

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG024-01	A01	HQP008970	NM_198834	ACACA
QG024-01	A02	HQP013061	NM_004035	ACOX1
QG024-01	A03	HQP003812	NM_000025	ADRB3
QG024-01	A04	HQP004154	NM_001129	AEBP1
QG024-01	A05	HQP054002	NM_005163	AKT1
QG024-01	A06	HQP004995	NM_001626	AKT2
QG024-01	A07	HQP000001	NM_005465	AKT3
QG024-01	A08	HQP007968	NM_001145	ANG
QG024-01	A09	HQP009832	NM_001654	ARAF
QG024-01	A10	HQP016238	NM_138578	BCL2L1
QG024-01	A11	HQP017733	NM_004333	BRAF
QG024-01	A12	HQP000590	NM_006367	CAP1
QG024-01	B01	HQP021430	NM_005188	CBL
QG024-01	B02	HQP000615	NM_004364	CEBPA
QG024-01	B03	HQP004426	NM_001381	DOK1
QG024-01	B04	HQP021953	NM_003974	DOK2
QG024-01	B05	HQP019396	NM_024872	DOK3
QG024-01	B06	HQP001227	NM_007026	DUSP14
QG024-01	B07	HQP004650	NM_001414	EIF2B1
QG024-01	B08	HQP004676	NM_004095	EIF4EBP1
QG024-01	B09	HQP004974	NM_001983	ERCC1
QG024-01	B10	HQP005134	NM_004104	FASN
QG024-01	B11	HQP005224	NM_000507	FBP1
QG024-01	B12	HQP006188	NM_005252	FOS
QG024-01	C01	HQP006426	NM_004958	FRAP1
QG024-01	C02	HQP054010	NM_006654	FRS2
QG024-01	C03	HQP000954	NM_006653	FRS3
QG024-01	C04	HQP006508	NM_000151	G6PC
QG024-01	C05	HQP007239	NM_000162	GCK
QG024-01	C06	HQP007812	NM_005276	GPD1
QG024-01	C07	HQP008291	NM_002086	GRB2
QG024-01	C08	HQP008468	NM_019884	GSK3A
QG024-01	C09	HQP054075	NM_002093	GSK3B
QG024-01	C10	HQP008843	NM_000189	HK2
QG024-01	C11	HQP009036	NM_005343	HRAS
QG024-01	C12	HQP009523	NM_000875	IGF1R
QG024-01	D01	HQP009529	NM_000612	IGF2
QG024-01	D02	HQP009539	NM_000596	IGFBP1
QG024-01	D03	HQP009749	NM_000207	INS
QG024-01	D04	HQP009764	NM_000208	INSR
QG024-01	D05	HQP009788	NM_005544	IRS1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG024-01	D06	HQP020912	NM_003604	IRS4
QG024-01	D07	HQP009853	NM_002228	JUN
QG024-01	D08	HQP010577	NM_000527	LDLR
QG024-01	D09	HQP010581	NM_000230	LEP
QG024-01	D10	HQP014907	NM_002755	MAP2K1
QG024-01	D11	HQP014848	NM_002745	MAPK1
QG024-01	D12	HQP011697	NM_006153	NCK1
QG024-01	E01	HQP011866	NM_000625	NOS2A
QG024-01	E02	HQP011874	NM_000905	NPY
QG024-01	E03	HQP012354	NM_004563	PCK2
QG024-01	E04	HQP012979	NM_002613	PDPK1
QG024-01	E05	HQP013150	NM_006218	PIK3CA
QG024-01	E06	HQP013155	NM_181504	PIK3R1
QG024-01	E07	HQP013158	NM_005027	PIK3R2
QG024-01	E08	HQP013185	NM_002654	PKM2
QG024-01	E09	HQP013634	NM_015869	PPARG
QG024-01	E10	HQP013954	NM_002708	PPP1CA
QG024-01	E11	HQP014751	NM_002739	PRKCG
QG024-01	E12	HQP014771	NM_002740	PRKCI
QG024-01	F01	HQP015024	NM_000948	PRL
QG024-01	F02	HQP015828	NM_002827	PTPN1
QG024-01	F03	HQP015921	NM_002840	PTPRF
QG024-01	F04	HQP016088	NM_002880	RAF1
QG024-01	F05	HQP015160	NM_020415	RETN
QG024-01	F06	HQP016502	NM_006270	RRAS
QG024-01	F07	HQP005458	NM_012250	RRAS2
QG024-01	F08	HQP012154	NM_000602	SERPINE1
QG024-01	F09	HQP017080	NM_003029	SHC1
QG024-01	F10	HQP017350	NM_006516	SLC2A1
QG024-01	F11	HQP053960	NM_001042	SLC2A4
QG024-01	F12	HQP017625	NM_005633	SOS1
QG024-01	G01	HQP018042	NM_003235	TG
QG024-01	G02	HQP018402	NM_021833	UCP1
QG024-01	G03	HQP018481	NM_003376	VEGFA
QG024-01	G04	HQP008973	NM_198838	ACACA
QG024-01	G05	HQP009026	NM_000194	HPRT1
QG024-01	G06	HQP012353	NM_001018073	PCK2
QG024-01	G07	HQP012980	NM_031268	PDPK1
QG024-01	G08	HQP014829	NM_001033581	PRKCZ
QG024-01	G09	HQP017081	NM_183001	SHC1
QG024-01	G10	HQP000695	NM_001034954	SORBS1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG024-01	G11	HQP000623	NM_005194	CEBPB
QG024-01	G12	HQP006558	NM_002039	GAB1
QG024-01	H01	HGDC		
QG024-01	H02	HGDC		
QG024-01	H03	HQP006940	NM_002046	GAPDH
QG024-01	H04	HQP016381	NM_001101	ACTB
QG024-01	H05	HQP015171	NM_004048	B2M
QG024-01	H06	HQP006171	NM_012423	RPL13A
QG024-01	H07	HQP009026	NM_000194	HPRT1
QG024-01	H08	HQP054253	NR_003286	RN18S1
QG024-01	H09	RT		
QG024-01	H10	RT		
QG024-01	H11	PCR		
QG024-01	H12	PCR		