

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG074-01	A01	HQP009670	NM_000600	IL6
QG074-01	A02	HQP018141	NM_000594	TNF
QG074-01	A03	HQP011547	NM_005957	MTHFR
QG074-01	A04	HQP013100	NM_000927	ABCB1
QG074-01	A05	HQP009641	NM_000576	IL1B
QG074-01	A06	HQP018342	NM_001071	TYMS
QG074-01	A07	HQP011810	NM_020529	NFKBIA
QG074-01	A08	HQP011807	NM_003998	NFKB1
QG074-01	A09	HQP009685	NM_000572	IL10
QG074-01	A10	HQP009672	NM_000565	IL6R
QG074-01	A11	HQP003809	NM_000769	CYP2C19
QG074-01	A12	HQP018564	NM_005432	XRCC3
QG074-01	B01	HQP018175	NM_000546	TP53
QG074-01	B02	HQP013473	NM_000446	PON1
QG074-01	B03	HQP011555	NM_002454	MTRR
QG074-01	B04	HQP011554	NM_000254	MTR
QG074-01	B05	HQP009781	NM_002460	IRF4
QG074-01	B06	HQP009675	NM_175767	IL6ST
QG074-01	B07	HQP009664	NM_000418	IL4R
QG074-01	B08	HQP009662	NM_000589	IL4
QG074-01	B09	HQP009645	NM_000577	IL1RN
QG074-01	B10	HQP009640	NM_000575	IL1A
QG074-01	B11	HQP009523	NM_000875	IGF1R
QG074-01	B12	HQP005280	NM_021642	FCGR2A
QG074-01	C01	HQP004948	NM_000120	EPHX1
QG074-01	C02	HQP003772	NM_000499	CYP1A1
QG074-01	C03	HQP003499	NM_001037631	CTLA4
QG074-01	C04	HQP000970	NM_006663	PPP1R13L
QG074-01	C05	HQP004317	NM_000903	NQO1
QG074-01	C06	HQP054047	BC008403	HLA-DRB1
QG074-01	C07	HQP023467	NM_000662	NAT1
QG074-01	C08	HQP023168	NM_004356	CD81
QG074-01	C09	HQP023155	NM_000626	CD79B
QG074-01	C10	HQP022955	NM_001250	CD40
QG074-01	C11	HQP022699	NM_006139	CD28
QG074-01	C12	HQP022337	NM_014207	CD5
QG074-01	D01	HQP022316	NM_000616	CD4
QG074-01	D02	HQP021834	NM_005092	TNFSF18
QG074-01	D03	HQP021604	NM_003879	CFLAR
QG074-01	D04	HQP021557	NM_003844	TNFRSF10A
QG074-01	D05	HQP021553	NM_003842	TNFRSF10B

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG074-01	D06	HQP021526	NM_003824	FADD
QG074-01	D07	HQP021522	NM_003820	TNFRSF14
QG074-01	D08	HQP021502	NM_003810	TNFSF10
QG074-01	D09	HQP021462	NM_003789	TRADD
QG074-01	D10	HQP020297	NM_004346	CASP3
QG074-01	D11	HQP019982	NM_001017388	TLR10
QG074-01	D12	HQP019768	NM_130445	COL18A1
QG074-01	E01	HQP018966	NM_001080124	CASP8
QG074-01	E02	HQP018568	NM_021141	XRCC5
QG074-01	E03	HQP018565	NM_003401	XRCC4
QG074-01	E04	HQP018475	NM_001025366	VEGFA
QG074-01	E05	HQP018466	NM_080682	VCAM1
QG074-01	E06	HQP018237	NM_004620	TRAF6
QG074-01	E07	HQP018236	NM_001033910	TRAF5
QG074-01	E08	HQP018233	NM_021138	TRAF2
QG074-01	E09	HQP018149	NM_001066	TNFRSF1B
QG074-01	E10	HQP018113	NM_003263	TLR1
QG074-01	E11	HQP016487	NM_001022	RPS19
QG074-01	E12	HQP016214	NM_006509	RELB
QG074-01	F01	HQP016213	NM_021975	RELA
QG074-01	F02	HQP016208	NM_002908	REL
QG074-01	F03	HQP015598	NM_000963	PTGS2
QG074-01	F04	HQP015194	NM_002791	PSMA6
QG074-01	F05	HQP013633	NM_005037	PPARG
QG074-01	F06	HQP012651	NM_016639	TNFRSF12A
QG074-01	F07	HQP012100	NM_002562	P2RX7
QG074-01	F08	HQP011813	NM_004556	NFKBIE
QG074-01	F09	HQP011808	NM_001077493	NFKB2
QG074-01	F10	HQP011325	NM_019899	ABCC1
QG074-01	F11	HQP011309	NM_000250	MPO
QG074-01	F12	HQP010613	NM_002312	LIG4
QG074-01	G01	HQP009851	NM_000215	JAK3
QG074-01	G02	HQP009788	NM_005544	IRS1
QG074-01	G03	HQP009718	NM_001562	IL18
QG074-01	G04	HQP009716	NM_001561	TNFRSF9
QG074-01	G05	HQP009708	NM_000585	IL15
QG074-01	G06	HQP009697	NM_002188	IL13
QG074-01	G07	HQP009696	NM_001559	IL12RB2
QG074-01	G08	HQP009694	NM_005535	IL12RB1
QG074-01	G09	HQP009693	NM_002187	IL12B
QG074-01	G10	HQP009692	NM_000882	IL12A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG074-01	G11	HQP009687	NM_000628	IL10RB
QG074-01	G12	HQP009681	NM_001557	IL8RB
QG074-01	H01	HGDC		
QG074-01	H02	HGDC		
QG074-01	H03	HQP006940	NM_002046	GAPDH
QG074-01	H04	HQP016381	NM_001101	ACTB
QG074-01	H05	HQP015171	NM_004048	B2M
QG074-01	H06	HQP006171	NM_012423	RPL13A
QG074-01	H07	HQP009026	NM_000194	HPRT1
QG074-01	H08	HQP054253	NR_003286	RN18S1
QG074-01	H09	RT		
QG074-01	H10	RT		
QG074-01	H11	PCR		
QG074-01	H12	PCR		