

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG045-01	A01	HQP018291	NM_133379	TTN
QG045-01	A02	HQP012736	NM_016240	SCARA3
QG045-01	A03	HQP008289	NM_003996	GPX5
QG045-01	A04	HQP009026	NM_000194	HPRT1
QG045-01	A05	HQP000708	NM_006440	TXNRD2
QG045-01	A06	HQP018336	NM_003330	TXNRD1
QG045-01	A07	HQP020551	NM_032243	TXNDC2
QG045-01	A08	HQP018289	NM_003319	TTN
QG045-01	A09	HQP018214	NM_000547	TPO
QG045-01	A10	HQP000598	NM_006374	STK25
QG045-01	A11	HQP003061	NM_080725	SRXN1
QG045-01	A12	HQP017616	NM_000636	SOD2
QG045-01	B01	HQP017615	NM_000454	SOD1
QG045-01	B02	HQP005604	NM_012237	SIRT2
QG045-01	B03	HQP016980	NM_003019	SFTPD
QG045-01	B04	HQP016819	NM_005410	SEPP1
QG045-01	B05	HQP054043	NM_203472	SELS
QG045-01	B06	HQP018781	NM_012293	PXDN
QG045-01	B07	HQP015598	NM_000963	PTGS2
QG045-01	B08	HQP015596	NM_000962	PTGS1
QG045-01	B09	HQP015732	NM_020820	PREX1
QG045-01	B10	HQP022952	NM_004905	PRDX6
QG045-01	B11	HQP000655	NM_006406	PRDX4
QG045-01	B12	HQP001067	NM_006793	PRDX3
QG045-01	C01	HQP018000	NM_005809	PRDX2
QG045-01	C02	HQP012152	NM_002574	PRDX1
QG045-01	C03	HQP006987	NM_015553	PIP3-E
QG045-01	C04	HQP022118	NM_020992	PDLIM1
QG045-01	C05	HQP023391	NM_005109	OXSRI
QG045-01	C06	HQP019025	NM_024505	NOX5
QG045-01	C07	HQP011866	NM_000625	NOS2A
QG045-01	C08	HQP020335	NM_003551	NME5
QG045-01	C09	HQP011693	NM_000433	NCF2
QG045-01	C10	HQP017417	NM_000265	NCF1
QG045-01	C11	HQP011539	NM_005954	MT3
QG045-01	C12	HQP011523	NM_012331	MSRA
QG045-01	D01	HQP011314	NM_002437	MPV17
QG045-01	D02	HQP011309	NM_000250	MPO
QG045-01	D03	HQP011210	NM_004528	MGST3
QG045-01	D04	HQP011077	NM_000242	MBL2
QG045-01	D05	HQP010136	NM_006121	KRT1

QG045-01	D06	HQP008490	NM_001513	GSTZ1
QG045-01	D07	HQP008474	NM_000178	GSS
QG045-01	D08	HQP008473	NM_000637	GSR
QG045-01	D09	HQP006685	NM_182701	GPX6
QG045-01	D10	HQP008288	NM_001509	GPX5
QG045-01	D11	HQP008285	NM_002085	GPX4
QG045-01	D12	HQP008282	NM_002084	GPX3
QG045-01	E01	HQP008281	NM_002083	GPX2
QG045-01	E02	HQP008279	NM_000581	GPX1
QG045-01	E03	HQP004150	NM_153002	GPR156
QG045-01	E04	HQP005712	NM_021953	FOXMI
QG045-01	E05	HQP020089	NM_000502	EPX
QG045-01	E06	HQP004950	NM_001979	EPHX2
QG045-01	E07	HQP004498	NM_004417	DUSP1
QG045-01	E08	HQP012142	NM_014080	DUOX2
QG045-01	E09	HQP054035	NM_175940	DUOX1
QG045-01	E10	HQP004308	NM_014762	DHCR24
QG045-01	E11	HQP002957	NM_001013742	DGKK
QG045-01	E12	HQP001679	NM_134268	CYGB
QG045-01	F01	HQP018772	NM_007158	CSDE1
QG045-01	F02	HQP016626	NM_002985	CCL5
QG045-01	F03	HQP017619	NM_004052	BNIP3
QG045-01	F04	HQP011770	NM_004045	ATOX1
QG045-01	F05	HQP009556	NM_000041	APOE
QG045-01	F06	HQP008905	NM_001159	AOX1
QG045-01	F07	HQP000272	NM_021146	ANGPTL7
QG045-01	F08	HQP006356	NM_000697	ALOX12
QG045-01	F09	HQP005047	NM_000477	ALB
QG045-01	F10	HQP010851	NM_006151	LPO
QG045-01	F11	HQP017618	NM_003102	SOD3
QG045-01	F12	HQP020946	NM_001752	CAT
QG045-01	G01	HQP023439	NM_005125	CCS
QG045-01	G02	HQP003737	NM_000101	CYBA
QG045-01	G03	HQP008506	NM_001518	GTF2I
QG045-01	G04	HQP023011	NM_004923	MTL5
QG045-01	G05	HQP001521	NM_007254	PNKP
QG045-01	G06	HQP006759	NM_181652	PRDX5
QG045-01	G07	HQP000461	NM_006093	PRG3
QG045-01	G08	HQP054040	NM_183079	PRNP
QG045-01	G09	HQP002882	NM_144651	PXDNL
QG045-01	G10	HQP022982	NM_014245	RNF7
QG045-01	G11	HQP012737	NM_182826	SCARA3
QG045-01	G12	HQP000143	NM_016276	SGK2

QG045-01	H01	HGDC		
QG045-01	H02	HGDC		
QG045-01	H03	HQP006940	NM_002046	GAPDH
QG045-01	H04	HQP016381	NM_001101	ACTB
QG045-01	H05	HQP015171	NM_004048	B2M
QG045-01	H06	HQP006171	NM_012423	RPL13A
QG045-01	H07	HQP009026	NM_000194	HPRT1
QG045-01	H08	HQP054253	NR_003286	RN18S1
QG045-01	H09	RT		
QG045-01	H10	RT		
QG045-01	H11	PCR		
QG045-01	H12	PCR		