

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG066-01	A01	HQP018175	NM_000546	TP53
QG066-01	A02	HQP054047	BC008403	HLA-DRB1
QG066-01	A03	HQP018141	NM_000594	TNF
QG066-01	A04	HQP011547	NM_005957	MTHFR
QG066-01	A05	HQP009685	NM_000572	IL10
QG066-01	A06	HQP018562	NM_006297	XRCC1
QG066-01	A07	HQP008849	NM_002116	HLA-A
QG066-01	A08	HQP003772	NM_000499	CYP1A1
QG066-01	A09	HQP009671	NM_000639	FASLG
QG066-01	A10	HQP003499	NM_001037631	CTLA4
QG066-01	A11	HQP018044	NM_000660	TGFB1
QG066-01	A12	HQP017899	NM_000593	TAP1
QG066-01	B01	HQP009645	NM_000577	IL1RN
QG066-01	B02	HQP017900	NM_000544	TAP2
QG066-01	B03	HQP016204	NM_053056	CCND1
QG066-01	B04	HQP009670	NM_000600	IL6
QG066-01	B05	HQP009641	NM_000576	IL1B
QG066-01	B06	HQP009693	NM_002187	IL12B
QG066-01	B07	HQP004605	NM_005228	EGFR
QG066-01	B08	HQP003817	NM_000773	CYP2E1
QG066-01	B09	HQP012021	NM_002542	OGG1
QG066-01	B10	HQP004317	NM_000903	NQO1
QG066-01	B11	HQP022699	NM_006139	CD28
QG066-01	B12	HQP018564	NM_005432	XRCC3
QG066-01	C01	HQP011135	NM_002392	MDM2
QG066-01	C02	HQP009718	NM_001562	IL18
QG066-01	C03	HQP008554	NM_012092	ICOS
QG066-01	C04	HQP004948	NM_000120	EPHX1
QG066-01	C05	HQP001136	NM_000015	NAT2
QG066-01	C06	HQP022500	NM_001040280	CD83
QG066-01	C07	HQP018475	NM_001025366	VEGFA
QG066-01	C08	HQP017616	NM_000636	SOD2
QG066-01	C09	HQP015598	NM_000963	PTGS2
QG066-01	C10	HQP015311	NM_002800	PSMB9
QG066-01	C11	HQP015292	NM_148919	PSMB8
QG066-01	C12	HQP011866	NM_000625	NOS2A
QG066-01	D01	HQP011554	NM_000254	MTR
QG066-01	D02	HQP011309	NM_000250	MPO
QG066-01	D03	HQP011255	NM_002421	MMP1
QG066-01	D04	HQP009778	NM_002198	IRF1
QG066-01	D05	HQP009692	NM_000882	IL12A

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG066-01	D06	HQP009061	NM_001641	APEX1
QG066-01	D07	HQP002210	NM_000579	CCR5
QG066-01	D08	HQP018180	NM_005427	TP73
QG066-01	D09	HQP010096	NM_006737	KIR3DL2
QG066-01	D10	HQP054686	NM_001123396	LOC729230
QG066-01	D11	HQP054057	BC071181	TGFBR1
QG066-01	D12	HQP054030	NM_130398	EXO1
QG066-01	E01	HQP023466	NM_004360	CDH1
QG066-01	E02	HQP023354	NM_001254	CDC6
QG066-01	E03	HQP022747	NM_175862	CD86
QG066-01	E04	HQP022722	NM_005191	CD80
QG066-01	E05	HQP021603	NM_003878	GGH
QG066-01	E06	HQP018556	NM_004628	XPC
QG066-01	E07	HQP017903	NM_172208	TAPBP
QG066-01	E08	HQP017764	NM_007315	STAT1
QG066-01	E09	HQP017759	NM_021978	ST14
QG066-01	E10	HQP017753	NM_000059	BRCA2
QG066-01	E11	HQP017498	NM_003062	SLIT3
QG066-01	E12	HQP017259	NM_006747	SIPA1
QG066-01	F01	HQP016670	NM_001033886	CXCL12
QG066-01	F02	HQP016224	NM_002913	RFC1
QG066-01	F03	HQP015946	NM_000465	BARD1
QG066-01	F04	HQP014161	NM_002716	PPP2R1B
QG066-01	F05	HQP013312	NM_000534	PMS1
QG066-01	F06	HQP013150	NM_006218	PIK3CA
QG066-01	F07	HQP012755	NM_001037872	REV1
QG066-01	F08	HQP012420	NM_002592	PCNA
QG066-01	F09	HQP011868	NM_000603	NOS3
QG066-01	F10	HQP011687	NM_002485	NBN
QG066-01	F11	HQP011555	NM_002454	MTRR
QG066-01	F12	HQP011263	NM_004994	MMP9
QG066-01	G01	HQP011256	NM_004530	MMP2
QG066-01	G02	HQP010966	NM_005904	SMAD7
QG066-01	G03	HQP010613	NM_002312	LIG4
QG066-01	G04	HQP010095	NM_013289	KIR3DL1
QG066-01	G05	HQP010089	NM_002255	KIR2DL4
QG066-01	G06	HQP010086	NM_015868	KIR2DL3
QG066-01	G07	HQP010083	NM_014218	KIR2DL1
QG066-01	G08	HQP010070	NM_002253	KDR
QG066-01	G09	HQP009851	NM_000215	JAK3
QG066-01	G10	HQP009780	NM_001571	IRF3

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG066-01	G11	HQP009681	NM_001557	IL8RB
QG066-01	G12	HQP009679	NM_000634	IL8RA
QG066-01	H01	HGDC		
QG066-01	H02	HGDC		
QG066-01	H03	HQP006940	NM_002046	GAPDH
QG066-01	H04	HQP016381	NM_001101	ACTB
QG066-01	H05	HQP015171	NM_004048	B2M
QG066-01	H06	HQP006171	NM_012423	RPL13A
QG066-01	H07	HQP009026	NM_000194	HPRT1
QG066-01	H08	HQP054253	NR_003286	RN18S1
QG066-01	H09	RT		
QG066-01	H10	RT		
QG066-01	H11	PCR		
QG066-01	H12	PCR		