

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-01	A01	HQP010133	NM_004985	KRAS
QG065-01	A02	HQP011547	NM_005957	MTHFR
QG065-01	A03	HQP011235	NM_000249	MLH1
QG065-01	A04	HQP011491	NM_000251	MSH2
QG065-01	A05	HQP018175	NM_000546	TP53
QG065-01	A06	HQP017733	NM_004333	BRAF
QG065-01	A07	HQP009024	NM_000038	APC
QG065-01	A08	HQP018342	NM_001071	TYMS
QG065-01	A09	HQP008493	NM_000179	MSH6
QG065-01	A10	HQP001136	NM_000015	NAT2
QG065-01	A11	HQP018562	NM_006297	XRCC1
QG065-01	A12	HQP018474	NM_000376	VDR
QG065-01	B01	HQP003772	NM_000499	CYP1A1
QG065-01	B02	HQP004976	NM_000400	ERCC2
QG065-01	B03	HQP023467	NM_000662	NAT1
QG065-01	B04	HQP016204	NM_053056	CCND1
QG065-01	B05	HQP015598	NM_000963	PTGS2
QG065-01	B06	HQP013150	NM_006218	PIK3CA
QG065-01	B07	HQP011554	NM_000254	MTR
QG065-01	B08	HQP013100	NM_000927	ABCB1
QG065-01	B09	HQP004975	NM_202001	ERCC1
QG065-01	B10	HQP003817	NM_000773	CYP2E1
QG065-01	B11	HQP023466	NM_004360	CDH1
QG065-01	B12	HQP004948	NM_000120	EPHX1
QG065-01	C01	HQP004317	NM_000903	NQO1
QG065-01	C02	HQP009670	NM_000600	IL6
QG065-01	C03	HQP003811	NM_000771	CYP2C9
QG065-01	C04	HQP012021	NM_002542	OGG1
QG065-01	C05	HQP018475	NM_001025366	VEGFA
QG065-01	C06	HQP009518	NM_000618	IGF1
QG065-01	C07	HQP004605	NM_005228	EGFR
QG065-01	C08	HQP018044	NM_000660	TGFB1
QG065-01	C09	HQP013633	NM_005037	PPARG
QG065-01	C10	HQP011555	NM_002454	MTRR
QG065-01	C11	HQP010966	NM_005904	SMAD7
QG065-01	C12	HQP018141	NM_000594	TNF
QG065-01	D01	HQP017811	NM_177536	SULT1A1
QG065-01	D02	HQP005087	NM_000690	ALDH2
QG065-01	D03	HQP004439	NM_000110	DPYD
QG065-01	D04	HQP003775	NM_000104	CYP1B1
QG065-01	D05	HQP001396	NM_001005735	CHEK2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-01	D06	HQP054057	BC071181	TGFBR1
QG065-01	D07	HQP018564	NM_005432	XRCC3
QG065-01	D08	HQP016801	NM_022162	NOD2
QG065-01	D09	HQP011255	NM_002421	MMP1
QG065-01	D10	HQP009556	NM_000041	APOE
QG065-01	D11	HQP013581	NM_019077	UGT1A7
QG065-01	D12	HQP013352	NM_000535	PMS2
QG065-01	E01	HQP009544	NM_000598	IGFBP3
QG065-01	E02	HQP011135	NM_002392	MDM2
QG065-01	E03	HQP009685	NM_000572	IL10
QG065-01	E04	HQP008757	NM_000410	HFE
QG065-01	E05	HQP021215	NM_033103	RHPN2
QG065-01	E06	HQP013583	NM_205862	UGT1A6
QG065-01	E07	HQP009788	NM_005544	IRS1
QG065-01	E08	HQP009678	NM_000584	IL8
QG065-01	E09	HQP009641	NM_000576	IL1B
QG065-01	E10	HQP008483	NM_000849	GSTM3
QG065-01	E11	HQP003814	NM_000106	CYP2D6
QG065-01	E12	HQP002413	NM_000669	ADH1C
QG065-01	F01	HQP016418	NM_004327	BCR
QG065-01	F02	HQP015596	NM_000962	PTGS1
QG065-01	F03	HQP011807	NM_003998	NFKB1
QG065-01	F04	HQP011263	NM_004994	MMP9
QG065-01	F05	HQP011257	NM_002422	MMP3
QG065-01	F06	HQP004599	NM_001963	EGF
QG065-01	F07	HQP003539	NM_001904	CTNNB1
QG065-01	F08	HQP000370	NM_058195	CDKN2A
QG065-01	F09	HQP004985	NM_000123	ERCC5
QG065-01	F10	HQP020899	NM_000388	CASR
QG065-01	F11	HQP011256	NM_004530	MMP2
QG065-01	F12	HQP004998	NM_000125	ESR1
QG065-01	G01	HQP002331	NM_000668	ADH1B
QG065-01	G02	HQP018180	NM_005427	TP73
QG065-01	G03	HQP013597	NM_021027	UGT1A9
QG065-01	G04	HQP018556	NM_004628	XPC
QG065-01	G05	HQP017753	NM_000059	BRCA2
QG065-01	G06	HQP009662	NM_000589	IL4
QG065-01	G07	HQP009645	NM_000577	IL1RN
QG065-01	G08	HQP009061	NM_001641	APEX1
QG065-01	G09	HQP009036	NM_005343	HRAS
QG065-01	G10	HQP008832	NM_181054	HIF1A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-01	G11	HQP007385	NM_000515	GH1
QG065-01	G12	HQP005280	NM_021642	FCGR2A
QG065-01	H01	HGDC		
QG065-01	H02	HGDC		
QG065-01	H03	HQP006940	NM_002046	GAPDH
QG065-01	H04	HQP016381	NM_001101	ACTB
QG065-01	H05	HQP015171	NM_004048	B2M
QG065-01	H06	HQP006171	NM_012423	RPL13A
QG065-01	H07	HQP009026	NM_000194	HPRT1
QG065-01	H08	HQP054253	NR_003286	RN18S1
QG065-01	H09	RT		
QG065-01	H10	RT		
QG065-01	H11	PCR		
QG065-01	H12	PCR		
QG065-02	A01	HQP003499	NM_001037631	CTLA4
QG065-02	A02	HQP003120	NM_001618	PARP1
QG065-02	A03	HQP011544	NM_002452	NUDT1
QG065-02	A04	HQP004414	NM_006892	DNMT3B
QG065-02	A05	HQP054030	NM_130398	EXO1
QG065-02	A06	HQP053910	NM_130851	BMP4
QG065-02	A07	HQP022625	NM_004797	ADIPOQ
QG065-02	A08	HQP021518	NM_000071	CBS
QG065-02	A09	HQP021417	NM_003756	EIF3S3
QG065-02	A10	HQP018966	NM_001080124	CASP8
QG065-02	A11	HQP017616	NM_000636	SOD2
QG065-02	A12	HQP015535	NM_000314	PTEN
QG065-02	B01	HQP012013	NM_002539	ODC1
QG065-02	B02	HQP011868	NM_000603	NOS3
QG065-02	B03	HQP011492	NM_002439	MSH3
QG065-02	B04	HQP010584	NM_002303	LEPR
QG065-02	B05	HQP010581	NM_000230	LEP
QG065-02	B06	HQP010570	NM_002299	LCT
QG065-02	B07	HQP006356	NM_000697	ALOX12
QG065-02	B08	HQP005002	NM_001437	ESR2
QG065-02	B09	HQP004081	NM_000789	ACE
QG065-02	B10	HQP003904	NM_000103	CYP19A1
QG065-02	B11	HQP003888	NM_000102	CYP17A1
QG065-02	B12	HQP003841	NM_000777	CYP3A5
QG065-02	C01	HQP003809	NM_000769	CYP2C19
QG065-02	C02	HQP002671	NM_000754	COMT
QG065-02	C03	HQP021586	NM_198061	CES2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-02	C04	HQP020114	NM_003502	AXIN1
QG065-02	C05	HQP018565	NM_003401	XRCC4
QG065-02	C06	HQP017979	NM_000355	TCN2
QG065-02	C07	HQP017482	NM_194255	SLC19A1
QG065-02	C08	HQP017462	NM_000452	SLC10A2
QG065-02	C09	HQP016819	NM_005410	SEPP1
QG065-02	C10	HQP016670	NM_001033886	CXCL12
QG065-02	C11	HQP015160	NM_020415	RETN
QG065-02	C12	HQP013204	NM_002658	PLAU
QG065-02	D01	HQP012154	NM_000602	SERPINE1
QG065-02	D02	HQP011914	NM_002524	NRAS
QG065-02	D03	HQP011866	NM_000625	NOS2A
QG065-02	D04	HQP011687	NM_002485	NBN
QG065-02	D05	HQP010961	NM_005359	SMAD4
QG065-02	D06	HQP009801	NM_000044	AR
QG065-02	D07	HQP009664	NM_000418	IL4R
QG065-02	D08	HQP009523	NM_000875	IGF1R
QG065-02	D09	HQP006359	NM_000698	ALOX5
QG065-02	D10	HQP006129	NM_001014986	FOLH1
QG065-02	D11	HQP005058	NM_000130	F5
QG065-02	D12	HQP004991	NM_001014431	AKT1
QG065-02	E01	HQP004986	NM_000124	ERCC6
QG065-02	E02	HQP004984	NM_005236	ERCC4
QG065-02	E03	HQP004658	NM_001621	AHR
QG065-02	E04	HQP004446	NM_000795	DRD2
QG065-02	E05	HQP022821	NM_000072	CD36
QG065-02	E06	HQP022490	NM_000591	CD14
QG065-02	E07	HQP021906	NM_001239	CCNH
QG065-02	E08	HQP020116	NM_004655	AXIN2
QG065-02	E09	HQP018568	NM_021141	XRCC5
QG065-02	E10	HQP018563	NM_005431	XRCC2
QG065-02	E11	HQP018425	NM_000373	UMPS
QG065-02	E12	HQP016874	NM_003013	SFRP2
QG065-02	F01	HQP016844	NM_000057	BLM
QG065-02	F02	HQP016744	NM_000450	SELE
QG065-02	F03	HQP016526	NM_002957	RXRA
QG065-02	F04	HQP016076	NM_002874	RAD23B
QG065-02	F05	HQP015925	NM_002843	PTPRJ
QG065-02	F06	HQP015897	NM_006264	PTPN13
QG065-02	F07	HQP013627	NM_006238	PPARD
QG065-02	F08	HQP013528	NM_006235	POU2AF1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-02	F09	HQP013422	NM_002690	POLB
QG065-02	F10	HQP013312	NM_000534	PMS1
QG065-02	F11	HQP013194	NM_000300	PLA2G2A
QG065-02	F12	HQP013099	NM_000926	PGR
QG065-02	G01	HQP012996	NM_016362	GHRL
QG065-02	G02	HQP011800	NM_006164	NFE2L2
QG065-02	G03	HQP011597	NM_002467	MYC
QG065-02	G04	HQP011325	NM_019899	ABCC1
QG065-02	G05	HQP011097	NM_005912	MC4R
QG065-02	G06	HQP009818	NM_000212	ITGB3
QG065-02	G07	HQP009764	NM_000208	INSR
QG065-02	G08	HQP009749	NM_000207	INS
QG065-02	G09	HQP009671	NM_000639	FASLG
QG065-02	G10	HQP009640	NM_000575	IL1A
QG065-02	G11	HQP009066	NM_002153	HSD17B2
QG065-02	G12	HQP009022	NM_000860	HPGD
QG065-02	H01	HGDC		
QG065-02	H02	HGDC		
QG065-02	H03	HQP006940	NM_002046	GAPDH
QG065-02	H04	HQP016381	NM_001101	ACTB
QG065-02	H05	HQP015171	NM_004048	B2M
QG065-02	H06	HQP006171	NM_012423	RPL13A
QG065-02	H07	HQP009026	NM_000194	HPRT1
QG065-02	H08	HQP054253	NR_003286	RN18S1
QG065-02	H09	RT		
QG065-02	H10	RT		
QG065-02	H11	PCR		
QG065-02	H12	PCR		
QG065-03	A01	HQP008285	NM_002085	GPX4
QG065-03	A02	HQP006431	NM_001463	FRZB
QG065-03	A03	HQP006426	NM_004958	FRAP1
QG065-03	A04	HQP005959	NM_001002294	FMO3
QG065-03	A05	HQP004940	NM_017449	EPHB2
QG065-03	A06	HQP004309	NM_000791	DHFR
QG065-03	A07	HQP004053	NM_005215	DCC
QG065-03	A08	HQP003923	NM_000785	CYP27B1
QG065-03	A09	HQP003917	NM_000782	CYP24A1
QG065-03	A10	HQP003810	NM_000770	CYP2C8
QG065-03	A11	HQP001282	NM_001274	CHEK1
QG065-03	A12	HQP000970	NM_006663	PPP1R13L
QG065-03	B01	HQP000285	NM_001799	CDK7

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-03	B02	HQP000132	NM_000022	ADA
QG065-03	B03	HQP011601	NM_005376	MYCL1
QG065-03	B04	HQP003859	NM_000780	CYP7A1
QG065-03	B05	HQP054043	NM_203472	SELS
QG065-03	B06	HQP053962	NM_001045	SLC6A4
QG065-03	B07	HQP023464	NM_012115	CASP8AP2
QG065-03	B08	HQP023414	NM_005847	SLC23A1
QG065-03	B09	HQP023203	NM_001785	CDA
QG065-03	B10	HQP023134	NM_014707	HDAC9
QG065-03	B11	HQP022853	NM_004864	GDF15
QG065-03	B12	HQP022804	NM_004849	ATG5
QG065-03	C01	HQP022747	NM_175862	CD86
QG065-03	C02	HQP022699	NM_006139	CD28
QG065-03	C03	HQP022491	NM_001616	ACVR2A
QG065-03	C04	HQP022361	NM_004747	DLG5
QG065-03	C05	HQP022168	NM_004707	ATG12
QG065-03	C06	HQP021752	NM_003939	BTRC
QG065-03	C07	HQP021683	NM_021090	MTMR3
QG065-03	C08	HQP021630	NM_003889	NR1I2
QG065-03	C09	HQP021610	NM_003882	WISP1
QG065-03	C10	HQP021603	NM_003878	GGH
QG065-03	C11	HQP021557	NM_003844	TNFRSF10A
QG065-03	C12	HQP021553	NM_003842	TNFRSF10B
QG065-03	D01	HQP021492	NM_003804	RIPK1
QG065-03	D02	HQP021459	NM_003786	ABCC3
QG065-03	D03	HQP021313	NM_001753	CAV1
QG065-03	D04	HQP021276	NM_003672	CDC14A
QG065-03	D05	HQP021250	NM_033389	SSH2
QG065-03	D06	HQP020982	NM_001039970	PTPN5
QG065-03	D07	HQP020596	NM_001042544	LTBP4
QG065-03	D08	HQP020481	NM_033338	CASP7
QG065-03	D09	HQP020427	NM_001226	CASP6
QG065-03	D10	HQP020297	NM_004346	CASP3
QG065-03	D11	HQP019829	NM_030782	CLPTM1L
QG065-03	D12	HQP019462	NM_024945	RMI1
QG065-03	E01	HQP019284	NM_024757	EHMT1
QG065-03	E02	HQP019172	NM_024642	GALNT12
QG065-03	E03	HQP019051	NM_007088	CALB2
QG065-03	E04	HQP018802	NM_001008540	CXCR4
QG065-03	E05	HQP018755	NM_001007257	PRDM2
QG065-03	E06	HQP018659	NM_001215	CA6

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-03	E07	HQP018545	NM_000553	WRN
QG065-03	E08	HQP018540	NM_004185	WNT2B
QG065-03	E09	HQP018485	NM_000551	VHL
QG065-03	E10	HQP018418	NM_001074	UGT2B7
QG065-03	E11	HQP018337	NM_182729	TXNRD1
QG065-03	E12	HQP018288	NM_003318	TTK
QG065-03	F01	HQP018265	NM_000548	TSC2
QG065-03	F02	HQP018263	NM_000368	TSC1
QG065-03	F03	HQP018176	NM_005657	TP53BP1
QG065-03	F04	HQP018104	NM_005077	TLE1
QG065-03	F05	HQP017958	NM_003202	TCF7
QG065-03	F06	HQP017954	NM_000545	TCF1
QG065-03	F07	HQP017535	NM_001204	BMPR2
QG065-03	F08	HQP017496	NM_003060	SLC22A5
QG065-03	F09	HQP017495	NM_003059	SLC22A4
QG065-03	F10	HQP016830	NM_003010	MAP2K4
QG065-03	F11	HQP016746	NM_003005	SELP
QG065-03	F12	HQP016652	NM_002996	CX3CL1
QG065-03	G01	HQP016634	NM_001713	BHMT
QG065-03	G02	HQP016621	NM_002982	CCL2
QG065-03	G03	HQP016620	NM_002981	CCL1
QG065-03	G04	HQP016471	NM_003161	RPS6KB1
QG065-03	G05	HQP016468	NM_001006932	RPS6KA2
QG065-03	G06	HQP016466	NM_001006665	RPS6KA1
QG065-03	G07	HQP016303	NM_021133	RNASEL
QG065-03	G08	HQP016224	NM_002913	RFC1
QG065-03	G09	HQP016168	NM_021629	GNB4
QG065-03	G10	HQP016131	NM_000321	RB1
QG065-03	G11	HQP016087	NM_134424	RAD52
QG065-03	G12	HQP015946	NM_000465	BARD1
QG065-03	H01	HGDC		
QG065-03	H02	HGDC		
QG065-03	H03	HQP006940	NM_002046	GAPDH
QG065-03	H04	HQP016381	NM_001101	ACTB
QG065-03	H05	HQP015171	NM_004048	B2M
QG065-03	H06	HQP006171	NM_012423	RPL13A
QG065-03	H07	HQP009026	NM_000194	HPRT1
QG065-03	H08	HQP054253	NR_003286	RN18S1
QG065-03	H09	RT		
QG065-03	H10	RT		
QG065-03	H11	PCR		



Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-03	H12	PCR		
QG065-04	A01	HQP015939	NM_002850	PTPRS
QG065-04	A02	HQP015563	NM_000958	PTGER4
QG065-04	A03	HQP015544	NM_000956	PTGER2
QG065-04	A04	HQP015530	NM_000264	PTCH1
QG065-04	A05	HQP015311	NM_002800	PSMB9
QG065-04	A06	HQP015190	NM_020165	RAD18
QG065-04	A07	HQP015087	NM_019845	RPRM
QG065-04	A08	HQP015032	NM_000311	PRNP
QG065-04	A09	HQP014691	NM_001008744	TDP1
QG065-04	A10	HQP014530	NM_006251	PRKAA1
QG065-04	A11	HQP014500	NM_017633	FAM46A
QG065-04	A12	HQP014362	NM_018407	LAPTM4B
QG065-04	B01	HQP014052	NM_018036	C14orf103
QG065-04	B02	HQP014024	NM_002711	PPP1R3A
QG065-04	B03	HQP013734	NM_017672	TRPM7
QG065-04	B04	HQP013596	NM_001184	ATR
QG065-04	B05	HQP013473	NM_000446	PON1
QG065-04	B06	HQP013467	NM_000939	POMC
QG065-04	B07	HQP013463	NM_018984	SSH1
QG065-04	B08	HQP013155	NM_181504	PIK3R1
QG065-04	B09	HQP013122	NM_000295	SERPINA1
QG065-04	B10	HQP013026	NM_002615	SERPINF1
QG065-04	B11	HQP012993	NM_016327	UPB1
QG065-04	B12	HQP012972	NM_016231	NLK
QG065-04	C01	HQP012723	NM_016203	PRKAG2
QG065-04	C02	HQP012675	NM_015891	CDC40
QG065-04	C03	HQP012420	NM_002592	PCNA
QG065-04	C04	HQP012393	NM_016009	SH3GLB1
QG065-04	C05	HQP012386	NM_015999	ADIPOR1
QG065-04	C06	HQP012161	NM_020525	IL22
QG065-04	C07	HQP012131	NM_015713	RRM2B
QG065-04	C08	HQP012121	NM_002569	FURIN
QG065-04	C09	HQP011932	NM_002526	NT5E
QG065-04	C10	HQP011920	NM_005173	ATP2A3
QG065-04	C11	HQP011909	NM_001681	ATP2A2
QG065-04	C12	HQP011853	NM_000269	NME1
QG065-04	D01	HQP011810	NM_020529	NFKBIA
QG065-04	D02	HQP011727	NM_004044	ATIC
QG065-04	D03	HQP011549	NM_000252	MTM1
QG065-04	D04	HQP011309	NM_000250	MPO



Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-04	D05	HQP011266	NM_002426	MMP12
QG065-04	D06	HQP011211	NM_000246	CIITA
QG065-04	D07	HQP011208	NM_020300	MGST1
QG065-04	D08	HQP011047	NM_005911	MAT2A
QG065-04	D09	HQP011006	NM_005908	MANBA
QG065-04	D10	HQP010960	NM_005902	SMAD3
QG065-04	D11	HQP010959	NM_001003652	SMAD2
QG065-04	D12	HQP010924	NM_001668	ARNT
QG065-04	E01	HQP010912	NM_206943	LTBP1
QG065-04	E02	HQP010626	NM_000236	LIPC
QG065-04	E03	HQP010613	NM_002312	LIG4
QG065-04	E04	HQP010612	NM_013975	LIG3
QG065-04	E05	HQP010609	NM_000234	LIG1
QG065-04	E06	HQP010070	NM_002253	KDR
QG065-04	E07	HQP009718	NM_001562	IL18
QG065-04	E08	HQP009693	NM_002187	IL12B
QG065-04	E09	HQP009692	NM_000882	IL12A
QG065-04	E10	HQP009639	NM_001556	IKBKB
QG065-04	E11	HQP009539	NM_000596	IGFBP1
QG065-04	E12	HQP009334	NM_198498	C11orf53
QG065-04	F01	HQP009101	NM_001168	BIRC5
QG065-04	F02	HQP009065	NM_000197	HSD17B3
QG065-04	F03	HQP008891	NM_000859	HMGCR
QG065-04	F04	HQP008482	NM_000848	GSTM2
QG065-04	F05	HQP008479	NM_001512	GSTA4
QG065-04	F06	HQP008476	NM_000846	GSTA2
QG065-04	F07	HQP008434	NM_013264	DDX25
QG065-04	F08	HQP008402	NM_001020825	NR3C1
QG065-04	F09	HQP008282	NM_002084	GPX3
QG065-04	F10	HQP008281	NM_002083	GPX2
QG065-04	F11	HQP008198	NM_173681	ATG9B
QG065-04	F12	HQP007755	NM_000516	GNAS
QG065-04	G01	HQP007703	NM_000168	GLI3
QG065-04	G02	HQP007526	NM_001018057	DKK3
QG065-04	G03	HQP007507	NM_000036	AMPD1
QG065-04	G04	HQP007213	NM_001634	AMD1
QG065-04	G05	HQP007110	NM_000819	GART
QG065-04	G06	HQP007058	NM_015617	PYGO1
QG065-04	G07	HQP006736	NM_012342	BAMBI
QG065-04	G08	HQP006715	NM_012415	RAD54B
QG065-04	G09	HQP006261	NM_014324	AMACR

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-04	G10	HQP006228	NM_001018078	FPGS
QG065-04	G11	HQP006144	NM_000802	FOLR1
QG065-04	G12	HQP005879	NM_002019	FLT1
QG065-04	H01	HGDC		
QG065-04	H02	HGDC		
QG065-04	H03	HQP006940	NM_002046	GAPDH
QG065-04	H04	HQP016381	NM_001101	ACTB
QG065-04	H05	HQP015171	NM_004048	B2M
QG065-04	H06	HQP006171	NM_012423	RPL13A
QG065-04	H07	HQP009026	NM_000194	HPRT1
QG065-04	H08	HQP054253	NR_003286	RN18S1
QG065-04	H09	RT		
QG065-04	H10	RT		
QG065-04	H11	PCR		
QG065-04	H12	PCR		