

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG040-01	A01	HQP013618	NM_032644	PPARA
QG040-01	A02	HQP001030	NM_006696	BRD8
QG040-01	A03	HQP009802	NM_001011645	AR
QG040-01	A04	HQP009026	NM_000194	HPRT1
QG040-01	A05	HQP018228	NM_001032287	NR2C1
QG040-01	A06	HQP018474	NM_000376	VDR
QG040-01	A07	HQP022535	NM_016213	TRIP4
QG040-01	A08	HQP018079	NM_000461	THRB
QG040-01	A09	HQP000044	NM_005481	THRAP5
QG040-01	A10	HQP023303	NM_014815	THRAP4
QG040-01	A11	HQP023420	NM_005121	THRAP1
QG040-01	A12	HQP018077	NM_003250	THRA
QG040-01	B01	HQP016527	NM_021976	RXRΒ
QG040-01	B02	HQP016526	NM_002957	RXRA
QG040-01	B03	HQP009574	NM_005349	RBPJ
QG040-01	B04	HQP016118	NM_000966	RARG
QG040-01	B05	HQP016116	NM_000965	RARB
QG040-01	B06	HQP016114	NM_000964	RARA
QG040-01	B07	HQP015346	NM_002804	PSMC3
QG040-01	B08	HQP002750	NM_133263	PPARGC1B
QG040-01	B09	HQP001016	NM_013261	PPARGC1A
QG040-01	B10	HQP013634	NM_015869	PPARG
QG040-01	B11	HQP013627	NM_006238	PPARD
QG040-01	B12	HQP013635	NM_004774	PPARBP
QG040-01	C01	HQP054001	NM_005036	PPARA
QG040-01	C02	HQP021621	NM_003884	PCAF
QG040-01	C03	HQP020042	NM_003489	NRIP1
QG040-01	C04	HQP008900	NM_002135	NR4A1
QG040-01	C05	HQP011251	NM_000901	NR3C2
QG040-01	C06	HQP008401	NM_000176	NR3C1
QG040-01	C07	HQP004966	NM_005234	NR2F6
QG040-01	C08	HQP018031	NM_021005	NR2F2
QG040-01	C09	HQP018230	NM_003298	NR2C2
QG040-01	C10	HQP021631	NM_022002	NR1I2
QG040-01	C11	HQP023437	NM_005123	NR1H4
QG040-01	C12	HQP000086	NM_005693	NR1H3
QG040-01	D01	HQP018431	NM_007121	NR1H2
QG040-01	D02	HQP023440	NM_005126	NR1D2
QG040-01	D03	HQP022925	NM_021724	NR1D1
QG040-01	D04	HQP020670	NM_021969	NR0B2
QG040-01	D05	HQP011875	NM_024408	NOTCH2

QG040-01	D06	HQP011864	NM_007363	NONO
QG040-01	D07	HQP053985	NM_002502	NFKB2
QG040-01	D08	HQP022978	NM_006311	NCOR1
QG040-01	D09	HQP023063	NM_024831	TGS1
QG040-01	D10	HQP019661	NM_005437	NCOA4
QG040-01	D11	HQP000603	NM_006540	NCOA2
QG040-01	D12	HQP022096	NM_004689	MTA1
QG040-01	E01	HQP008386	NM_014166	MED4
QG040-01	E02	HQP023419	NM_005120	MED12
QG040-01	E03	HQP000630	NM_006388	HTATIP
QG040-01	E04	HQP012862	NM_016596	HDAC7A
QG040-01	E05	HQP000022	NM_006044	HDAC6
QG040-01	E06	HQP000024	NM_005474	HDAC5
QG040-01	E07	HQP023167	NM_006037	HDAC4
QG040-01	E08	HQP021612	NM_003883	HDAC3
QG040-01	E09	HQP008746	NM_001527	HDAC2
QG040-01	E10	HQP008745	NM_004964	HDAC1
QG040-01	E11	HQP005005	NM_001438	ESRRG
QG040-01	E12	HQP005003	NM_004451	ESRRA
QG040-01	F01	HQP005002	NM_001437	ESR2
QG040-01	F02	HQP004998	NM_000125	ESR1
QG040-01	F03	HQP004144	NM_004396	DDX5
QG040-01	F04	HQP022755	NM_004268	CRSP6
QG040-01	F05	HQP022465	NM_004229	CRSP2
QG040-01	F06	HQP002921	NM_004380	CREBBP
QG040-01	F07	HQP022524	NM_004236	COPS2
QG040-01	F08	HQP010924	NM_001668	ARNT
QG040-01	F09	HQP009801	NM_000044	AR
QG040-01	F10	HQP004658	NM_001621	AHR
QG040-01	F11	HQP004563	NM_000475	NR0B1
QG040-01	F12	HQP016374	NM_134260	RORA
QG040-01	G01	HQP016529	NM_006917	RXRG
QG040-01	G02	HQP001032	NM_183359	BRD8
QG040-01	G03	HQP005004	NM_004452	ESRRB
QG040-01	G04	HQP054033	NM_145904	HMGA1
QG040-01	G05	HQP008912	NM_178849	HNF4A
QG040-01	G06	HQP015364	NM_002805	PSMC5
QG040-01	G07	HQP021389	NM_003743	NCOA1
QG040-01	G08	HQP020041	NM_181659	NCOA3
QG040-01	G09	HQP023436	NM_005122	NR1I3
QG040-01	G10	HQP018229	NM_003297	NR2C1
QG040-01	G11	HQP000004	NM_014249	NR2E3
QG040-01	G12	HQP018030	NM_005654	NR2F1

QG040-01	H01	HGDC		
QG040-01	H02	HGDC		
QG040-01	H03	HQP006940	NM_002046	GAPDH
QG040-01	H04	HQP016381	NM_001101	ACTB
QG040-01	H05	HQP015171	NM_004048	B2M
QG040-01	H06	HQP006171	NM_012423	RPL13A
QG040-01	H07	HQP009026	NM_000194	HPRT1
QG040-01	H08	HQP054253	NR_003286	RN18S1
QG040-01	H09	RT		
QG040-01	H10	RT		
QG040-01	H11	PCR		
QG040-01	H12	PCR		