

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG069-01	A01	HQP018485	NM_000551	VHL
QG069-01	A02	HQP009685	NM_000572	IL10
QG069-01	A03	HQP018475	NM_001025366	VEGFA
QG069-01	A04	HQP018474	NM_000376	VDR
QG069-01	A05	HQP013100	NM_000927	ABCB1
QG069-01	A06	HQP011255	NM_002421	MMP1
QG069-01	A07	HQP018562	NM_006297	XRCC1
QG069-01	A08	HQP018556	NM_004628	XPC
QG069-01	A09	HQP018141	NM_000594	TNF
QG069-01	A10	HQP011547	NM_005957	MTHFR
QG069-01	A11	HQP008832	NM_181054	HIF1A
QG069-01	A12	HQP004976	NM_000400	ERCC2
QG069-01	B01	HQP018564	NM_005432	XRCC3
QG069-01	B02	HQP018546	NM_000378	WT1
QG069-01	B03	HQP017906	NM_004178	TARBP2
QG069-01	B04	HQP016689	NM_003000	SDHB
QG069-01	B05	HQP015667	NM_020750	XPO5
QG069-01	B06	HQP012169	NM_015721	GEMIN4
QG069-01	B07	HQP011257	NM_002422	MMP3
QG069-01	B08	HQP009662	NM_000589	IL4
QG069-01	B09	HQP007287	NM_012199	EIF2C1
QG069-01	B10	HQP005446	NM_000143	FH
QG069-01	B11	HQP003841	NM_000777	CYP3A5
QG069-01	B12	HQP003499	NM_001037631	CTLA4
QG069-01	C01	HQP001136	NM_000015	NAT2
QG069-01	C02	HQP004985	NM_000123	ERCC5
QG069-01	C03	HQP054057	BC071181	TGFBR1
QG069-01	C04	HQP054047	BC008403	HLA-DRB1
QG069-01	C05	HQP023466	NM_004360	CDH1
QG069-01	C06	HQP022833	NM_005505	SCARB1
QG069-01	C07	HQP022764	NM_004832	GSTO1
QG069-01	C08	HQP020472	NM_032141	CCDC55
QG069-01	C09	HQP019141	NM_024608	NEIL1
QG069-01	C10	HQP018966	NM_001080124	CASP8
QG069-01	C11	HQP018895	NM_001080432	FTO
QG069-01	C12	HQP018668	NM_001216	CA9
QG069-01	D01	HQP018342	NM_001071	TYMS
QG069-01	D02	HQP018175	NM_000546	TP53
QG069-01	D03	HQP018044	NM_000660	TGFB1
QG069-01	D04	HQP017835	NM_003174	SVIL
QG069-01	D05	HQP017753	NM_000059	BRCA2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG069-01	D06	HQP016694	NM_001035511	SDHC
QG069-01	D07	HQP016626	NM_002985	CCL5
QG069-01	D08	HQP016621	NM_002982	CCL2
QG069-01	D09	HQP016526	NM_002957	RXRA
QG069-01	D10	HQP016281	NM_002931	RING1
QG069-01	D11	HQP016224	NM_002913	RFC1
QG069-01	D12	HQP016212	NM_000657	BCL2
QG069-01	E01	HQP016204	NM_053056	CCND1
QG069-01	E02	HQP015832	NM_017893	SEMA4G
QG069-01	E03	HQP014841	NM_019099	C1orf183
QG069-01	E04	HQP014766	NM_018462	C3orf10
QG069-01	E05	HQP014292	NM_018315	FBXW7
QG069-01	E06	HQP013633	NM_005037	PPARG
QG069-01	E07	HQP012062	NM_001008505	OPRM1
QG069-01	E08	HQP011602	NM_005378	MYCN
QG069-01	E09	HQP011554	NM_000254	MTR
QG069-01	E10	HQP011491	NM_000251	MSH2
QG069-01	E11	HQP011235	NM_000249	MLH1
QG069-01	E12	HQP011135	NM_002392	MDM2
QG069-01	F01	HQP011110	NM_005915	MCM6
QG069-01	F02	HQP010966	NM_005904	SMAD7
QG069-01	F03	HQP010133	NM_004985	KRAS
QG069-01	F04	HQP010070	NM_002253	KDR
QG069-01	F05	HQP009818	NM_000212	ITGB3
QG069-01	F06	HQP009801	NM_000044	AR
QG069-01	F07	HQP009671	NM_000639	FASLG
QG069-01	F08	HQP009670	NM_000600	IL6
QG069-01	F09	HQP009664	NM_000418	IL4R
QG069-01	F10	HQP009640	NM_000575	IL1A
QG069-01	F11	HQP009556	NM_000041	APOE
QG069-01	F12	HQP009544	NM_000598	IGFBP3
QG069-01	G01	HQP008849	NM_002116	HLA-A
QG069-01	G02	HQP008483	NM_000849	GSTM3
QG069-01	G03	HQP008107	NM_020345	NKIRAS1
QG069-01	G04	HQP007755	NM_000516	GNAS
QG069-01	G05	HQP007558	NM_012154	EIF2C2
QG069-01	G06	HQP007527	NM_014421	DKK2
QG069-01	G07	HQP007526	NM_001018057	DKK3
QG069-01	G08	HQP007525	NM_014420	DKK4
QG069-01	G09	HQP006425	NM_001140	ALOX15
QG069-01	G10	HQP006152	NM_015345	DAAM2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG069-01	G11	HQP005434	NM_000142	FGFR3
QG069-01	G12	HQP004998	NM_000125	ESR1
QG069-01	H01	HGDC		
QG069-01	H02	HGDC		
QG069-01	H03	HQP006940	NM_002046	GAPDH
QG069-01	H04	HQP016381	NM_001101	ACTB
QG069-01	H05	HQP015171	NM_004048	B2M
QG069-01	H06	HQP006171	NM_012423	RPL13A
QG069-01	H07	HQP009026	NM_000194	HPRT1
QG069-01	H08	HQP054253	NR_003286	RN18S1
QG069-01	H09	RT		
QG069-01	H10	RT		
QG069-01	H11	PCR		
QG069-01	H12	PCR		