

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-01	A01	HQP018175	NM_000546	TP53
QG062-01	A02	HQP009351	NM_005896	IDH1
QG062-01	A03	HQP004599	NM_001963	EGF
QG062-01	A04	HQP009353	NM_002168	IDH2
QG062-01	A05	HQP011135	NM_002392	MDM2
QG062-01	A06	HQP004605	NM_005228	EGFR
QG062-01	A07	HQP000370	NM_058195	CDKN2A
QG062-01	A08	HQP018562	NM_006297	XRCC1
QG062-01	A09	HQP018564	NM_005432	XRCC3
QG062-01	A10	HQP015535	NM_000314	PTEN
QG062-01	A11	HQP008849	NM_002116	HLA-A
QG062-01	A12	HQP054047	BC008403	HLA-DRB1
QG062-01	B01	HQP013006	NM_016434	RTEL1
QG062-01	B02	HQP011547	NM_005957	MTHFR
QG062-01	B03	HQP004976	NM_000400	ERCC2
QG062-01	B04	HQP013473	NM_000446	PON1
QG062-01	B05	HQP009697	NM_002188	IL13
QG062-01	B06	HQP009664	NM_000418	IL4R
QG062-01	B07	HQP018966	NM_001080124	CASP8
QG062-01	B08	HQP018568	NM_021141	XRCC5
QG062-01	B09	HQP018565	NM_003401	XRCC4
QG062-01	B10	HQP017615	NM_000454	SOD1
QG062-01	B11	HQP011807	NM_003998	NFKB1
QG062-01	B12	HQP011693	NM_000433	NCF2
QG062-01	C01	HQP011687	NM_002485	NBN
QG062-01	C02	HQP011554	NM_000254	MTR
QG062-01	C03	HQP010613	NM_002312	LIG4
QG062-01	C04	HQP009670	NM_000600	IL6
QG062-01	C05	HQP009662	NM_000589	IL4
QG062-01	C06	HQP008483	NM_000849	GSTM3
QG062-01	C07	HQP005843	NM_015157	PHLDB1
QG062-01	C08	HQP001136	NM_000015	NAT2
QG062-01	C09	HQP005011	NM_000031	ALAD
QG062-01	C10	HQP004317	NM_000903	NQO1
QG062-01	C11	HQP020707	NM_003579	RAD54L
QG062-01	C12	HQP018563	NM_005431	XRCC2
QG062-01	D01	HQP018552	NM_000379	XDH
QG062-01	D02	HQP017979	NM_000355	TCN2
QG062-01	D03	HQP017764	NM_007315	STAT1
QG062-01	D04	HQP017733	NM_004333	BRAF
QG062-01	D05	HQP017616	NM_000636	SOD2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-01	D06	HQP016746	NM_003005	SELP
QG062-01	D07	HQP016621	NM_002982	CCL2
QG062-01	D08	HQP016204	NM_053056	CCND1
QG062-01	D09	HQP016087	NM_134424	RAD52
QG062-01	D10	HQP016064	NM_018890	RAC1
QG062-01	D11	HQP015598	NM_000963	PTGS2
QG062-01	D12	HQP013100	NM_000927	ABCB1
QG062-01	E01	HQP011868	NM_000603	NOS3
QG062-01	E02	HQP011866	NM_000625	NOS2A
QG062-01	E03	HQP011778	NM_000268	NF2
QG062-01	E04	HQP011694	NM_000631	NCF4
QG062-01	E05	HQP011309	NM_000250	MPO
QG062-01	E06	HQP011257	NM_002422	MMP3
QG062-01	E07	HQP011255	NM_002421	MMP1
QG062-01	E08	HQP010609	NM_000234	LIG1
QG062-01	E09	HQP010133	NM_004985	KRAS
QG062-01	E10	HQP010095	NM_013289	KIR3DL1
QG062-01	E11	HQP010086	NM_015868	KIR2DL3
QG062-01	E12	HQP010083	NM_014218	KIR2DL1
QG062-01	F01	HQP009061	NM_001641	APEX1
QG062-01	F02	HQP008673	NM_015711	GLTSCR1
QG062-01	F03	HQP006547	NM_001469	XRCC6
QG062-01	F04	HQP006359	NM_000698	ALOX5
QG062-01	F05	HQP005280	NM_021642	FCGR2A
QG062-01	F06	HQP004975	NM_202001	ERCC1
QG062-01	F07	HQP003772	NM_000499	CYP1A1
QG062-01	F08	HQP003539	NM_001904	CTNNB1
QG062-01	F09	HQP003120	NM_001618	PARP1
QG062-01	F10	HQP001408	NM_007199	IRAK3
QG062-01	F11	HQP000245	NM_000075	CDK4
QG062-01	F12	HQP017618	NM_003102	SOD3
QG062-01	G01	HQP011865	NM_000620	NOS1
QG062-01	G02	HQP010851	NM_006151	LPO
QG062-01	G03	HQP054326	NM_001099287	ICHTHYIN
QG062-01	G04	HQP054030	NM_130398	EXO1
QG062-01	G05	HQP023467	NM_000662	NAT1
QG062-01	G06	HQP023466	NM_004360	CDH1
QG062-01	G07	HQP023300	NM_014813	LRIG2
QG062-01	G08	HQP023012	NM_006536	CLCA2
QG062-01	G09	HQP022998	NM_004917	KLK4
QG062-01	G10	HQP022584	NM_004787	SLIT2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-01	G11	HQP022490	NM_000591	CD14
QG062-01	G12	HQP022406	NM_004756	NUMBL
QG062-01	H01	HGDC		
QG062-01	H02	HGDC		
QG062-01	H03	HQP006940	NM_002046	GAPDH
QG062-01	H04	HQP016381	NM_001101	ACTB
QG062-01	H05	HQP015171	NM_004048	B2M
QG062-01	H06	HQP006171	NM_012423	RPL13A
QG062-01	H07	HQP009026	NM_000194	HPRT1
QG062-01	H08	HQP054253	NR_003286	RN18S1
QG062-01	H09	RT		
QG062-01	H10	RT		
QG062-01	H11	PCR		
QG062-01	H12	PCR		
QG062-02	A01	HQP022361	NM_004747	DLG5
QG062-02	A02	HQP022323	NM_004217	AURKB
QG062-02	A03	HQP021906	NM_001239	CCNH
QG062-02	A04	HQP021889	NM_003955	SOCS3
QG062-02	A05	HQP021834	NM_005092	TNFSF18
QG062-02	A06	HQP021636	NM_032454	STK19
QG062-02	A07	HQP021602	NM_003877	SOCS2
QG062-02	A08	HQP021518	NM_000071	CBS
QG062-02	A09	HQP021484	NM_033274	ADAM19
QG062-02	A10	HQP021399	NM_003745	SOCS1
QG062-02	A11	HQP021260	NM_033035	TSLP
QG062-02	A12	HQP021225	NM_006769	LMO4
QG062-02	B01	HQP021130	NM_003632	CNTNAP1
QG062-02	B02	HQP020906	NM_032622	LNX1
QG062-02	B03	HQP020499	NM_032169	ACAD11
QG062-02	B04	HQP019982	NM_001017388	TLR10
QG062-02	B05	HQP019941	NM_030931	DEFB126
QG062-02	B06	HQP019268	NM_006304	SHFM1
QG062-02	B07	HQP018854	NM_024051	C7orf24
QG062-02	B08	HQP018556	NM_004628	XPC
QG062-02	B09	HQP018545	NM_000553	WRN
QG062-02	B10	HQP018485	NM_000551	VHL
QG062-02	B11	HQP018475	NM_001025366	VEGFA
QG062-02	B12	HQP018466	NM_080682	VCAM1
QG062-02	C01	HQP018411	NM_003357	SCGB1A1
QG062-02	C02	HQP018391	NM_000606	C8G
QG062-02	C03	HQP018380	NM_000066	C8B

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-02	C04	HQP018353	NM_000587	C7
QG062-02	C05	HQP018340	NM_003331	TYK2
QG062-02	C06	HQP018326	NM_001736	C5AR1
QG062-02	C07	HQP018276	NM_000716	C4BPB
QG062-02	C08	HQP018259	NM_000715	C4BPA
QG062-02	C09	HQP018226	NM_000063	C2
QG062-02	C10	HQP018170	NM_172369	C1QC
QG062-02	C11	HQP018141	NM_000594	TNF
QG062-02	C12	HQP018115	NM_003265	TLR3
QG062-02	D01	HQP018113	NM_003263	TLR1
QG062-02	D02	HQP018106	NM_001212	C1QBP
QG062-02	D03	HQP018044	NM_000660	TGFB1
QG062-02	D04	HQP017811	NM_177536	SULT1A1
QG062-02	D05	HQP017770	NM_003151	STAT4
QG062-02	D06	HQP017767	NM_003150	STAT3
QG062-02	D07	HQP017766	NM_005419	STAT2
QG062-02	D08	HQP017753	NM_000059	BRCA2
QG062-02	D09	HQP017744	NM_001050	SSTR2
QG062-02	D10	HQP017702	NM_001725	BPI
QG062-02	D11	HQP017673	NM_000582	SPP1
QG062-02	D12	HQP017625	NM_005633	SOS1
QG062-02	E01	HQP017549	NM_030807	SLC2A11
QG062-02	E02	HQP017526	NM_003073	SMARCB1
QG062-02	E03	HQP016982	NM_001033855	DCLRE1C
QG062-02	E04	HQP016801	NM_022162	NOD2
QG062-02	E05	HQP016747	NM_003006	SELPLG
QG062-02	E06	HQP016744	NM_000450	SELE
QG062-02	E07	HQP016738	NM_003004	SECTM1
QG062-02	E08	HQP016644	NM_002991	CCL24
QG062-02	E09	HQP016639	NM_004591	CCL20
QG062-02	E10	HQP016637	NM_002988	CCL18
QG062-02	E11	HQP016630	NM_005408	CCL13
QG062-02	E12	HQP016629	NM_002986	CCL11
QG062-02	F01	HQP016626	NM_002985	CCL5
QG062-02	F02	HQP016620	NM_002981	CCL1
QG062-02	F03	HQP016581	NM_006919	SERPINB3
QG062-02	F04	HQP016565	NM_001710	CFB
QG062-02	F05	HQP016500	NM_001032295	SERPING1
QG062-02	F06	HQP016224	NM_002913	RFC1
QG062-02	F07	HQP016170	NM_021632	ZNF350
QG062-02	F08	HQP016131	NM_000321	RB1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-02	F09	HQP016108	NM_000055	BCHE
QG062-02	F10	HQP016088	NM_002880	RAF1
QG062-02	F11	HQP016082	NM_133509	RAD51L1
QG062-02	F12	HQP016076	NM_002874	RAD23B
QG062-02	G01	HQP015946	NM_000465	BARD1
QG062-02	G02	HQP015912	NM_001040712	PTPRD
QG062-02	G03	HQP015625	NM_020706	SFRS15
QG062-02	G04	HQP015537	NM_000953	PTGDR
QG062-02	G05	HQP015248	NM_001080452	GPR108
QG062-02	G06	HQP015240	NM_020162	DHX33
QG062-02	G07	HQP015102	NM_001012965	KLK6
QG062-02	G08	HQP015096	NM_005046	KLK7
QG062-02	G09	HQP014909	NM_030662	MAP2K2
QG062-02	G10	HQP014907	NM_002755	MAP2K1
QG062-02	G11	HQP014732	NM_018416	FOXJ2
QG062-02	G12	HQP014543	NM_017797	BTBD2
QG062-02	H01	HGDC		
QG062-02	H02	HGDC		
QG062-02	H03	HQP006940	NM_002046	GAPDH
QG062-02	H04	HQP016381	NM_001101	ACTB
QG062-02	H05	HQP015171	NM_004048	B2M
QG062-02	H06	HQP006171	NM_012423	RPL13A
QG062-02	H07	HQP009026	NM_000194	HPRT1
QG062-02	H08	HQP054253	NR_003286	RN18S1
QG062-02	H09	RT		
QG062-02	H10	RT		
QG062-02	H11	PCR		
QG062-02	H12	PCR		
QG062-03	A01	HQP014462	NM_017509	KLK15
QG062-03	A02	HQP013496	NM_019009	TOLLIP
QG062-03	A03	HQP013422	NM_002690	POLB
QG062-03	A04	HQP013388	NM_017442	TLR9
QG062-03	A05	HQP013150	NM_006218	PIK3CA
QG062-03	A06	HQP013133	NM_005025	SERPINI1
QG062-03	A07	HQP013071	NM_006212	PFKFB2
QG062-03	A08	HQP012755	NM_001037872	REV1
QG062-03	A09	HQP012584	NM_016546	C1RL
QG062-03	A10	HQP012441	NM_016123	IRAK4
QG062-03	A11	HQP012420	NM_002592	PCNA
QG062-03	A12	HQP012166	NM_022047	DEF6
QG062-03	B01	HQP012155	NM_002575	SERPINB2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-03	B02	HQP012154	NM_000602	SERPINE1
QG062-03	B03	HQP012062	NM_001008505	OPRM1
QG062-03	B04	HQP011938	NM_001007792	NTRK1
QG062-03	B05	HQP011810	NM_020529	NFKBIA
QG062-03	B06	HQP011728	NM_004550	NDUFS2
QG062-03	B07	HQP011663	NM_005967	NAB2
QG062-03	B08	HQP011620	NM_000488	SERPINC1
QG062-03	B09	HQP011603	NM_002468	MYD88
QG062-03	B10	HQP011555	NM_002454	MTRR
QG062-03	B11	HQP011491	NM_000251	MSH2
QG062-03	B12	HQP011320	NM_005590	MRE11A
QG062-03	C01	HQP011263	NM_004994	MMP9
QG062-03	C02	HQP011219	NM_002415	MIF
QG062-03	C03	HQP011171	NM_000244	MEN1
QG062-03	C04	HQP011113	NM_002389	CD46
QG062-03	C05	HQP011085	NM_001025081	MBP
QG062-03	C06	HQP010996	NM_022438	MAL
QG062-03	C07	HQP010932	NM_005582	CD180
QG062-03	C08	HQP010690	NM_002315	LMO1
QG062-03	C09	HQP010660	NM_005570	LMAN1
QG062-03	C10	HQP010560	NM_004139	LBP
QG062-03	C11	HQP010104	NM_000892	KLKB1
QG062-03	C12	HQP010100	NM_002257	KLK1
QG062-03	D01	HQP010092	NM_012313	KIR2DS3
QG062-03	D02	HQP010070	NM_002253	KDR
QG062-03	D03	HQP009851	NM_000215	JAK3
QG062-03	D04	HQP009776	NM_001570	IRAK2
QG062-03	D05	HQP009693	NM_002187	IL12B
QG062-03	D06	HQP009681	NM_001557	IL8RB
QG062-03	D07	HQP009679	NM_000634	IL8RA
QG062-03	D08	HQP009678	NM_000584	IL8
QG062-03	D09	HQP009556	NM_000041	APOE
QG062-03	D10	HQP009544	NM_000598	IGFBP3
QG062-03	D11	HQP009543	NM_182625	FLJ40869
QG062-03	D12	HQP009529	NM_000612	IGF2
QG062-03	E01	HQP009523	NM_000875	IGF1R
QG062-03	E02	HQP009518	NM_000618	IGF1
QG062-03	E03	HQP009201	NM_001039132	ICAM4
QG062-03	E04	HQP009139	NM_001643	APOA2
QG062-03	E05	HQP009085	NM_002155	HSPA6
QG062-03	E06	HQP009036	NM_005343	HRAS

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-03	E07	HQP009024	NM_000038	APC
QG062-03	E08	HQP008846	NM_001607	ACAA1
QG062-03	E09	HQP008808	NM_021155	CD209
QG062-03	E10	HQP008801	NM_001010931	HGF
QG062-03	E11	HQP008198	NM_173681	ATG9B
QG062-03	E12	HQP007767	NM_002075	GNB3
QG062-03	F01	HQP007722	NM_014905	GLS
QG062-03	F02	HQP007100	NM_015670	SENP3
QG062-03	F03	HQP006715	NM_012415	RAD54B
QG062-03	F04	HQP006427	NM_001039130	ALOX15B
QG062-03	F05	HQP006425	NM_001140	ALOX15
QG062-03	F06	HQP006424	NM_145659	IL27
QG062-03	F07	HQP006396	NM_153289	DEFB119
QG062-03	F08	HQP006375	NM_001629	ALOX5AP
QG062-03	F09	HQP006356	NM_000697	ALOX12
QG062-03	F10	HQP006296	NM_015364	LY96
QG062-03	F11	HQP006131	NM_014294	TRAM1
QG062-03	F12	HQP005959	NM_001002294	FMO3
QG062-03	G01	HQP005939	NM_002021	FMO1
QG062-03	G02	HQP005589	NM_012072	CD93
QG062-03	G03	HQP005403	NM_002006	FGF2
QG062-03	G04	HQP005242	NM_002002	FCER2
QG062-03	G05	HQP005240	NM_004106	FCER1G
QG062-03	G06	HQP005238	NM_002001	FCER1A
QG062-03	G07	HQP005100	NM_000691	ALDH3A1
QG062-03	G08	HQP005058	NM_000130	F5
QG062-03	G09	HQP004984	NM_005236	ERCC4
QG062-03	G10	HQP004971	NM_001982	ERBB3
QG062-03	G11	HQP004948	NM_000120	EPHX1
QG062-03	G12	HQP004445	NM_000794	DRD1
QG062-03	H01	HGDC		
QG062-03	H02	HGDC		
QG062-03	H03	HQP006940	NM_002046	GAPDH
QG062-03	H04	HQP016381	NM_001101	ACTB
QG062-03	H05	HQP015171	NM_004048	B2M
QG062-03	H06	HQP006171	NM_012423	RPL13A
QG062-03	H07	HQP009026	NM_000194	HPRT1
QG062-03	H08	HQP054253	NR_003286	RN18S1
QG062-03	H09	RT		
QG062-03	H10	RT		
QG062-03	H11	PCR		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG062-03	H12	PCR		