

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG023-01	A01	HQP021191	NM_033034	TRIM5
QG023-01	A02	HQP011784	NM_006162	NFATC1
QG023-01	A03	HQP005231	NM_133273	FCAR
QG023-01	A04	HQP022974	NM_001001390	CD44
QG023-01	A05	HQP020550	NM_033358	CASP8
QG023-01	A06	HQP009026	NM_000194	HPRT1
QG023-01	A07	HQP004752	NM_001006666	APOBEC3F
QG023-01	A08	HQP018570	NM_003403	YY1
QG023-01	A09	HQP018561	NM_003400	XPO1
QG023-01	A10	HQP007577	NM_013245	VPS4A
QG023-01	A11	HQP018269	NM_006292	TSG101
QG023-01	A12	HQP021502	NM_003810	TNFSF10
QG023-01	B01	HQP018149	NM_001066	TNFRSF1B
QG023-01	B02	HQP018141	NM_000594	TNF
QG023-01	B03	HQP018044	NM_000660	TGFB1
QG023-01	B04	HQP018029	NM_005653	TFCP2
QG023-01	B05	HQP017767	NM_003150	STAT3
QG023-01	B06	HQP017764	NM_007315	STAT1
QG023-01	B07	HQP017526	NM_003073	SMARCB1
QG023-01	B08	HQP017501	NM_003064	SLPI
QG023-01	B09	HQP011620	NM_000488	SERPINC1
QG023-01	B10	HQP013122	NM_000295	SERPINA1
QG023-01	B11	HQP016745	NM_000655	SELL
QG023-01	B12	HQP016172	NM_005611	RBL2
QG023-01	C01	HQP005093	NM_004103	PTK2B
QG023-01	C02	HQP012152	NM_002574	PRDX1
QG023-01	C03	HQP013694	NM_021130	PPIA
QG023-01	C04	HQP011810	NM_020529	NFKBIA
QG023-01	C05	HQP011788	NM_172390	NFATC1
QG023-01	C06	HQP011077	NM_000242	MBL2
QG023-01	C07	HQP010915	NM_002342	LTBR
QG023-01	C08	HQP010112	NM_002262	KLRD1
QG023-01	C09	HQP009779	NM_002199	IRF2
QG023-01	C10	HQP009778	NM_002198	IRF1
QG023-01	C11	HQP009678	NM_000584	IL8
QG023-01	C12	HQP009649	NM_000586	IL2
QG023-01	D01	HQP009641	NM_000576	IL1B
QG023-01	D02	HQP009693	NM_002187	IL12B
QG023-01	D03	HQP009685	NM_000572	IL10
QG023-01	D04	HQP009467	NM_000619	IFNG
QG023-01	D05	HQP009463	NM_002176	IFNB1

QG023-01	D06	HQP009419	NM_024013	IFNA1
QG023-01	D07	HQP007681	NM_014500	HTATSF1
QG023-01	D08	HQP008739	NM_002110	HCK
QG023-01	D09	HQP004125	NM_001924	GADD45A
QG023-01	D10	HQP006188	NM_005252	FOS
QG023-01	D11	HQP004897	NM_001429	EP300
QG023-01	D12	HQP004689	NM_001972	ELA2
QG023-01	E01	HQP018803	NM_003467	CXCR4
QG023-01	E02	HQP016669	NM_000609	CXCL12
QG023-01	E03	HQP016652	NM_002996	CX3CL1
QG023-01	E04	HQP002921	NM_004380	CREBBP
QG023-01	E05	HQP001117	NM_006833	COPS6
QG023-01	E06	HQP000623	NM_005194	CEBPB
QG023-01	E07	HQP000331	NM_000389	CDKN1A
QG023-01	E08	HQP000321	NM_001261	CDK9
QG023-01	E09	HQP000285	NM_001799	CDK7
QG023-01	E10	HQP023094	NM_001781	CD69
QG023-01	E11	HQP022972	NM_000610	CD44
QG023-01	E12	HQP022316	NM_000616	CD4
QG023-01	F01	HQP022295	NM_000734	CD247
QG023-01	F02	HQP008808	NM_021155	CD209
QG023-01	F03	HQP002210	NM_000579	CCR5
QG023-01	F04	HQP002209	NM_005508	CCR4
QG023-01	F05	HQP002201	NM_000648	CCR2
QG023-01	F06	HQP016628	NM_005623	CCL8
QG023-01	F07	HQP016626	NM_002985	CCL5
QG023-01	F08	HQP016625	NM_002984	CCL4
QG023-01	F09	HQP016621	NM_002982	CCL2
QG023-01	F10	HQP006127	NM_012117	CBX5
QG023-01	F11	HQP020297	NM_004346	CASP3
QG023-01	F12	HQP021753	NM_033637	BTRC
QG023-01	G01	HQP016211	NM_000633	BCL2
QG023-01	G02	HQP015964	NM_004324	BAX
QG023-01	G03	HQP015538	NM_004322	BAD
QG023-01	G04	HQP016317	NM_021822	APOBEC3G
QG023-01	G05	HQP004753	NM_145298	APOBEC3F
QG023-01	G06	HQP054029	NM_080649	APEX1
QG023-01	G07	HQP020548	NM_001228	CASP8
QG023-01	G08	HQP002207	NM_001837	CCR3
QG023-01	G09	HQP021575	NM_003860	BANF1
QG023-01	G10	HQP023129	NM_004355	CD74
QG023-01	G11	HQP002890	NM_001877	CR2
QG023-01	G12	HQP005227	NM_002000	FCAR

QG023-01	H01	HGDC		
QG023-01	H02	HGDC		
QG023-01	H03	HQP006940	NM_002046	GAPDH
QG023-01	H04	HQP016381	NM_001101	ACTB
QG023-01	H05	HQP015171	NM_004048	B2M
QG023-01	H06	HQP006171	NM_012423	RPL13A
QG023-01	H07	HQP009026	NM_000194	HPRT1
QG023-01	H08	HQP054253	NR_003286	RN18S1
QG023-01	H09	RT		
QG023-01	H10	RT		
QG023-01	H11	PCR		
QG023-01	H12	PCR		