

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG046-01	A01	HQP016478	NM_001015051	RUNX2
QG046-01	A02	HQP006023	NM_054034	FN1
QG046-01	A03	HQP005430	NM_023107	FGFR1
QG046-01	A04	HQP009026	NM_000194	HPRT1
QG046-01	A05	HQP018481	NM_003376	VEGFA
QG046-01	A06	HQP018474	NM_000376	VDR
QG046-01	A07	HQP053975	NM_001078	VCAM1
QG046-01	A08	HQP018328	NM_000474	TWIST1
QG046-01	A09	HQP018320	NM_020127	TUFT1
QG046-01	A10	HQP018141	NM_000594	TNF
QG046-01	A11	HQP018051	NM_004612	TGFBR1
QG046-01	A12	HQP018048	NM_003239	TGFB3
QG046-01	B01	HQP018047	NM_003238	TGFB2
QG046-01	B02	HQP018044	NM_000660	TGFB1
QG046-01	B03	HQP054013	NM_012143	TFIP11
QG046-01	B04	HQP017635	NM_000346	SOX9
QG046-01	B05	HQP010961	NM_005359	SMAD4
QG046-01	B06	HQP010960	NM_005902	SMAD3
QG046-01	B07	HQP054007	NM_005901	SMAD2
QG046-01	B08	HQP054006	NM_005900	SMAD1
QG046-01	B09	HQP021470	NM_001235	SERPINH1
QG046-01	B10	HQP022833	NM_005505	SCARB1
QG046-01	B11	HQP021334	NM_004348	RUNX2
QG046-01	B12	HQP013108	NM_000444	PHEX
QG046-01	C01	HQP012847	NM_002607	PDGFA
QG046-01	C02	HQP011807	NM_003998	NFKB1
QG046-01	C03	HQP011263	NM_004994	MMP9
QG046-01	C04	HQP011262	NM_002424	MMP8
QG046-01	C05	HQP011264	NM_002425	MMP10
QG046-01	C06	HQP022913	NM_004897	MINPP1
QG046-01	C07	HQP009810	NM_002211	ITGB1
QG046-01	C08	HQP009807	NM_000632	ITGAM
QG046-01	C09	HQP009794	NM_002203	ITGA2
QG046-01	C10	HQP009793	NM_181501	ITGA1
QG046-01	C11	HQP009529	NM_000612	IGF2
QG046-01	C12	HQP009523	NM_000875	IGF1R
QG046-01	D01	HQP009518	NM_000618	IGF1
QG046-01	D02	HQP009184	NM_000201	ICAM1
QG046-01	D03	HQP006022	NM_002026	FN1
QG046-01	D04	HQP005879	NM_002019	FLT1
QG046-01	D05	HQP005437	NM_000141	FGFR2

QG046-01	D06	HQP005403	NM_002006	FGF2
QG046-01	D07	HQP005400	NM_000800	FGF1
QG046-01	D08	HQP004605	NM_005228	EGFR
QG046-01	D09	HQP004599	NM_001963	EGF
QG046-01	D10	HQP004478	NM_014208	DSPP
QG046-01	D11	HQP004375	NM_004407	DMP1
QG046-01	D12	HQP003626	NM_000396	CTSK
QG046-01	E01	HQP003173	NM_000759	CSF3
QG046-01	E02	HQP003159	NM_000758	CSF2
QG046-01	E03	HQP002670	NM_000095	COMP
QG046-01	E04	HQP002541	NM_000093	COL5A1
QG046-01	E05	HQP002478	NM_000090	COL3A1
QG046-01	E06	HQP002473	NM_001844	COL2A1
QG046-01	E07	HQP002462	NM_000088	COL1A1
QG046-01	E08	HQP002644	NM_001855	COL15A1
QG046-01	E09	HQP018426	NM_021110	COL14A1
QG046-01	E10	HQP002612	NM_004370	COL12A1
QG046-01	E11	HQP002602	NM_080629	COL11A1
QG046-01	E12	HQP000131	NM_001797	CDH11
QG046-01	F01	HQP022821	NM_000072	CD36
QG046-01	F02	HQP019452	NM_001742	CALCR
QG046-01	F03	HQP017433	NM_021073	BMP5
QG046-01	F04	HQP053910	NM_130851	BMP4
QG046-01	F05	HQP017357	NM_001201	BMP3
QG046-01	F06	HQP017333	NM_001200	BMP2
QG046-01	F07	HQP017282	NM_006129	BMP1
QG046-01	F08	HQP016610	NM_001711	BGN
QG046-01	F09	HQP016599	NM_199173	BGLAP
QG046-01	F10	HQP008829	NM_001154	ANXA5
QG046-01	F11	HQP006840	NM_016519	AMBN
QG046-01	F12	HQP006440	NM_000478	ALPL
QG046-01	G01	HQP011526	NM_002448	MSX1
QG046-01	G02	HQP002613	NM_080645	COL12A1
QG046-01	G03	HQP018054	NM_003242	TGFBR2
QG046-01	G04	HQP018482	NM_003377	VEGFB
QG046-01	G05	HQP004606	NM_201282	EGFR
QG046-01	G06	HQP002596	NM_000493	COL10A1
QG046-01	G07	HQP002466	NM_000089	COL1A2
QG046-01	G08	HQP002508	NM_000091	COL4A3
QG046-01	G09	HQP000152	NM_031889	ENAM
QG046-01	G10	HQP005404	NM_005247	FGF3
QG046-01	G11	HQP005427	NM_015850	FGFR1
QG046-01	G12	HQP007326	NM_004962	GDF10

QG046-01	H01	HGDC		
QG046-01	H02	HGDC		
QG046-01	H03	HQP006940	NM_002046	GAPDH
QG046-01	H04	HQP016381	NM_001101	ACTB
QG046-01	H05	HQP015171	NM_004048	B2M
QG046-01	H06	HQP006171	NM_012423	RPL13A
QG046-01	H07	HQP009026	NM_000194	HPRT1
QG046-01	H08	HQP054253	NR_003286	RN18S1
QG046-01	H09	RT		
QG046-01	H10	RT		
QG046-01	H11	PCR		
QG046-01	H12	PCR		