

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG020-01	A01	HQP018148	NM_001065	TNFRSF1A
QG020-01	A02	HQP012049	NM_002546	TNFRSF11B
QG020-01	A03	HQP021550	NM_003839	TNFRSF11A
QG020-01	A04	HQP021557	NM_003844	TNFRSF10A
QG020-01	A05	HQP018145	NM_006290	TNFAIP3
QG020-01	A06	HQP018141	NM_000594	TNF
QG020-01	A07	HQP018064	NM_000360	TH
QG020-01	A08	HQP018051	NM_004612	TGFBR1
QG020-01	A09	HQP018047	NM_003238	TGFB2
QG020-01	A10	HQP018044	NM_000660	TGFB1
QG020-01	A11	HQP018043	NM_003236	TGFA
QG020-01	A12	HQP018032	NM_007111	TFDP1
QG020-01	B01	HQP014691	NM_001008744	TDP1
QG020-01	B02	HQP017996	NM_003211	TDG
QG020-01	B03	HQP017984	NM_001008897	TCP1
QG020-01	B04	HQP017899	NM_000593	TAP1
QG020-01	B05	HQP017784	NM_005420	SULT1E1
QG020-01	B06	HQP017810	NM_001055	SULT1A1
QG020-01	B07	HQP017771	NM_003152	STAT5A
QG020-01	B08	HQP017698	NM_000348	SRD5A2
QG020-01	B09	HQP017616	NM_000636	SOD2
QG020-01	B10	HQP017615	NM_000454	SOD1
QG020-01	B11	HQP013340	NM_017435	SLCO1C1
QG020-01	B12	HQP017350	NM_006516	SLC2A1
QG020-01	C01	HQP001134	NM_005094	SLC27A4
QG020-01	C02	HQP010011	NM_198580	SLC27A1
QG020-01	C03	HQP017462	NM_000452	SLC10A2
QG020-01	C04	HQP017461	NM_003049	SLC10A1
QG020-01	C05	HQP006080	NM_012238	SIRT1
QG020-01	C06	HQP017098	NM_000193	SHH
QG020-01	C07	HQP012154	NM_000602	SERPINE1
QG020-01	C08	HQP016689	NM_003000	SDHB
QG020-01	C09	HQP016613	NM_001007098	SCP2
QG020-01	C10	HQP016583	NM_005063	SCD
QG020-01	C11	HQP016555	NM_000331	SAA1
QG020-01	C12	HQP016527	NM_021976	RXRΒ
QG020-01	D01	HQP016526	NM_002957	RXRA
QG020-01	D02	HQP016214	NM_006509	RELB
QG020-01	D03	HQP016131	NM_000321	RB1
QG020-01	D04	HQP016114	NM_000964	RARA
QG020-01	D05	HQP016089	NM_000448	RAG1

QG020-01	D06	HQP016088	NM_002880	RAF1
QG020-01	D07	HQP016087	NM_134424	RAD52
QG020-01	D08	HQP016077	NM_002875	RAD51
QG020-01	D09	HQP000145	NM_005732	RAD50
QG020-01	D10	HQP016075	NM_005053	RAD23A
QG020-01	D11	HQP016074	NM_006265	RAD21
QG020-01	D12	HQP015190	NM_020165	RAD18
QG020-01	E01	HQP015947	NM_002853	RAD1
QG020-01	E02	HQP016063	NM_006908	RAC1
QG020-01	E03	HQP015598	NM_000963	PTGS2
QG020-01	E04	HQP015596	NM_000962	PTGS1
QG020-01	E05	HQP015535	NM_000314	PTEN
QG020-01	E06	HQP015530	NM_000264	PTCH1
QG020-01	E07	HQP012152	NM_002574	PRDX1
QG020-01	E08	HQP001016	NM_013261	PPARGC1A
QG020-01	E09	HQP013633	NM_005037	PPARG
QG020-01	E10	HQP013627	NM_006238	PPARD
QG020-01	E11	HQP013504	NM_000941	POR
QG020-01	E12	HQP013493	NM_000940	PON3
QG020-01	F01	HQP007689	NM_013274	POLL
QG020-01	F02	HQP012725	NM_016218	POLK
QG020-01	F03	HQP001399	NM_007195	POLI
QG020-01	F04	HQP013422	NM_002690	POLB
QG020-01	F05	HQP013352	NM_000535	PMS2
QG020-01	F06	HQP013280	NM_006227	PLTP
QG020-01	F07	HQP013194	NM_000300	PLA2G2A
QG020-01	F08	HQP013181	NM_000298	PKLR
QG020-01	F09	HQP013099	NM_000926	PGR
QG020-01	F10	HQP016753	NM_022121	PERP
QG020-01	F11	HQP012420	NM_002592	PCNA
QG020-01	F12	HQP012078	NM_000607	ORM1
QG020-01	G01	HQP011914	NM_002524	NRAS
QG020-01	G02	HQP021630	NM_003889	NR1I2
QG020-01	G03	HQP004317	NM_000903	NQO1
QG020-01	G04	HQP011868	NM_000603	NOS3
QG020-01	G05	HQP011866	NM_000625	NOS2A
QG020-01	G06	HQP011860	NM_006169	NNMT
QG020-01	G07	HQP011807	NM_003998	NFKB1
QG020-01	G08	HQP011774	NM_000267	NF1
QG020-01	G09	HQP011730	NM_007103	NDUFV1
QG020-01	G10	HQP011731	NM_002495	NDUFS4
QG020-01	G11	HQP011726	NM_005006	NDUFS1
QG020-01	G12	HQP001136	NM_000015	NAT2

QG020-01	H01	HGDC		
QG020-01	H02	HGDC		
QG020-01	H03	HQP006940	NM_002046	GAPDH
QG020-01	H04	HQP016381	NM_001101	ACTB
QG020-01	H05	HQP015171	NM_004048	B2M
QG020-01	H06	HQP006171	NM_012423	RPL13A
QG020-01	H07	HQP009026	NM_000194	HPRT1
QG020-01	H08	HQP054253	NR_003286	RN18S1
QG020-01	H09	RT		
QG020-01	H10	RT		
QG020-01	H11	PCR		
QG020-01	H12	PCR		
QG020-02	A01	HQP023467	NM_000662	NAT1
QG020-02	A02	HQP011597	NM_002467	MYC
QG020-02	A03	HQP023412	NM_005115	MVP
QG020-02	A04	HQP011553	NM_000253	MTTP
QG020-02	A05	HQP011539	NM_005954	MT3
QG020-02	A06	HQP011538	NM_005953	MT2A
QG020-02	A07	HQP008493	NM_000179	MSH6
QG020-02	A08	HQP011491	NM_000251	MSH2
QG020-02	A09	HQP011309	NM_000250	MPO
QG020-02	A10	HQP011304	NM_001015052	MPG
QG020-02	A11	HQP011263	NM_004994	MMP9
QG020-02	A12	HQP011256	NM_004530	MMP2
QG020-02	B01	HQP011241	NM_005933	MLL
QG020-02	B02	HQP011235	NM_000249	MLH1
QG020-02	B03	HQP011219	NM_002415	MIF
QG020-02	B04	HQP011208	NM_020300	MGST1
QG020-02	B05	HQP011181	NM_000245	MET
QG020-02	B06	HQP011135	NM_002392	MDM2
QG020-02	B07	HQP011104	NM_021960	MCL1
QG020-02	B08	HQP011077	NM_000242	MBL2
QG020-02	B09	HQP021736	NM_003925	MBD4
QG020-02	B10	HQP011071	NM_005368	MB
QG020-02	B11	HQP014886	NM_002750	MAPK8
QG020-02	B12	HQP014848	NM_002745	MAPK1
QG020-02	C01	HQP011008	NM_000898	MAOB
QG020-02	C02	HQP011007	NM_000240	MAOA
QG020-02	C03	HQP010915	NM_002342	LTBR
QG020-02	C04	HQP010907	NM_000595	LTA
QG020-02	C05	HQP010613	NM_002312	LIG4
QG020-02	C06	HQP010609	NM_000234	LIG1
QG020-02	C07	HQP010577	NM_000527	LDLR

QG020-02	C08	HQP010560	NM_004139	LBP
QG020-02	C09	HQP009854	NM_002229	JUNB
QG020-02	C10	HQP009853	NM_002228	JUN
QG020-02	C11	HQP009845	NM_002225	IVD
QG020-02	C12	HQP009815	NM_000211	ITGB2
QG020-02	D01	HQP009806	NM_002209	ITGAL
QG020-02	D02	HQP009749	NM_000207	INS
QG020-02	D03	HQP009670	NM_000600	IL6
QG020-02	D04	HQP009641	NM_000576	IL1B
QG020-02	D05	HQP009640	NM_000575	IL1A
QG020-02	D06	HQP009718	NM_001562	IL18
QG020-02	D07	HQP009555	NM_002178	IGFBP6
QG020-02	D08	HQP009544	NM_000598	IGFBP3
QG020-02	D09	HQP009541	NM_000597	IGFBP2
QG020-02	D10	HQP009539	NM_000596	IGFBP1
QG020-02	D11	HQP009532	NM_000876	IGF2R
QG020-02	D12	HQP009523	NM_000875	IGF1R
QG020-02	E01	HQP009518	NM_000618	IGF1
QG020-02	E02	HQP000633	NM_006389	HYOU1
QG020-02	E03	HQP009138	NM_004507	HUS1
QG020-02	E04	HQP009098	NM_002156	HSPD1
QG020-02	E05	HQP009090	NM_001541	HSPB2
QG020-02	E06	HQP009089	NM_001540	HSPB1
QG020-02	E07	HQP009088	NM_004134	HSPA9
QG020-02	E08	HQP009083	NM_005347	HSPA5
QG020-02	E09	HQP009080	NM_021979	HSPA2
QG020-02	E10	HQP009079	NM_005527	HSPA1L
QG020-02	E11	HQP009078	NM_005346	HSPA1B
QG020-02	E12	HQP009077	NM_005345	HSPA1A
QG020-02	F01	HQP009092	NM_001017963	HSP90AA1
QG020-02	F02	HQP009068	NM_005526	HSF1
QG020-02	F03	HQP009035	NM_000613	HPX
QG020-02	F04	HQP009026	NM_000194	HPRT1
QG020-02	F05	HQP008898	NM_002133	HMOX1
QG020-02	F06	HQP008831	NM_001530	HIF1A
QG020-02	F07	HQP008800	NM_000601	HGF
QG020-02	F08	HQP004493	NM_001945	HBEGF
QG020-02	F09	HQP013447	NM_017545	HAO1
QG020-02	F10	HQP008502	NM_005316	GTF2H1
QG020-02	F11	HQP008479	NM_001512	GSTA4
QG020-02	F12	HQP008473	NM_000637	GSR
QG020-02	G01	HQP008469	NM_002093	GSK3B
QG020-02	G02	HQP008288	NM_001509	GPX5

QG020-02	G03	HQP008281	NM_002083	GPX2
QG020-02	G04	HQP008279	NM_000581	GPX1
QG020-02	G05	HQP007813	NM_000408	GPD2
QG020-02	G06	HQP006930	NM_002045	GAP43
QG020-02	G07	HQP004125	NM_001924	GADD45A
QG020-02	G08	HQP011894	NM_001680	FXVD2
QG020-02	G09	HQP005759	NM_001455	FOXO3A
QG020-02	G10	HQP005712	NM_021953	FOXM1
QG020-02	G11	HQP006188	NM_005252	FOS
QG020-02	G12	HQP005968	NM_002022	FMO4
QG020-02	H01	HGDC		
QG020-02	H02	HGDC		
QG020-02	H03	HQP006940	NM_002046	GAPDH
QG020-02	H04	HQP016381	NM_001101	ACTB
QG020-02	H05	HQP015171	NM_004048	B2M
QG020-02	H06	HQP006171	NM_012423	RPL13A
QG020-02	H07	HQP009026	NM_000194	HPRT1
QG020-02	H08	HQP054253	NR_003286	RN18S1
QG020-02	H09	RT		
QG020-02	H10	RT		
QG020-02	H11	PCR		
QG020-02	H12	PCR		
QG020-03	A01	HQP005959	NM_001002294	FMO3
QG020-03	A02	HQP005939	NM_002021	FMO1
QG020-03	A03	HQP005437	NM_000141	FGFR2
QG020-03	A04	HQP005403	NM_002006	FGF2
QG020-03	A05	HQP005400	NM_000800	FGF1
QG020-03	A06	HQP005391	NM_004111	FEN1
QG020-03	A07	HQP005387	NM_004110	FDXR
QG020-03	A08	HQP009671	NM_000639	FASLG
QG020-03	A09	HQP009651	NM_000043	FAS
QG020-03	A10	HQP005083	NM_000136	FANCC
QG020-03	A11	HQP005081	NM_000135	FANCA
QG020-03	A12	HQP021526	NM_003824	FADD
QG020-03	B01	HQP005080	NM_001446	FABP7
QG020-03	B02	HQP005073	NM_001443	FABP1
QG020-03	B03	HQP005001	NM_001040275	ESR2
QG020-03	B04	HQP004998	NM_000125	ESR1
QG020-03	B05	HQP004983	NM_000122	ERCC3
QG020-03	B06	HQP004976	NM_000400	ERCC2
QG020-03	B07	HQP004974	NM_001983	ERCC1
QG020-03	B08	HQP004970	NM_001005915	ERBB3
QG020-03	B09	HQP004968	NM_001005862	ERBB2

QG020-03	B10	HQP020089	NM_000502	EPX
QG020-03	B11	HQP004948	NM_000120	EPHX1
QG020-03	B12	HQP004749	NM_005229	ELK1
QG020-03	C01	HQP004612	NM_001964	EGR1
QG020-03	C02	HQP004605	NM_005228	EGFR
QG020-03	C03	HQP004599	NM_001963	EGF
QG020-03	C04	HQP004564	NM_000115	EDNRB
QG020-03	C05	HQP004557	NM_001955	EDN1
QG020-03	C06	HQP004528	NM_001950	E2F4
QG020-03	C07	HQP004524	NM_005225	E2F1
QG020-03	C08	HQP004439	NM_000110	DPYD
QG020-03	C09	HQP009103	NM_006145	DNAJB1
QG020-03	C10	HQP009076	NM_001539	DNAJA1
QG020-03	C11	HQP015116	NM_019887	DIABLO
QG020-03	C12	HQP004309	NM_000791	DHFR
QG020-03	D01	HQP004127	NM_004083	DDIT3
QG020-03	D02	HQP004115	NM_000790	DDC
QG020-03	D03	HQP004114	NM_000107	DDB2
QG020-03	D04	HQP016982	NM_001033855	DCLRE1C
QG020-03	D05	HQP003975	NM_004394	DAP
QG020-03	D06	HQP003865	NM_004391	CYP8B1
QG020-03	D07	HQP022723	NM_004820	CYP7B1
QG020-03	D08	HQP003859	NM_000780	CYP7A1
QG020-03	D09	HQP003924	NM_000786	CYP51A1
QG020-03	D10	HQP003855	NM_000779	CYP4B1
QG020-03	D11	HQP003847	NM_000778	CYP4A11
QG020-03	D12	HQP000977	NM_006668	CYP46A1
QG020-03	E01	HQP003841	NM_000777	CYP3A5
QG020-03	E02	HQP003836	NM_017460	CYP3A4
QG020-03	E03	HQP008529	NM_030622	CYP2S1
QG020-03	E04	HQP003818	NM_000774	CYP2F1
QG020-03	E05	HQP003817	NM_000773	CYP2E1
QG020-03	E06	HQP003814	NM_000106	CYP2D6
QG020-03	E07	HQP003811	NM_000771	CYP2C9
QG020-03	E08	HQP003808	NM_000767	CYP2B6
QG020-03	E09	HQP003804	NM_000766	CYP2A13
QG020-03	E10	HQP003923	NM_000785	CYP27B1
QG020-03	E11	HQP003922	NM_000784	CYP27A1
QG020-03	E12	HQP003919	NM_000783	CYP26A1
QG020-03	F01	HQP003917	NM_000782	CYP24A1
QG020-03	F02	HQP003775	NM_000104	CYP1B1
QG020-03	F03	HQP003774	NM_000761	CYP1A2
QG020-03	F04	HQP003772	NM_000499	CYP1A1

QG020-03	F05	HQP003904	NM_000103	CYP19A1
QG020-03	F06	HQP003888	NM_000102	CYP17A1
QG020-03	F07	HQP003871	NM_000781	CYP11A1
QG020-03	F08	HQP004315	NM_000398	CYB5R3
QG020-03	F09	HQP009746	NM_001565	CXCL10
QG020-03	F10	HQP003572	NM_001905	CTPS
QG020-03	F11	HQP003539	NM_001904	CTNNB1
QG020-03	F12	HQP003218	NM_004383	CSK
QG020-03	G01	HQP003159	NM_000758	CSF2
QG020-03	G02	HQP002867	NM_001031847	CPT1A
QG020-03	G03	HQP002820	NM_000096	CP
QG020-03	G04	HQP002671	NM_000754	COMT
QG020-03	G05	HQP002462	NM_000088	COL1A1
QG020-03	G06	HQP002037	NM_001831	CLU
QG020-03	G07	HQP001396	NM_001005735	CHEK2
QG020-03	G08	HQP001182	NM_020549	CHAT
QG020-03	G09	HQP000815	NM_001025194	CES1
QG020-03	G10	HQP000369	NM_000077	CDKN2A
QG020-03	G11	HQP000342	NM_004064	CDKN1B
QG020-03	G12	HQP000331	NM_000389	CDKN1A
QG020-03	H01	HGDC		
QG020-03	H02	HGDC		
QG020-03	H03	HQP006940	NM_002046	GAPDH
QG020-03	H04	HQP016381	NM_001101	ACTB
QG020-03	H05	HQP015171	NM_004048	B2M
QG020-03	H06	HQP006171	NM_012423	RPL13A
QG020-03	H07	HQP009026	NM_000194	HPRT1
QG020-03	H08	HQP054253	NR_003286	RN18S1
QG020-03	H09	RT		
QG020-03	H10	RT		
QG020-03	H11	PCR		
QG020-03	H12	PCR		
QG020-04	A01	HQP000245	NM_000075	CDK4
QG020-04	A02	HQP000225	NM_001798	CDK2
QG020-04	A03	HQP023354	NM_001254	CDC6
QG020-04	A04	HQP001310	NM_007065	CDC37
QG020-04	A05	HQP023409	NM_001790	CDC25C
QG020-04	A06	HQP023399	NM_004358	CDC25B
QG020-04	A07	HQP023385	NM_001789	CDC25A
QG020-04	A08	HQP023365	NM_001255	CDC20
QG020-04	A09	HQP022955	NM_001250	CD40
QG020-04	A10	HQP022821	NM_000072	CD36
QG020-04	A11	HQP021882	NM_004354	CCNG2

QG020-04	A12	HQP021819	NM_001238	CCNE1
QG020-04	B01	HQP021754	NM_001759	CCND2
QG020-04	B02	HQP016204	NM_053056	CCND1
QG020-04	B03	HQP021734	NM_001013399	CCNC
QG020-04	B04	HQP021727	NM_031966	CCNB1
QG020-04	B05	HQP021701	NM_001237	CCNA2
QG020-04	B06	HQP016625	NM_002984	CCL4
QG020-04	B07	HQP020648	NM_001229	CASP9
QG020-04	B08	HQP018966	NM_001080124	CASP8
QG020-04	B09	HQP020297	NM_004346	CASP3
QG020-04	B10	HQP017753	NM_000059	BRCA2
QG020-04	B11	HQP017713	NM_007294	BRCA1
QG020-04	B12	HQP017642	NM_032515	BOK
QG020-04	C01	HQP017631	NM_004331	BNIP3L
QG020-04	C02	HQP017619	NM_004052	BNIP3
QG020-04	C03	HQP016844	NM_000057	BLM
QG020-04	C04	HQP009099	NM_001012270	BIRC5
QG020-04	C05	HQP016653	NM_001196	BID
QG020-04	C06	HQP016418	NM_004327	BCR
QG020-04	C07	HQP016328	NM_001706	BCL6
QG020-04	C08	HQP016237	NM_001191	BCL2L1
QG020-04	C09	HQP016211	NM_000633	BCL2
QG020-04	C10	HQP016108	NM_000055	BCHE
QG020-04	C11	HQP015964	NM_004324	BAX
QG020-04	C12	HQP015917	NM_001188	BAK1
QG020-04	D01	HQP015574	NM_004323	BAG1
QG020-04	D02	HQP015538	NM_004322	BAD
QG020-04	D03	HQP011898	NM_004320	ATP2A1
QG020-04	D04	HQP011736	NM_000051	ATM
QG020-04	D05	HQP010924	NM_001668	ARNT
QG020-04	D06	HQP009801	NM_000044	AR
QG020-04	D07	HQP009578	NM_000484	APP
QG020-04	D08	HQP009556	NM_000041	APOE
QG020-04	D09	HQP009447	NM_000483	APOC2
QG020-04	D10	HQP009125	NM_000039	APOA1
QG020-04	D11	HQP009061	NM_001641	APEX1
QG020-04	D12	HQP009024	NM_000038	APC
QG020-04	E01	HQP008934	NM_001160	APAF1
QG020-04	E02	HQP021378	NM_003734	AOC3
QG020-04	E03	HQP008829	NM_001154	ANXA5
QG020-04	E04	HQP006359	NM_000698	ALOX5
QG020-04	E05	HQP005075	NM_000689	ALDH1A1
QG020-04	E06	HQP004991	NM_001014431	AKT1

QG020-04	E07	HQP004658	NM_001621	AHR
QG020-04	E08	HQP002234	NM_001122	ADFP
QG020-04	E09	HQP010411	NM_000019	ACAT1
QG020-04	E10	HQP007422	NM_001091	ABP1
QG020-04	E11	HQP006954	NM_005157	ABL1
QG020-04	E12	HQP016876	NM_022437	ABCG8
QG020-04	F01	HQP016875	NM_022436	ABCG5
QG020-04	F02	HQP022745	NM_004827	ABCG2
QG020-04	F03	HQP015971	NM_002858	ABCD3
QG020-04	F04	HQP005426	NM_005164	ABCD2
QG020-04	F05	HQP005065	NM_000033	ABCD1
QG020-04	F06	HQP009816	NM_001079528	ABCC6
QG020-04	F07	HQP000074	NM_001023587	ABCC5
QG020-04	F08	HQP021459	NM_003786	ABCC3
QG020-04	F09	HQP002260	NM_000392	ABCC2
QG020-04	F10	HQP011322	NM_004996	ABCC1
QG020-04	F11	HQP013101	NM_000443	ABCB4
QG020-04	F12	HQP021388	NM_003742	ABCB11
QG020-04	G01	HQP013100	NM_000927	ABCB1
QG020-04	G02	HQP004727	NM_005502	ABCA1
QG020-04	G03	HQP006558	NM_002039	GAB1
QG020-04	G04	HQP011604	NM_015675	GADD45B
QG020-04	G05	HQP007231	NM_000159	GCDH
QG020-04	G06	HQP011720	NM_004146	NDUFB7
QG020-04	G07	HQP011937	NM_002528	NTHL1
QG020-04	G08	HQP022046	NM_004203	PKMYT1
QG020-04	G09	HQP013296	NM_021127	PMAIP1
QG020-04	G10	HQP009626	NM_000485	APRT
QG020-04	G11	HQP022987	NM_004915	ABCG1
QG020-04	G12	HQP009833	NM_001609	ACADSB
QG020-04	H01	HGDC		
QG020-04	H02	HGDC		
QG020-04	H03	HQP006940	NM_002046	GAPDH
QG020-04	H04	HQP016381	NM_001101	ACTB
QG020-04	H05	HQP015171	NM_004048	B2M
QG020-04	H06	HQP006171	NM_012423	RPL13A
QG020-04	H07	HQP009026	NM_000194	HPRT1
QG020-04	H08	HQP054253	NR_003286	RN18S1
QG020-04	H09	RT		
QG020-04	H10	RT		
QG020-04	H11	PCR		
QG020-04	H12	PCR		