

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG027-01	A01	HQP002833	NM_000675	ADORA2A
QG027-01	A02	HQP018304	NM_001735	C5
QG027-01	A03	HQP018362	NM_000562	C8A
QG027-01	A04	HQP020046	NM_004345	CAMP
QG027-01	A05	HQP020208	NM_033292	CASP1
QG027-01	A06	HQP020332	NM_001225	CASP4
QG027-01	A07	HQP016621	NM_002982	CCL2
QG027-01	A08	HQP022490	NM_000591	CD14
QG027-01	A09	HQP022129	NM_001766	CD1D
QG027-01	A10	HQP003948	NM_000574	CD55
QG027-01	A11	HQP001708	NM_001278	CHUK
QG027-01	A12	HQP019828	NM_130386	COLEC12
QG027-01	B01	HQP002994	NM_000567	CRP
QG027-01	B02	HQP018803	NM_003467	CXCR4
QG027-01	B03	HQP003741	NM_000397	CYBB
QG027-01	B04	HQP004193	NM_004942	DEFB4
QG027-01	B05	HQP004353	NM_004406	DMBT1
QG027-01	B06	HQP006022	NM_002026	FN1
QG027-01	B07	HQP008898	NM_002133	HMOX1
QG027-01	B08	HQP009419	NM_024013	IFNA1
QG027-01	B09	HQP009463	NM_002176	IFNB1
QG027-01	B10	HQP009469	NM_000416	IFNGR1
QG027-01	B11	HQP009472	NM_005534	IFNGR2
QG027-01	B12	HQP009639	NM_001556	IKKBK
QG027-01	C01	HQP009685	NM_000572	IL10
QG027-01	C02	HQP009696	NM_001559	IL12RB2
QG027-01	C03	HQP009640	NM_000575	IL1A
QG027-01	C04	HQP009641	NM_000576	IL1B
QG027-01	C05	HQP007289	NM_012275	IL1F5
QG027-01	C06	HQP007573	NM_014440	IL1F6
QG027-01	C07	HQP007567	NM_173178	IL1F8
QG027-01	C08	HQP015063	NM_019618	IL1F9
QG027-01	C09	HQP009642	NM_000877	IL1R1
QG027-01	C10	HQP018800	NM_004633	IL1R2
QG027-01	C11	HQP009643	NM_002182	IL1RAP
QG027-01	C12	HQP007197	NM_017416	IL1RAPL2
QG027-01	D01	HQP021569	NM_003854	IL1RL2
QG027-01	D02	HQP009645	NM_000577	IL1RN
QG027-01	D03	HQP009670	NM_000600	IL6
QG027-01	D04	HQP009773	NM_001569	IRAK1
QG027-01	D05	HQP009776	NM_001570	IRAK2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG027-01	D06	HQP009778	NM_002198	IRF1
QG027-01	D07	HQP010489	NM_002289	LALBA
QG027-01	D08	HQP010560	NM_004139	LBP
QG027-01	D09	HQP010919	NM_002343	LTF
QG027-01	D10	HQP006296	NM_015364	LY96
QG027-01	D11	HQP010937	NM_000239	LYZ
QG027-01	D12	HQP003133	NM_001315	MAPK14
QG027-01	E01	HQP014886	NM_002750	MAPK8
QG027-01	E02	HQP011219	NM_002415	MIF
QG027-01	E03	HQP011603	NM_002468	MYD88
QG027-01	E04	HQP011694	NM_000631	NCF4
QG027-01	E05	HQP011807	NM_003998	NFKB1
QG027-01	E06	HQP053985	NM_002502	NFKB2
QG027-01	E07	HQP011810	NM_020529	NFKBIA
QG027-01	E08	HQP015999	NM_021209	NLRC4
QG027-01	E09	HQP011866	NM_000625	NOS2A
QG027-01	E10	HQP001682	NM_052890	PGLYRP2
QG027-01	E11	HQP013662	NM_002704	PPBP
QG027-01	E12	HQP015041	NM_000312	PROC
QG027-01	F01	HQP015524	NM_000952	PTAFR
QG027-01	F02	HQP016551	NM_005621	S100A12
QG027-01	F03	HQP013122	NM_000295	SERPINA1
QG027-01	F04	HQP012154	NM_000602	SERPINE1
QG027-01	F05	HQP016980	NM_003019	SFTPD
QG027-01	F06	HQP018044	NM_000660	TGFB1
QG027-01	F07	HQP018113	NM_003263	TLR1
QG027-01	F08	HQP054024	NM_030956	TLR10
QG027-01	F09	HQP018114	NM_003264	TLR2
QG027-01	F10	HQP018115	NM_003265	TLR3
QG027-01	F11	HQP018116	NM_138554	TLR4
QG027-01	F12	HQP000415	NM_006068	TLR6
QG027-01	G01	HQP012625	NM_138636	TLR8
QG027-01	G02	HQP013388	NM_017442	TLR9
QG027-01	G03	HQP018141	NM_000594	TNF
QG027-01	G04	HQP018148	NM_001065	TNFRSF1A
QG027-01	G05	HQP013496	NM_019009	TOLLIP
QG027-01	G06	HQP018237	NM_004620	TRAF6
QG027-01	G07	HQP020210	NM_033294	CASP1
QG027-01	G08	HQP009026	NM_000194	HPRT1
QG027-01	G09	HQP006023	NM_054034	FN1
QG027-01	G10	HQP009644	NM_134470	IL1RAP

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG027-01	G11	HQP002207	NM_001837	CCR3
QG027-01	G12	HQP001683	NM_052891	PGLYRP3
QG027-01	H01	HGDC		
QG027-01	H02	HGDC		
QG027-01	H03	HQP006940	NM_002046	GAPDH
QG027-01	H04	HQP016381	NM_001101	ACTB
QG027-01	H05	HQP015171	NM_004048	B2M
QG027-01	H06	HQP006171	NM_012423	RPL13A
QG027-01	H07	HQP009026	NM_000194	HPRT1
QG027-01	H08	HQP054253	NR_003286	RN18S1
QG027-01	H09	RT		
QG027-01	H10	RT		
QG027-01	H11	PCR		
QG027-01	H12	PCR		