

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG088-01	A01	HQP020596	NM_001042544	LTBP4
QG088-01	A02	HQP008818	NM_004495	NRG1
QG088-01	A03	HQP003150	NM_172210	CSF1
QG088-01	A04	HQP013043	NM_017424	CECR1
QG088-01	A05	HQP009026	NM_000194	HPRT1
QG088-01	A06	HQP018483	NM_005429	VEGFC
QG088-01	A07	HQP018481	NM_003376	VEGFA
QG088-01	A08	HQP018074	NM_000460	THPO
QG088-01	A09	HQP018044	NM_000660	TGFB1
QG088-01	A10	HQP017997	NM_003212	TDGF1
QG088-01	A11	HQP017673	NM_000582	SPP1
QG088-01	A12	HQP014905	NM_016205	PDGFC
QG088-01	B01	HQP054039	NM_182981	OSGIN1
QG088-01	B02	HQP011933	NM_002527	NTF3
QG088-01	B03	HQP022888	NM_013982	NRG2
QG088-01	B04	HQP008820	NM_013957	NRG1
QG088-01	B05	HQP011861	NM_018055	NODAL
QG088-01	B06	HQP011827	NM_002506	NGFB
QG088-01	B07	HQP010607	NM_002309	LIF
QG088-01	B08	HQP018049	NM_003240	LEFTY2
QG088-01	B09	HQP000772	NM_020997	LEFTY1
QG088-01	B10	HQP009847	NM_002226	JAG2
QG088-01	B11	HQP004470	NM_000214	JAG1
QG088-01	B12	HQP009744	NM_002193	INHBB
QG088-01	C01	HQP009743	NM_002192	INHBA
QG088-01	C02	HQP009742	NM_002191	INHA
QG088-01	C03	HQP009662	NM_000589	IL4
QG088-01	C04	HQP009649	NM_000586	IL2

QG088-01	C05	HQP009641	NM_000576	IL1B
QG088-01	C06	HQP009640	NM_000575	IL1A
QG088-01	C07	HQP009718	NM_001562	IL18
QG088-01	C08	HQP009693	NM_002187	IL12B
QG088-01	C09	HQP009685	NM_000572	IL10
QG088-01	C10	HQP009529	NM_000612	IGF2
QG088-01	C11	HQP009518	NM_000618	IGF1
QG088-01	C12	HQP004493	NM_001945	HBEGF
QG088-01	D01	HQP007814	NM_000175	GPI
QG088-01	D02	HQP007346	NM_000514	GDNF
QG088-01	D03	HQP007323	NM_005259	GDF8
QG088-01	D04	HQP000275	NM_005811	GDF11
QG088-01	D05	HQP005451	NM_004469	FIGF
QG088-01	D06	HQP005416	NM_002010	FGF9
QG088-01	D07	HQP005411	NM_002009	FGF7
QG088-01	D08	HQP005408	NM_004464	FGF5
QG088-01	D09	HQP005403	NM_002006	FGF2
QG088-01	D10	HQP021583	NM_003867	FGF17
QG088-01	D11	HQP005424	NM_004115	FGF14
QG088-01	D12	HQP005419	NM_004112	FGF11
QG088-01	E01	HQP005400	NM_000800	FGF1
QG088-01	E02	HQP004978	NM_001432	EREG
QG088-01	E03	HQP004538	NM_001953	ECGF1
QG088-01	E04	HQP005612	NM_012242	DKK1
QG088-01	E05	HQP008456	NM_001511	CXCL1
QG088-01	E06	HQP003173	NM_000759	CSF3
QG088-01	E07	HQP003159	NM_000758	CSF2
QG088-01	E08	HQP003149	NM_000757	CSF1
QG088-01	E09	HQP001978	NM_001828	CLC
QG088-01	E10	HQP013044	NM_177405	CECR1

QG088-01	E11	HQP017477	NM_001720	BMP8B
QG088-01	E12	HQP017467	NM_001719	BMP7
QG088-01	F01	HQP017433	NM_021073	BMP5
QG088-01	F02	HQP053910	NM_130851	BMP4
QG088-01	F03	HQP017357	NM_001201	BMP3
QG088-01	F04	HQP017333	NM_001200	BMP2
QG088-01	F05	HQP007656	NM_014482	BMP10
QG088-01	F06	HQP017282	NM_006129	BMP1
QG088-01	F07	HQP005409	NM_033143	FGF5
QG088-01	F08	HQP011700	NM_000266	NDP
QG088-01	F09	HQP015033	NM_004158	PSPN
QG088-01	F10	HQP018154	NM_003283	TNNT1
QG088-01	F11	HQP007394	NM_000479	AMH
QG088-01	F12	HQP013010	NM_016442	ARTS-1
QG088-01	G01	HQP016545	NM_001709	BDNF
QG088-01	G02	HQP013089	NM_002632	PGF
QG088-01	G03	HQP005422	NM_004114	FGF13
QG088-01	G04	HQP023416	NM_005117	FGF19
QG088-01	G05	HQP007429	NM_020637	FGF22
QG088-01	G06	HQP019743	NM_020638	FGF23
QG088-01	G07	HQP007326	NM_004962	GDF10
QG088-01	G08	HQP009688	NM_000641	IL11
QG088-01	G09	HQP009660	NM_000588	IL3
QG088-01	G10	HQP020598	NM_003573	LTBP4
QG088-01	G11	HQP053984	NM_002391	MDK
QG088-01	G12	HQP000852	NM_001010848	NRG3
QG088-01	H01	HGDC		
QG088-01	H02	HGDC		
QG088-01	H03	HQP006940	NM_002046	GAPDH
QG088-01	H04	HQP016381	NM_001101	ACTB

QG088-01	H05	HQP015171	NM_004048	B2M
QG088-01	H06	HQP006171	NM_012423	RPL13A
QG088-01	H07	HQP009026	NM_000194	HPRT1
QG088-01	H08	HQP054253	NR_003286	RN18S1
QG088-01	H09	RT		
QG088-01	H10	RT		
QG088-01	H11	PCR		
QG088-01	H12	PCR		