

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG072-01	A01	HQP018175	NM_000546	TP53
QG072-01	A02	HQP018141	NM_000594	TNF
QG072-01	A03	HQP008757	NM_000410	HFE
QG072-01	A04	HQP011547	NM_005957	MTHFR
QG072-01	A05	HQP009685	NM_000572	IL10
QG072-01	A06	HQP009641	NM_000576	IL1B
QG072-01	A07	HQP001136	NM_000015	NAT2
QG072-01	A08	HQP018562	NM_006297	XRCC1
QG072-01	A09	HQP018044	NM_000660	TGFB1
QG072-01	A10	HQP013581	NM_019077	UGT1A7
QG072-01	A11	HQP003817	NM_000773	CYP2E1
QG072-01	A12	HQP003772	NM_000499	CYP1A1
QG072-01	B01	HQP054047	BC008403	HLA-DRB1
QG072-01	B02	HQP009670	NM_000600	IL6
QG072-01	B03	HQP011263	NM_004994	MMP9
QG072-01	B04	HQP011135	NM_002392	MDM2
QG072-01	B05	HQP009718	NM_001562	IL18
QG072-01	B06	HQP005087	NM_000690	ALDH2
QG072-01	B07	HQP004948	NM_000120	EPHX1
QG072-01	B08	HQP004599	NM_001963	EGF
QG072-01	B09	HQP023467	NM_000662	NAT1
QG072-01	B10	HQP018556	NM_004628	XPC
QG072-01	B11	HQP017616	NM_000636	SOD2
QG072-01	B12	HQP016670	NM_001033886	CXCL12
QG072-01	C01	HQP016204	NM_053056	CCND1
QG072-01	C02	HQP011257	NM_002422	MMP3
QG072-01	C03	HQP011255	NM_002421	MMP1
QG072-01	C04	HQP009801	NM_000044	AR
QG072-01	C05	HQP009692	NM_000882	IL12A
QG072-01	C06	HQP009645	NM_000577	IL1RN
QG072-01	C07	HQP004605	NM_005228	EGFR
QG072-01	C08	HQP002671	NM_000754	COMT
QG072-01	C09	HQP013597	NM_021027	UGT1A9
QG072-01	C10	HQP023354	NM_001254	CDC6
QG072-01	C11	HQP018802	NM_001008540	CXCR4
QG072-01	C12	HQP018475	NM_001025366	VEGFA
QG072-01	D01	HQP018342	NM_001071	TYMS
QG072-01	D02	HQP011810	NM_020529	NFKBIA
QG072-01	D03	HQP011807	NM_003998	NFKB1
QG072-01	D04	HQP011309	NM_000250	MPO
QG072-01	D05	HQP011256	NM_004530	MMP2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG072-01	D06	HQP010133	NM_004985	KRAS
QG072-01	D07	HQP009662	NM_000589	IL4
QG072-01	D08	HQP009518	NM_000618	IGF1
QG072-01	D09	HQP009458	NM_000629	IFNAR1
QG072-01	D10	HQP008483	NM_000849	GSTM3
QG072-01	D11	HQP004976	NM_000400	ERCC2
QG072-01	D12	HQP003888	NM_000102	CYP17A1
QG072-01	E01	HQP003814	NM_000106	CYP2D6
QG072-01	E02	HQP003809	NM_000769	CYP2C19
QG072-01	E03	HQP003775	NM_000104	CYP1B1
QG072-01	E04	HQP003499	NM_001037631	CTLA4
QG072-01	E05	HQP002210	NM_000579	CCR5
QG072-01	E06	HQP012021	NM_002542	OGG1
QG072-01	E07	HQP004985	NM_000123	ERCC5
QG072-01	E08	HQP004414	NM_006892	DNMT3B
QG072-01	E09	HQP004317	NM_000903	NQO1
QG072-01	E10	HQP054684	NM_001033	RRM1
QG072-01	E11	HQP053978	NM_001300	KLF6
QG072-01	E12	HQP053916	NM_001076	UGT2B15
QG072-01	F01	HQP023466	NM_004360	CDH1
QG072-01	F02	HQP023284	NM_014805	EPM2AIP1
QG072-01	F03	HQP023237	NM_014779	TSC22D2
QG072-01	F04	HQP023168	NM_004356	CD81
QG072-01	F05	HQP023134	NM_014707	HDAC9
QG072-01	F06	HQP023005	NM_001778	CD48
QG072-01	F07	HQP022764	NM_004832	GSTO1
QG072-01	F08	HQP022722	NM_005191	CD80
QG072-01	F09	HQP022680	NM_004810	GRAP2
QG072-01	F10	HQP022575	NM_130785	TPTE2
QG072-01	F11	HQP022238	NM_004720	EDG4
QG072-01	F12	HQP022063	NM_001037334	USP14
QG072-01	G01	HQP021995	NM_005443	PAPSS1
QG072-01	G02	HQP021865	NM_005679	TAF1C
QG072-01	G03	HQP021754	NM_001759	CCND2
QG072-01	G04	HQP021752	NM_003939	BTRC
QG072-01	G05	HQP021612	NM_003883	HDAC3
QG072-01	G06	HQP020843	NM_032562	PLA2G12B
QG072-01	G07	HQP020394	NM_032019	HDAC10
QG072-01	G08	HQP020331	NM_001013836	MAD1L1
QG072-01	G09	HQP019650	NM_005436	CCDC6
QG072-01	G10	HQP019452	NM_001742	CALCR

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG072-01	G11	HQP018768	NM_001954	DDR1
QG072-01	G12	HQP018564	NM_005432	XRCC3
QG072-01	H01	HGDC		
QG072-01	H02	HGDC		
QG072-01	H03	HQP006940	NM_002046	GAPDH
QG072-01	H04	HQP016381	NM_001101	ACTB
QG072-01	H05	HQP015171	NM_004048	B2M
QG072-01	H06	HQP006171	NM_012423	RPL13A
QG072-01	H07	HQP009026	NM_000194	HPRT1
QG072-01	H08	HQP054253	NR_003286	RN18S1
QG072-01	H09	RT		
QG072-01	H10	RT		
QG072-01	H11	PCR		
QG072-01	H12	PCR		
QG072-02	A01	HQP018474	NM_000376	VDR
QG072-02	A02	HQP018418	NM_001074	UGT2B7
QG072-02	A03	HQP018276	NM_000716	C4BPB
QG072-02	A04	HQP018246	NM_007118	TRIO
QG072-02	A05	HQP018237	NM_004620	TRAF6
QG072-02	A06	HQP018127	NM_003273	TM7SF2
QG072-02	A07	HQP018045	NM_001042454	TGFB111
QG072-02	A08	HQP018016	NM_005652	TERF2
QG072-02	A09	HQP018014	NM_003218	TERF1
QG072-02	A10	HQP017767	NM_003150	STAT3
QG072-02	A11	HQP017764	NM_007315	STAT1
QG072-02	A12	HQP017733	NM_004333	BRAF
QG072-02	B01	HQP017705	NM_004599	SREBF2
QG072-02	B02	HQP017701	NM_005989	AKR1D1
QG072-02	B03	HQP017653	NM_003118	SPARC
QG072-02	B04	HQP017647	NM_003113	SP100
QG072-02	B05	HQP017498	NM_003062	SLIT3
QG072-02	B06	HQP017124	NM_022743	SMYD3
QG072-02	B07	HQP016694	NM_001035511	SDHC
QG072-02	B08	HQP016626	NM_002985	CCL5
QG072-02	B09	HQP016621	NM_002982	CCL2
QG072-02	B10	HQP016561	NM_005622	ACSM3
QG072-02	B11	HQP016284	NM_012421	RLF
QG072-02	B12	HQP016225	NM_002914	RFC2
QG072-02	C01	HQP016212	NM_000657	BCL2
QG072-02	C02	HQP016131	NM_000321	RB1
QG072-02	C03	HQP016087	NM_134424	RAD52

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG072-02	C04	HQP015598	NM_000963	PTGS2
QG072-02	C05	HQP015530	NM_000264	PTCH1
QG072-02	C06	HQP015311	NM_002800	PSMB9
QG072-02	C07	HQP015059	NM_000313	PROS1
QG072-02	C08	HQP014847	NM_006259	PRKG2
QG072-02	C09	HQP013682	NM_017589	BTG4
QG072-02	C10	HQP013616	NM_019093	UGT1A3
QG072-02	C11	HQP013614	NM_007120	UGT1A4
QG072-02	C12	HQP013583	NM_205862	UGT1A6
QG072-02	D01	HQP013579	NM_019075	UGT1A10
QG072-02	D02	HQP013493	NM_000940	PON3
QG072-02	D03	HQP013459	NM_005035	POLRMT
QG072-02	D04	HQP013416	NM_001018111	PODXL
QG072-02	D05	HQP013257	NM_000301	PLG
QG072-02	D06	HQP013240	NM_002661	PLCG2
QG072-02	D07	HQP013181	NM_000298	PKLR
QG072-02	D08	HQP013122	NM_000295	SERPINA1
QG072-02	D09	HQP013100	NM_000927	ABCB1
QG072-02	D10	HQP013088	NM_002631	PGD
QG072-02	D11	HQP012155	NM_002575	SERPINB2
QG072-02	D12	HQP012154	NM_000602	SERPINE1
QG072-02	E01	HQP011866	NM_000625	NOS2A
QG072-02	E02	HQP011808	NM_001077493	NFKB2
QG072-02	E03	HQP011597	NM_002467	MYC
QG072-02	E04	HQP011582	NM_002462	MX1
QG072-02	E05	HQP011555	NM_002454	MTRR
QG072-02	E06	HQP011537	NM_005952	MT1X
QG072-02	E07	HQP011491	NM_000251	MSH2
QG072-02	E08	HQP011325	NM_019899	ABCC1
QG072-02	E09	HQP011266	NM_002426	MMP12
QG072-02	E10	HQP011262	NM_002424	MMP8
QG072-02	E11	HQP011235	NM_000249	MLH1
QG072-02	E12	HQP011211	NM_000246	CIITA
QG072-02	F01	HQP010913	NM_000428	LTBP2
QG072-02	F02	HQP010626	NM_000236	LIPC
QG072-02	F03	HQP010099	NM_000222	KIT
QG072-02	F04	HQP010095	NM_013289	KIR3DL1
QG072-02	F05	HQP010092	NM_012313	KIR2DS3
QG072-02	F06	HQP010089	NM_002255	KIR2DL4
QG072-02	F07	HQP010086	NM_015868	KIR2DL3
QG072-02	F08	HQP010083	NM_014218	KIR2DL1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG072-02	F09	HQP010070	NM_002253	KDR
QG072-02	F10	HQP010005	NM_002239	KCNJ3
QG072-02	F11	HQP009849	NM_002227	JAK1
QG072-02	F12	HQP009808	NM_002210	ITGAV
QG072-02	G01	HQP009803	NM_002207	ITGA9
QG072-02	G02	HQP009749	NM_000207	INS
QG072-02	G03	HQP009746	NM_001565	CXCL10
QG072-02	G04	HQP009726	NM_005536	IMPA1
QG072-02	G05	HQP009694	NM_005535	IL12RB1
QG072-02	G06	HQP009678	NM_000584	IL8
QG072-02	G07	HQP009671	NM_000639	FASLG
QG072-02	G08	HQP009658	NM_000878	IL2RB
QG072-02	G09	HQP009649	NM_000586	IL2
QG072-02	G10	HQP009640	NM_000575	IL1A
QG072-02	G11	HQP009555	NM_002178	IGFBP6
QG072-02	G12	HQP009544	NM_000598	IGFBP3
QG072-02	H01	HGDC		
QG072-02	H02	HGDC		
QG072-02	H03	HQP006940	NM_002046	GAPDH
QG072-02	H04	HQP016381	NM_001101	ACTB
QG072-02	H05	HQP015171	NM_004048	B2M
QG072-02	H06	HQP006171	NM_012423	RPL13A
QG072-02	H07	HQP009026	NM_000194	HPRT1
QG072-02	H08	HQP054253	NR_003286	RN18S1
QG072-02	H09	RT		
QG072-02	H10	RT		
QG072-02	H11	PCR		
QG072-02	H12	PCR		