

Plate	Position	Catalog No. of	Accession No. of	Symbol
		Primer	Gene	
QG093-01	A01	HQP008818	NM_004495	NRG1
QG093-01	A02	HQP005430	NM_023107	FGFR1
QG093-01	A03	HQP009026	NM_000194	HPRT1
QG093-01	A04	HQP009653	NM_152872	FAS
QG093-01	A05	HQP018250	NM_016157	TRO
QG093-01	A06	HQP018175	NM_000546	TP53
QG093-01	A07	HQP018044	NM_000660	TGFB1
QG093-01	A08	HQP018043	NM_003236	TGFA
QG093-01	A09	HQP054009	NM_006070	TFG
QG093-01	A10	HQP017867	NM_001058	TACR1
QG093-01	A11	HQP017770	NM_003151	STAT4
QG093-01	A12	HQP017767	NM_003150	STAT3
QG093-01	B01	HQP017766	NM_005419	STAT2
QG093-01	B02	HQP017764	NM_007315	STAT1
QG093-01	B03	HQP015544	NM_000956	PTGER2
QG093-01	B04	HQP011945	NM_006180	NTRK2
QG093-01	B05	HQP011933	NM_002527	NTF3
QG093-01	B06	HQP003284	NM_138573	NRG4
QG093-01	B07	HQP022888	NM_013982	NRG2
QG093-01	B08	HQP008820	NM_013957	NRG1
QG093-01	B09	HQP021631	NM_022002	NR112
QG093-01	B10	HQP011907	NM_000910	NPY2R
QG093-01	B11	HQP011906	NM_000909	NPY1R
QG093-01	B12	HQP011874	NM_000905	NPY
QG093-01	C01	HQP021348	NM_003717	NPFF
QG093-01	C02	HQP011828	NM_002507	NGFR
QG093-01	C03	HQP011827	NM_002506	NGFB
QG093-01	C04	HQP011774	NM_000267	NF1
QG093-01	C05	HQP011760	NM_006157	NELL1
QG093-01	C06	HQP011597	NM_002467	MYC
QG093-01	C07	HQP011539	NM_005954	MT3
QG093-01	C08	HQP011151	NM_002397	MEF2C
QG093-01	C09	HQP010608	NM_002310	LIFR
QG093-01	C10	HQP010607	NM_002309	LIF
QG093-01	C11	HQP009674	NM_002184	IL6ST
QG093-01	C12	HQP009672	NM_000565	IL6R
QG093-01	D01	HQP009670	NM_000600	IL6
QG093-01	D02	HQP009642	NM_000877	IL1R1
QG093-01	D03	HQP009641	NM_000576	IL1B
QG093-01	D04	HQP009686	NM_001558	IL10RA
QG093-01	D05	HQP009685	NM_000572	IL10

QG093-01	D06	HQP009089	NM_001540	HSPB1
QG093-01	D07	HQP008464	NM_005314	GRPR
QG093-01	D08	HQP001011	NM_053036	NPFFR2
QG093-01	D09	HQP022878	NM_004877	GMFG
QG093-01	D10	HQP007740	NM_004124	GMFB
QG093-01	D11	HQP007376	NM_001496	GFRA3
QG093-01	D12	HQP007374	NM_001495	GFRA2
QG093-01	E01	HQP007346	NM_000514	GDNF
QG093-01	E02	HQP006820	NM_001480	GALR1
QG093-01	E03	HQP006451	NM_004960	FUS
QG093-01	E04	HQP000954	NM_006653	FRS3
QG093-01	E05	HQP054010	NM_006654	FRS2
QG093-01	E06	HQP006188	NM_005252	FOS
QG093-01	E07	HQP005416	NM_002010	FGF9
QG093-01	E08	HQP005403	NM_002006	FGF2
QG093-01	E09	HQP009651	NM_000043	FAS
QG093-01	E10	HQP018803	NM_003467	CXCR4
QG093-01	E11	HQP003691	NM_001337	CX3CR1
QG093-01	E12	HQP002967	NM_001882	CRHBP
QG093-01	F01	HQP002427	NM_001842	CNTFR
QG093-01	F02	HQP002424	NM_000614	CNTF
QG093-01	F03	HQP022955	NM_001250	CD40
QG093-01	F04	HQP021647	NM_000730	CCKAR
QG093-01	F05	HQP021445	NM_004352	CBLN1
QG093-01	F06	HQP016211	NM_000633	BCL2
QG093-01	F07	HQP015964	NM_004324	BAX
QG093-01	F08	HQP001980	NM_001118	ADCYAP1R1
QG093-01	F09	HQP002971	NM_004382	CRHR1
QG093-01	F10	HQP011094	NM_000529	MC2R
QG093-01	F11	HQP011934	NM_006179	NTF5
QG093-01	F12	HQP011959	NM_002531	NTSR1
QG093-01	G01	HQP015033	NM_004158	PSPN
QG093-01	G02	HQP018400	NM_003353	UCN
QG093-01	G03	HQP016545	NM_001709	BDNF
QG093-01	G04	HQP002962	NM_000756	CRH
QG093-01	G05	HQP005427	NM_015850	FGFR1
QG093-01	G06	HQP008741	NM_001524	HCRT
QG093-01	G07	HQP054012	NM_006986	MAGED1
QG093-01	G08	HQP007435	NM_014380	NGFRAP1
QG093-01	G09	HQP011940	NM_002529	NTRK1
QG093-01	G10	HQP013298	NM_006228	PNOC
QG093-01	G11	HQP014388	NM_005972	PPYR1
QG093-01	G12	HQP018249	NM_001039705	TRO

QG093-01	H01	HGDC		
QG093-01	H02	HGDC		
QG093-01	H03	HQP006940	NM_002046	GAPDH
QG093-01	H04	HQP016381	NM_001101	ACTB
QG093-01	H05	HQP015171	NM_004048	B2M
QG093-01	H06	HQP006171	NM_012423	RPL13A
QG093-01	H07	HQP009026	NM_000194	HPRT1
QG093-01	H08	HQP054253	NR_003286	RN18S1
QG093-01	H09	RT		
QG093-01	H10	RT		
QG093-01	H11	PCR		
QG093-01	H12	PCR		