

Plate	Position	Catalog No. of	Accession No. of	Symbol
		Primer	Gene	
QG006-01	A01	HQP009026	NM_000194	HPRT1
QG006-01	A02	HQP018357	NM_003334	UBE1
QG006-01	A03	HQP018175	NM_000546	TP53
QG006-01	A04	HQP018033	NM_006286	TFDP2
QG006-01	A05	HQP018032	NM_007111	TFDP1
QG006-01	A06	HQP053990	NM_003352	SUMO1
QG006-01	A07	HQP017300	NM_005983	SKP2
QG006-01	A08	HQP008623	NM_013376	SERTAD1
QG006-01	A09	HQP016391	NM_002947	RPA3
QG006-01	A10	HQP016172	NM_005611	RBL2
QG006-01	A11	HQP016160	NM_002895	RBL1
QG006-01	A12	HQP016155	NM_002894	RBBP8
QG006-01	B01	HQP016131	NM_000321	RB1
QG006-01	B02	HQP016077	NM_002875	RAD51
QG006-01	B03	HQP016070	NM_002873	RAD17
QG006-01	B04	HQP015947	NM_002853	RAD1
QG006-01	B05	HQP054038	NM_182649	PCNA
QG006-01	B06	HQP011687	NM_002485	NBN
QG006-01	B07	HQP011320	NM_005590	MRE11A
QG006-01	B08	HQP011278	NM_002431	MNAT1
QG006-01	B09	HQP011232	NM_002417	MKI67
QG006-01	B10	HQP011109	NM_006739	MCM5
QG006-01	B11	HQP011108	NM_005914	MCM4
QG006-01	B12	HQP011107	NM_002388	MCM3
QG006-01	C01	HQP011106	NM_004526	MCM2
QG006-01	C02	HQP000552	NM_006341	MAD2L2
QG006-01	C03	HQP010957	NM_002358	MAD2L1
QG006-01	C04	HQP010125	NM_002266	KPNA2
QG006-01	C05	HQP023139	NM_014708	KNTC1
QG006-01	C06	HQP009138	NM_004507	HUS1
QG006-01	C07	HQP012794	NM_016426	GTSE1
QG006-01	C08	HQP008502	NM_005316	GTF2H1
QG006-01	C09	HQP004125	NM_001924	GADD45A
QG006-01	C10	HQP004528	NM_001950	E2F4
QG006-01	C11	HQP004406	NM_004945	DNM2
QG006-01	C12	HQP004162	NM_004399	DDX11
QG006-01	D01	HQP020782	NM_003590	CUL3
QG006-01	D02	HQP020788	NM_003591	CUL2
QG006-01	D03	HQP020798	NM_003592	CUL1
QG006-01	D04	HQP001894	NM_001827	CKS2
QG006-01	D05	HQP001884	NM_001826	CKS1B

QG006-01	D06	HQP001282	NM_001274	CHEK1
QG006-01	D07	HQP000418	NM_005192	CDKN3
QG006-01	D08	HQP000369	NM_000077	CDKN2A
QG006-01	D09	HQP000342	NM_004064	CDKN1B
QG006-01	D10	HQP000331	NM_000389	CDKN1A
QG006-01	D11	HQP000285	NM_001799	CDK7
QG006-01	D12	HQP000274	NM_001259	CDK6
QG006-01	E01	HQP012938	NM_016408	CDK5RAP1
QG006-01	E02	HQP021622	NM_003885	CDK5R1
QG006-01	E03	HQP000245	NM_000075	CDK4
QG006-01	E04	HQP000225	NM_001798	CDK2
QG006-01	E05	HQP023443	NM_004359	CDC34
QG006-01	E06	HQP023365	NM_001255	CDC20
QG006-01	E07	HQP023261	NM_001786	CDC2
QG006-01	E08	HQP053993	NM_003903	CDC16
QG006-01	E09	HQP021906	NM_001239	CCNH
QG006-01	E10	HQP021882	NM_004354	CCNG2
QG006-01	E11	HQP021857	NM_004060	CCNG1
QG006-01	E12	HQP021842	NM_001761	CCNF
QG006-01	F01	HQP021819	NM_001238	CCNE1
QG006-01	F02	HQP021754	NM_001759	CCND2
QG006-01	F03	HQP016204	NM_053056	CCND1
QG006-01	F04	HQP022141	NM_004701	CCNB2
QG006-01	F05	HQP021727	NM_031966	CCNB1
QG006-01	F06	HQP017753	NM_000059	BRCA2
QG006-01	F07	HQP017713	NM_007294	BRCA1
QG006-01	F08	HQP009101	NM_001168	BIRC5
QG006-01	F09	HQP016211	NM_000633	BCL2
QG006-01	F10	HQP015124	NM_016567	BCCIP
QG006-01	F11	HQP015964	NM_004324	BAX
QG006-01	F12	HQP013596	NM_001184	ATR
QG006-01	G01	HQP011736	NM_000051	ATM
QG006-01	G02	HQP008560	NM_013366	ANAPC2
QG006-01	G03	HQP006954	NM_005157	ABL1
QG006-01	G04	HQP022027	NM_004675	DIRAS3
QG006-01	G05	HQP008614	NM_013367	ANAPC4
QG006-01	G06	HQP021735	NM_005190	CCNC
QG006-01	G07	HQP021991	NM_001241	CCNT2
QG006-01	G08	HQP000312	NM_001260	CDK8
QG006-01	G09	HQP000382	NM_004936	CDKN2B
QG006-01	G10	HQP001397	NM_007194	CHEK2
QG006-01	G11	HQP017714	NM_007297	BRCA1
QG006-01	G12	HQP011737	NM_138292	ATM

QG006-01	H01	HGDC		
QG006-01	H02	HGDC		
QG006-01	H03	HQP006940	NM_002046	GAPDH
QG006-01	H04	HQP016381	NM_001101	ACTB
QG006-01	H05	HQP015171	NM_004048	B2M
QG006-01	H06	HQP006171	NM_012423	RPL13A
QG006-01	H07	HQP009026	NM_000194	HPRT1
QG006-01	H08	HQP054253	NR_003286	RN18S1
QG006-01	H09	RT		
QG006-01	H10	RT		
QG006-01	H11	PCR		
QG006-01	H12	PCR		