

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG047-01	A01	HQP000472	NM_001111	ADAR
QG047-01	A02	HQP054002	NM_005163	AKT1
QG047-01	A03	HQP004995	NM_001626	AKT2
QG047-01	A04	HQP000001	NM_005465	AKT3
QG047-01	A05	HQP009024	NM_000038	APC
QG047-01	A06	HQP015538	NM_004322	BAD
QG047-01	A07	HQP017988	NM_000061	BTK
QG047-01	A08	HQP020648	NM_001229	CASP9
QG047-01	A09	HQP016204	NM_053056	CCND1
QG047-01	A10	HQP022490	NM_000591	CD14
QG047-01	A11	HQP053981	NM_001791	CDC42
QG047-01	A12	HQP000342	NM_004064	CDKN1B
QG047-01	B01	HQP001708	NM_001278	CHUK
QG047-01	B02	HQP003277	NM_001895	CSNK2A1
QG047-01	B03	HQP003539	NM_001904	CTNNB1
QG047-01	B04	HQP014948	NM_002759	EIF2AK2
QG047-01	B05	HQP004674	NM_001417	EIF4B
QG047-01	B06	HQP004675	NM_001968	EIF4E
QG047-01	B07	HQP004676	NM_004095	EIF4EBP1
QG047-01	B08	HQP004749	NM_005229	ELK1
QG047-01	B09	HQP009671	NM_000639	FASLG
QG047-01	B10	HQP006188	NM_005252	FOS
QG047-01	B11	HQP005747	NM_002015	FOXO1A
QG047-01	B12	HQP005759	NM_001455	FOXO3A
QG047-01	C01	HQP006426	NM_004958	FRAP1
QG047-01	C02	HQP007411	NM_000165	GJA1
QG047-01	C03	HQP008296	NM_005311	GRB10
QG047-01	C04	HQP008291	NM_002086	GRB2
QG047-01	C05	HQP054075	NM_002093	GSK3B
QG047-01	C06	HQP009036	NM_005343	HRAS
QG047-01	C07	HQP009089	NM_001540	HSPB1
QG047-01	C08	HQP009518	NM_000618	IGF1
QG047-01	C09	HQP009523	NM_000875	IGF1R
QG047-01	C10	HQP053997	NM_004517	ILK
QG047-01	C11	HQP009773	NM_001569	IRAK1
QG047-01	C12	HQP009788	NM_005544	IRS1
QG047-01	D01	HQP009810	NM_002211	ITGB1
QG047-01	D02	HQP009853	NM_002228	JUN
QG047-01	D03	HQP014907	NM_002755	MAP2K1
QG047-01	D04	HQP014848	NM_002745	MAPK1
QG047-01	D05	HQP003133	NM_001315	MAPK14

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG047-01	D06	HQP014855	NM_002746	MAPK3
QG047-01	D07	HQP014886	NM_002750	MAPK8
QG047-01	D08	HQP011542	NM_001018025	MTCP1
QG047-01	D09	HQP011603	NM_002468	MYD88
QG047-01	D10	HQP011807	NM_003998	NFKB1
QG047-01	D11	HQP011810	NM_020529	NFKBIA
QG047-01	D12	HQP012156	NM_002576	PAK1
QG047-01	E01	HQP012866	NM_006206	PDGFRA
QG047-01	E02	HQP012923	NM_002610	PDK1
QG047-01	E03	HQP012931	NM_002611	PDK2
QG047-01	E04	HQP012979	NM_002613	PDPK1
QG047-01	E05	HQP013150	NM_006218	PIK3CA
QG047-01	E06	HQP013154	NM_002649	PIK3CG
QG047-01	E07	HQP013155	NM_181504	PIK3R1
QG047-01	E08	HQP013158	NM_005027	PIK3R2
QG047-01	E09	HQP014706	NM_002737	PRKCA
QG047-01	E10	HQP014718	NM_002738	PRKCB1
QG047-01	E11	HQP015535	NM_000314	PTEN
QG047-01	E12	HQP015639	NM_005607	PTK2
QG047-01	F01	HQP015878	NM_002834	PTPN11
QG047-01	F02	HQP016063	NM_006908	RAC1
QG047-01	F03	HQP016088	NM_002880	RAF1
QG047-01	F04	HQP016125	NM_002890	RASA1
QG047-01	F05	HQP016172	NM_005611	RBL2
QG047-01	F06	HQP016276	NM_005614	RHEB
QG047-01	F07	HQP010229	NM_001664	RHOA
QG047-01	F08	HQP016471	NM_003161	RPS6KB1
QG047-01	F09	HQP017080	NM_003029	SHC1
QG047-01	F10	HQP017625	NM_005633	SOS1
QG047-01	F11	HQP019839	NM_021966	TCL1A
QG047-01	F12	HQP001674	NM_001039661	TIRAP
QG047-01	G01	HQP018116	NM_138554	TLR4
QG047-01	G02	HQP013496	NM_019009	TOLLIP
QG047-01	G03	HQP018263	NM_000368	TSC1
QG047-01	G04	HQP018265	NM_000548	TSC2
QG047-01	G05	HQP021768	NM_003941	WASL
QG047-01	G06	HQP018575	NM_003405	YWHAH
QG047-01	G07	HQP000471	NM_001025107	ADAR
QG047-01	G08	HQP009026	NM_000194	HPRT1
QG047-01	G09	HQP020649	NM_032996	CASP9
QG047-01	G10	HQP003278	NM_177560	CSNK2A1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG047-01	G11	HQP004679	NM_004953	EIF4G1
QG047-01	G12	HQP014829	NM_001033581	PRKCZ
QG047-01	H01	HGDC		
QG047-01	H02	HGDC		
QG047-01	H03	HQP006940	NM_002046	GAPDH
QG047-01	H04	HQP016381	NM_001101	ACTB
QG047-01	H05	HQP015171	NM_004048	B2M
QG047-01	H06	HQP006171	NM_012423	RPL13A
QG047-01	H07	HQP009026	NM_000194	HPRT1
QG047-01	H08	HQP054253	NR_003286	RN18S1
QG047-01	H09	RT		
QG047-01	H10	RT		
QG047-01	H11	PCR		
QG047-01	H12	PCR		