

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG007-01	A01	HQP013100	NM_000927	ABCB1
QG007-01	A02	HQP011322	NM_004996	ABCC1
QG007-01	A03	HQP002260	NM_000392	ABCC2
QG007-01	A04	HQP021459	NM_003786	ABCC3
QG007-01	A05	HQP000075	NM_005688	ABCC5
QG007-01	A06	HQP009817	NM_001171	ABCC6
QG007-01	A07	HQP022745	NM_004827	ABCG2
QG007-01	A08	HQP004658	NM_001621	AHR
QG007-01	A09	HQP001961	NM_001283	AP1S1
QG007-01	A10	HQP009024	NM_000038	APC
QG007-01	A11	HQP009801	NM_000044	AR
QG007-01	A12	HQP010924	NM_001668	ARNT
QG007-01	B01	HQP011736	NM_000051	ATM
QG007-01	B02	HQP015964	NM_004324	BAX
QG007-01	B03	HQP016211	NM_000633	BCL2
QG007-01	B04	HQP016238	NM_138578	BCL2L1
QG007-01	B05	HQP016897	NM_000386	BLMH
QG007-01	B06	HQP017713	NM_007294	BRCA1
QG007-01	B07	HQP017753	NM_000059	BRCA2
QG007-01	B08	HQP016204	NM_053056	CCND1
QG007-01	B09	HQP021819	NM_001238	CCNE1
QG007-01	B10	HQP000225	NM_001798	CDK2
QG007-01	B11	HQP000245	NM_000075	CDK4
QG007-01	B12	HQP000331	NM_000389	CDKN1A
QG007-01	C01	HQP000342	NM_004064	CDKN1B
QG007-01	C02	HQP000369	NM_000077	CDKN2A
QG007-01	C03	HQP000408	NM_001800	CDKN2D
QG007-01	C04	HQP019829	NM_030782	CLPTM1L
QG007-01	C05	HQP003772	NM_000499	CYP1A1
QG007-01	C06	HQP003774	NM_000761	CYP1A2
QG007-01	C07	HQP003808	NM_000767	CYP2B6
QG007-01	C08	HQP003809	NM_000769	CYP2C19
QG007-01	C09	HQP003810	NM_000770	CYP2C8
QG007-01	C10	HQP003811	NM_000771	CYP2C9
QG007-01	C11	HQP003814	NM_000106	CYP2D6
QG007-01	C12	HQP003817	NM_000773	CYP2E1
QG007-01	D01	HQP003836	NM_017460	CYP3A4
QG007-01	D02	HQP003841	NM_000777	CYP3A5
QG007-01	D03	HQP004309	NM_000791	DHFR
QG007-01	D04	HQP004605	NM_005228	EGFR
QG007-01	D05	HQP004749	NM_005229	ELK1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG007-01	D06	HQP004948	NM_000120	EPHX1
QG007-01	D07	HQP004969	NM_004448	ERBB2
QG007-01	D08	HQP004971	NM_001982	ERBB3
QG007-01	D09	HQP004983	NM_000122	ERCC3
QG007-01	D10	HQP004998	NM_000125	ESR1
QG007-01	D11	HQP005002	NM_001437	ESR2
QG007-01	D12	HQP005403	NM_002006	FGF2
QG007-01	E01	HQP006188	NM_005252	FOS
QG007-01	E02	HQP008468	NM_019884	GSK3A
QG007-01	E03	HQP008487	NM_000852	GSTP1
QG007-01	E04	HQP008831	NM_001530	HIF1A
QG007-01	E05	HQP009523	NM_000875	IGF1R
QG007-01	E06	HQP009532	NM_000876	IGF2R
QG007-01	E07	HQP011181	NM_000245	MET
QG007-01	E08	HQP011491	NM_000251	MSH2
QG007-01	E09	HQP054016	NM_017458	MVP
QG007-01	E10	HQP011597	NM_002467	MYC
QG007-01	E11	HQP001136	NM_000015	NAT2
QG007-01	E12	HQP011807	NM_003998	NFKB1
QG007-01	F01	HQP053985	NM_002502	NFKB2
QG007-01	F02	HQP011812	NM_002503	NFKBIB
QG007-01	F03	HQP011813	NM_004556	NFKBIE
QG007-01	F04	HQP054001	NM_005036	PPARA
QG007-01	F05	HQP013627	NM_006238	PPARD
QG007-01	F06	HQP013634	NM_015869	PPARG
QG007-01	F07	HQP016114	NM_000964	RARA
QG007-01	F08	HQP016116	NM_000965	RARB
QG007-01	F09	HQP016118	NM_000966	RARG
QG007-01	F10	HQP016131	NM_000321	RB1
QG007-01	F11	HQP016214	NM_006509	RELB
QG007-01	F12	HQP016526	NM_002957	RXRA
QG007-01	G01	HQP016527	NM_021976	RXRB
QG007-01	G02	HQP017615	NM_000454	SOD1
QG007-01	G03	HQP017784	NM_005420	SULT1E1
QG007-01	G04	HQP021550	NM_003839	TNFRSF11A
QG007-01	G05	HQP018171	NM_003286	TOP1
QG007-01	G06	HQP018172	NM_001067	TOP2A
QG007-01	G07	HQP018173	NM_001068	TOP2B
QG007-01	G08	HQP018175	NM_000546	TP53
QG007-01	G09	HQP018556	NM_004628	XPC
QG007-01	G10	HQP009802	NM_001011645	AR

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG007-01	G11	HQP009026	NM_000194	HPRT1
QG007-01	G12	HQP000074	NM_001023587	ABCC5
QG007-01	H01	HGDC		
QG007-01	H02	HGDC		
QG007-01	H03	HQP006940	NM_002046	GAPDH
QG007-01	H04	HQP016381	NM_001101	ACTB
QG007-01	H05	HQP015171	NM_004048	B2M
QG007-01	H06	HQP006171	NM_012423	RPL13A
QG007-01	H07	HQP009026	NM_000194	HPRT1
QG007-01	H08	HQP054253	NR_003286	RN18S1
QG007-01	H09	RT		
QG007-01	H10	RT		
QG007-01	H11	PCR		
QG007-01	H12	PCR		