

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG038-01	A01	HQP016064	NM_018890	RAC1
QG038-01	A02	HQP009026	NM_000194	HPRT1
QG038-01	A03	HQP018175	NM_000546	TP53
QG038-01	A04	HQP010961	NM_005359	SMAD4
QG038-01	A05	HQP007802	NM_006142	SFN
QG038-01	A06	HQP016131	NM_000321	RB1
QG038-01	A07	HQP016088	NM_002880	RAF1
QG038-01	A08	HQP016063	NM_006908	RAC1
QG038-01	A09	HQP022952	NM_004905	PRDX6
QG038-01	A10	HQP012156	NM_002576	PAK1
QG038-01	A11	HQP011914	NM_002524	NRAS
QG038-01	A12	HQP011796	NM_004554	NFATC4
QG038-01	B01	HQP011597	NM_002467	MYC
QG038-01	B02	HQP011299	NM_005372	MOS
QG038-01	B03	HQP021299	NM_003684	MKNK1
QG038-01	B04	HQP011151	NM_002397	MEF2C
QG038-01	B05	HQP018812	NM_004635	MAPKAPK3
QG038-01	B06	HQP014896	NM_002752	MAPK9
QG038-01	B07	HQP006190	NM_012324	MAPK8IP2
QG038-01	B08	HQP014886	NM_002750	MAPK8
QG038-01	B09	HQP014876	NM_002748	MAPK6
QG038-01	B10	HQP014855	NM_002746	MAPK3
QG038-01	B11	HQP003133	NM_001315	MAPK14
QG038-01	B12	HQP014906	NM_002754	MAPK13
QG038-01	C01	HQP014895	NM_002751	MAPK11
QG038-01	C02	HQP014900	NM_002753	MAPK10
QG038-01	C03	HQP014848	NM_002745	MAPK1
QG038-01	C04	HQP011166	NM_005922	MAP3K4
QG038-01	C05	HQP011164	NM_002401	MAP3K3
QG038-01	C06	HQP000881	NM_006609	MAP3K2
QG038-01	C07	HQP053966	NM_005921	MAP3K1
QG038-01	C08	HQP014926	NM_145185	MAP2K7
QG038-01	C09	HQP014919	NM_002758	MAP2K6
QG038-01	C10	HQP014917	NM_002757	MAP2K5
QG038-01	C11	HQP016830	NM_003010	MAP2K4
QG038-01	C12	HQP014914	NM_002756	MAP2K3
QG038-01	D01	HQP014909	NM_030662	MAP2K2
QG038-01	D02	HQP021392	NM_021970	MAP2K1IP1
QG038-01	D03	HQP014907	NM_002755	MAP2K1
QG038-01	D04	HQP021615	NM_014238	KSR1
QG038-01	D05	HQP009853	NM_002228	JUN

QG038-01	D06	HQP009089	NM_001540	HSPB1
QG038-01	D07	HQP009083	NM_005347	HSPA5
QG038-01	D08	HQP009036	NM_005343	HRAS
QG038-01	D09	HQP008291	NM_002086	GRB2
QG038-01	D10	HQP006188	NM_005252	FOS
QG038-01	D11	HQP005015	NM_005239	ETS2
QG038-01	D12	HQP005014	NM_005238	ETS1
QG038-01	E01	HQP004749	NM_005229	ELK1
QG038-01	E02	HQP004612	NM_001964	EGR1
QG038-01	E03	HQP004605	NM_005228	EGFR
QG038-01	E04	HQP004524	NM_005225	E2F1
QG038-01	E05	HQP021547	NM_003836	DLK1
QG038-01	E06	HQP002921	NM_004380	CREBBP
QG038-01	E07	HQP002907	NM_004379	CREB1
QG038-01	E08	HQP002462	NM_000088	COL1A1
QG038-01	E09	HQP001708	NM_001278	CHUK
QG038-01	E10	HQP000408	NM_001800	CDKN2D
QG038-01	E11	HQP000369	NM_000077	CDKN2A
QG038-01	E12	HQP000356	NM_000076	CDKN1C
QG038-01	F01	HQP000342	NM_004064	CDKN1B
QG038-01	F02	HQP000331	NM_000389	CDKN1A
QG038-01	F03	HQP000274	NM_001259	CDK6
QG038-01	F04	HQP000245	NM_000075	CDK4
QG038-01	F05	HQP000225	NM_001798	CDK2
QG038-01	F06	HQP053981	NM_001791	CDC42
QG038-01	F07	HQP021819	NM_001238	CCNE1
QG038-01	F08	HQP021757	NM_001760	CCND3
QG038-01	F09	HQP021754	NM_001759	CCND2
QG038-01	F10	HQP016204	NM_053056	CCND1
QG038-01	F11	HQP022141	NM_004701	CCNB2
QG038-01	F12	HQP021727	NM_031966	CCNB1
QG038-01	G01	HQP021701	NM_001237	CCNA2
QG038-01	G02	HQP021692	NM_003914	CCNA1
QG038-01	G03	HQP017733	NM_004333	BRAF
QG038-01	G04	HQP002912	NM_001880	ATF2
QG038-01	G05	HQP009832	NM_001654	ARAF
QG038-01	G06	HQP000382	NM_004936	CDKN2B
QG038-01	G07	HQP053904	NM_078626	CDKN2C
QG038-01	G08	HQP001372	NM_007181	MAP4K1
QG038-01	G09	HQP004606	NM_201282	EGFR
QG038-01	G10	HQP014877	NM_002749	MAPK7
QG038-01	G11	HQP022425	NM_004759	MAPKAPK2
QG038-01	G12	HQP011059	NM_002382	MAX

QG038-01	H01	HGDC		
QG038-01	H02	HGDC		
QG038-01	H03	HQP006940	NM_002046	GAPDH
QG038-01	H04	HQP016381	NM_001101	ACTB
QG038-01	H05	HQP015171	NM_004048	B2M
QG038-01	H06	HQP006171	NM_012423	RPL13A
QG038-01	H07	HQP009026	NM_000194	HPRT1
QG038-01	H08	HQP054253	NR_003286	RN18S1
QG038-01	H09	RT		
QG038-01	H10	RT		
QG038-01	H11	PCR		
QG038-01	H12	PCR		