

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG002-01	A01	HQP008801	NM_001010931	HGF
QG002-01	A02	HQP019768	NM_130445	COL18A1
QG002-01	A03	HQP008804	NM_001010934	HGF
QG002-01	A04	HQP021591	NM_201264	NRP2
QG002-01	A05	HQP009026	NM_000194	HPRT1
QG002-01	A06	HQP018483	NM_005429	VEGFC
QG002-01	A07	HQP018481	NM_003376	VEGFA
QG002-01	A08	HQP018144	NM_006291	TNFAIP2
QG002-01	A09	HQP018141	NM_000594	TNF
QG002-01	A10	HQP018094	NM_000362	TIMP3
QG002-01	A11	HQP018093	NM_003255	TIMP2
QG002-01	A12	HQP018092	NM_003254	TIMP1
QG002-01	B01	HQP018069	NM_003247	THBS2
QG002-01	B02	HQP018068	NM_003246	THBS1
QG002-01	B03	HQP018051	NM_004612	TGFBR1
QG002-01	B04	HQP018047	NM_003238	TGFB2
QG002-01	B05	HQP018044	NM_000660	TGFB1
QG002-01	B06	HQP018043	NM_003236	TGFA
QG002-01	B07	HQP018012	NM_000459	TEK
QG002-01	B08	HQP005820	NM_015136	STAB1
QG002-01	B09	HQP021658	NM_021972	SPHK1
QG002-01	B10	HQP013026	NM_002615	SERPINF1
QG002-01	B11	HQP015596	NM_000962	PTGS1
QG002-01	B12	HQP016353	NM_021935	PROK2
QG002-01	C01	HQP015418	NM_020405	PLXDC1
QG002-01	C02	HQP013257	NM_000301	PLG
QG002-01	C03	HQP013204	NM_002658	PLAU
QG002-01	C04	HQP013015	NM_000442	PECAM1
QG002-01	C05	HQP012847	NM_002607	PDGFA
QG002-01	C06	HQP021589	NM_003872	NRP2
QG002-01	C07	HQP021597	NM_003873	NRP1
QG002-01	C08	HQP011263	NM_004994	MMP9
QG002-01	C09	HQP010581	NM_000230	LEP
QG002-01	C10	HQP010070	NM_002253	KDR
QG002-01	C11	HQP004470	NM_000214	JAG1
QG002-01	C12	HQP009818	NM_000212	ITGB3
QG002-01	D01	HQP009808	NM_002210	ITGAV
QG002-01	D02	HQP009678	NM_000584	IL8
QG002-01	D03	HQP009670	NM_000600	IL6
QG002-01	D04	HQP009641	NM_000576	IL1B
QG002-01	D05	HQP009518	NM_000618	IGF1

QG002-01	D06	HQP009467	NM_000619	IFNG
QG002-01	D07	HQP009463	NM_002176	IFNB1
QG002-01	D08	HQP009419	NM_024013	IFNA1
QG002-01	D09	HQP009282	NM_002167	ID3
QG002-01	D10	HQP009266	NM_002165	ID1
QG002-01	D11	HQP000974	NM_006665	HPSE
QG002-01	D12	HQP008831	NM_001530	HIF1A
QG002-01	E01	HQP008800	NM_000601	HGF
QG002-01	E02	HQP022788	NM_021973	HAND2
QG002-01	E03	HQP005879	NM_002019	FLT1
QG002-01	E04	HQP005451	NM_004469	FIGF
QG002-01	E05	HQP005434	NM_000142	FGFR3
QG002-01	E06	HQP005403	NM_002006	FGF2
QG002-01	E07	HQP005400	NM_000800	FGF1
QG002-01	E08	HQP004978	NM_001432	EREG
QG002-01	E09	HQP004945	NM_004444	EPHB4
QG002-01	E10	HQP004856	NM_000118	ENG
QG002-01	E11	HQP004599	NM_001963	EGF
QG002-01	E12	HQP004597	NM_004093	EFNB2
QG002-01	F01	HQP004591	NM_004952	EFNA3
QG002-01	F02	HQP004589	NM_182685	EFNA1
QG002-01	F03	HQP004554	NM_001400	EDG1
QG002-01	F04	HQP004538	NM_001953	ECGF1
QG002-01	F05	HQP011220	NM_002416	CXCL9
QG002-01	F06	HQP016648	NM_002993	CXCL6
QG002-01	F07	HQP016650	NM_002994	CXCL5
QG002-01	F08	HQP008459	NM_002090	CXCL3
QG002-01	F09	HQP009746	NM_001565	CXCL10
QG002-01	F10	HQP008456	NM_001511	CXCL1
QG002-01	F11	HQP000052	NM_001795	CDH5
QG002-01	F12	HQP016621	NM_002982	CCL2
QG002-01	G01	HQP016629	NM_002986	CCL11
QG002-01	G02	HQP015745	NM_001702	BAI1
QG002-01	G03	HQP008414	NM_001150	ANPEP
QG002-01	G04	HQP012432	NM_001039667	ANGPTL4
QG002-01	G05	HQP007675	NM_014495	ANGPTL3
QG002-01	G06	HQP008202	NM_001147	ANGPT2
QG002-01	G07	HQP054002	NM_005163	AKT1
QG002-01	G08	HQP008097	NM_001146	ANGPT1
QG002-01	G09	HQP019767	NM_030582	COL18A1
QG002-01	G10	HQP002508	NM_000091	COL4A3
QG002-01	G11	HQP013057	NM_002619	PF4
QG002-01	G12	HQP001209	NM_007015	LECT1

QG002-01	H01	HGDC		
QG002-01	H02	HGDC		
QG002-01	H03	HQP006940	NM_002046	GAPDH
QG002-01	H04	HQP016381	NM_001101	ACTB
QG002-01	H05	HQP015171	NM_004048	B2M
QG002-01	H06	HQP006171	NM_012423	RPL13A
QG002-01	H07	HQP009026	NM_000194	HPRT1
QG002-01	H08	HQP054253	NR_003286	RN18S1
QG002-01	H09	RT		
QG002-01	H10	RT		
QG002-01	H11	PCR		
QG002-01	H12	PCR		