

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG012-01	A01	HQP018565	NM_003401	XRCC4
QG012-01	A02	HQP018564	NM_005432	XRCC3
QG012-01	A03	HQP018563	NM_005431	XRCC2
QG012-01	A04	HQP018562	NM_006297	XRCC1
QG012-01	A05	HQP018556	NM_004628	XPC
QG012-01	A06	HQP015272	NM_020196	XAB2
QG012-01	A07	HQP008445	NM_001048201	UHRF1
QG012-01	A08	HQP018176	NM_005657	TP53BP1
QG012-01	A09	HQP018175	NM_000546	TP53
QG012-01	A10	HQP014691	NM_001008744	TDP1
QG012-01	A11	HQP017996	NM_003211	TDG
QG012-01	A12	HQP006236	NM_014311	SMUG1
QG012-01	B01	HQP016223	NM_002912	REV3L
QG012-01	B02	HQP012755	NM_001037872	REV1
QG012-01	B03	HQP016087	NM_134424	RAD52
QG012-01	B04	HQP016079	NM_002876	RAD51C
QG012-01	B05	HQP016077	NM_002875	RAD51
QG012-01	B06	HQP016076	NM_002874	RAD23B
QG012-01	B07	HQP016075	NM_005053	RAD23A
QG012-01	B08	HQP016074	NM_006265	RAD21
QG012-01	B09	HQP015190	NM_020165	RAD18
QG012-01	B10	HQP016070	NM_002873	RAD17
QG012-01	B11	HQP015947	NM_002853	RAD1
QG012-01	B12	HQP007689	NM_013274	POLL
QG012-01	C01	HQP012725	NM_016218	POLK
QG012-01	C02	HQP001399	NM_007195	POLI
QG012-01	C03	HQP013425	NM_006231	POLE
QG012-01	C04	HQP013423	NM_002691	POLD1
QG012-01	C05	HQP013422	NM_002690	POLB
QG012-01	C06	HQP013352	NM_000535	PMS2
QG012-01	C07	HQP012420	NM_002592	PCNA
QG012-01	C08	HQP014222	NM_018248	NEIL3
QG012-01	C09	HQP006460	NM_145043	NEIL2
QG012-01	C10	HQP019141	NM_024608	NEIL1
QG012-01	C11	HQP008493	NM_000179	MSH6
QG012-01	C12	HQP011492	NM_002439	MSH3
QG012-01	D01	HQP011491	NM_000251	MSH2
QG012-01	D02	HQP011320	NM_005590	MRE11A
QG012-01	D03	HQP011304	NM_001015052	MPG
QG012-01	D04	HQP011241	NM_005933	MLL
QG012-01	D05	HQP011235	NM_000249	MLH1

QG012-01	D06	HQP021736	NM_003925	MBD4
QG012-01	D07	HQP010613	NM_002312	LIG4
QG012-01	D08	HQP010609	NM_000234	LIG1
QG012-01	D09	HQP009138	NM_004507	HUS1
QG012-01	D10	HQP008505	NM_001517	GTF2H4
QG012-01	D11	HQP008504	NM_001516	GTF2H3
QG012-01	D12	HQP008502	NM_005316	GTF2H1
QG012-01	E01	HQP001042	NM_006705	GADD45G
QG012-01	E02	HQP004125	NM_001924	GADD45A
QG012-01	E03	HQP005391	NM_004111	FEN1
QG012-01	E04	HQP005083	NM_000136	FANCC
QG012-01	E05	HQP005081	NM_000135	FANCA
QG012-01	E06	HQP004984	NM_005236	ERCC4
QG012-01	E07	HQP004983	NM_000122	ERCC3
QG012-01	E08	HQP004976	NM_000400	ERCC2
QG012-01	E09	HQP004974	NM_001983	ERCC1
QG012-01	E10	HQP004114	NM_000107	DDB2
QG012-01	E11	HQP004111	NM_001923	DDB1
QG012-01	E12	HQP016982	NM_001033855	DCLRE1C
QG012-01	F01	HQP017213	NM_022836	DCLRE1B
QG012-01	F02	HQP023381	NM_014881	DCLRE1A
QG012-01	F03	HQP017753	NM_000059	BRCA2
QG012-01	F04	HQP017713	NM_007294	BRCA1
QG012-01	F05	HQP016844	NM_000057	BLM
QG012-01	F06	HQP013636	NM_000489	ATRX
QG012-01	F07	HQP011736	NM_000051	ATM
QG012-01	F08	HQP007655	NM_014481	APEX2
QG012-01	F09	HQP009061	NM_001641	APEX1
QG012-01	F10	HQP011937	NM_002528	NTHL1
QG012-01	F11	HQP005099	NM_004629	FANCG
QG012-01	F12	HQP008703	NM_002105	H2AFX
QG012-01	G01	HQP018427	NM_003362	UNG
QG012-01	G02	HQP018555	NM_000380	XPA
QG012-01	G03	HQP011494	NM_002441	MSH5
QG012-01	G04	HQP011576	NM_001048171	MUTYH
QG012-01	G05	HQP011686	NM_001024688	NBN
QG012-01	G06	HQP000048	NM_001042618	PARP2
QG012-01	G07	HQP001521	NM_007254	PNKP
QG012-01	G08	HQP013429	NM_006502	POLH
QG012-01	G09	HQP014837	NM_001081640	PRKDC
QG012-01	G10	HQP011493	NM_002440	MSH4
QG012-01	G11	HQP016230	NM_007370	RFC5
QG012-01	G12	HQP001503	NM_016381	TREX1

QG012-01	H01	HGDC		
QG012-01	H02	HGDC		
QG012-01	H03	HQP006940	NM_002046	GAPDH
QG012-01	H04	HQP016381	NM_001101	ACTB
QG012-01	H05	HQP015171	NM_004048	B2M
QG012-01	H06	HQP006171	NM_012423	RPL13A
QG012-01	H07	HQP009026	NM_000194	HPRT1
QG012-01	H08	HQP054253	NR_003286	RN18S1
QG012-01	H09	RT		
QG012-01	H10	RT		
QG012-01	H11	PCR		
QG012-01	H12	PCR		