

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG032-01	A01	HQP013694	NM_021130	PPIA
QG032-01	A02	HQP008650	NM_013444	UBQLN2
QG032-01	A03	HQP016381	NM_001101	ACTB
QG032-01	A04	HQP015171	NM_004048	B2M
QG032-01	A05	HQP008515	NM_013314	BLNK
QG032-01	A06	HQP002198	NM_001295	CCR1
QG032-01	A07	HQP022722	NM_005191	CD80
QG032-01	A08	HQP003159	NM_000758	CSF2
QG032-01	A09	HQP003173	NM_000759	CSF3
QG032-01	A10	HQP016649	NM_005409	CXCL11
QG032-01	A11	HQP008458	NM_002089	CXCL2
QG032-01	A12	HQP008459	NM_002090	CXCL3
QG032-01	B01	HQP011220	NM_002416	CXCL9
QG032-01	B02	HQP008583	NM_000181	GUSB
QG032-01	B03	HQP009026	NM_000194	HPRT1
QG032-01	B04	HQP000670	NM_006417	IFI44
QG032-01	B05	HQP009416	NM_001031683	IFIT3
QG032-01	B06	HQP009685	NM_000572	IL10
QG032-01	B07	HQP009708	NM_000585	IL15
QG032-01	B08	HQP009641	NM_000576	IL1B
QG032-01	B09	HQP012859	NM_016584	IL23A
QG032-01	B10	HQP009785	NM_001572	IRF7
QG032-01	B11	HQP009789	NM_002201	ISG20
QG032-01	B12	HQP001164	NM_006863	LILRA1
QG032-01	C01	HQP002725	NM_005204	MAP3K8
QG032-01	C02	HQP011257	NM_002422	MMP3
QG032-01	C03	HQP011263	NM_004994	MMP9
QG032-01	C04	HQP011582	NM_002462	MX1
QG032-01	C05	HQP011983	NM_001032409	OAS1
QG032-01	C06	HQP011988	NM_001032731	OAS2
QG032-01	C07	HQP012857	NM_033016	PDGFB
QG032-01	C08	HQP006171	NM_012423	RPL13A
QG032-01	C09	HQP017764	NM_007315	STAT1
QG032-01	C10	HQP018113	NM_003263	TLR1
QG032-01	C11	HQP018115	NM_003265	TLR3
QG032-01	C12	HQP012591	NM_016562	TLR7
QG032-01	D01	HQP001496	NM_017414	USP18
QG032-01	D02	HQP006359	NM_000698	ALOX5
QG032-01	D03	HQP019227	NM_024701	ASB13
QG032-01	D04	HQP001840	NM_138456	BATF2
QG032-01	D05	HQP018239	NM_004054	C3AR1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG032-01	D06	HQP022548	NM_004244	CD163
QG032-01	D07	HQP023056	NM_000611	CD59
QG032-01	D08	HQP022747	NM_175862	CD86
QG032-01	D09	HQP008456	NM_001511	CXCL1
QG032-01	D10	HQP006239	NM_014314	DDX58
QG032-01	D11	HQP014498	NM_017631	FLJ20035
QG032-01	D12	HQP004173	NM_004084	DEFA1
QG032-01	E01	HQP018936	NM_024119	LGP2
QG032-01	E02	HQP004712	NM_172373	ELF1
QG032-01	E03	HQP004735	NM_001421	ELF4
QG032-01	E04	HQP005022	NM_001987	ETV6
QG032-01	E05	HQP005251	NM_000566	FCGR1A
QG032-01	E06	HQP006244	NM_001005738	FPRL1
QG032-01	E07	HQP001794	NM_052941	GBP4
QG032-01	E08	HQP021850	NM_177977	HAP1
QG032-01	E09	HQP008843	NM_000189	HK2
QG032-01	E10	HQP009097	NM_007355	HSP90AB1
QG032-01	E11	HQP010070	NM_002253	KDR
QG032-01	E12	HQP010563	NM_002296	LBR
QG032-01	F01	HQP022587	NM_004789	LHX2
QG032-01	F02	HQP022648	NM_006033	LIPG
QG032-01	F03	HQP011149	NM_005587	MEF2A
QG032-01	F04	HQP011687	NM_002485	NBN
QG032-01	F05	HQP000603	NM_006540	NCOA2
QG032-01	F06	HQP011864	NM_007363	NONO
QG032-01	F07	HQP012662	NM_005018	PDCD1
QG032-01	F08	HQP013072	NM_004566	PFKFB3
QG032-01	F09	HQP015292	NM_148919	PSMB8
QG032-01	F10	HQP019923	NM_030916	PVRL4
QG032-01	F11	HQP003517	NM_152501	PYHIN1
QG032-01	F12	HQP017647	NM_003113	SP100
QG032-01	G01	HQP023244	NM_006038	SPATA2
QG032-01	G02	HQP022852	NM_004863	SPTLC2
QG032-01	G03	HQP017928	NM_003194	TBP
QG032-01	G04	HQP018041	NM_003234	TFRC
QG032-01	G05	HQP018116	NM_138554	TLR4
QG032-01	G06	HQP023251	NM_033220	TRIM14
QG032-01	G07	HQP023408	NM_006313	USP15
QG032-01	G08	HQP013661	NM_199139	XAF1
QG032-01	G09	HQP019822	NM_030776	ZBP1
QG032-01	G10	HQP020411	NM_032036	FAM14A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG032-01	G11	HQP015598	NM_000963	PTGS2
QG032-01	G12	HQP006424	NM_145659	IL27
QG032-01	H01	HGDC		
QG032-01	H02	HGDC		
QG032-01	H03	HQP006940	NM_002046	GAPDH
QG032-01	H04	HQP016381	NM_001101	ACTB
QG032-01	H05	HQP015171	NM_004048	B2M
QG032-01	H06	HQP006171	NM_012423	RPL13A
QG032-01	H07	HQP009026	NM_000194	HPRT1
QG032-01	H08	HQP054253	NR_003286	RN18S1
QG032-01	H09	RT		
QG032-01	H10	RT		
QG032-01	H11	PCR		
QG032-01	H12	PCR		