

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG098-01	A01	HQP022754	NM_152942	TNFRSF8
QG098-01	A02	HQP017180	NM_173217	ST6GAL1
QG098-01	A03	HQP009653	NM_152872	FAS
QG098-01	A04	HQP009026	NM_000194	HPRT1
QG098-01	A05	HQP018504	NM_000552	VWF
QG098-01	A06	HQP053975	NM_001078	VCAM1
QG098-01	A07	HQP018221	NM_003294	TPSAB1
QG098-01	A08	HQP022753	NM_001243	TNFRSF8
QG098-01	A09	HQP018012	NM_000459	TEK
QG098-01	A10	HQP010942	NM_002354	TACSTD1
QG098-01	A11	HQP017179	NM_003032	ST6GAL1
QG098-01	A12	HQP016746	NM_003005	SELP
QG098-01	B01	HQP016544	NM_002964	S100A8
QG098-01	B02	HQP015160	NM_020415	RETN
QG098-01	B03	HQP013015	NM_000442	PECAM1
QG098-01	B04	HQP011932	NM_002526	NT5E
QG098-01	B05	HQP011868	NM_000603	NOS3
QG098-01	B06	HQP022608	NM_153604	MYOCD
QG098-01	B07	HQP011614	NM_002473	MYH9
QG098-01	B08	HQP011615	NM_005964	MYH10
QG098-01	B09	HQP022527	NM_021950	MS4A1
QG098-01	B10	HQP010141	NM_000424	KRT5
QG098-01	B11	HQP010197	NM_000224	KRT18
QG098-01	B12	HQP010112	NM_002262	KLRD1
QG098-01	C01	HQP010106	NM_002258	KLRB1
QG098-01	C02	HQP009794	NM_002203	ITGA2
QG098-01	C03	HQP009793	NM_181501	ITGA1
QG098-01	C04	HQP009650	NM_000417	IL2RA
QG098-01	C05	HQP018800	NM_004633	IL1R2
QG098-01	C06	HQP009694	NM_005535	IL12RB1
QG098-01	C07	HQP009187	NM_000873	ICAM2
QG098-01	C08	HQP008866	NM_019111	HLA-DRA
QG098-01	C09	HQP008849	NM_002116	HLA-A
QG098-01	C10	HQP005251	NM_000566	FCGR1A
QG098-01	C11	HQP005242	NM_002002	FCER2
QG098-01	C12	HQP005238	NM_002001	FCER1A
QG098-01	D01	HQP009651	NM_000043	FAS
QG098-01	D02	HQP004856	NM_000118	ENG
QG098-01	D03	HQP004434	NM_001935	DPP4
QG098-01	D04	HQP003500	NM_005214	CTLA4
QG098-01	D05	HQP003158	NM_005211	CSF1R

QG098-01	D06	HQP002462	NM_000088	COL1A1
QG098-01	D07	HQP022817	NM_004854	CHST10
QG098-01	D08	HQP000279	NM_005816	CD96
QG098-01	D09	HQP022438	NM_004931	CD8B
QG098-01	D10	HQP022746	NM_006889	CD86
QG098-01	D11	HQP022722	NM_005191	CD80
QG098-01	D12	HQP023155	NM_000626	CD79B
QG098-01	E01	HQP023117	NM_001782	CD72
QG098-01	E02	HQP023108	NM_001252	CD70
QG098-01	E03	HQP022399	NM_006137	CD7
QG098-01	E04	HQP023094	NM_001781	CD69
QG098-01	E05	HQP022384	NM_006725	CD6
QG098-01	E06	HQP022337	NM_014207	CD5
QG098-01	E07	HQP022962	NM_000074	CD40LG
QG098-01	E08	HQP022955	NM_001250	CD40
QG098-01	E09	HQP022316	NM_000616	CD4
QG098-01	E10	HQP022256	NM_000073	CD3G
QG098-01	E11	HQP022212	NM_000732	CD3D
QG098-01	E12	HQP022870	NM_001775	CD38
QG098-01	F01	HQP022699	NM_006139	CD28
QG098-01	F02	HQP022295	NM_000734	CD247
QG098-01	F03	HQP013003	NM_016382	CD244
QG098-01	F04	HQP022579	NM_013230	CD24
QG098-01	F05	HQP022561	NM_001771	CD22
QG098-01	F06	HQP008808	NM_021155	CD209
QG098-01	F07	HQP022190	NM_001767	CD2
QG098-01	F08	HQP022129	NM_001766	CD1D
QG098-01	F09	HQP022111	NM_001765	CD1C
QG098-01	F10	HQP022070	NM_001763	CD1A
QG098-01	F11	HQP022504	NM_001770	CD19
QG098-01	F12	HQP022548	NM_004244	CD163
QG098-01	G01	HQP001290	NM_007053	CD160
QG098-01	G02	HQP018326	NM_001736	C5AR1
QG098-01	G03	HQP023157	NM_021602	CD79B
QG098-01	G04	HQP022856	NM_001774	CD37
QG098-01	G05	HQP022418	NM_001768	CD8A
QG098-01	G06	HQP023129	NM_004355	CD74
QG098-01	G07	HQP023144	NM_001783	CD79A
QG098-01	G08	HQP022501	NM_004233	CD83
QG098-01	G09	HQP002466	NM_000089	COL1A2
QG098-01	G10	HQP002890	NM_001877	CR2
QG098-01	G11	HQP009796	NM_002204	ITGA3
QG098-01	G12	HQP010107	NM_002259	KLRC1

QG098-01	H01	HGDC		
QG098-01	H02	HGDC		
QG098-01	H03	HQP006940	NM_002046	GAPDH
QG098-01	H04	HQP016381	NM_001101	ACTB
QG098-01	H05	HQP015171	NM_004048	B2M
QG098-01	H06	HQP006171	NM_012423	RPL13A
QG098-01	H07	HQP009026	NM_000194	HPRT1
QG098-01	H08	HQP054253	NR_003286	RN18S1
QG098-01	H09	RT		
QG098-01	H10	RT		
QG098-01	H11	PCR		
QG098-01	H12	PCR		