

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG041-01	A01	HQP018291	NM_133379	TTN
QG041-01	A02	HQP012736	NM_016240	SCARA3
QG041-01	A03	HQP008289	NM_003996	GPX5
QG041-01	A04	HQP009026	NM_000194	HPRT1
QG041-01	A05	HQP018481	NM_003376	VEGFA
QG041-01	A06	HQP000708	NM_006440	TXNRD2
QG041-01	A07	HQP018289	NM_003319	TTN
QG041-01	A08	HQP000043	NM_005480	TROAP
QG041-01	A09	HQP018214	NM_000547	TPO
QG041-01	A10	HQP003061	NM_080725	SRXN1
QG041-01	A11	HQP017616	NM_000636	SOD2
QG041-01	A12	HQP017615	NM_000454	SOD1
QG041-01	B01	HQP005604	NM_012237	SIRT2
QG041-01	B02	HQP016819	NM_005410	SEPP1
QG041-01	B03	HQP054043	NM_203472	SELS
QG041-01	B04	HQP014706	NM_002737	PRKCA
QG041-01	B05	HQP014672	NM_002735	PRKAR1B
QG041-01	B06	HQP015732	NM_020820	PREX1
QG041-01	B07	HQP022952	NM_004905	PRDX6
QG041-01	B08	HQP018000	NM_005809	PRDX2
QG041-01	B09	HQP014309	NM_000944	PPP3CA
QG041-01	B10	HQP004317	NM_000903	NQO1
QG041-01	B11	HQP019025	NM_024505	NOX5
QG041-01	B12	HQP011868	NM_000603	NOS3
QG041-01	C01	HQP011866	NM_000625	NOS2A
QG041-01	C02	HQP020335	NM_003551	NME5
QG041-01	C03	HQP011693	NM_000433	NCF2
QG041-01	C04	HQP017417	NM_000265	NCF1
QG041-01	C05	HQP011587	NM_005375	MYB
QG041-01	C06	HQP011539	NM_005954	MT3
QG041-01	C07	HQP011523	NM_012331	MSRA
QG041-01	C08	HQP011309	NM_000250	MPO
QG041-01	C09	HQP011077	NM_000242	MBL2
QG041-01	C10	HQP010136	NM_006121	KRT1
QG041-01	C11	HQP009853	NM_002228	JUN
QG041-01	C12	HQP009749	NM_000207	INS
QG041-01	D01	HQP009678	NM_000584	IL8
QG041-01	D02	HQP009685	NM_000572	IL10
QG041-01	D03	HQP009097	NM_007355	HSP90AB1
QG041-01	D04	HQP008474	NM_000178	GSS
QG041-01	D05	HQP008380	NM_000836	GRIN2D

QG041-01	D06	HQP006685	NM_182701	GPX6
QG041-01	D07	HQP008288	NM_001509	GPX5
QG041-01	D08	HQP008285	NM_002085	GPX4
QG041-01	D09	HQP008282	NM_002084	GPX3
QG041-01	D10	HQP008281	NM_002083	GPX2
QG041-01	D11	HQP008279	NM_000581	GPX1
QG041-01	D12	HQP004150	NM_153002	GPR156
QG041-01	E01	HQP007574	NM_000169	GLA
QG041-01	E02	HQP007235	NM_000161	GCH1
QG041-01	E03	HQP005712	NM_021953	FOXMI
QG041-01	E04	HQP020089	NM_000502	EPX
QG041-01	E05	HQP004605	NM_005228	EGFR
QG041-01	E06	HQP004498	NM_004417	DUSP1
QG041-01	E07	HQP012142	NM_014080	DUOX2
QG041-01	E08	HQP054035	NM_175940	DUOX1
QG041-01	E09	HQP004342	NM_001365	DLG4
QG041-01	E10	HQP006223	NM_013974	DDAH2
QG041-01	E11	HQP001679	NM_134268	CYGB
QG041-01	E12	HQP018772	NM_007158	CSDE1
QG041-01	F01	HQP000331	NM_000389	CDKN1A
QG041-01	F02	HQP021692	NM_003914	CCNA1
QG041-01	F03	HQP021198	NM_003656	CAMK1
QG041-01	F04	HQP011770	NM_004045	ATOX1
QG041-01	F05	HQP010138	NM_001172	ARG2
QG041-01	F06	HQP009556	NM_000041	APOE
QG041-01	F07	HQP006356	NM_000697	ALOX12
QG041-01	F08	HQP054002	NM_005163	AKT1
QG041-01	F09	HQP007238	NM_005258	GCHFR
QG041-01	F10	HQP010851	NM_006151	LPO
QG041-01	F11	HQP017618	NM_003102	SOD3
QG041-01	F12	HQP020946	NM_001752	CAT
QG041-01	G01	HQP023439	NM_005125	CCS
QG041-01	G02	HQP003737	NM_000101	CYBA
QG041-01	G03	HQP053992	NM_003746	DYNLL1
QG041-01	G04	HQP023011	NM_004923	MTL5
QG041-01	G05	HQP001521	NM_007254	PNKP
QG041-01	G06	HQP006759	NM_181652	PRDX5
QG041-01	G07	HQP000461	NM_006093	PRG3
QG041-01	G08	HQP054040	NM_183079	PRNP
QG041-01	G09	HQP022982	NM_014245	RNF7
QG041-01	G10	HQP012737	NM_182826	SCARA3
QG041-01	G11	HQP021265	NM_033129	SCRT2
QG041-01	G12	HQP000143	NM_016276	SGK2

QG041-01	H01	HGDC		
QG041-01	H02	HGDC		
QG041-01	H03	HQP006940	NM_002046	GAPDH
QG041-01	H04	HQP016381	NM_001101	ACTB
QG041-01	H05	HQP015171	NM_004048	B2M
QG041-01	H06	HQP006171	NM_012423	RPL13A
QG041-01	H07	HQP009026	NM_000194	HPRT1
QG041-01	H08	HQP054253	NR_003286	RN18S1
QG041-01	H09	RT		
QG041-01	H10	RT		
QG041-01	H11	PCR		
QG041-01	H12	PCR		