

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG044-01	A01	HQP001713	NM_001013031	SORCS1
QG044-01	A02	HQP008367	NM_000832	GRIN1
QG044-01	A03	HQP009026	NM_000194	HPRT1
QG044-01	A04	HQP001182	NM_020549	CHAT
QG044-01	A05	HQP018195	NM_004179	TPH1
QG044-01	A06	HQP017862	NM_001057	TACR2
QG044-01	A07	HQP017867	NM_001058	TACR1
QG044-01	A08	HQP017745	NM_001051	SSTR3
QG044-01	A09	HQP017744	NM_001050	SSTR2
QG044-01	A10	HQP015688	NM_020777	SORCS2
QG044-01	A11	HQP011907	NM_000910	NPY2R
QG044-01	A12	HQP011906	NM_000909	NPY1R
QG044-01	B01	HQP015244	NM_020167	NMUR2
QG044-01	B02	HQP000392	NM_006056	NMUR1
QG044-01	B03	HQP011007	NM_000240	MAOA
QG044-01	B04	HQP022253	NM_006028	HTR3B
QG044-01	B05	HQP009123	NM_000869	HTR3A
QG044-01	B06	HQP009120	NM_000621	HTR2A
QG044-01	B07	HQP009116	NM_000863	HTR1B
QG044-01	B08	HQP008743	NM_001526	HCRTR2
QG044-01	B09	HQP008464	NM_005314	GRPR
QG044-01	B10	HQP008302	NM_000827	GRIA1
QG044-01	B11	HQP001011	NM_053036	NPFFR2
QG044-01	B12	HQP008082	NM_005297	MCHR1
QG044-01	C01	HQP007235	NM_000161	GCH1
QG044-01	C02	HQP020973	NM_003614	GALR3
QG044-01	C03	HQP006820	NM_001480	GALR1
QG044-01	C04	HQP006683	NM_000817	GAD1
QG044-01	C05	HQP006666	NM_002042	GABRR1
QG044-01	C06	HQP014797	NM_018558	GABRQ
QG044-01	C07	HQP006660	NM_014211	GABRP
QG044-01	C08	HQP006651	NM_000816	GABRG2
QG044-01	C09	HQP006647	NM_173536	GABRG1
QG044-01	C10	HQP006624	NM_000813	GABRB2
QG044-01	C11	HQP006620	NM_000812	GABRB1
QG044-01	C12	HQP006601	NM_000807	GABRA2
QG044-01	D01	HQP006599	NM_000806	GABRA1
QG044-01	D02	HQP004446	NM_000795	DRD2
QG044-01	D03	HQP004445	NM_000794	DRD1
QG044-01	D04	HQP002671	NM_000754	COMT
QG044-01	D05	HQP001678	NM_005199	CHRNA

QG044-01	D06	HQP001673	NM_000080	CHRNE
QG044-01	D07	HQP001667	NM_000750	CHRN4
QG044-01	D08	HQP001644	NM_000746	CHRNA7
QG044-01	D09	HQP001643	NM_000745	CHRNA5
QG044-01	D10	HQP001637	NM_000744	CHRNA4
QG044-01	D11	HQP001632	NM_000743	CHRNA3
QG044-01	D12	HQP001614	NM_000079	CHRNA1
QG044-01	E01	HQP001562	NM_000740	CHRM3
QG044-01	E02	HQP001535	NM_000739	CHRM2
QG044-01	E03	HQP001526	NM_000738	CHRM1
QG044-01	E04	HQP054020	NM_020985	CHAT
QG044-01	E05	HQP021662	NM_176875	CCKBR
QG044-01	E06	HQP021647	NM_000730	CCKAR
QG044-01	E07	HQP017802	NM_001727	BRS3
QG044-01	E08	HQP020523	NM_003568	ANXA9
QG044-01	E09	HQP011345	NM_000665	ACHE
QG044-01	E10	HQP004553	NM_000663	ABAT
QG044-01	E11	HQP006638	NM_000815	GABRD
QG044-01	E12	HQP007238	NM_005258	GCHFR
QG044-01	F01	HQP007712	NM_002063	GLRA2
QG044-01	F02	HQP020492	NM_198179	GPR103
QG044-01	F03	HQP016788	NM_022146	NPFFR1
QG044-01	F04	HQP002517	NM_144773	PROKR2
QG044-01	F05	HQP008368	NM_007327	GRIN1
QG044-01	F06	HQP007913	NM_004248	PRLHR
QG044-01	F07	HQP001714	NM_052918	SORCS1
QG044-01	F08	HQP017746	NM_001052	SSTR4
QG044-01	F09	HQP018081	NM_000714	TSPO
QG044-01	F10	HQP021761	NM_004198	CHRNA6
QG044-01	F11	HQP001010	NM_004885	NPFFR2
QG044-01	F12	HQP001658	NM_000748	CHRN2
QG044-01	G01	HQP001668	NM_000751	CHRN1
QG044-01	G02	HQP004448	NM_000796	DRD3
QG044-01	G03	HQP006604	NM_000808	GABRA3
QG044-01	G04	HQP054022	NM_021990	GABRE
QG044-01	G05	HQP006672	NM_002043	GABRR2
QG044-01	G06	HQP007709	NM_000171	GLRA1
QG044-01	G07	HQP009124	NM_213621	HTR3A
QG044-01	G08	HQP006652	NM_198903	GABRG2
QG044-01	G09	HQP011851	NM_002511	NMBR
QG044-01	G10	HQP010839	NM_005169	PHOX2A
QG044-01	G11	HQP014388	NM_005972	PPYR1
QG044-01	G12	HQP016312	NM_021815	SLC5A7

QG044-01	H01	HGDC		
QG044-01	H02	HGDC		
QG044-01	H03	HQP006940	NM_002046	GAPDH
QG044-01	H04	HQP016381	NM_001101	ACTB
QG044-01	H05	HQP015171	NM_004048	B2M
QG044-01	H06	HQP006171	NM_012423	RPL13A
QG044-01	H07	HQP009026	NM_000194	HPRT1
QG044-01	H08	HQP054253	NR_003286	RN18S1
QG044-01	H09	RT		
QG044-01	H10	RT		
QG044-01	H11	PCR		
QG044-01	H12	PCR		