

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG019-01	A01	HQP006023	NM_054034	FN1
QG019-01	A02	HQP005401	NM_033136	FGF1
QG019-01	A03	HQP019768	NM_130445	COL18A1
QG019-01	A04	HQP008804	NM_001010934	HGF
QG019-01	A05	HQP008801	NM_001010931	HGF
QG019-01	A06	HQP009026	NM_000194	HPRT1
QG019-01	A07	HQP018481	NM_003376	VEGFA
QG019-01	A08	HQP018153	NM_000363	TNNI3
QG019-01	A09	HQP018152	NM_003282	TNNI2
QG019-01	A10	HQP018141	NM_000594	TNF
QG019-01	A11	HQP018094	NM_000362	TIMP3
QG019-01	A12	HQP018093	NM_003255	TIMP2
QG019-01	B01	HQP018092	NM_003254	TIMP1
QG019-01	B02	HQP018091	NM_005424	TIE1
QG019-01	B03	HQP018068	NM_003246	THBS1
QG019-01	B04	HQP018044	NM_000660	TGFB1
QG019-01	B05	HQP018043	NM_003236	TGFA
QG019-01	B06	HQP005820	NM_015136	STAB1
QG019-01	B07	HQP001140	NM_006846	SPINK5
QG019-01	B08	HQP013026	NM_002615	SERPINF1
QG019-01	B09	HQP012154	NM_000602	SERPINE1
QG019-01	B10	HQP011620	NM_000488	SERPINC1
QG019-01	B11	HQP021347	NM_001754	RUNX1
QG019-01	B12	HQP016332	NM_002939	RNH1
QG019-01	C01	HQP015024	NM_000948	PRL
QG019-01	C02	HQP013662	NM_002704	PPBP
QG019-01	C03	HQP013257	NM_000301	PLG
QG019-01	C04	HQP019651	NM_025208	PDGFD
QG019-01	C05	HQP012856	NM_002608	PDGFB
QG019-01	C06	HQP010581	NM_000230	LEP
QG019-01	C07	HQP009637	NM_001648	KLK3
QG019-01	C08	HQP011205	NM_003994	KITLG
QG019-01	C09	HQP009678	NM_000584	IL8
QG019-01	C10	HQP009670	NM_000600	IL6
QG019-01	C11	HQP001495	NM_052872	IL17F
QG019-01	C12	HQP009693	NM_002187	IL12B
QG019-01	D01	HQP009692	NM_000882	IL12A
QG019-01	D02	HQP009685	NM_000572	IL10
QG019-01	D03	HQP009467	NM_000619	IFNG
QG019-01	D04	HQP009463	NM_002176	IFNB1
QG019-01	D05	HQP009419	NM_024013	IFNA1

QG019-01	D06	HQP008800	NM_000601	HGF
QG019-01	D07	HQP008462	NM_002091	GRP
QG019-01	D08	HQP006022	NM_002026	FN1
QG019-01	D09	HQP005451	NM_004469	FIGF
QG019-01	D10	HQP023445	NM_005130	FGFBP1
QG019-01	D11	HQP005403	NM_002006	FGF2
QG019-01	D12	HQP005400	NM_000800	FGF1
QG019-01	E01	HQP004978	NM_001432	EREG
QG019-01	E02	HQP004538	NM_001953	ECGF1
QG019-01	E03	HQP011220	NM_002416	CXCL9
QG019-01	E04	HQP016648	NM_002993	CXCL6
QG019-01	E05	HQP016650	NM_002994	CXCL5
QG019-01	E06	HQP008459	NM_002090	CXCL3
QG019-01	E07	HQP008458	NM_002089	CXCL2
QG019-01	E08	HQP022893	NM_004887	CXCL14
QG019-01	E09	HQP000672	NM_006419	CXCL13
QG019-01	E10	HQP016669	NM_000609	CXCL12
QG019-01	E11	HQP016649	NM_005409	CXCL11
QG019-01	E12	HQP009746	NM_001565	CXCL10
QG019-01	F01	HQP003173	NM_000759	CSF3
QG019-01	F02	HQP001309	NM_001275	CHGA
QG019-01	F03	HQP023056	NM_000611	CD59
QG019-01	F04	HQP003948	NM_000574	CD55
QG019-01	F05	HQP016621	NM_002982	CCL2
QG019-01	F06	HQP054027	NM_032965	CCL15
QG019-01	F07	HQP017983	NM_001731	BTG1
QG019-01	F08	HQP017333	NM_001200	BMP2
QG019-01	F09	HQP015745	NM_001702	BAI1
QG019-01	F10	HQP022013	NM_004673	ANGPTL1
QG019-01	F11	HQP008202	NM_001147	ANGPT2
QG019-01	F12	HQP007968	NM_001145	ANG
QG019-01	G01	HQP003781	NM_133265	AMOT
QG019-01	G02	HQP014059	NM_018046	AGGF1
QG019-01	G03	HQP013057	NM_002619	PF4
QG019-01	G04	HQP013089	NM_002632	PGF
QG019-01	G05	HQP008097	NM_001146	ANGPT1
QG019-01	G06	HQP020717	NM_032414	PROK1
QG019-01	G07	HQP002508	NM_000091	COL4A3
QG019-01	G08	HQP005422	NM_004114	FGF13
QG019-01	G09	HQP008330	NM_002087	GRN
QG019-01	G10	HQP053984	NM_002391	MDK
QG019-01	G11	HQP011897	NM_002521	NPPB
QG019-01	G12	HQP010327	NM_004040	RHOB

QG019-01	H01	HGDC		
QG019-01	H02	HGDC		
QG019-01	H03	HQP006940	NM_002046	GAPDH
QG019-01	H04	HQP016381	NM_001101	ACTB
QG019-01	H05	HQP015171	NM_004048	B2M
QG019-01	H06	HQP006171	NM_012423	RPL13A
QG019-01	H07	HQP009026	NM_000194	HPRT1
QG019-01	H08	HQP054253	NR_003286	RN18S1
QG019-01	H09	RT		
QG019-01	H10	RT		
QG019-01	H11	PCR		
QG019-01	H12	PCR		