

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG031-01	A01	HQP053905	NM_001079	ZAP70
QG031-01	A02	HQP007848	NM_001024644	XCR1
QG031-01	A03	HQP018466	NM_080682	VCAM1
QG031-01	A04	HQP018461	NM_005428	VAV1
QG031-01	A05	HQP018237	NM_004620	TRAF6
QG031-01	A06	HQP018234	NM_003300	TRAF3
QG031-01	A07	HQP018233	NM_021138	TRAF2
QG031-01	A08	HQP018232	NM_005658	TRAF1
QG031-01	A09	HQP013496	NM_019009	TOLLIP
QG031-01	A10	HQP000821	NM_006573	TNFSF13B
QG031-01	A11	HQP021502	NM_003810	TNFSF10
QG031-01	A12	HQP018149	NM_001066	TNFRSF1B
QG031-01	B01	HQP016367	NM_001192	TNFRSF17
QG031-01	B02	HQP006148	NM_012452	TNFRSF13B
QG031-01	B03	HQP021550	NM_003839	TNFRSF11A
QG031-01	B04	HQP018147	NM_007115	TNFAIP6
QG031-01	B05	HQP018141	NM_000594	TNF
QG031-01	B06	HQP013388	NM_017442	TLR9
QG031-01	B07	HQP012591	NM_016562	TLR7
QG031-01	B08	HQP018116	NM_138554	TLR4
QG031-01	B09	HQP018114	NM_003264	TLR2
QG031-01	B10	HQP001674	NM_001039661	TIRAP
QG031-01	B11	HQP018083	NM_006288	THY1
QG031-01	B12	HQP018051	NM_004612	TGFBR1
QG031-01	C01	HQP018044	NM_000660	TGFB1
QG031-01	C02	HQP018043	NM_003236	TGFA
QG031-01	C03	HQP018041	NM_003234	TFRC
QG031-01	C04	HQP017958	NM_003202	TCF7
QG031-01	C05	HQP008682	NM_013351	TBX21
QG031-01	C06	HQP017902	NM_003190	TAPBP
QG031-01	C07	HQP017899	NM_000593	TAP1
QG031-01	C08	HQP017895	NM_003189	TAL1
QG031-01	C09	HQP017845	NM_003177	SYK
QG031-01	C10	HQP017775	NM_003153	STAT6
QG031-01	C11	HQP017774	NM_012448	STAT5B
QG031-01	C12	HQP017771	NM_003152	STAT5A
QG031-01	D01	HQP017767	NM_003150	STAT3
QG031-01	D02	HQP017764	NM_007315	STAT1
QG031-01	D03	HQP017696	NM_005417	SRC
QG031-01	D04	HQP017673	NM_000582	SPP1
QG031-01	D05	HQP017615	NM_000454	SOD1

QG031-01	D06	HQP021889	NM_003955	SOCS3
QG031-01	D07	HQP021399	NM_003745	SOCS1
QG031-01	D08	HQP017303	NM_003037	SLAMF1
QG031-01	D09	HQP016150	NM_021805	SIGIRR
QG031-01	D10	HQP016747	NM_003006	SELPLG
QG031-01	D11	HQP016745	NM_000655	SELL
QG031-01	D12	HQP016744	NM_000450	SELE
QG031-01	E01	HQP016378	NM_001001523	RORC
QG031-01	E02	HQP016213	NM_021975	RELA
QG031-01	E03	HQP016208	NM_002908	REL
QG031-01	E04	HQP009574	NM_005349	RBPJ
QG031-01	E05	HQP016089	NM_000448	RAG1
QG031-01	E06	HQP015908	NM_002838	PTPRC
QG031-01	E07	HQP007113	NM_012411	PTPN22
QG031-01	E08	HQP015598	NM_000963	PTGS2
QG031-01	E09	HQP015524	NM_000952	PTAFR
QG031-01	E10	HQP015311	NM_002800	PSMB9
QG031-01	E11	HQP015291	NM_004159	PSMB8
QG031-01	E12	HQP021305	NM_003690	PRKRA
QG031-01	F01	HQP014731	NM_006254	PRKCD
QG031-01	F02	HQP016740	NM_001198	PRDM1
QG031-01	F03	HQP013633	NM_005037	PPARG
QG031-01	F04	HQP013205	NM_001005376	PLAUR
QG031-01	F05	HQP013204	NM_002658	PLAU
QG031-01	F06	HQP013015	NM_000442	PECAM1
QG031-01	F07	HQP012889	NM_002609	PDGFRB
QG031-01	F08	HQP012856	NM_002608	PDGFB
QG031-01	F09	HQP012662	NM_005018	PDCD1
QG031-01	F10	HQP012212	NM_016734	PAX5
QG031-01	F11	HQP011988	NM_001032731	OAS2
QG031-01	F12	HQP011873	NM_017617	NOTCH1
QG031-01	G01	HQP011866	NM_000625	NOS2A
QG031-01	G02	HQP016801	NM_022162	NOD2
QG031-01	G03	HQP014556	NM_017852	NLRP2
QG031-01	G04	HQP011810	NM_020529	NFKBIA
QG031-01	G05	HQP011808	NM_001077493	NFKB2
QG031-01	G06	HQP011807	NM_003998	NFKB1
QG031-01	G07	HQP011789	NM_012340	NFATC2
QG031-01	G08	HQP011784	NM_006162	NFATC1
QG031-01	G09	HQP022750	NM_004829	NCR1
QG031-01	G10	HQP011694	NM_000631	NCF4
QG031-01	G11	HQP011603	NM_002468	MYD88
QG031-01	G12	HQP011582	NM_002462	MX1

QG031-01	H01	HGDC		
QG031-01	H02	HGDC		
QG031-01	H03	HQP006940	NM_002046	GAPDH
QG031-01	H04	HQP016381	NM_001101	ACTB
QG031-01	H05	HQP015171	NM_004048	B2M
QG031-01	H06	HQP006171	NM_012423	RPL13A
QG031-01	H07	HQP009026	NM_000194	HPRT1
QG031-01	H08	HQP054253	NR_003286	RN18S1
QG031-01	H09	RT		
QG031-01	H10	RT		
QG031-01	H11	PCR		
QG031-01	H12	PCR		
QG031-02	A01	HQP011547	NM_005957	MTHFR
QG031-02	A02	HQP022527	NM_021950	MS4A1
QG031-02	A03	HQP011263	NM_004994	MMP9
QG031-02	A04	HQP011256	NM_004530	MMP2
QG031-02	A05	HQP011267	NM_002427	MMP13
QG031-02	A06	HQP011219	NM_002415	MIF
QG031-02	A07	HQP011154	NM_000243	MEFV
QG031-02	A08	HQP011077	NM_000242	MBL2
QG031-02	A09	HQP000882	NM_006610	MASP2
QG031-02	A10	HQP014886	NM_002750	MAPK8
QG031-02	A11	HQP014854	NM_001040056	MAPK3
QG031-02	A12	HQP003133	NM_001315	MAPK14
QG031-02	B01	HQP014848	NM_002745	MAPK1
QG031-02	B02	HQP017891	NM_003188	MAP3K7
QG031-02	B03	HQP001017	NM_006785	MALT1
QG031-02	B04	HQP006296	NM_015364	LY96
QG031-02	B05	HQP010933	NM_002349	LY75
QG031-02	B06	HQP010919	NM_002343	LTF
QG031-02	B07	HQP010915	NM_002342	LTBR
QG031-02	B08	HQP010907	NM_000595	LTA
QG031-02	B09	HQP010608	NM_002310	LIFR
QG031-02	B10	HQP010607	NM_002309	LIF
QG031-02	B11	HQP010582	NM_001003679	LEPR
QG031-02	B12	HQP010581	NM_000230	LEP
QG031-02	C01	HQP012480	NM_016269	LEF1
QG031-02	C02	HQP010569	NM_005565	LCP2
QG031-02	C03	HQP010565	NM_001042771	LCK
QG031-02	C04	HQP010560	NM_004139	LBP
QG031-02	C05	HQP007461	NM_001014987	LAT
QG031-02	C06	HQP010136	NM_006121	KRT1
QG031-02	C07	HQP005584	NM_007360	KLRK1

QG031-02	C08	HQP010112	NM_002262	KLRD1
QG031-02	C09	HQP000884	NM_006611	KLRA1
QG031-02	C10	HQP011204	NM_000899	KITLG
QG031-02	C11	HQP010099	NM_000222	KIT
QG031-02	C12	HQP009851	NM_000215	JAK3
QG031-02	D01	HQP009850	NM_004972	JAK2
QG031-02	D02	HQP009849	NM_002227	JAK1
QG031-02	D03	HQP009815	NM_000211	ITGB2
QG031-02	D04	HQP009810	NM_002211	ITGB1
QG031-02	D05	HQP009809	NM_000887	ITGAX
QG031-02	D06	HQP009807	NM_000632	ITGAM
QG031-02	D07	HQP009806	NM_002209	ITGAL
QG031-02	D08	HQP009798	NM_000885	ITGA4
QG031-02	D09	HQP009795	NM_000419	ITGA2B
QG031-02	D10	HQP009793	NM_181501	ITGA1
QG031-02	D11	HQP009251	NM_002163	IRF8
QG031-02	D12	HQP009780	NM_001571	IRF3
QG031-02	E01	HQP009778	NM_002198	IRF1
QG031-02	E02	HQP009776	NM_001570	IRAK2
QG031-02	E03	HQP009771	NM_001025242	IRAK1
QG031-02	E04	HQP009754	NM_001017915	INPP5D
QG031-02	E05	HQP009736	NM_002164	INDO
QG031-02	E06	HQP009681	NM_001557	IL8RB
QG031-02	E07	HQP009679	NM_000634	IL8RA
QG031-02	E08	HQP009678	NM_000584	IL8
QG031-02	E09	HQP009677	NM_002185	IL7R
QG031-02	E10	HQP009676	NM_000880	IL7
QG031-02	E11	HQP009674	NM_002184	IL6ST
QG031-02	E12	HQP009672	NM_000565	IL6R
QG031-02	F01	HQP009670	NM_000600	IL6
QG031-02	F02	HQP009666	NM_000879	IL5
QG031-02	F03	HQP009664	NM_000418	IL4R
QG031-02	F04	HQP009662	NM_000589	IL4
QG031-02	F05	HQP009659	NM_000206	IL2RG
QG031-02	F06	HQP009658	NM_000878	IL2RB
QG031-02	F07	HQP009650	NM_000417	IL2RA
QG031-02	F08	HQP022792	NM_004843	IL27RA
QG031-02	F09	HQP003495	NM_144701	IL23R
QG031-02	F10	HQP012859	NM_016584	IL23A
QG031-02	F11	HQP012160	NM_021798	IL21R
QG031-02	F12	HQP016102	NM_021803	IL21
QG031-02	G01	HQP009649	NM_000586	IL2
QG031-02	G02	HQP009645	NM_000577	IL1RN

QG031-02	G03	HQP022241	NM_003856	IL1RL1
QG031-02	G04	HQP009643	NM_002182	IL1RAP
QG031-02	G05	HQP018800	NM_004633	IL1R2
QG031-02	G06	HQP009642	NM_000877	IL1R1
QG031-02	G07	HQP009641	NM_000576	IL1B
QG031-02	G08	HQP009640	NM_000575	IL1A
QG031-02	G09	HQP009718	NM_001562	IL18
QG031-02	G10	HQP001495	NM_052872	IL17F
QG031-02	G11	HQP009717	NM_002190	IL17A
QG031-02	G12	HQP009708	NM_000585	IL15
QG031-02	H01	HGDC		
QG031-02	H02	HGDC		
QG031-02	H03	HQP006940	NM_002046	GAPDH
QG031-02	H04	HQP016381	NM_001101	ACTB
QG031-02	H05	HQP015171	NM_004048	B2M
QG031-02	H06	HQP006171	NM_012423	RPL13A
QG031-02	H07	HQP009026	NM_000194	HPRT1
QG031-02	H08	HQP054253	NR_003286	RN18S1
QG031-02	H09	RT		
QG031-02	H10	RT		
QG031-02	H11	PCR		
QG031-02	H12	PCR		
QG031-03	A01	HQP009700	NM_001560	IL13RA1
QG031-03	A02	HQP009697	NM_002188	IL13
QG031-03	A03	HQP009694	NM_005535	IL12RB1
QG031-03	A04	HQP009693	NM_002187	IL12B
QG031-03	A05	HQP009692	NM_000882	IL12A
QG031-03	A06	HQP009690	NM_004512	IL11RA
QG031-03	A07	HQP009686	NM_001558	IL10RA
QG031-03	A08	HQP009685	NM_000572	IL10
QG031-03	A09	HQP021140	NM_003639	IKBKG
QG031-03	A10	HQP009544	NM_000598	IGFBP3
QG031-03	A11	HQP009467	NM_000619	IFNG
QG031-03	A12	HQP009463	NM_002176	IFNB1
QG031-03	B01	HQP009460	NM_000874	IFNAR2
QG031-03	B02	HQP009458	NM_000629	IFNAR1
QG031-03	B03	HQP009430	NM_021057	IFNA7
QG031-03	B04	HQP021142	NM_003641	IFITM1
QG031-03	B05	HQP009403	NM_001547	IFIT2
QG031-03	B06	HQP009393	NM_005533	IFI35
QG031-03	B07	HQP005977	NM_015259	ICOSLG
QG031-03	B08	HQP008554	NM_012092	ICOS
QG031-03	B09	HQP009187	NM_000873	ICAM2

QG031-03	B10	HQP009184	NM_000201	ICAM1
QG031-03	B11	HQP009077	NM_005345	HSPA1A
QG031-03	B12	HQP018231	NM_003299	HSP90B1
QG031-03	C01	HQP016162	NM_021624	HRH4
QG031-03	C02	HQP008883	NM_002128	HMGB1
QG031-03	C03	HQP008867	NM_002124	HLA-DRB1
QG031-03	C04	HQP008866	NM_019111	HLA-DRA
QG031-03	C05	HQP008861	NM_002122	HLA-DQA1
QG031-03	C06	HQP008851	NM_002117	HLA-C
QG031-03	C07	HQP008849	NM_002116	HLA-A
QG031-03	C08	HQP008757	NM_000410	HFE
QG031-03	C09	HQP008690	NM_004131	GZMB
QG031-03	C10	HQP008689	NM_006144	GZMA
QG031-03	C11	HQP007165	NM_001002295	GATA3
QG031-03	C12	HQP006484	NM_002037	FYN
QG031-03	D01	HQP006426	NM_004958	FRAP1
QG031-03	D02	HQP012269	NM_014009	FOXP3
QG031-03	D03	HQP006188	NM_005252	FOS
QG031-03	D04	HQP006022	NM_002026	FN1
QG031-03	D05	HQP005323	NM_004107	FCGRT
QG031-03	D06	HQP005308	NM_000569	FCGR3A
QG031-03	D07	HQP005290	NM_001002273	FCGR2B
QG031-03	D08	HQP005280	NM_021642	FCGR2A
QG031-03	D09	HQP005251	NM_000566	FCGR1A
QG031-03	D10	HQP005242	NM_002002	FCER2
QG031-03	D11	HQP005238	NM_002001	FCER1A
QG031-03	D12	HQP009671	NM_000639	FASLG
QG031-03	E01	HQP009651	NM_000043	FAS
QG031-03	E02	HQP005068	NM_000505	F12
QG031-03	E03	HQP022883	NM_001776	ENTPD1
QG031-03	E04	HQP004605	NM_005228	EGFR
QG031-03	E05	HQP004599	NM_001963	EGF
QG031-03	E06	HQP004557	NM_001955	EDN1
QG031-03	E07	HQP004536	NM_004951	EBI2
QG031-03	E08	HQP004534	NM_024007	EBF1
QG031-03	E09	HQP004434	NM_001935	DPP4
QG031-03	E10	HQP015116	NM_019887	DIABLO
QG031-03	E11	HQP004193	NM_004942	DEFB4
QG031-03	E12	HQP004191	NM_005218	DEFB1
QG031-03	F01	HQP004173	NM_004084	DEFA1
QG031-03	F02	HQP003992	NM_001350	DAXX
QG031-03	F03	HQP018802	NM_001008540	CXCR4
QG031-03	F04	HQP007900	NM_001504	CXCR3

QG031-03	F05	HQP011220	NM_002416	CXCL9
QG031-03	F06	HQP008459	NM_002090	CXCL3
QG031-03	F07	HQP008458	NM_002089	CXCL2
QG031-03	F08	HQP016669	NM_000609	CXCL12
QG031-03	F09	HQP009746	NM_001565	CXCL10
QG031-03	F10	HQP003691	NM_001337	CX3CR1
QG031-03	F11	HQP016652	NM_002996	CX3CL1
QG031-03	F12	HQP003612	NM_001911	CTSG
QG031-03	G01	HQP003599	NM_001909	CTSD
QG031-03	G02	HQP003596	NM_001908	CTSB
QG031-03	G03	HQP003499	NM_001037631	CTLA4
QG031-03	G04	HQP003185	NM_000760	CSF3R
QG031-03	G05	HQP003173	NM_000759	CSF3
QG031-03	G06	HQP003159	NM_000758	CSF2
QG031-03	G07	HQP003158	NM_005211	CSF1R
QG031-03	G08	HQP003149	NM_000757	CSF1
QG031-03	G09	HQP002994	NM_000567	CRP
QG031-03	G10	HQP002889	NM_001006658	CR2
QG031-03	G11	HQP002242	NM_004072	CMKLR1
QG031-03	G12	HQP002037	NM_001831	CLU
QG031-03	H01	HGDC		
QG031-03	H02	HGDC		
QG031-03	H03	HQP006940	NM_002046	GAPDH
QG031-03	H04	HQP016381	NM_001101	ACTB
QG031-03	H05	HQP015171	NM_004048	B2M
QG031-03	H06	HQP006171	NM_012423	RPL13A
QG031-03	H07	HQP009026	NM_000194	HPRT1
QG031-03	H08	HQP054253	NR_003286	RN18S1
QG031-03	H09	RT		
QG031-03	H10	RT		
QG031-03	H11	PCR		
QG031-03	H12	PCR		
QG031-04	A01	HQP006263	NM_013252	CLEC5A
QG031-04	A02	HQP011211	NM_000246	CIITA
QG031-04	A03	HQP001708	NM_001278	CHUK
QG031-04	A04	HQP008755	NM_000186	CFH
QG031-04	A05	HQP004198	NM_001928	CFD
QG031-04	A06	HQP016565	NM_001710	CFB
QG031-04	A07	HQP016623	NM_001024912	CEACAM1
QG031-04	A08	HQP000331	NM_000389	CDKN1A
QG031-04	A09	HQP000052	NM_001795	CDH5
QG031-04	A10	HQP011212	NM_002414	CD99
QG031-04	A11	HQP022474	NM_001769	CD9

QG031-04	A12	HQP022746	NM_006889	CD86
QG031-04	B01	HQP022722	NM_005191	CD80
QG031-04	B02	HQP023094	NM_001781	CD69
QG031-04	B03	HQP022384	NM_006725	CD6
QG031-04	B04	HQP023056	NM_000611	CD59
QG031-04	B05	HQP003948	NM_000574	CD55
QG031-04	B06	HQP022337	NM_014207	CD5
QG031-04	B07	HQP022972	NM_000610	CD44
QG031-04	B08	HQP022962	NM_000074	CD40LG
QG031-04	B09	HQP022955	NM_001250	CD40
QG031-04	B10	HQP022316	NM_000616	CD4
QG031-04	B11	HQP000971	NM_012099	CD3EAP
QG031-04	B12	HQP022236	NM_000733	CD3E
QG031-04	C01	HQP022821	NM_000072	CD36
QG031-04	C02	HQP022812	NM_001025109	CD34
QG031-04	C03	HQP006266	NM_012120	CD2AP
QG031-04	C04	HQP022699	NM_006139	CD28
QG031-04	C05	HQP022295	NM_000734	CD247
QG031-04	C06	HQP013003	NM_016382	CD244
QG031-04	C07	HQP022579	NM_013230	CD24
QG031-04	C08	HQP022561	NM_001771	CD22
QG031-04	C09	HQP008808	NM_021155	CD209
QG031-04	C10	HQP022190	NM_001767	CD2
QG031-04	C11	HQP022129	NM_001766	CD1D
QG031-04	C12	HQP022504	NM_001770	CD19
QG031-04	D01	HQP010932	NM_005582	CD180
QG031-04	D02	HQP022490	NM_000591	CD14
QG031-04	D03	HQP002222	NM_005201	CCR8
QG031-04	D04	HQP002217	NM_001838	CCR7
QG031-04	D05	HQP002212	NM_004367	CCR6
QG031-04	D06	HQP002200	NM_000647	CCR2
QG031-04	D07	HQP016626	NM_002985	CCL5
QG031-04	D08	HQP016625	NM_002984	CCL4
QG031-04	D09	HQP053971	NM_021006	CCL3L1
QG031-04	D10	HQP016641	NM_002990	CCL22
QG031-04	D11	HQP016639	NM_004591	CCL20
QG031-04	D12	HQP016638	NM_006274	CCL19
QG031-04	E01	HQP016630	NM_005408	CCL13
QG031-04	E02	HQP016629	NM_002986	CCL11
QG031-04	E03	HQP002229	NM_001296	CCBP2
QG031-04	E04	HQP021430	NM_005188	CBL
QG031-04	E05	HQP018966	NM_001080124	CASP8
QG031-04	E06	HQP020297	NM_004346	CASP3



QG031-04	E07	HQP020249	NM_032982	CASP2
QG031-04	E08	HQP020207	NM_001223	CASP1
QG031-04	E09	HQP020718	NM_032415	CARD11
QG031-04	E10	HQP020046	NM_004345	CAMP
QG031-04	E11	HQP019841	NM_004343	CALR
QG031-04	E12	HQP018414	NM_001737	C9
QG031-04	F01	HQP018362	NM_000562	C8A
QG031-04	F02	HQP018344	NM_000065	C6
QG031-04	F03	HQP018304	NM_001735	C5
QG031-04	F04	HQP018248	NM_007293	C4A
QG031-04	F05	HQP018238	NM_000064	C3
QG031-04	F06	HQP018226	NM_000063	C2
QG031-04	F07	HQP018206	NM_001734	C1S
QG031-04	F08	HQP018146	NM_015991	C1QA
QG031-04	F09	HQP017988	NM_000061	BTK
QG031-04	F10	HQP017834	NM_004334	BST1
QG031-04	F11	HQP008515	NM_013314	BLNK
QG031-04	F12	HQP009084	NM_001165	BIRC3
QG031-04	G01	HQP016653	NM_001196	BID
QG031-04	G02	HQP016328	NM_001706	BCL6
QG031-04	G03	HQP016287	NM_005178	BCL3
QG031-04	G04	HQP016237	NM_001191	BCL2L1
QG031-04	G05	HQP016211	NM_000633	BCL2
QG031-04	G06	HQP000166	NM_005745	BCAP31
QG031-04	G07	HQP015964	NM_004324	BAX
QG031-04	G08	HQP015538	NM_004322	BAD
QG031-04	G09	HQP015171	NM_004048	B2M
QG031-04	G10	HQP011736	NM_000051	ATM
QG031-04	G11	HQP011667	NM_005171	ATF1
QG031-04	G12	HQP004991	NM_001014431	AKT1
QG031-04	H01	HGDC		
QG031-04	H02	HGDC		
QG031-04	H03	HQP006940	NM_002046	GAPDH
QG031-04	H04	HQP016381	NM_001101	ACTB
QG031-04	H05	HQP015171	NM_004048	B2M
QG031-04	H06	HQP006171	NM_012423	RPL13A
QG031-04	H07	HQP009026	NM_000194	HPRT1
QG031-04	H08	HQP054253	NR_003286	RN18S1
QG031-04	H09	RT		
QG031-04	H10	RT		
QG031-04	H11	PCR		
QG031-04	H12	PCR		