

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG053-01	A01	HQP006954	NM_005157	ABL1
QG053-01	A02	HQP004658	NM_001621	AHR
QG053-01	A03	HQP004995	NM_001626	AKT2
QG053-01	A04	HQP000460	NM_014885	ANAPC10
QG053-01	A05	HQP011736	NM_000051	ATM
QG053-01	A06	HQP017713	NM_007294	BRCA1
QG053-01	A07	HQP021701	NM_001237	CCNA2
QG053-01	A08	HQP021727	NM_031966	CCNB1
QG053-01	A09	HQP022141	NM_004701	CCNB2
QG053-01	A10	HQP016204	NM_053056	CCND1
QG053-01	A11	HQP021754	NM_001759	CCND2
QG053-01	A12	HQP021757	NM_001760	CCND3
QG053-01	B01	HQP021819	NM_001238	CCNE1
QG053-01	B02	HQP021857	NM_004060	CCNG1
QG053-01	B03	HQP021882	NM_004354	CCNG2
QG053-01	B04	HQP021906	NM_001239	CCNH
QG053-01	B05	HQP001120	NM_006835	CCNI
QG053-01	B06	HQP021572	NM_003858	CCNK
QG053-01	B07	HQP021276	NM_003672	CDC14A
QG053-01	B08	HQP021664	NM_001078645	CDC16
QG053-01	B09	HQP023261	NM_001786	CDC2
QG053-01	B10	HQP023365	NM_001255	CDC20
QG053-01	B11	HQP021443	NM_004661	CDC23
QG053-01	B12	HQP023385	NM_001789	CDC25A
QG053-01	C01	HQP023399	NM_004358	CDC25B
QG053-01	C02	HQP023409	NM_001790	CDC25C
QG053-01	C03	HQP023421	NM_001256	CDC27
QG053-01	C04	HQP021349	NM_003718	CDC2L5
QG053-01	C05	HQP001310	NM_007065	CDC37
QG053-01	C06	HQP020120	NM_003504	CDC45L
QG053-01	C07	HQP023354	NM_001254	CDC6
QG053-01	C08	HQP000225	NM_001798	CDK2
QG053-01	C09	HQP000245	NM_000075	CDK4
QG053-01	C10	HQP000274	NM_001259	CDK6
QG053-01	C11	HQP000285	NM_001799	CDK7
QG053-01	C12	HQP000321	NM_001261	CDK9
QG053-01	D01	HQP021574	NM_004196	CDKL1
QG053-01	D02	HQP000331	NM_000389	CDKN1A
QG053-01	D03	HQP000342	NM_004064	CDKN1B
QG053-01	D04	HQP000369	NM_000077	CDKN2A
QG053-01	D05	HQP000396	NM_001262	CDKN2C

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG053-01	D06	HQP000408	NM_001800	CDKN2D
QG053-01	D07	HQP000418	NM_005192	CDKN3
QG053-01	D08	HQP000842	NM_004344	CETN2
QG053-01	D09	HQP001282	NM_001274	CHEK1
QG053-01	D10	HQP001396	NM_001005735	CHEK2
QG053-01	D11	HQP020798	NM_003592	CUL1
QG053-01	D12	HQP004127	NM_004083	DDIT3
QG053-01	E01	HQP001314	NM_007068	DMC1
QG053-01	E02	HQP004498	NM_004417	DUSP1
QG053-01	E03	HQP004524	NM_005225	E2F1
QG053-01	E04	HQP004526	NM_004091	E2F2
QG053-01	E05	HQP004527	NM_001949	E2F3
QG053-01	E06	HQP004528	NM_001950	E2F4
QG053-01	E07	HQP004125	NM_001924	GADD45A
QG053-01	E08	HQP008745	NM_004964	HDAC1
QG053-01	E09	HQP009036	NM_005343	HRAS
QG053-01	E10	HQP009853	NM_002228	JUN
QG053-01	E11	HQP009854	NM_002229	JUNB
QG053-01	E12	HQP000552	NM_006341	MAD2L2
QG053-01	F01	HQP014848	NM_002745	MAPK1
QG053-01	F02	HQP011106	NM_004526	MCM2
QG053-01	F03	HQP011107	NM_002388	MCM3
QG053-01	F04	HQP011110	NM_005915	MCM6
QG053-01	F05	HQP011135	NM_002392	MDM2
QG053-01	F06	HQP022945	NM_016195	MPHOSPH1
QG053-01	F07	HQP011597	NM_002467	MYC
QG053-01	F08	HQP011914	NM_002524	NRAS
QG053-01	F09	HQP001189	NM_006999	POLS
QG053-01	F10	HQP015535	NM_000314	PTEN
QG053-01	F11	HQP016070	NM_002873	RAD17
QG053-01	F12	HQP016074	NM_006265	RAD21
QG053-01	G01	HQP016088	NM_002880	RAF1
QG053-01	G02	HQP016131	NM_000321	RB1
QG053-01	G03	HQP017887	NM_006284	TAF10
QG053-01	G04	HQP018032	NM_007111	TFDP1
QG053-01	G05	HQP018033	NM_006286	TFDP2
QG053-01	G06	HQP018175	NM_000546	TP53
QG053-01	G07	HQP008445	NM_001048201	UHRF1
QG053-01	G08	HQP018519	NM_003390	WEE1
QG053-01	G09	HQP018546	NM_000378	WT1
QG053-01	G10	HQP006205	NM_006732	FOSB

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG053-01	G11	HQP012070	NM_004153	ORC1L
QG053-01	G12	HQP022046	NM_004203	PKMYT1
QG053-01	H01	HGDC		
QG053-01	H02	HGDC		
QG053-01	H03	HQP006940	NM_002046	GAPDH
QG053-01	H04	HQP016381	NM_001101	ACTB
QG053-01	H05	HQP015171	NM_004048	B2M
QG053-01	H06	HQP006171	NM_012423	RPL13A
QG053-01	H07	HQP009026	NM_000194	HPRT1
QG053-01	H08	HQP054253	NR_003286	RN18S1
QG053-01	H09	RT		
QG053-01	H10	RT		
QG053-01	H11	PCR		
QG053-01	H12	PCR		