

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG025-01	A01	HQP010996	NM_022438	MAL
QG025-01	A02	HQP015909	NM_080921	PTPRC
QG025-01	A03	HQP009675	NM_175767	IL6ST
QG025-01	A04	HQP023134	NM_014707	HDAC9
QG025-01	A05	HQP003150	NM_172210	CSF1
QG025-01	A06	HQP009026	NM_000194	HPRT1
QG025-01	A07	HQP022974	NM_001001390	CD44
QG025-01	A08	HQP021771	NM_033131	WNT3A
QG025-01	A09	HQP018481	NM_003376	VEGFA
QG025-01	A10	HQP018461	NM_005428	VAV1
QG025-01	A11	HQP000139	NM_006778	TRIM10
QG025-01	A12	HQP018116	NM_138554	TLR4
QG025-01	B01	HQP018115	NM_003265	TLR3
QG025-01	B02	HQP018012	NM_000459	TEK
QG025-01	B03	HQP017895	NM_003189	TAL1
QG025-01	B04	HQP015757	NM_020860	STIM2
QG025-01	B05	HQP017767	NM_003150	STAT3
QG025-01	B06	HQP017764	NM_007315	STAT1
QG025-01	B07	HQP017673	NM_000582	SPP1
QG025-01	B08	HQP054032	NM_144949	SOCS5
QG025-01	B09	HQP022694	NM_022754	SFXN1
QG025-01	B10	HQP021347	NM_001754	RUNX1
QG025-01	B11	HQP009574	NM_005349	RBPJ
QG025-01	B12	HQP015908	NM_002838	PTPRC
QG025-01	C01	HQP013015	NM_000442	PECAM1
QG025-01	C02	HQP012212	NM_016734	PAX5
QG025-01	C03	HQP011875	NM_024408	NOTCH2
QG025-01	C04	HQP011873	NM_017617	NOTCH1
QG025-01	C05	HQP011866	NM_000625	NOS2A
QG025-01	C06	HQP011263	NM_004994	MMP9
QG025-01	C07	HQP010995	NM_002371	MAL
QG025-01	C08	HQP010868	NM_006152	LRMP
QG025-01	C09	HQP012480	NM_016269	LEF1
QG025-01	C10	HQP011205	NM_003994	KITLG
QG025-01	C11	HQP010099	NM_000222	KIT
QG025-01	C12	HQP010070	NM_002253	KDR
QG025-01	D01	HQP009847	NM_002226	JAG2
QG025-01	D02	HQP004470	NM_000214	JAG1
QG025-01	D03	HQP009743	NM_002192	INHBA
QG025-01	D04	HQP009742	NM_002191	INHA
QG025-01	D05	HQP009674	NM_002184	IL6ST

QG025-01	D06	HQP002747	NM_139017	IL31RA
QG025-01	D07	HQP012157	NM_018724	IL20
QG025-01	D08	HQP009649	NM_000586	IL2
QG025-01	D09	HQP009640	NM_000575	IL1A
QG025-01	D10	HQP009693	NM_002187	IL12B
QG025-01	D11	HQP009685	NM_000572	IL10
QG025-01	D12	HQP012862	NM_016596	HDAC7A
QG025-01	E01	HQP000024	NM_005474	HDAC5
QG025-01	E02	HQP023167	NM_006037	HDAC4
QG025-01	E03	HQP007152	NM_032638	GATA2
QG025-01	E04	HQP007146	NM_002049	GATA1
QG025-01	E05	HQP012676	NM_015892	GALNAC4S-6S T
QG025-01	E06	HQP020133	NM_003505	FZD1
QG025-01	E07	HQP005022	NM_001987	ETV6
QG025-01	E08	HQP005014	NM_005238	ETS1
QG025-01	E09	HQP008111	NM_005618	DLL1
QG025-01	E10	HQP003159	NM_000758	CSF2
QG025-01	E11	HQP003149	NM_000757	CSF1
QG025-01	E12	HQP000656	NM_001806	CEBPG
QG025-01	F01	HQP022746	NM_006889	CD86
QG025-01	F02	HQP022722	NM_005191	CD80
QG025-01	F03	HQP022972	NM_000610	CD44
QG025-01	F04	HQP022316	NM_000616	CD4
QG025-01	F05	HQP022256	NM_000073	CD3G
QG025-01	F06	HQP022212	NM_000732	CD3D
QG025-01	F07	HQP022667	NM_001242	CD27
QG025-01	F08	HQP022190	NM_001767	CD2
QG025-01	F09	HQP022129	NM_001766	CD1D
QG025-01	F10	HQP022490	NM_000591	CD14
QG025-01	F11	HQP002198	NM_001295	CCR1
QG025-01	F12	HQP008515	NM_013314	BLNK
QG025-01	G01	HQP022017	NM_004674	ASH2L
QG025-01	G02	HQP009024	NM_000038	APC
QG025-01	G03	HQP000648	NM_001805	CEBPE
QG025-01	G04	HQP010701	NM_005574	LMO2
QG025-01	G05	HQP008097	NM_001146	ANGPT1
QG025-01	G06	HQP021407	NM_001755	CBFB
QG025-01	G07	HQP021521	NM_006016	CD164
QG025-01	G08	HQP022813	NM_001773	CD34
QG025-01	G09	HQP005902	NM_001459	FLT3LG
QG025-01	G10	HQP020928	NM_032664	FUT10
QG025-01	G11	HQP023138	NM_178425	HDAC9

QG025-01	G12	HQP009688	NM_000641	IL11
QG025-01	H01	HGDC		
QG025-01	H02	HGDC		
QG025-01	H03	HQP006940	NM_002046	GAPDH
QG025-01	H04	HQP016381	NM_001101	ACTB
QG025-01	H05	HQP015171	NM_004048	B2M
QG025-01	H06	HQP006171	NM_012423	RPL13A
QG025-01	H07	HQP009026	NM_000194	HPRT1
QG025-01	H08	HQP054253	NR_003286	RN18S1
QG025-01	H09	RT		
QG025-01	H10	RT		
QG025-01	H11	PCR		
QG025-01	H12	PCR		