

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG082-01	A01	HQP018120	NM_004615	TSPAN7
QG082-01	A02	HQP022531	NM_004240	TRIP10
QG082-01	A03	HQP018175	NM_000546	TP53
QG082-01	A04	HQP018093	NM_003255	TIMP2
QG082-01	A05	HQP018092	NM_003254	TIMP1
QG082-01	A06	HQP018063	NM_003245	TGM3
QG082-01	A07	HQP018061	NM_004613	TGM2
QG082-01	A08	HQP018060	NM_000359	TGM1
QG082-01	A09	HQP017928	NM_003194	TBP
QG082-01	A10	HQP017640	NM_138473	SP1
QG082-01	A11	HQP017318	NM_004171	SLC1A2
QG082-01	A12	HQP017005	NM_005627	SGK
QG082-01	B01	HQP023308	NM_012432	SETDB1
QG082-01	B02	HQP008383	NM_014159	SETD2
QG082-01	B03	HQP016600	NM_174934	SCN4B
QG082-01	B04	HQP021524	NM_003821	RIPK2
QG082-01	B05	HQP016125	NM_002890	RASA1
QG082-01	B06	HQP000241	NM_005789	PSME3
QG082-01	B07	HQP006702	NM_001031698	PRPF40B
QG082-01	B08	HQP015032	NM_000311	PRNP
QG082-01	B09	HQP020918	NM_032639	PLEKHA8
QG082-01	B10	HQP000968	NM_006661	PDE10A
QG082-01	B11	HQP011874	NM_000905	NPY
QG082-01	B12	HQP022978	NM_006311	NCOR1
QG082-01	C01	HQP011547	NM_005957	MTHFR
QG082-01	C02	HQP011008	NM_000898	MAOB
QG082-01	C03	HQP010581	NM_000230	LEP
QG082-01	C04	HQP021287	NM_003679	KMO
QG082-01	C05	HQP010007	NM_004981	KCNJ4
QG082-01	C06	HQP010003	NM_000891	KCNJ2
QG082-01	C07	HQP021838	NM_001024660	KALRN
QG082-01	C08	HQP014035	NM_018010	IFT57
QG082-01	C09	HQP008834	NM_005339	HIP2
QG082-01	C10	HQP008833	NM_005338	HIP1
QG082-01	C11	HQP008744	NM_002111	HD
QG082-01	C12	HQP008371	NM_000833	GRIN2A
QG082-01	D01	HQP008367	NM_000832	GRIN1
QG082-01	D02	HQP008291	NM_002086	GRB2
QG082-01	D03	HQP008324	NM_014030	GIT1
QG082-01	D04	HQP012996	NM_016362	GHRL
QG082-01	D05	HQP007346	NM_000514	GDNF

QG082-01	D06	HQP006940	NM_002046	GAPDH
QG082-01	D07	HQP006683	NM_000817	GAD1
QG082-01	D08	HQP006426	NM_004958	FRAP1
QG082-01	D09	HQP006188	NM_005252	FOS
QG082-01	D10	HQP005403	NM_002006	FGF2
QG082-01	D11	HQP004897	NM_001429	EP300
QG082-01	D12	HQP004612	NM_001964	EGR1
QG082-01	E01	HQP009074	NM_001039550	DNAJB2
QG082-01	E02	HQP004336	NM_000108	DLD
QG082-01	E03	HQP004096	NM_004082	DCTN1
QG082-01	E04	HQP002920	NM_001079846	CREBBP
QG082-01	E05	HQP002128	NM_004859	CLTC
QG082-01	E06	HQP000615	NM_004364	CEBPA
QG082-01	E07	HQP000261	NM_004935	CDK5
QG082-01	E08	HQP021518	NM_000071	CBS
QG082-01	E09	HQP018966	NM_001080124	CASP8
QG082-01	E10	HQP020427	NM_001226	CASP6
QG082-01	E11	HQP020297	NM_004346	CASP3
QG082-01	E12	HQP020207	NM_001223	CASP1
QG082-01	F01	HQP006132	NM_014296	CAPN7
QG082-01	F02	HQP018286	NM_004055	CAPN5
QG082-01	F03	HQP001299	NM_021251	CAPN10
QG082-01	F04	HQP020065	NM_005186	CAPN1
QG082-01	F05	HQP013039	NM_017422	CALML5
QG082-01	F06	HQP019834	NM_005185	CALML3
QG082-01	F07	HQP019811	NM_005184	CALM3
QG082-01	F08	HQP019707	NM_001743	CALM2
QG082-01	F09	HQP019580	NM_006888	CALM1
QG082-01	F10	HQP019024	NM_004929	CALB1
QG082-01	F11	HQP015964	NM_004324	BAX
QG082-01	F12	HQP016578	NM_000332	ATXN1
QG082-01	G01	HQP002912	NM_001880	ATF2
QG082-01	G02	HQP004991	NM_001014431	AKT1
QG082-01	G03	HQP002833	NM_000675	ADORA2A
QG082-01	G04	HQP006134	NM_014298	QPRT
QG082-01	G05	HQP001045	NM_001040006	TCERG1
QG082-01	G06	HQP002110	NM_001076677	CLTA
QG082-01	G07	HQP004203	NM_001004285	DFFB
QG082-01	G08	HQP008703	NM_002105	H2AFX
QG082-01	G09	HQP021847	NM_001079870	HAP1
QG082-01	G10	HQP015551	NM_020655	JPH3
QG082-01	G11	HQP053965	NM_001105579	LOC646658
QG082-01	G12	HQP020511	NM_032192	PPP1R1B

QG082-01	H01	HGDC		
QG082-01	H02	HGDC		
QG082-01	H03	HQP006940	NM_002046	GAPDH
QG082-01	H04	HQP016381	NM_001101	ACTB
QG082-01	H05	HQP015171	NM_004048	B2M
QG082-01	H06	HQP006171	NM_012423	RPL13A
QG082-01	H07	HQP009026	NM_000194	HPRT1
QG082-01	H08	HQP054253	NR_003286	RN18S1
QG082-01	H09	RT		
QG082-01	H10	RT		
QG082-01	H11	PCR		
QG082-01	H12	PCR		