

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG015-01	A01	HQP054002	NM_005163	AKT1
QG015-01	A02	HQP004995	NM_001626	AKT2
QG015-01	A03	HQP000001	NM_005465	AKT3
QG015-01	A04	HQP009832	NM_001654	ARAF
QG015-01	A05	HQP011667	NM_005171	ATF1
QG015-01	A06	HQP002912	NM_001880	ATF2
QG015-01	A07	HQP015538	NM_004322	BAD
QG015-01	A08	HQP016211	NM_000633	BCL2
QG015-01	A09	HQP017733	NM_004333	BRAF
QG015-01	A10	HQP020297	NM_004346	CASP3
QG015-01	A11	HQP020648	NM_001229	CASP9
QG015-01	A12	HQP021430	NM_005188	CBL
QG015-01	B01	HQP016204	NM_053056	CCND1
QG015-01	B02	HQP001708	NM_001278	CHUK
QG015-01	B03	HQP002462	NM_000088	COL1A1
QG015-01	B04	HQP002907	NM_004379	CREB1
QG015-01	B05	HQP003277	NM_001895	CSNK2A1
QG015-01	B06	HQP004498	NM_004417	DUSP1
QG015-01	B07	HQP004504	NM_001946	DUSP6
QG015-01	B08	HQP004599	NM_001963	EGF
QG015-01	B09	HQP004605	NM_005228	EGFR
QG015-01	B10	HQP004612	NM_001964	EGR1
QG015-01	B11	HQP004675	NM_001968	EIF4E
QG015-01	B12	HQP004749	NM_005229	ELK1
QG015-01	C01	HQP004961	NM_004447	EPS8
QG015-01	C02	HQP009671	NM_000639	FASLG
QG015-01	C03	HQP006022	NM_002026	FN1
QG015-01	C04	HQP006188	NM_005252	FOS
QG015-01	C05	HQP005759	NM_001455	FOXO3A
QG015-01	C06	HQP008291	NM_002086	GRB2
QG015-01	C07	HQP008468	NM_019884	GSK3A
QG015-01	C08	HQP054075	NM_002093	GSK3B
QG015-01	C09	HQP004493	NM_001945	HBEGF
QG015-01	C10	HQP009036	NM_005343	HRAS
QG015-01	C11	HQP009639	NM_001556	IKBKB
QG015-01	C12	HQP009649	NM_000586	IL2
QG015-01	D01	HQP009849	NM_002227	JAK1
QG015-01	D02	HQP009853	NM_002228	JUN
QG015-01	D03	HQP010907	NM_000595	LTA
QG015-01	D04	HQP014907	NM_002755	MAP2K1
QG015-01	D05	HQP016830	NM_003010	MAP2K4

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG015-01	D06	HQP014926	NM_145185	MAP2K7
QG015-01	D07	HQP000881	NM_006609	MAP3K2
QG015-01	D08	HQP014848	NM_002745	MAPK1
QG015-01	D09	HQP014900	NM_002753	MAPK10
QG015-01	D10	HQP014855	NM_002746	MAPK3
QG015-01	D11	HQP014886	NM_002750	MAPK8
QG015-01	D12	HQP014896	NM_002752	MAPK9
QG015-01	E01	HQP021299	NM_003684	MKNK1
QG015-01	E02	HQP011258	NM_002423	MMP7
QG015-01	E03	HQP011792	NM_004555	NFATC3
QG015-01	E04	HQP011807	NM_003998	NFKB1
QG015-01	E05	HQP011914	NM_002524	NRAS
QG015-01	E06	HQP006291	NM_012346	NUP62
QG015-01	E07	HQP012847	NM_002607	PDGFA
QG015-01	E08	HQP012856	NM_002608	PDGFB
QG015-01	E09	HQP012866	NM_006206	PDGFRA
QG015-01	E10	HQP012979	NM_002613	PDPK1
QG015-01	E11	HQP013150	NM_006218	PIK3CA
QG015-01	E12	HQP013155	NM_181504	PIK3R1
QG015-01	F01	HQP013158	NM_005027	PIK3R2
QG015-01	F02	HQP013201	NM_000930	PLAT
QG015-01	F03	HQP013238	NM_002660	PLCG1
QG015-01	F04	HQP014115	NM_002715	PPP2CA
QG015-01	F05	HQP014706	NM_002737	PRKCA
QG015-01	F06	HQP015535	NM_000314	PTEN
QG015-01	F07	HQP016088	NM_002880	RAF1
QG015-01	F08	HQP053987	NM_002884	RAP1A
QG015-01	F09	HQP016125	NM_002890	RASA1
QG015-01	F10	HQP010229	NM_001664	RHOA
QG015-01	F11	HQP022404	NM_004755	RPS6KA5
QG015-01	F12	HQP016471	NM_003161	RPS6KB1
QG015-01	G01	HQP017080	NM_003029	SHC1
QG015-01	G02	HQP017696	NM_005417	SRC
QG015-01	G03	HQP017764	NM_007315	STAT1
QG015-01	G04	HQP017767	NM_003150	STAT3
QG015-01	G05	HQP017771	NM_003152	STAT5A
QG015-01	G06	HQP018175	NM_000546	TP53
QG015-01	G07	HQP020649	NM_032996	CASP9
QG015-01	G08	HQP009026	NM_000194	HPRT1
QG015-01	G09	HQP003278	NM_177560	CSNK2A1
QG015-01	G10	HQP004505	NM_022652	DUSP6

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG015-01	G11	HQP020713	NM_001004722	NCK2
QG015-01	G12	HQP016126	NM_022650	RASA1
QG015-01	H01	HGDC		
QG015-01	H02	HGDC		
QG015-01	H03	HQP006940	NM_002046	GAPDH
QG015-01	H04	HQP016381	NM_001101	ACTB
QG015-01	H05	HQP015171	NM_004048	B2M
QG015-01	H06	HQP006171	NM_012423	RPL13A
QG015-01	H07	HQP009026	NM_000194	HPRT1
QG015-01	H08	HQP054253	NR_003286	RN18S1
QG015-01	H09	RT		
QG015-01	H10	RT		
QG015-01	H11	PCR		
QG015-01	H12	PCR		