

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG079-01	A01	HQP009641	NM_000576	IL1B
QG079-01	A02	HQP009645	NM_000577	IL1RN
QG079-01	A03	HQP018141	NM_000594	TNF
QG079-01	A04	HQP011547	NM_005957	MTHFR
QG079-01	A05	HQP009685	NM_000572	IL10
QG079-01	A06	HQP018175	NM_000546	TP53
QG079-01	A07	HQP009678	NM_000584	IL8
QG079-01	A08	HQP023466	NM_004360	CDH1
QG079-01	A09	HQP018562	NM_006297	XRCC1
QG079-01	A10	HQP003817	NM_000773	CYP2E1
QG079-01	A11	HQP018342	NM_001071	TYMS
QG079-01	A12	HQP004976	NM_000400	ERCC2
QG079-01	B01	HQP003772	NM_000499	CYP1A1
QG079-01	B02	HQP018044	NM_000660	TGFB1
QG079-01	B03	HQP009670	NM_000600	IL6
QG079-01	B04	HQP012021	NM_002542	OGG1
QG079-01	B05	HQP015598	NM_000963	PTGS2
QG079-01	B06	HQP009662	NM_000589	IL4
QG079-01	B07	HQP018475	NM_001025366	VEGFA
QG079-01	B08	HQP004975	NM_202001	ERCC1
QG079-01	B09	HQP054047	BC008403	HLA-DRB1
QG079-01	B10	HQP018564	NM_005432	XRCC3
QG079-01	B11	HQP001136	NM_000015	NAT2
QG079-01	B12	HQP011866	NM_000625	NOS2A
QG079-01	C01	HQP011235	NM_000249	MLH1
QG079-01	C02	HQP009640	NM_000575	IL1A
QG079-01	C03	HQP022490	NM_000591	CD14
QG079-01	C04	HQP016204	NM_053056	CCND1
QG079-01	C05	HQP013100	NM_000927	ABCB1
QG079-01	C06	HQP016801	NM_022162	NOD2
QG079-01	C07	HQP011491	NM_000251	MSH2
QG079-01	C08	HQP010133	NM_004985	KRAS
QG079-01	C09	HQP004081	NM_000789	ACE
QG079-01	C10	HQP013633	NM_005037	PPARG
QG079-01	C11	HQP013087	NM_002630	PGC
QG079-01	C12	HQP011554	NM_000254	MTR
QG079-01	D01	HQP011256	NM_004530	MMP2
QG079-01	D02	HQP011135	NM_002392	MDM2
QG079-01	D03	HQP009692	NM_000882	IL12A
QG079-01	D04	HQP009671	NM_000639	FASLG
QG079-01	D05	HQP009664	NM_000418	IL4R

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG079-01	D06	HQP009649	NM_000586	IL2
QG079-01	D07	HQP004439	NM_000110	DPYD
QG079-01	D08	HQP011601	NM_005376	MYCL1
QG079-01	D09	HQP018565	NM_003401	XRCC4
QG079-01	D10	HQP018556	NM_004628	XPC
QG079-01	D11	HQP015596	NM_000962	PTGS1
QG079-01	D12	HQP011555	NM_002454	MTRR
QG079-01	E01	HQP011309	NM_000250	MPO
QG079-01	E02	HQP011263	NM_004994	MMP9
QG079-01	E03	HQP009061	NM_001641	APEX1
QG079-01	E04	HQP004605	NM_005228	EGFR
QG079-01	E05	HQP004599	NM_001963	EGF
QG079-01	E06	HQP000370	NM_058195	CDKN2A
QG079-01	E07	HQP004414	NM_006892	DNMT3B
QG079-01	E08	HQP053978	NM_001300	KLF6
QG079-01	E09	HQP023467	NM_000662	NAT1
QG079-01	E10	HQP018563	NM_005431	XRCC2
QG079-01	E11	HQP017811	NM_177536	SULT1A1
QG079-01	E12	HQP017616	NM_000636	SOD2
QG079-01	F01	HQP017482	NM_194255	SLC19A1
QG079-01	F02	HQP016626	NM_002985	CCL5
QG079-01	F03	HQP013388	NM_017442	TLR9
QG079-01	F04	HQP011687	NM_002485	NBN
QG079-01	F05	HQP011559	NM_001018016	MUC1
QG079-01	F06	HQP010099	NM_000222	KIT
QG079-01	F07	HQP009718	NM_001562	IL18
QG079-01	F08	HQP009717	NM_002190	IL17A
QG079-01	F09	HQP009544	NM_000598	IGFBP3
QG079-01	F10	HQP009469	NM_000416	IFNGR1
QG079-01	F11	HQP009218	NM_000384	APOB
QG079-01	F12	HQP008849	NM_002116	HLA-A
QG079-01	G01	HQP008483	NM_000849	GSTM3
QG079-01	G02	HQP005087	NM_000690	ALDH2
QG079-01	G03	HQP004948	NM_000120	EPHX1
QG079-01	G04	HQP003814	NM_000106	CYP2D6
QG079-01	G05	HQP003171	NM_006437	PARP4
QG079-01	G06	HQP003120	NM_001618	PARP1
QG079-01	G07	HQP002413	NM_000669	ADH1C
QG079-01	G08	HQP001495	NM_052872	IL17F
QG079-01	G09	HQP054686	NM_001123396	LOC729230
QG079-01	G10	HQP054043	NM_203472	SELS

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG079-01	G11	HQP054030	NM_130398	EXO1
QG079-01	G12	HQP053959	NM_001040	SHBG
QG079-01	H01	HGDC		
QG079-01	H02	HGDC		
QG079-01	H03	HQP006940	NM_002046	GAPDH
QG079-01	H04	HQP016381	NM_001101	ACTB
QG079-01	H05	HQP015171	NM_004048	B2M
QG079-01	H06	HQP006171	NM_012423	RPL13A
QG079-01	H07	HQP009026	NM_000194	HPRT1
QG079-01	H08	HQP054253	NR_003286	RN18S1
QG079-01	H09	RT		
QG079-01	H10	RT		
QG079-01	H11	PCR		
QG079-01	H12	PCR		
QG079-02	A01	HQP023414	NM_005847	SLC23A1
QG079-02	A02	HQP023203	NM_001785	CDA
QG079-02	A03	HQP021752	NM_003939	BTRC
QG079-02	A04	HQP021683	NM_021090	MTMR3
QG079-02	A05	HQP021557	NM_003844	TNFRSF10A
QG079-02	A06	HQP021518	NM_000071	CBS
QG079-02	A07	HQP021276	NM_003672	CDC14A
QG079-02	A08	HQP021250	NM_033389	SSH2
QG079-02	A09	HQP020984	NM_032782	HAVCR2
QG079-02	A10	HQP020481	NM_033338	CASP7
QG079-02	A11	HQP020427	NM_001226	CASP6
QG079-02	A12	HQP020297	NM_004346	CASP3
QG079-02	B01	HQP019768	NM_130445	COL18A1
QG079-02	B02	HQP018425	NM_000373	UMPS
QG079-02	B03	HQP018176	NM_005657	TP53BP1
QG079-02	B04	HQP017979	NM_000355	TCN2
QG079-02	B05	HQP017783	NM_006948	STCH
QG079-02	B06	HQP017753	NM_000059	BRCA2
QG079-02	B07	HQP016641	NM_002990	CCL22
QG079-02	B08	HQP016084	NM_014226	RAGE
QG079-02	B09	HQP015828	NM_002827	PTPN1
QG079-02	B10	HQP015565	NM_020661	AICDA
QG079-02	B11	HQP015535	NM_000314	PTEN
QG079-02	B12	HQP015171	NM_004048	B2M
QG079-02	C01	HQP014760	NM_006255	PRKCH
QG079-02	C02	HQP014362	NM_018407	LAPTM4B
QG079-02	C03	HQP013581	NM_019077	UGT1A7

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG079-02	C04	HQP013463	NM_018984	SSH1
QG079-02	C05	HQP013422	NM_002690	POLB
QG079-02	C06	HQP013150	NM_006218	PIK3CA
QG079-02	C07	HQP012501	NM_016341	PLCE1
QG079-02	C08	HQP012448	NM_016161	A4GNT
QG079-02	C09	HQP012393	NM_016009	SH3GLB1
QG079-02	C10	HQP012154	NM_000602	SERPINE1
QG079-02	C11	HQP011810	NM_020529	NFKBIA
QG079-02	C12	HQP011807	NM_003998	NFKB1
QG079-02	D01	HQP011800	NM_006164	NFE2L2
QG079-02	D02	HQP011597	NM_002467	MYC
QG079-02	D03	HQP011549	NM_000252	MTM1
QG079-02	D04	HQP011266	NM_002426	MMP12
QG079-02	D05	HQP011255	NM_002421	MMP1
QG079-02	D06	HQP011219	NM_002415	MIF
QG079-02	D07	HQP010961	NM_005359	SMAD4
QG079-02	D08	HQP010847	NM_000237	LPL
QG079-02	D09	HQP010127	NM_000045	ARG1
QG079-02	D10	HQP009749	NM_000207	INS
QG079-02	D11	HQP009693	NM_002187	IL12B
QG079-02	D12	HQP009681	NM_001557	IL8RB
QG079-02	E01	HQP009679	NM_000634	IL8RA
QG079-02	E02	HQP009677	NM_002185	IL7R
QG079-02	E03	HQP009666	NM_000879	IL5
QG079-02	E04	HQP009539	NM_000596	IGFBP1
QG079-02	E05	HQP009529	NM_000612	IGF2
QG079-02	E06	HQP009523	NM_000875	IGF1R
QG079-02	E07	HQP009472	NM_005534	IFNGR2
QG079-02	E08	HQP009101	NM_001168	BIRC5
QG079-02	E09	HQP009064	NM_000413	HSD17B1
QG079-02	E10	HQP009036	NM_005343	HRAS
QG079-02	E11	HQP009024	NM_000038	APC
QG079-02	E12	HQP008832	NM_181054	HIF1A
QG079-02	F01	HQP008494	NM_015859	GTF2A1
QG079-02	F02	HQP008493	NM_000179	MSH6
QG079-02	F03	HQP008469	NM_002093	GSK3B
QG079-02	F04	HQP008282	NM_002084	GPX3
QG079-02	F05	HQP007755	NM_000516	GNAS
QG079-02	F06	HQP006296	NM_015364	LY96
QG079-02	F07	HQP006144	NM_000802	FOLR1
QG079-02	F08	HQP006129	NM_001014986	FOLH1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG079-02	F09	HQP005280	NM_021642	FCGR2A
QG079-02	F10	HQP005058	NM_000130	F5
QG079-02	F11	HQP005052	NM_000506	F2
QG079-02	F12	HQP004537	NM_001397	ECE1
QG079-02	G01	HQP004525	NM_000686	AGTR2
QG079-02	G02	HQP004494	NM_000029	AGT
QG079-02	G03	HQP004411	NM_022552	DNMT3A
QG079-02	G04	HQP003904	NM_000103	CYP19A1
QG079-02	G05	HQP003888	NM_000102	CYP17A1
QG079-02	G06	HQP003809	NM_000769	CYP2C19
QG079-02	G07	HQP003775	NM_000104	CYP1B1
QG079-02	G08	HQP003539	NM_001904	CTNNB1
QG079-02	G09	HQP003499	NM_001037631	CTLA4
QG079-02	G10	HQP003495	NM_144701	IL23R
QG079-02	G11	HQP003299	NM_004385	VCAN
QG079-02	G12	HQP002671	NM_000754	COMT
QG079-02	H01	HGDC		
QG079-02	H02	HGDC		
QG079-02	H03	HQP006940	NM_002046	GAPDH
QG079-02	H04	HQP016381	NM_001101	ACTB
QG079-02	H05	HQP015171	NM_004048	B2M
QG079-02	H06	HQP006171	NM_012423	RPL13A
QG079-02	H07	HQP009026	NM_000194	HPRT1
QG079-02	H08	HQP054253	NR_003286	RN18S1
QG079-02	H09	RT		
QG079-02	H10	RT		
QG079-02	H11	PCR		
QG079-02	H12	PCR		