

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG076-01	A01	HQP010133	NM_004985	KRAS
QG076-01	A02	HQP017753	NM_000059	BRCA2
QG076-01	A03	HQP000370	NM_058195	CDKN2A
QG076-01	A04	HQP018175	NM_000546	TP53
QG076-01	A05	HQP011547	NM_005957	MTHFR
QG076-01	A06	HQP018562	NM_006297	XRCC1
QG076-01	A07	HQP017666	NM_003122	SPINK1
QG076-01	A08	HQP004605	NM_005228	EGFR
QG076-01	A09	HQP023467	NM_000662	NAT1
QG076-01	A10	HQP015083	NM_002769	PRSS1
QG076-01	A11	HQP001136	NM_000015	NAT2
QG076-01	A12	HQP012021	NM_002542	OGG1
QG076-01	B01	HQP018141	NM_000594	TNF
QG076-01	B02	HQP011171	NM_000244	MEN1
QG076-01	B03	HQP003772	NM_000499	CYP1A1
QG076-01	B04	HQP020707	NM_003579	RAD54L
QG076-01	B05	HQP018564	NM_005432	XRCC3
QG076-01	B06	HQP018556	NM_004628	XPC
QG076-01	B07	HQP017616	NM_000636	SOD2
QG076-01	B08	HQP013596	NM_001184	ATR
QG076-01	B09	HQP011491	NM_000251	MSH2
QG076-01	B10	HQP011235	NM_000249	MLH1
QG076-01	B11	HQP010961	NM_005359	SMAD4
QG076-01	B12	HQP009670	NM_000600	IL6
QG076-01	C01	HQP009024	NM_000038	APC
QG076-01	C02	HQP008493	NM_000179	MSH6
QG076-01	C03	HQP008362	NM_020469	ABO
QG076-01	C04	HQP004976	NM_000400	ERCC2
QG076-01	C05	HQP001396	NM_001005735	CHEK2
QG076-01	C06	HQP054684	NM_001033	RRM1
QG076-01	C07	HQP053999	NM_004955	SLC29A1
QG076-01	C08	HQP023466	NM_004360	CDH1
QG076-01	C09	HQP023203	NM_001785	CDA
QG076-01	C10	HQP022202	NM_004213	SLC28A1
QG076-01	C11	HQP018563	NM_005431	XRCC2
QG076-01	C12	HQP018342	NM_001071	TYMS
QG076-01	D01	HQP017794	NM_000455	STK11
QG076-01	D02	HQP017733	NM_004333	BRAF
QG076-01	D03	HQP015598	NM_000963	PTGS2
QG076-01	D04	HQP015084	NM_002770	PRSS2
QG076-01	D05	HQP013422	NM_002690	POLB

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG076-01	D06	HQP013150	NM_006218	PIK3CA
QG076-01	D07	HQP011555	NM_002454	MTRR
QG076-01	D08	HQP011320	NM_005590	MRE11A
QG076-01	D09	HQP011135	NM_002392	MDM2
QG076-01	D10	HQP009641	NM_000576	IL1B
QG076-01	D11	HQP009529	NM_000612	IGF2
QG076-01	D12	HQP009518	NM_000618	IGF1
QG076-01	E01	HQP009061	NM_001641	APEX1
QG076-01	E02	HQP005676	NM_016081	PALLD
QG076-01	E03	HQP005083	NM_000136	FANCC
QG076-01	E04	HQP004078	NM_001012732	DCTD
QG076-01	E05	HQP004066	NM_000788	DCK
QG076-01	E06	HQP003775	NM_000104	CYP1B1
QG076-01	E07	HQP001282	NM_001274	CHEK1
QG076-01	E08	HQP018180	NM_005427	TP73
QG076-01	E09	HQP054689	NM_194294	INDOL1
QG076-01	E10	HQP054054	BC039243	FGFR2
QG076-01	E11	HQP054045	BC004257	RET
QG076-01	E12	HQP054030	NM_130398	EXO1
QG076-01	F01	HQP054003	NM_005265	GGT1
QG076-01	F02	HQP023038	NM_014641	MDC1
QG076-01	F03	HQP022197	NM_004212	SLC28A2
QG076-01	F04	HQP021459	NM_003786	ABCC3
QG076-01	F05	HQP021347	NM_001754	RUNX1
QG076-01	F06	HQP020924	NM_181672	OGT
QG076-01	F07	HQP020912	NM_003604	IRS4
QG076-01	F08	HQP019829	NM_030782	CLPTM1L
QG076-01	F09	HQP019268	NM_006304	SHFM1
QG076-01	F10	HQP019129	NM_024596	MCPH1
QG076-01	F11	HQP018966	NM_001080124	CASP8
QG076-01	F12	HQP018546	NM_000378	WT1
QG076-01	G01	HQP018485	NM_000551	VHL
QG076-01	G02	HQP018447	NM_001017415	USP1
QG076-01	G03	HQP018271	NM_000369	TSHR
QG076-01	G04	HQP018176	NM_005657	TP53BP1
QG076-01	G05	HQP018044	NM_000660	TGFB1
QG076-01	G06	HQP018036	NM_005423	TFF2
QG076-01	G07	HQP017954	NM_000545	TCF1
QG076-01	G08	HQP017898	NM_001730	KLF5
QG076-01	G09	HQP017811	NM_177536	SULT1A1
QG076-01	G10	HQP017563	NM_005631	SMO

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG076-01	G11	HQP017526	NM_003073	SMARCB1
QG076-01	G12	HQP016847	NM_022362	MMS19L
QG076-01	H01	HGDC		
QG076-01	H02	HGDC		
QG076-01	H03	HQP006940	NM_002046	GAPDH
QG076-01	H04	HQP016381	NM_001101	ACTB
QG076-01	H05	HQP015171	NM_004048	B2M
QG076-01	H06	HQP006171	NM_012423	RPL13A
QG076-01	H07	HQP009026	NM_000194	HPRT1
QG076-01	H08	HQP054253	NR_003286	RN18S1
QG076-01	H09	RT		
QG076-01	H10	RT		
QG076-01	H11	PCR		
QG076-01	H12	PCR		
QG076-02	A01	HQP016830	NM_003010	MAP2K4
QG076-02	A02	HQP016763	NM_022127	SLC28A3
QG076-02	A03	HQP016744	NM_000450	SELE
QG076-02	A04	HQP016645	NM_001003398	BICD1
QG076-02	A05	HQP016626	NM_002985	CCL5
QG076-02	A06	HQP016303	NM_021133	RNASEL
QG076-02	A07	HQP016155	NM_002894	RBBP8
QG076-02	A08	HQP016131	NM_000321	RB1
QG076-02	A09	HQP015535	NM_000314	PTEN
QG076-02	A10	HQP015530	NM_000264	PTCH1
QG076-02	A11	HQP014292	NM_018315	FBXW7
QG076-02	A12	HQP013633	NM_005037	PPARG
QG076-02	B01	HQP013581	NM_019077	UGT1A7
QG076-02	B02	HQP013312	NM_000534	PMS1
QG076-02	B03	HQP013155	NM_181504	PIK3R1
QG076-02	B04	HQP013100	NM_000927	ABCB1
QG076-02	B05	HQP012504	NM_182944	NIN
QG076-02	B06	HQP011946	NM_001007156	NTRK3
QG076-02	B07	HQP011914	NM_002524	NRAS
QG076-02	B08	HQP011807	NM_003998	NFKB1
QG076-02	B09	HQP011778	NM_000268	NF2
QG076-02	B10	HQP011774	NM_000267	NF1
QG076-02	B11	HQP011687	NM_002485	NBN
QG076-02	B12	HQP011554	NM_000254	MTR
QG076-02	C01	HQP011492	NM_002439	MSH3
QG076-02	C02	HQP011325	NM_019899	ABCC1
QG076-02	C03	HQP011309	NM_000250	MPO

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG076-02	C04	HQP011308	NM_005373	MPL
QG076-02	C05	HQP011181	NM_000245	MET
QG076-02	C06	HQP010960	NM_005902	SMAD3
QG076-02	C07	HQP010959	NM_001003652	SMAD2
QG076-02	C08	HQP010613	NM_002312	LIG4
QG076-02	C09	HQP010612	NM_013975	LIG3
QG076-02	C10	HQP010099	NM_000222	KIT
QG076-02	C11	HQP009850	NM_004972	JAK2
QG076-02	C12	HQP009788	NM_005544	IRS1
QG076-02	D01	HQP009671	NM_000639	FASLG
QG076-02	D02	HQP009664	NM_000418	IL4R
QG076-02	D03	HQP009662	NM_000589	IL4
QG076-02	D04	HQP009551	NM_000599	IGFBP5
QG076-02	D05	HQP009544	NM_000598	IGFBP3
QG076-02	D06	HQP009539	NM_000596	IGFBP1
QG076-02	D07	HQP009523	NM_000875	IGF1R
QG076-02	D08	HQP009351	NM_005896	IDH1
QG076-02	D09	HQP009101	NM_001168	BIRC5
QG076-02	D10	HQP009036	NM_005343	HRAS
QG076-02	D11	HQP008849	NM_002116	HLA-A
QG076-02	D12	HQP008479	NM_001512	GSTA4
QG076-02	E01	HQP007814	NM_000175	GPI
QG076-02	E02	HQP007755	NM_000516	GNAS
QG076-02	E03	HQP007366	NM_002056	GFPT1
QG076-02	E04	HQP007239	NM_000162	GCK
QG076-02	E05	HQP006715	NM_012415	RAD54B
QG076-02	E06	HQP006438	NM_205860	NR5A2
QG076-02	E07	HQP005890	NM_004119	FLT3
QG076-02	E08	HQP005434	NM_000142	FGFR3
QG076-02	E09	HQP005098	NM_022725	FANCF
QG076-02	E10	HQP005087	NM_000690	ALDH2
QG076-02	E11	HQP005084	NM_001018115	FANCD2
QG076-02	E12	HQP005081	NM_000135	FANCA
QG076-02	F01	HQP004984	NM_005236	ERCC4
QG076-02	F02	HQP004924	NM_005233	EPHA3
QG076-02	F03	HQP004599	NM_001963	EGF
QG076-02	F04	HQP004439	NM_000110	DPYD
QG076-02	F05	HQP004435	NM_001039350	DPP6
QG076-02	F06	HQP003888	NM_000102	CYP17A1
QG076-02	F07	HQP003817	NM_000773	CYP2E1
QG076-02	F08	HQP003804	NM_000766	CYP2A13

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG076-02	F09	HQP003539	NM_001904	CTNNB1
QG076-02	F10	HQP003158	NM_005211	CSF1R
QG076-02	F11	HQP002413	NM_000669	ADH1C
QG076-02	F12	HQP002331	NM_000668	ADH1B
QG076-02	G01	HQP002210	NM_000579	CCR5
QG076-02	G02	HQP001507	NM_007249	KLF12
QG076-02	G03	HQP001504	NM_032166	TREX1
QG076-02	G04	HQP001285	NM_007050	PTPRT
QG076-02	G05	HQP001228	NM_007027	TOPBP1
QG076-02	G06	HQP000659	NM_006409	ARPC1A
QG076-02	G07	HQP000615	NM_004364	CEBPA
QG076-02	G08	HQP000245	NM_000075	CDK4
QG076-02	G09	HQP000145	NM_005732	RAD50
QG076-02	G10	HQP000075	NM_005688	ABCC5
QG076-02	G11	HQP013597	NM_021027	UGT1A9
QG076-02	G12	HQP008556	NM_001522	GUCY2F
QG076-02	H01	HGDC		
QG076-02	H02	HGDC		
QG076-02	H03	HQP006940	NM_002046	GAPDH
QG076-02	H04	HQP016381	NM_001101	ACTB
QG076-02	H05	HQP015171	NM_004048	B2M
QG076-02	H06	HQP006171	NM_012423	RPL13A
QG076-02	H07	HQP009026	NM_000194	HPRT1
QG076-02	H08	HQP054253	NR_003286	RN18S1
QG076-02	H09	RT		
QG076-02	H10	RT		
QG076-02	H11	PCR		
QG076-02	H12	PCR		