

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-01	A01	HQP018175	NM_000546	TP53
QG078-01	A02	HQP000370	NM_058195	CDKN2A
QG078-01	A03	HQP004976	NM_000400	ERCC2
QG078-01	A04	HQP018562	NM_006297	XRCC1
QG078-01	A05	HQP003772	NM_000499	CYP1A1
QG078-01	A06	HQP017733	NM_004333	BRAF
QG078-01	A07	HQP009685	NM_000572	IL10
QG078-01	A08	HQP005087	NM_000690	ALDH2
QG078-01	A09	HQP018474	NM_000376	VDR
QG078-01	A10	HQP011547	NM_005957	MTHFR
QG078-01	A11	HQP003817	NM_000773	CYP2E1
QG078-01	A12	HQP002331	NM_000668	ADH1B
QG078-01	B01	HQP004605	NM_005228	EGFR
QG078-01	B02	HQP054047	BC008403	HLA-DRB1
QG078-01	B03	HQP018564	NM_005432	XRCC3
QG078-01	B04	HQP016204	NM_053056	CCND1
QG078-01	B05	HQP017753	NM_000059	BRCA2
QG078-01	B06	HQP015598	NM_000963	PTGS2
QG078-01	B07	HQP004599	NM_001963	EGF
QG078-01	B08	HQP018556	NM_004628	XPC
QG078-01	B09	HQP011135	NM_002392	MDM2
QG078-01	B10	HQP010133	NM_004985	KRAS
QG078-01	B11	HQP018141	NM_000594	TNF
QG078-01	B12	HQP009641	NM_000576	IL1B
QG078-01	C01	HQP002413	NM_000669	ADH1C
QG078-01	C02	HQP018044	NM_000660	TGFB1
QG078-01	C03	HQP012008	NM_000275	OCA2
QG078-01	C04	HQP009670	NM_000600	IL6
QG078-01	C05	HQP001136	NM_000015	NAT2
QG078-01	C06	HQP000245	NM_000075	CDK4
QG078-01	C07	HQP015530	NM_000264	PTCH1
QG078-01	C08	HQP011914	NM_002524	NRAS
QG078-01	C09	HQP009678	NM_000584	IL8
QG078-01	C10	HQP018475	NM_001025366	VEGFA
QG078-01	C11	HQP018343	NM_000372	TYR
QG078-01	C12	HQP011303	NM_001672	ASIP
QG078-01	D01	HQP011257	NM_002422	MMP3
QG078-01	D02	HQP010099	NM_000222	KIT
QG078-01	D03	HQP009671	NM_000639	FASLG
QG078-01	D04	HQP004985	NM_000123	ERCC5
QG078-01	D05	HQP016418	NM_004327	BCR

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-01	D06	HQP009662	NM_000589	IL4
QG078-01	D07	HQP008849	NM_002116	HLA-A
QG078-01	D08	HQP003775	NM_000104	CYP1B1
QG078-01	D09	HQP003499	NM_001037631	CTLA4
QG078-01	D10	HQP001396	NM_001005735	CHEK2
QG078-01	D11	HQP012021	NM_002542	OGG1
QG078-01	D12	HQP004317	NM_000903	NQO1
QG078-01	E01	HQP018563	NM_005431	XRCC2
QG078-01	E02	HQP018349	NM_000550	TYRP1
QG078-01	E03	HQP012457	NM_001012509	SLC45A2
QG078-01	E04	HQP011554	NM_000254	MTR
QG078-01	E05	HQP011263	NM_004994	MMP9
QG078-01	E06	HQP011255	NM_002421	MMP1
QG078-01	E07	HQP009781	NM_002460	IRF4
QG078-01	E08	HQP007755	NM_000516	GNAS
QG078-01	E09	HQP004984	NM_005236	ERCC4
QG078-01	E10	HQP004975	NM_202001	ERCC1
QG078-01	E11	HQP003814	NM_000106	CYP2D6
QG078-01	E12	HQP019829	NM_030782	CLPTM1L
QG078-01	F01	HQP018176	NM_005657	TP53BP1
QG078-01	F02	HQP016212	NM_000657	BCL2
QG078-01	F03	HQP013100	NM_000927	ABCB1
QG078-01	F04	HQP011687	NM_002485	NBN
QG078-01	F05	HQP011540	NM_002451	MTAP
QG078-01	F06	HQP011256	NM_004530	MMP2
QG078-01	F07	HQP009718	NM_001562	IL18
QG078-01	F08	HQP009645	NM_000577	IL1RN
QG078-01	F09	HQP009061	NM_001641	APEX1
QG078-01	F10	HQP008483	NM_000849	GSTM3
QG078-01	F11	HQP005440	NM_022963	FGFR4
QG078-01	F12	HQP004948	NM_000120	EPHX1
QG078-01	G01	HQP004081	NM_000789	ACE
QG078-01	G02	HQP018180	NM_005427	TP73
QG078-01	G03	HQP023467	NM_000662	NAT1
QG078-01	G04	HQP020803	NM_139321	ATRN
QG078-01	G05	HQP020418	NM_001004426	PLA2G6
QG078-01	G06	HQP018966	NM_001080124	CASP8
QG078-01	G07	HQP018565	NM_003401	XRCC4
QG078-01	G08	HQP018149	NM_001066	TNFRSF1B
QG078-01	G09	HQP017636	NM_006941	SOX10
QG078-01	G10	HQP017616	NM_000636	SOD2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-01	G11	HQP016670	NM_001033886	CXCL12
QG078-01	G12	HQP015535	NM_000314	PTEN
QG078-01	H01	HGDC		
QG078-01	H02	HGDC		
QG078-01	H03	HQP006940	NM_002046	GAPDH
QG078-01	H04	HQP016381	NM_001101	ACTB
QG078-01	H05	HQP015171	NM_004048	B2M
QG078-01	H06	HQP006171	NM_012423	RPL13A
QG078-01	H07	HQP009026	NM_000194	HPRT1
QG078-01	H08	HQP054253	NR_003286	RN18S1
QG078-01	H09	RT		
QG078-01	H10	RT		
QG078-01	H11	PCR		
QG078-01	H12	PCR		
QG078-02	A01	HQP013633	NM_005037	PPARG
QG078-02	A02	HQP013467	NM_000939	POMC
QG078-02	A03	HQP012501	NM_016341	PLCE1
QG078-02	A04	HQP011807	NM_003998	NFKB1
QG078-02	A05	HQP011597	NM_002467	MYC
QG078-02	A06	HQP011555	NM_002454	MTRR
QG078-02	A07	HQP011491	NM_000251	MSH2
QG078-02	A08	HQP011309	NM_000250	MPO
QG078-02	A09	HQP011235	NM_000249	MLH1
QG078-02	A10	HQP011223	NM_000248	MITF
QG078-02	A11	HQP010613	NM_002312	LIG4
QG078-02	A12	HQP009672	NM_000565	IL6R
QG078-02	B01	HQP009649	NM_000586	IL2
QG078-02	B02	HQP004564	NM_000115	EDNRB
QG078-02	B03	HQP003841	NM_000777	CYP3A5
QG078-02	B04	HQP003804	NM_000766	CYP2A13
QG078-02	B05	HQP002192	NM_153646	SLC24A4
QG078-02	B06	HQP004414	NM_006892	DNMT3B
QG078-02	B07	HQP054684	NM_001033	RRM1
QG078-02	B08	HQP054054	BC039243	FGFR2
QG078-02	B09	HQP053910	NM_130851	BMP4
QG078-02	B10	HQP023466	NM_004360	CDH1
QG078-02	B11	HQP023394	NM_005462	MAGEC1
QG078-02	B12	HQP023341	NM_138448	ACYP2
QG078-02	C01	HQP023203	NM_001785	CDA
QG078-02	C02	HQP022699	NM_006139	CD28
QG078-02	C03	HQP022630	NM_001031689	PLAA

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-02	C04	HQP022500	NM_001040280	CD83
QG078-02	C05	HQP021829	NM_014240	LIMD1
QG078-02	C06	HQP021597	NM_003873	NRP1
QG078-02	C07	HQP021542	NM_003834	RGS11
QG078-02	C08	HQP021518	NM_000071	CBS
QG078-02	C09	HQP021252	NM_003664	AP3B1
QG078-02	C10	HQP021205	NM_033345	RGS8
QG078-02	C11	HQP021125	NM_001005476	PKP4
QG078-02	C12	HQP021023	NM_003617	RGS5
QG078-02	D01	HQP020297	NM_004346	CASP3
QG078-02	D02	HQP019568	NM_025113	C13orf18
QG078-02	D03	HQP019462	NM_024945	RMI1
QG078-02	D04	HQP019310	NM_024780	TMC5
QG078-02	D05	HQP019268	NM_006304	SHFM1
QG078-02	D06	HQP019189	NM_001006636	GTDC1
QG078-02	D07	HQP019141	NM_024608	NEIL1
QG078-02	D08	HQP018802	NM_001008540	CXCR4
QG078-02	D09	HQP018568	NM_021141	XRCC5
QG078-02	D10	HQP018552	NM_000379	XDH
QG078-02	D11	HQP018372	NM_152653	UBE2E2
QG078-02	D12	HQP017897	NM_006755	TALDO1
QG078-02	E01	HQP017811	NM_177536	SULT1A1
QG078-02	E02	HQP017794	NM_000455	STK11
QG078-02	E03	HQP017673	NM_000582	SPP1
QG078-02	E04	HQP017238	NM_006928	SILV
QG078-02	E05	HQP017124	NM_022743	SMYD3
QG078-02	E06	HQP017001	NM_000337	SGCD
QG078-02	E07	HQP016332	NM_002939	RNH1
QG078-02	E08	HQP016295	NM_002935	RNASE3
QG078-02	E09	HQP016269	NM_002928	RGS16
QG078-02	E10	HQP016268	NM_002927	RGS13
QG078-02	E11	HQP016263	NM_001005339	RGS10
QG078-02	E12	HQP016262	NM_002924	RGS7
QG078-02	F01	HQP016259	NM_005613	RGS4
QG078-02	F02	HQP016253	NM_017790	RGS3
QG078-02	F03	HQP016252	NM_002923	RGS2
QG078-02	F04	HQP016131	NM_000321	RB1
QG078-02	F05	HQP015997	NM_021205	RHOU
QG078-02	F06	HQP015897	NM_006264	PTPN13
QG078-02	F07	HQP015845	NM_020964	KIAA1632
QG078-02	F08	HQP015807	NM_020932	MAGEE1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-02	F09	HQP015771	NM_020884	MYH7B
QG078-02	F10	HQP015756	NM_020859	SHROOM3
QG078-02	F11	HQP015669	NM_020752	GPR158
QG078-02	F12	HQP015589	NM_020682	AS3MT
QG078-02	G01	HQP015179	NM_020124	IFNK
QG078-02	G02	HQP015124	NM_016567	BCCIP
QG078-02	G03	HQP014851	NM_018847	KLHL9
QG078-02	G04	HQP014839	NM_018715	RCC2
QG078-02	G05	HQP014116	NM_001011722	ARHGEF10L
QG078-02	G06	HQP013150	NM_006218	PIK3CA
QG078-02	G07	HQP013147	NM_002647	PIK3C3
QG078-02	G08	HQP013144	NM_002645	PIK3C2A
QG078-02	G09	HQP012796	NM_016448	DTL
QG078-02	G10	HQP012755	NM_001037872	REV1
QG078-02	G11	HQP012605	NM_016583	PLUNC
QG078-02	G12	HQP012591	NM_016562	TLR7
QG078-02	H01	HGDC		
QG078-02	H02	HGDC		
QG078-02	H03	HQP006940	NM_002046	GAPDH
QG078-02	H04	HQP016381	NM_001101	ACTB
QG078-02	H05	HQP015171	NM_004048	B2M
QG078-02	H06	HQP006171	NM_012423	RPL13A
QG078-02	H07	HQP009026	NM_000194	HPRT1
QG078-02	H08	HQP054253	NR_003286	RN18S1
QG078-02	H09	RT		
QG078-02	H10	RT		
QG078-02	H11	PCR		
QG078-02	H12	PCR		
QG078-03	A01	HQP012210	NM_181461	PAX3
QG078-03	A02	HQP011926	NM_002197	ACO1
QG078-03	A03	HQP011866	NM_000625	NOS2A
QG078-03	A04	HQP011810	NM_020529	NFKBIA
QG078-03	A05	HQP011800	NM_006164	NFE2L2
QG078-03	A06	HQP011648	NM_000260	MYO7A
QG078-03	A07	HQP011609	NM_002470	MYH3
QG078-03	A08	HQP011262	NM_002424	MMP8
QG078-03	A09	HQP010979	NM_001011548	MAGEA4
QG078-03	A10	HQP010977	NM_004988	MAGEA1
QG078-03	A11	HQP010961	NM_005359	SMAD4
QG078-03	A12	HQP010584	NM_002303	LEPR
QG078-03	B01	HQP010581	NM_000230	LEP

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-03	B02	HQP010141	NM_000424	KRT5
QG078-03	B03	HQP010086	NM_015868	KIR2DL3
QG078-03	B04	HQP009818	NM_000212	ITGB3
QG078-03	B05	HQP009693	NM_002187	IL12B
QG078-03	B06	HQP009692	NM_000882	IL12A
QG078-03	B07	HQP009664	NM_000418	IL4R
QG078-03	B08	HQP009640	NM_000575	IL1A
QG078-03	B09	HQP009558	NM_001553	IGFBP7
QG078-03	B10	HQP009529	NM_000612	IGF2
QG078-03	B11	HQP009518	NM_000618	IGF1
QG078-03	B12	HQP009463	NM_002176	IFNB1
QG078-03	C01	HQP009446	NM_002173	IFNA16
QG078-03	C02	HQP009427	NM_002169	IFNA5
QG078-03	C03	HQP009422	NM_000605	IFNA2
QG078-03	C04	HQP009353	NM_002168	IDH2
QG078-03	C05	HQP009351	NM_005896	IDH1
QG078-03	C06	HQP009181	NM_176891	IFNE1
QG078-03	C07	HQP009101	NM_001168	BIRC5
QG078-03	C08	HQP009027	NM_000195	HPS1
QG078-03	C09	HQP009019	NM_005143	HP
QG078-03	C10	HQP008934	NM_001160	APAF1
QG078-03	C11	HQP008832	NM_181054	HIF1A
QG078-03	C12	HQP008757	NM_000410	HFE
QG078-03	D01	HQP008669	NM_020804	PACSIN1
QG078-03	D02	HQP008554	NM_012092	ICOS
QG078-03	D03	HQP008479	NM_001512	GSTA4
QG078-03	D04	HQP008436	NM_000838	GRM1
QG078-03	D05	HQP008402	NM_001020825	NR3C1
QG078-03	D06	HQP008231	NM_001004125	TUSC1
QG078-03	D07	HQP008118	NM_173653	SLC9A9
QG078-03	D08	HQP007821	NM_019844	SLCO1B3
QG078-03	D09	HQP007803	NM_000173	GP1BA
QG078-03	D10	HQP007767	NM_002075	GNB3
QG078-03	D11	HQP007743	NM_002067	GNA11
QG078-03	D12	HQP007346	NM_000514	GDNF
QG078-03	E01	HQP007098	NM_015668	RGS22
QG078-03	E02	HQP006547	NM_001469	XRCC6
QG078-03	E03	HQP006460	NM_145043	NEIL2
QG078-03	E04	HQP006441	NM_000350	ABCA4
QG078-03	E05	HQP006227	NM_012387	PADI4
QG078-03	E06	HQP006038	NM_015316	PPP1R13B

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-03	E07	HQP006010	NM_033071	SYNE1
QG078-03	E08	HQP005965	NM_015246	MGRN1
QG078-03	E09	HQP005925	NM_015210	KIAA0802
QG078-03	E10	HQP005899	NM_182734	PLCB1
QG078-03	E11	HQP005879	NM_002019	FLT1
QG078-03	E12	HQP005736	NM_015061	JMJD2C
QG078-03	F01	HQP005434	NM_000142	FGFR3
QG078-03	F02	HQP005093	NM_004103	PTK2B
QG078-03	F03	HQP005058	NM_000130	F5
QG078-03	F04	HQP005052	NM_000506	F2
QG078-03	F05	HQP004986	NM_000124	ERCC6
QG078-03	F06	HQP004494	NM_000029	AGT
QG078-03	F07	HQP003888	NM_000102	CYP17A1
QG078-03	F08	HQP003171	NM_006437	PARP4
QG078-03	F09	HQP003120	NM_001618	PARP1
QG078-03	F10	HQP002701	NM_000673	ADH7
QG078-03	F11	HQP002534	NM_080476	PIGU
QG078-03	F12	HQP002301	NM_000667	ADH1A
QG078-03	G01	HQP002210	NM_000579	CCR5
QG078-03	G02	HQP001887	NM_133445	GRIN3A
QG078-03	G03	HQP001712	NM_052917	GALNT13
QG078-03	G04	HQP001682	NM_052890	PGLYRP2
QG078-03	G05	HQP001643	NM_000745	CHRNA5
QG078-03	G06	HQP001399	NM_007195	POLI
QG078-03	G07	HQP000771	NM_006480	RGS14
QG078-03	G08	HQP000396	NM_001262	CDKN2C
QG078-03	G09	HQP018876	NM_024071	ZFYVE21
QG078-03	G10	HQP016965	NM_130782	RGS18
QG078-03	G11	HQP016265	NM_002926	RGS12
QG078-03	G12	HQP012055	NM_016190	CRNN
QG078-03	H01	HGDC		
QG078-03	H02	HGDC		
QG078-03	H03	HQP006940	NM_002046	GAPDH
QG078-03	H04	HQP016381	NM_001101	ACTB
QG078-03	H05	HQP015171	NM_004048	B2M
QG078-03	H06	HQP006171	NM_012423	RPL13A
QG078-03	H07	HQP009026	NM_000194	HPRT1
QG078-03	H08	HQP054253	NR_003286	RN18S1
QG078-03	H09	RT		
QG078-03	H10	RT		
QG078-03	H11	PCR		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG078-03	H12	PCR		