

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG014-01	A01	HQP018437	NM_003366	UQCRC2
QG014-01	A02	HQP018436	NM_003365	UQCRC1
QG014-01	A03	HQP000708	NM_006440	TXNRD2
QG014-01	A04	HQP018336	NM_003330	TXNRD1
QG014-01	A05	HQP019885	NM_030755	TXNDC1
QG014-01	A06	HQP006764	NM_012473	TXN2
QG014-01	A07	HQP018335	NM_003329	TXN
QG014-01	A08	HQP016694	NM_001035511	SDHC
QG014-01	A09	HQP016689	NM_003000	SDHB
QG014-01	A10	HQP016682	NM_004168	SDHA
QG014-01	A11	HQP015586	NM_000961	PTGIS
QG014-01	A12	HQP001016	NM_013261	PPARGC1A
QG014-01	B01	HQP013504	NM_000941	POR
QG014-01	B02	HQP012109	NM_000918	P4HB
QG014-01	B03	HQP000086	NM_005693	NR1H3
QG014-01	B04	HQP011858	NM_000904	NQO2
QG014-01	B05	HQP004317	NM_000903	NQO1
QG014-01	B06	HQP011868	NM_000603	NOS3
QG014-01	B07	HQP011866	NM_000625	NOS2A
QG014-01	B08	HQP011735	NM_021074	NDUFV2
QG014-01	B09	HQP011730	NM_007103	NDUFV1
QG014-01	B10	HQP011731	NM_002495	NDUFS4
QG014-01	B11	HQP011726	NM_005006	NDUFS1
QG014-01	B12	HQP011721	NM_005004	NDUFB8
QG014-01	C01	HQP011706	NM_002490	NDUFA6
QG014-01	C02	HQP011008	NM_000898	MAOB
QG014-01	C03	HQP011007	NM_000240	MAOA
QG014-01	C04	HQP009845	NM_002225	IVD
QG014-01	C05	HQP013447	NM_017545	HAO1
QG014-01	C06	HQP008473	NM_000637	GSR
QG014-01	C07	HQP007813	NM_000408	GPD2
QG014-01	C08	HQP005968	NM_002022	FMO4
QG014-01	C09	HQP005959	NM_001002294	FMO3
QG014-01	C10	HQP005951	NM_001460	FMO2
QG014-01	C11	HQP005939	NM_002021	FMO1
QG014-01	C12	HQP005387	NM_004110	FDXR
QG014-01	D01	HQP003865	NM_004391	CYP8B1
QG014-01	D02	HQP022723	NM_004820	CYP7B1
QG014-01	D03	HQP003859	NM_000780	CYP7A1
QG014-01	D04	HQP003924	NM_000786	CYP51A1
QG014-01	D05	HQP010910	NM_000896	CYP4F3

QG014-01	D06	HQP017537	NM_023944	CYP4F12
QG014-01	D07	HQP003855	NM_000779	CYP4B1
QG014-01	D08	HQP003847	NM_000778	CYP4A11
QG014-01	D09	HQP000977	NM_006668	CYP46A1
QG014-01	D10	HQP003841	NM_000777	CYP3A5
QG014-01	D11	HQP017181	NM_022820	CYP3A43
QG014-01	D12	HQP003836	NM_017460	CYP3A4
QG014-01	E01	HQP012613	NM_016593	CYP39A1
QG014-01	E02	HQP008529	NM_030622	CYP2S1
QG014-01	E03	HQP002083	NM_024514	CYP2R1
QG014-01	E04	HQP003823	NM_000775	CYP2J2
QG014-01	E05	HQP003818	NM_000774	CYP2F1
QG014-01	E06	HQP003817	NM_000773	CYP2E1
QG014-01	E07	HQP003814	NM_000106	CYP2D6
QG014-01	E08	HQP003811	NM_000771	CYP2C9
QG014-01	E09	HQP003809	NM_000769	CYP2C19
QG014-01	E10	HQP003808	NM_000767	CYP2B6
QG014-01	E11	HQP003787	NM_000762	CYP2A6
QG014-01	E12	HQP003804	NM_000766	CYP2A13
QG014-01	F01	HQP003923	NM_000785	CYP27B1
QG014-01	F02	HQP003922	NM_000784	CYP27A1
QG014-01	F03	HQP003919	NM_000783	CYP26A1
QG014-01	F04	HQP003917	NM_000782	CYP24A1
QG014-01	F05	HQP003908	NM_000500	CYP21A2
QG014-01	F06	HQP003775	NM_000104	CYP1B1
QG014-01	F07	HQP003774	NM_000761	CYP1A2
QG014-01	F08	HQP003772	NM_000499	CYP1A1
QG014-01	F09	HQP003904	NM_000103	CYP19A1
QG014-01	F10	HQP003888	NM_000102	CYP17A1
QG014-01	F11	HQP003871	NM_000781	CYP11A1
QG014-01	F12	HQP004315	NM_000398	CYP5B3
QG014-01	G01	HQP017020	NM_000712	BLVRA
QG014-01	G02	HQP006359	NM_000698	ALOX5
QG014-01	G03	HQP006356	NM_000697	ALOX12
QG014-01	G04	HQP013061	NM_004035	ACOX1
QG014-01	G05	HQP010079	NM_000018	ACADVL
QG014-01	G06	HQP009706	NM_000017	ACADS
QG014-01	G07	HQP009284	NM_001608	ACADL
QG014-01	G08	HQP007422	NM_001091	ABP1
QG014-01	G09	HQP003750	NM_001916	CYC1
QG014-01	G10	HQP003878	NM_000497	CYP11B1
QG014-01	G11	HQP009569	NM_000016	ACADM
QG014-01	G12	HQP009833	NM_001609	ACADSB

QG014-01	H01	HGDC		
QG014-01	H02	HGDC		
QG014-01	H03	HQP006940	NM_002046	GAPDH
QG014-01	H04	HQP016381	NM_001101	ACTB
QG014-01	H05	HQP015171	NM_004048	B2M
QG014-01	H06	HQP006171	NM_012423	RPL13A
QG014-01	H07	HQP009026	NM_000194	HPRT1
QG014-01	H08	HQP054253	NR_003286	RN18S1
QG014-01	H09	RT		
QG014-01	H10	RT		
QG014-01	H11	PCR		
QG014-01	H12	PCR		