

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG060-01	A01	HQP021608	NM_003881	WISP2
QG060-01	A02	HQP021610	NM_003882	WISP1
QG060-01	A03	HQP018483	NM_005429	VEGFC
QG060-01	A04	HQP018475	NM_001025366	VEGFA
QG060-01	A05	HQP003299	NM_004385	VCAN
QG060-01	A06	HQP018328	NM_000474	TWIST1
QG060-01	A07	HQP018104	NM_005077	TLE1
QG060-01	A08	HQP018090	NM_003253	TIAM1
QG060-01	A09	HQP018047	NM_003238	TGFB2
QG060-01	A10	HQP017962	NM_030756	TCF7L2
QG060-01	A11	HQP020164	NM_031283	TCF7L1
QG060-01	A12	HQP017958	NM_003202	TCF7
QG060-01	B01	HQP017951	NM_003199	TCF4
QG060-01	B02	HQP017957	NM_003200	TCF3
QG060-01	B03	HQP017955	NM_000458	TCF2
QG060-01	B04	HQP017954	NM_000545	TCF1
QG060-01	B05	HQP019196	NM_024665	TBL1XR1
QG060-01	B06	HQP017927	NM_005647	TBL1X
QG060-01	B07	HQP017857	NM_003181	T
QG060-01	B08	HQP017635	NM_000346	SOX9
QG060-01	B09	HQP015461	NM_020436	SALL4
QG060-01	B10	HQP016478	NM_001015051	RUNX2
QG060-01	B11	HQP015997	NM_021205	RHOA
QG060-01	B12	HQP022362	NM_004219	PTTG1
QG060-01	C01	HQP015598	NM_000963	PTGS2
QG060-01	C02	HQP013627	NM_006238	PPARD
QG060-01	C03	HQP000764	NM_006475	POSTN
QG060-01	C04	HQP013205	NM_001005376	PLAUR
QG060-01	C05	HQP011915	NM_001037132	NRCAM
QG060-01	C06	HQP011866	NM_000625	NOS2A
QG060-01	C07	HQP019390	NM_024865	NANOG
QG060-01	C08	HQP007208	NM_012333	MYCBP
QG060-01	C09	HQP011597	NM_002467	MYC
QG060-01	C10	HQP011263	NM_004994	MMP9
QG060-01	C11	HQP011258	NM_002423	MMP7
QG060-01	C12	HQP011257	NM_002422	MMP3
QG060-01	D01	HQP011256	NM_004530	MMP2
QG060-01	D02	HQP011181	NM_000245	MET
QG060-01	D03	HQP010877	NM_002336	LRP6
QG060-01	D04	HQP012480	NM_016269	LEF1
QG060-01	D05	HQP010390	NM_000425	L1CAM

QG060-01	D06	HQP009853	NM_002228	JUN
QG060-01	D07	HQP004470	NM_000214	JAG1
QG060-01	D08	HQP009791	NM_002202	ISL1
QG060-01	D09	HQP009670	NM_000600	IL6
QG060-01	D10	HQP009529	NM_000612	IGF2
QG060-01	D11	HQP009518	NM_000618	IGF1
QG060-01	D12	HQP009273	NM_002166	ID2
QG060-01	E01	HQP007411	NM_000165	GJA1
QG060-01	E02	HQP007233	NM_002054	GCG
QG060-01	E03	HQP020140	NM_031866	FZD8
QG060-01	E04	HQP020136	NM_003507	FZD7
QG060-01	E05	HQP020808	NM_003593	FOXN1
QG060-01	E06	HQP019708	NM_005438	FOSL1
QG060-01	E07	HQP005416	NM_002010	FGF9
QG060-01	E08	HQP005405	NM_002007	FGF4
QG060-01	E09	HQP007198	NM_019851	FGF20
QG060-01	E10	HQP004942	NM_004443	EPHB3
QG060-01	E11	HQP004939	NM_004442	EPHB2
QG060-01	E12	HQP004605	NM_005228	EGFR
QG060-01	F01	HQP004596	NM_004429	EFNB1
QG060-01	F02	HQP004557	NM_001955	EDN1
QG060-01	F03	HQP004544	NM_001005609	EDA
QG060-01	F04	HQP008111	NM_005618	DLL1
QG060-01	F05	HQP005612	NM_012242	DKK1
QG060-01	F06	HQP003539	NM_001904	CTNNB1
QG060-01	F07	HQP022026	NM_021101	CLDN1
QG060-01	F08	HQP022972	NM_000610	CD44
QG060-01	F09	HQP021752	NM_003939	BTRC
QG060-01	F10	HQP017396	NM_001202	BMP4
QG060-01	F11	HQP009099	NM_001012270	BIRC5
QG060-01	F12	HQP016599	NM_199173	BGLAP
QG060-01	G01	HQP020116	NM_004655	AXIN2
QG060-01	G02	HQP004183	NM_001130	AES
QG060-01	G03	HQP011762	NM_005172	ATOH1
QG060-01	G04	HQP000534	NM_001804	CDX1
QG060-01	G05	HQP000567	NM_005193	CDX4
QG060-01	G06	HQP000564	NM_006350	FST
QG060-01	G07	HQP011773	NM_006161	NEUROG1
QG060-01	G08	HQP023466	NM_004360	CDH1
QG060-01	G09	HQP009792	NM_005545	ISLR
QG060-01	G10	HQP021321	NM_003701	TNFSF11
QG060-01	G11	HQP008470	NM_000177	GSN
QG060-01	G12	HQP018993	NM_024336	IRX3

QG060-01	H01	HGDC		
QG060-01	H02	HGDC		
QG060-01	H03	HQP006940	NM_002046	GAPDH
QG060-01	H04	HQP016381	NM_001101	ACTB
QG060-01	H05	HQP015171	NM_004048	B2M
QG060-01	H06	HQP006171	NM_012423	RPL13A
QG060-01	H07	HQP009026	NM_000194	HPRT1
QG060-01	H08	HQP054253	NR_003286	RN18S1
QG060-01	H09	RT		
QG060-01	H10	RT		
QG060-01	H11	PCR		
QG060-01	H12	PCR		