

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG029-01	A01	HQP009467	NM_000619	IFNG
QG029-01	A02	HQP009685	NM_000572	IL10
QG029-01	A03	HQP009649	NM_000586	IL2
QG029-01	A04	HQP008866	NM_019111	HLA-DRA
QG029-01	A05	HQP009670	NM_000600	IL6
QG029-01	A06	HQP021502	NM_003810	TNFSF10
QG029-01	A07	HQP009641	NM_000576	IL1B
QG029-01	A08	HQP017767	NM_003150	STAT3
QG029-01	A09	HQP009693	NM_002187	IL12B
QG029-01	A10	HQP018044	NM_000660	TGFB1
QG029-01	A11	HQP017899	NM_000593	TAP1
QG029-01	A12	HQP011220	NM_002416	CXCL9
QG029-01	B01	HQP011582	NM_002462	MX1
QG029-01	B02	HQP018340	NM_003331	TYK2
QG029-01	B03	HQP009850	NM_004972	JAK2
QG029-01	B04	HQP005052	NM_000506	F2
QG029-01	B05	HQP009746	NM_001565	CXCL10
QG029-01	B06	HQP009463	NM_002176	IFNB1
QG029-01	B07	HQP021271	NM_016166	PIAS1
QG029-01	B08	HQP011597	NM_002467	MYC
QG029-01	B09	HQP014948	NM_002759	EIF2AK2
QG029-01	B10	HQP008745	NM_004964	HDAC1
QG029-01	B11	HQP023021	NM_014002	IKBKE
QG029-01	B12	HQP003420	NM_182919	TICAM1
QG029-01	C01	HQP018175	NM_000546	TP53
QG029-01	C02	HQP011211	NM_000246	CIITA
QG029-01	C03	HQP016131	NM_000321	RB1
QG029-01	C04	HQP009469	NM_000416	IFNGR1
QG029-01	C05	HQP009664	NM_000418	IL4R
QG029-01	C06	HQP009687	NM_000628	IL10RB
QG029-01	C07	HQP009408	NM_001548	IFIT1
QG029-01	C08	HQP009780	NM_001571	IRF3
QG029-01	C09	HQP007221	NM_002053	GBP1
QG029-01	C10	HQP009251	NM_002163	IRF8
QG029-01	C11	HQP009674	NM_002184	IL6ST
QG029-01	C12	HQP009778	NM_002198	IRF1
QG029-01	D01	HQP009779	NM_002199	IRF2
QG029-01	D02	HQP009849	NM_002227	JAK1
QG029-01	D03	HQP009781	NM_002460	IRF4
QG029-01	D04	HQP011603	NM_002468	MYD88
QG029-01	D05	HQP017729	NM_003141	TRIM21

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG029-01	D06	HQP017770	NM_003151	STAT4
QG029-01	D07	HQP017766	NM_005419	STAT2
QG029-01	D08	HQP009379	NM_005531	IFI16
QG029-01	D09	HQP009472	NM_005534	IFNGR2
QG029-01	D10	HQP009784	NM_006147	IRF6
QG029-01	D11	HQP000520	NM_006332	IFI30
QG029-01	D12	HQP000670	NM_006417	IFI44
QG029-01	E01	HQP017764	NM_007315	STAT1
QG029-01	E02	HQP012591	NM_016562	TLR7
QG029-01	E03	HQP015659	NM_020746	VISA
QG029-01	E04	HQP016303	NM_021133	RNASEL
QG029-01	E05	HQP016808	NM_022168	IFIH1
QG029-01	E06	HQP009651	NM_000043	FAS
QG029-01	E07	HQP017673	NM_000582	SPP1
QG029-01	E08	HQP009458	NM_000629	IFNAR1
QG029-01	E09	HQP009460	NM_000874	IFNAR2
QG029-01	E10	HQP000471	NM_001025107	ADAR
QG029-01	E11	HQP009416	NM_001031683	IFIT3
QG029-01	E12	HQP011983	NM_001032409	OAS1
QG029-01	F01	HQP011988	NM_001032731	OAS2
QG029-01	F02	HQP004683	NM_001042559	EIF4G2
QG029-01	F03	HQP009703	NM_001077397	IRF2BP2
QG029-01	F04	HQP017662	NM_001080547	SPI1
QG029-01	F05	HQP020207	NM_001223	CASP1
QG029-01	F06	HQP009403	NM_001547	IFIT2
QG029-01	F07	HQP009785	NM_001572	IRF7
QG029-01	F08	HQP008849	NM_002116	HLA-A
QG029-01	F09	HQP009427	NM_002169	IFNA5
QG029-01	F10	HQP009446	NM_002173	IFNA16
QG029-01	F11	HQP009782	NM_002200	IRF5
QG029-01	F12	HQP009789	NM_002201	ISG20
QG029-01	G01	HQP015311	NM_002800	PSMB9
QG029-01	G02	HQP021142	NM_003641	IFITM1
QG029-01	G03	HQP021305	NM_003690	PRKRA
QG029-01	G04	HQP021399	NM_003745	SOCS1
QG029-01	G05	HQP014969	NM_004705	PRKRIR
QG029-01	G06	HQP023013	NM_005101	ISG15
QG029-01	G07	HQP016649	NM_005409	CXCL11
QG029-01	G08	HQP009385	NM_005532	IFI27
QG029-01	G09	HQP009393	NM_005533	IFI35
QG029-01	G10	HQP015499	NM_006263	PSME1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG029-01	G11	HQP007080	NM_015649	IRF2BP1
QG029-01	G12	HQP001496	NM_017414	USP18
QG029-01	H01	HGDC		
QG029-01	H02	HGDC		
QG029-01	H03	HQP006940	NM_002046	GAPDH
QG029-01	H04	HQP016381	NM_001101	ACTB
QG029-01	H05	HQP015171	NM_004048	B2M
QG029-01	H06	HQP006171	NM_012423	RPL13A
QG029-01	H07	HQP009026	NM_000194	HPRT1
QG029-01	H08	HQP054253	NR_003286	RN18S1
QG029-01	H09	RT		
QG029-01	H10	RT		
QG029-01	H11	PCR		
QG029-01	H12	PCR		