQP021889	NM_003955	SOCS3
QG073-02	B11	HQP021492	NM_003804	RIPK1
QG073-02	B12	HQP020481	NM_033338	CASP7
QG073-02	C01	HQP020427	NM_001226	CASP6
QG073-02	C02	HQP020384	NM_004347	CASP5
QG073-02	C03	HQP020332	NM_001225	CASP4

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-02	C04	HQP020207	NM_001223	CASP1
QG073-02	C05	HQP019982	NM_001017388	TLR10
QG073-02	C06	HQP018565	NM_003401	XRCC4
QG073-02	C07	HQP018552	NM_000379	XDH
QG073-02	C08	HQP018380	NM_000066	C8B
QG073-02	C09	HQP018353	NM_000587	C7
QG073-02	C10	HQP018343	NM_000372	TYR
QG073-02	C11	HQP018326	NM_001736	C5AR1
QG073-02	C12	HQP018276	NM_000716	C4BPB
QG073-02	D01	HQP018259	NM_000715	C4BPA
QG073-02	D02	HQP018226	NM_000063	C2
QG073-02	D03	HQP018170	NM_172369	C1QC
QG073-02	D04	HQP018149	NM_001066	TNFRSF1B
QG073-02	D05	HQP017979	NM_000355	TCN2
QG073-02	D06	HQP017978	NM_001062	TCN1
QG073-02	D07	HQP017770	NM_003151	STAT4
QG073-02	D08	HQP017764	NM_007315	STAT1
QG073-02	D09	HQP016844	NM_000057	BLM
QG073-02	D10	HQP016744	NM_000450	SELE
QG073-02	D11	HQP016634	NM_001713	BHMT
QG073-02	D12	HQP016621	NM_002982	CCL2
QG073-02	E01	HQP016565	NM_001710	CFB
QG073-02	E02	HQP016500	NM_001032295	SERPING1
QG073-02	E03	HQP016260	NM_004050	BCL2L2
QG073-02	E04	HQP016064	NM_018890	RAC1
QG073-02	E05	HQP015917	NM_001188	BAK1
QG073-02	E06	HQP015538	NM_004322	BAD
QG073-02	E07	HQP015024	NM_000948	PRL
QG073-02	E08	HQP013473	NM_000446	PON1
QG073-02	E09	HQP012996	NM_016362	GHRL
QG073-02	E10	HQP012584	NM_016546	C1RL
QG073-02	E11	HQP012154	NM_000602	SERPINE1
QG073-02	E12	HQP011874	NM_000905	NPY
QG073-02	F01	HQP011812	NM_002503	NFKBIB
QG073-02	F02	HQP011810	NM_020529	NFKBIA
QG073-02	F03	HQP011694	NM_000631	NCF4
QG073-02	F04	HQP011693	NM_000433	NCF2
QG073-02	F05	HQP011603	NM_002468	MYD88
QG073-02	F06	HQP011256	NM_004530	MMP2
QG073-02	F07	HQP011219	NM_002415	MIF
QG073-02	F08	HQP011113	NM_002389	CD46

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-02	F09	HQP010133	NM_004985	KRAS
QG073-02	F10	HQP010095	NM_013289	KIR3DL1
QG073-02	F11	HQP010092	NM_012313	KIR2DS3
QG073-02	F12	HQP010086	NM_015868	KIR2DL3
QG073-02	G01	HQP010083	NM_014218	KIR2DL1
QG073-02	G02	HQP009851	NM_000215	JAK3
QG073-02	G03	HQP009708	NM_000585	IL15
QG073-02	G04	HQP009681	NM_001557	IL8RB
QG073-02	G05	HQP009677	NM_002185	IL7R
QG073-02	G06	HQP009472	NM_005534	IFNGR2
QG073-02	G07	HQP009139	NM_001643	APOA2
QG073-02	G08	HQP008402	NM_001020825	NR3C1
QG073-02	G09	HQP008285	NM_002085	GPX4
QG073-02	G10	HQP008198	NM_173681	ATG9B
QG073-02	G11	HQP005589	NM_012072	CD93
QG073-02	G12	HQP005242	NM_002002	FCER2
QG073-02	H01	HGDC		
QG073-02	H02	HGDC		
QG073-02	H03	HQP006940	NM_002046	GAPDH
QG073-02	H04	HQP016381	NM_001101	ACTB
QG073-02	H05	HQP015171	NM_004048	B2M
QG073-02	H06	HQP006171	NM_012423	RPL13A
QG073-02	H07	HQP009026	NM_000194	HPRT1
QG073-02	H08	HQP054253	NR_003286	RN18S1
QG073-02	H09	RT		
QG073-02	H10	RT		
QG073-02	H11	PCR		
QG073-02	H12	PCR		
QG073-03	A01	HQP005052	NM_000506	F2
QG073-03	A02	HQP004998	NM_000125	ESR1
QG073-03	A03	HQP004986	NM_000124	ERCC6
QG073-03	A04	HQP004975	NM_202001	ERCC1
QG073-03	A05	HQP004557	NM_001955	EDN1
QG073-03	A06	HQP003808	NM_000767	CYP2B6
QG073-03	A07	HQP003791	NM_000024	ADRB2
QG073-03	A08	HQP003120	NM_001618	PARP1
QG073-03	A09	HQP002671	NM_000754	COMT
QG073-03	A10	HQP000415	NM_006068	TLR6
QG073-03	A11	HQP018159	NM_000491	C1QB
QG073-03	A12	HQP017618	NM_003102	SOD3
QG073-03	B01	HQP011865	NM_000620	NOS1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-03	B02	HQP000027	NM_020396	BCL2L10
QG073-03	B03	HQP054688	NM_032453	ZNF527
QG073-03	B04	HQP054326	NM_001099287	ICHTHYIN
QG073-03	B05	HQP054057	BC071181	TGFBR1
QG073-03	B06	HQP053998	NM_004873	BAG5
QG073-03	B07	HQP053959	NM_001040	SHBG
QG073-03	B08	HQP023464	NM_012115	CASP8AP2
QG073-03	B09	HQP023414	NM_005847	SLC23A1
QG073-03	B10	HQP023354	NM_001254	CDC6
QG073-03	B11	HQP023203	NM_001785	CDA
QG073-03	B12	HQP023186	NM_014739	BCLAF1
QG073-03	C01	HQP023095	NM_012291	ESPL1
QG073-03	C02	HQP023012	NM_006536	CLCA2
QG073-03	C03	HQP022998	NM_004917	KLK4
QG073-03	C04	HQP022884	NM_004881	TP53I3
QG073-03	C05	HQP022872	NM_004281	BAG3
QG073-03	C06	HQP022764	NM_004832	GSTO1
QG073-03	C07	HQP022722	NM_005191	CD80
QG073-03	C08	HQP022625	NM_004797	ADIPOQ
QG073-03	C09	HQP022361	NM_004747	DLG5
QG073-03	C10	HQP022337	NM_014207	CD5
QG073-03	C11	HQP021834	NM_005092	TNFSF18
QG073-03	C12	HQP021737	NM_003927	MBD2
QG073-03	D01	HQP021725	NM_003921	BCL10
QG073-03	D02	HQP021636	NM_032454	STK19
QG073-03	D03	HQP021604	NM_003879	CFLAR
QG073-03	D04	HQP021602	NM_003877	SOCS2
QG073-03	D05	HQP021557	NM_003844	TNFRSF10A
QG073-03	D06	HQP021524	NM_003821	RIPK2
QG073-03	D07	HQP021522	NM_003820	TNFRSF14
QG073-03	D08	HQP021484	NM_033274	ADAM19
QG073-03	D09	HQP021462	NM_003789	TRADD
QG073-03	D10	HQP021420	NM_001756	SERPINA6
QG073-03	D11	HQP021399	NM_003745	SOCS1
QG073-03	D12	HQP021347	NM_001754	RUNX1
QG073-03	E01	HQP021260	NM_033035	TSLP
QG073-03	E02	HQP021130	NM_003632	CNTNAP1
QG073-03	E03	HQP021129	NM_003631	PARG
QG073-03	E04	HQP020808	NM_003593	FOXN1
QG073-03	E05	HQP020707	NM_003579	RAD54L
QG073-03	E06	HQP020499	NM_032169	ACAD11

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-03	E07	HQP020390	NM_032016	STARD3NL
QG073-03	E08	HQP020245	NM_001040668	BCL2L12
QG073-03	E09	HQP020040	NM_006534	NCOA3
QG073-03	E10	HQP019941	NM_030931	DEFB126
QG073-03	E11	HQP019873	NM_030787	CFHR5
QG073-03	E12	HQP019021	NM_030766	BCL2L14
QG073-03	F01	HQP018983	NM_004639	BAT3
QG073-03	F02	HQP018854	NM_024051	C7orf24
QG073-03	F03	HQP018802	NM_001008540	CXCR4
QG073-03	F04	HQP018768	NM_001954	DDR1
QG073-03	F05	HQP018760	NM_003463	PTP4A1
QG073-03	F06	HQP018556	NM_004628	XPC
QG073-03	F07	HQP018431	NM_007121	NR1H2
QG073-03	F08	HQP018411	NM_003357	SCGB1A1
QG073-03	F09	HQP018391	NM_000606	C8G
QG073-03	F10	HQP018340	NM_003331	TYK2
QG073-03	F11	HQP018237	NM_004620	TRAF6
QG073-03	F12	HQP018236	NM_001033910	TRAF5
QG073-03	G01	HQP018233	NM_021138	TRAF2
QG073-03	G02	HQP018172	NM_001067	TOP2A
QG073-03	G03	HQP018118	NM_003268	TLR5
QG073-03	G04	HQP018115	NM_003265	TLR3
QG073-03	G05	HQP018106	NM_001212	C1QBP
QG073-03	G06	HQP017866	NM_003183	ADAM17
QG073-03	G07	HQP017766	NM_005419	STAT2
QG073-03	G08	HQP017762	NM_000349	STAR
QG073-03	G09	HQP017702	NM_001725	BPI
QG073-03	G10	HQP017697	NM_001047	SRD5A1
QG073-03	G11	HQP017619	NM_004052	BNIP3
QG073-03	G12	HQP017549	NM_030807	SLC2A11
QG073-03	H01	HGDC		
QG073-03	H02	HGDC		
QG073-03	H03	HQP006940	NM_002046	GAPDH
QG073-03	H04	HQP016381	NM_001101	ACTB
QG073-03	H05	HQP015171	NM_004048	B2M
QG073-03	H06	HQP006171	NM_012423	RPL13A
QG073-03	H07	HQP009026	NM_000194	HPRT1
QG073-03	H08	HQP054253	NR_003286	RN18S1
QG073-03	H09	RT		
QG073-03	H10	RT		
QG073-03	H11	PCR		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-03	H12	PCR		
QG073-04	A01	HQP017469	NM_022444	SLC13A1
QG073-04	A02	HQP017117	NM_022735	ACBD3
QG073-04	A03	HQP016747	NM_003006	SELPLG
QG073-04	A04	HQP016746	NM_003005	SELP
QG073-04	A05	HQP016738	NM_003004	SECTM1
QG073-04	A06	HQP016644	NM_002991	CCL24
QG073-04	A07	HQP016639	NM_004591	CCL20
QG073-04	A08	HQP016637	NM_002988	CCL18
QG073-04	A09	HQP016630	NM_005408	CCL13
QG073-04	A10	HQP016629	NM_002986	CCL11
QG073-04	A11	HQP016620	NM_002981	CCL1
QG073-04	A12	HQP016581	NM_006919	SERPINB3
QG073-04	B01	HQP016344	NM_001024808	BCL7A
QG073-04	B02	HQP016208	NM_002908	REL
QG073-04	B03	HQP016131	NM_000321	RB1
QG073-04	B04	HQP016125	NM_002890	RASA1
QG073-04	B05	HQP016076	NM_002874	RAD23B
QG073-04	B06	HQP015828	NM_002827	PTPN1
QG073-04	B07	HQP015625	NM_020706	SFRS15
QG073-04	B08	HQP015596	NM_000962	PTGS1
QG073-04	B09	HQP015565	NM_020661	AICDA
QG073-04	B10	HQP015537	NM_000953	PTGDR
QG073-04	B11	HQP015248	NM_001080452	GPR108
QG073-04	B12	HQP015240	NM_020162	DHX33
QG073-04	C01	HQP015102	NM_001012965	KLK6
QG073-04	C02	HQP015096	NM_005046	KLK7
QG073-04	C03	HQP015061	NM_019619	PARD3
QG073-04	C04	HQP014732	NM_018416	FOXJ2
QG073-04	C05	HQP014462	NM_017509	KLK15
QG073-04	C06	HQP013991	NM_017944	USP47
QG073-04	C07	HQP013957	NM_001611	ACP5
QG073-04	C08	HQP013696	NM_017628	KIAA1546
QG073-04	C09	HQP013633	NM_005037	PPARG
QG073-04	C10	HQP013496	NM_019009	TOLLIP
QG073-04	C11	HQP013467	NM_000939	POMC
QG073-04	C12	HQP013422	NM_002690	POLB
QG073-04	D01	HQP013152	NM_002648	PIM1
QG073-04	D02	HQP013150	NM_006218	PIK3CA
QG073-04	D03	HQP013133	NM_005025	SERPINI1
QG073-04	D04	HQP013099	NM_000926	PGR

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-04	D05	HQP013071	NM_006212	PFKFB2
QG073-04	D06	HQP012441	NM_016123	IRAK4
QG073-04	D07	HQP012436	NM_001040443	PHF11
QG073-04	D08	HQP012212	NM_016734	PAX5
QG073-04	D09	HQP012211	NM_006193	PAX4
QG073-04	D10	HQP012166	NM_022047	DEF6
QG073-04	D11	HQP012155	NM_002575	SERPINB2
QG073-04	D12	HQP012008	NM_000275	OCA2
QG073-04	E01	HQP011914	NM_002524	NRAS
QG073-04	E02	HQP011885	NM_002518	NPAS2
QG073-04	E03	HQP011860	NM_006169	NNMT
QG073-04	E04	HQP011728	NM_004550	NDUFS2
QG073-04	E05	HQP011663	NM_005967	NAB2
QG073-04	E06	HQP011620	NM_000488	SERPINC1
QG073-04	E07	HQP011540	NM_002451	MTAP
QG073-04	E08	HQP011325	NM_019899	ABCC1
QG073-04	E09	HQP011264	NM_002425	MMP10
QG073-04	E10	HQP011263	NM_004994	MMP9
QG073-04	E11	HQP011257	NM_002422	MMP3
QG073-04	E12	HQP011255	NM_002421	MMP1
QG073-04	F01	HQP011235	NM_000249	MLH1
QG073-04	F02	HQP011223	NM_000248	MITF
QG073-04	F03	HQP011097	NM_005912	MC4R
QG073-04	F04	HQP011085	NM_001025081	MBP
QG073-04	F05	HQP010996	NM_022438	MAL
QG073-04	F06	HQP010932	NM_005582	CD180
QG073-04	F07	HQP010878	NM_002335	LRP5
QG073-04	F08	HQP010847	NM_000237	LPL
QG073-04	F09	HQP010660	NM_005570	LMAN1
QG073-04	F10	HQP010626	NM_000236	LIPC
QG073-04	F11	HQP010612	NM_013975	LIG3
QG073-04	F12	HQP010609	NM_000234	LIG1
QG073-04	G01	HQP010560	NM_004139	LBP
QG073-04	G02	HQP010497	NM_000426	LAMA2
QG073-04	G03	HQP010104	NM_000892	KLKB1
QG073-04	G04	HQP010100	NM_002257	KLK1
QG073-04	G05	HQP009849	NM_002227	JAK1
QG073-04	G06	HQP009837	NM_033453	ITPA
QG073-04	G07	HQP009801	NM_000044	AR
QG073-04	G08	HQP009776	NM_001570	IRAK2
QG073-04	G09	HQP009745	NM_005538	INHBC

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-04	G10	HQP009718	NM_001562	IL18
QG073-04	G11	HQP009710	NM_002189	IL15RA
QG073-04	G12	HQP009696	NM_001559	IL12RB2
QG073-04	H01	HGDC		
QG073-04	H02	HGDC		
QG073-04	H03	HQP006940	NM_002046	GAPDH
QG073-04	H04	HQP016381	NM_001101	ACTB
QG073-04	H05	HQP015171	NM_004048	B2M
QG073-04	H06	HQP006171	NM_012423	RPL13A
QG073-04	H07	HQP009026	NM_000194	HPRT1
QG073-04	H08	HQP054253	NR_003286	RN18S1
QG073-04	H09	RT		
QG073-04	H10	RT		
QG073-04	H11	PCR		
QG073-04	H12	PCR		
QG073-05	A01	HQP009694	NM_005535	IL12RB1
QG073-05	A02	HQP009679	NM_000634	IL8RA
QG073-05	A03	HQP009650	NM_000417	IL2RA
QG073-05	A04	HQP009639	NM_001556	IKBKB
QG073-05	A05	HQP009544	NM_000598	IGFBP3
QG073-05	A06	HQP009539	NM_000596	IGFBP1
QG073-05	A07	HQP009529	NM_000612	IGF2
QG073-05	A08	HQP009523	NM_000875	IGF1R
QG073-05	A09	HQP009518	NM_000618	IGF1
QG073-05	A10	HQP009469	NM_000416	IFNGR1
QG073-05	A11	HQP009351	NM_005896	IDH1
QG073-05	A12	HQP009218	NM_000384	APOB
QG073-05	B01	HQP009201	NM_001039132	ICAM4
QG073-05	B02	HQP009171	NM_000482	APOA4
QG073-05	B03	HQP009085	NM_002155	HSPA6
QG073-05	B04	HQP009066	NM_002153	HSD17B2
QG073-05	B05	HQP009064	NM_000413	HSD17B1
QG073-05	B06	HQP009061	NM_001641	APEX1
QG073-05	B07	HQP009060	NM_000198	HSD3B2
QG073-05	B08	HQP009059	NM_000862	HSD3B1
QG073-05	B09	HQP009019	NM_005143	HP
QG073-05	B10	HQP008893	NM_005518	HMGCS2
QG073-05	B11	HQP008892	NM_002130	HMGCS1
QG073-05	B12	HQP008846	NM_001607	ACAA1
QG073-05	C01	HQP008808	NM_021155	CD209
QG073-05	C02	HQP008801	NM_001010931	HGF

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-05	C03	HQP008620	NM_013371	IL19
QG073-05	C04	HQP008554	NM_012092	ICOS
QG073-05	C05	HQP008498	NM_005513	GTF2E1
QG073-05	C06	HQP008302	NM_000827	GRIA1
QG073-05	C07	HQP008282	NM_002084	GPX3
QG073-05	C08	HQP008281	NM_002083	GPX2
QG073-05	C09	HQP007821	NM_019844	SLCO1B3
QG073-05	C10	HQP007722	NM_014905	GLS
QG073-05	C11	HQP007385	NM_000515	GH1
QG073-05	C12	HQP007100	NM_015670	SENP3
QG073-05	D01	HQP006427	NM_001039130	ALOX15B
QG073-05	D02	HQP006425	NM_001140	ALOX15
QG073-05	D03	HQP006396	NM_153289	DEFB119
QG073-05	D04	HQP006375	NM_001629	ALOX5AP
QG073-05	D05	HQP006359	NM_000698	ALOX5
QG073-05	D06	HQP006356	NM_000697	ALOX12
QG073-05	D07	HQP006350	NM_015367	BCL2L13
QG073-05	D08	HQP006296	NM_015364	LY96
QG073-05	D09	HQP006245	NM_014317	PDSS1
QG073-05	D10	HQP006233	NM_012114	CASP14
QG073-05	D11	HQP006202	NM_012276	LILRA4
QG073-05	D12	HQP006131	NM_014294	TRAM1
QG073-05	E01	HQP006080	NM_012238	SIRT1
QG073-05	E02	HQP005879	NM_002019	FLT1
QG073-05	E03	HQP005403	NM_002006	FGF2
QG073-05	E04	HQP005369	NM_001010873	BZRPL1
QG073-05	E05	HQP005357	NM_004462	FDFT1
QG073-05	E06	HQP005292	NM_001002275	FCGR2B
QG073-05	E07	HQP005240	NM_004106	FCER1G
QG073-05	E08	HQP005239	NM_000139	MS4A2
QG073-05	E09	HQP005238	NM_002001	FCER1A
QG073-05	E10	HQP005022	NM_001987	ETV6
QG073-05	E11	HQP005002	NM_001437	ESR2
QG073-05	E12	HQP004991	NM_001014431	AKT1
QG073-05	F01	HQP004984	NM_005236	ERCC4
QG073-05	F02	HQP004983	NM_000122	ERCC3
QG073-05	F03	HQP004950	NM_001979	EPHX2
QG073-05	F04	HQP004724	NM_001623	AIF1
QG073-05	F05	HQP004451	NM_000798	DRD5
QG073-05	F06	HQP004446	NM_000795	DRD2
QG073-05	F07	HQP004242	NM_194320	ZNF169

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG073-05	F08	HQP004198	NM_001928	CFD
QG073-05	F09	HQP004186	NM_021010	DEFA5
QG073-05	F10	HQP004182	NM_001925	DEFA4
QG073-05	F11	HQP004081	NM_000789	ACE
QG073-05	F12	HQP004066	NM_000788	DCK
QG073-05	G01	HQP003992	NM_001350	DAXX
QG073-05	G02	HQP003904	NM_000103	CYP19A1
QG073-05	G03	HQP003812	NM_000025	ADRB3
QG073-05	G04	HQP003809	NM_000769	CYP2C19
QG073-05	G05	HQP003539	NM_001904	CTNNB1
QG073-05	G06	HQP003420	NM_182919	TICAM1
QG073-05	G07	HQP003393	NM_144685	HIPK4
QG073-05	G08	HQP003173	NM_000759	CSF3
QG073-05	G09	HQP003084	NM_021117	CRY2
QG073-05	G10	HQP003069	NM_139074	DEFB127
QG073-05	G11	HQP002881	NM_000651	CR1
QG073-05	G12	HQP002873	NM_000098	CPT2
QG073-05	H01	HGDC		
QG073-05	H02	HGDC		
QG073-05	H03	HQP006940	NM_002046	GAPDH
QG073-05	H04	HQP016381	NM_001101	ACTB
QG073-05	H05	HQP015171	NM_004048	B2M
QG073-05	H06	HQP006171	NM_012423	RPL13A
QG073-05	H07	HQP009026	NM_000194	HPRT1
QG073-05	H08	HQP054253	NR_003286	RN18S1
QG073-05	H09	RT		
QG073-05	H10	RT		
QG073-05	H11	PCR		
QG073-05	H12	PCR		