

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG026-01	A01	HQP017693	NM_178313	SPTBN1
QG026-01	A02	HQP013618	NM_032644	PPARA
QG026-01	A03	HQP009675	NM_175767	IL6ST
QG026-01	A04	HQP009026	NM_000194	HPRT1
QG026-01	A05	HQP020210	NM_033294	CASP1
QG026-01	A06	HQP018481	NM_003376	VEGFA
QG026-01	A07	HQP018403	NM_003355	UCP2
QG026-01	A08	HQP018298	NM_006000	TUBA4A
QG026-01	A09	HQP018280	NM_003312	TST
QG026-01	A10	HQP018064	NM_000360	TH
QG026-01	A11	HQP017692	NM_003128	SPTBN1
QG026-01	A12	HQP017587	NM_003089	SNRP70
QG026-01	B01	HQP053960	NM_001042	SLC2A4
QG026-01	B02	HQP017350	NM_006516	SLC2A1
QG026-01	B03	HQP000072	NM_005500	SAE1
QG026-01	B04	HQP053973	NM_001011	RPS7
QG026-01	B05	HQP016454	NM_002952	RPS2
QG026-01	B06	HQP016435	NM_000994	RPL32
QG026-01	B07	HQP016432	NM_000991	RPL28
QG026-01	B08	HQP016114	NM_000964	RARA
QG026-01	B09	HQP015241	NM_002795	PSMB3
QG026-01	B10	HQP014530	NM_006251	PRKAA1
QG026-01	B11	HQP054001	NM_005036	PPARA
QG026-01	B12	HQP021796	NM_001084	PLOD3
QG026-01	C01	HQP013204	NM_002658	PLAU
QG026-01	C02	HQP006987	NM_015553	PIP3-E
QG026-01	C03	HQP008954	NM_001161	NUDT2
QG026-01	C04	HQP011874	NM_000905	NPY
QG026-01	C05	HQP011873	NM_017617	NOTCH1
QG026-01	C06	HQP011866	NM_000625	NOS2A
QG026-01	C07	HQP011593	NM_002466	MYBL2
QG026-01	C08	HQP011539	NM_005954	MT3
QG026-01	C09	HQP011005	NM_000528	MAN2B1
QG026-01	C10	HQP010581	NM_000230	LEP
QG026-01	C11	HQP010570	NM_002299	LCT
QG026-01	C12	HQP010099	NM_000222	KIT
QG026-01	D01	HQP021301	NM_003685	KHSRP
QG026-01	D02	HQP021588	NM_003870	IQGAP1
QG026-01	D03	HQP009674	NM_002184	IL6ST
QG026-01	D04	HQP009670	NM_000600	IL6
QG026-01	D05	HQP009640	NM_000575	IL1A

QG026-01	D06	HQP009539	NM_000596	IGFBP1
QG026-01	D07	HQP009529	NM_000612	IGF2
QG026-01	D08	HQP000633	NM_006389	HYOU1
QG026-01	D09	HQP000630	NM_006388	HTATIP
QG026-01	D10	HQP008898	NM_002133	HMOX1
QG026-01	D11	HQP008843	NM_000189	HK2
QG026-01	D12	HQP014564	NM_017902	HIF1AN
QG026-01	E01	HQP008831	NM_001530	HIF1A
QG026-01	E02	HQP008729	NM_000518	HBB
QG026-01	E03	HQP008279	NM_000581	GPX1
QG026-01	E04	HQP007814	NM_000175	GPI
QG026-01	E05	HQP007743	NM_002067	GNA11
QG026-01	E06	HQP004903	NM_001430	EPAS1
QG026-01	E07	HQP004897	NM_001429	EP300
QG026-01	E08	HQP004860	NM_001428	ENO1
QG026-01	E09	HQP004570	NM_001402	EEF1A1
QG026-01	E10	HQP004537	NM_001397	ECE1
QG026-01	E11	HQP000649	NM_006400	DCTN2
QG026-01	E12	HQP003982	NM_001348	DAPK3
QG026-01	F01	HQP001679	NM_134268	CYGB
QG026-01	F02	HQP003388	NM_000100	CSTB
QG026-01	F03	HQP002921	NM_004380	CREBBP
QG026-01	F04	HQP002462	NM_000088	COL1A1
QG026-01	F05	HQP001309	NM_001275	CHGA
QG026-01	F06	HQP053981	NM_001791	CDC42
QG026-01	F07	HQP020208	NM_033292	CASP1
QG026-01	F08	HQP018641	NM_001738	CA1
QG026-01	F09	HQP009101	NM_001168	BIRC5
QG026-01	F10	HQP021270	NM_003670	BHLHB2
QG026-01	F11	HQP015964	NM_004324	BAX
QG026-01	F12	HQP023361	NM_014862	ARNT2
QG026-01	G01	HQP012432	NM_001039667	ANGPTL4
QG026-01	G02	HQP005957	NM_015239	AGTPBP1
QG026-01	G03	HQP002765	NM_001124	ADM
QG026-01	G04	HQP000663	NM_006412	AGPAT2
QG026-01	G05	HQP004443	NM_001938	DR1
QG026-01	G06	HQP004956	NM_000799	EPO
QG026-01	G07	HQP007660	NM_014484	MOCS3
QG026-01	G08	HQP020079	NM_003491	ARD1A
QG026-01	G09	HQP020946	NM_001752	CAT
QG026-01	G10	HQP000642	NM_006396	SSSCA1
QG026-01	G11	HQP016928	NM_152794	HIF3A
QG026-01	G12	HQP017106	NM_006849	PDIA2

QG026-01	H01	HGDC		
QG026-01	H02	HGDC		
QG026-01	H03	HQP006940	NM_002046	GAPDH
QG026-01	H04	HQP016381	NM_001101	ACTB
QG026-01	H05	HQP015171	NM_004048	B2M
QG026-01	H06	HQP006171	NM_012423	RPL13A
QG026-01	H07	HQP009026	NM_000194	HPRT1
QG026-01	H08	HQP054253	NR_003286	RN18S1
QG026-01	H09	RT		
QG026-01	H10	RT		
QG026-01	H11	PCR		
QG026-01	H12	PCR		