

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG004-01	A01	HQP004749	NM_005229	ELK1
QG004-01	A02	HQP009122	NM_000868	HTR2C
QG004-01	A03	HQP000623	NM_005194	CEBPB
QG004-01	A04	HQP009544	NM_000598	IGFBP3
QG004-01	A05	HQP013181	NM_000298	PKLR
QG004-01	A06	HQP009749	NM_000207	INS
QG004-01	A07	HQP018116	NM_138554	TLR4
QG004-01	A08	HQP012369	NM_015973	GAL
QG004-01	A09	HQP014547	NM_006253	PRKAB1
QG004-01	A10	HQP018114	NM_003264	TLR2
QG004-01	A11	HQP015986	NM_002862	PYGB
QG004-01	A12	HQP005934	NM_001013839	EXOC7
QG004-01	B01	HQP018743	NM_000721	CACNA1E
QG004-01	B02	HQP018723	NM_000719	CACNA1C
QG004-01	B03	HQP007230	NM_000583	GC
QG004-01	B04	HQP009147	NM_000415	IAPP
QG004-01	B05	HQP013115	NM_000293	PHKB
QG004-01	B06	HQP010847	NM_000237	LPL
QG004-01	B07	HQP007395	NM_000163	GHR
QG004-01	B08	HQP001136	NM_000015	NAT2
QG004-01	B09	HQP021438	NM_170662	CBLB
QG004-01	B10	HQP018402	NM_021833	UCP1
QG004-01	B11	HQP007396	NM_021081	GHRH
QG004-01	B12	HQP015244	NM_020167	NMUR2
QG004-01	C01	HQP009467	NM_000619	IFNG
QG004-01	C02	HQP008257	NM_001004056	GRK4
QG004-01	C03	HQP002424	NM_000614	CNTF
QG004-01	C04	HQP017962	NM_030756	TCF7L2
QG004-01	C05	HQP013077	NM_000289	PFKM
QG004-01	C06	HQP004676	NM_004095	EIF4EBP1
QG004-01	C07	HQP012219	NM_000280	PAX6
QG004-01	C08	HQP007239	NM_000162	GCK
QG004-01	C09	HQP015828	NM_002827	PTPN1
QG004-01	C10	HQP013151	NM_006219	PIK3CB
QG004-01	C11	HQP014907	NM_002755	MAP2K1
QG004-01	C12	HQP022870	NM_001775	CD38
QG004-01	D01	HQP004494	NM_000029	AGT
QG004-01	D02	HQP008970	NM_198834	ACACA
QG004-01	D03	HQP010011	NM_198580	SLC27A1
QG004-01	D04	HQP013155	NM_181504	PIK3R1
QG004-01	D05	HQP021662	NM_176875	CCKBR

QG004-01	D06	HQP012371	NM_032951	MLXIPL
QG004-01	D07	HQP020344	NM_031917	ANGPTL6
QG004-01	D08	HQP019539	NM_025072	PTGES2
QG004-01	D09	HQP020670	NM_021969	NR0B2
QG004-01	D10	HQP008906	NM_021784	FOXA2
QG004-01	D11	HQP001299	NM_021251	CAPN10
QG004-01	D12	HQP015857	NM_021158	TRIB3
QG004-01	E01	HQP015160	NM_020415	RETN
QG004-01	E02	HQP018498	NM_018727	TRPV1
QG004-01	E03	HQP001722	NM_018030	OSBPL1A
QG004-01	E04	HQP008273	NM_017572	MKNK2
QG004-01	E05	HQP013039	NM_017422	CALML5
QG004-01	E06	HQP012996	NM_016362	GHRL
QG004-01	E07	HQP002408	NM_016083	CNR1
QG004-01	E08	HQP006982	NM_015550	OSBPL3
QG004-01	E09	HQP023059	NM_014654	SDC3
QG004-01	E10	HQP006699	NM_012435	SHC2
QG004-01	E11	HQP006101	NM_012249	RHOQ
QG004-01	E12	HQP021442	NM_012079	DGAT1
QG004-01	F01	HQP018241	NM_007117	TRH
QG004-01	F02	HQP016063	NM_006908	RAC1
QG004-01	F03	HQP000957	NM_006656	NEU3
QG004-01	F04	HQP014731	NM_006254	PRKCD
QG004-01	F05	HQP014530	NM_006251	PRKAA1
QG004-01	F06	HQP013280	NM_006227	PLTP
QG004-01	F07	HQP011846	NM_006168	NKX6-1
QG004-01	F08	HQP018520	NM_006005	WFS1
QG004-01	F09	HQP000167	NM_005746	PBEF1
QG004-01	F10	HQP016276	NM_005614	RHEB
QG004-01	F11	HQP009062	NM_005525	HSD11B1
QG004-01	F12	HQP000001	NM_005465	AKT3
QG004-01	G01	HQP010627	NM_005357	LIPE
QG004-01	G02	HQP009036	NM_005343	HRAS
QG004-01	G03	HQP008082	NM_005297	MCHR1
QG004-01	G04	HQP019834	NM_005185	CALML3
QG004-01	G05	HQP011920	NM_005173	ATP2A3
QG004-01	G06	HQP013633	NM_005037	PPARG
QG004-01	G07	HQP009893	NM_004975	KCNB1
QG004-01	G08	HQP022625	NM_004797	ADIPOQ
QG004-01	G09	HQP008904	NM_004496	FOXA1
QG004-01	G10	HQP015538	NM_004322	BAD
QG004-01	G11	HQP022914	NM_004285	H6PD
QG004-01	G12	HQP016501	NM_004165	RRAD

QG004-01	H01	HGDC		
QG004-01	H02	HGDC		
QG004-01	H03	HQP006940	NM_002046	GAPDH
QG004-01	H04	HQP016381	NM_001101	ACTB
QG004-01	H05	HQP015171	NM_004048	B2M
QG004-01	H06	HQP006171	NM_012423	RPL13A
QG004-01	H07	HQP009026	NM_000194	HPRT1
QG004-01	H08	HQP054253	NR_003286	RN18S1
QG004-01	H09	RT		
QG004-01	H10	RT		
QG004-01	H11	PCR		
QG004-01	H12	PCR		
QG004-02	A01	HQP005093	NM_004103	PTK2B
QG004-02	A02	HQP004677	NM_004096	EIF4EBP2
QG004-02	A03	HQP021889	NM_003955	SOCS3
QG004-02	A04	HQP021602	NM_003877	SOCS2
QG004-02	A05	HQP021399	NM_003745	SOCS1
QG004-02	A06	HQP020912	NM_003604	IRS4
QG004-02	A07	HQP018490	NM_003381	VIP
QG004-02	A08	HQP018404	NM_003356	UCP3
QG004-02	A09	HQP018295	NM_003320	TUB
QG004-02	A10	HQP017080	NM_003029	SHC1
QG004-02	A11	HQP015921	NM_002840	PTPRF
QG004-02	A12	HQP015839	NM_002828	PTPN2
QG004-02	B01	HQP014900	NM_002753	MAPK10
QG004-02	B02	HQP014896	NM_002752	MAPK9
QG004-02	B03	HQP014886	NM_002750	MAPK8
QG004-02	B04	HQP014848	NM_002745	MAPK1
QG004-02	B05	HQP014718	NM_002738	PRKCB1
QG004-02	B06	HQP014650	NM_002734	PRKAR1A
QG004-02	B07	HQP014627	NM_002733	PRKAG1
QG004-02	B08	HQP014575	NM_002730	PRKACA
QG004-02	B09	HQP014024	NM_002711	PPP1R3A
QG004-02	B10	HQP013970	NM_002709	PPP1CB
QG004-02	B11	HQP013185	NM_002654	PKM2
QG004-02	B12	HQP011219	NM_002415	MIF
QG004-02	C01	HQP008894	NM_002131	HMGA1
QG004-02	C02	HQP008291	NM_002086	GRB2
QG004-02	C03	HQP007767	NM_002075	GNB3
QG004-02	C04	HQP004557	NM_001955	EDN1
QG004-02	C05	HQP000679	NM_001807	CEL
QG004-02	C06	HQP021313	NM_001753	CAV1
QG004-02	C07	HQP019707	NM_001743	CALM2

QG004-02	C08	HQP011909	NM_001681	ATP2A2
QG004-02	C09	HQP011303	NM_001672	ASIP
QG004-02	C10	HQP021320	NM_001234	CAV3
QG004-02	C11	HQP004453	NM_001138	AGRP
QG004-02	C12	HQP002765	NM_001124	ADM
QG004-02	D01	HQP011640	NM_001080779	MYO1C
QG004-02	D02	HQP018895	NM_001080432	FTO
QG004-02	D03	HQP005680	NM_001080421	UNC13A
QG004-02	D04	HQP017744	NM_001050	SSTR2
QG004-02	D05	HQP053963	NM_001048	SST
QG004-02	D06	HQP053960	NM_001042	SLC2A4
QG004-02	D07	HQP012722	NM_001040633	PRKAG2
QG004-02	D08	HQP012960	NM_001040092	ENPP2
QG004-02	D09	HQP023083	NM_001040059	CD68
QG004-02	D10	HQP023454	NM_001039802	CDC42
QG004-02	D11	HQP012432	NM_001039667	ANGPTL4
QG004-02	D12	HQP000695	NM_001034954	SORBS1
QG004-02	E01	HQP019178	NM_001033952	CALCA
QG004-02	E02	HQP014829	NM_001033581	PRKCZ
QG004-02	E03	HQP016623	NM_001024912	CEACAM1
QG004-02	E04	HQP012353	NM_001018073	PCK2
QG004-02	E05	HQP009754	NM_001017915	INPP5D
QG004-02	E06	HQP004991	NM_001014431	AKT1
QG004-02	E07	HQP053957	NM_001010	RPS6
QG004-02	E08	HQP013953	NM_001008709	PPP1CA
QG004-02	E09	HQP016472	NM_001007071	RPS6KB2
QG004-02	E10	HQP017703	NM_001005291	SREBF1
QG004-02	E11	HQP010582	NM_001003679	LEPR
QG004-02	E12	HQP013074	NM_001002021	PFKL
QG004-02	F01	HQP013617	NM_001001928	PPARA
QG004-02	F02	HQP018256	NM_001001188	TRPM2
QG004-02	F03	HQP015539	NM_000954	PTGDS
QG004-02	F04	HQP013467	NM_000939	POMC
QG004-02	F05	HQP011906	NM_000909	NPY1R
QG004-02	F06	HQP011874	NM_000905	NPY
QG004-02	F07	HQP004446	NM_000795	DRD2
QG004-02	F08	HQP021637	NM_000729	CCK
QG004-02	F09	HQP002200	NM_000647	CCR2
QG004-02	F10	HQP011866	NM_000625	NOS2A
QG004-02	F11	HQP009518	NM_000618	IGF1
QG004-02	F12	HQP009529	NM_000612	IGF2
QG004-02	G01	HQP022972	NM_000610	CD44
QG004-02	G02	HQP011868	NM_000603	NOS3

QG004-02	G03	HQP009539	NM_000596	IGFBP1
QG004-02	G04	HQP018265	NM_000548	TSC2
QG004-02	G05	HQP017954	NM_000545	TCF1
QG004-02	G06	HQP010015	NM_000525	KCNJ11
QG004-02	G07	HQP007385	NM_000515	GH1
QG004-02	G08	HQP008908	NM_000457	HNF4A
QG004-02	G09	HQP017794	NM_000455	STK11
QG004-02	G10	HQP008757	NM_000410	HFE
QG004-02	G11	HQP005406	NM_000382	ALDH3A2
QG004-02	G12	HQP018474	NM_000376	VDR
QG004-02	H01	HGDC		
QG004-02	H02	HGDC		
QG004-02	H03	HQP006940	NM_002046	GAPDH
QG004-02	H04	HQP016381	NM_001101	ACTB
QG004-02	H05	HQP015171	NM_004048	B2M
QG004-02	H06	HQP006171	NM_012423	RPL13A
QG004-02	H07	HQP009026	NM_000194	HPRT1
QG004-02	H08	HQP054253	NR_003286	RN18S1
QG004-02	H09	RT		
QG004-02	H10	RT		
QG004-02	H11	PCR		
QG004-02	H12	PCR		
QG004-03	A01	HQP018067	NM_000361	THBD
QG004-03	A02	HQP009764	NM_000208	INSR
QG004-03	A03	HQP008401	NM_000176	NR3C1
QG004-03	A04	HQP005074	NM_000134	FABP2
QG004-03	A05	HQP003904	NM_000103	CYP19A1
QG004-03	A06	HQP022821	NM_000072	CD36
QG004-03	A07	HQP009651	NM_000043	FAS
QG004-03	A08	HQP009556	NM_000041	APOE
QG004-03	A09	HQP009125	NM_000039	APOA1
QG004-03	A10	HQP003812	NM_000025	ADRB3
QG004-03	A11	HQP003791	NM_000024	ADRB2
QG004-03	A12	HQP010646	NM_203349	SHC4
QG004-03	B01	HQP006461	NM_178509	STXBP4
QG004-03	B02	HQP005328	NM_175061	JAZF1
QG004-03	B03	HQP004229	NM_173851	SLC30A8
QG004-03	B04	HQP003977	NM_173527	REM2
QG004-03	B05	HQP005830	NM_145693	LPIN1
QG004-03	B06	HQP020845	NM_032564	DGAT2
QG004-03	B07	HQP015989	NM_031885	BBS2
QG004-03	B08	HQP015979	NM_024649	BBS1
QG004-03	B09	HQP018897	NM_024090	ELOVL6

QG004-03	B10	HQP017235	NM_022896	LPIN3
QG004-03	B11	HQP008662	NM_021957	GYS2
QG004-03	B12	HQP000726	NM_020979	SH2B2
QG004-03	C01	HQP015796	NM_020918	GPAM
QG004-03	C02	HQP015674	NM_020761	KIAA1303
QG004-03	C03	HQP017348	NM_018979	WNK1
QG004-03	C04	HQP020036	NM_018848	MKKS
QG004-03	C05	HQP013698	NM_017636	TRPM4
QG004-03	C06	HQP014460	NM_017508	SOX6
QG004-03	C07	HQP013237	NM_016848	SHC3
QG004-03	C08	HQP012444	NM_016133	INSIG2
QG004-03	C09	HQP018790	NM_015120	ALMS1
QG004-03	C10	HQP023050	NM_014646	LPIN2
QG004-03	C11	HQP007675	NM_014495	ANGPTL3
QG004-03	C12	HQP001016	NM_013261	PPARGC1A
QG004-03	D01	HQP006080	NM_012238	SIRT1
QG004-03	D02	HQP001291	NM_007054	KIF3A
QG004-03	D03	HQP017626	NM_006939	SOS2
QG004-03	D04	HQP019580	NM_006888	CALM1
QG004-03	D05	HQP000998	NM_006681	NMU
QG004-03	D06	HQP000601	NM_006377	UNC13B
QG004-03	D07	HQP014804	NM_006257	PRKCQ
QG004-03	D08	HQP014541	NM_006252	PRKAA2
QG004-03	D09	HQP013150	NM_006218	PIK3CA
QG004-03	D10	HQP013128	NM_006216	SERPINE2
QG004-03	D11	HQP013118	NM_006213	PHKG1
QG004-03	D12	HQP011097	NM_005912	MC4R
QG004-03	E01	HQP000263	NM_005803	FLOT1
QG004-03	E02	HQP017625	NM_005633	SOS1
QG004-03	E03	HQP015988	NM_005609	PYGM
QG004-03	E04	HQP009788	NM_005544	IRS1
QG004-03	E05	HQP004727	NM_005502	ABCA1
QG004-03	E06	HQP017696	NM_005417	SRC
QG004-03	E07	HQP008298	NM_005312	RAPGEF1
QG004-03	E08	HQP007724	NM_005271	GLUD1
QG004-03	E09	HQP002989	NM_005207	CRKL
QG004-03	E10	HQP021430	NM_005188	CBL
QG004-03	E11	HQP019811	NM_005184	CALM3
QG004-03	E12	HQP017252	NM_005068	SIM1
QG004-03	F01	HQP014987	NM_005044	PRKX
QG004-03	F02	HQP013158	NM_005027	PIK3R2
QG004-03	F03	HQP013153	NM_005026	PIK3CD
QG004-03	F04	HQP022927	NM_004898	CLOCK

QG004-03	F05	HQP022800	NM_004846	EIF4E2
QG004-03	F06	HQP017800	NM_004603	STX1A
QG004-03	F07	HQP016470	NM_004586	RPS6KA3
QG004-03	F08	HQP016049	NM_004578	RAB4A
QG004-03	F09	HQP010077	NM_004521	KIF5B
QG004-03	F10	HQP005860	NM_004475	FLOT2
QG004-03	F11	HQP022970	NM_004291	CARTPT
QG004-03	F12	HQP022531	NM_004240	TRIP10
QG004-03	G01	HQP008600	NM_004130	GYG1
QG004-03	G02	HQP007406	NM_004123	GIP
QG004-03	G03	HQP005134	NM_004104	FASN
QG004-03	G04	HQP021527	NM_003825	SNAP23
QG004-03	G05	HQP021299	NM_003684	MKNK1
QG004-03	G06	HQP021127	NM_003629	PIK3R3
QG004-03	G07	HQP020830	NM_003597	KLF11
QG004-03	G08	HQP018403	NM_003355	UCP2
QG004-03	G09	HQP017845	NM_003177	SYK
QG004-03	G10	HQP017574	NM_003081	SNAP25
QG004-03	G11	HQP015873	NM_002833	PTPN9
QG004-03	G12	HQP015062	NM_002763	PROX1
QG004-03	H01	HGDC		
QG004-03	H02	HGDC		
QG004-03	H03	HQP006940	NM_002046	GAPDH
QG004-03	H04	HQP016381	NM_001101	ACTB
QG004-03	H05	HQP015171	NM_004048	B2M
QG004-03	H06	HQP006171	NM_012423	RPL13A
QG004-03	H07	HQP009026	NM_000194	HPRT1
QG004-03	H08	HQP054253	NR_003286	RN18S1
QG004-03	H09	RT		
QG004-03	H10	RT		
QG004-03	H11	PCR		
QG004-03	H12	PCR		
QG004-04	A01	HQP014672	NM_002735	PRKAR1B
QG004-04	A02	HQP013261	NM_002666	PLIN
QG004-04	A03	HQP013113	NM_002637	PHKA1
QG004-04	A04	HQP012979	NM_002613	PDPK1
QG004-04	A05	HQP012576	NM_002594	PCSK2
QG004-04	A06	HQP012342	NM_002591	PCK1
QG004-04	A07	HQP012080	NM_002556	OSBP
QG004-04	A08	HQP011914	NM_002524	NRAS
QG004-04	A09	HQP009808	NM_002210	ITGAV
QG004-04	A10	HQP008653	NM_002103	GYS1
QG004-04	A11	HQP007233	NM_002054	GCG

QG004-04	A12	HQP005747	NM_002015	FOXO1A
QG004-04	B01	HQP004675	NM_001968	EIF4E
QG004-04	B02	HQP021420	NM_001756	SERPINA6
QG004-04	B03	HQP004995	NM_001626	AKT2
QG004-04	B04	HQP003816	NM_001619	ADRBK1
QG004-04	B05	HQP009756	NM_001567	INPPL1
QG004-04	B06	HQP007245	NM_001486	GCKR
QG004-04	B07	HQP004978	NM_001432	EREG
QG004-04	B08	HQP009073	NM_001093	ACACB
QG004-04	B09	HQP015024	NM_000948	PRL
QG004-04	B10	HQP012717	NM_000922	PDE3B
QG004-04	B11	HQP011907	NM_000910	NPY2R
QG004-04	B12	HQP009523	NM_000875	IGF1R
QG004-04	C01	HQP004081	NM_000789	ACE
QG004-04	C02	HQP003817	NM_000773	CYP2E1
QG004-04	C03	HQP021647	NM_000730	CCKAR
QG004-04	C04	HQP018502	NM_000638	VTN
QG004-04	C05	HQP009120	NM_000621	HTR2A
QG004-04	C06	HQP012154	NM_000602	SERPINE1
QG004-04	C07	HQP010577	NM_000527	LDLR
QG004-04	C08	HQP009171	NM_000482	APOA4
QG004-04	C09	HQP016744	NM_000450	SELE
QG004-04	C10	HQP013473	NM_000446	PON1
QG004-04	C11	HQP012539	NM_000439	PCSK1
QG004-04	C12	HQP009218	NM_000384	APOB
QG004-04	D01	HQP018263	NM_000368	TSC1
QG004-04	D02	HQP015535	NM_000314	PTEN
QG004-04	D03	HQP013114	NM_000292	PHKA2
QG004-04	D04	HQP010626	NM_000236	LIPC
QG004-04	D05	HQP010581	NM_000230	LEP
QG004-04	D06	HQP009818	NM_000212	ITGB3
QG004-04	D07	HQP007408	NM_000164	GIPR
QG004-04	D08	HQP006508	NM_000151	G6PC
QG004-04	D09	HQP005047	NM_000477	ALB
QG004-04	D10	HQP014854	NM_001040056	MAPK3
QG004-04	D11	HQP013627	NM_006238	PPARD
QG004-04	D12	HQP011807	NM_003998	NFKB1
QG004-04	E01	HQP021550	NM_003839	TNFRSF11A
QG004-04	E02	HQP008469	NM_002093	GSK3B
QG004-04	E03	HQP009639	NM_001556	IKKB
QG004-04	E04	HQP005001	NM_001040275	ESR2
QG004-04	E05	HQP004998	NM_000125	ESR1
QG004-04	E06	HQP008810	NM_014598	SOCS7



QG004-04	E07	HQP006077	NM_012240	SIRT4
QG004-04	E08	HQP012957	NM_006208	ENPP1
QG004-04	E09	HQP023437	NM_005123	NR1H4
QG004-04	E10	HQP006426	NM_004958	FRAP1
QG004-04	E11	HQP017733	NM_004333	BRAF
QG004-04	E12	HQP014687	NM_004157	PRKAR2A
QG004-04	F01	HQP021548	NM_003837	FBP2
QG004-04	F02	HQP015987	NM_002863	PYGL
QG004-04	F03	HQP014798	NM_002742	PRKD1
QG004-04	F04	HQP005072	NM_001442	FABP4
QG004-04	F05	HQP018020	NM_001063	TF
QG004-04	F06	HQP002820	NM_000096	CP
QG004-04	F07	HQP016471	NM_003161	RPS6KB1
QG004-04	F08	HQP014909	NM_030662	MAP2K2
QG004-04	F09	HQP019090	NM_024551	ADIPOR2
QG004-04	F10	HQP012386	NM_015999	ADIPOR1
QG004-04	F11	HQP006038	NM_015316	PPP1R13B
QG004-04	F12	HQP014760	NM_006255	PRKCH
QG004-04	G01	HQP011547	NM_005957	MTHFR
QG004-04	G02	HQP000039	NM_005476	GNE
QG004-04	G03	HQP016630	NM_005408	CCL13
QG004-04	G04	HQP012695	NM_005020	PDE1C
QG004-04	G05	HQP009850	NM_004972	JAK2
QG004-04	G06	HQP016088	NM_002880	RAF1
QG004-04	G07	HQP014706	NM_002737	PRKCA
QG004-04	G08	HQP013078	NM_002627	PFKP
QG004-04	G09	HQP007708	NM_002062	GLP1R
QG004-04	G10	HQP002847	NM_001873	CPE
QG004-04	G11	HQP009832	NM_001654	ARAF
QG004-04	G12	HQP006375	NM_001629	ALOX5AP
QG004-04	H01	HGDC		
QG004-04	H02	HGDC		
QG004-04	H03	HQP006940	NM_002046	GAPDH
QG004-04	H04	HQP016381	NM_001101	ACTB
QG004-04	H05	HQP015171	NM_004048	B2M
QG004-04	H06	HQP006171	NM_012423	RPL13A
QG004-04	H07	HQP009026	NM_000194	HPRT1
QG004-04	H08	HQP054253	NR_003286	RN18S1
QG004-04	H09	RT		
QG004-04	H10	RT		
QG004-04	H11	PCR		
QG004-04	H12	PCR		