

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-01	A01	HQP017753	NM_000059	BRCA2
QG075-01	A02	HQP018175	NM_000546	TP53
QG075-01	A03	HQP013099	NM_000926	PGR
QG075-01	A04	HQP018475	NM_001025366	VEGFA
QG075-01	A05	HQP003772	NM_000499	CYP1A1
QG075-01	A06	HQP010133	NM_004985	KRAS
QG075-01	A07	HQP009801	NM_000044	AR
QG075-01	A08	HQP000370	NM_058195	CDKN2A
QG075-01	A09	HQP018474	NM_000376	VDR
QG075-01	A10	HQP017733	NM_004333	BRAF
QG075-01	A11	HQP001396	NM_001005735	CHEK2
QG075-01	A12	HQP016131	NM_000321	RB1
QG075-01	B01	HQP004976	NM_000400	ERCC2
QG075-01	B02	HQP004975	NM_202001	ERCC1
QG075-01	B03	HQP002671	NM_000754	COMT
QG075-01	B04	HQP013100	NM_000927	ABCB1
QG075-01	B05	HQP011547	NM_005957	MTHFR
QG075-01	B06	HQP011255	NM_002421	MMP1
QG075-01	B07	HQP003775	NM_000104	CYP1B1
QG075-01	B08	HQP021819	NM_001238	CCNE1
QG075-01	B09	HQP021754	NM_001759	CCND2
QG075-01	B10	HQP016204	NM_053056	CCND1
QG075-01	B11	HQP011687	NM_002485	NBN
QG075-01	B12	HQP011257	NM_002422	MMP3
QG075-01	C01	HQP011235	NM_000249	MLH1
QG075-01	C02	HQP009670	NM_000600	IL6
QG075-01	C03	HQP003841	NM_000777	CYP3A5
QG075-01	C04	HQP000245	NM_000075	CDK4
QG075-01	C05	HQP018562	NM_006297	XRCC1
QG075-01	C06	HQP017616	NM_000636	SOD2
QG075-01	C07	HQP016087	NM_134424	RAD52
QG075-01	C08	HQP015535	NM_000314	PTEN
QG075-01	C09	HQP013150	NM_006218	PIK3CA
QG075-01	C10	HQP011491	NM_000251	MSH2
QG075-01	C11	HQP009641	NM_000576	IL1B
QG075-01	C12	HQP008493	NM_000179	MSH6
QG075-01	D01	HQP006895	NM_000155	GALT
QG075-01	D02	HQP005002	NM_001437	ESR2
QG075-01	D03	HQP004998	NM_000125	ESR1
QG075-01	D04	HQP003888	NM_000102	CYP17A1
QG075-01	D05	HQP000408	NM_001800	CDKN2D

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-01	D06	HQP000396	NM_001262	CDKN2C
QG075-01	D07	HQP000225	NM_001798	CDK2
QG075-01	D08	HQP054054	BC039243	FGFR2
QG075-01	D09	HQP054047	BC008403	HLA-DRB1
QG075-01	D10	HQP018966	NM_001080124	CASP8
QG075-01	D11	HQP018564	NM_005432	XRCC3
QG075-01	D12	HQP018563	NM_005431	XRCC2
QG075-01	E01	HQP017698	NM_000348	SRD5A2
QG075-01	E02	HQP015598	NM_000963	PTGS2
QG075-01	E03	HQP013352	NM_000535	PMS2
QG075-01	E04	HQP013312	NM_000534	PMS1
QG075-01	E05	HQP013104	NM_002634	PHB
QG075-01	E06	HQP011492	NM_002439	MSH3
QG075-01	E07	HQP011320	NM_005590	MRE11A
QG075-01	E08	HQP011135	NM_002392	MDM2
QG075-01	E09	HQP009718	NM_001562	IL18
QG075-01	E10	HQP009685	NM_000572	IL10
QG075-01	E11	HQP009645	NM_000577	IL1RN
QG075-01	E12	HQP009640	NM_000575	IL1A
QG075-01	F01	HQP009544	NM_000598	IGFBP3
QG075-01	F02	HQP009518	NM_000618	IGF1
QG075-01	F03	HQP006435	NM_000145	FSHR
QG075-01	F04	HQP004599	NM_001963	EGF
QG075-01	F05	HQP004411	NM_022552	DNMT3A
QG075-01	F06	HQP003904	NM_000103	CYP19A1
QG075-01	F07	HQP003810	NM_000770	CYP2C8
QG075-01	F08	HQP000274	NM_001259	CDK6
QG075-01	F09	HQP053959	NM_001040	SHBG
QG075-01	F10	HQP023466	NM_004360	CDH1
QG075-01	F11	HQP022093	NM_004688	NMI
QG075-01	F12	HQP020040	NM_006534	NCOA3
QG075-01	G01	HQP018141	NM_000594	TNF
QG075-01	G02	HQP018044	NM_000660	TGFB1
QG075-01	G03	HQP017811	NM_177536	SULT1A1
QG075-01	G04	HQP016170	NM_021632	ZNF350
QG075-01	G05	HQP016155	NM_002894	RBBP8
QG075-01	G06	HQP016079	NM_002876	RAD51C
QG075-01	G07	HQP015946	NM_000465	BARD1
QG075-01	G08	HQP015596	NM_000962	PTGS1
QG075-01	G09	HQP013473	NM_000446	PON1
QG075-01	G10	HQP013204	NM_002658	PLAU

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-01	G11	HQP012154	NM_000602	SERPINE1
QG075-01	G12	HQP011810	NM_020529	NFKBIA
QG075-01	H01	HGDC		
QG075-01	H02	HGDC		
QG075-01	H03	HQP006940	NM_002046	GAPDH
QG075-01	H04	HQP016381	NM_001101	ACTB
QG075-01	H05	HQP015171	NM_004048	B2M
QG075-01	H06	HQP006171	NM_012423	RPL13A
QG075-01	H07	HQP009026	NM_000194	HPRT1
QG075-01	H08	HQP054253	NR_003286	RN18S1
QG075-01	H09	RT		
QG075-01	H10	RT		
QG075-01	H11	PCR		
QG075-01	H12	PCR		
QG075-02	A01	HQP011559	NM_001018016	MUC1
QG075-02	A02	HQP011309	NM_000250	MPO
QG075-02	A03	HQP011266	NM_002426	MMP12
QG075-02	A04	HQP011263	NM_004994	MMP9
QG075-02	A05	HQP011256	NM_004530	MMP2
QG075-02	A06	HQP010613	NM_002312	LIG4
QG075-02	A07	HQP009818	NM_000212	ITGB3
QG075-02	A08	HQP009749	NM_000207	INS
QG075-02	A09	HQP009642	NM_000877	IL1R1
QG075-02	A10	HQP009551	NM_000599	IGFBP5
QG075-02	A11	HQP009539	NM_000596	IGFBP1
QG075-02	A12	HQP009523	NM_000875	IGF1R
QG075-02	B01	HQP009024	NM_000038	APC
QG075-02	B02	HQP008849	NM_002116	HLA-A
QG075-02	B03	HQP006839	NM_020474	GALNT1
QG075-02	B04	HQP005403	NM_002006	FGF2
QG075-02	B05	HQP004948	NM_000120	EPHX1
QG075-02	B06	HQP004605	NM_005228	EGFR
QG075-02	B07	HQP003817	NM_000773	CYP2E1
QG075-02	B08	HQP003811	NM_000771	CYP2C9
QG075-02	B09	HQP003808	NM_000767	CYP2B6
QG075-02	B10	HQP001136	NM_000015	NAT2
QG075-02	B11	HQP000145	NM_005732	RAD50
QG075-02	B12	HQP015106	NM_001077500	KLK10
QG075-02	C01	HQP013203	NM_033011	PLAT
QG075-02	C02	HQP004985	NM_000123	ERCC5
QG075-02	C03	HQP054691	NM_001159542	POU5F1P1

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-02	C04	HQP054057	BC071181	TGFBR1
QG075-02	C05	HQP054045	BC004257	RET
QG075-02	C06	HQP054030	NM_130398	EXO1
QG075-02	C07	HQP053966	NM_005921	MAP3K1
QG075-02	C08	HQP053910	NM_130851	BMP4
QG075-02	C09	HQP023399	NM_004358	CDC25B
QG075-02	C10	HQP023385	NM_001789	CDC25A
QG075-02	C11	HQP023365	NM_001255	CDC20
QG075-02	C12	HQP023132	NM_014705	DOCK4
QG075-02	D01	HQP023013	NM_005101	ISG15
QG075-02	D02	HQP022978	NM_006311	NCOR1
QG075-02	D03	HQP022893	NM_004887	CXCL14
QG075-02	D04	HQP022884	NM_004881	TP53I3
QG075-02	D05	HQP022677	NM_024690	MUC16
QG075-02	D06	HQP022491	NM_001616	ACVR2A
QG075-02	D07	HQP022362	NM_004219	PTTG1
QG075-02	D08	HQP022143	NM_057749	CCNE2
QG075-02	D09	HQP022141	NM_004701	CCNB2
QG075-02	D10	HQP022071	NM_001105	ACVR1
QG075-02	D11	HQP021882	NM_004354	CCNG2
QG075-02	D12	HQP021857	NM_004060	CCNG1
QG075-02	E01	HQP021763	NM_004199	P4HA2
QG075-02	E02	HQP021727	NM_031966	CCNB1
QG075-02	E03	HQP021646	NM_003896	ST3GAL5
QG075-02	E04	HQP021621	NM_003884	PCAF
QG075-02	E05	HQP021610	NM_003882	WISP1
QG075-02	E06	HQP021604	NM_003879	CFLAR
QG075-02	E07	HQP021557	NM_003844	TNFRSF10A
QG075-02	E08	HQP021502	NM_003810	TNFSF10
QG075-02	E09	HQP021456	NM_001040147	SERPINB7
QG075-02	E10	HQP021385	NM_003739	AKR1C3
QG075-02	E11	HQP021368	NM_001075099	SKAP1
QG075-02	E12	HQP021329	NM_003707	RUVBL1
QG075-02	F01	HQP021305	NM_003690	PRKRA
QG075-02	F02	HQP021142	NM_003641	IFITM1
QG075-02	F03	HQP020999	NM_032797	AIFM2
QG075-02	F04	HQP020924	NM_181672	OGT
QG075-02	F05	HQP020798	NM_003592	CUL1
QG075-02	F06	HQP020384	NM_004347	CASP5
QG075-02	F07	HQP020117	NM_004656	BAP1
QG075-02	F08	HQP020116	NM_004655	AXIN2

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-02	F09	HQP020042	NM_003489	NRIP1
QG075-02	F10	HQP019539	NM_025072	PTGES2
QG075-02	F11	HQP019178	NM_001033952	CALCA
QG075-02	F12	HQP018760	NM_003463	PTP4A1
QG075-02	G01	HQP018755	NM_001007257	PRDM2
QG075-02	G02	HQP018568	NM_021141	XRCC5
QG075-02	G03	HQP018565	NM_003401	XRCC4
QG075-02	G04	HQP018556	NM_004628	XPC
QG075-02	G05	HQP018538	NM_003394	WNT10B
QG075-02	G06	HQP018531	NM_003392	WNT5A
QG075-02	G07	HQP018529	NM_003391	WNT2
QG075-02	G08	HQP018488	NM_003379	VIL2
QG075-02	G09	HQP018483	NM_005429	VEGFC
QG075-02	G10	HQP018342	NM_001071	TYMS
QG075-02	G11	HQP018340	NM_003331	TYK2
QG075-02	G12	HQP018328	NM_000474	TWIST1
QG075-02	H01	HGDC		
QG075-02	H02	HGDC		
QG075-02	H03	HQP006940	NM_002046	GAPDH
QG075-02	H04	HQP016381	NM_001101	ACTB
QG075-02	H05	HQP015171	NM_004048	B2M
QG075-02	H06	HQP006171	NM_012423	RPL13A
QG075-02	H07	HQP009026	NM_000194	HPRT1
QG075-02	H08	HQP054253	NR_003286	RN18S1
QG075-02	H09	RT		
QG075-02	H10	RT		
QG075-02	H11	PCR		
QG075-02	H12	PCR		
QG075-03	A01	HQP018149	NM_001066	TNFRSF1B
QG075-03	A02	HQP018145	NM_006290	TNFAIP3
QG075-03	A03	HQP018144	NM_006291	TNFAIP2
QG075-03	A04	HQP018085	NM_005655	KLF10
QG075-03	A05	HQP018072	NM_003248	THBS4
QG075-03	A06	HQP018047	NM_003238	TGFB2
QG075-03	A07	HQP018033	NM_006286	TFDP2
QG075-03	A08	HQP018032	NM_007111	TFDP1
QG075-03	A09	HQP017999	NM_004336	BUB1
QG075-03	A10	HQP017969	NM_005649	ZNF354A
QG075-03	A11	HQP017938	NM_001061	TBXAS1
QG075-03	A12	HQP017844	NM_004335	BST2
QG075-03	B01	HQP017767	NM_003150	STAT3

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-03	B02	HQP017764	NM_007315	STAT1
QG075-03	B03	HQP017729	NM_003141	TRIM21
QG075-03	B04	HQP017697	NM_001047	SRD5A1
QG075-03	B05	HQP017673	NM_000582	SPP1
QG075-03	B06	HQP017653	NM_003118	SPARC
QG075-03	B07	HQP017635	NM_000346	SOX9
QG075-03	B08	HQP017623	NM_003104	SORD
QG075-03	B09	HQP017615	NM_000454	SOD1
QG075-03	B10	HQP017300	NM_005983	SKP2
QG075-03	B11	HQP016961	NM_001012398	AKTIP
QG075-03	B12	HQP016801	NM_022162	NOD2
QG075-03	C01	HQP016648	NM_002993	CXCL6
QG075-03	C02	HQP016630	NM_005408	CCL13
QG075-03	C03	HQP016629	NM_002986	CCL11
QG075-03	C04	HQP016572	NM_002970	SAT1
QG075-03	C05	HQP016543	NM_002963	S100A7
QG075-03	C06	HQP016537	NM_002959	SORT1
QG075-03	C07	HQP016500	NM_001032295	SERPING1
QG075-03	C08	HQP016426	NM_021134	MRPL23
QG075-03	C09	HQP016418	NM_004327	BCR
QG075-03	C10	HQP016252	NM_002923	RGS2
QG075-03	C11	HQP016212	NM_000657	BCL2
QG075-03	C12	HQP016172	NM_005611	RBL2
QG075-03	D01	HQP016160	NM_002895	RBL1
QG075-03	D02	HQP015828	NM_002827	PTPN1
QG075-03	D03	HQP015610	NM_002820	PTHLH
QG075-03	D04	HQP015586	NM_000961	PTGIS
QG075-03	D05	HQP015555	NM_198714	PTGER3
QG075-03	D06	HQP015540	NM_000955	PTGER1
QG075-03	D07	HQP015537	NM_000953	PTGDR
QG075-03	D08	HQP015496	NM_001186	BACH1
QG075-03	D09	HQP015184	NM_152278	TCEAL7
QG075-03	D10	HQP014969	NM_004705	PRKRIR
QG075-03	D11	HQP014948	NM_002759	EIF2AK2
QG075-03	D12	HQP014848	NM_002745	MAPK1
QG075-03	E01	HQP014787	NM_018479	ECHDC1
QG075-03	E02	HQP013633	NM_005037	PPARG
QG075-03	E03	HQP013627	NM_006238	PPARD
QG075-03	E04	HQP013596	NM_001184	ATR
QG075-03	E05	HQP013583	NM_205862	UGT1A6
QG075-03	E06	HQP013454	NM_005034	POLR2K

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-03	E07	HQP013416	NM_001018111	PODXL
QG075-03	E08	HQP013271	NM_182943	PLOD2
QG075-03	E09	HQP013269	NM_000302	PLOD1
QG075-03	E10	HQP013262	NM_005030	PLK1
QG075-03	E11	HQP013205	NM_001005376	PLAUR
QG075-03	E12	HQP012998	NM_016373	WWOX
QG075-03	F01	HQP012986	NM_016290	UIMC1
QG075-03	F02	HQP012889	NM_002609	PDGFRB
QG075-03	F03	HQP012857	NM_033016	PDGFB
QG075-03	F04	HQP012770	NM_016371	HSD17B7
QG075-03	F05	HQP012694	NM_015986	CRLF3
QG075-03	F06	HQP012651	NM_016639	TNFRSF12A
QG075-03	F07	HQP012447	NM_016142	HSD17B12
QG075-03	F08	HQP012420	NM_002592	PCNA
QG075-03	F09	HQP012155	NM_002575	SERPINB2
QG075-03	F10	HQP011983	NM_001032409	OAS1
QG075-03	F11	HQP011955	NM_001014796	DDR2
QG075-03	F12	HQP011932	NM_002526	NT5E
QG075-03	G01	HQP011876	NM_000435	NOTCH3
QG075-03	G02	HQP011868	NM_000603	NOS3
QG075-03	G03	HQP011866	NM_000625	NOS2A
QG075-03	G04	HQP011852	NM_001678	ATP1B2
QG075-03	G05	HQP011812	NM_002503	NFKBIB
QG075-03	G06	HQP011808	NM_001077493	NFKB2
QG075-03	G07	HQP011597	NM_002467	MYC
QG075-03	G08	HQP011582	NM_002462	MX1
QG075-03	G09	HQP011325	NM_019899	ABCC1
QG075-03	G10	HQP011268	NM_004995	MMP14
QG075-03	G11	HQP011220	NM_002416	CXCL9
QG075-03	G12	HQP011187	NM_002404	MFAP4
QG075-03	H01	HGDC		
QG075-03	H02	HGDC		
QG075-03	H03	HQP006940	NM_002046	GAPDH
QG075-03	H04	HQP016381	NM_001101	ACTB
QG075-03	H05	HQP015171	NM_004048	B2M
QG075-03	H06	HQP006171	NM_012423	RPL13A
QG075-03	H07	HQP009026	NM_000194	HPRT1
QG075-03	H08	HQP054253	NR_003286	RN18S1
QG075-03	H09	RT		
QG075-03	H10	RT		
QG075-03	H11	PCR		

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-03	H12	PCR		
QG075-04	A01	HQP011179	NM_002402	MEST
QG075-04	A02	HQP011140	NM_002393	MDM4
QG075-04	A03	HQP011110	NM_005915	MCM6
QG075-04	A04	HQP011106	NM_004526	MCM2
QG075-04	A05	HQP011024	NM_016835	MAPT
QG075-04	A06	HQP010996	NM_022438	MAL
QG075-04	A07	HQP010960	NM_005902	SMAD3
QG075-04	A08	HQP010942	NM_002354	TACSTD1
QG075-04	A09	HQP010912	NM_206943	LTBP1
QG075-04	A10	HQP010612	NM_013975	LIG3
QG075-04	A11	HQP010609	NM_000234	LIG1
QG075-04	A12	HQP010390	NM_000425	L1CAM
QG075-04	B01	HQP010125	NM_002266	KPNA2
QG075-04	B02	HQP009941	NM_001657	AREG
QG075-04	B03	HQP009853	NM_002228	JUN
QG075-04	B04	HQP009810	NM_002211	ITGB1
QG075-04	B05	HQP009808	NM_002210	ITGAV
QG075-04	B06	HQP009785	NM_001572	IRF7
QG075-04	B07	HQP009774	NM_000210	ITGA6
QG075-04	B08	HQP009764	NM_000208	INSR
QG075-04	B09	HQP009745	NM_005538	INHBC
QG075-04	B10	HQP009744	NM_002193	INHBB
QG075-04	B11	HQP009681	NM_001557	IL8RB
QG075-04	B12	HQP009678	NM_000584	IL8
QG075-04	C01	HQP009675	NM_175767	IL6ST
QG075-04	C02	HQP009671	NM_000639	FASLG
QG075-04	C03	HQP009555	NM_002178	IGFBP6
QG075-04	C04	HQP009548	NM_001552	IGFBP4
QG075-04	C05	HQP009541	NM_000597	IGFBP2
QG075-04	C06	HQP009529	NM_000612	IGF2
QG075-04	C07	HQP009408	NM_001548	IFIT1
QG075-04	C08	HQP009393	NM_005533	IFI35
QG075-04	C09	HQP009379	NM_005531	IFI16
QG075-04	C10	HQP009273	NM_002166	ID2
QG075-04	C11	HQP009101	NM_001168	BIRC5
QG075-04	C12	HQP009067	NM_000414	HSD17B4
QG075-04	D01	HQP009064	NM_000413	HSD17B1
QG075-04	D02	HQP009061	NM_001641	APEX1
QG075-04	D03	HQP009036	NM_005343	HRAS
QG075-04	D04	HQP008832	NM_181054	HIF1A



Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-04	D05	HQP008751	NM_018063	HELLS
QG075-04	D06	HQP008746	NM_001527	HDAC2
QG075-04	D07	HQP008745	NM_004964	HDAC1
QG075-04	D08	HQP008693	NM_005318	H1FO
QG075-04	D09	HQP008616	NM_013369	DNMT3L
QG075-04	D10	HQP008483	NM_000849	GSTM3
QG075-04	D11	HQP008459	NM_002090	CXCL3
QG075-04	D12	HQP008346	NM_014059	C13orf15
QG075-04	E01	HQP008285	NM_002085	GPX4
QG075-04	E02	HQP007781	NM_000406	GNRHR
QG075-04	E03	HQP007421	NM_020547	AMHR2
QG075-04	E04	HQP007382	NM_012449	STEAP1
QG075-04	E05	HQP006992	NM_015557	CHD5
QG075-04	E06	HQP006547	NM_001469	XRCC6
QG075-04	E07	HQP006446	NM_000147	FUCA1
QG075-04	E08	HQP006432	NM_000510	FSHB
QG075-04	E09	HQP006427	NM_001039130	ALOX15B
QG075-04	E10	HQP006368	NM_015368	PANX1
QG075-04	E11	HQP006323	NM_031412	GABARAPL1
QG075-04	E12	HQP006188	NM_005252	FOS
QG075-04	F01	HQP006023	NM_054034	FN1
QG075-04	F02	HQP006010	NM_033071	SYNE1
QG075-04	F03	HQP005663	NM_005250	FOXL1
QG075-04	F04	HQP005612	NM_012242	DKK1
QG075-04	F05	HQP005100	NM_000691	ALDH3A1
QG075-04	F06	HQP005098	NM_022725	FANCF
QG075-04	F07	HQP005084	NM_001018115	FANCD2
QG075-04	F08	HQP005083	NM_000136	FANCC
QG075-04	F09	HQP005081	NM_000135	FANCA
QG075-04	F10	HQP005075	NM_000689	ALDH1A1
QG075-04	F11	HQP005069	NM_000129	F13A1
QG075-04	F12	HQP005058	NM_000130	F5
QG075-04	G01	HQP005052	NM_000506	F2
QG075-04	G02	HQP005003	NM_004451	ESRRA
QG075-04	G03	HQP004995	NM_001626	AKT2
QG075-04	G04	HQP004991	NM_001014431	AKT1
QG075-04	G05	HQP004986	NM_000124	ERCC6
QG075-04	G06	HQP004984	NM_005236	ERCC4
QG075-04	G07	HQP004983	NM_000122	ERCC3
QG075-04	G08	HQP004971	NM_001982	ERBB3
QG075-04	G09	HQP004897	NM_001429	EP300

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG075-04	G10	HQP004677	NM_004096	EIF4EBP2
QG075-04	G11	HQP004613	NM_000399	EGR2
QG075-04	G12	HQP004528	NM_001950	E2F4
QG075-04	H01	HGDC		
QG075-04	H02	HGDC		
QG075-04	H03	HQP006940	NM_002046	GAPDH
QG075-04	H04	HQP016381	NM_001101	ACTB
QG075-04	H05	HQP015171	NM_004048	B2M
QG075-04	H06	HQP006171	NM_012423	RPL13A
QG075-04	H07	HQP009026	NM_000194	HPRT1
QG075-04	H08	HQP054253	NR_003286	RN18S1
QG075-04	H09	RT		
QG075-04	H10	RT		
QG075-04	H11	PCR		
QG075-04	H12	PCR		