

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG022-01	A01	HQP003236	NM_001025105	CSNK1A1
QG022-01	A02	HQP021752	NM_003939	BTRC
QG022-01	A03	HQP009026	NM_000194	HPRT1
QG022-01	A04	HQP018596	NM_007129	ZIC2
QG022-01	A05	HQP018543	NM_003396	WNT9B
QG022-01	A06	HQP018542	NM_003395	WNT9A
QG022-01	A07	HQP018534	NM_058238	WNT7B
QG022-01	A08	HQP018533	NM_004625	WNT7A
QG022-01	A09	HQP054026	NM_032642	WNT5B
QG022-01	A10	HQP018531	NM_003392	WNT5A
QG022-01	A11	HQP013446	NM_030761	WNT4
QG022-01	A12	HQP021771	NM_033131	WNT3A
QG022-01	B01	HQP018530	NM_030753	WNT3
QG022-01	B02	HQP018540	NM_004185	WNT2B
QG022-01	B03	HQP018529	NM_003391	WNT2
QG022-01	B04	HQP018539	NM_004626	WNT11
QG022-01	B05	HQP018538	NM_003394	WNT10B
QG022-01	B06	HQP019670	NM_025216	WNT10A
QG022-01	B07	HQP018528	NM_005430	WNT1
QG022-01	B08	HQP001390	NM_007191	WIF1
QG022-01	B09	HQP012958	NM_016169	SUFU
QG022-01	B10	HQP007549	NM_015690	STK36
QG022-01	B11	HQP017563	NM_005631	SMO
QG022-01	B12	HQP017150	NM_003031	SIAH1
QG022-01	C01	HQP017098	NM_000193	SHH
QG022-01	C02	HQP016865	NM_003012	SFRP1
QG022-01	C03	HQP054041	NM_183227	RAB23
QG022-01	C04	HQP009877	NM_001034842	PTCHD3
QG022-01	C05	HQP002968	NM_173495	PTCHD1
QG022-01	C06	HQP021384	NM_003738	PTCH2
QG022-01	C07	HQP015530	NM_000264	PTCH1
QG022-01	C08	HQP014987	NM_005044	PRKX
QG022-01	C09	HQP014594	NM_002732	PRKACG
QG022-01	C10	HQP014575	NM_002730	PRKACA
QG022-01	C11	HQP012085	NM_021728	OTX2
QG022-01	C12	HQP023201	NM_014751	MTSS1
QG022-01	D01	HQP014848	NM_002745	MAPK1
QG022-01	D02	HQP010871	NM_004525	LRP2
QG022-01	D03	HQP003354	NM_001002914	KCTD11
QG022-01	D04	HQP009632	NM_002181	IHH
QG022-01	D05	HQP012391	NM_016004	IFT52

QG022-01	D06	HQP016957	NM_022475	HHIP
QG022-01	D07	HQP014640	NM_018194	HHAT
QG022-01	D08	HQP054075	NM_002093	GSK3B
QG022-01	D09	HQP007703	NM_000168	GLI3
QG022-01	D10	HQP007702	NM_005270	GLI2
QG022-01	D11	HQP007701	NM_005269	GLI1
QG022-01	D12	HQP006344	NM_012181	FKBP8
QG022-01	E01	HQP005434	NM_000142	FGFR3
QG022-01	E02	HQP005416	NM_002010	FGF9
QG022-01	E03	HQP005960	NM_012300	FBXW11
QG022-01	E04	HQP003539	NM_001904	CTNNB1
QG022-01	E05	HQP003257	NM_001319	CSNK1G2
QG022-01	E06	HQP013348	NM_022048	CSNK1G1
QG022-01	E07	HQP003249	NM_001894	CSNK1E
QG022-01	E08	HQP003240	NM_001893	CSNK1D
QG022-01	E09	HQP002148	NM_145203	CSNK1A1L
QG022-01	E10	HQP003237	NM_001892	CSNK1A1
QG022-01	E11	HQP012543	NM_016441	CRIM1
QG022-01	E12	HQP012262	NM_016952	CDON
QG022-01	F01	HQP019423	NM_024899	CEP76
QG022-01	F02	HQP008588	NM_013326	C18orf8
QG022-01	F03	HQP021753	NM_033637	BTRC
QG022-01	F04	HQP017477	NM_001720	BMP8B
QG022-01	F05	HQP009619	NM_181809	BMP8A
QG022-01	F06	HQP017467	NM_001719	BMP7
QG022-01	F07	HQP017433	NM_021073	BMP5
QG022-01	F08	HQP053910	NM_130851	BMP4
QG022-01	F09	HQP017333	NM_001200	BMP2
QG022-01	F10	HQP007313	NM_013372	GREM1
QG022-01	F11	HQP018535	NM_058244	WNT8A
QG022-01	F12	HQP018536	NM_003393	WNT8B
QG022-01	G01	HQP018595	NM_003412	ZIC1
QG022-01	G02	HQP005962	NM_033645	FBXW11
QG022-01	G03	HQP011480	NM_001013732	RP3-402H5.2
QG022-01	G04	HQP012238	NM_021044	DHH
QG022-01	G05	HQP012699	NM_057168	WNT16
QG022-01	G06	HQP005704	NM_004473	FOXE1
QG022-01	G07	HQP007115	NM_002048	GAS1
QG022-01	G08	HQP008559	NM_013389	NPC1L1
QG022-01	G09	HQP018532	NM_006522	WNT6
QG022-01	G10	HQP014584	NM_182948	PRKACB
QG022-01	G11	HQP015021	NM_002760	PRKY
QG022-01	G12	HQP054019	NM_020780	PTCHD2

QG022-01	H01	HGDC		
QG022-01	H02	HGDC		
QG022-01	H03	HQP006940	NM_002046	GAPDH
QG022-01	H04	HQP016381	NM_001101	ACTB
QG022-01	H05	HQP015171	NM_004048	B2M
QG022-01	H06	HQP006171	NM_012423	RPL13A
QG022-01	H07	HQP009026	NM_000194	HPRT1
QG022-01	H08	HQP054253	NR_003286	RN18S1
QG022-01	H09	RT		
QG022-01	H10	RT		
QG022-01	H11	PCR		
QG022-01	H12	PCR		