

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG095-01	A01	HQP002611	NM_145259	ACVR1C
QG095-01	A02	HQP002467	NM_175866	UHMK1
QG095-01	A03	HQP002148	NM_145203	CSNK1A1L
QG095-01	A04	HQP002101	NM_198578	LRRK2
QG095-01	A05	HQP002012	NM_152309	PIK3AP1
QG095-01	A06	HQP001700	NM_052902	STK11IP
QG095-01	A07	HQP001522	NM_033550	TP53RK
QG095-01	A08	HQP022604	NM_033115	MGC16169
QG095-01	A09	HQP022363	NM_001003786	LYK5
QG095-01	A10	HQP022258	NM_182493	MLCK
QG095-01	A11	HQP022249	NM_033116	NEK9
QG095-01	A12	HQP022169	NM_033276	XRCC6BP1
QG095-01	B01	HQP022059	NM_052853	ADCK2
QG095-01	B02	HQP021196	NM_033118	MYLK2
QG095-01	B03	HQP021043	NM_032844	MASTL
QG095-01	B04	HQP020831	NM_032538	TTBK1
QG095-01	B05	HQP020737	NM_032435	KIAA1804
QG095-01	B06	HQP020729	NM_032430	BRSK1
QG095-01	B07	HQP020544	NM_032237	FLJ23356
QG095-01	B08	HQP020412	NM_032037	TSSK6
QG095-01	B09	HQP020401	NM_032028	TSSK1B
QG095-01	B10	HQP020391	NM_032017	STK40
QG095-01	B11	HQP019977	NM_030952	NUAK2
QG095-01	B12	HQP019947	NM_052841	TSSK3
QG095-01	C01	HQP019919	NM_030911	CDADC1
QG095-01	C02	HQP019655	NM_030626	LRRC27
QG095-01	C03	HQP019596	NM_025144	ALPK1
QG095-01	C04	HQP019401	NM_024876	ADCK4
QG095-01	C05	HQP019183	NM_024652	LRRK1
QG095-01	C06	HQP018898	NM_024091	FASTKD3
QG095-01	C07	HQP018852	NM_024046	CAMKV
QG095-01	C08	HQP017508	NM_030906	STK33
QG095-01	C09	HQP017387	NM_006648	WNK2
QG095-01	C10	HQP017385	NM_001002838	WNK3
QG095-01	C11	HQP017384	NM_032387	WNK4
QG095-01	C12	HQP017348	NM_018979	WNK1
QG095-01	D01	HQP017313	NM_139158	ALS2CR7
QG095-01	D02	HQP017170	NM_001031715	IQCH
QG095-01	D03	HQP016818	NM_022344	C17orf75
QG095-01	D04	HQP015862	NM_031417	MARK4
QG095-01	D05	HQP015857	NM_021158	TRIB3

QG095-01	D06	HQP015705	NM_020791	TAOK1
QG095-01	D07	HQP015689	NM_020778	ALPK3
QG095-01	D08	HQP015587	NM_001048218	SCYL1
QG095-01	D09	HQP015572	NM_020666	CLK4
QG095-01	D10	HQP015467	NM_020439	CAMK1G
QG095-01	D11	HQP015440	NM_020423	SCYL3
QG095-01	D12	HQP015408	NM_020397	CAMK1D
QG095-01	E01	HQP015320	NM_020247	CABC1
QG095-01	E02	HQP015124	NM_016567	BCCIP
QG095-01	E03	HQP015017	NM_031414	STK31
QG095-01	E04	HQP014794	NM_018492	PBK
QG095-01	E05	HQP014622	NM_018110	DOK4
QG095-01	E06	HQP014588	NM_017988	SCYL2
QG095-01	E07	HQP014489	NM_017593	BMP2K
QG095-01	E08	HQP014471	NM_017525	CDC42BPG
QG095-01	E09	HQP014403	NM_018571	ALS2CR2
QG095-01	E10	HQP014366	NM_018423	STYK1
QG095-01	E11	HQP014360	NM_018401	STK32B
QG095-01	E12	HQP014201	NM_018216	PANK4
QG095-01	F01	HQP013865	NM_017811	UBE2R2
QG095-01	F02	HQP013842	NM_017771	PXK
QG095-01	F03	HQP013788	NM_017719	SNRK
QG095-01	F04	HQP013734	NM_017672	TRPM7
QG095-01	F05	HQP013496	NM_019009	TOLLIP
QG095-01	F06	HQP013385	NM_020639	RIPK4
QG095-01	F07	HQP013348	NM_022048	CSNK1G1
QG095-01	F08	HQP013028	NM_016653	ZAK
QG095-01	F09	HQP013022	NM_001042452	RP6-213H19.1
QG095-01	F10	HQP013013	NM_016507	CRKRS
QG095-01	F11	HQP012972	NM_016231	NLK
QG095-01	F12	HQP012722	NM_001040633	PRKAG2
QG095-01	G01	HQP012666	NM_016281	TAOK3
QG095-01	G02	HQP012572	NM_016508	CDKL3
QG095-01	G03	HQP012543	NM_016441	CRIM1
QG095-01	G04	HQP012541	NM_001025778	VRK3
QG095-01	G05	HQP012441	NM_016123	IRAK4
QG095-01	G06	HQP012375	NM_015978	TNNI3K
QG095-01	G07	HQP012241	NM_016388	TRAT1
QG095-01	G08	HQP008817	NM_014602	PIK3R4
QG095-01	G09	HQP008771	NM_014586	HUNK
QG095-01	G10	HQP008685	NM_001024666	SH3KBP1
QG095-01	G11	HQP008669	NM_020804	PACSIN1
QG095-01	G12	HQP008633	NM_013392	NRBP1

QG095-01	H01	HGDC		
QG095-01	H02	HGDC		
QG095-01	H03	HQP006940	NM_002046	GAPDH
QG095-01	H04	HQP016381	NM_001101	ACTB
QG095-01	H05	HQP015171	NM_004048	B2M
QG095-01	H06	HQP006171	NM_012423	RPL13A
QG095-01	H07	HQP009026	NM_000194	HPRT1
QG095-01	H08	HQP054253	NR_003286	RN18S1
QG095-01	H09	RT		
QG095-01	H10	RT		
QG095-01	H11	PCR		
QG095-01	H12	PCR		
QG095-02	A01	HQP008609	NM_013355	PKN3
QG095-02	A02	HQP008578	NM_013302	EEF2K
QG095-02	A03	HQP008428	NM_013254	TBK1
QG095-02	A04	HQP008358	NM_022740	HIPK2
QG095-02	A05	HQP007693	NM_013233	STK39
QG095-02	A06	HQP007676	NM_014496	RPS6KA6
QG095-02	A07	HQP007549	NM_015690	STK36
QG095-02	A08	HQP007510	NM_014413	EIF2AK1
QG095-02	A09	HQP007372	NM_012424	RPS6KC1
QG095-02	A10	HQP007305	NM_014370	SRPK3
QG095-02	A11	HQP007288	NM_014572	LATS2
QG095-02	A12	HQP007224	NM_014365	HSPB8
QG095-02	B01	HQP006710	NM_015375	RIPK5
QG095-02	B02	HQP006319	NM_005813	PRKD3
QG095-02	B03	HQP006291	NM_012346	NUP62
QG095-02	B04	HQP006274	NM_053006	TSSK2
QG095-02	B05	HQP006265	NM_014326	DAPK2
QG095-02	B06	HQP006135	NM_014299	BRD4
QG095-02	B07	HQP006056	NM_025164	KIAA0999
QG095-02	B08	HQP005897	NM_015191	SNF1LK2
QG095-02	B09	HQP005832	NM_015148	PASK
QG095-02	B10	HQP005791	NM_015112	MAST2
QG095-02	B11	HQP005773	NM_015093	MAP3K7IP2
QG095-02	B12	HQP005703	NM_015092	SMG1
QG095-02	C01	HQP005699	NM_015028	TNIK
QG095-02	C02	HQP005665	NM_015000	STK38L
QG095-02	C03	HQP005522	NM_014929	FASTKD2
QG095-02	C04	HQP005508	NM_014920	ICK
QG095-02	C05	HQP005505	NM_014916	LMTK2
QG095-02	C06	HQP005501	NM_014911	AAK1
QG095-02	C07	HQP001608	NM_007284	TWF2

QG095-02	C08	HQP001580	NM_007271	STK38
QG095-02	C09	HQP001409	NM_006738	AKAP13
QG095-02	C10	HQP001408	NM_007199	IRAK3
QG095-02	C11	HQP001396	NM_001005735	CHEK2
QG095-02	C12	HQP001390	NM_007191	WIF1
QG095-02	D01	HQP001370	NM_006575	MAP4K5
QG095-02	D02	HQP001184	NM_006875	PIM2
QG095-02	D03	HQP001177	NM_006871	RIPK3
QG095-02	D04	HQP001055	NM_006712	FASTK
QG095-02	D05	HQP000920	NM_014397	NEK6
QG095-02	D06	HQP000904	NM_006622	PLK2
QG095-02	D07	HQP000881	NM_006609	MAP3K2
QG095-02	D08	HQP000868	NM_014264	PLK4
QG095-02	D09	HQP000781	NM_006549	CAMKK2
QG095-02	D10	HQP000776	NM_006545	TUSC4
QG095-02	D11	HQP000718	NM_033266	ERN2
QG095-02	D12	HQP000598	NM_006374	STK25
QG095-02	E01	HQP000555	NM_006343	MERTK
QG095-02	E02	HQP000503	NM_007170	TESK2
QG095-02	E03	HQP000419	NM_001007565	TFG
QG095-02	E04	HQP000365	NM_001014831	PAK4
QG095-02	E05	HQP000361	NM_005881	BCKDK
QG095-02	E06	HQP000276	NM_025195	TRIB1
QG095-02	E07	HQP000231	NM_001010938	TNK2
QG095-02	E08	HQP000149	NM_001048200	HIPK3
QG095-02	E09	HQP000113	NM_005713	COL4A3BP
QG095-02	E10	HQP000001	NM_005465	AKT3
QG095-02	E11	HQP023402	NM_005113	GOLGA5
QG095-02	E12	HQP023391	NM_005109	OXSR1
QG095-02	F01	HQP023331	NM_014840	NUAK1
QG095-02	F02	HQP023312	NM_012290	TLK1
QG095-02	F03	HQP023256	NM_014791	MELK
QG095-02	F04	HQP023153	NM_014720	SLK
QG095-02	F05	HQP023102	NM_014683	ULK2
QG095-02	F06	HQP023021	NM_014002	IKBKE
QG095-02	F07	HQP022933	NM_006035	CDC42BPB
QG095-02	F08	HQP022805	NM_004850	ROCK2
QG095-02	F09	HQP022771	NM_004836	EIF2AK3
QG095-02	F10	HQP022766	NM_004834	MAP4K4
QG095-02	F11	HQP022570	NM_004783	TAOK2
QG095-02	F12	HQP022428	NM_004760	STK17A
QG095-02	G01	HQP022427	NM_004226	STK17B
QG095-02	G02	HQP022404	NM_004755	RPS6KA5

QG095-02	G03	HQP022380	NM_004749	TBRG4
QG095-02	G04	HQP022323	NM_004217	AURKB
QG095-02	G05	HQP022304	NM_004734	DCLK1
QG095-02	G06	HQP022250	NM_004721	MAP3K13
QG095-02	G07	HQP022099	NM_004690	LATS1
QG095-02	G08	HQP021888	NM_003954	MAP3K14
QG095-02	G09	HQP021841	NM_003948	CDKL2
QG095-02	G10	HQP021838	NM_001024660	KALRN
QG095-02	G11	HQP021742	NM_003930	SKAP2
QG095-02	G12	HQP021688	NM_003913	PRPF4B
QG095-02	H01	HGDC		
QG095-02	H02	HGDC		
QG095-02	H03	HQP006940	NM_002046	GAPDH
QG095-02	H04	HQP016381	NM_001101	ACTB
QG095-02	H05	HQP015171	NM_004048	B2M
QG095-02	H06	HQP006171	NM_012423	RPL13A
QG095-02	H07	HQP009026	NM_000194	HPRT1
QG095-02	H08	HQP054253	NR_003286	RN18S1
QG095-02	H09	RT		
QG095-02	H10	RT		
QG095-02	H11	PCR		
QG095-02	H12	PCR		
QG095-03	A01	HQP021681	NM_003909	CPNE3
QG095-03	A02	HQP021622	NM_003885	CDK5R1
QG095-03	A03	HQP021615	NM_014238	KSR1
QG095-03	A04	HQP021589	NM_003872	NRP2
QG095-03	A05	HQP021574	NM_004196	CDKL1
QG095-03	A06	HQP021534	NM_003831	RIOK3
QG095-03	A07	HQP021524	NM_003821	RIPK2
QG095-03	A08	HQP021492	NM_003804	RIPK1
QG095-03	A09	HQP021368	NM_001075099	SKAP1
QG095-03	A10	HQP021349	NM_003718	CDC2L5
QG095-03	A11	HQP021306	NM_001008910	STK16
QG095-03	A12	HQP021305	NM_003690	PRKRA
QG095-03	B01	HQP021303	NM_003688	CASK
QG095-03	B02	HQP021299	NM_003684	MKNK1
QG095-03	B03	HQP021198	NM_003656	CAMK1
QG095-03	B04	HQP021140	NM_003639	IKBK1
QG095-03	B05	HQP021033	NM_003618	MAP4K3
QG095-03	B06	HQP020935	NM_003607	CDC42BPA
QG095-03	B07	HQP020755	NM_003585	DOC2B
QG095-03	B08	HQP020743	NM_003583	DYRK2
QG095-03	B09	HQP020734	NM_001004023	DYRK3

QG095-03	B10	HQP020596	NM_001042544	LTBP4
QG095-03	B11	HQP020476	NM_003565	ULK1
QG095-03	B12	HQP019579	NM_007371	BRD3
QG095-03	C01	HQP018812	NM_004635	MAPKAPK3
QG095-03	C02	HQP018750	NM_006301	MAP3K12
QG095-03	C03	HQP018569	NM_005433	YES1
QG095-03	C04	HQP018519	NM_003390	WEE1
QG095-03	C05	HQP018500	NM_006296	VRK2
QG095-03	C06	HQP018346	NM_006293	TYRO3
QG095-03	C07	HQP018340	NM_003331	TYK2
QG095-03	C08	HQP018334	NM_003328	TXK
QG095-03	C09	HQP018289	NM_003319	TTN
QG095-03	C10	HQP018288	NM_003318	TTK
QG095-03	C11	HQP018246	NM_007118	TRIO
QG095-03	C12	HQP018091	NM_005424	TIE1
QG095-03	D01	HQP018055	NM_003243	TGFBR3
QG095-03	D02	HQP018051	NM_004612	TGFBR1
QG095-03	D03	HQP018047	NM_003238	TGFB2
QG095-03	D04	HQP018043	NM_003236	TGFA
QG095-03	D05	HQP018019	NM_006285	TESK1
QG095-03	D06	HQP018012	NM_000459	TEK
QG095-03	D07	HQP018008	NM_003215	TEC
QG095-03	D08	HQP017891	NM_003188	MAP3K7
QG095-03	D09	HQP017873	NM_004606	TAF1
QG095-03	D10	HQP017845	NM_003177	SYK
QG095-03	D11	HQP017795	NM_001015878	AURKC
QG095-03	D12	HQP017794	NM_000455	STK11
QG095-03	E01	HQP017793	NM_005990	STK10
QG095-03	E02	HQP017791	NM_003600	AURKA
QG095-03	E03	HQP017789	NM_006282	STK4
QG095-03	E04	HQP017787	NM_003157	NEK4
QG095-03	E05	HQP017696	NM_005417	SRC
QG095-03	E06	HQP017005	NM_005627	SGK
QG095-03	E07	HQP016830	NM_003010	MAP2K4
QG095-03	E08	HQP016625	NM_002984	CCL4
QG095-03	E09	HQP016621	NM_002982	CCL2
QG095-03	E10	HQP016530	NM_001005861	RYK
QG095-03	E11	HQP016472	NM_001007071	RPS6KB2
QG095-03	E12	HQP016471	NM_003161	RPS6KB1
QG095-03	F01	HQP016470	NM_004586	RPS6KA3
QG095-03	F02	HQP016468	NM_001006932	RPS6KA2
QG095-03	F03	HQP016466	NM_001006665	RPS6KA1
QG095-03	F04	HQP016380	NM_002944	ROS1

QG095-03	F05	HQP016371	NM_005406	ROCK1
QG095-03	F06	HQP016309	NM_005104	BRD2
QG095-03	F07	HQP016303	NM_021133	RNASEL
QG095-03	F08	HQP016233	NM_006510	TRIM27
QG095-03	F09	HQP016088	NM_002880	RAF1
QG095-03	F10	HQP016084	NM_014226	RAGE
QG095-03	F11	HQP015701	NM_002821	PTK7
QG095-03	F12	HQP015692	NM_005975	PTK6
QG095-03	G01	HQP015639	NM_005607	PTK2
QG095-03	G02	HQP015174	NM_006742	PSKH1
QG095-03	G03	HQP015082	NM_001008530	LGMN
QG095-03	G04	HQP014987	NM_005044	PRKX
QG095-03	G05	HQP014948	NM_002759	EIF2AK2
QG095-03	G06	HQP014926	NM_145185	MAP2K7
QG095-03	G07	HQP014919	NM_002758	MAP2K6
QG095-03	G08	HQP014917	NM_002757	MAP2K5
QG095-03	G09	HQP014914	NM_002756	MAP2K3
QG095-03	G10	HQP014909	NM_030662	MAP2K2
QG095-03	G11	HQP014907	NM_002755	MAP2K1
QG095-03	G12	HQP014906	NM_002754	MAPK13
QG095-03	H01	HGDC		
QG095-03	H02	HGDC		
QG095-03	H03	HQP006940	NM_002046	GAPDH
QG095-03	H04	HQP016381	NM_001101	ACTB
QG095-03	H05	HQP015171	NM_004048	B2M
QG095-03	H06	HQP006171	NM_012423	RPL13A
QG095-03	H07	HQP009026	NM_000194	HPRT1
QG095-03	H08	HQP054253	NR_003286	RN18S1
QG095-03	H09	RT		
QG095-03	H10	RT		
QG095-03	H11	PCR		
QG095-03	H12	PCR		
QG095-04	A01	HQP014900	NM_002753	MAPK10
QG095-04	A02	HQP014896	NM_002752	MAPK9
QG095-04	A03	HQP014895	NM_002751	MAPK11
QG095-04	A04	HQP014886	NM_002750	MAPK8
QG095-04	A05	HQP014876	NM_002748	MAPK6
QG095-04	A06	HQP014865	NM_002747	MAPK4
QG095-04	A07	HQP014854	NM_001040056	MAPK3
QG095-04	A08	HQP014848	NM_002745	MAPK1
QG095-04	A09	HQP014847	NM_006259	PRKG2
QG095-04	A10	HQP014844	NM_006258	PRKG1
QG095-04	A11	HQP014829	NM_001033581	PRKCZ

QG095-04	A12	HQP014813	NM_001001329	PRKCSH
QG095-04	B01	HQP014804	NM_006257	PRKCQ
QG095-04	B02	HQP014798	NM_002742	PRKD1
QG095-04	B03	HQP014791	NM_006256	PKN2
QG095-04	B04	HQP014771	NM_002740	PRKCI
QG095-04	B05	HQP014760	NM_006255	PRKCH
QG095-04	B06	HQP014751	NM_002739	PRKCG
QG095-04	B07	HQP014731	NM_006254	PRKCD
QG095-04	B08	HQP014718	NM_002738	PRKCB1
QG095-04	B09	HQP014706	NM_002737	PRKCA
QG095-04	B10	HQP014687	NM_004157	PRKAR2A
QG095-04	B11	HQP014627	NM_002733	PRKAG1
QG095-04	B12	HQP014594	NM_002732	PRKACG
QG095-04	C01	HQP014575	NM_002730	PRKACA
QG095-04	C02	HQP014541	NM_006252	PRKAA2
QG095-04	C03	HQP014530	NM_006251	PRKAA1
QG095-04	C04	HQP014321	NM_002720	PPP4C
QG095-04	C05	HQP013350	NM_001001998	EXOSC10
QG095-04	C06	HQP013262	NM_005030	PLK1
QG095-04	C07	HQP013152	NM_002648	PIM1
QG095-04	C08	HQP013119	NM_000294	PHKG2
QG095-04	C09	HQP013118	NM_006213	PHKG1
QG095-04	C10	HQP013115	NM_000293	PHKB
QG095-04	C11	HQP013114	NM_000292	PHKA2
QG095-04	C12	HQP013113	NM_002637	PHKA1
QG095-04	D01	HQP013082	NM_012395	PFTK1
QG095-04	D02	HQP012979	NM_002613	PDPK1
QG095-04	D03	HQP012953	NM_002612	PDK4
QG095-04	D04	HQP012942	NM_005391	PDK3
QG095-04	D05	HQP012931	NM_002611	PDK2
QG095-04	D06	HQP012923	NM_002610	PDK1
QG095-04	D07	HQP012889	NM_002609	PDGFRB
QG095-04	D08	HQP012871	NM_006207	PDGFRL
QG095-04	D09	HQP012866	NM_006206	PDGFRA
QG095-04	D10	HQP012596	NM_002595	PCTK2
QG095-04	D11	HQP012585	NM_006201	PCTK1
QG095-04	D12	HQP012174	NM_002578	PAK3
QG095-04	E01	HQP012170	NM_002577	PAK2
QG095-04	E02	HQP012156	NM_002576	PAK1
QG095-04	E03	HQP011955	NM_001014796	DDR2
QG095-04	E04	HQP011954	NM_004560	ROR2
QG095-04	E05	HQP011946	NM_001007156	NTRK3
QG095-04	E06	HQP011938	NM_001007792	NTRK1



QG095-04	E07	HQP011902	NM_003995	NPR2
QG095-04	E08	HQP011765	NM_002498	NEK3
QG095-04	E09	HQP011764	NM_002497	NEK2
QG095-04	E10	HQP011763	NM_012224	NEK1
QG095-04	E11	HQP011760	NM_006157	NELL1
QG095-04	E12	HQP011574	NM_005592	MUSK
QG095-04	F01	HQP011525	NM_002447	MST1R
QG095-04	F02	HQP011299	NM_005372	MOS
QG095-04	F03	HQP011240	NM_002419	MAP3K11
QG095-04	F04	HQP011237	NM_002446	MAP3K10
QG095-04	F05	HQP011236	NM_033141	MAP3K9
QG095-04	F06	HQP011181	NM_000245	MET
QG095-04	F07	HQP011166	NM_005922	MAP3K4
QG095-04	F08	HQP011164	NM_002401	MAP3K3
QG095-04	F09	HQP053966	NM_005921	MAP3K1
QG095-04	F10	HQP011048	NM_002378	MATK
QG095-04	F11	HQP010994	NM_005906	MAK
QG095-04	F12	HQP010935	NM_002350	LYN
QG095-04	G01	HQP010920	NM_002344	LTK
QG095-04	G02	HQP010911	NM_000627	LTBP1
QG095-04	G03	HQP010620	NM_001031801	LIMK2
QG095-04	G04	HQP010619	NM_002314	LIMK1
QG095-04	G05	HQP010565	NM_001042771	LCK
QG095-04	G06	HQP010099	NM_000222	KIT
QG095-04	G07	HQP010070	NM_002253	KDR
QG095-04	G08	HQP009851	NM_000215	JAK3
QG095-04	G09	HQP009850	NM_004972	JAK2
QG095-04	G10	HQP009849	NM_002227	JAK1
QG095-04	G11	HQP009835	NM_005546	ITK
QG095-04	G12	HQP009776	NM_001570	IRAK2
QG095-04	H01	HGDC		
QG095-04	H02	HGDC		
QG095-04	H03	HQP006940	NM_002046	GAPDH
QG095-04	H04	HQP016381	NM_001101	ACTB
QG095-04	H05	HQP015171	NM_004048	B2M
QG095-04	H06	HQP006171	NM_012423	RPL13A
QG095-04	H07	HQP009026	NM_000194	HPRT1
QG095-04	H08	HQP054253	NR_003286	RN18S1
QG095-04	H09	RT		
QG095-04	H10	RT		
QG095-04	H11	PCR		
QG095-04	H12	PCR		
QG095-05	A01	HQP009771	NM_001025242	IRAK1

QG095-05	A02	HQP009766	NM_014215	INSRR
QG095-05	A03	HQP009764	NM_000208	INSR
QG095-05	A04	HQP009639	NM_001556	IKBKB
QG095-05	A05	HQP009523	NM_000875	IGF1R
QG095-05	A06	HQP009138	NM_004507	HUS1
QG095-05	A07	HQP008739	NM_002110	HCK
QG095-05	A08	HQP008683	NM_000180	GUCY2D
QG095-05	A09	HQP008552	NM_004963	GUCY2C
QG095-05	A10	HQP008502	NM_005316	GTF2H1
QG095-05	A11	HQP054075	NM_002093	GSK3B
QG095-05	A12	HQP008468	NM_019884	GSK3A
QG095-05	B01	HQP008273	NM_017572	MKNK2
QG095-05	B02	HQP008260	NM_005308	GRK5
QG095-05	B03	HQP008257	NM_001004056	GRK4
QG095-05	B04	HQP006741	NM_005255	GAK
QG095-05	B05	HQP006484	NM_002037	FYN
QG095-05	B06	HQP006377	NM_002031	FRK
QG095-05	B07	HQP005915	NM_002020	FLT4
QG095-05	B08	HQP005890	NM_004119	FLT3
QG095-05	B09	HQP005879	NM_002019	FLT1
QG095-05	B10	HQP005444	NM_001042729	FGR
QG095-05	B11	HQP005439	NM_002011	FGFR4
QG095-05	B12	HQP005437	NM_000141	FGFR2
QG095-05	C01	HQP005434	NM_000142	FGFR3
QG095-05	C02	HQP005395	NM_002005	FES
QG095-05	C03	HQP005394	NM_005246	FER
QG095-05	C04	HQP005093	NM_004103	PTK2B
QG095-05	C05	HQP004992	NM_001433	ERN1
QG095-05	C06	HQP004970	NM_001005915	ERBB3
QG095-05	C07	HQP004968	NM_001005862	ERBB2
QG095-05	C08	HQP004947	NM_004445	EPHB6
QG095-05	C09	HQP004945	NM_004444	EPHB4
QG095-05	C10	HQP004942	NM_004443	EPHB3
QG095-05	C11	HQP004939	NM_004442	EPHB2
QG095-05	C12	HQP004930	NM_004440	EPHA7
QG095-05	D01	HQP004926	NM_004438	EPHA4
QG095-05	D02	HQP004924	NM_005233	EPHA3
QG095-05	D03	HQP004922	NM_005232	EPHA1
QG095-05	D04	HQP004657	NM_004431	EPHA2
QG095-05	D05	HQP004605	NM_005228	EGFR
QG095-05	D06	HQP004598	NM_001406	EFNB3
QG095-05	D07	HQP004592	NM_005227	EFNA4
QG095-05	D08	HQP004591	NM_004952	EFNA3

QG095-05	D09	HQP003982	NM_001348	DAPK3
QG095-05	D10	HQP003979	NM_004938	DAPK1
QG095-05	D11	HQP003285	NM_001896	CSNK2A2
QG095-05	D12	HQP003277	NM_001895	CSNK2A1
QG095-05	E01	HQP003257	NM_001319	CSNK1G2
QG095-05	E02	HQP003249	NM_001894	CSNK1E
QG095-05	E03	HQP003240	NM_001893	CSNK1D
QG095-05	E04	HQP003236	NM_001025105	CSNK1A1
QG095-05	E05	HQP003218	NM_004383	CSK
QG095-05	E06	HQP003158	NM_005211	CSF1R
QG095-05	E07	HQP003133	NM_001315	MAPK14
QG095-05	E08	HQP002989	NM_005207	CRKL
QG095-05	E09	HQP002725	NM_005204	MAP3K8
QG095-05	E10	HQP002379	NM_004073	PLK3
QG095-05	E11	HQP001894	NM_001827	CKS2
QG095-05	E12	HQP001884	NM_001826	CKS1B
QG095-05	F01	HQP001708	NM_001278	CHUK
QG095-05	F02	HQP001282	NM_001274	CHEK1
QG095-05	F03	HQP000331	NM_000389	CDKN1A
QG095-05	F04	HQP000321	NM_001261	CDK9
QG095-05	F05	HQP000285	NM_001799	CDK7
QG095-05	F06	HQP000274	NM_001259	CDK6
QG095-05	F07	HQP000261	NM_004935	CDK5
QG095-05	F08	HQP000245	NM_000075	CDK4
QG095-05	F09	HQP000225	NM_001798	CDK2
QG095-05	F10	HQP023261	NM_001786	CDC2
QG095-05	F11	HQP021906	NM_001239	CCNH
QG095-05	F12	HQP021757	NM_001760	CCND3
QG095-05	G01	HQP019986	NM_001221	CAMK2D
QG095-05	G02	HQP019959	NM_001220	CAMK2B
QG095-05	G03	HQP019917	NM_015981	CAMK2A
QG095-05	G04	HQP019874	NM_001744	CAMK4
QG095-05	G05	HQP019707	NM_001743	CALM2
QG095-05	G06	HQP018928	NM_004341	CAD
QG095-05	G07	HQP018768	NM_001954	DDR1
QG095-05	G08	HQP018022	NM_001211	BUB1B
QG095-05	G09	HQP017999	NM_004336	BUB1
QG095-05	G10	HQP017988	NM_000061	BTK
QG095-05	G11	HQP017733	NM_004333	BRAF
QG095-05	G12	HQP017566	NM_001721	BMX
QG095-05	H01	HGDC		
QG095-05	H02	HGDC		
QG095-05	H03	HQP006940	NM_002046	GAPDH

QG095-05	H04	HQP016381	NM_001101	ACTB
QG095-05	H05	HQP015171	NM_004048	B2M
QG095-05	H06	HQP006171	NM_012423	RPL13A
QG095-05	H07	HQP009026	NM_000194	HPRT1
QG095-05	H08	HQP054253	NR_003286	RN18S1
QG095-05	H09	RT		
QG095-05	H10	RT		
QG095-05	H11	PCR		
QG095-05	H12	PCR		
QG095-06	A01	HQP017535	NM_001204	BMPR2
QG095-06	A02	HQP017500	NM_001203	BMPR1B
QG095-06	A03	HQP017489	NM_004329	BMPR1A
QG095-06	A04	HQP016783	NM_001715	BLK
QG095-06	A05	HQP016418	NM_004327	BCR
QG095-06	A06	HQP016204	NM_053056	CCND1
QG095-06	A07	HQP014815	NM_001699	AXL
QG095-06	A08	HQP013596	NM_001184	ATR
QG095-06	A09	HQP011736	NM_000051	ATM
QG095-06	A10	HQP009832	NM_001654	ARAF
QG095-06	A11	HQP007421	NM_020547	AMHR2
QG095-06	A12	HQP006353	NM_004304	ALK
QG095-06	B01	HQP004995	NM_001626	AKT2
QG095-06	B02	HQP004991	NM_001014431	AKT1
QG095-06	B03	HQP003848	NM_005160	ADRBK2
QG095-06	B04	HQP003816	NM_001619	ADRBK1
QG095-06	B05	HQP022834	NM_000020	ACVRL1
QG095-06	B06	HQP022668	NM_001106	ACVR2B
QG095-06	B07	HQP022491	NM_001616	ACVR2A
QG095-06	B08	HQP022297	NM_004302	ACVR1B
QG095-06	B09	HQP022071	NM_001105	ACVR1
QG095-06	B10	HQP006954	NM_005157	ABL1
QG095-06	B11	HQP004933	NM_004441	EPHB1
QG095-06	B12	HQP014583	NM_002731	PRKACB
QG095-06	C01	HQP017725	NM_182691	SRPK2
QG095-06	C02	HQP020636	NM_001032296	STK24
QG095-06	C03	HQP021896	NM_003957	BRSK2
QG095-06	C04	HQP022182	NM_138370	LOC91461
QG095-06	C05	HQP021807	NM_001001874	TPD52L3
QG095-06	C06	HQP020433	NM_052843	OBSCN
QG095-06	C07	HQP020387	NM_031965	GSG2
QG095-06	C08	HQP014698	NM_018343	RIOK2
QG095-06	C09	HQP007785	NM_005158	ABL2
QG095-06	C10	HQP015008	NM_031272	TEX14

QG095-06	C11	HQP020019	NM_001222	CAMK2G
QG095-06	C12	HQP023275	NM_001787	CDC2L1
QG095-06	D01	HQP000234	NM_001258	CDK3
QG095-06	D02	HQP000312	NM_001260	CDK8
QG095-06	D03	HQP000382	NM_004936	CDKN2B
QG095-06	D04	HQP001845	NM_001825	CKMT2
QG095-06	D05	HQP002051	NM_001024646	CLK1
QG095-06	D06	HQP002068	NM_001292	CLK3
QG095-06	D07	HQP003259	NM_001031812	CSNK1G3
QG095-06	D08	HQP003289	NM_001320	CSNK2B
QG095-06	D09	HQP004379	NM_001081560	DMPK
QG095-06	D10	HQP016322	NM_021826	FASTKD5
QG095-06	D11	HQP022007	NM_004672	MAP3K6
QG095-06	D12	HQP004928	NM_004439	EPHA5
QG095-06	E01	HQP004931	NM_001006943	EPHA8
QG095-06	E02	HQP016685	NM_022076	DUSP21
QG095-06	E03	HQP005472	NM_000801	FKBP1A
QG095-06	E04	HQP001147	NM_006852	TLK2
QG095-06	E05	HQP001371	NM_001042600	MAP4K1
QG095-06	E06	HQP011031	NM_002376	MARK3
QG095-06	E07	HQP011168	NM_005923	MAP3K5
QG095-06	E08	HQP018931	NM_001006617	MAPKAP1
QG095-06	E09	HQP011941	NM_001007097	NTRK2
QG095-06	E10	HQP006208	NM_001039803	CCRK
QG095-06	E11	HQP012609	NM_002596	PCTK3
QG095-06	E12	HQP014742	NM_005400	PRKCE
QG095-06	F01	HQP014782	NM_002741	PKN1
QG095-06	F02	HQP014837	NM_001081640	PRKDC
QG095-06	F03	HQP014877	NM_002749	MAPK7
QG095-06	F04	HQP019328	NM_024800	NEK11
QG095-06	F05	HQP006316	NM_001033578	SGK3
QG095-06	F06	HQP016055	NM_004579	MAP4K2
QG095-06	F07	HQP016220	NM_020630	RET
QG095-06	F08	HQP019524	NM_001018046	YSK4
QG095-06	F09	HQP006801	NM_001079880	PRKD2
QG095-06	F10	HQP020310	NM_031480	RIOK1
QG095-06	F11	HQP008313	NM_021643	TRIB2
QG095-06	F12	HQP017788	NM_006281	STK3
QG095-06	G01	HQP017792	NM_001037343	CDKL5
QG095-06	G02	HQP018053	NM_001024847	TGFBR2
QG095-06	G03	HQP018499	NM_003384	VRK1
QG095-06	G04	HQP018578	NM_207519	ZAP70
QG095-06	G05	HQP021267	NM_003668	MAPKAPK5

QG095-06	G06	HQP012135	NM_001024937	MINK1
QG095-06	G07	HQP021457	NM_003985	TNK1
QG095-06	G08	HQP021558	NM_003845	DYRK4
QG095-06	G09	HQP021595	NM_001024628	NRP1
QG095-06	G10	HQP013335	NM_017433	MYO3A
QG095-06	G11	HQP021799	NM_001006944	RPS6KA4
QG095-06	G12	HQP053901	NM_017886	ULK4
QG095-06	H01	HGDC		
QG095-06	H02	HGDC		
QG095-06	H03	HQP006940	NM_002046	GAPDH
QG095-06	H04	HQP016381	NM_001101	ACTB
QG095-06	H05	HQP015171	NM_004048	B2M
QG095-06	H06	HQP006171	NM_012423	RPL13A
QG095-06	H07	HQP009026	NM_000194	HPRT1
QG095-06	H08	HQP054253	NR_003286	RN18S1
QG095-06	H09	RT		
QG095-06	H10	RT		
QG095-06	H11	PCR		
QG095-06	H12	PCR		