

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG048-01	A01	HQP016328	NM_001706	BCL6
QG048-01	A02	HQP018238	NM_000064	C3
QG048-01	A03	HQP018239	NM_004054	C3AR1
QG048-01	A04	HQP018248	NM_007293	C4A
QG048-01	A05	HQP016629	NM_002986	CCL11
QG048-01	A06	HQP016630	NM_005408	CCL13
QG048-01	A07	HQP016635	NM_004590	CCL16
QG048-01	A08	HQP016636	NM_002987	CCL17
QG048-01	A09	HQP016638	NM_006274	CCL19
QG048-01	A10	HQP016621	NM_002982	CCL2
QG048-01	A11	HQP016640	NM_002989	CCL21
QG048-01	A12	HQP016641	NM_002990	CCL22
QG048-01	B01	HQP016642	NM_005064	CCL23
QG048-01	B02	HQP016644	NM_002991	CCL24
QG048-01	B03	HQP016622	NM_002983	CCL3
QG048-01	B04	HQP016625	NM_002984	CCL4
QG048-01	B05	HQP016626	NM_002985	CCL5
QG048-01	B06	HQP016627	NM_006273	CCL7
QG048-01	B07	HQP016628	NM_005623	CCL8
QG048-01	B08	HQP002198	NM_001295	CCR1
QG048-01	B09	HQP002201	NM_000648	CCR2
QG048-01	B10	HQP002209	NM_005508	CCR4
QG048-01	B11	HQP002217	NM_001838	CCR7
QG048-01	B12	HQP022955	NM_001250	CD40
QG048-01	C01	HQP022962	NM_000074	CD40LG
QG048-01	C02	HQP002994	NM_000567	CRP
QG048-01	C03	HQP003149	NM_000757	CSF1
QG048-01	C04	HQP008456	NM_001511	CXCL1
QG048-01	C05	HQP009746	NM_001565	CXCL10
QG048-01	C06	HQP008458	NM_002089	CXCL2
QG048-01	C07	HQP008459	NM_002090	CXCL3
QG048-01	C08	HQP016650	NM_002994	CXCL5
QG048-01	C09	HQP016648	NM_002993	CXCL6
QG048-01	C10	HQP011220	NM_002416	CXCL9
QG048-01	C11	HQP018803	NM_003467	CXCR4
QG048-01	C12	HQP009671	NM_000639	FASLG
QG048-01	D01	HQP006188	NM_005252	FOS
QG048-01	D02	HQP023167	NM_006037	HDAC4
QG048-01	D03	HQP009467	NM_000619	IFNG
QG048-01	D04	HQP009685	NM_000572	IL10
QG048-01	D05	HQP009687	NM_000628	IL10RB

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG048-01	D06	HQP009718	NM_001562	IL18
QG048-01	D07	HQP021568	NM_003853	IL18RAP
QG048-01	D08	HQP009640	NM_000575	IL1A
QG048-01	D09	HQP009641	NM_000576	IL1B
QG048-01	D10	HQP009642	NM_000877	IL1R1
QG048-01	D11	HQP009643	NM_002182	IL1RAP
QG048-01	D12	HQP009645	NM_000577	IL1RN
QG048-01	E01	HQP012161	NM_020525	IL22
QG048-01	E02	HQP001881	NM_052962	IL22RA2
QG048-01	E03	HQP012859	NM_016584	IL23A
QG048-01	E04	HQP003495	NM_144701	IL23R
QG048-01	E05	HQP009670	NM_000600	IL6
QG048-01	E06	HQP009672	NM_000565	IL6R
QG048-01	E07	HQP009678	NM_000584	IL8
QG048-01	E08	HQP009679	NM_000634	IL8RA
QG048-01	E09	HQP009681	NM_001557	IL8RB
QG048-01	E10	HQP009815	NM_000211	ITGB2
QG048-01	E11	HQP010114	NM_000893	KNG1
QG048-01	E12	HQP010907	NM_000595	LTA
QG048-01	F01	HQP006296	NM_015364	LY96
QG048-01	F02	HQP011603	NM_002468	MYD88
QG048-01	F03	HQP011792	NM_004555	NFATC3
QG048-01	F04	HQP011807	NM_003998	NFKB1
QG048-01	F05	HQP011866	NM_000625	NOS2A
QG048-01	F06	HQP008401	NM_000176	NR3C1
QG048-01	F07	HQP021524	NM_003821	RIPK2
QG048-01	F08	HQP001674	NM_001039661	TIRAP
QG048-01	F09	HQP018113	NM_003263	TLR1
QG048-01	F10	HQP018114	NM_003264	TLR2
QG048-01	F11	HQP018115	NM_003265	TLR3
QG048-01	F12	HQP018116	NM_138554	TLR4
QG048-01	G01	HQP018118	NM_003268	TLR5
QG048-01	G02	HQP000415	NM_006068	TLR6
QG048-01	G03	HQP012591	NM_016562	TLR7
QG048-01	G04	HQP018141	NM_000594	TNF
QG048-01	G05	HQP021496	NM_003807	TNFSF14
QG048-01	G06	HQP013496	NM_019009	TOLLIP
QG048-01	G07	HQP003150	NM_172210	CSF1
QG048-01	G08	HQP009026	NM_000194	HPRT1
QG048-01	G09	HQP009644	NM_134470	IL1RAP
QG048-01	G10	HQP009673	NM_181359	IL6R

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG048-01	G11	HQP016643	NM_145898	CCL23
QG048-01	G12	HQP002207	NM_001837	CCR3
QG048-01	H01	HGDC		
QG048-01	H02	HGDC		
QG048-01	H03	HQP006940	NM_002046	GAPDH
QG048-01	H04	HQP016381	NM_001101	ACTB
QG048-01	H05	HQP015171	NM_004048	B2M
QG048-01	H06	HQP006171	NM_012423	RPL13A
QG048-01	H07	HQP009026	NM_000194	HPRT1
QG048-01	H08	HQP054253	NR_003286	RN18S1
QG048-01	H09	RT		
QG048-01	H10	RT		
QG048-01	H11	PCR		
QG048-01	H12	PCR		