

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG033-01	A01	HQP018214	NM_000547	TPO
QG033-01	A02	HQP018068	NM_003246	THBS1
QG033-01	A03	HQP017673	NM_000582	SPP1
QG033-01	A04	HQP003078	NM_001040022	SIRPA
QG033-01	A05	HQP012154	NM_000602	SERPINE1
QG033-01	A06	HQP015059	NM_000313	PROS1
QG033-01	A07	HQP013257	NM_000301	PLG
QG033-01	A08	HQP013204	NM_002658	PLAU
QG033-01	A09	HQP013201	NM_000930	PLAT
QG033-01	A10	HQP023437	NM_005123	NR1H4
QG033-01	A11	HQP011268	NM_004995	MMP14
QG033-01	A12	HQP011077	NM_000242	MBL2
QG033-01	B01	HQP000882	NM_006610	MASP2
QG033-01	B02	HQP010870	NM_002332	LRP1
QG033-01	B03	HQP010136	NM_006121	KRT1
QG033-01	B04	HQP010100	NM_002257	KLK1
QG033-01	B05	HQP009818	NM_000212	ITGB3
QG033-01	B06	HQP009815	NM_000211	ITGB2
QG033-01	B07	HQP009809	NM_000887	ITGAX
QG033-01	B08	HQP009808	NM_002210	ITGAV
QG033-01	B09	HQP009807	NM_000632	ITGAM
QG033-01	B10	HQP009430	NM_021057	IFNA7
QG033-01	B11	HQP009048	NM_000412	HRG
QG033-01	B12	HQP007230	NM_000583	GC
QG033-01	C01	HQP005339	NM_002003	FCN1
QG033-01	C02	HQP002994	NM_000567	CRP
QG033-01	C03	HQP002889	NM_001006658	CR2
QG033-01	C04	HQP002857	NM_001308	CPN1
QG033-01	C05	HQP002037	NM_001831	CLU
QG033-01	C06	HQP013060	NM_002621	CFP
QG033-01	C07	HQP019873	NM_030787	CFHR5
QG033-01	C08	HQP001001	NM_006684	CFHR4
QG033-01	C09	HQP008755	NM_000186	CFH
QG033-01	C10	HQP004198	NM_001928	CFD
QG033-01	C11	HQP016565	NM_001710	CFB
QG033-01	C12	HQP023178	NM_001025160	CD97
QG033-01	D01	HQP005589	NM_012072	CD93
QG033-01	D02	HQP023056	NM_000611	CD59
QG033-01	D03	HQP003948	NM_000574	CD55
QG033-01	D04	HQP011113	NM_002389	CD46
QG033-01	D05	HQP022972	NM_000610	CD44

QG033-01	D06	HQP008808	NM_021155	CD209
QG033-01	D07	HQP017515	NM_001002259	CAPRIN2
QG033-01	D08	HQP018414	NM_001737	C9
QG033-01	D09	HQP018391	NM_000606	C8G
QG033-01	D10	HQP018380	NM_000066	C8B
QG033-01	D11	HQP018362	NM_000562	C8A
QG033-01	D12	HQP018353	NM_000587	C7
QG033-01	E01	HQP018344	NM_000065	C6
QG033-01	E02	HQP018326	NM_001736	C5AR1
QG033-01	E03	HQP018304	NM_001735	C5
QG033-01	E04	HQP018276	NM_000716	C4BPB
QG033-01	E05	HQP018259	NM_000715	C4BPA
QG033-01	E06	HQP018248	NM_007293	C4A
QG033-01	E07	HQP018239	NM_004054	C3AR1
QG033-01	E08	HQP018238	NM_000064	C3
QG033-01	E09	HQP018226	NM_000063	C2
QG033-01	E10	HQP018206	NM_001734	C1S
QG033-01	E11	HQP012584	NM_016546	C1RL
QG033-01	E12	HQP018179	NM_001733	C1R
QG033-01	F01	HQP009207	NM_001008223	C1QL4
QG033-01	F02	HQP010407	NM_001010908	C1QL3
QG033-01	F03	HQP001007	NM_006688	C1QL1
QG033-01	F04	HQP018170	NM_172369	C1QC
QG033-01	F05	HQP018106	NM_001212	C1QBP
QG033-01	F06	HQP018146	NM_015991	C1QA
QG033-01	F07	HQP008346	NM_014059	C13orf15
QG033-01	F08	HQP016610	NM_001711	BGN
QG033-01	F09	HQP009031	NM_001639	APCS
QG033-01	F10	HQP004137	NM_182528	C1QL2
QG033-01	F11	HQP017007	NM_033225	CSMD1
QG033-01	F12	HQP009173	NM_004967	IBSP
QG033-01	G01	HQP015091	NM_001031849	MASP1
QG033-01	G02	HQP008768	NM_002113	CFHR1
QG033-01	G03	HQP008770	NM_005666	CFHR2
QG033-01	G04	HQP001002	NM_021023	CFHR3
QG033-01	G05	HQP009376	NM_000204	CFI
QG033-01	G06	HQP002836	NM_001872	CPB2
QG033-01	G07	HQP002880	NM_000573	CR1
QG033-01	G08	HQP014437	NM_001083116	PRF1
QG033-01	G09	HQP004071	NM_001920	DCN
QG033-01	G10	HQP010280	NM_001014980	FAM132A
QG033-01	G11	HQP005991	NM_002023	FMOD
QG033-01	G12	HQP007592	NM_018485	GPR77

QG033-01	H01	HGDC		
QG033-01	H02	HGDC		
QG033-01	H03	HQP006940	NM_002046	GAPDH
QG033-01	H04	HQP016381	NM_001101	ACTB
QG033-01	H05	HQP015171	NM_004048	B2M
QG033-01	H06	HQP006171	NM_012423	RPL13A
QG033-01	H07	HQP009026	NM_000194	HPRT1
QG033-01	H08	HQP054253	NR_003286	RN18S1
QG033-01	H09	RT		
QG033-01	H10	RT		
QG033-01	H11	PCR		
QG033-01	H12	PCR		