

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG039-01	A01	HQP010053	NM_172109	KCNQ2
QG039-01	A02	HQP010042	NM_170782	KCNN3
QG039-01	A03	HQP002003	NM_021735	CLCN6
QG039-01	A04	HQP018779	NM_199248	CACNB1
QG039-01	A05	HQP009026	NM_000194	HPRT1
QG039-01	A06	HQP053962	NM_001045	SLC6A4
QG039-01	A07	HQP017431	NM_014229	SLC6A11
QG039-01	A08	HQP017549	NM_030807	SLC2A11
QG039-01	A09	HQP017350	NM_006516	SLC2A1
QG039-01	A10	HQP017318	NM_004171	SLC1A2
QG039-01	A11	HQP017312	NM_004170	SLC1A1
QG039-01	A12	HQP017478	NM_003053	SLC18A1
QG039-01	B01	HQP016605	NM_002977	SCN9A
QG039-01	B02	HQP016603	NM_002976	SCN7A
QG039-01	B03	HQP016601	NM_000335	SCN5A
QG039-01	B04	HQP016598	NM_000334	SCN4A
QG039-01	B05	HQP016597	NM_006922	SCN3A
QG039-01	B06	HQP016594	NM_004588	SCN2B
QG039-01	B07	HQP053902	NM_001037	SCN1B
QG039-01	B08	HQP016589	NM_006920	SCN1A
QG039-01	B09	HQP010069	NM_002252	KCNS3
QG039-01	B10	HQP010063	NM_020697	KCNS2
QG039-01	B11	HQP010058	NM_002251	KCNS1
QG039-01	B12	HQP022139	NM_004700	KCNQ4
QG039-01	C01	HQP010054	NM_004519	KCNQ3
QG039-01	C02	HQP010049	NM_004518	KCNQ2
QG039-01	C03	HQP010046	NM_000218	KCNQ1
QG039-01	C04	HQP010043	NM_002250	KCNN4
QG039-01	C05	HQP010041	NM_002249	KCNN3
QG039-01	C06	HQP010039	NM_021614	KCNN2
QG039-01	C07	HQP010035	NM_004137	KCNMB1
QG039-01	C08	HQP010010	NM_002240	KCNJ6
QG039-01	C09	HQP010009	NM_000890	KCNJ5
QG039-01	C10	HQP010007	NM_004981	KCNJ4
QG039-01	C11	HQP010005	NM_002239	KCNJ3
QG039-01	C12	HQP010015	NM_000525	KCNJ11
QG039-01	D01	HQP010001	NM_000220	KCNJ1
QG039-01	D02	HQP009998	NM_000238	KCNH2
QG039-01	D03	HQP009990	NM_002238	KCNH1
QG039-01	D04	HQP009980	NM_002237	KCNG1
QG039-01	D05	HQP009969	NM_000219	KCNE1

QG039-01	D06	HQP009949	NM_012281	KCND2
QG039-01	D07	HQP009947	NM_004979	KCND1
QG039-01	D08	HQP009900	NM_004976	KCNC1
QG039-01	D09	HQP022515	NM_004770	KCNB2
QG039-01	D10	HQP009893	NM_004975	KCNB1
QG039-01	D11	HQP009873	NM_002234	KCNA5
QG039-01	D12	HQP009869	NM_002233	KCNA4
QG039-01	E01	HQP009868	NM_002232	KCNA3
QG039-01	E02	HQP002044	NM_001289	CLIC2
QG039-01	E03	HQP002039	NM_001288	CLIC1
QG039-01	E04	HQP002010	NM_001287	CLCN7
QG039-01	E05	HQP002002	NM_001286	CLCN6
QG039-01	E06	HQP002001	NM_000084	CLCN5
QG039-01	E07	HQP001985	NM_001830	CLCN4
QG039-01	E08	HQP001982	NM_004366	CLCN2
QG039-01	E09	HQP001981	NM_000083	CLCN1
QG039-01	E10	HQP000948	NM_000492	CFTR
QG039-01	E11	HQP018799	NM_000725	CACNB3
QG039-01	E12	HQP018782	NM_000724	CACNB2
QG039-01	F01	HQP018777	NM_000723	CACNB1
QG039-01	F02	HQP018758	NM_000069	CACNA1S
QG039-01	F03	HQP018723	NM_000719	CACNA1C
QG039-01	F04	HQP018714	NM_000718	CACNA1B
QG039-01	F05	HQP012018	NM_000704	ATP4A
QG039-01	F06	HQP011841	NM_001677	ATP1B1
QG039-01	F07	HQP011776	NM_000701	ATP1A1
QG039-01	F08	HQP010975	NM_001094	ACCN1
QG039-01	F09	HQP018704	NM_000068	CACNA1A
QG039-01	F10	HQP001979	NM_001285	CLCA1
QG039-01	F11	HQP054021	NM_021007	SCN2A
QG039-01	F12	HQP010040	NM_170775	KCNN2
QG039-01	G01	HQP011145	NM_020039	ACCN2
QG039-01	G02	HQP011899	NM_173201	ATP2A1
QG039-01	G03	HQP017320	NM_004172	SLC1A3
QG039-01	G04	HQP006684	NM_013445	GAD1
QG039-01	G05	HQP000439	NM_006078	CACNG2
QG039-01	G06	HQP001983	NM_001829	CLCN3
QG039-01	G07	HQP009864	NM_000217	KCNA1
QG039-01	G08	HQP009865	NM_004974	KCNA2
QG039-01	G09	HQP001013	NM_016540	GPR83
QG039-01	G10	HQP009973	NM_002236	KCNF1
QG039-01	G11	HQP010037	NM_002248	KCNN1
QG039-01	G12	HQP017479	NM_003054	SLC18A2

QG039-01	H01	HGDC		
QG039-01	H02	HGDC		
QG039-01	H03	HQP006940	NM_002046	GAPDH
QG039-01	H04	HQP016381	NM_001101	ACTB
QG039-01	H05	HQP015171	NM_004048	B2M
QG039-01	H06	HQP006171	NM_012423	RPL13A
QG039-01	H07	HQP009026	NM_000194	HPRT1
QG039-01	H08	HQP054253	NR_003286	RN18S1
QG039-01	H09	RT		
QG039-01	H10	RT		
QG039-01	H11	PCR		
QG039-01	H12	PCR		