

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG052-01	A01	HQP015565	NM_020661	AICDA
QG052-01	A02	HQP009139	NM_001643	APOA2
QG052-01	A03	HQP008515	NM_013314	BLNK
QG052-01	A04	HQP016959	NM_001716	BLR1
QG052-01	A05	HQP016622	NM_002983	CCL3
QG052-01	A06	HQP016625	NM_002984	CCL4
QG052-01	A07	HQP002198	NM_001295	CCR1
QG052-01	A08	HQP002201	NM_000648	CCR2
QG052-01	A09	HQP002209	NM_005508	CCR4
QG052-01	A10	HQP002210	NM_000579	CCR5
QG052-01	A11	HQP022190	NM_001767	CD2
QG052-01	A12	HQP022667	NM_001242	CD27
QG052-01	B01	HQP022699	NM_006139	CD28
QG052-01	B02	HQP022212	NM_000732	CD3D
QG052-01	B03	HQP022236	NM_000733	CD3E
QG052-01	B04	HQP022256	NM_000073	CD3G
QG052-01	B05	HQP022316	NM_000616	CD4
QG052-01	B06	HQP022955	NM_001250	CD40
QG052-01	B07	HQP022962	NM_000074	CD40LG
QG052-01	B08	HQP022337	NM_014207	CD5
QG052-01	B09	HQP022399	NM_006137	CD7
QG052-01	B10	HQP022722	NM_005191	CD80
QG052-01	B11	HQP023168	NM_004356	CD81
QG052-01	B12	HQP022746	NM_006889	CD86
QG052-01	C01	HQP022438	NM_004931	CD8B
QG052-01	C02	HQP003159	NM_000758	CSF2
QG052-01	C03	HQP016652	NM_002996	CX3CL1
QG052-01	C04	HQP007900	NM_001504	CXCR3
QG052-01	C05	HQP018803	NM_003467	CXCR4
QG052-01	C06	HQP009651	NM_000043	FAS
QG052-01	C07	HQP009671	NM_000639	FASLG
QG052-01	C08	HQP012676	NM_015892	GALNAC4S-6ST
QG052-01	C09	HQP023167	NM_006037	HDAC4
QG052-01	C10	HQP000024	NM_005474	HDAC5
QG052-01	C11	HQP012862	NM_016596	HDAC7A
QG052-01	C12	HQP008866	NM_019111	HLA-DRA
QG052-01	D01	HQP005977	NM_015259	ICOSLG
QG052-01	D02	HQP009463	NM_002176	IFNB1
QG052-01	D03	HQP009467	NM_000619	IFNG
QG052-01	D04	HQP009469	NM_000416	IFNGR1
QG052-01	D05	HQP009472	NM_005534	IFNGR2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG052-01	D06	HQP009505	NM_001551	IGBP1
QG052-01	D07	HQP009685	NM_000572	IL10
QG052-01	D08	HQP009692	NM_000882	IL12A
QG052-01	D09	HQP009693	NM_002187	IL12B
QG052-01	D10	HQP009694	NM_005535	IL12RB1
QG052-01	D11	HQP009696	NM_001559	IL12RB2
QG052-01	D12	HQP009697	NM_002188	IL13
QG052-01	E01	HQP009718	NM_001562	IL18
QG052-01	E02	HQP009649	NM_000586	IL2
QG052-01	E03	HQP009650	NM_000417	IL2RA
QG052-01	E04	HQP009662	NM_000589	IL4
QG052-01	E05	HQP009664	NM_000418	IL4R
QG052-01	E06	HQP009666	NM_000879	IL5
QG052-01	E07	HQP009678	NM_000584	IL8
QG052-01	E08	HQP009742	NM_002191	INHA
QG052-01	E09	HQP009743	NM_002192	INHBA
QG052-01	E10	HQP009781	NM_002460	IRF4
QG052-01	E11	HQP005972	NM_015254	KIF13B
QG052-01	E12	HQP053978	NM_001300	KLF6
QG052-01	F01	HQP022527	NM_021950	MS4A1
QG052-01	F02	HQP011697	NM_006153	NCK1
QG052-01	F03	HQP011866	NM_000625	NOS2A
QG052-01	F04	HQP015027	NM_000949	PRLR
QG052-01	F05	HQP015956	NM_002855	PVRL1
QG052-01	F06	HQP016251	NM_002922	RGS1
QG052-01	F07	HQP016980	NM_003019	SFTPD
QG052-01	F08	HQP054032	NM_144949	SOCS5
QG052-01	F09	HQP018044	NM_000660	TGFB1
QG052-01	F10	HQP018113	NM_003263	TLR1
QG052-01	F11	HQP018114	NM_003264	TLR2
QG052-01	F12	HQP018116	NM_138554	TLR4
QG052-01	G01	HQP000415	NM_006068	TLR6
QG052-01	G02	HQP013388	NM_017442	TLR9
QG052-01	G03	HQP021496	NM_003807	TNFSF14
QG052-01	G04	HQP013496	NM_019009	TOLLIP
QG052-01	G05	HQP017052	NM_197954	CLEC7A
QG052-01	G06	HQP009026	NM_000194	HPRT1
QG052-01	G07	HQP009653	NM_152872	FAS
QG052-01	G08	HQP023134	NM_014707	HDAC9
QG052-01	G09	HQP020713	NM_001004722	NCK2
QG052-01	G10	HQP022418	NM_001768	CD8A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG052-01	G11	HQP009688	NM_000641	IL11
QG052-01	G12	HQP021570	NM_003855	IL18R1
QG052-01	H01	HGDC		
QG052-01	H02	HGDC		
QG052-01	H03	HQP006940	NM_002046	GAPDH
QG052-01	H04	HQP016381	NM_001101	ACTB
QG052-01	H05	HQP015171	NM_004048	B2M
QG052-01	H06	HQP006171	NM_012423	RPL13A
QG052-01	H07	HQP009026	NM_000194	HPRT1
QG052-01	H08	HQP054253	NR_003286	RN18S1
QG052-01	H09	RT		
QG052-01	H10	RT		
QG052-01	H11	PCR		
QG052-01	H12	PCR		