

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG049-01	A01	HQP005430	NM_023107	FGFR1
QG049-01	A02	HQP004513	NM_181870	DVL1
QG049-01	A03	HQP002581	NM_078485	COL9A1
QG049-01	A04	HQP022974	NM_001001390	CD44
QG049-01	A05	HQP009026	NM_000194	HPRT1
QG049-01	A06	HQP000471	NM_001025107	ADAR
QG049-01	A07	HQP018528	NM_005430	WNT1
QG049-01	A08	HQP017627	NM_005986	SOX1
QG049-01	A09	HQP013634	NM_015869	PPARG
QG049-01	A10	HQP013627	NM_006238	PPARD
QG049-01	A11	HQP000343	NM_005866	OPRS1
QG049-01	A12	HQP011875	NM_024408	NOTCH2
QG049-01	B01	HQP020510	NM_032188	MYST1
QG049-01	B02	HQP010168	NM_002275	KRT15
QG049-01	B03	HQP009088	NM_004134	HSPA9
QG049-01	B04	HQP008746	NM_001527	HDAC2
QG049-01	B05	HQP007477	NM_004004	GJB2
QG049-01	B06	HQP007316	NM_016204	GDF2
QG049-01	B07	HQP007253	NM_021078	GCN5L2
QG049-01	B08	HQP020133	NM_003505	FZD1
QG049-01	B09	HQP000042	NM_005479	FRAT1
QG049-01	B10	HQP004897	NM_001429	EP300
QG049-01	B11	HQP004512	NM_004421	DVL1
QG049-01	B12	HQP001642	NM_020892	DTX2
QG049-01	C01	HQP004495	NM_004416	DTX1
QG049-01	C02	HQP000828	NM_016941	DLL3
QG049-01	C03	HQP008111	NM_005618	DLL1
QG049-01	C04	HQP003515	NM_001903	CTNNA1
QG049-01	C05	HQP000018	NM_001792	CDH2
QG049-01	C06	HQP053981	NM_001791	CDC42
QG049-01	C07	HQP023261	NM_001786	CDC2
QG049-01	C08	HQP022438	NM_004931	CD8B
QG049-01	C09	HQP022212	NM_000732	CD3D
QG049-01	C10	HQP021819	NM_001238	CCNE1
QG049-01	C11	HQP021701	NM_001237	CCNA2
QG049-01	C12	HQP021753	NM_033637	BTRC
QG049-01	D01	HQP017357	NM_001201	BMP3
QG049-01	D02	HQP017282	NM_006129	BMP1
QG049-01	D03	HQP016599	NM_199173	BGLAP
QG049-01	D04	HQP020114	NM_003502	AXIN1
QG049-01	D05	HQP006433	NM_001631	ALPI

QG049-01	D06	HQP005087	NM_000690	ALDH2
QG049-01	D07	HQP005075	NM_000689	ALDH1A1
QG049-01	D08	HQP000472	NM_001111	ADAR
QG049-01	D09	HQP018117	NM_005159	ACTC1
QG049-01	D10	HQP018519	NM_003390	WEE1
QG049-01	D11	HQP018489	NM_003380	VIM
QG049-01	D12	HQP018475	NM_001025366	VEGFA
QG049-01	E01	HQP018474	NM_000376	VDR
QG049-01	E02	HQP018466	NM_080682	VCAM1
QG049-01	E03	HQP018388	NM_000462	UBE3A
QG049-01	E04	HQP000446	NM_006086	TUBB3
QG049-01	E05	HQP023250	NM_014788	TRIM14
QG049-01	E06	HQP021361	NM_003722	TP73L
QG049-01	E07	HQP018175	NM_000546	TP53
QG049-01	E08	HQP021502	NM_003810	TNFSF10
QG049-01	E09	HQP018148	NM_001065	TNFRSF1A
QG049-01	E10	HQP018141	NM_000594	TNF
QG049-01	E11	HQP009140	NM_002160	TNC
QG049-01	E12	HQP018116	NM_138554	TLR4
QG049-01	F01	HQP018092	NM_003254	TIMP1
QG049-01	F02	HQP018083	NM_006288	THY1
QG049-01	F03	HQP018074	NM_000460	THPO
QG049-01	F04	HQP018047	NM_003238	TGFB2
QG049-01	F05	HQP018044	NM_000660	TGFB1
QG049-01	F06	HQP018041	NM_003234	TFRC
QG049-01	F07	HQP019038	NM_007162	TFEB
QG049-01	F08	HQP018027	NM_003222	TFAP2C
QG049-01	F09	HQP018023	NM_001032280	TFAP2A
QG049-01	F10	HQP018012	NM_000459	TEK
QG049-01	F11	HQP020838	NM_003598	TEAD2
QG049-01	F12	HQP017997	NM_003212	TDGF1
QG049-01	G01	HQP019839	NM_021966	TCL1A
QG049-01	G02	HQP020164	NM_031283	TCF7L1
QG049-01	G03	HQP017957	NM_003200	TCF3
QG049-01	G04	HQP017945	NM_003196	TCEA3
QG049-01	G05	HQP017895	NM_003189	TAL1
QG049-01	G06	HQP017877	NM_001001522	TAGLN
QG049-01	G07	HQP000554	NM_006342	TACC3
QG049-01	G08	HQP017857	NM_003181	T
QG049-01	G09	HQP017850	NM_003179	SYP
QG049-01	G10	HQP017810	NM_001055	SULT1A1
QG049-01	G11	HQP017774	NM_012448	STAT5B
QG049-01	G12	HQP017771	NM_003152	STAT5A

QG049-01	H01	HGDC		
QG049-01	H02	HGDC		
QG049-01	H03	HQP006940	NM_002046	GAPDH
QG049-01	H04	HQP016381	NM_001101	ACTB
QG049-01	H05	HQP015171	NM_004048	B2M
QG049-01	H06	HQP006171	NM_012423	RPL13A
QG049-01	H07	HQP009026	NM_000194	HPRT1
QG049-01	H08	HQP054253	NR_003286	RN18S1
QG049-01	H09	RT		
QG049-01	H10	RT		
QG049-01	H11	PCR		
QG049-01	H12	PCR		
QG049-02	A01	HQP017767	NM_003150	STAT3
QG049-02	A02	HQP053963	NM_001048	SST
QG049-02	A03	HQP017703	NM_001005291	SREBF1
QG049-02	A04	HQP017673	NM_000582	SPP1
QG049-02	A05	HQP017670	NM_001030288	SPN
QG049-02	A06	HQP017662	NM_001080547	SPI1
QG049-02	A07	HQP017635	NM_000346	SOX9
QG049-02	A08	HQP014460	NM_017508	SOX6
QG049-02	A09	HQP017628	NM_003106	SOX2
QG049-02	A10	HQP016906	NM_022454	SOX17
QG049-02	A11	HQP021889	NM_003955	SOCS3
QG049-02	A12	HQP017512	NM_003072	SMARCA4
QG049-02	B01	HQP017353	NM_006931	SLC2A3
QG049-02	B02	HQP012076	NM_002555	SLC22A18
QG049-02	B03	HQP017303	NM_003037	SLAMF1
QG049-02	B04	HQP002151	NM_001040153	SLAIN1
QG049-02	B05	HQP017259	NM_006747	SIPA1
QG049-02	B06	HQP017102	NM_004169	SHMT1
QG049-02	B07	HQP017098	NM_000193	SHH
QG049-02	B08	HQP017080	NM_003029	SHC1
QG049-02	B09	HQP000726	NM_020979	SH2B2
QG049-02	B10	HQP016874	NM_003013	SFRP2
QG049-02	B11	HQP016865	NM_003012	SFRP1
QG049-02	B12	HQP016744	NM_000450	SELE
QG049-02	C01	HQP019428	NM_001037582	SCD5
QG049-02	C02	HQP015461	NM_020436	SALL4
QG049-02	C03	HQP016564	NM_002968	SALL1
QG049-02	C04	HQP016553	NM_006272	S100B
QG049-02	C05	HQP016542	NM_014624	S100A6
QG049-02	C06	HQP016540	NM_002961	S100A4
QG049-02	C07	HQP016478	NM_001015051	RUNX2

QG049-02	C08	HQP022404	NM_004755	RPS6KA5
QG049-02	C09	HQP016371	NM_005406	ROCK1
QG049-02	C10	HQP016307	NM_007212	RNF2
QG049-02	C11	HQP016131	NM_000321	RB1
QG049-02	C12	HQP016118	NM_000966	RARG
QG049-02	D01	HQP016063	NM_006908	RAC1
QG049-02	D02	HQP016047	NM_002867	RAB3B
QG049-02	D03	HQP006039	NM_015317	PUM2
QG049-02	D04	HQP015921	NM_002840	PTPRF
QG049-02	D05	HQP015908	NM_002838	PTPRC
QG049-02	D06	HQP015878	NM_002834	PTPN11
QG049-02	D07	HQP015535	NM_000314	PTEN
QG049-02	D08	HQP015530	NM_000264	PTCH1
QG049-02	D09	HQP015123	NM_000021	PSEN1
QG049-02	D10	HQP003586	NM_144707	PROM2
QG049-02	D11	HQP021613	NM_006017	PROM1
QG049-02	D12	HQP013717	NM_000943	PPIC
QG049-02	E01	HQP013633	NM_005037	PPARG
QG049-02	E02	HQP013600	NM_002701	POU5F1
QG049-02	E03	HQP013416	NM_001018111	PODXL
QG049-02	E04	HQP013201	NM_000930	PLAT
QG049-02	E05	HQP014077	NM_018068	PIWIL2
QG049-02	E06	HQP012632	NM_016621	PHF21A
QG049-02	E07	HQP013015	NM_000442	PECAM1
QG049-02	E08	HQP012856	NM_002608	PDGFB
QG049-02	E09	HQP018681	NM_007144	PCGF2
QG049-02	E10	HQP012219	NM_000280	PAX6
QG049-02	E11	HQP003171	NM_006437	PARP4
QG049-02	E12	HQP012140	NM_000430	PAFAH1B1
QG049-02	F01	HQP012100	NM_002562	P2RX7
QG049-02	F02	HQP012085	NM_021728	OTX2
QG049-02	F03	HQP021395	NM_001005743	NUMB
QG049-02	F04	HQP011968	NM_006186	NR4A2
QG049-02	F05	HQP011873	NM_017617	NOTCH1
QG049-02	F06	HQP011861	NM_018055	NODAL
QG049-02	F07	HQP016801	NM_022162	NOD2
QG049-02	F08	HQP021210	NM_033120	NKD2
QG049-02	F09	HQP011807	NM_003998	NFKB1
QG049-02	F10	HQP000899	NM_006617	NES
QG049-02	F11	HQP019390	NM_024865	NANOG
QG049-02	F12	HQP011605	NM_005593	MYF5
QG049-02	G01	HQP011597	NM_002467	MYC
QG049-02	G02	HQP011555	NM_002454	MTRR

QG049-02	G03	HQP011547	NM_005957	MTHFR
QG049-02	G04	HQP011491	NM_000251	MSH2
QG049-02	G05	HQP022527	NM_021950	MS4A1
QG049-02	G06	HQP011308	NM_005373	MPL
QG049-02	G07	HQP011263	NM_004994	MMP9
QG049-02	G08	HQP011254	NM_000902	MME
QG049-02	G09	HQP011223	NM_000248	MITF
QG049-02	G10	HQP011179	NM_002402	MEST
QG049-02	G11	HQP011171	NM_000244	MEN1
QG049-02	G12	HQP023256	NM_014791	MELK
QG049-02	H01	HGDC		
QG049-02	H02	HGDC		
QG049-02	H03	HQP006940	NM_002046	GAPDH
QG049-02	H04	HQP016381	NM_001101	ACTB
QG049-02	H05	HQP015171	NM_004048	B2M
QG049-02	H06	HQP006171	NM_012423	RPL13A
QG049-02	H07	HQP009026	NM_000194	HPRT1
QG049-02	H08	HQP054253	NR_003286	RN18S1
QG049-02	H09	RT		
QG049-02	H10	RT		
QG049-02	H11	PCR		
QG049-02	H12	PCR		
QG049-03	A01	HQP021938	NM_139279	MCFD2
QG049-03	A02	HQP011077	NM_000242	MBL2
QG049-03	A03	HQP007713	NM_013283	MAT2B
QG049-03	A04	HQP011047	NM_005911	MAT2A
QG049-03	A05	HQP003133	NM_001315	MAPK14
QG049-03	A06	HQP053966	NM_005921	MAP3K1
QG049-03	A07	HQP014907	NM_002755	MAP2K1
QG049-03	A08	HQP014450	NM_018717	MAML3
QG049-03	A09	HQP010928	NM_002346	LY6E
QG049-03	A10	HQP015246	NM_020169	LXN
QG049-03	A11	HQP006969	NM_015541	LRIG1
QG049-03	A12	HQP010608	NM_002310	LIFR
QG049-03	B01	HQP010607	NM_002309	LIF
QG049-03	B02	HQP022587	NM_004789	LHX2
QG049-03	B03	HQP010581	NM_000230	LEP
QG049-03	B04	HQP018049	NM_003240	LEFTY2
QG049-03	B05	HQP008000	NM_005559	LAMA1
QG049-03	B06	HQP010233	NM_002276	KRT19
QG049-03	B07	HQP010186	NM_000422	KRT17
QG049-03	B08	HQP011204	NM_000899	KITLG
QG049-03	B09	HQP010099	NM_000222	KIT

QG049-03	B10	HQP010070	NM_002253	KDR
QG049-03	B11	HQP016008	NM_021219	JAM2
QG049-03	B12	HQP009850	NM_004972	JAK2
QG049-03	C01	HQP009847	NM_002226	JAG2
QG049-03	C02	HQP004470	NM_000214	JAG1
QG049-03	C03	HQP009815	NM_000211	ITGB2
QG049-03	C04	HQP009810	NM_002211	ITGB1
QG049-03	C05	HQP009807	NM_000632	ITGAM
QG049-03	C06	HQP009774	NM_000210	ITGA6
QG049-03	C07	HQP009795	NM_000419	ITGA2B
QG049-03	C08	HQP009794	NM_002203	ITGA2
QG049-03	C09	HQP009793	NM_181501	ITGA1
QG049-03	C10	HQP009791	NM_002202	ISL1
QG049-03	C11	HQP009744	NM_002193	INHBB
QG049-03	C12	HQP009676	NM_000880	IL7
QG049-03	D01	HQP009674	NM_002184	IL6ST
QG049-03	D02	HQP009670	NM_000600	IL6
QG049-03	D03	HQP009666	NM_000879	IL5
QG049-03	D04	HQP009662	NM_000589	IL4
QG049-03	D05	HQP009649	NM_000586	IL2
QG049-03	D06	HQP009708	NM_000585	IL15
QG049-03	D07	HQP009693	NM_002187	IL12B
QG049-03	D08	HQP009692	NM_000882	IL12A
QG049-03	D09	HQP009685	NM_000572	IL10
QG049-03	D10	HQP009532	NM_000876	IGF2R
QG049-03	D11	HQP009529	NM_000612	IGF2
QG049-03	D12	HQP009518	NM_000618	IGF1
QG049-03	E01	HQP009467	NM_000619	IFNG
QG049-03	E02	HQP009184	NM_000201	ICAM1
QG049-03	E03	HQP009107	NM_005529	HSPG2
QG049-03	E04	HQP009089	NM_001540	HSPB1
QG049-03	E05	HQP009077	NM_005345	HSPA1A
QG049-03	E06	HQP018231	NM_003299	HSP90B1
QG049-03	E07	HQP008990	NM_024015	HOXB4
QG049-03	E08	HQP008980	NM_152739	HOXA9
QG049-03	E09	HQP008981	NM_018951	HOXA10
QG049-03	E10	HQP008894	NM_002131	HMGA1
QG049-03	E11	HQP008831	NM_001530	HIF1A
QG049-03	E12	HQP008800	NM_000601	HGF
QG049-03	F01	HQP013609	NM_019089	HES2
QG049-03	F02	HQP009057	NM_005524	HES1
QG049-03	F03	HQP023134	NM_014707	HDAC9
QG049-03	F04	HQP008745	NM_004964	HDAC1

QG049-03	F05	HQP008469	NM_002093	GSK3B
QG049-03	F06	HQP007599	NM_018960	GNMT
QG049-03	F07	HQP007762	NM_000172	GNAT1
QG049-03	F08	HQP007701	NM_005269	GLI1
QG049-03	F09	HQP007467	NM_000166	GJB1
QG049-03	F10	HQP007411	NM_000165	GJA1
QG049-03	F11	HQP007385	NM_000515	GH1
QG049-03	F12	HQP007346	NM_000514	GDNF
QG049-03	G01	HQP000275	NM_005811	GDF11
QG049-03	G02	HQP021145	NM_003643	GCM1
QG049-03	G03	HQP007196	NM_005257	GATA6
QG049-03	G04	HQP007183	NM_002052	GATA4
QG049-03	G05	HQP007165	NM_001002295	GATA3
QG049-03	G06	HQP007152	NM_032638	GATA2
QG049-03	G07	HQP006628	NM_000814	GABRB3
QG049-03	G08	HQP006580	NM_002040	GABPA
QG049-03	G09	HQP006515	NM_000402	G6PD
QG049-03	G10	HQP020136	NM_003507	FZD7
QG049-03	G11	HQP006455	NM_002033	FUT4
QG049-03	G12	HQP006426	NM_004958	FRAP1
QG049-03	H01	HGDC		
QG049-03	H02	HGDC		
QG049-03	H03	HQP006940	NM_002046	GAPDH
QG049-03	H04	HQP016381	NM_001101	ACTB
QG049-03	H05	HQP015171	NM_004048	B2M
QG049-03	H06	HQP006171	NM_012423	RPL13A
QG049-03	H07	HQP009026	NM_000194	HPRT1
QG049-03	H08	HQP054253	NR_003286	RN18S1
QG049-03	H09	RT		
QG049-03	H10	RT		
QG049-03	H11	PCR		
QG049-03	H12	PCR		
QG049-04	A01	HQP006231	NM_002029	FPR1
QG049-04	A02	HQP012269	NM_014009	FOXP3
QG049-04	A03	HQP005759	NM_001455	FOXO3A
QG049-04	A04	HQP005663	NM_005250	FOXL1
QG049-04	A05	HQP005335	NM_001037165	FOXK1
QG049-04	A06	HQP008906	NM_021784	FOXA2
QG049-04	A07	HQP006188	NM_005252	FOS
QG049-04	A08	HQP006144	NM_000802	FOLR1
QG049-04	A09	HQP006022	NM_002026	FN1
QG049-04	A10	HQP005890	NM_004119	FLT3
QG049-04	A11	HQP005879	NM_002019	FLT1

QG049-04	A12	HQP005792	NM_002017	FLI1
QG049-04	B01	HQP004797	NM_144606	FLCN
QG049-04	B02	HQP005434	NM_000142	FGFR3
QG049-04	B03	HQP005411	NM_002009	FGF7
QG049-04	B04	HQP005408	NM_004464	FGF5
QG049-04	B05	HQP005405	NM_002007	FGF4
QG049-04	B06	HQP005403	NM_002006	FGF2
QG049-04	B07	HQP005400	NM_000800	FGF1
QG049-04	B08	HQP005196	NM_001004019	FBLN2
QG049-04	B09	HQP005106	NM_001996	FBLN1
QG049-04	B10	HQP009671	NM_000639	FASLG
QG049-04	B11	HQP009651	NM_000043	FAS
QG049-04	B12	HQP005052	NM_000506	F2
QG049-04	C01	HQP005050	NM_004456	EZH2
QG049-04	C02	HQP005020	NM_004454	ETV5
QG049-04	C03	HQP004998	NM_000125	ESR1
QG049-04	C04	HQP004939	NM_004442	EPHB2
QG049-04	C05	HQP004657	NM_004431	EPHA2
QG049-04	C06	HQP004903	NM_001430	EPAS1
QG049-04	C07	HQP020132	NM_005442	EOMES
QG049-04	C08	HQP004818	NM_001423	EMP1
QG049-04	C09	HQP018897	NM_024090	ELOVL6
QG049-04	C10	HQP004616	NM_004430	EGR3
QG049-04	C11	HQP004605	NM_005228	EGFR
QG049-04	C12	HQP004524	NM_005225	E2F1
QG049-04	D01	HQP014176	NM_018189	DPPA4
QG049-04	D02	HQP003652	NM_138815	DPPA2
QG049-04	D03	HQP008616	NM_013369	DNMT3L
QG049-04	D04	HQP004411	NM_022552	DNMT3A
QG049-04	D05	HQP023200	NM_014750	DLG7
QG049-04	D06	HQP007526	NM_001018057	DKK3
QG049-04	D07	HQP004334	NM_001363	DKC1
QG049-04	D08	HQP004196	NM_001927	DES
QG049-04	D09	HQP003859	NM_000780	CYP7A1
QG049-04	D10	HQP003919	NM_000783	CYP26A1
QG049-04	D11	HQP018802	NM_001008540	CXCR4
QG049-04	D12	HQP016669	NM_000609	CXCL12
QG049-04	E01	HQP003539	NM_001904	CTNNB1
QG049-04	E02	HQP003185	NM_000760	CSF3R
QG049-04	E03	HQP003173	NM_000759	CSF3
QG049-04	E04	HQP003159	NM_000758	CSF2
QG049-04	E05	HQP003158	NM_005211	CSF1R
QG049-04	E06	HQP003149	NM_000757	CSF1



QG049-04	E07	HQP002867	NM_001031847	CPT1A
QG049-04	E08	HQP002478	NM_000090	COL3A1
QG049-04	E09	HQP002473	NM_001844	COL2A1
QG049-04	E10	HQP002462	NM_000088	COL1A1
QG049-04	E11	HQP004198	NM_001928	CFD
QG049-04	E12	HQP000615	NM_004364	CEBPA
QG049-04	F01	HQP000553	NM_001265	CDX2
QG049-04	F02	HQP000396	NM_001262	CDKN2C
QG049-04	F03	HQP000369	NM_000077	CDKN2A
QG049-04	F04	HQP000342	NM_004064	CDKN1B
QG049-04	F05	HQP000331	NM_000389	CDKN1A
QG049-04	F06	HQP000052	NM_001795	CDH5
QG049-04	F07	HQP023454	NM_001039802	CDC42
QG049-04	F08	HQP000279	NM_005816	CD96
QG049-04	F09	HQP022474	NM_001769	CD9
QG049-04	F10	HQP023083	NM_001040059	CD68
QG049-04	F11	HQP022972	NM_000610	CD44
QG049-04	F12	HQP022316	NM_000616	CD4
QG049-04	G01	HQP022812	NM_001025109	CD34
QG049-04	G02	HQP022504	NM_001770	CD19
QG049-04	G03	HQP021754	NM_001759	CCND2
QG049-04	G04	HQP016204	NM_053056	CCND1
QG049-04	G05	HQP016626	NM_002985	CCL5
QG049-04	G06	HQP021518	NM_000071	CBS
QG049-04	G07	HQP020297	NM_004346	CASP3
QG049-04	G08	HQP006321	NM_014333	CADM1
QG049-04	G09	HQP018071	NM_004053	BYSL
QG049-04	G10	HQP017713	NM_007294	BRCA1
QG049-04	G11	HQP017535	NM_001204	BMPR2
QG049-04	G12	HQP017396	NM_001202	BMP4
QG049-04	H01	HGDC		
QG049-04	H02	HGDC		
QG049-04	H03	HQP006940	NM_002046	GAPDH
QG049-04	H04	HQP016381	NM_001101	ACTB
QG049-04	H05	HQP015171	NM_004048	B2M
QG049-04	H06	HQP006171	NM_012423	RPL13A
QG049-04	H07	HQP009026	NM_000194	HPRT1
QG049-04	H08	HQP054253	NR_003286	RN18S1
QG049-04	H09	RT		
QG049-04	H10	RT		
QG049-04	H11	PCR		
QG049-04	H12	PCR		
QG049-05	A01	HQP017333	NM_001200	BMP2

QG049-05	A02	HQP017234	NM_005180	BMI1
QG049-05	A03	HQP006329	NM_017614	BHMT2
QG049-05	A04	HQP016418	NM_004327	BCR
QG049-05	A05	HQP016211	NM_000633	BCL2
QG049-05	A06	HQP010924	NM_001668	ARNT
QG049-05	A07	HQP009578	NM_000484	APP
QG049-05	A08	HQP009024	NM_000038	APC
QG049-05	A09	HQP008829	NM_001154	ANXA5
QG049-05	A10	HQP008414	NM_001150	ANPEP
QG049-05	A11	HQP006440	NM_000478	ALPL
QG049-05	A12	HQP004991	NM_001014431	AKT1
QG049-05	B01	HQP004494	NM_000029	AGT
QG049-05	B02	HQP004351	NM_001134	AFP
QG049-05	B03	HQP022625	NM_004797	ADIPOQ
QG049-05	B04	HQP004081	NM_000789	ACE
QG049-05	B05	HQP004388	NM_001135	ACAN
QG049-05	B06	HQP022745	NM_004827	ABCG2
QG049-05	B07	HQP013100	NM_000927	ABCB1
QG049-05	B08	HQP000567	NM_005193	CDX4
QG049-05	B09	HQP004956	NM_000799	EPO
QG049-05	B10	HQP000564	NM_006350	FST
QG049-05	B11	HQP007229	NM_001485	GBX2
QG049-05	B12	HQP009680	NM_000590	IL9
QG049-05	C01	HQP015279	NM_020203	MEPE
QG049-05	C02	HQP011654	NM_002478	MYOD1
QG049-05	C03	HQP004563	NM_000475	NR0B1
QG049-05	C04	HQP018119	NM_003269	NR2E1
QG049-05	C05	HQP002580	NM_001851	COL9A1
QG049-05	C06	HQP011526	NM_002448	MSX1
QG049-05	C07	HQP016727	NM_024019	NEUROG2
QG049-05	C08	HQP001925	NM_001117	ADCYAP1
QG049-05	C09	HQP008716	NM_001002857	ANXA2
QG049-05	C10	HQP020946	NM_001752	CAT
QG049-05	C11	HQP002245	NM_181657	LTB4R
QG049-05	C12	HQP002852	NM_001307	CLDN7
QG049-05	D01	HQP003933	NM_001343	DAB2
QG049-05	D02	HQP004407	NM_001379	DNMT1
QG049-05	D03	HQP004414	NM_006892	DNMT3B
QG049-05	D04	HQP009699	NM_199286	DPPA3
QG049-05	D05	HQP009293	NM_001025290	DPPA5
QG049-05	D06	HQP005422	NM_004114	FGF13
QG049-05	D07	HQP005404	NM_005247	FGF3
QG049-05	D08	HQP005427	NM_015850	FGFR1

QG049-05	D09	HQP005902	NM_001459	FLT3LG
QG049-05	D10	HQP007441	NM_012183	FOXD3
QG049-05	D11	HQP022926	NM_020634	GDF3
QG049-05	D12	HQP007430	NM_021954	GJA3
QG049-05	E01	HQP008607	NM_002099	GYPA
QG049-05	E02	HQP008703	NM_002105	H2AFX
QG049-05	E03	HQP022724	NM_004821	HAND1
QG049-05	E04	HQP008992	NM_018952	HOXB6
QG049-05	E05	HQP009688	NM_000641	IL11
QG049-05	E06	HQP009660	NM_000588	IL3
QG049-05	E07	HQP000436	NM_016270	KLF2
QG049-05	E08	HQP011688	NM_000615	NCAM1
QG049-05	E09	HQP011771	NM_002500	NEUROD1
QG049-05	E10	HQP011823	NM_002505	NFYA
QG049-05	E11	HQP007259	NM_001489	NR6A1
QG049-05	E12	HQP013880	NM_001079839	OCIAD1
QG049-05	F01	HQP009769	NM_000209	PDX1
QG049-05	F02	HQP001583	NM_007273	PHB2
QG049-05	F03	HQP004566	NM_004426	PHC1
QG049-05	F04	HQP013174	NM_000325	PITX2
QG049-05	F05	HQP022448	NM_004764	PIWIL1
QG049-05	F06	HQP019461	NM_005672	PSCA
QG049-05	F07	HQP023091	NM_001020658	PUM1
QG049-05	F08	HQP021346	NM_001001890	RUNX1
QG049-05	F09	HQP016575	NM_001017369	SC4MOL
QG049-05	F10	HQP017557	NM_000344	SMN1
QG049-05	F11	HQP001231	NM_007029	STMN2
QG049-05	F12	HQP018392	NM_001005781	SUMO1
QG049-05	G01	HQP006162	NM_015355	SUZ12
QG049-05	G02	HQP018017	NM_198253	TERT
QG049-05	G03	HQP020063	NM_001039590	USP9X
QG049-05	G04	HQP011253	NM_005170	ASCL2
QG049-05	G05	HQP023466	NM_004360	CDH1
QG049-05	G06	HQP012238	NM_021044	DHH
QG049-05	G07	HQP001313	NM_007067	MYST2
QG049-05	G08	HQP022418	NM_001768	CD8A
QG049-05	G09	HQP012245	NM_016948	PARD6A
QG049-05	G10	HQP017280	NM_001199	BMP1
QG049-05	G11	HQP011689	NM_001076682	NCAM1
QG049-05	G12	HQP021189	NM_003655	CBX4
QG049-05	H01	HGDC		
QG049-05	H02	HGDC		
QG049-05	H03	HQP006940	NM_002046	GAPDH

QG049-05	H04	HQP016381	NM_001101	ACTB
QG049-05	H05	HQP015171	NM_004048	B2M
QG049-05	H06	HQP006171	NM_012423	RPL13A
QG049-05	H07	HQP009026	NM_000194	HPRT1
QG049-05	H08	HQP054253	NR_003286	RN18S1
QG049-05	H09	RT		
QG049-05	H10	RT		
QG049-05	H11	PCR		
QG049-05	H12	PCR		