

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG084-01	A01	HQP018381	NM_003347	UBE2L3
QG084-01	A02	HQP018357	NM_003334	UBE1
QG084-01	A03	HQP018175	NM_000546	TP53
QG084-01	A04	HQP018064	NM_000360	TH
QG084-01	A05	HQP018061	NM_004613	TGM2
QG084-01	A06	HQP017740	NM_006280	SSR4
QG084-01	A07	HQP017615	NM_000454	SOD1
QG084-01	A08	HQP023002	NM_005460	SNCAIP
QG084-01	A09	HQP017582	NM_000345	SNCA
QG084-01	A10	HQP017318	NM_004171	SLC1A2
QG084-01	A11	HQP016125	NM_002890	RASA1
QG084-01	A12	HQP014773	NM_172341	PSENN
QG084-01	B01	HQP015123	NM_000021	PSEN1
QG084-01	B02	HQP015032	NM_000311	PRNP
QG084-01	B03	HQP014309	NM_000944	PPP3CA
QG084-01	B04	HQP020918	NM_032639	PLEKHA8
QG084-01	B05	HQP001549	NM_007262	PARK7
QG084-01	B06	HQP012193	NM_004562	PARK2
QG084-01	B07	HQP011800	NM_006164	NFE2L2
QG084-01	B08	HQP011751	NM_005382	NEFM
QG084-01	B09	HQP011761	NM_006158	NEFL
QG084-01	B10	HQP011759	NM_021076	NEFH
QG084-01	B11	HQP006054	NM_015331	NCSTN
QG084-01	B12	HQP022978	NM_006311	NCOR1
QG084-01	C01	HQP011254	NM_000902	MME
QG084-01	C02	HQP022342	NM_001033057	MAGI1
QG084-01	C03	HQP010870	NM_002332	LRP1
QG084-01	C04	HQP010847	NM_000237	LPL
QG084-01	C05	HQP010527	NM_002291	LAMB1
QG084-01	C06	HQP008000	NM_005559	LAMA1
QG084-01	C07	HQP009863	NM_005548	KARS
QG084-01	C08	HQP020315	NM_031483	ITCH
QG084-01	C09	HQP009764	NM_000208	INSR
QG084-01	C10	HQP009749	NM_000207	INS
QG084-01	C11	HQP014035	NM_018010	IFT57
QG084-01	C12	HQP009098	NM_002156	HSPD1
QG084-01	D01	HQP009083	NM_005347	HSPA5
QG084-01	D02	HQP008834	NM_005339	HIP2
QG084-01	D03	HQP008833	NM_005338	HIP1
QG084-01	D04	HQP000022	NM_006044	HDAC6
QG084-01	D05	HQP008744	NM_002111	HD

QG084-01	D06	HQP008469	NM_002093	GSK3B
QG084-01	D07	HQP008291	NM_002086	GRB2
QG084-01	D08	HQP008279	NM_000581	GPX1
QG084-01	D09	HQP008221	NM_005302	GPR37
QG084-01	D10	HQP007360	NM_002055	GFAP
QG084-01	D11	HQP006940	NM_002046	GAPDH
QG084-01	D12	HQP004096	NM_004082	DCTN1
QG084-01	E01	HQP002920	NM_001079846	CREBBP
QG084-01	E02	HQP002128	NM_004859	CLTC
QG084-01	E03	HQP021518	NM_000071	CBS
QG084-01	E04	HQP018966	NM_001080124	CASP8
QG084-01	E05	HQP020427	NM_001226	CASP6
QG084-01	E06	HQP020297	NM_004346	CASP3
QG084-01	E07	HQP020207	NM_001223	CASP1
QG084-01	E08	HQP013039	NM_017422	CALML5
QG084-01	E09	HQP019811	NM_005184	CALM3
QG084-01	E10	HQP019707	NM_001743	CALM2
QG084-01	E11	HQP019580	NM_006888	CALM1
QG084-01	E12	HQP016237	NM_001191	BCL2L1
QG084-01	F01	HQP016211	NM_000633	BCL2
QG084-01	F02	HQP015964	NM_004324	BAX
QG084-01	F03	HQP000549	NM_006340	BAIAP2
QG084-01	F04	HQP015538	NM_004322	BAD
QG084-01	F05	HQP006760	NM_012105	BACE2
QG084-01	F06	HQP006277	NM_012104	BACE1
QG084-01	F07	HQP004456	NM_001007026	ATN1
QG084-01	F08	HQP021666	NM_001018159	APPBP1
QG084-01	F09	HQP009578	NM_000484	APP
QG084-01	F10	HQP009556	NM_000041	APOE
QG084-01	F11	HQP012404	NM_001077628	APH1A
QG084-01	F12	HQP009007	NM_001164	APBB1
QG084-01	G01	HQP015797	NM_020919	ALS2
QG084-01	G02	HQP008678	NM_000014	A2M
QG084-01	G03	HQP013394	NM_002688	SEPT5
QG084-01	G04	HQP011748	NM_001042681	RERE
QG084-01	G05	HQP016545	NM_001709	BDNF
QG084-01	G06	HQP020946	NM_001752	CAT
QG084-01	G07	HQP010543	NM_001012321	RPSA
QG084-01	G08	HQP018375	NM_003343	UBE2G2
QG084-01	G09	HQP021847	NM_001079870	HAP1
QG084-01	G10	HQP009349	NM_004969	IDE
QG084-01	G11	HQP010532	NM_002293	LAMC1
QG084-01	G12	HQP011022	NM_005910	MAPT

QG084-01	H01	HGDC		
QG084-01	H02	HGDC		
QG084-01	H03	HQP006940	NM_002046	GAPDH
QG084-01	H04	HQP016381	NM_001101	ACTB
QG084-01	H05	HQP015171	NM_004048	B2M
QG084-01	H06	HQP006171	NM_012423	RPL13A
QG084-01	H07	HQP009026	NM_000194	HPRT1
QG084-01	H08	HQP054253	NR_003286	RN18S1
QG084-01	H09	RT		
QG084-01	H10	RT		
QG084-01	H11	PCR		
QG084-01	H12	PCR		