

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG061-01	A01	HQP004605	NM_005228	EGFR
QG061-01	A02	HQP010133	NM_004985	KRAS
QG061-01	A03	HQP018175	NM_000546	TP53
QG061-01	A04	HQP011547	NM_005957	MTHFR
QG061-01	A05	HQP009024	NM_000038	APC
QG061-01	A06	HQP017733	NM_004333	BRAF
QG061-01	A07	HQP018562	NM_006297	XRCC1
QG061-01	A08	HQP004976	NM_000400	ERCC2
QG061-01	A09	HQP009641	NM_000576	IL1B
QG061-01	A10	HQP015598	NM_000963	PTGS2
QG061-01	A11	HQP003772	NM_000499	CYP1A1
QG061-01	A12	HQP018342	NM_001071	TYMS
QG061-01	B01	HQP012021	NM_002542	OGG1
QG061-01	B02	HQP018474	NM_000376	VDR
QG061-01	B03	HQP009645	NM_000577	IL1RN
QG061-01	B04	HQP009685	NM_000572	IL10
QG061-01	B05	HQP001136	NM_000015	NAT2
QG061-01	B06	HQP018564	NM_005432	XRCC3
QG061-01	B07	HQP011491	NM_000251	MSH2
QG061-01	B08	HQP011235	NM_000249	MLH1
QG061-01	B09	HQP009678	NM_000584	IL8
QG061-01	B10	HQP018141	NM_000594	TNF
QG061-01	B11	HQP018044	NM_000660	TGFB1
QG061-01	B12	HQP017753	NM_000059	BRCA2
QG061-01	C01	HQP013633	NM_005037	PPARG
QG061-01	C02	HQP013150	NM_006218	PIK3CA
QG061-01	C03	HQP011554	NM_000254	MTR
QG061-01	C04	HQP009670	NM_000600	IL6
QG061-01	C05	HQP009518	NM_000618	IGF1
QG061-01	C06	HQP004975	NM_202001	ERCC1
QG061-01	C07	HQP004317	NM_000903	NQO1
QG061-01	C08	HQP018556	NM_004628	XPC
QG061-01	C09	HQP018475	NM_001025366	VEGFA
QG061-01	C10	HQP015083	NM_002769	PRSS1
QG061-01	C11	HQP013100	NM_000927	ABCB1
QG061-01	C12	HQP010961	NM_005359	SMAD4
QG061-01	D01	HQP009544	NM_000598	IGFBP3
QG061-01	D02	HQP009523	NM_000875	IGF1R
QG061-01	D03	HQP009036	NM_005343	HRAS
QG061-01	D04	HQP004599	NM_001963	EGF
QG061-01	D05	HQP003817	NM_000773	CYP2E1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG061-01	D06	HQP000370	NM_058195	CDKN2A
QG061-01	D07	HQP023467	NM_000662	NAT1
QG061-01	D08	HQP021961	NM_003977	AIP
QG061-01	D09	HQP018176	NM_005657	TP53BP1
QG061-01	D10	HQP011135	NM_002392	MDM2
QG061-01	D11	HQP009671	NM_000639	FASLG
QG061-01	D12	HQP009662	NM_000589	IL4
QG061-01	E01	HQP009529	NM_000612	IGF2
QG061-01	E02	HQP009061	NM_001641	APEX1
QG061-01	E03	HQP008757	NM_000410	HFE
QG061-01	E04	HQP008493	NM_000179	MSH6
QG061-01	E05	HQP008402	NM_001020825	NR3C1
QG061-01	E06	HQP004948	NM_000120	EPHX1
QG061-01	E07	HQP003904	NM_000103	CYP19A1
QG061-01	E08	HQP003814	NM_000106	CYP2D6
QG061-01	E09	HQP001643	NM_000745	CHRNA5
QG061-01	E10	HQP020481	NM_033338	CASP7
QG061-01	E11	HQP020427	NM_001226	CASP6
QG061-01	E12	HQP020297	NM_004346	CASP3
QG061-01	F01	HQP018563	NM_005431	XRCC2
QG061-01	F02	HQP017794	NM_000455	STK11
QG061-01	F03	HQP016204	NM_053056	CCND1
QG061-01	F04	HQP015596	NM_000962	PTGS1
QG061-01	F05	HQP015535	NM_000314	PTEN
QG061-01	F06	HQP015084	NM_002770	PRSS2
QG061-01	F07	HQP012013	NM_002539	ODC1
QG061-01	F08	HQP011914	NM_002524	NRAS
QG061-01	F09	HQP011866	NM_000625	NOS2A
QG061-01	F10	HQP011492	NM_002439	MSH3
QG061-01	F11	HQP010584	NM_002303	LEPR
QG061-01	F12	HQP009801	NM_000044	AR
QG061-01	G01	HQP009664	NM_000418	IL4R
QG061-01	G02	HQP009556	NM_000041	APOE
QG061-01	G03	HQP007767	NM_002075	GNB3
QG061-01	G04	HQP007755	NM_000516	GNAS
QG061-01	G05	HQP007385	NM_000515	GH1
QG061-01	G06	HQP005087	NM_000690	ALDH2
QG061-01	G07	HQP004991	NM_001014431	AKT1
QG061-01	G08	HQP004446	NM_000795	DRD2
QG061-01	G09	HQP003888	NM_000102	CYP17A1
QG061-01	G10	HQP003811	NM_000771	CYP2C9

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG061-01	G11	HQP003775	NM_000104	CYP1B1
QG061-01	G12	HQP002413	NM_000669	ADH1C
QG061-01	H01	HGDC		
QG061-01	H02	HGDC		
QG061-01	H03	HQP006940	NM_002046	GAPDH
QG061-01	H04	HQP016381	NM_001101	ACTB
QG061-01	H05	HQP015171	NM_004048	B2M
QG061-01	H06	HQP006171	NM_012423	RPL13A
QG061-01	H07	HQP009026	NM_000194	HPRT1
QG061-01	H08	HQP054253	NR_003286	RN18S1
QG061-01	H09	RT		
QG061-01	H10	RT		
QG061-01	H11	PCR		
QG061-01	H12	PCR		
QG061-02	A01	HQP001396	NM_001005735	CHEK2
QG061-02	A02	HQP018180	NM_005427	TP73
QG061-02	A03	HQP011544	NM_002452	NUDT1
QG061-02	A04	HQP004414	NM_006892	DNMT3B
QG061-02	A05	HQP054684	NM_001033	RRM1
QG061-02	A06	HQP054057	BC071181	TGFBR1
QG061-02	A07	HQP054047	BC008403	HLA-DRB1
QG061-02	A08	HQP054045	BC004257	RET
QG061-02	A09	HQP054030	NM_130398	EXO1
QG061-02	A10	HQP053916	NM_001076	UGT2B15
QG061-02	A11	HQP023466	NM_004360	CDH1
QG061-02	A12	HQP023414	NM_005847	SLC23A1
QG061-02	B01	HQP023203	NM_001785	CDA
QG061-02	B02	HQP023038	NM_014641	MDC1
QG061-02	B03	HQP022500	NM_001040280	CD83
QG061-02	B04	HQP022490	NM_000591	CD14
QG061-02	B05	HQP021597	NM_003873	NRP1
QG061-02	B06	HQP021518	NM_000071	CBS
QG061-02	B07	HQP021459	NM_003786	ABCC3
QG061-02	B08	HQP021351	NM_001029851	PDE8B
QG061-02	B09	HQP020912	NM_003604	IRS4
QG061-02	B10	HQP020384	NM_004347	CASP5
QG061-02	B11	HQP020332	NM_001225	CASP4
QG061-02	B12	HQP020207	NM_001223	CASP1
QG061-02	C01	HQP020116	NM_004655	AXIN2
QG061-02	C02	HQP019829	NM_030782	CLPTM1L
QG061-02	C03	HQP019268	NM_006304	SHFM1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG061-02	C04	HQP019141	NM_024608	NEIL1
QG061-02	C05	HQP019129	NM_024596	MCPH1
QG061-02	C06	HQP018983	NM_004639	BAT3
QG061-02	C07	HQP018966	NM_001080124	CASP8
QG061-02	C08	HQP018568	NM_021141	XRCC5
QG061-02	C09	HQP018565	NM_003401	XRCC4
QG061-02	C10	HQP018447	NM_001017415	USP1
QG061-02	C11	HQP018425	NM_000373	UMPS
QG061-02	C12	HQP018418	NM_001074	UGT2B7
QG061-02	D01	HQP018337	NM_182729	TXNRD1
QG061-02	D02	HQP017979	NM_000355	TCN2
QG061-02	D03	HQP017616	NM_000636	SOD2
QG061-02	D04	HQP017482	NM_194255	SLC19A1
QG061-02	D05	HQP017462	NM_000452	SLC10A2
QG061-02	D06	HQP016847	NM_022362	MMS19L
QG061-02	D07	HQP016819	NM_005410	SEPP1
QG061-02	D08	HQP016801	NM_022162	NOD2
QG061-02	D09	HQP016744	NM_000450	SELE
QG061-02	D10	HQP016526	NM_002957	RXRA
QG061-02	D11	HQP016155	NM_002894	RBBP8
QG061-02	D12	HQP016125	NM_002890	RASA1
QG061-02	E01	HQP015563	NM_000958	PTGER4
QG061-02	E02	HQP015544	NM_000956	PTGER2
QG061-02	E03	HQP015530	NM_000264	PTCH1
QG061-02	E04	HQP014650	NM_002734	PRKAR1A
QG061-02	E05	HQP014243	NM_018272	CASC1
QG061-02	E06	HQP014222	NM_018248	NEIL3
QG061-02	E07	HQP013734	NM_017672	TRPM7
QG061-02	E08	HQP013616	NM_019093	UGT1A3
QG061-02	E09	HQP013614	NM_007120	UGT1A4
QG061-02	E10	HQP013596	NM_001184	ATR
QG061-02	E11	HQP013583	NM_205862	UGT1A6
QG061-02	E12	HQP013579	NM_019075	UGT1A10
QG061-02	F01	HQP013388	NM_017442	TLR9
QG061-02	F02	HQP013312	NM_000534	PMS1
QG061-02	F03	HQP012979	NM_002613	PDPK1
QG061-02	F04	HQP012501	NM_016341	PLCE1
QG061-02	F05	HQP011810	NM_020529	NFKBIA
QG061-02	F06	HQP011807	NM_003998	NFKB1
QG061-02	F07	HQP011800	NM_006164	NFE2L2
QG061-02	F08	HQP011687	NM_002485	NBN

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG061-02	F09	HQP011555	NM_002454	MTRR
QG061-02	F10	HQP011325	NM_019899	ABCC1
QG061-02	F11	HQP011320	NM_005590	MRE11A
QG061-02	F12	HQP011309	NM_000250	MPO
QG061-02	G01	HQP011266	NM_002426	MMP12
QG061-02	G02	HQP011257	NM_002422	MMP3
QG061-02	G03	HQP011256	NM_004530	MMP2
QG061-02	G04	HQP011255	NM_002421	MMP1
QG061-02	G05	HQP011171	NM_000244	MEN1
QG061-02	G06	HQP010868	NM_006152	LRMP
QG061-02	G07	HQP010613	NM_002312	LIG4
QG061-02	G08	HQP009788	NM_005544	IRS1
QG061-02	G09	HQP009718	NM_001562	IL18
QG061-02	G10	HQP009693	NM_002187	IL12B
QG061-02	G11	HQP009692	NM_000882	IL12A
QG061-02	G12	HQP009640	NM_000575	IL1A
QG061-02	H01	HGDC		
QG061-02	H02	HGDC		
QG061-02	H03	HQP006940	NM_002046	GAPDH
QG061-02	H04	HQP016381	NM_001101	ACTB
QG061-02	H05	HQP015171	NM_004048	B2M
QG061-02	H06	HQP006171	NM_012423	RPL13A
QG061-02	H07	HQP009026	NM_000194	HPRT1
QG061-02	H08	HQP054253	NR_003286	RN18S1
QG061-02	H09	RT		
QG061-02	H10	RT		
QG061-02	H11	PCR		
QG061-02	H12	PCR		