

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG097-01	A01	HQP001281	NM_007048	BTN3A1
QG097-01	A02	HQP000996	NM_006678	CD300C
QG097-01	A03	HQP000971	NM_012099	CD3EAP
QG097-01	A04	HQP000821	NM_006573	TNFSF13B
QG097-01	A05	HQP000811	NM_006566	CD226
QG097-01	A06	HQP000808	NM_006564	CXCR6
QG097-01	A07	HQP000652	NM_006404	PROCR
QG097-01	A08	HQP000613	NM_006378	SEMA4D
QG097-01	A09	HQP000415	NM_006068	TLR6
QG097-01	A10	HQP000411	NM_014257	CLEC4M
QG097-01	A11	HQP000279	NM_005816	CD96
QG097-01	A12	HQP023346	NM_006039	MRC2
QG097-01	B01	HQP022750	NM_004829	NCR1
QG097-01	B02	HQP022745	NM_004827	ABCG2
QG097-01	B03	HQP022665	NM_004258	IGSF2
QG097-01	B04	HQP022548	NM_004244	CD163
QG097-01	B05	HQP022500	NM_001040280	CD83
QG097-01	B06	HQP021613	NM_006017	PROM1
QG097-01	B07	HQP021599	NM_003874	CD84
QG097-01	B08	HQP021557	NM_003844	TNFRSF10A
QG097-01	B09	HQP021553	NM_003842	TNFRSF10B
QG097-01	B10	HQP021552	NM_003841	TNFRSF10C
QG097-01	B11	HQP021551	NM_003840	TNFRSF10D
QG097-01	B12	HQP021550	NM_003839	TNFRSF11A
QG097-01	C01	HQP021533	NM_003830	SIGLEC5
QG097-01	C02	HQP021502	NM_003810	TNFSF10
QG097-01	C03	HQP021496	NM_003807	TNFSF14
QG097-01	C04	HQP021142	NM_003641	IFITM1
QG097-01	C05	HQP020964	NM_003612	SEMA7A
QG097-01	C06	HQP020141	NM_003508	FZD9
QG097-01	C07	HQP020134	NM_012193	FZD4
QG097-01	C08	HQP018802	NM_001008540	CXCR4
QG097-01	C09	HQP018466	NM_080682	VCAM1
QG097-01	C10	HQP018329	NM_003326	TNFSF4
QG097-01	C11	HQP018120	NM_004615	TSPAN7
QG097-01	C12	HQP018116	NM_138554	TLR4
QG097-01	D01	HQP018115	NM_003265	TLR3
QG097-01	D02	HQP018114	NM_003264	TLR2
QG097-01	D03	HQP018113	NM_003263	TLR1
QG097-01	D04	HQP018083	NM_006288	THY1
QG097-01	D05	HQP018067	NM_000361	THBD

QG097-01	D06	HQP018041	NM_003234	TFRC
QG097-01	D07	HQP017670	NM_001030288	SPN
QG097-01	D08	HQP017365	NM_000342	SLC4A1
QG097-01	D09	HQP017303	NM_003037	SLAMF1
QG097-01	D10	HQP017179	NM_003032	ST6GAL1
QG097-01	D11	HQP016747	NM_003006	SELPLG
QG097-01	D12	HQP016746	NM_003005	SELP
QG097-01	E01	HQP016745	NM_000655	SELL
QG097-01	E02	HQP016659	NM_001006946	SDC1
QG097-01	E03	HQP016270	NM_000324	RHAG
QG097-01	E04	HQP015962	NM_001042724	PVRL2
QG097-01	E05	HQP015956	NM_002855	PVRL1
QG097-01	E06	HQP015954	NM_006505	PVR
QG097-01	E07	HQP015925	NM_002843	PTPRJ
QG097-01	E08	HQP015908	NM_002838	PTPRC
QG097-01	E09	HQP015570	NM_020440	PTGFRN
QG097-01	E10	HQP015032	NM_000311	PRNP
QG097-01	E11	HQP013205	NM_001005376	PLAUR
QG097-01	E12	HQP013100	NM_000927	ABCB1
QG097-01	F01	HQP013015	NM_000442	PECAM1
QG097-01	F02	HQP012662	NM_005018	PDCD1
QG097-01	F03	HQP011932	NM_002526	NT5E
QG097-01	F04	HQP011828	NM_002507	NGFR
QG097-01	F05	HQP011559	NM_001018016	MUC1
QG097-01	F06	HQP011525	NM_002447	MST1R
QG097-01	F07	HQP011319	NM_002438	MRC1
QG097-01	F08	HQP011308	NM_005373	MPL
QG097-01	F09	HQP011301	NM_001004196	CD200
QG097-01	F10	HQP011254	NM_000902	MME
QG097-01	F11	HQP011212	NM_002414	CD99
QG097-01	F12	HQP011189	NM_005929	MFI2
QG097-01	G01	HQP011113	NM_002389	CD46
QG097-01	G02	HQP011099	NM_006500	MCAM
QG097-01	G03	HQP010942	NM_002354	TACSTD1
QG097-01	G04	HQP010933	NM_002349	LY75
QG097-01	G05	HQP010932	NM_005582	CD180
QG097-01	G06	HQP010930	NM_001033667	LY9
QG097-01	G07	HQP010922	NM_001013257	BCAM
QG097-01	G08	HQP010915	NM_002342	LTBR
QG097-01	G09	HQP010870	NM_002332	LRP1
QG097-01	G10	HQP010608	NM_002310	LIFR
QG097-01	G11	HQP010582	NM_001003679	LEPR
QG097-01	G12	HQP010478	NM_002288	LAIR2

QG097-01	H01	HGDC		
QG097-01	H02	HGDC		
QG097-01	H03	HQP006940	NM_002046	GAPDH
QG097-01	H04	HQP016381	NM_001101	ACTB
QG097-01	H05	HQP015171	NM_004048	B2M
QG097-01	H06	HQP006171	NM_012423	RPL13A
QG097-01	H07	HQP009026	NM_000194	HPRT1
QG097-01	H08	HQP054253	NR_003286	RN18S1
QG097-01	H09	RT		
QG097-01	H10	RT		
QG097-01	H11	PCR		
QG097-01	H12	PCR		
QG097-02	A01	HQP010467	NM_002287	LAIR1
QG097-02	A02	HQP010390	NM_000425	L1CAM
QG097-02	A03	HQP010112	NM_002262	KLRD1
QG097-02	A04	HQP010106	NM_002258	KLRB1
QG097-02	A05	HQP010099	NM_000222	KIT
QG097-02	A06	HQP010095	NM_013289	KIR3DL1
QG097-02	A07	HQP010083	NM_014218	KIR2DL1
QG097-02	A08	HQP010071	NM_000420	KEL
QG097-02	A09	HQP010070	NM_002253	KDR
QG097-02	A10	HQP009819	NM_000213	ITGB4
QG097-02	A11	HQP009818	NM_000212	ITGB3
QG097-02	A12	HQP009815	NM_000211	ITGB2
QG097-02	B01	HQP009810	NM_002211	ITGB1
QG097-02	B02	HQP009808	NM_002210	ITGAV
QG097-02	B03	HQP009805	NM_002208	ITGAE
QG097-02	B04	HQP009804	NM_005353	ITGAD
QG097-02	B05	HQP009795	NM_000419	ITGA2B
QG097-02	B06	HQP009789	NM_002201	ISG20
QG097-02	B07	HQP009764	NM_000208	INSR
QG097-02	B08	HQP009716	NM_001561	TNFRSF9
QG097-02	B09	HQP009694	NM_005535	IL12RB1
QG097-02	B10	HQP009683	NM_002186	IL9R
QG097-02	B11	HQP009681	NM_001557	IL8RB
QG097-02	B12	HQP009679	NM_000634	IL8RA
QG097-02	C01	HQP009677	NM_002185	IL7R
QG097-02	C02	HQP009674	NM_002184	IL6ST
QG097-02	C03	HQP009672	NM_000565	IL6R
QG097-02	C04	HQP009667	NM_000564	IL5RA
QG097-02	C05	HQP009664	NM_000418	IL4R
QG097-02	C06	HQP009661	NM_002183	IL3RA
QG097-02	C07	HQP009659	NM_000206	IL2RG

QG097-02	C08	HQP009658	NM_000878	IL2RB
QG097-02	C09	HQP009650	NM_000417	IL2RA
QG097-02	C10	HQP009532	NM_000876	IGF2R
QG097-02	C11	HQP009523	NM_000875	IGF1R
QG097-02	C12	HQP009469	NM_000416	IFNGR1
QG097-02	D01	HQP009201	NM_001039132	ICAM4
QG097-02	D02	HQP009187	NM_000873	ICAM2
QG097-02	D03	HQP009184	NM_000201	ICAM1
QG097-02	D04	HQP008896	NM_012484	HMMR
QG097-02	D05	HQP008634	NM_002101	GYPC
QG097-02	D06	HQP007900	NM_001504	CXCR3
QG097-02	D07	HQP007803	NM_000173	GP1BA
QG097-02	D08	HQP006479	NM_002036	DARC
QG097-02	D09	HQP006455	NM_002033	FUT4
QG097-02	D10	HQP005890	NM_004119	FLT3
QG097-02	D11	HQP005439	NM_002011	FGFR4
QG097-02	D12	HQP005437	NM_000141	FGFR2
QG097-02	E01	HQP005434	NM_000142	FGFR3
QG097-02	E02	HQP005308	NM_000569	FCGR3A
QG097-02	E03	HQP005290	NM_001002273	FCGR2B
QG097-02	E04	HQP005280	NM_021642	FCGR2A
QG097-02	E05	HQP005251	NM_000566	FCGR1A
QG097-02	E06	HQP005242	NM_002002	FCER2
QG097-02	E07	HQP005057	NM_001993	F3
QG097-02	E08	HQP004968	NM_001005862	ERBB2
QG097-02	E09	HQP004867	NM_001977	ENPEP
QG097-02	E10	HQP004856	NM_000118	ENG
QG097-02	E11	HQP004434	NM_001935	DPP4
QG097-02	E12	HQP004081	NM_000789	ACE
QG097-02	F01	HQP003948	NM_000574	CD55
QG097-02	F02	HQP003499	NM_001037631	CTLA4
QG097-02	F03	HQP003185	NM_000760	CSF3R
QG097-02	F04	HQP003170	NM_000395	CSF2RB
QG097-02	F05	HQP003158	NM_005211	CSF1R
QG097-02	F06	HQP002889	NM_001006658	CR2
QG097-02	F07	HQP002217	NM_001838	CCR7
QG097-02	F08	HQP002212	NM_004367	CCR6
QG097-02	F09	HQP002210	NM_000579	CCR5
QG097-02	F10	HQP002198	NM_001295	CCR1
QG097-02	F11	HQP001014	NM_001816	CEACAM8
QG097-02	F12	HQP000593	NM_004363	CEACAM5
QG097-02	G01	HQP000525	NM_001803	CD52
QG097-02	G02	HQP000052	NM_001795	CDH5

QG097-02	G03	HQP000018	NM_001792	CDH2
QG097-02	G04	HQP023192	NM_001039490	CD151
QG097-02	G05	HQP023178	NM_001025160	CD97
QG097-02	G06	HQP023168	NM_004356	CD81
QG097-02	G07	HQP023155	NM_000626	CD79B
QG097-02	G08	HQP023117	NM_001782	CD72
QG097-02	G09	HQP023108	NM_001252	CD70
QG097-02	G10	HQP023094	NM_001781	CD69
QG097-02	G11	HQP023083	NM_001040059	CD68
QG097-02	G12	HQP023069	NM_001040034	CD63
QG097-02	H01	HGDC		
QG097-02	H02	HGDC		
QG097-02	H03	HQP006940	NM_002046	GAPDH
QG097-02	H04	HQP016381	NM_001101	ACTB
QG097-02	H05	HQP015171	NM_004048	B2M
QG097-02	H06	HQP006171	NM_012423	RPL13A
QG097-02	H07	HQP009026	NM_000194	HPRT1
QG097-02	H08	HQP054253	NR_003286	RN18S1
QG097-02	H09	RT		
QG097-02	H10	RT		
QG097-02	H11	PCR		
QG097-02	H12	PCR		
QG097-03	A01	HQP023056	NM_000611	CD59
QG097-03	A02	HQP023046	NM_001779	CD58
QG097-03	A03	HQP023005	NM_001778	CD48
QG097-03	A04	HQP022972	NM_000610	CD44
QG097-03	A05	HQP022962	NM_000074	CD40LG
QG097-03	A06	HQP022955	NM_001250	CD40
QG097-03	A07	HQP022934	NM_001249	ENTPD5
QG097-03	A08	HQP022921	NM_001248	ENTPD3
QG097-03	A09	HQP022912	NM_001247	ENTPD6
QG097-03	A10	HQP022894	NM_001246	ENTPD2
QG097-03	A11	HQP022883	NM_001776	ENTPD1
QG097-03	A12	HQP022870	NM_001775	CD38
QG097-03	B01	HQP022855	NM_001040031	CD37
QG097-03	B02	HQP022846	NM_005506	SCARB2
QG097-03	B03	HQP022821	NM_000072	CD36
QG097-03	B04	HQP022812	NM_001025109	CD34
QG097-03	B05	HQP022769	NM_001244	TNFSF8
QG097-03	B06	HQP022753	NM_001243	TNFRSF8
QG097-03	B07	HQP022746	NM_006889	CD86
QG097-03	B08	HQP022722	NM_005191	CD80
QG097-03	B09	HQP022699	NM_006139	CD28

QG097-03	B10	HQP022667	NM_001242	CD27
QG097-03	B11	HQP022561	NM_001771	CD22
QG097-03	B12	HQP022504	NM_001770	CD19
QG097-03	C01	HQP022490	NM_000591	CD14
QG097-03	C02	HQP022474	NM_001769	CD9
QG097-03	C03	HQP022438	NM_004931	CD8B
QG097-03	C04	HQP022399	NM_006137	CD7
QG097-03	C05	HQP022384	NM_006725	CD6
QG097-03	C06	HQP022357	NM_005894	CD5L
QG097-03	C07	HQP022337	NM_014207	CD5
QG097-03	C08	HQP022316	NM_000616	CD4
QG097-03	C09	HQP022295	NM_000734	CD247
QG097-03	C10	HQP022256	NM_000073	CD3G
QG097-03	C11	HQP022236	NM_000733	CD3E
QG097-03	C12	HQP022212	NM_000732	CD3D
QG097-03	D01	HQP022190	NM_001767	CD2
QG097-03	D02	HQP022159	NM_001042583	CD1E
QG097-03	D03	HQP022129	NM_001766	CD1D
QG097-03	D04	HQP022111	NM_001765	CD1C
QG097-03	D05	HQP022090	NM_001764	CD1B
QG097-03	D06	HQP022070	NM_001763	CD1A
QG097-03	D07	HQP018768	NM_001954	DDR1
QG097-03	D08	HQP018326	NM_001736	C5AR1
QG097-03	D09	HQP017844	NM_004335	BST2
QG097-03	D10	HQP017834	NM_004334	BST1
QG097-03	D11	HQP017820	NM_001728	BSG
QG097-03	D12	HQP017489	NM_004329	BMPR1A
QG097-03	E01	HQP016959	NM_001716	BLR1
QG097-03	E02	HQP016623	NM_001024912	CEACAM1
QG097-03	E03	HQP016367	NM_001192	TNFRSF17
QG097-03	E04	HQP011863	NM_001679	ATP1B3
QG097-03	E05	HQP011153	NM_021071	ART4
QG097-03	E06	HQP009671	NM_000639	FASLG
QG097-03	E07	HQP009651	NM_000043	FAS
QG097-03	E08	HQP008414	NM_001150	ANPEP
QG097-03	E09	HQP006353	NM_004304	ALK
QG097-03	E10	HQP004470	NM_000214	JAG1
QG097-03	E11	HQP011127	NM_004314	ART1
QG097-03	E12	HQP002207	NM_001837	CCR3
QG097-03	F01	HQP009192	NM_002162	ICAM3
QG097-03	F02	HQP010087	NM_001080770	KIR2DL4
QG097-03	F03	HQP011520	NM_002445	MSR1
QG097-03	F04	HQP000246	NM_001109	ADAM8

QG097-03	F05	HQP010462	NM_002286	LAG3
QG097-03	F06	HQP022541	NM_001031666	MS4A3
QG097-03	F07	HQP021167	NM_001082618	CD33
QG097-03	F08	HQP022797	NM_001245	SIGLEC6
QG097-03	F09	HQP021321	NM_003701	TNFSF11
QG097-03	F10	HQP023019	NM_000560	CD53
QG097-03	F11	HQP023127	NM_001025158	CD74
QG097-03	F12	HQP023144	NM_001783	CD79A
QG097-03	G01	HQP023466	NM_004360	CDH1
QG097-03	G02	HQP000972	NM_001815	CEACAM3
QG097-03	G03	HQP002880	NM_000573	CR1
QG097-03	G04	HQP003163	NM_006140	CSF2RA
QG097-03	G05	HQP005227	NM_002000	FCAR
QG097-03	G06	HQP005427	NM_015850	FGFR1
QG097-03	G07	HQP006454	NM_000149	FUT3
QG097-03	G08	HQP007378	NM_001032364	GGT1
QG097-03	G09	HQP007809	NM_004488	GP5
QG097-03	G10	HQP023380	NM_014880	CD302
QG097-03	G11	HQP009860	NM_001024844	CD82
QG097-03	G12	HQP000353	NM_001080978	LILRB2
QG097-03	H01	HGDC		
QG097-03	H02	HGDC		
QG097-03	H03	HQP006940	NM_002046	GAPDH
QG097-03	H04	HQP016381	NM_001101	ACTB
QG097-03	H05	HQP015171	NM_004048	B2M
QG097-03	H06	HQP006171	NM_012423	RPL13A
QG097-03	H07	HQP009026	NM_000194	HPRT1
QG097-03	H08	HQP054253	NR_003286	RN18S1
QG097-03	H09	RT		
QG097-03	H10	RT		
QG097-03	H11	PCR		
QG097-03	H12	PCR		