

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG080-01	A01	HQP017733	NM_004333	BRAF
QG080-01	A02	HQP054045	BC004257	RET
QG080-01	A03	HQP016418	NM_004327	BCR
QG080-01	A04	HQP011914	NM_002524	NRAS
QG080-01	A05	HQP010133	NM_004985	KRAS
QG080-01	A06	HQP013150	NM_006218	PIK3CA
QG080-01	A07	HQP009036	NM_005343	HRAS
QG080-01	A08	HQP018564	NM_005432	XRCC3
QG080-01	A09	HQP018562	NM_006297	XRCC1
QG080-01	A10	HQP018175	NM_000546	TP53
QG080-01	A11	HQP004991	NM_001014431	AKT1
QG080-01	A12	HQP054047	BC008403	HLA-DRB1
QG080-01	B01	HQP018271	NM_000369	TSHR
QG080-01	B02	HQP018097	NM_003317	TITF1
QG080-01	B03	HQP018077	NM_003250	THRA
QG080-01	B04	HQP015925	NM_002843	PTPRJ
QG080-01	B05	HQP011235	NM_000249	MLH1
QG080-01	B06	HQP001136	NM_000015	NAT2
QG080-01	B07	HQP005704	NM_004473	FOXE1
QG080-01	B08	HQP022764	NM_004832	GSTO1
QG080-01	B09	HQP019650	NM_005436	CCDC6
QG080-01	B10	HQP018565	NM_003401	XRCC4
QG080-01	B11	HQP018475	NM_001025366	VEGFA
QG080-01	B12	HQP018042	NM_003235	TG
QG080-01	C01	HQP016088	NM_002880	RAF1
QG080-01	C02	HQP011938	NM_001007792	NTRK1
QG080-01	C03	HQP011597	NM_002467	MYC
QG080-01	C04	HQP011492	NM_002439	MSH3
QG080-01	C05	HQP010613	NM_002312	LIG4
QG080-01	C06	HQP007767	NM_002075	GNB3
QG080-01	C07	HQP004605	NM_005228	EGFR
QG080-01	C08	HQP003772	NM_000499	CYP1A1
QG080-01	C09	HQP016706	NM_003002	SDHD
QG080-01	C10	HQP007370	NM_005264	GFRA1
QG080-01	C11	HQP054052	BC015035	FGFR1
QG080-01	C12	HQP054035	NM_175940	DUOX1
QG080-01	D01	HQP023420	NM_005121	THRAP1
QG080-01	D02	HQP022741	NM_170751	CDYL
QG080-01	D03	HQP022530	NM_004239	TRIP11
QG080-01	D04	HQP022528	NM_004238	TRIP12
QG080-01	D05	HQP022428	NM_004760	STK17A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG080-01	D06	HQP022427	NM_004226	STK17B
QG080-01	D07	HQP021961	NM_003977	AIP
QG080-01	D08	HQP020440	NM_032119	GPR98
QG080-01	D09	HQP019661	NM_005437	NCOA4
QG080-01	D10	HQP018802	NM_001008540	CXCR4
QG080-01	D11	HQP018794	NM_003466	PAX8
QG080-01	D12	HQP018750	NM_006301	MAP3K12
QG080-01	E01	HQP018568	NM_021141	XRCC5
QG080-01	E02	HQP018563	NM_005431	XRCC2
QG080-01	E03	HQP018546	NM_000378	WT1
QG080-01	E04	HQP018504	NM_000552	VWF
QG080-01	E05	HQP018502	NM_000638	VTN
QG080-01	E06	HQP018474	NM_000376	VDR
QG080-01	E07	HQP018402	NM_021833	UCP1
QG080-01	E08	HQP018340	NM_003331	TYK2
QG080-01	E09	HQP018237	NM_004620	TRAF6
QG080-01	E10	HQP018214	NM_000547	TPO
QG080-01	E11	HQP017965	NM_003205	TCF12
QG080-01	E12	HQP017950	NM_003198	TCEB3
QG080-01	F01	HQP017767	NM_003150	STAT3
QG080-01	F02	HQP017766	NM_005419	STAT2
QG080-01	F03	HQP017764	NM_007315	STAT1
QG080-01	F04	HQP017696	NM_005417	SRC
QG080-01	F05	HQP017563	NM_005631	SMO
QG080-01	F06	HQP017163	NM_005067	SIAH2
QG080-01	F07	HQP016743	NM_005065	SEL1L
QG080-01	F08	HQP016694	NM_001035511	SDHC
QG080-01	F09	HQP016689	NM_003000	SDHB
QG080-01	F10	HQP016626	NM_002985	CCL5
QG080-01	F11	HQP016549	NM_002966	S100A10
QG080-01	F12	HQP016212	NM_000657	BCL2
QG080-01	G01	HQP016125	NM_002890	RASA1
QG080-01	G02	HQP016103	NM_001010935	RAP1A
QG080-01	G03	HQP016087	NM_134424	RAD52
QG080-01	G04	HQP015610	NM_002820	PTHLH
QG080-01	G05	HQP015538	NM_004322	BAD
QG080-01	G06	HQP015535	NM_000314	PTEN
QG080-01	G07	HQP015530	NM_000264	PTCH1
QG080-01	G08	HQP014848	NM_002745	MAPK1
QG080-01	G09	HQP014707	NM_018371	ChGn
QG080-01	G10	HQP014650	NM_002734	PRKAR1A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG080-01	G11	HQP013633	NM_005037	PPARG
QG080-01	G12	HQP013352	NM_000535	PMS2
QG080-01	H01	HGDC		
QG080-01	H02	HGDC		
QG080-01	H03	HQP006940	NM_002046	GAPDH
QG080-01	H04	HQP016381	NM_001101	ACTB
QG080-01	H05	HQP015171	NM_004048	B2M
QG080-01	H06	HQP006171	NM_012423	RPL13A
QG080-01	H07	HQP009026	NM_000194	HPRT1
QG080-01	H08	HQP054253	NR_003286	RN18S1
QG080-01	H09	RT		
QG080-01	H10	RT		
QG080-01	H11	PCR		
QG080-01	H12	PCR		