

Plate	Position	Catalog No. of	Accession No. of	Symbol
		Primer	Gene	
QG092-01	A01	HQP021631	NM_022002	NR1I2
QG092-01	A02	HQP018272	NM_001018036	TSHR
QG092-01	A03	HQP016379	NM_005060	RORC
QG092-01	A04	HQP016375	NM_134261	RORA
QG092-01	A05	HQP016117	NM_016152	RARB
QG092-01	A06	HQP013634	NM_015869	PPARG
QG092-01	A07	HQP008402	NM_001020825	NR3C1
QG092-01	A08	HQP006438	NM_205860	NR5A2
QG092-01	A09	HQP005002	NM_001437	ESR2
QG092-01	A10	HQP009802	NM_001011645	AR
QG092-01	A11	HQP001030	NM_006696	BRD8
QG092-01	A12	HQP000508	NM_006320	PGRMC2
QG092-01	B01	HQP000086	NM_005693	NR1H3
QG092-01	B02	HQP023440	NM_005126	NR1D2
QG092-01	B03	HQP023437	NM_005123	NR1H4
QG092-01	B04	HQP022925	NM_021724	NR1D1
QG092-01	B05	HQP021630	NM_003889	NR1I2
QG092-01	B06	HQP020670	NM_021969	NR0B2
QG092-01	B07	HQP019536	NM_006981	NR4A3
QG092-01	B08	HQP018474	NM_000376	VDR
QG092-01	B09	HQP018431	NM_007121	NR1H2
QG092-01	B10	HQP018271	NM_000369	TSHR
QG092-01	B11	HQP018242	NM_003301	TRHR
QG092-01	B12	HQP018241	NM_007117	TRH
QG092-01	C01	HQP018230	NM_003298	NR2C2
QG092-01	C02	HQP018228	NM_001032287	NR2C1
QG092-01	C03	HQP018079	NM_000461	THRB
QG092-01	C04	HQP018077	NM_003250	THRA
QG092-01	C05	HQP018031	NM_021005	NR2F2
QG092-01	C06	HQP016527	NM_021976	RXRB
QG092-01	C07	HQP016526	NM_002957	RXRA
QG092-01	C08	HQP016378	NM_001001523	RORC
QG092-01	C09	HQP016377	NM_006914	RORB
QG092-01	C10	HQP016118	NM_000966	RARG
QG092-01	C11	HQP016116	NM_000965	RARB
QG092-01	C12	HQP016114	NM_000964	RARA
QG092-01	D01	HQP015552	NM_000957	PTGER3
QG092-01	D02	HQP013633	NM_005037	PPARG
QG092-01	D03	HQP013627	NM_006238	PPARD
QG092-01	D04	HQP013617	NM_001001928	PPARA
QG092-01	D05	HQP013099	NM_000926	PGR

QG092-01	D06	HQP011968	NM_006186	NR4A2
QG092-01	D07	HQP011891	NM_001037738	NPM1
QG092-01	D08	HQP011251	NM_000901	NR3C2
QG092-01	D09	HQP008914	NM_004133	HNF4G
QG092-01	D10	HQP008908	NM_000457	HNF4A
QG092-01	D11	HQP008900	NM_002135	NR4A1
QG092-01	D12	HQP008401	NM_000176	NR3C1
QG092-01	E01	HQP007781	NM_000406	GNRHR
QG092-01	E02	HQP007395	NM_000163	GHR
QG092-01	E03	HQP005005	NM_001438	ESRRG
QG092-01	E04	HQP005003	NM_004451	ESRRA
QG092-01	E05	HQP005001	NM_001040275	ESR2
QG092-01	E06	HQP004998	NM_000125	ESR1
QG092-01	E07	HQP004966	NM_005234	NR2F6
QG092-01	E08	HQP004829	NM_001974	EMR1
QG092-01	E09	HQP009801	NM_000044	AR
QG092-01	E10	HQP004658	NM_001621	AHR
QG092-01	E11	HQP004563	NM_000475	NR0B1
QG092-01	E12	HQP002971	NM_004382	CRHR1
QG092-01	F01	HQP007399	NM_004122	GHSR
QG092-01	F02	HQP016373	NM_002943	RORA
QG092-01	F03	HQP018119	NM_003269	NR2E1
QG092-01	F04	HQP016529	NM_006917	RXRG
QG092-01	F05	HQP005004	NM_004452	ESRRB
QG092-01	F06	HQP006437	NM_003822	NR5A2
QG092-01	F07	HQP054190	NM_001130101	NR1H3
QG092-01	F08	HQP007259	NM_001489	NR6A1
QG092-01	F09	HQP016528	NM_001009598	RXRG
QG092-01	F10	HQP018030	NM_005654	NR2F1
QG092-01	F11	HQP023422	NM_001077469	NR1I3
QG092-01	F12	HQP000004	NM_014249	NR2E3
QG092-01	G01	HQP000005	NM_016346	NR2E3
QG092-01	G02	HQP054187	NM_001134285	ESRRG
QG092-01	G03	HQP007398	NM_001009824	GHRHR
QG092-01	G04	HQP007400	NM_198407	GHSR
QG092-01	G05	HQP007782	NM_001012763	GNRHR
QG092-01	G06	HQP008909	NM_001030003	HNF4A
QG092-01	G07	HQP011892	NM_002520	NPM1
QG092-01	G08	HQP023423	NM_001077470	NR1I3
QG092-01	G09	HQP054189	NM_001145425	NR1D2
QG092-01	G10	HQP018078	NM_199334	THRA
QG092-01	G11	HQP018229	NM_003297	NR2C1
QG092-01	G12	HQP019537	NM_173199	NR4A3

QG092-01	H01	HGDC		
QG092-01	H02	HGDC		
QG092-01	H03	HQP006940	NM_002046	GAPDH
QG092-01	H04	HQP016381	NM_001101	ACTB
QG092-01	H05	HQP015171	NM_004048	B2M
QG092-01	H06	HQP006171	NM_012423	RPL13A
QG092-01	H07	HQP009026	NM_000194	HPRT1
QG092-01	H08	HQP054253	NR_003286	RN18S1
QG092-01	H09	RT		
QG092-01	H10	RT		
QG092-01	H11	PCR		
QG092-01	H12	PCR		