

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG009-01	A01	HQP012353	NM_001018073	PCK2
QG009-01	A02	HQP009026	NM_000194	HPRT1
QG009-01	A03	HQP018490	NM_003381	VIP
QG009-01	A04	HQP018141	NM_000594	TNF
QG009-01	A05	HQP018068	NM_003246	THBS1
QG009-01	A06	HQP018064	NM_000360	TH
QG009-01	A07	HQP018048	NM_003239	TGFB3
QG009-01	A08	HQP017867	NM_001058	TACR1
QG009-01	A09	HQP017767	NM_003150	STAT3
QG009-01	A10	HQP017744	NM_001050	SSTR2
QG009-01	A11	HQP053963	NM_001048	SST
QG009-01	A12	HQP017706	NM_003131	SRF
QG009-01	B01	HQP017616	NM_000636	SOD2
QG009-01	B02	HQP017478	NM_003053	SLC18A1
QG009-01	B03	HQP017005	NM_005627	SGK
QG009-01	B04	HQP018805	NM_003469	SCG2
QG009-01	B05	HQP016542	NM_014624	S100A6
QG009-01	B06	HQP016551	NM_005621	S100A12
QG009-01	B07	HQP016131	NM_000321	RB1
QG009-01	B08	HQP015598	NM_000963	PTGS2
QG009-01	B09	HQP015024	NM_000948	PRL
QG009-01	B10	HQP014650	NM_002734	PRKAR1A
QG009-01	B11	HQP014115	NM_002715	PPP2CA
QG009-01	B12	HQP006298	NM_014330	PPP1R15A
QG009-01	C01	HQP013528	NM_006235	POU2AF1
QG009-01	C02	HQP013268	NM_002667	PLN
QG009-01	C03	HQP013201	NM_000930	PLAT
QG009-01	C04	HQP054038	NM_182649	PCNA
QG009-01	C05	HQP012354	NM_004563	PCK2
QG009-01	C06	HQP011968	NM_006186	NR4A2
QG009-01	C07	HQP011874	NM_000905	NPY
QG009-01	C08	HQP011866	NM_000625	NOS2A
QG009-01	C09	HQP011774	NM_000267	NF1
QG009-01	C10	HQP011219	NM_002415	MIF
QG009-01	C11	HQP010969	NM_005360	MAF
QG009-01	C12	HQP010571	NM_005566	LDHA
QG009-01	D01	HQP009873	NM_002234	KCNA5
QG009-01	D02	HQP009854	NM_002229	JUNB
QG009-01	D03	HQP009743	NM_002192	INHBA
QG009-01	D04	HQP009670	NM_000600	IL6
QG009-01	D05	HQP009649	NM_000586	IL2

QG009-01	D06	HQP009083	NM_005347	HSPA5
QG009-01	D07	HQP009081	NM_002154	HSPA4
QG009-01	D08	HQP008843	NM_000189	HK2
QG009-01	D09	HQP007408	NM_000164	GIPR
QG009-01	D10	HQP007353	NM_005261	GEM
QG009-01	D11	HQP007233	NM_002054	GCG
QG009-01	D12	HQP006188	NM_005252	FOS
QG009-01	E01	HQP004864	NM_001975	ENO2
QG009-01	E02	HQP004613	NM_000399	EGR2
QG009-01	E03	HQP004612	NM_001964	EGR1
QG009-01	E04	HQP004498	NM_004417	DUSP1
QG009-01	E05	HQP004127	NM_004083	DDIT3
QG009-01	E06	HQP003479	NM_001330	CTF1
QG009-01	E07	HQP002950	NM_183011	CREM
QG009-01	E08	HQP002907	NM_004379	CREB1
QG009-01	E09	HQP002384	NM_001299	CNN1
QG009-01	E10	HQP001309	NM_001275	CHGA
QG009-01	E11	HQP000956	NM_000735	CGA
QG009-01	E12	HQP000261	NM_004935	CDK5
QG009-01	F01	HQP016204	NM_053056	CCND1
QG009-01	F02	HQP021692	NM_003914	CCNA1
QG009-01	F03	HQP019841	NM_004343	CALR
QG009-01	F04	HQP019580	NM_006888	CALM1
QG009-01	F05	HQP019024	NM_004929	CALB1
QG009-01	F06	HQP017713	NM_007294	BRCA1
QG009-01	F07	HQP016211	NM_000633	BCL2
QG009-01	F08	HQP053980	NM_001674	ATF3
QG009-01	F09	HQP009941	NM_001657	AREG
QG009-01	F10	HQP007213	NM_001634	AMD1
QG009-01	F11	HQP004658	NM_001621	AHR
QG009-01	F12	HQP003754	NM_000684	ADRB1
QG009-01	G01	HQP005410	NM_020996	FGF6
QG009-01	G02	HQP006205	NM_006732	FOSB
QG009-01	G03	HQP013033	NM_006211	PENK
QG009-01	G04	HQP013296	NM_021127	PMAIP1
QG009-01	G05	HQP013519	NM_000306	POU1F1
QG009-01	G06	HQP016545	NM_001709	BDNF
QG009-01	G07	HQP019049	NM_001740	CALB2
QG009-01	G08	HQP000382	NM_004936	CDKN2B
QG009-01	G09	HQP009855	NM_005354	JUND
QG009-01	G10	HQP011688	NM_000615	NCAM1
QG009-01	G11	HQP013046	NM_002616	PER1
QG009-01	G12	HQP018467	NM_003373	VCL

QG009-01	H01	HGDC		
QG009-01	H02	HGDC		
QG009-01	H03	HQP006940	NM_002046	GAPDH
QG009-01	H04	HQP016381	NM_001101	ACTB
QG009-01	H05	HQP015171	NM_004048	B2M
QG009-01	H06	HQP006171	NM_012423	RPL13A
QG009-01	H07	HQP009026	NM_000194	HPRT1
QG009-01	H08	HQP054253	NR_003286	RN18S1
QG009-01	H09	RT		
QG009-01	H10	RT		
QG009-01	H11	PCR		
QG009-01	H12	PCR		