

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG008-01	A01	HQP004535	NM_005161	AGTRL1
QG008-01	A02	HQP009651	NM_000043	FAS
QG008-01	A03	HQP016367	NM_001192	TNFRSF17
QG008-01	A04	HQP016959	NM_001716	BLR1
QG008-01	A05	HQP022667	NM_001242	CD27
QG008-01	A06	HQP022753	NM_001243	TNFRSF8
QG008-01	A07	HQP022955	NM_001250	CD40
QG008-01	A08	HQP001708	NM_001278	CHUK
QG008-01	A09	HQP002198	NM_001295	CCR1
QG008-01	A10	HQP002210	NM_000579	CCR5
QG008-01	A11	HQP002212	NM_004367	CCR6
QG008-01	A12	HQP002217	NM_001838	CCR7
QG008-01	B01	HQP002222	NM_005201	CCR8
QG008-01	B02	HQP002242	NM_004072	CMKLR1
QG008-01	B03	HQP003691	NM_001337	CX3CR1
QG008-01	B04	HQP004324	NM_006729	DIAPH2
QG008-01	B05	HQP004605	NM_005228	EGFR
QG008-01	B06	HQP004958	NM_000121	EPOR
QG008-01	B07	HQP006479	NM_002036	DARC
QG008-01	B08	HQP007833	NM_016602	CCR10
QG008-01	B09	HQP007900	NM_001504	CXCR3
QG008-01	B10	HQP008143	NM_005299	GPR31
QG008-01	B11	HQP009458	NM_000629	IFNAR1
QG008-01	B12	HQP009460	NM_000874	IFNAR2
QG008-01	C01	HQP009469	NM_000416	IFNGR1
QG008-01	C02	HQP009472	NM_005534	IFNGR2
QG008-01	C03	HQP009523	NM_000875	IGF1R
QG008-01	C04	HQP009532	NM_000876	IGF2R
QG008-01	C05	HQP009642	NM_000877	IL1R1
QG008-01	C06	HQP009643	NM_002182	IL1RAP
QG008-01	C07	HQP009650	NM_000417	IL2RA
QG008-01	C08	HQP009658	NM_000878	IL2RB
QG008-01	C09	HQP009659	NM_000206	IL2RG
QG008-01	C10	HQP009661	NM_002183	IL3RA
QG008-01	C11	HQP009664	NM_000418	IL4R
QG008-01	C12	HQP009667	NM_000564	IL5RA
QG008-01	D01	HQP009672	NM_000565	IL6R
QG008-01	D02	HQP009674	NM_002184	IL6ST
QG008-01	D03	HQP009677	NM_002185	IL7R
QG008-01	D04	HQP009679	NM_000634	IL8RA
QG008-01	D05	HQP009681	NM_001557	IL8RB

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG008-01	D06	HQP009683	NM_002186	IL9R
QG008-01	D07	HQP009686	NM_001558	IL10RA
QG008-01	D08	HQP009687	NM_000628	IL10RB
QG008-01	D09	HQP009690	NM_004512	IL11RA
QG008-01	D10	HQP009694	NM_005535	IL12RB1
QG008-01	D11	HQP009696	NM_001559	IL12RB2
QG008-01	D12	HQP009700	NM_001560	IL13RA1
QG008-01	E01	HQP009702	NM_000640	IL13RA2
QG008-01	E02	HQP009710	NM_002189	IL15RA
QG008-01	E03	HQP009716	NM_001561	TNFRSF9
QG008-01	E04	HQP012049	NM_002546	TNFRSF11B
QG008-01	E05	HQP012866	NM_006206	PDGFRA
QG008-01	E06	HQP012871	NM_006207	PDGFRL
QG008-01	E07	HQP012889	NM_002609	PDGFRB
QG008-01	E08	HQP018114	NM_003264	TLR2
QG008-01	E09	HQP018116	NM_138554	TLR4
QG008-01	E10	HQP018148	NM_001065	TNFRSF1A
QG008-01	E11	HQP018149	NM_001066	TNFRSF1B
QG008-01	E12	HQP018800	NM_004633	IL1R2
QG008-01	F01	HQP018802	NM_001008540	CXCR4
QG008-01	F02	HQP021399	NM_003745	SOCS1
QG008-01	F03	HQP021522	NM_003820	TNFRSF14
QG008-01	F04	HQP021536	NM_004195	TNFRSF18
QG008-01	F05	HQP021550	NM_003839	TNFRSF11A
QG008-01	F06	HQP021551	NM_003840	TNFRSF10D
QG008-01	F07	HQP021552	NM_003841	TNFRSF10C
QG008-01	F08	HQP021553	NM_003842	TNFRSF10B
QG008-01	F09	HQP021557	NM_003844	TNFRSF10A
QG008-01	F10	HQP021568	NM_003853	IL18RAP
QG008-01	F11	HQP021569	NM_003854	IL1RL2
QG008-01	F12	HQP021602	NM_003877	SOCS2
QG008-01	G01	HQP021889	NM_003955	SOCS3
QG008-01	G02	HQP021922	NM_003965	CCRL2
QG008-01	G03	HQP022241	NM_003856	IL1RL1
QG008-01	G04	HQP022792	NM_004843	IL27RA
QG008-01	G05	HQP023037	NM_014011	SOCS5
QG008-01	G06	HQP000186	NM_005755	EBI3
QG008-01	G07	HQP000808	NM_006564	CXCR6
QG008-01	G08	HQP000939	NM_006641	CCR9
QG008-01	G09	HQP002207	NM_001837	CCR3
QG008-01	G10	HQP002245	NM_181657	LTB4R

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG008-01	G11	HQP005427	NM_015850	FGFR1
QG008-01	G12	HQP018332	NM_003327	TNFRSF4
QG008-01	H01	HGDC		
QG008-01	H02	HGDC		
QG008-01	H03	HQP006940	NM_002046	GAPDH
QG008-01	H04	HQP016381	NM_001101	ACTB
QG008-01	H05	HQP015171	NM_004048	B2M
QG008-01	H06	HQP006171	NM_012423	RPL13A
QG008-01	H07	HQP009026	NM_000194	HPRT1
QG008-01	H08	HQP054253	NR_003286	RN18S1
QG008-01	H09	RT		
QG008-01	H10	RT		
QG008-01	H11	PCR		
QG008-01	H12	PCR		