

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG003-01	A01	HQP016367	NM_001192	TNFRSF17
QG003-01	A02	HQP018114	NM_003264	TLR2
QG003-01	A03	HQP018074	NM_000460	THPO
QG003-01	A04	HQP018044	NM_000660	TGFB1
QG003-01	A05	HQP017984	NM_001008897	TCP1
QG003-01	A06	HQP017902	NM_003190	TAPBP
QG003-01	A07	HQP017900	NM_000544	TAP2
QG003-01	A08	HQP017899	NM_000593	TAP1
QG003-01	A09	HQP017845	NM_003177	SYK
QG003-01	A10	HQP017764	NM_007315	STAT1
QG003-01	A11	HQP021399	NM_003745	SOCS1
QG003-01	A12	HQP010229	NM_001664	RHOA
QG003-01	B01	HQP016247	NM_000538	RFXAP
QG003-01	B02	HQP021359	NM_003721	RFXANK
QG003-01	B03	HQP016246	NM_000449	RFX5
QG003-01	B04	HQP016063	NM_006908	RAC1
QG003-01	B05	HQP015512	NM_002818	PSME2
QG003-01	B06	HQP015499	NM_006263	PSME1
QG003-01	B07	HQP015311	NM_002800	PSMB9
QG003-01	B08	HQP015291	NM_004159	PSMB8
QG003-01	B09	HQP014731	NM_006254	PRKCD
QG003-01	B10	HQP008463	NM_005313	PDIA3
QG003-01	B11	HQP002334	NM_130771	OSCAR
QG003-01	B12	HQP022857	NM_006310	NPEPPS
QG003-01	C01	HQP003133	NM_001315	MAPK14
QG003-01	C02	HQP010933	NM_002349	LY75
QG003-01	C03	HQP015082	NM_001008530	LGMN
QG003-01	C04	HQP012337	NM_015907	LAP3
QG003-01	C05	HQP010540	NM_002294	LAMP2
QG003-01	C06	HQP009251	NM_002163	IRF8
QG003-01	C07	HQP009779	NM_002199	IRF2
QG003-01	C08	HQP009467	NM_000619	IFNG
QG003-01	C09	HQP000520	NM_006332	IFI30
QG003-01	C10	HQP005977	NM_015259	ICOSLG
QG003-01	C11	HQP008554	NM_012092	ICOS
QG003-01	C12	HQP009184	NM_000201	ICAM1
QG003-01	D01	HQP009083	NM_005347	HSPA5
QG003-01	D02	HQP009077	NM_005345	HSPA1A
QG003-01	D03	HQP018231	NM_003299	HSP90B1
QG003-01	D04	HQP009097	NM_007355	HSP90AB1
QG003-01	D05	HQP009092	NM_001017963	HSP90AA1

QG003-01	D06	HQP008874	NM_005516	HLA-E
QG003-01	D07	HQP008867	NM_002124	HLA-DRB1
QG003-01	D08	HQP008866	NM_019111	HLA-DRA
QG003-01	D09	HQP008861	NM_002122	HLA-DQA1
QG003-01	D10	HQP008860	NM_002121	HLA-DPB1
QG003-01	D11	HQP008859	NM_033554	HLA-DPA1
QG003-01	D12	HQP008858	NM_002120	HLA-DOB
QG003-01	E01	HQP008857	NM_002119	HLA-DOA
QG003-01	E02	HQP008853	NM_002118	HLA-DMB
QG003-01	E03	HQP008852	NM_006120	HLA-DMA
QG003-01	E04	HQP008851	NM_002117	HLA-C
QG003-01	E05	HQP008849	NM_002116	HLA-A
QG003-01	E06	HQP008757	NM_000410	HFE
QG003-01	E07	HQP005290	NM_001002273	FCGR2B
QG003-01	E08	HQP004968	NM_001005862	ERBB2
QG003-01	E09	HQP003668	NM_004079	CTSS
QG003-01	E10	HQP003605	NM_001910	CTSE
QG003-01	E11	HQP003599	NM_001909	CTSD
QG003-01	E12	HQP003596	NM_001908	CTSB
QG003-01	F01	HQP002907	NM_004379	CREB1
QG003-01	F02	HQP011211	NM_000246	CIITA
QG003-01	F03	HQP023454	NM_001039802	CDC42
QG003-01	F04	HQP022746	NM_006889	CD86
QG003-01	F05	HQP011113	NM_002389	CD46
QG003-01	F06	HQP022955	NM_001250	CD40
QG003-01	F07	HQP008808	NM_021155	CD209
QG003-01	F08	HQP012139	NM_015717	CD207
QG003-01	F09	HQP022129	NM_001766	CD1D
QG003-01	F10	HQP019917	NM_015981	CAMK2A
QG003-01	F11	HQP019841	NM_004343	CALR
QG003-01	F12	HQP016897	NM_000386	BLMH
QG003-01	G01	HQP015171	NM_004048	B2M
QG003-01	G02	HQP013009	NM_001040458	ARTS-1
QG003-01	G03	HQP007413	NM_012095	AP3M1
QG003-01	G04	HQP021252	NM_003664	AP3B1
QG003-01	G05	HQP008865	NM_182549	HLA-DQB2
QG003-01	G06	HQP008876	NM_002127	HLA-G
QG003-01	G07	HQP008878	NM_001531	MR1
QG003-01	G08	HQP008850	NM_005514	HLA-B
QG003-01	G09	HQP008875	NM_018950	HLA-F
QG003-01	G10	HQP020051	NM_001024649	CANX
QG003-01	G11	HQP023127	NM_001025158	CD74
QG003-01	G12	HQP013299	NM_002675	PML

QG003-01	H01	HGDC		
QG003-01	H02	HGDC		
QG003-01	H03	HQP006940	NM_002046	GAPDH
QG003-01	H04	HQP016381	NM_001101	ACTB
QG003-01	H05	HQP015171	NM_004048	B2M
QG003-01	H06	HQP006171	NM_012423	RPL13A
QG003-01	H07	HQP009026	NM_000194	HPRT1
QG003-01	H08	HQP054253	NR_003286	RN18S1
QG003-01	H09	RT		
QG003-01	H10	RT		
QG003-01	H11	PCR		
QG003-01	H12	PCR		