

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG042-01	A01	HQP021591	NM_201264	NRP2
QG042-01	A02	HQP008818	NM_004495	NRG1
QG042-01	A03	HQP011915	NM_001037132	NRCAM
QG042-01	A04	HQP009026	NM_000194	HPRT1
QG042-01	A05	HQP018575	NM_003405	YWHAH
QG042-01	A06	HQP018481	NM_003376	VEGFA
QG042-01	A07	HQP017767	NM_003150	STAT3
QG042-01	A08	HQP022584	NM_004787	SLIT2
QG042-01	A09	HQP017098	NM_000193	SHH
QG042-01	A10	HQP000613	NM_006378	SEMA4D
QG042-01	A11	HQP016553	NM_006272	S100B
QG042-01	A12	HQP016542	NM_014624	S100A6
QG042-01	B01	HQP015432	NM_007008	RTN4
QG042-01	B02	HQP016063	NM_006908	RAC1
QG042-01	B03	HQP013574	NM_006236	POU3F3
QG042-01	B04	HQP012219	NM_000280	PAX6
QG042-01	B05	HQP012212	NM_016734	PAX5
QG042-01	B06	HQP012210	NM_181461	PAX3
QG042-01	B07	HQP020816	NM_032521	PARD6B
QG042-01	B08	HQP015061	NM_019619	PARD3
QG042-01	B09	HQP012140	NM_000430	PAFAH1B1
QG042-01	B10	HQP000223	NM_014253	ODZ1
QG042-01	B11	HQP022735	NM_004822	NTN1
QG042-01	B12	HQP021589	NM_003872	NRP2
QG042-01	C01	HQP021597	NM_003873	NRP1
QG042-01	C02	HQP008820	NM_013957	NRG1
QG042-01	C03	HQP011904	NM_002522	NPTX1
QG042-01	C04	HQP011875	NM_024408	NOTCH2
QG042-01	C05	HQP011241	NM_005933	MLL
QG042-01	C06	HQP011151	NM_002397	MEF2C
QG042-01	C07	HQP009743	NM_002192	INHBA
QG042-01	C08	HQP006147	NM_012259	HEY2
QG042-01	C09	HQP009057	NM_005524	HES1
QG042-01	C10	HQP012862	NM_016596	HDAC7A
QG042-01	C11	HQP023167	NM_006037	HDAC4
QG042-01	C12	HQP007814	NM_000175	GPI
QG042-01	D01	HQP007752	NM_020988	GNAO1
QG042-01	D02	HQP007346	NM_000514	GDNF
QG042-01	D03	HQP005403	NM_002006	FGF2
QG042-01	D04	HQP023016	NM_005103	FEZ1
QG042-01	D05	HQP004969	NM_004448	ERBB2

QG042-01	D06	HQP004897	NM_001429	EP300
QG042-01	D07	HQP004599	NM_001963	EGF
QG042-01	D08	HQP004596	NM_004429	EFNB1
QG042-01	D09	HQP004516	NM_004423	DVL3
QG042-01	D10	HQP004446	NM_000795	DRD2
QG042-01	D11	HQP004445	NM_000794	DRD1
QG042-01	D12	HQP008111	NM_005618	DLL1
QG042-01	E01	HQP004342	NM_001365	DLG4
QG042-01	E02	HQP008456	NM_001511	CXCL1
QG042-01	E03	HQP001535	NM_000739	CHRM2
QG042-01	E04	HQP019636	NM_176096	CDK5RAP3
QG042-01	E05	HQP012938	NM_016408	CDK5RAP1
QG042-01	E06	HQP021622	NM_003885	CDK5R1
QG042-01	E07	HQP017477	NM_001720	BMP8B
QG042-01	E08	HQP053910	NM_130851	BMP4
QG042-01	E09	HQP017333	NM_001200	BMP2
QG042-01	E10	HQP015745	NM_001702	BAI1
QG042-01	E11	HQP023361	NM_014862	ARNT2
QG042-01	E12	HQP009556	NM_000041	APOE
QG042-01	F01	HQP009007	NM_001164	APBB1
QG042-01	F02	HQP006353	NM_004304	ALK
QG042-01	F03	HQP002833	NM_000675	ADORA2A
QG042-01	F04	HQP002799	NM_000674	ADORA1
QG042-01	F05	HQP011345	NM_000665	ACHE
QG042-01	F06	HQP011244	NM_004316	ASCL1
QG042-01	F07	HQP022320	NM_005448	BMP15
QG042-01	F08	HQP008368	NM_007327	GRIN1
QG042-01	F09	HQP011700	NM_000266	NDP
QG042-01	F10	HQP008772	NM_014587	SOX8
QG042-01	F11	HQP015433	NM_020532	RTN4
QG042-01	F12	HQP016545	NM_001709	BDNF
QG042-01	G01	HQP018166	NM_003285	TNR
QG042-01	G02	HQP005422	NM_004114	FGF13
QG042-01	G03	HQP006122	NM_012258	HEY1
QG042-01	G04	HQP009660	NM_000588	IL3
QG042-01	G05	HQP053984	NM_002391	MDK
QG042-01	G06	HQP011699	NM_002487	NDN
QG042-01	G07	HQP011771	NM_002500	NEUROD1
QG042-01	G08	HQP054071	NM_005450	NOG
QG042-01	G09	HQP011917	NM_005010	NRCAM
QG042-01	G10	HQP013585	NM_006237	POU4F1
QG042-01	G11	HQP015778	NM_002825	PTN
QG042-01	G12	HQP016368	NM_002941	ROBO1

QG042-01	H01	HGDC		
QG042-01	H02	HGDC		
QG042-01	H03	HQP006940	NM_002046	GAPDH
QG042-01	H04	HQP016381	NM_001101	ACTB
QG042-01	H05	HQP015171	NM_004048	B2M
QG042-01	H06	HQP006171	NM_012423	RPL13A
QG042-01	H07	HQP009026	NM_000194	HPRT1
QG042-01	H08	HQP054253	NR_003286	RN18S1
QG042-01	H09	RT		
QG042-01	H10	RT		
QG042-01	H11	PCR		
QG042-01	H12	PCR		