

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG091-01	A01	HQP004278	NM_133329	KCNG3
QG091-01	A02	HQP004260	NM_130770	HTR3C
QG091-01	A03	HQP004022	NM_145068	TRPV3
QG091-01	A04	HQP003549	NM_014406	CESK1
QG091-01	A05	HQP003354	NM_001002914	KCTD11
QG091-01	A06	HQP003207	NM_032735	BEST3
QG091-01	A07	HQP003059	NM_017662	TRPM6
QG091-01	A08	HQP003051	NM_183240	TMEM37
QG091-01	A09	HQP002662	NM_144633	KCNH8
QG091-01	A10	HQP002617	NM_152387	KCTD18
QG091-01	A11	HQP001927	NM_053054	CATSPER1
QG091-01	A12	HQP001887	NM_133445	GRIN3A
QG091-01	B01	HQP001877	NM_052961	SLC26A8
QG091-01	B02	HQP001498	NM_052874	STX1B2
QG091-01	B03	HQP022461	NM_138392	SHKBP1
QG091-01	B04	HQP021784	NM_139320	CHRFAM7A
QG091-01	B05	HQP021034	NM_032834	ALG10
QG091-01	B06	HQP020684	NM_001040107	HVCN1
QG091-01	B07	HQP020622	NM_032320	BTBD10
QG091-01	B08	HQP020377	NM_031954	KCTD10
QG091-01	B09	HQP020330	NM_032115	KCNK16
QG091-01	B10	HQP019824	NM_030779	KCNH6
QG091-01	B11	HQP019488	NM_001007470	TRPM3
QG091-01	B12	HQP018882	NM_024076	KCTD15
QG091-01	C01	HQP018880	NM_024074	TMEM38A
QG091-01	C02	HQP017522	NM_023930	KCTD14
QG091-01	C03	HQP016163	NM_021625	TRPV4
QG091-01	C04	HQP015783	NM_020897	HCN3
QG091-01	C05	HQP015679	NM_020768	KCTD16
QG091-01	C06	HQP015489	NM_020533	MCOLN1
QG091-01	C07	HQP015357	NM_020402	CHRNA10
QG091-01	C08	HQP015244	NM_020167	NMUR2
QG091-01	C09	HQP015199	NM_020122	KCMF1
QG091-01	C10	HQP015143	NM_022054	KCNK13
QG091-01	C11	HQP015065	NM_019841	TRPV5
QG091-01	C12	HQP014797	NM_018558	GABRQ
QG091-01	D01	HQP014717	NM_018398	CACNA2D3
QG091-01	D02	HQP014423	NM_018646	TRPV6
QG091-01	D03	HQP014269	NM_018298	MCOLN3
QG091-01	D04	HQP014107	NM_018112	TMEM38B
QG091-01	D05	HQP013745	NM_017682	BEST2

QG091-01	D06	HQP013734	NM_017672	TRPM7
QG091-01	D07	HQP013698	NM_017636	TRPM4
QG091-01	D08	HQP013697	NM_017634	KCTD9
QG091-01	D09	HQP013471	NM_018992	KCTD5
QG091-01	D10	HQP013412	NM_021161	KCNK10
QG091-01	D11	HQP013386	NM_053277	CLIC6
QG091-01	D12	HQP013318	NM_173160	FXYD4
QG091-01	E01	HQP013317	NM_014164	FXYD5
QG091-01	E02	HQP013316	NM_022003	FXYD6
QG091-01	E03	HQP013315	NM_022006	FXYD7
QG091-01	E04	HQP013255	NM_016929	CLIC5
QG091-01	E05	HQP013242	NM_017901	TPCN1
QG091-01	E06	HQP012707	NM_016113	TRPV2
QG091-01	E07	HQP012616	NM_016601	KCNK9
QG091-01	E08	HQP012439	NM_016121	KCTD3
QG091-01	E09	HQP012381	NM_015993	PLLP
QG091-01	E10	HQP008827	NM_014604	TAX1BP3
QG091-01	E11	HQP008787	NM_014591	KCNIP2
QG091-01	E12	HQP008553	NM_014555	TRPM5
QG091-01	F01	HQP007691	NM_014505	KCNMB4
QG091-01	F02	HQP007540	NM_139318	KCNH5
QG091-01	F03	HQP007497	NM_014407	KCNMB3
QG091-01	F04	HQP007496	NM_014405	CACNG4
QG091-01	F05	HQP007459	NM_014386	PKD2L2
QG091-01	F06	HQP007453	NM_007052	NOX1
QG091-01	F07	HQP007447	NM_001001485	ATP2C1
QG091-01	F08	HQP007069	NM_015638	TRPC4AP
QG091-01	F09	HQP006320	NM_080671	KCNE4
QG091-01	F10	HQP006307	NM_014331	SLC7A11
QG091-01	F11	HQP006287	NM_012282	KCNE1L
QG091-01	F12	HQP006160	NM_015353	KCTD2
QG091-01	G01	HQP005868	NM_015166	MLC1
QG091-01	G02	HQP005805	NM_001048210	CLCC1
QG091-01	G03	HQP001511	NM_014139	SCN11A
QG091-01	G04	HQP000579	NM_001042537	SLC9A6
QG091-01	G05	HQP000540	NM_006114	TOMM40
QG091-01	G06	HQP000420	NM_006071	PKDREJ
QG091-01	G07	HQP000081	NM_005691	ABCC9
QG091-01	G08	HQP000040	NM_005477	HCN4
QG091-01	G09	HQP000016	NM_005472	KCNE3
QG091-01	G10	HQP023462	NM_172201	KCNE2
QG091-01	G11	HQP023012	NM_006536	CLCA2
QG091-01	G12	HQP022739	NM_004823	KCNK6

QG091-01	H01	HGDC		
QG091-01	H02	HGDC		
QG091-01	H03	HQP006940	NM_002046	GAPDH
QG091-01	H04	HQP016381	NM_001101	ACTB
QG091-01	H05	HQP015171	NM_004048	B2M
QG091-01	H06	HQP006171	NM_012423	RPL13A
QG091-01	H07	HQP009026	NM_000194	HPRT1
QG091-01	H08	HQP054253	NR_003286	RN18S1
QG091-01	H09	RT		
QG091-01	H10	RT		
QG091-01	H11	PCR		
QG091-01	H12	PCR		
QG091-02	A01	HQP022515	NM_004770	KCNB2
QG091-02	A02	HQP022407	NM_001005505	CACNA2D2
QG091-02	A03	HQP022253	NM_006028	HTR3B
QG091-02	A04	HQP022139	NM_004700	KCNQ4
QG091-02	A05	HQP022123	NM_005446	P2RXL1
QG091-02	A06	HQP022088	NM_004685	MTMR6
QG091-02	A07	HQP021891	NM_004669	CLIC3
QG091-02	A08	HQP021818	NM_007332	TRPA1
QG091-02	A09	HQP021709	NM_018896	CACNA1G
QG091-02	A10	HQP021703	NM_001003406	CACNA1I
QG091-02	A11	HQP021386	NM_003740	KCNK5
QG091-02	A12	HQP021220	NM_003661	APOL1
QG091-02	B01	HQP019709	NM_003478	CUL5
QG091-02	B02	HQP019470	NM_001042543	GLRA3
QG091-02	B03	HQP018767	NM_057176	BSND
QG091-02	B04	HQP018498	NM_018727	TRPV1
QG091-02	B05	HQP018496	NM_004183	BEST1
QG091-02	B06	HQP018472	NM_005662	VDAC3
QG091-02	B07	HQP018470	NM_003374	VDAC1
QG091-02	B08	HQP018256	NM_001001188	TRPM2
QG091-02	B09	HQP018255	NM_004621	TRPC6
QG091-02	B10	HQP018253	NM_016179	TRPC4
QG091-02	B11	HQP018252	NM_003305	TRPC3
QG091-02	B12	HQP018251	NM_003304	TRPC1
QG091-02	C01	HQP018223	NM_199259	TPTE
QG091-02	C02	HQP018143	NM_021137	TNFAIP1
QG091-02	C03	HQP017900	NM_000544	TAP2
QG091-02	C04	HQP017899	NM_000593	TAP1
QG091-02	C05	HQP017460	NM_004594	SLC9A5
QG091-02	C06	HQP017459	NM_004174	SLC9A3
QG091-02	C07	HQP017457	NM_003048	SLC9A2

QG091-02	C08	HQP017456	NM_003047	SLC9A1
QG091-02	C09	HQP016608	NM_000336	SCNN1B
QG091-02	C10	HQP016605	NM_002977	SCN9A
QG091-02	C11	HQP016603	NM_002976	SCN7A
QG091-02	C12	HQP016601	NM_000335	SCN5A
QG091-02	D01	HQP016600	NM_174934	SCN4B
QG091-02	D02	HQP016598	NM_000334	SCN4A
QG091-02	D03	HQP016594	NM_004588	SCN2B
QG091-02	D04	HQP016592	NM_001040142	SCN2A
QG091-02	D05	HQP016591	NM_199037	SCN1B
QG091-02	D06	HQP016589	NM_006920	SCN1A
QG091-02	D07	HQP016535	NM_001036	RYR3
QG091-02	D08	HQP016532	NM_000540	RYR1
QG091-02	D09	HQP013264	NM_005971	FXYD3
QG091-02	D10	HQP013180	NM_000297	PKD2
QG091-02	D11	HQP012100	NM_002562	P2RX7
QG091-02	D12	HQP012095	NM_002559	P2RX3
QG091-02	E01	HQP012094	NM_002558	P2RX1
QG091-02	E02	HQP011252	NM_002420	TRPM1
QG091-02	E03	HQP010069	NM_002252	KCNS3
QG091-02	E04	HQP010063	NM_020697	KCNS2
QG091-02	E05	HQP010058	NM_002251	KCNS1
QG091-02	E06	HQP010054	NM_004519	KCNQ3
QG091-02	E07	HQP010049	NM_004518	KCNQ2
QG091-02	E08	HQP010046	NM_000218	KCNQ1
QG091-02	E09	HQP010043	NM_002250	KCNN4
QG091-02	E10	HQP010041	NM_002249	KCNN3
QG091-02	E11	HQP010039	NM_021614	KCNN2
QG091-02	E12	HQP010035	NM_004137	KCNMB1
QG091-02	F01	HQP010033	NM_001014797	KCNMA1
QG091-02	F02	HQP010027	NM_001017424	KCNK2
QG091-02	F03	HQP010024	NM_002245	KCNK1
QG091-02	F04	HQP010022	NM_002243	KCNJ15
QG091-02	F05	HQP010018	NM_002242	KCNJ13
QG091-02	F06	HQP010015	NM_000525	KCNJ11
QG091-02	F07	HQP010012	NM_004982	KCNJ8
QG091-02	F08	HQP010010	NM_002240	KCNJ6
QG091-02	F09	HQP010009	NM_000890	KCNJ5
QG091-02	F10	HQP010007	NM_004981	KCNJ4
QG091-02	F11	HQP010005	NM_002239	KCNJ3
QG091-02	F12	HQP010003	NM_000891	KCNJ2
QG091-02	G01	HQP010001	NM_000220	KCNJ1
QG091-02	G02	HQP009998	NM_000238	KCNH2

QG091-02	G03	HQP009990	NM_002238	KCNH1
QG091-02	G04	HQP009980	NM_002237	KCNG1
QG091-02	G05	HQP009969	NM_000219	KCNE1
QG091-02	G06	HQP009959	NM_004980	KCND3
QG091-02	G07	HQP009949	NM_012281	KCND2
QG091-02	G08	HQP009947	NM_004979	KCND1
QG091-02	G09	HQP009938	NM_001039574	KCNC4
QG091-02	G10	HQP009906	NM_139136	KCNC2
QG091-02	G11	HQP009900	NM_004976	KCNC1
QG091-02	G12	HQP009893	NM_004975	KCNB1
QG091-02	H01	HGDC		
QG091-02	H02	HGDC		
QG091-02	H03	HQP006940	NM_002046	GAPDH
QG091-02	H04	HQP016381	NM_001101	ACTB
QG091-02	H05	HQP015171	NM_004048	B2M
QG091-02	H06	HQP006171	NM_012423	RPL13A
QG091-02	H07	HQP009026	NM_000194	HPRT1
QG091-02	H08	HQP054253	NR_003286	RN18S1
QG091-02	H09	RT		
QG091-02	H10	RT		
QG091-02	H11	PCR		
QG091-02	H12	PCR		
QG091-03	A01	HQP009873	NM_002234	KCNA5
QG091-03	A02	HQP009869	NM_002233	KCNA4
QG091-03	A03	HQP009868	NM_002232	KCNA3
QG091-03	A04	HQP009844	NM_002224	ITPR3
QG091-03	A05	HQP009843	NM_002223	ITPR2
QG091-03	A06	HQP009842	NM_002222	ITPR1
QG091-03	A07	HQP009123	NM_000869	HTR3A
QG091-03	A08	HQP008380	NM_000836	GRIN2D
QG091-03	A09	HQP008376	NM_000835	GRIN2C
QG091-03	A10	HQP008371	NM_000833	GRIN2A
QG091-03	A11	HQP008367	NM_000832	GRIN1
QG091-03	A12	HQP008365	NM_002088	GRIK5
QG091-03	B01	HQP008363	NM_014619	GRIK4
QG091-03	B02	HQP008353	NM_021956	GRIK2
QG091-03	B03	HQP008340	NM_000830	GRIK1
QG091-03	B04	HQP008321	NM_001510	GRID2
QG091-03	B05	HQP008312	NM_017551	GRID1
QG091-03	B06	HQP008306	NM_000828	GRIA3
QG091-03	B07	HQP008303	NM_000826	GRIA2
QG091-03	B08	HQP008302	NM_000827	GRIA1
QG091-03	B09	HQP007719	NM_000824	GLRB

QG091-03	B10	HQP006666	NM_002042	GABRR1
QG091-03	B11	HQP006660	NM_014211	GABRP
QG091-03	B12	HQP006651	NM_000816	GABRG2
QG091-03	C01	HQP006647	NM_173536	GABRG1
QG091-03	C02	HQP006628	NM_000814	GABRB3
QG091-03	C03	HQP006624	NM_000813	GABRB2
QG091-03	C04	HQP006620	NM_000812	GABRB1
QG091-03	C05	HQP006613	NM_000810	GABRA5
QG091-03	C06	HQP006611	NM_000809	GABRA4
QG091-03	C07	HQP006601	NM_000807	GABRA2
QG091-03	C08	HQP006599	NM_000806	GABRA1
QG091-03	C09	HQP005239	NM_000139	MS4A2
QG091-03	C10	HQP003741	NM_000397	CYBB
QG091-03	C11	HQP002364	NM_001037329	CNGA4
QG091-03	C12	HQP002344	NM_005140	CNGA2
QG091-03	D01	HQP002330	NM_000087	CNGA1
QG091-03	D02	HQP002319	NM_001297	CNGB1
QG091-03	D03	HQP002099	NM_001293	CLNS1A
QG091-03	D04	HQP002044	NM_001289	CLIC2
QG091-03	D05	HQP002039	NM_001288	CLIC1
QG091-03	D06	HQP002021	NM_000085	CLCNKB
QG091-03	D07	HQP002010	NM_001287	CLCN7
QG091-03	D08	HQP002002	NM_001286	CLCN6
QG091-03	D09	HQP002001	NM_000084	CLCN5
QG091-03	D10	HQP001985	NM_001830	CLCN4
QG091-03	D11	HQP001982	NM_004366	CLCN2
QG091-03	D12	HQP001981	NM_000083	CLCN1
QG091-03	E01	HQP001678	NM_005199	CHRNA7
QG091-03	E02	HQP001673	NM_000080	CHRNE
QG091-03	E03	HQP001667	NM_000750	CHRNA4
QG091-03	E04	HQP001663	NM_000749	CHRNA3
QG091-03	E05	HQP001644	NM_000746	CHRNA7
QG091-03	E06	HQP001643	NM_000745	CHRNA5
QG091-03	E07	HQP001637	NM_000744	CHRNA4
QG091-03	E08	HQP001632	NM_000743	CHRNA3
QG091-03	E09	HQP001614	NM_000079	CHRNA1
QG091-03	E10	HQP001535	NM_000739	CHRM2
QG091-03	E11	HQP000948	NM_000492	CFTR
QG091-03	E12	HQP018799	NM_000725	CACNB3
QG091-03	F01	HQP018782	NM_000724	CACNB2
QG091-03	F02	HQP018777	NM_000723	CACNB1
QG091-03	F03	HQP018775	NM_000722	CACNA2D1
QG091-03	F04	HQP018758	NM_000069	CACNA1S

QG091-03	F05	HQP018752	NM_005183	CACNA1F
QG091-03	F06	HQP018743	NM_000721	CACNA1E
QG091-03	F07	HQP018734	NM_000720	CACNA1D
QG091-03	F08	HQP018723	NM_000719	CACNA1C
QG091-03	F09	HQP016388	NM_001194	HCN2
QG091-03	F10	HQP013383	NM_000053	ATP7B
QG091-03	F11	HQP012018	NM_000704	ATP4A
QG091-03	F12	HQP011991	NM_001001396	ATP2B4
QG091-03	G01	HQP011920	NM_005173	ATP2A3
QG091-03	G02	HQP011909	NM_001681	ATP2A2
QG091-03	G03	HQP011898	NM_004320	ATP2A1
QG091-03	G04	HQP011894	NM_001680	FXD2
QG091-03	G05	HQP011863	NM_001679	ATP1B3
QG091-03	G06	HQP011852	NM_001678	ATP1B2
QG091-03	G07	HQP011806	NM_152296	ATP1A3
QG091-03	G08	HQP011776	NM_000701	ATP1A1
QG091-03	G09	HQP009816	NM_001079528	ABCC6
QG091-03	G10	HQP009790	NM_020980	AQP9
QG091-03	G11	HQP009734	NM_001650	AQP4
QG091-03	G12	HQP009705	NM_000486	AQP2
QG091-03	H01	HGDC		
QG091-03	H02	HGDC		
QG091-03	H03	HQP006940	NM_002046	GAPDH
QG091-03	H04	HQP016381	NM_001101	ACTB
QG091-03	H05	HQP015171	NM_004048	B2M
QG091-03	H06	HQP006171	NM_012423	RPL13A
QG091-03	H07	HQP009026	NM_000194	HPRT1
QG091-03	H08	HQP054253	NR_003286	RN18S1
QG091-03	H09	RT		
QG091-03	H10	RT		
QG091-03	H11	PCR		
QG091-03	H12	PCR		
QG091-04	A01	HQP009682	NM_001649	SHROOM2
QG091-04	A02	HQP011144	NM_001095	ACCN2
QG091-04	A03	HQP010975	NM_001094	ACCN1
QG091-04	A04	HQP018704	NM_000068	CACNA1A
QG091-04	A05	HQP001597	NM_012125	CHRM5
QG091-04	A06	HQP001979	NM_001285	CLCA1
QG091-04	A07	HQP006638	NM_000815	GABRD
QG091-04	A08	HQP007712	NM_002063	GLRA2
QG091-04	A09	HQP016606	NM_006514	SCN10A
QG091-04	A10	HQP000115	NM_005714	KCNK7
QG091-04	A11	HQP000438	NM_006539	CACNG3

QG091-04	A12	HQP006085	NM_012284	KCNH3
QG091-04	B01	HQP007153	NM_012283	KCNG2
QG091-04	B02	HQP008796	NM_001034837	KCNIP1
QG091-04	B03	HQP013036	NM_017419	ACCN5
QG091-04	B04	HQP013644	NM_019098	CNGB3
QG091-04	B05	HQP014485	NM_017581	CHRNA9
QG091-04	B06	HQP019676	NM_001035003	KCNIP4
QG091-04	B07	HQP020992	NM_032790	TMEM142A
QG091-04	B08	HQP021783	NM_031460	KCNK17
QG091-04	B09	HQP021860	NM_033272	KCNH7
QG091-04	B10	HQP001888	NM_138690	GRIN3B
QG091-04	B11	HQP008854	NM_001156	ANXA7
QG091-04	B12	HQP011840	NM_001001787	ATP1B1
QG091-04	C01	HQP011935	NM_001001323	ATP2B1
QG091-04	C02	HQP012031	NM_000705	ATP4B
QG091-04	C03	HQP013334	NM_000052	ATP7A
QG091-04	C04	HQP015561	NM_001005367	TTYH1
QG091-04	C05	HQP016144	NM_031897	CACNG6
QG091-04	C06	HQP001581	NM_000741	CHRM4
QG091-04	C07	HQP019211	NM_024681	KCTD17
QG091-04	C08	HQP001658	NM_000748	CHRN2
QG091-04	C09	HQP001668	NM_000751	CHRND
QG091-04	C10	HQP001983	NM_001829	CLCN3
QG091-04	C11	HQP002013	NM_001042704	CLCNKA
QG091-04	C12	HQP002350	NM_001079878	CNGA3
QG091-04	D01	HQP006604	NM_000808	GABRA3
QG091-04	D02	HQP006617	NM_000811	GABRA6
QG091-04	D03	HQP006642	NM_004961	GABRE
QG091-04	D04	HQP006657	NM_033223	GABRG3
QG091-04	D05	HQP006672	NM_002043	GABRR2
QG091-04	D06	HQP007709	NM_000171	GLRA1
QG091-04	D07	HQP008309	NM_000829	GRIA4
QG091-04	D08	HQP008361	NM_000831	GRIK3
QG091-04	D09	HQP009864	NM_000217	KCNA1
QG091-04	D10	HQP009865	NM_004974	KCNA2
QG091-04	D11	HQP019721	NM_025250	TTYH3
QG091-04	D12	HQP009884	NM_031886	KCNA7
QG091-04	E01	HQP020406	NM_032034	SLC4A11
QG091-04	E02	HQP009926	NM_004977	KCNC3
QG091-04	E03	HQP009973	NM_002236	KCNF1
QG091-04	E04	HQP010013	NM_004983	KCNJ9
QG091-04	E05	HQP010014	NM_002241	KCNJ10
QG091-04	E06	HQP010021	NM_013348	KCNJ14



QG091-04	E07	HQP010023	NM_018658	KCNJ16
QG091-04	E08	HQP010030	NM_002246	KCNK3
QG091-04	E09	HQP010037	NM_002248	KCNN1
QG091-04	E10	HQP012096	NM_002560	P2RX4
QG091-04	E11	HQP012097	NM_002561	P2RX5
QG091-04	E12	HQP013263	NM_005031	FXYD1
QG091-04	F01	HQP016595	NM_001081676	SCN3A
QG091-04	F02	HQP018254	NM_012471	TRPC5
QG091-04	F03	HQP018471	NM_003375	VDAC2
QG091-04	F04	HQP018820	NM_003471	KCNAB1
QG091-04	F05	HQP021136	NM_003636	KCNAB2
QG091-04	F06	HQP021761	NM_004198	CHRNA6
QG091-04	F07	HQP021920	NM_016112	PKD2L1
QG091-04	F08	HQP022290	NM_004732	KCNAB3
QG091-04	F09	HQP023004	NM_004921	CLCA3
QG091-04	F10	HQP000302	NM_005832	KCNMB2
QG091-04	F11	HQP000439	NM_006078	CACNG2
QG091-04	F12	HQP005460	NM_012128	CLCA4
QG091-04	G01	HQP005619	NM_012226	P2RX2
QG091-04	G02	HQP006084	NM_012285	KCNH4
QG091-04	G03	HQP006899	NM_013943	CLIC4
QG091-04	G04	HQP007433	NM_014379	KCNV1
QG091-04	G05	HQP007494	NM_014404	CACNG5
QG091-04	G06	HQP001773	NM_052832	SLC26A7
QG091-04	G07	HQP012213	NM_033310	KCNK4
QG091-04	G08	HQP013222	NM_024047	NUDT9
QG091-04	G09	HQP013462	NM_018983	NOLA1
QG091-04	G10	HQP001780	NM_138444	KCTD12
QG091-04	G11	HQP015090	NM_019842	KCNQ5
QG091-04	G12	HQP015144	NM_022055	KCNK12
QG091-04	H01	HGDC		
QG091-04	H02	HGDC		
QG091-04	H03	HQP006940	NM_002046	GAPDH
QG091-04	H04	HQP016381	NM_001101	ACTB
QG091-04	H05	HQP015171	NM_004048	B2M
QG091-04	H06	HQP006171	NM_012423	RPL13A
QG091-04	H07	HQP009026	NM_000194	HPRT1
QG091-04	H08	HQP054253	NR_003286	RN18S1
QG091-04	H09	RT		
QG091-04	H10	RT		
QG091-04	H11	PCR		
QG091-04	H12	PCR		