

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG083-01	A01	HQP014423	NM_018646	TRPV6
QG083-01	A02	HQP018256	NM_001001188	TRPM2
QG083-01	A03	HQP022528	NM_004238	TRIP12
QG083-01	A04	HQP007081	NM_015650	TRAF3IP1
QG083-01	A05	HQP018232	NM_005658	TRAF1
QG083-01	A06	HQP002121	NM_173353	TPH2
QG083-01	A07	HQP018195	NM_004179	TPH1
QG083-01	A08	HQP018141	NM_000594	TNF
QG083-01	A09	HQP018104	NM_005077	TLE1
QG083-01	A10	HQP018077	NM_003250	THRA
QG083-01	A11	HQP018056	NM_003244	TGIF1
QG083-01	A12	HQP018044	NM_000660	TGFB1
QG083-01	B01	HQP018026	NM_003221	TFAP2B
QG083-01	B02	HQP018023	NM_001032280	TFAP2A
QG083-01	B03	HQP018020	NM_001063	TF
QG083-01	B04	HQP017958	NM_003202	TCF7
QG083-01	B05	HQP017951	NM_003199	TCF4
QG083-01	B06	HQP020054	NM_003490	SYN3
QG083-01	B07	HQP017841	NM_005638	SYBL1
QG083-01	B08	HQP017810	NM_001055	SULT1A1
QG083-01	B09	HQP017762	NM_000349	STAR
QG083-01	B10	HQP019846	NM_006011	ST8SIA2
QG083-01	B11	HQP053963	NM_001048	SST
QG083-01	B12	HQP015850	NM_020971	SPTBN4
QG083-01	C01	HQP017688	NM_003127	SPTAN1
QG083-01	C02	HQP017676	NM_003124	SPR
QG083-01	C03	HQP017636	NM_006941	SOX10
QG083-01	C04	HQP017615	NM_000454	SOD1
QG083-01	C05	HQP022567	NM_004782	SNAP29
QG083-01	C06	HQP017556	NM_003079	SMARCE1
QG083-01	C07	HQP053962	NM_001045	SLC6A4
QG083-01	C08	HQP017397	NM_001043	SLC6A2
QG083-01	C09	HQP017388	NM_006933	SLC5A3
QG083-01	C10	HQP017478	NM_003053	SLC18A1
QG083-01	C11	HQP017471	NM_005073	SLC15A1
QG083-01	C12	HQP015830	NM_001029880	SFMBT2
QG083-01	D01	HQP018805	NM_003469	SCG2
QG083-01	D02	HQP016572	NM_002970	SAT1
QG083-01	D03	HQP016548	NM_002965	S100A9
QG083-01	D04	HQP016478	NM_001015051	RUNX2
QG083-01	D05	HQP015432	NM_007008	RTN4

QG083-01	D06	HQP016307	NM_007212	RNF2
QG083-01	D07	HQP021492	NM_003804	RIPK1
QG083-01	D08	HQP016259	NM_005613	RGS4
QG083-01	D09	HQP015094	NM_005045	RELN
QG083-01	D10	HQP016213	NM_021975	RELA
QG083-01	D11	HQP016107	NM_002885	RAP1GAP
QG083-01	D12	HQP000062	NM_005493	RANBP9
QG083-01	E01	HQP018819	NM_004637	RAB7A
QG083-01	E02	HQP015948	NM_005859	PURA
QG083-01	E03	HQP015598	NM_000963	PTGS2
QG083-01	E04	HQP015444	NM_002812	PSMD8
QG083-01	E05	HQP015369	NM_002806	PSMC6
QG083-01	E06	HQP006319	NM_005813	PRKD3
QG083-01	E07	HQP014829	NM_001033581	PRKCZ
QG083-01	E08	HQP014804	NM_006257	PRKCQ
QG083-01	E09	HQP014771	NM_002740	PRKCI
QG083-01	E10	HQP014751	NM_002739	PRKCG
QG083-01	E11	HQP014731	NM_006254	PRKCD
QG083-01	E12	HQP014718	NM_002738	PRKCB1
QG083-01	F01	HQP014706	NM_002737	PRKCA
QG083-01	F02	HQP014575	NM_002730	PRKACA
QG083-01	F03	HQP014431	NM_002726	PREP
QG083-01	F04	HQP014262	NM_002719	PPP2R5C
QG083-01	F05	HQP014202	NM_020416	PPP2R2C
QG083-01	F06	HQP000535	NM_006112	PPIE
QG083-01	F07	HQP021090	NM_015053	PPFIA4
QG083-01	F08	HQP013428	NM_002693	POLG
QG083-01	F09	HQP015379	NM_020353	PLSCR4
QG083-01	F10	HQP013272	NM_000533	PLP1
QG083-01	F11	HQP012381	NM_015993	PLLP
QG083-01	F12	HQP013238	NM_002660	PLCG1
QG083-01	G01	HQP021328	NM_003706	PLA2G4C
QG083-01	G02	HQP013195	NM_024420	PLA2G4A
QG083-01	G03	HQP012134	NM_015715	PLA2G3
QG083-01	G04	HQP013194	NM_000300	PLA2G2A
QG083-01	G05	HQP013193	NM_000928	PLA2G1B
QG083-01	G06	HQP006336	NM_012399	PITPNB
QG083-01	G07	HQP013171	NM_005028	PIP5K2A
QG083-01	G08	HQP013161	NM_002651	PIK4CB
QG083-01	G09	HQP013159	NM_002650	PIK4CA
QG083-01	G10	HQP013147	NM_002647	PIK3C3
QG083-01	G11	HQP013093	NM_000291	PGK1
QG083-01	G12	HQP013080	NM_002628	PFN2

QG083-01	H01	HGDC		
QG083-01	H02	HGDC		
QG083-01	H03	HQP006940	NM_002046	GAPDH
QG083-01	H04	HQP016381	NM_001101	ACTB
QG083-01	H05	HQP015171	NM_004048	B2M
QG083-01	H06	HQP006171	NM_012423	RPL13A
QG083-01	H07	HQP009026	NM_000194	HPRT1
QG083-01	H08	HQP054253	NR_003286	RN18S1
QG083-01	H09	RT		
QG083-01	H10	RT		
QG083-01	H11	PCR		
QG083-01	H12	PCR		
QG083-02	A01	HQP021640	NM_016831	PER3
QG083-02	A02	HQP012750	NM_006203	PDE4D
QG083-02	A03	HQP012474	NM_006031	PCNT
QG083-02	A04	HQP018681	NM_007144	PCGF2
QG083-02	A05	HQP012212	NM_016734	PAX5
QG083-02	A06	HQP012197	NM_002583	PAWR
QG083-02	A07	HQP003120	NM_001618	PARP1
QG083-02	A08	HQP012140	NM_000430	PAFAH1B1
QG083-02	A09	HQP012100	NM_002562	P2RX7
QG083-02	A10	HQP012085	NM_021728	OTX2
QG083-02	A11	HQP011938	NM_001007792	NTRK1
QG083-02	A12	HQP011933	NM_002527	NTF3
QG083-02	B01	HQP008818	NM_004495	NRG1
QG083-02	B02	HQP008401	NM_000176	NR3C1
QG083-02	B03	HQP022925	NM_021724	NR1D1
QG083-02	B04	HQP011874	NM_000905	NPY
QG083-02	B05	HQP011885	NM_002518	NPAS2
QG083-02	B06	HQP011868	NM_000603	NOS3
QG083-02	B07	HQP023119	NM_014697	NOS1AP
QG083-02	B08	HQP011811	NM_001001716	NFKBIB
QG083-02	B09	HQP011808	NM_001077493	NFKB2
QG083-02	B10	HQP011807	NM_003998	NFKB1
QG083-02	B11	HQP011735	NM_021074	NDUFV2
QG083-02	B12	HQP011730	NM_007103	NDUFV1
QG083-02	C01	HQP011734	NM_002496	NDUFS8
QG083-02	C02	HQP009875	NM_024407	NDUFS7
QG083-02	C03	HQP019903	NM_001025579	NDEL1
QG083-02	C04	HQP021529	NM_003826	NAPG
QG083-02	C05	HQP011597	NM_002467	MYC
QG083-02	C06	HQP011555	NM_002454	MTRR
QG083-02	C07	HQP011554	NM_000254	MTR

QG083-02	C08	HQP011547	NM_005957	MTHFR
QG083-02	C09	HQP011281	NM_182935	MOBP
QG083-02	C10	HQP005868	NM_015166	MLC1
QG083-02	C11	HQP000555	NM_006343	MERTK
QG083-02	C12	HQP011146	NM_002396	ME2
QG083-02	D01	HQP011098	NM_005913	MC5R
QG083-02	D02	HQP011085	NM_001025081	MBP
QG083-02	D03	HQP014854	NM_001040056	MAPK3
QG083-02	D04	HQP014848	NM_002745	MAPK1
QG083-02	D05	HQP011010	NM_002373	MAP1A
QG083-02	D06	HQP011007	NM_000240	MAOA
QG083-02	D07	HQP010995	NM_002371	MAL
QG083-02	D08	HQP010907	NM_000595	LTA
QG083-02	D09	HQP021587	NM_004664	LIN7A
QG083-02	D10	HQP012480	NM_016269	LEF1
QG083-02	D11	HQP006061	NM_015340	LARS2
QG083-02	D12	HQP010049	NM_004518	KCNQ2
QG083-02	E01	HQP010041	NM_002249	KCNN3
QG083-02	E02	HQP010024	NM_002245	KCNK1
QG083-02	E03	HQP000373	NM_005886	KATNB1
QG083-02	E04	HQP020437	NM_001014380	KATNAL1
QG083-02	E05	HQP021838	NM_001024660	KALRN
QG083-02	E06	HQP009853	NM_002228	JUN
QG083-02	E07	HQP017035	NM_001001132	ITSN1
QG083-02	E08	HQP009842	NM_002222	ITPR1
QG083-02	E09	HQP012769	NM_016368	ISYNA1
QG083-02	E10	HQP009747	NM_002194	INPP1
QG083-02	E11	HQP009727	NM_014214	IMPA2
QG083-02	E12	HQP009726	NM_005536	IMPA1
QG083-02	F01	HQP009683	NM_002186	IL9R
QG083-02	F02	HQP009670	NM_000600	IL6
QG083-02	F03	HQP009658	NM_000878	IL2RB
QG083-02	F04	HQP009645	NM_000577	IL1RN
QG083-02	F05	HQP009641	NM_000576	IL1B
QG083-02	F06	HQP009718	NM_001562	IL18
QG083-02	F07	HQP009541	NM_000597	IGFBP2
QG083-02	F08	HQP009518	NM_000618	IGF1
QG083-02	F09	HQP009123	NM_000869	HTR3A
QG083-02	F10	HQP009122	NM_000868	HTR2C
QG083-02	F11	HQP009120	NM_000621	HTR2A
QG083-02	F12	HQP009117	NM_000864	HTR1D
QG083-02	G01	HQP009116	NM_000863	HTR1B
QG083-02	G02	HQP009083	NM_005347	HSPA5

QG083-02	G03	HQP018231	NM_003299	HSP90B1
QG083-02	G04	HQP008866	NM_019111	HLA-DRA
QG083-02	G05	HQP021900	NM_003959	HIP1R
QG083-02	G06	HQP008745	NM_004964	HDAC1
QG083-02	G07	HQP008690	NM_004131	GZMB
QG083-02	G08	HQP008469	NM_002093	GSK3B
QG083-02	G09	HQP008449	NM_000841	GRM4
QG083-02	G10	HQP008448	NM_000840	GRM3
QG083-02	G11	HQP008380	NM_000836	GRIN2D
QG083-02	G12	HQP008371	NM_000833	GRIN2A
QG083-02	H01	HGDC		
QG083-02	H02	HGDC		
QG083-02	H03	HQP006940	NM_002046	GAPDH
QG083-02	H04	HQP016381	NM_001101	ACTB
QG083-02	H05	HQP015171	NM_004048	B2M
QG083-02	H06	HQP006171	NM_012423	RPL13A
QG083-02	H07	HQP009026	NM_000194	HPRT1
QG083-02	H08	HQP054253	NR_003286	RN18S1
QG083-02	H09	RT		
QG083-02	H10	RT		
QG083-02	H11	PCR		
QG083-02	H12	PCR		
QG083-03	A01	HQP008367	NM_000832	GRIN1
QG083-03	A02	HQP008363	NM_014619	GRIK4
QG083-03	A03	HQP008302	NM_000827	GRIA1
QG083-03	A04	HQP008281	NM_002083	GPX2
QG083-03	A05	HQP008279	NM_000581	GPX1
QG083-03	A06	HQP005436	NM_004466	GPC5
QG083-03	A07	HQP007788	NM_004486	GOLGA2
QG083-03	A08	HQP013589	NM_053004	GNB1L
QG083-03	A09	HQP007755	NM_000516	GNAS
QG083-03	A10	HQP007360	NM_002055	GFAP
QG083-03	A11	HQP007235	NM_000161	GCH1
QG083-03	A12	HQP006750	NM_000153	GALC
QG083-03	B01	HQP006683	NM_000817	GAD1
QG083-03	B02	HQP006651	NM_000816	GABRG2
QG083-03	B03	HQP006613	NM_000810	GABRA5
QG083-03	B04	HQP006599	NM_000806	GABRA1
QG083-03	B05	HQP006515	NM_000402	G6PD
QG083-03	B06	HQP019238	NM_017412	FZD3
QG083-03	B07	HQP019708	NM_005438	FOSL1
QG083-03	B08	HQP006188	NM_005252	FOS
QG083-03	B09	HQP005563	NM_004117	FKBP5

QG083-03	B10	HQP023016	NM_005103	FEZ1
QG083-03	B11	HQP005357	NM_004462	FDFT1
QG083-03	B12	HQP005141	NM_005245	FAT
QG083-03	C01	HQP009671	NM_000639	FASLG
QG083-03	C02	HQP009651	NM_000043	FAS
QG083-03	C03	HQP005050	NM_004456	EZH2
QG083-03	C04	HQP005001	NM_001040275	ESR2
QG083-03	C05	HQP004998	NM_000125	ESR1
QG083-03	C06	HQP004970	NM_001005915	ERBB3
QG083-03	C07	HQP004918	NM_001978	EPB49
QG083-03	C08	HQP004897	NM_001429	EP300
QG083-03	C09	HQP012960	NM_001040092	ENPP2
QG083-03	C10	HQP004842	NM_004098	EMX2
QG083-03	C11	HQP021417	NM_003756	EIF3S3
QG083-03	C12	HQP021680	NM_003908	EIF2S2
QG083-03	D01	HQP004643	NM_004094	EIF2S1
QG083-03	D02	HQP004650	NM_001414	EIF2B1
QG083-03	D03	HQP004605	NM_005228	EGFR
QG083-03	D04	HQP021480	NM_003797	EED
QG083-03	D05	HQP004555	NM_001401	EDG2
QG083-03	D06	HQP004504	NM_001946	DUSP6
QG083-03	D07	HQP020444	NM_032122	DTNBP1
QG083-03	D08	HQP004451	NM_000798	DRD5
QG083-03	D09	HQP004450	NM_000797	DRD4
QG083-03	D10	HQP004446	NM_000795	DRD2
QG083-03	D11	HQP004445	NM_000794	DRD1
QG083-03	D12	HQP004441	NM_001386	DPYSL2
QG083-03	E01	HQP006013	NM_015296	DOCK9
QG083-03	E02	HQP009103	NM_006145	DNAJB1
QG083-03	E03	HQP004342	NM_001365	DLG4
QG083-03	E04	HQP007578	NM_001012957	DISC1
QG083-03	E05	HQP001244	NM_022105	DIDO1
QG083-03	E06	HQP003959	NM_152910	DGKH
QG083-03	E07	HQP004032	NM_001352	DBP
QG083-03	E08	HQP003970	NM_001917	DAO
QG083-03	E09	HQP008458	NM_002089	CXCL2
QG083-03	E10	HQP016652	NM_002996	CX3CL1
QG083-03	E11	HQP005986	NM_015267	CUTL2
QG083-03	E12	HQP003464	NM_001012614	CTBP1
QG083-03	F01	HQP006044	NM_015321	CRTC1
QG083-03	F02	HQP002967	NM_001882	CRHBP
QG083-03	F03	HQP002920	NM_001079846	CREBBP
QG083-03	F04	HQP002907	NM_004379	CREB1

QG083-03	F05	HQP002873	NM_000098	CPT2
QG083-03	F06	HQP002788	NM_004374	COX6C
QG083-03	F07	HQP022636	NM_004255	COX5A
QG083-03	F08	HQP002671	NM_000754	COMT
QG083-03	F09	HQP002400	NM_033133	CNP
QG083-03	F10	HQP022927	NM_004898	CLOCK
QG083-03	F11	HQP012083	NM_005602	CLDN11
QG083-03	F12	HQP001829	NM_001824	CKM
QG083-03	G01	HQP001708	NM_001278	CHUK
QG083-03	G02	HQP001644	NM_000746	CHRNA7
QG083-03	G03	HQP001614	NM_000079	CHRNA1
QG083-03	G04	HQP001535	NM_000739	CHRM2
QG083-03	G05	HQP019569	NM_025114	CEP290
QG083-03	G06	HQP021757	NM_001760	CCND3
QG083-03	G07	HQP021754	NM_001759	CCND2
QG083-03	G08	HQP016204	NM_053056	CCND1
QG083-03	G09	HQP053971	NM_021006	CCL3L1
QG083-03	G10	HQP016630	NM_005408	CCL13
QG083-03	G11	HQP021647	NM_000730	CCKAR
QG083-03	G12	HQP020920	NM_005189	CBX2
QG083-03	H01	HGDC		
QG083-03	H02	HGDC		
QG083-03	H03	HQP006940	NM_002046	GAPDH
QG083-03	H04	HQP016381	NM_001101	ACTB
QG083-03	H05	HQP015171	NM_004048	B2M
QG083-03	H06	HQP006171	NM_012423	RPL13A
QG083-03	H07	HQP009026	NM_000194	HPRT1
QG083-03	H08	HQP054253	NR_003286	RN18S1
QG083-03	H09	RT		
QG083-03	H10	RT		
QG083-03	H11	PCR		
QG083-03	H12	PCR		
QG083-04	A01	HQP018966	NM_001080124	CASP8
QG083-04	A02	HQP020427	NM_001226	CASP6
QG083-04	A03	HQP020249	NM_032982	CASP2
QG083-04	A04	HQP000781	NM_006549	CAMKK2
QG083-04	A05	HQP019841	NM_004343	CALR
QG083-04	A06	HQP000445	NM_006085	BPNT1
QG083-04	A07	HQP017234	NM_005180	BMI1
QG083-04	A08	HQP016653	NM_001196	BID
QG083-04	A09	HQP016418	NM_004327	BCR
QG083-04	A10	HQP016211	NM_000633	BCL2
QG083-04	A11	HQP015964	NM_004324	BAX

QG083-04	A12	HQP015917	NM_001188	BAK1
QG083-04	B01	HQP014386	NM_000707	AVPR1B
QG083-04	B02	HQP016578	NM_000332	ATXN1
QG083-04	B03	HQP013091	NM_001003696	ATP5J
QG083-04	B04	HQP011909	NM_001681	ATP2A2
QG083-04	B05	HQP011806	NM_152296	ATP1A3
QG083-04	B06	HQP011736	NM_000051	ATM
QG083-04	B07	HQP014637	NM_018179	ATF7IP
QG083-04	B08	HQP005471	NM_012068	ATF5
QG083-04	B09	HQP011497	NM_000049	ASPA
QG083-04	B10	HQP009801	NM_000044	AR
QG083-04	B11	HQP006348	NM_030882	APOL2
QG083-04	B12	HQP009556	NM_000041	APOE
QG083-04	C01	HQP008934	NM_001160	APAF1
QG083-04	C02	HQP013821	NM_017747	ANKHD1
QG083-04	C03	HQP006359	NM_000698	ALOX5
QG083-04	C04	HQP006356	NM_000697	ALOX12
QG083-04	C05	HQP019261	NM_001077690	ALG9
QG083-04	C06	HQP004991	NM_001014431	AKT1
QG083-04	C07	HQP017203	NM_031279	AGXT2L1
QG083-04	C08	HQP003928	NM_001126	ADSS
QG083-04	C09	HQP003848	NM_005160	ADRBK2
QG083-04	C10	HQP001836	NM_001116	ADCY9
QG083-04	C11	HQP021691	NM_001103	ACTN2
QG083-04	C12	HQP004081	NM_000789	ACE
QG083-04	D01	HQP012188	NM_002580	REG3A
QG083-04	D02	HQP022179	NM_004711	SYNGR1
QG083-04	D03	HQP017910	NM_005992	TBX1
QG083-04	D04	HQP021724	NM_003920	TIMELESS
QG083-04	D05	HQP011291	NM_001008228	MOG
QG083-04	D06	HQP000268	NM_005806	OLIG2
QG083-04	D07	HQP013033	NM_006211	PENK
QG083-04	D08	HQP014583	NM_002731	PRKACB
QG083-04	D09	HQP021643	NM_003895	SYNJ1
QG083-04	D10	HQP018154	NM_003283	TNNT1
QG083-04	D11	HQP022987	NM_004915	ABCG1
QG083-04	D12	HQP000604	NM_001033049	ADARB1
QG083-04	E01	HQP001925	NM_001117	ADCYAP1
QG083-04	E02	HQP003417	NM_000679	ADRA1B
QG083-04	E03	HQP021431	NM_005090	PLA2G4B
QG083-04	E04	HQP012286	NM_001001973	ATP5C1
QG083-04	E05	HQP013049	NM_001002258	ATP5G3
QG083-04	E06	HQP016545	NM_001709	BDNF

QG083-04	E07	HQP015425	NM_020412	CHMP1B
QG083-04	E08	HQP007944	NM_178034	PLA2G4D
QG083-04	E09	HQP006178	NM_013246	CLCF1
QG083-04	E10	HQP002962	NM_000756	CRH
QG083-04	E11	HQP004331	NM_000793	DIO2
QG083-04	E12	HQP004407	NM_001379	DNMT1
QG083-04	F01	HQP004448	NM_000796	DRD3
QG083-04	F02	HQP013311	NM_000304	PMP22
QG083-04	F03	HQP006604	NM_000808	GABRA3
QG083-04	F04	HQP008361	NM_000831	GRIK3
QG083-04	F05	HQP009025	NM_020995	HPR
QG083-04	F06	HQP006867	NM_025015	HSPA12A
QG083-04	F07	HQP009115	NM_000524	HTR1A
QG083-04	F08	HQP009126	NM_000870	HTR4
QG083-04	F09	HQP001124	NM_006839	IMMT
QG083-04	F10	HQP015090	NM_019842	KCNQ5
QG083-04	H05	HQP009026	NM_000194	HPRT0
QG083-04	H04	HQP054253	NR_003285	RN18S0
QG083-04	H03	RT		
QG083-04	H02	RT		
QG083-04	H01	PCR		
QG083-04	H00	PCR		
QG083-04	G05	HQP005996	NM_015277	NEDD4L
QG083-04	G06	HQP011941	NM_001007097	NTRK2
QG083-04	G07	HQP001343	NM_019094	NUDT4
QG083-04	G08	HQP001890	NM_138983	OLIG1
QG083-04	G09	HQP012146	NM_002573	PAFAH1B3
QG083-04	G10	HQP000737	NM_001011513	PDLIM5
QG083-04	G11	HQP013046	NM_002616	PER1
QG083-04	G12	HQP021641	NM_003894	PER2
QG083-04	H01	HGDC		
QG083-04	H02	HGDC		
QG083-04	H03	HQP006940	NM_002046	GAPDH
QG083-04	H04	HQP016381	NM_001101	ACTB
QG083-04	H05	HQP015171	NM_004048	B2M
QG083-04	H06	HQP006171	NM_012423	RPL13A
QG083-04	H07	HQP009026	NM_000194	HPRT1
QG083-04	H08	HQP054253	NR_003286	RN18S1
QG083-04	H09	RT		
QG083-04	H10	RT		
QG083-04	H11	PCR		
QG083-04	H12	PCR		