

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG087-01	A01	HQP009816	NM_001079528	ABCC6
QG087-01	A02	HQP000074	NM_001023587	ABCC5
QG087-01	A03	HQP009026	NM_000194	HPRT1
QG087-01	A04	HQP018470	NM_003374	VDAC1
QG087-01	A05	HQP017900	NM_000544	TAP2
QG087-01	A06	HQP017899	NM_000593	TAP1
QG087-01	A07	HQP007819	NM_016354	SLCO4A1
QG087-01	A08	HQP007820	NM_013272	SLCO3A1
QG087-01	A09	HQP001541	NM_007256	SLCO2B1
QG087-01	A10	HQP017486	NM_005630	SLCO2A1
QG087-01	A11	HQP007821	NM_019844	SLCO1B3
QG087-01	A12	HQP000720	NM_006446	SLCO1B1
QG087-01	B01	HQP001306	NM_014270	SLC7A9
QG087-01	B02	HQP006094	NM_182728	SLC7A8
QG087-01	B03	HQP021984	NM_003982	SLC7A7
QG087-01	B04	HQP053994	NM_003983	SLC7A6
QG087-01	B05	HQP006307	NM_014331	SLC7A11
QG087-01	B06	HQP017389	NM_014227	SLC5A4
QG087-01	B07	HQP017368	NM_000343	SLC5A1
QG087-01	B08	HQP017363	NM_002394	SLC3A2
QG087-01	B09	HQP017356	NM_000341	SLC3A1
QG087-01	B10	HQP022451	NM_033518	SLC38A5
QG087-01	B11	HQP013453	NM_018976	SLC38A2
QG087-01	B12	HQP002694	NM_001859	SLC31A1
QG087-01	C01	HQP017353	NM_006931	SLC2A3
QG087-01	C02	HQP017351	NM_000340	SLC2A2
QG087-01	C03	HQP017350	NM_006516	SLC2A1
QG087-01	C04	HQP008931	NM_001532	SLC29A2

QG087-01	C05	HQP053999	NM_004955	SLC29A1
QG087-01	C06	HQP016763	NM_022127	SLC28A3
QG087-01	C07	HQP022197	NM_004212	SLC28A2
QG087-01	C08	HQP022202	NM_004213	SLC28A1
QG087-01	C09	HQP001672	NM_080866	SLC22A9
QG087-01	C10	HQP022635	NM_004254	SLC22A8
QG087-01	C11	HQP017492	NM_021977	SLC22A3
QG087-01	C12	HQP017493	NM_003058	SLC22A2
QG087-01	D01	HQP017490	NM_003057	SLC22A1
QG087-01	D02	HQP019711	NM_025243	SLC19A3
QG087-01	D03	HQP000669	NM_006996	SLC19A2
QG087-01	D04	HQP017482	NM_194255	SLC19A1
QG087-01	D05	HQP053996	NM_004207	SLC16A3
QG087-01	D06	HQP017474	NM_006517	SLC16A2
QG087-01	D07	HQP017473	NM_003051	SLC16A1
QG087-01	D08	HQP017472	NM_021082	SLC15A2
QG087-01	D09	HQP017471	NM_005073	SLC15A1
QG087-01	D10	HQP017462	NM_000452	SLC10A2
QG087-01	D11	HQP017461	NM_003049	SLC10A1
QG087-01	D12	HQP054016	NM_017458	MVP
QG087-01	E01	HQP013383	NM_000053	ATP7B
QG087-01	E02	HQP009790	NM_020980	AQP9
QG087-01	E03	HQP009689	NM_198098	AQP1
QG087-01	E04	HQP016876	NM_022437	ABCG8
QG087-01	E05	HQP022745	NM_004827	ABCG2
QG087-01	E06	HQP006358	NM_001090	ABCF1
QG087-01	E07	HQP015971	NM_002858	ABCD3
QG087-01	E08	HQP005065	NM_000033	ABCD1
QG087-01	E09	HQP009817	NM_001171	ABCC6
QG087-01	E10	HQP000075	NM_005688	ABCC5

QG087-01	E11	HQP021459	NM_003786	ABCC3
QG087-01	E12	HQP002260	NM_000392	ABCC2
QG087-01	F01	HQP022719	NM_033226	ABCC12
QG087-01	F02	HQP021179	NM_032583	ABCC11
QG087-01	F03	HQP021787	NM_033450	ABCC10
QG087-01	F04	HQP011322	NM_004996	ABCC1
QG087-01	F05	HQP000076	NM_005689	ABCB6
QG087-01	F06	HQP009299	NM_178559	ABCB5
QG087-01	F07	HQP013101	NM_000443	ABCB4
QG087-01	F08	HQP021388	NM_003742	ABCB11
QG087-01	F09	HQP013100	NM_000927	ABCB1
QG087-01	F10	HQP000427	NM_080283	ABCA9
QG087-01	F11	HQP006441	NM_000350	ABCA4
QG087-01	F12	HQP005199	NM_001089	ABCA3
QG087-01	G01	HQP003777	NM_152701	ABCA13
QG087-01	G02	HQP004727	NM_005502	ABCA1
QG087-01	G03	HQP007090	NM_173076	ABCA12
QG087-01	G04	HQP004999	NM_001606	ABCA2
QG087-01	G05	HQP015972	NM_005050	ABCD4
QG087-01	G06	HQP013139	NM_001694	ATP6V0C
QG087-01	G07	HQP013334	NM_000052	ATP7A
QG087-01	G08	HQP022588	NM_004790	SLC22A6
QG087-01	G09	HQP000986	NM_006672	SLC22A7
QG087-01	G10	HQP000210	NM_014251	SLC25A13
QG087-01	G11	HQP017487	NM_005075	SLCO1A2
QG087-01	G12	HQP018471	NM_003375	VDAC2
QG087-01	H01	HGDC		
QG087-01	H02	HGDC		
QG087-01	H03	HQP006940	NM_002046	GAPDH
QG087-01	H04	HQP016381	NM_001101	ACTB

QG087-01	H05	HQP015171	NM_004048	B2M
QG087-01	H06	HQP006171	NM_012423	RPL13A
QG087-01	H07	HQP009026	NM_000194	HPRT1
QG087-01	H08	HQP054253	NR_003286	RN18S1
QG087-01	H09	RT		
QG087-01	H10	RT		
QG087-01	H11	PCR		
QG087-01	H12	PCR		