

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG043-01	A01	HQP013633	NM_005037	PPARG
QG043-01	A02	HQP012980	NM_031268	PDPK1
QG043-01	A03	HQP009026	NM_000194	HPRT1
QG043-01	A04	HQP022974	NM_001001390	CD44
QG043-01	A05	HQP018596	NM_007129	ZIC2
QG043-01	A06	HQP018539	NM_004626	WNT11
QG043-01	A07	HQP021610	NM_003882	WISP1
QG043-01	A08	HQP018104	NM_005077	TLE1
QG043-01	A09	HQP018003	NM_021961	TEAD1
QG043-01	A10	HQP017243	NM_003035	STIL
QG043-01	A11	HQP010936	NM_002351	SH2D1A
QG043-01	A12	HQP021347	NM_001754	RUNX1
QG043-01	B01	HQP014773	NM_172341	PSENNEN
QG043-01	B02	HQP013634	NM_015869	PPARG
QG043-01	B03	HQP006158	NM_172236	POFUT1
QG043-01	B04	HQP012212	NM_016734	PAX5
QG043-01	B05	HQP011968	NM_006186	NR4A2
QG043-01	B06	HQP053985	NM_002502	NFKB2
QG043-01	B07	HQP011601	NM_005376	MYCL1
QG043-01	B08	HQP011258	NM_002423	MMP7
QG043-01	B09	HQP014926	NM_145185	MAP2K7
QG043-01	B10	HQP010809	NM_000427	LOR
QG043-01	B11	HQP010136	NM_006121	KRT1
QG043-01	B12	HQP009650	NM_000417	IL2RA
QG043-01	C01	HQP007583	NM_014443	IL17B
QG043-01	C02	HQP009467	NM_000619	IFNG
QG043-01	C03	HQP008990	NM_024015	HOXB4
QG043-01	C04	HQP054075	NM_002093	GSK3B
QG043-01	C05	HQP007701	NM_005269	GLI1
QG043-01	C06	HQP007223	NM_004120	GBP2
QG043-01	C07	HQP004969	NM_004448	ERBB2
QG043-01	C08	HQP053993	NM_003903	CDC16
QG043-01	C09	HQP022972	NM_000610	CD44
QG043-01	C10	HQP021819	NM_001238	CCNE1
QG043-01	C11	HQP021430	NM_005188	CBL
QG043-01	C12	HQP018044	NM_000660	TGFB1
QG043-01	D01	HQP017824	NM_003170	SUPT6H
QG043-01	D02	HQP012958	NM_016169	SUFU
QG043-01	D03	HQP017775	NM_003153	STAT6
QG043-01	D04	HQP017767	NM_003150	STAT3
QG043-01	D05	HQP005609	NM_012245	SNW1

QG043-01	D06	HQP017563	NM_005631	SMO
QG043-01	D07	HQP010959	NM_001003652	SMAD2
QG043-01	D08	HQP010958	NM_001003688	SMAD1
QG043-01	D09	HQP017098	NM_000193	SHH
QG043-01	D10	HQP016743	NM_005065	SEL1L
QG043-01	D11	HQP009574	NM_005349	RBPJ
QG043-01	D12	HQP015123	NM_000021	PSEN1
QG043-01	E01	HQP006157	NM_015352	POFUT1
QG043-01	E02	HQP012979	NM_002613	PDPK1
QG043-01	E03	HQP021621	NM_003884	PCAF
QG043-01	E04	HQP012219	NM_000280	PAX6
QG043-01	E05	HQP022406	NM_004756	NUMBL
QG043-01	E06	HQP021395	NM_001005743	NUMB
QG043-01	E07	HQP011876	NM_000435	NOTCH3
QG043-01	E08	HQP011875	NM_024408	NOTCH2
QG043-01	E09	HQP011873	NM_017617	NOTCH1
QG043-01	E10	HQP011808	NM_001077493	NFKB2
QG043-01	E11	HQP011807	NM_003998	NFKB1
QG043-01	E12	HQP006054	NM_015331	NCSTN
QG043-01	F01	HQP011597	NM_002467	MYC
QG043-01	F02	HQP011191	NM_002405	MFNG
QG043-01	F03	HQP014450	NM_018717	MAML3
QG043-01	F04	HQP009847	NM_002226	JAG2
QG043-01	F05	HQP004470	NM_000214	JAG1
QG043-01	F06	HQP009674	NM_002184	IL6ST
QG043-01	F07	HQP014725	NM_005144	HR
QG043-01	F08	HQP006121	NM_001040708	HEY1
QG043-01	F09	HQP009057	NM_005524	HES1
QG043-01	F10	HQP008746	NM_001527	HDAC2
QG043-01	F11	HQP008745	NM_004964	HDAC1
QG043-01	F12	HQP008469	NM_002093	GSK3B
QG043-01	G01	HQP007253	NM_021078	GCN5L2
QG043-01	G02	HQP007165	NM_001002295	GATA3
QG043-01	G03	HQP020136	NM_003507	FZD7
QG043-01	G04	HQP020135	NM_003506	FZD6
QG043-01	G05	HQP020134	NM_012193	FZD4
QG043-01	G06	HQP019238	NM_017412	FZD3
QG043-01	G07	HQP006492	NM_001466	FZD2
QG043-01	G08	HQP020133	NM_003505	FZD1
QG043-01	G09	HQP019708	NM_005438	FOSL1
QG043-01	G10	HQP006188	NM_005252	FOS
QG043-01	G11	HQP005451	NM_004469	FIGF
QG043-01	G12	HQP004968	NM_001005862	ERBB2

QG043-01	H01	HGDC		
QG043-01	H02	HGDC		
QG043-01	H03	HQP006940	NM_002046	GAPDH
QG043-01	H04	HQP016381	NM_001101	ACTB
QG043-01	H05	HQP015171	NM_004048	B2M
QG043-01	H06	HQP006171	NM_012423	RPL13A
QG043-01	H07	HQP009026	NM_000194	HPRT1
QG043-01	H08	HQP054253	NR_003286	RN18S1
QG043-01	H09	RT		
QG043-01	H10	RT		
QG043-01	H11	PCR		
QG043-01	H12	PCR		