

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG034-01	A01	HQP008678	NM_000014	A2M
QG034-01	A02	HQP000726	NM_020979	SH2B2
QG034-01	A03	HQP016238	NM_138578	BCL2L1
QG034-01	A04	HQP016204	NM_053056	CCND1
QG034-01	A05	HQP000331	NM_000389	CDKN1A
QG034-01	A06	HQP002988	NM_016823	CRK
QG034-01	A07	HQP002994	NM_000567	CRP
QG034-01	A08	HQP003158	NM_005211	CSF1R
QG034-01	A09	HQP003170	NM_000395	CSF2RB
QG034-01	A10	HQP011220	NM_002416	CXCL9
QG034-01	A11	HQP004605	NM_005228	EGFR
QG034-01	A12	HQP004958	NM_000121	EPOR
QG034-01	B01	HQP005052	NM_000506	F2
QG034-01	B02	HQP005053	NM_001992	F2R
QG034-01	B03	HQP009651	NM_000043	FAS
QG034-01	B04	HQP005238	NM_002001	FCER1A
QG034-01	B05	HQP005251	NM_000566	FCGR1A
QG034-01	B06	HQP023013	NM_005101	ISG15
QG034-01	B07	HQP007166	NM_002051	GATA3
QG034-01	B08	HQP007221	NM_002053	GBP1
QG034-01	B09	HQP007395	NM_000163	GHR
QG034-01	B10	HQP009458	NM_000629	IFNAR1
QG034-01	B11	HQP009467	NM_000619	IFNG
QG034-01	B12	HQP009469	NM_000416	IFNGR1
QG034-01	C01	HQP009686	NM_001558	IL10RA
QG034-01	C02	HQP012157	NM_018724	IL20
QG034-01	C03	HQP009650	NM_000417	IL2RA
QG034-01	C04	HQP009659	NM_000206	IL2RG
QG034-01	C05	HQP009662	NM_000589	IL4
QG034-01	C06	HQP009664	NM_000418	IL4R
QG034-01	C07	HQP009674	NM_002184	IL6ST
QG034-01	C08	HQP009764	NM_000208	INSR
QG034-01	C09	HQP009778	NM_002198	IRF1
QG034-01	C10	HQP009849	NM_002227	JAK1
QG034-01	C11	HQP009850	NM_004972	JAK2
QG034-01	C12	HQP009851	NM_000215	JAK3
QG034-01	D01	HQP009853	NM_002228	JUN
QG034-01	D02	HQP009854	NM_002229	JUNB
QG034-01	D03	HQP011257	NM_002422	MMP3
QG034-01	D04	HQP011308	NM_005373	MPL
QG034-01	D05	HQP011597	NM_002467	MYC

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG034-01	D06	HQP011807	NM_003998	NFKB1
QG034-01	D07	HQP011866	NM_000625	NOS2A
QG034-01	D08	HQP008401	NM_000176	NR3C1
QG034-01	D09	HQP011984	NM_002534	OAS1
QG034-01	D10	HQP012866	NM_006206	PDGFRA
QG034-01	D11	HQP021271	NM_016166	PIAS1
QG034-01	D12	HQP022005	NM_004671	PIAS2
QG034-01	E01	HQP014147	NM_014225	PPP2R1A
QG034-01	E02	HQP015027	NM_000949	PRLR
QG034-01	E03	HQP015828	NM_002827	PTPN1
QG034-01	E04	HQP015908	NM_002838	PTPRC
QG034-01	E05	HQP020528	NM_175077	SLA2
QG034-01	E06	HQP054006	NM_005900	SMAD1
QG034-01	E07	HQP054007	NM_005901	SMAD2
QG034-01	E08	HQP010960	NM_005902	SMAD3
QG034-01	E09	HQP010961	NM_005359	SMAD4
QG034-01	E10	HQP054008	NM_005903	SMAD5
QG034-01	E11	HQP021399	NM_003745	SOCS1
QG034-01	E12	HQP021602	NM_003877	SOCS2
QG034-01	F01	HQP021889	NM_003955	SOCS3
QG034-01	F02	HQP002181	NM_080867	SOCS4
QG034-01	F03	HQP054032	NM_144949	SOCS5
QG034-01	F04	HQP017640	NM_138473	SP1
QG034-01	F05	HQP017663	NM_003120	SPI1
QG034-01	F06	HQP017696	NM_005417	SRC
QG034-01	F07	HQP019637	NM_003473	STAM
QG034-01	F08	HQP017764	NM_007315	STAT1
QG034-01	F09	HQP017766	NM_005419	STAT2
QG034-01	F10	HQP017767	NM_003150	STAT3
QG034-01	F11	HQP017770	NM_003151	STAT4
QG034-01	F12	HQP017771	NM_003152	STAT5A
QG034-01	G01	HQP017774	NM_012448	STAT5B
QG034-01	G02	HQP017775	NM_003153	STAT6
QG034-01	G03	HQP018340	NM_003331	TYK2
QG034-01	G04	HQP018570	NM_003403	YY1
QG034-01	G05	HQP009026	NM_000194	HPRT1
QG034-01	G06	HQP009653	NM_152872	FAS
QG034-01	G07	HQP009675	NM_175767	IL6ST
QG034-01	G08	HQP015909	NM_080921	PTPRC
QG034-01	G09	HQP017769	NM_213662	STAT3
QG034-01	G10	HQP000623	NM_005194	CEBPB

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG034-01	G11	HQP012081	NM_020530	OSM
QG034-01	G12	HQP000334	NM_005861	STUB1
QG034-01	H01	HGDC		
QG034-01	H02	HGDC		
QG034-01	H03	HQP006940	NM_002046	GAPDH
QG034-01	H04	HQP016381	NM_001101	ACTB
QG034-01	H05	HQP015171	NM_004048	B2M
QG034-01	H06	HQP006171	NM_012423	RPL13A
QG034-01	H07	HQP009026	NM_000194	HPRT1
QG034-01	H08	HQP054253	NR_003286	RN18S1
QG034-01	H09	RT		
QG034-01	H10	RT		
QG034-01	H11	PCR		
QG034-01	H12	PCR		