

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG096-01	A01	HQP005152	NM_173555	TYSND1
QG096-01	A02	HQP011260	NM_001002036	ASTL
QG096-01	A03	HQP011040	NM_001001711	DDI1
QG096-01	A04	HQP009339	NM_183378	OVCH1
QG096-01	A05	HQP009338	NM_198185	OVCH2
QG096-01	A06	HQP009253	NM_183062	MPN2
QG096-01	A07	HQP006680	NM_152786	C9orf43
QG096-01	A08	HQP006518	NM_199227	MAP1D
QG096-01	A09	HQP006473	NM_173567	ABHD7
QG096-01	A10	HQP005110	NM_182488	USP12
QG096-01	A11	HQP004954	NM_152699	SENP5
QG096-01	A12	HQP004885	NM_145004	ADAM32
QG096-01	B01	HQP004884	NM_153692	HTRA4
QG096-01	B02	HQP004881	NM_198464	UNQ9391
QG096-01	B03	HQP004619	NM_144981	IMMP1L
QG096-01	B04	HQP004286	NM_133638	ADAMTS19
QG096-01	B05	HQP004273	NM_139057	ADAMTS17
QG096-01	B06	HQP004272	NM_139056	ADAMTS16
QG096-01	B07	HQP004271	NM_139055	ADAMTS15
QG096-01	B08	HQP004199	NM_153362	PRSS35
QG096-01	B09	HQP004021	NM_138328	RHBDL3
QG096-01	B10	HQP003993	NM_130901	OTUD7A
QG096-01	B11	HQP003916	NM_152586	USP54
QG096-01	B12	HQP003903	NM_201286	USP51
QG096-01	C01	HQP003631	NM_152792	SASP
QG096-01	C02	HQP003467	NM_152491	FLJ32569
QG096-01	C03	HQP003313	NM_173502	PRSS36
QG096-01	C04	HQP003141	NM_001025096	PSMA8
QG096-01	C05	HQP002849	NM_001001317	TRY1
QG096-01	C06	HQP002810	NM_001010853	ACY1L2
QG096-01	C07	HQP002808	NM_133493	CD109
QG096-01	C08	HQP002214	NM_152336	AGBL1
QG096-01	C09	HQP002146	NM_139015	UNQ1887
QG096-01	C10	HQP002050	NM_198148	CPXM2
QG096-01	C11	HQP002019	NM_147191	MMP21
QG096-01	C12	HQP001781	NM_145243	OMA1
QG096-01	D01	HQP001440	NM_033341	BIRC8
QG096-01	D02	HQP022682	NM_053044	HTRA3
QG096-01	D03	HQP022353	NM_144575	CAPN13
QG096-01	D04	HQP022234	NM_152271	LONRF1
QG096-01	D05	HQP022169	NM_033276	XRCC6BP1

QG096-01	D06	HQP021775	NM_033029	LMLN
QG096-01	D07	HQP021086	NM_032885	ATG4D
QG096-01	D08	HQP021058	NM_032859	ABHD13
QG096-01	D09	HQP021051	NM_032852	ATG4C
QG096-01	D10	HQP021002	NM_032802	SPPL2A
QG096-01	D11	HQP020987	NM_032785	AGBL4
QG096-01	D12	HQP020923	NM_032649	CNDP1
QG096-01	E01	HQP020865	NM_032582	USP32
QG096-01	E02	HQP020573	NM_032276	RHBDD1
QG096-01	E03	HQP020542	NM_001032730	USP48
QG096-01	E04	HQP020501	NM_032172	USP42
QG096-01	E05	HQP020484	NM_001042403	USP44
QG096-01	E06	HQP020402	NM_032549	IMMP2L
QG096-01	E07	HQP020371	NM_031948	PRSS27
QG096-01	E08	HQP020323	NM_031490	LONP2
QG096-01	E09	HQP020303	NM_031474	NRIP2
QG096-01	E10	HQP019983	NM_030957	ADAMTS10
QG096-01	E11	HQP019981	NM_030955	ADAMTS12
QG096-01	E12	HQP019526	NM_025054	VCPIP1
QG096-01	F01	HQP019346	NM_024817	THSD4
QG096-01	F02	HQP019313	NM_024783	AGBL2
QG096-01	F03	HQP019117	NM_024583	SCRN3
QG096-01	F04	HQP018955	NM_001032278	MMP28
QG096-01	F05	HQP017206	NM_022832	USP46
QG096-01	F06	HQP017018	NM_024164	TPSB2
QG096-01	F07	HQP016839	NM_022357	DPEP3
QG096-01	F08	HQP016833	NM_022353	OSGEPL1
QG096-01	F09	HQP016754	NM_022122	MMP27
QG096-01	F10	HQP016705	NM_022098	XPNPEP3
QG096-01	F11	HQP016665	NM_022060	ABHD4
QG096-01	F12	HQP016571	NM_033440	ELA2A
QG096-01	G01	HQP016354	NM_020318	PAPPA2
QG096-01	G02	HQP016338	NM_021928	SPCS3
QG096-01	G03	HQP016331	NM_001035507	AGBL5
QG096-01	G04	HQP016166	NM_021627	SENP2
QG096-01	G05	HQP016165	NM_021626	SCPEP1
QG096-01	G06	HQP016137	NM_021804	ACE2
QG096-01	G07	HQP015809	NM_020935	USP37
QG096-01	G08	HQP015773	NM_020886	USP28
QG096-01	G09	HQP015762	NM_001004360	DPP10
QG096-01	G10	HQP015746	NM_025090	USP36
QG096-01	G11	HQP015710	NM_020799	STAMBPL1
QG096-01	G12	HQP015643	NM_020726	NLN

QG096-01	H01	HGDC		
QG096-01	H02	HGDC		
QG096-01	H03	HQP006940	NM_002046	GAPDH
QG096-01	H04	HQP016381	NM_001101	ACTB
QG096-01	H05	HQP015171	NM_004048	B2M
QG096-01	H06	HQP006171	NM_012423	RPL13A
QG096-01	H07	HQP009026	NM_000194	HPRT1
QG096-01	H08	HQP054253	NR_003286	RN18S1
QG096-01	H09	RT		
QG096-01	H10	RT		
QG096-01	H11	PCR		
QG096-01	H12	PCR		
QG096-02	A01	HQP015571	NM_020665	TMEM27
QG096-02	A02	HQP015549	NM_001077203	SENP7
QG096-02	A03	HQP015431	NM_018226	RNPEPL1
QG096-02	A04	HQP015385	NM_020361	CPA6
QG096-02	A05	HQP015322	NM_182920	ADAMTS9
QG096-02	A06	HQP015282	NM_020205	OTUD7B
QG096-02	A07	HQP015212	NM_020134	DPYSL5
QG096-02	A08	HQP015151	NM_020645	NRIP3
QG096-02	A09	HQP014659	NM_018235	CNDP2
QG096-02	A10	HQP014544	NM_017807	OSGEP
QG096-02	A11	HQP014516	NM_017714	TASP1
QG096-02	A12	HQP014508	NM_017670	OTUB1
QG096-02	B01	HQP014494	NM_017602	OTUD5
QG096-02	B02	HQP014462	NM_017509	KLK15
QG096-02	B03	HQP014401	NM_018566	YOD1
QG096-02	B04	HQP014354	NM_018394	ABHD10
QG096-02	B05	HQP014328	NM_018359	C4orf20
QG096-02	B06	HQP014204	NM_018218	USP40
QG096-02	B07	HQP013991	NM_017944	USP47
QG096-02	B08	HQP013872	NM_017821	RHBDL2
QG096-02	B09	HQP013781	NM_017712	PGPEP1
QG096-02	B10	HQP013680	NM_017580	ZRANB1
QG096-02	B11	HQP013677	NM_017573	PCSK4
QG096-02	B12	HQP013547	NM_019050	USP53
QG096-02	C01	HQP013009	NM_001040458	ARTS-1
QG096-02	C02	HQP012692	NM_015984	UCHL5
QG096-02	C03	HQP012673	NM_015884	MBTPS2
QG096-02	C04	HQP012584	NM_016546	C1RL
QG096-02	C05	HQP012509	NM_016352	CPA4
QG096-02	C06	HQP012392	NM_016006	ABHD5
QG096-02	C07	HQP012337	NM_015907	LAP3

QG096-02	C08	HQP011329	NM_022046	KLK14
QG096-02	C09	HQP008625	NM_013379	DPP7
QG096-02	C10	HQP008550	NM_014554	SENP1
QG096-02	C11	HQP008516	NM_013396	USP25
QG096-02	C12	HQP008440	NM_013270	TSP50
QG096-02	D01	HQP008424	NM_013258	PYCARD
QG096-02	D02	HQP008345	NM_014058	TMPRSS11E
QG096-02	D03	HQP008333	NM_014041	SPCS1
QG096-02	D04	HQP007710	NM_013247	HTRA2
QG096-02	D05	HQP007649	NM_014479	ADAMDEC1
QG096-02	D06	HQP007631	NM_014464	TINAG
QG096-02	D07	HQP007100	NM_015670	SENP3
QG096-02	D08	HQP007024	NM_015596	KLK13
QG096-02	D09	HQP007008	NM_015571	SENP6
QG096-02	D10	HQP006828	NM_001001991	DKFZP586H2123
QG096-02	D11	HQP006799	NM_018561	USP49
QG096-02	D12	HQP006760	NM_012105	BACE2
QG096-02	E01	HQP006748	NM_001077491	KLK5
QG096-02	E02	HQP006726	NM_012413	QPCT
QG096-02	E03	HQP006312	NM_016192	TMEFF2
QG096-02	E04	HQP006277	NM_012104	BACE1
QG096-02	E05	HQP006233	NM_012114	CASP14
QG096-02	E06	HQP006204	NM_012100	DNPEP
QG096-02	E07	HQP006137	NM_014300	SEC11A
QG096-02	E08	HQP006132	NM_014296	CAPN7
QG096-02	E09	HQP005957	NM_015239	AGTPBP1
QG096-02	E10	HQP005864	NM_015160	PMPCA
QG096-02	E11	HQP005850	NM_013325	ATG4B
QG096-02	E12	HQP005828	NM_015143	METAP1
QG096-02	F01	HQP005745	NM_001040152	PEG10
QG096-02	F02	HQP005686	NM_015017	USP33
QG096-02	F03	HQP001582	NM_007272	CTRC
QG096-02	F04	HQP001496	NM_017414	USP18
QG096-02	F05	HQP001400	NM_007196	KLK8
QG096-02	F06	HQP001391	NM_007192	SUPT16H
QG096-02	F07	HQP001357	NM_014272	ADAMTS7
QG096-02	F08	HQP001299	NM_021251	CAPN10
QG096-02	F09	HQP001298	NM_007058	CAPN11
QG096-02	F10	HQP001260	NM_007173	PRSS23
QG096-02	F11	HQP001258	NM_007038	ADAMTS5
QG096-02	F12	HQP001257	NM_007037	ADAMTS8
QG096-02	G01	HQP001248	NM_014269	ADAM29
QG096-02	G02	HQP001247	NM_021794	ADAM30

QG096-02	G03	HQP001123	NM_006838	METAP2
QG096-02	G04	HQP001122	NM_006837	COPS5
QG096-02	G05	HQP001071	NM_006796	AFG3L2
QG096-02	G06	HQP001019	NM_006690	MMP24
QG096-02	G07	HQP001017	NM_006785	MALT1
QG096-02	G08	HQP000993	NM_006677	USP19
QG096-02	G09	HQP000992	NM_001008563	USP20
QG096-02	G10	HQP000984	NM_014265	ADAM28
QG096-02	G11	HQP000967	NM_006660	CLPX
QG096-02	G12	HQP000888	NM_006615	CAPN9
QG096-02	H01	HGDC		
QG096-02	H02	HGDC		
QG096-02	H03	HQP006940	NM_002046	GAPDH
QG096-02	H04	HQP016381	NM_001101	ACTB
QG096-02	H05	HQP015171	NM_004048	B2M
QG096-02	H06	HQP006171	NM_012423	RPL13A
QG096-02	H07	HQP009026	NM_000194	HPRT1
QG096-02	H08	HQP054253	NR_003286	RN18S1
QG096-02	H09	RT		
QG096-02	H10	RT		
QG096-02	H11	PCR		
QG096-02	H12	PCR		
QG096-03	A01	HQP000882	NM_006610	MASP2
QG096-03	A02	HQP000864	NM_014263	YME1L1
QG096-03	A03	HQP000846	NM_006590	USP39
QG096-03	A04	HQP000841	NM_006587	CORIN
QG096-03	A05	HQP000747	NM_006463	STAMPB
QG096-03	A06	HQP000723	NM_001001992	USP16
QG096-03	A07	HQP000640	NM_014889	PITRM1
QG096-03	A08	HQP000481	NM_006103	WFDC2
QG096-03	A09	HQP000480	NM_016134	PGCP
QG096-03	A10	HQP000341	NM_005865	PRSS16
QG096-03	A11	HQP000265	NM_005805	PSMD14
QG096-03	A12	HQP000168	NM_005747	ELA3A
QG096-03	B01	HQP000045	NM_005482	PIGK
QG096-03	B02	HQP000007	NM_005468	NAALADL1
QG096-03	B03	HQP000006	NM_005467	NAALAD2
QG096-03	B04	HQP023448	NM_001032279	RCE1
QG096-03	B05	HQP023411	NM_006537	USP3
QG096-03	B06	HQP023408	NM_006313	USP15
QG096-03	B07	HQP023370	NM_014871	USP52
QG096-03	B08	HQP023202	NM_014752	SPCS2
QG096-03	B09	HQP023140	NM_014709	USP34

QG096-03	B10	HQP023116	NM_014694	ADAMTSL2
QG096-03	B11	HQP023114	NM_001037324	ECE2
QG096-03	B12	HQP023095	NM_012291	ESPL1
QG096-03	C01	HQP022998	NM_004917	KLK4
QG096-03	C02	HQP022936	NM_001042385	PREPL
QG096-03	C03	HQP022857	NM_006310	NPEPPS
QG096-03	C04	HQP022847	NM_006988	ADAMTS1
QG096-03	C05	HQP022844	NM_014244	ADAMTS2
QG096-03	C06	HQP022843	NM_014243	ADAMTS3
QG096-03	C07	HQP022842	NM_005099	ADAMTS4
QG096-03	C08	HQP022806	NM_004851	NAPSA
QG096-03	C09	HQP022602	NM_004793	LONP1
QG096-03	C10	HQP022517	NM_004771	MMP20
QG096-03	C11	HQP022211	NM_004716	PCSK7
QG096-03	C12	HQP022076	NM_005154	USP8
QG096-03	D01	HQP022072	NM_005153	USP10
QG096-03	D02	HQP022065	NM_004505	USP6
QG096-03	D03	HQP022063	NM_001037334	USP14
QG096-03	D04	HQP021764	NM_003940	USP13
QG096-03	D05	HQP021700	NM_006025	P11
QG096-03	D06	HQP021604	NM_003879	CFLAR
QG096-03	D07	HQP021603	NM_003878	GGH
QG096-03	D08	HQP021517	NM_003817	ADAM7
QG096-03	D09	HQP021509	NM_003815	ADAM15
QG096-03	D10	HQP021507	NM_014237	ADAM18
QG096-03	D11	HQP021504	NM_003812	ADAM23
QG096-03	D12	HQP021493	NM_003805	CRADD
QG096-03	E01	HQP021483	NM_023038	ADAM19
QG096-03	E02	HQP021475	NM_003793	CTSF
QG096-03	E03	HQP021471	NM_003791	MBTPS1
QG096-03	E04	HQP021042	NM_003619	PRSS12
QG096-03	E05	HQP020117	NM_004656	BAP1
QG096-03	E06	HQP020088	NM_004654	USP9Y
QG096-03	E07	HQP020062	NM_004651	USP11
QG096-03	E08	HQP019770	NM_003481	USP5
QG096-03	E09	HQP018818	NM_003470	USP7
QG096-03	E10	HQP018560	NM_003399	XPNPEP2
QG096-03	E11	HQP018559	NM_020383	XPNPEP1
QG096-03	E12	HQP018447	NM_001017415	USP1
QG096-03	F01	HQP018437	NM_003366	UQCRC2
QG096-03	F02	HQP018436	NM_003365	UQCRC1
QG096-03	F03	HQP018429	NM_003363	USP4
QG096-03	F04	HQP018398	NM_006002	UCHL3

QG096-03	F05	HQP018397	NM_004181	UCHL1
QG096-03	F06	HQP018221	NM_003294	TPSAB1
QG096-03	F07	HQP018219	NM_003291	TPP2
QG096-03	F08	HQP018145	NM_006290	TNFAIP3
QG096-03	F09	HQP018136	NM_005656	TMPRSS2
QG096-03	F10	HQP018110	NM_012465	TLL2
QG096-03	F11	HQP018109	NM_012464	TLL1
QG096-03	F12	HQP018073	NM_003249	THOP1
QG096-03	G01	HQP018041	NM_003234	TFRC
QG096-03	G02	HQP017875	NM_003184	TAF2
QG096-03	G03	HQP017866	NM_003183	ADAM17
QG096-03	G04	HQP017759	NM_021978	ST14
QG096-03	G05	HQP017660	NM_003119	SPG7
QG096-03	G06	HQP017620	NM_005632	SOLH
QG096-03	G07	HQP017098	NM_000193	SHH
QG096-03	G08	HQP016333	NM_020216	RNPEP
QG096-03	G09	HQP016215	NM_000537	REN
QG096-03	G10	HQP016039	NM_002864	PZP
QG096-03	G11	HQP015430	NM_002811	PSMD7
QG096-03	G12	HQP015323	NM_002801	PSMB10
QG096-03	H01	HGDC		
QG096-03	H02	HGDC		
QG096-03	H03	HQP006940	NM_002046	GAPDH
QG096-03	H04	HQP016381	NM_001101	ACTB
QG096-03	H05	HQP015171	NM_004048	B2M
QG096-03	H06	HQP006171	NM_012423	RPL13A
QG096-03	H07	HQP009026	NM_000194	HPRT1
QG096-03	H08	HQP054253	NR_003286	RN18S1
QG096-03	H09	RT		
QG096-03	H10	RT		
QG096-03	H11	PCR		
QG096-03	H12	PCR		
QG096-04	A01	HQP015311	NM_002800	PSMB9
QG096-04	A02	HQP015291	NM_004159	PSMB8
QG096-04	A03	HQP015273	NM_002798	PSMB6
QG096-04	A04	HQP015262	NM_002797	PSMB5
QG096-04	A05	HQP015253	NM_002796	PSMB4
QG096-04	A06	HQP015241	NM_002795	PSMB3
QG096-04	A07	HQP015228	NM_002794	PSMB2
QG096-04	A08	HQP015219	NM_002793	PSMB1
QG096-04	A09	HQP015194	NM_002791	PSMA6
QG096-04	A10	HQP015192	NM_002789	PSMA4
QG096-04	A11	HQP015177	NM_002786	PSMA1

QG096-04	A12	HQP015123	NM_000021	PSEN1
QG096-04	B01	HQP015105	NM_002775	HTRA1
QG096-04	B02	HQP015100	NM_002773	PRSS8
QG096-04	B03	HQP015097	NM_002772	PRSS7
QG096-04	B04	HQP015096	NM_005046	KLK7
QG096-04	B05	HQP015084	NM_002770	PRSS2
QG096-04	B06	HQP015083	NM_002769	PRSS1
QG096-04	B07	HQP015082	NM_001008530	LGMN
QG096-04	B08	HQP015041	NM_000312	PROC
QG096-04	B09	HQP014431	NM_002726	PREP
QG096-04	B10	HQP013257	NM_000301	PLG
QG096-04	B11	HQP013204	NM_002658	PLAU
QG096-04	B12	HQP013201	NM_000930	PLAT
QG096-04	C01	HQP013125	NM_002639	SERPINB5
QG096-04	C02	HQP013122	NM_000295	SERPINA1
QG096-04	C03	HQP013108	NM_000444	PHEX
QG096-04	C04	HQP013087	NM_002630	PGC
QG096-04	C05	HQP013045	NM_000285	PEPD
QG096-04	C06	HQP012576	NM_002594	PCSK2
QG096-04	C07	HQP012566	NM_006200	PCSK5
QG096-04	C08	HQP012539	NM_000439	PCSK1
QG096-04	C09	HQP012332	NM_000624	SERPINA5
QG096-04	C10	HQP012189	NM_002581	PAPPA
QG096-04	C11	HQP012155	NM_002575	SERPINB2
QG096-04	C12	HQP012154	NM_000602	SERPINE1
QG096-04	D01	HQP012121	NM_002569	FURIN
QG096-04	D02	HQP011931	NM_006178	NSF
QG096-04	D03	HQP011918	NM_002525	NRD1
QG096-04	D04	HQP011272	NM_016155	MMP17
QG096-04	D05	HQP011270	NM_005941	MMP16
QG096-04	D06	HQP011269	NM_002428	MMP15
QG096-04	D07	HQP011268	NM_004995	MMP14
QG096-04	D08	HQP011267	NM_002427	MMP13
QG096-04	D09	HQP011266	NM_002426	MMP12
QG096-04	D10	HQP011265	NM_005940	MMP11
QG096-04	D11	HQP011264	NM_002425	MMP10
QG096-04	D12	HQP011263	NM_004994	MMP9
QG096-04	E01	HQP011262	NM_002424	MMP8
QG096-04	E02	HQP011258	NM_002423	MMP7
QG096-04	E03	HQP011257	NM_002422	MMP3
QG096-04	E04	HQP011255	NM_002421	MMP1
QG096-04	E05	HQP011254	NM_000902	MME
QG096-04	E06	HQP011222	NM_005932	MIPEP

QG096-04	E07	HQP011177	NM_005588	MEP1A
QG096-04	E08	HQP011129	NM_002390	ADAM11
QG096-04	E09	HQP010919	NM_002343	LTF
QG096-04	E10	HQP010906	NM_000895	LTA4H
QG096-04	E11	HQP010792	NM_005575	LNPEP
QG096-04	E12	HQP010104	NM_000892	KLKB1
QG096-04	F01	HQP010100	NM_002257	KLK1
QG096-04	F02	HQP010071	NM_000420	KEL
QG096-04	F03	HQP009632	NM_002181	IHH
QG096-04	F04	HQP009019	NM_005143	HP
QG096-04	F05	HQP008800	NM_000601	HGF
QG096-04	F06	HQP008711	NM_004132	HABP2
QG096-04	F07	HQP008691	NM_002104	GZMK
QG096-04	F08	HQP008690	NM_004131	GZMB
QG096-04	F09	HQP008689	NM_006144	GZMA
QG096-04	F10	HQP008675	NM_033423	GZMH
QG096-04	F11	HQP008463	NM_005313	PDIA3
QG096-04	F12	HQP006129	NM_001014986	FOLH1
QG096-04	G01	HQP005068	NM_000505	F12
QG096-04	G02	HQP005066	NM_000128	F11
QG096-04	G03	HQP005064	NM_000504	F10
QG096-04	G04	HQP005052	NM_000506	F2
QG096-04	G05	HQP004948	NM_000120	EPHX1
QG096-04	G06	HQP004867	NM_001977	ENPEP
QG096-04	G07	HQP004860	NM_001428	ENO1
QG096-04	G08	HQP004689	NM_001972	ELA2
QG096-04	G09	HQP004688	NM_001971	ELA1
QG096-04	G10	HQP004537	NM_001397	ECE1
QG096-04	G11	HQP004435	NM_001039350	DPP6
QG096-04	G12	HQP004434	NM_001935	DPP4
QG096-04	H01	HGDC		
QG096-04	H02	HGDC		
QG096-04	H03	HQP006940	NM_002046	GAPDH
QG096-04	H04	HQP016381	NM_001101	ACTB
QG096-04	H05	HQP015171	NM_004048	B2M
QG096-04	H06	HQP006171	NM_012423	RPL13A
QG096-04	H07	HQP009026	NM_000194	HPRT1
QG096-04	H08	HQP054253	NR_003286	RN18S1
QG096-04	H09	RT		
QG096-04	H10	RT		
QG096-04	H11	PCR		
QG096-04	H12	PCR		
QG096-05	A01	HQP004198	NM_001928	CFD

QG096-05	A02	HQP004081	NM_000789	ACE
QG096-05	A03	HQP003761	NM_001042355	CYLD
QG096-05	A04	HQP003668	NM_004079	CTSS
QG096-05	A05	HQP003658	NM_001334	CTSO
QG096-05	A06	HQP003630	NM_001912	CTSL1
QG096-05	A07	HQP003626	NM_000396	CTSK
QG096-05	A08	HQP003620	NM_004390	CTSH
QG096-05	A09	HQP003612	NM_001911	CTSG
QG096-05	A10	HQP003605	NM_001910	CTSE
QG096-05	A11	HQP003599	NM_001909	CTSD
QG096-05	A12	HQP003596	NM_001908	CTSB
QG096-05	B01	HQP003577	NM_001906	CTRB1
QG096-05	B02	HQP002857	NM_001308	CPN1
QG096-05	B03	HQP002855	NM_001005502	CPM
QG096-05	B04	HQP002847	NM_001873	CPE
QG096-05	B05	HQP002842	NM_001304	CPD
QG096-05	B06	HQP002835	NM_001871	CPB1
QG096-05	B07	HQP002832	NM_001870	CPA3
QG096-05	B08	HQP002821	NM_001868	CPA1
QG096-05	B09	HQP002142	NM_001836	CMA1
QG096-05	B10	HQP002078	NM_000391	TPP1
QG096-05	B11	HQP000894	NM_001814	CTSC
QG096-05	B12	HQP020709	NM_001230	CASP10
QG096-05	C01	HQP020648	NM_001229	CASP9
QG096-05	C02	HQP018966	NM_001080124	CASP8
QG096-05	C03	HQP020480	NM_001227	CASP7
QG096-05	C04	HQP020427	NM_001226	CASP6
QG096-05	C05	HQP020384	NM_004347	CASP5
QG096-05	C06	HQP020332	NM_001225	CASP4
QG096-05	C07	HQP020297	NM_004346	CASP3
QG096-05	C08	HQP020249	NM_032982	CASP2
QG096-05	C09	HQP020207	NM_001223	CASP1
QG096-05	C10	HQP020086	NM_014289	CAPN6
QG096-05	C11	HQP020071	NM_000070	CAPN3
QG096-05	C12	HQP020070	NM_001748	CAPN2
QG096-05	D01	HQP020065	NM_005186	CAPN1
QG096-05	D02	HQP018286	NM_004055	CAPN5
QG096-05	D03	HQP018238	NM_000064	C3
QG096-05	D04	HQP018226	NM_000063	C2
QG096-05	D05	HQP018206	NM_001734	C1S
QG096-05	D06	HQP018179	NM_001733	C1R
QG096-05	D07	HQP016897	NM_000386	BLMH
QG096-05	D08	HQP016565	NM_001710	CFB

QG096-05	D09	HQP009633	NM_001030047	KLK3
QG096-05	D10	HQP009091	NM_001167	BIRC4
QG096-05	D11	HQP009072	NM_001166	BIRC2
QG096-05	D12	HQP009056	NM_001640	APEH
QG096-05	E01	HQP008414	NM_001150	ANPEP
QG096-05	E02	HQP004378	NM_000027	AGA
QG096-05	E03	HQP004154	NM_001129	AEBP1
QG096-05	E04	HQP000372	NM_001110	ADAM10
QG096-05	E05	HQP022963	NM_000666	ACY1
QG096-05	E06	HQP012072	NM_001097	ACR
QG096-05	E07	HQP015156	NM_001700	AZU1
QG096-05	E08	HQP002566	NM_000094	COL7A1
QG096-05	E09	HQP005059	NM_000131	F7
QG096-05	E10	HQP005101	NM_004460	FAP
QG096-05	E11	HQP008692	NM_005317	GZMM
QG096-05	E12	HQP010101	NM_001002231	KLK2
QG096-05	F01	HQP015091	NM_001031849	MASP1
QG096-05	F02	HQP016196	NM_002900	RBP3
QG096-05	F03	HQP021904	NM_003961	RHBDL1
QG096-05	F04	HQP022693	NM_004262	TMPRSS11D
QG096-05	F05	HQP001148	NM_006853	KLK11
QG096-05	F06	HQP001358	NM_197941	ADAMTS6
QG096-05	F07	HQP007550	NM_015692	CPAMD8
QG096-05	F08	HQP011330	NM_019598	KLK12
QG096-05	F09	HQP019131	NM_001005498	RHBDF2
QG096-05	F10	HQP019504	NM_025003	ADAMTS20
QG096-05	F11	HQP019815	NM_030770	TMPRSS5
QG096-05	F12	HQP020337	NM_031907	USP26
QG096-05	G01	HQP021825	NM_173462	PAPLN
QG096-05	G02	HQP002727	NM_182502	TMPRSS11B
QG096-05	G03	HQP003053	NM_080722	ADAMTS14
QG096-05	G04	HQP003800	NM_133463	AMZ1
QG096-05	G05	HQP003867	NM_144966	FREM1
QG096-05	G06	HQP004119	NM_153609	TMPRSS6
QG096-05	G07	HQP006606	NM_174936	PCSK9
QG096-05	G08	HQP008030	NM_012315	KLK9
QG096-05	G09	HQP009051	NM_001042518	SERPINA9
QG096-05	G10	HQP009274	NM_182702	TESSP2
QG096-05	G11	HQP009439	NM_001042575	TMPRSS7
QG096-05	G12	HQP009712	NM_182973	TMPRSS9
QG096-05	H01	HGDC		
QG096-05	H02	HGDC		
QG096-05	H03	HQP006940	NM_002046	GAPDH

QG096-05	H04	HQP016381	NM_001101	ACTB
QG096-05	H05	HQP015171	NM_004048	B2M
QG096-05	H06	HQP006171	NM_012423	RPL13A
QG096-05	H07	HQP009026	NM_000194	HPRT1
QG096-05	H08	HQP054253	NR_003286	RN18S1
QG096-05	H09	RT		
QG096-05	H10	RT		
QG096-05	H11	PCR		
QG096-05	H12	PCR		
QG096-06	A01	HQP000246	NM_001109	ADAM8
QG096-06	A02	HQP004322	NM_001031805	AFG3L1
QG096-06	A03	HQP017280	NM_001199	BMP1
QG096-06	A04	HQP020082	NM_001003962	CAPNS1
QG096-06	A05	HQP002824	NM_001869	CPA2
QG096-06	A06	HQP002836	NM_001872	CPB2
QG096-06	A07	HQP002860	NM_001080513	CPN2
QG096-06	A08	HQP003587	NM_001907	CTRL
QG096-06	A09	HQP003635	NM_001333	CTSL2
QG096-06	A10	HQP003677	NM_001335	CTSW
QG096-06	A11	HQP003680	NM_001336	CTSZ
QG096-06	A12	HQP004430	NM_004413	DPEP1
QG096-06	B01	HQP020429	NM_001077263	TMPRSS13
QG096-06	B02	HQP007378	NM_001032364	GGT1
QG096-06	B03	HQP008811	NM_001528	HGFAC
QG096-06	B04	HQP009023	NM_002151	HPN
QG096-06	B05	HQP009349	NM_004969	IDE
QG096-06	B06	HQP009376	NM_000204	CFI
QG096-06	B07	HQP011178	NM_005925	MEP1B
QG096-06	B08	HQP054072	NM_004530	MMP2
QG096-06	B09	HQP020839	NM_032557	USP38
QG096-06	B10	HQP011524	NM_020998	MST1
QG096-06	B11	HQP012112	NM_006191	PA2G4
QG096-06	B12	HQP012122	NM_002570	PCSK6
QG096-06	C01	HQP012507	NM_001083314	CHMP1A
QG096-06	C02	HQP013084	NM_014224	PGA5
QG096-06	C03	HQP013686	NM_000308	CTSA
QG096-06	C04	HQP014413	NM_005040	PRCP
QG096-06	C05	HQP021022	NM_032823	C9orf3
QG096-06	C06	HQP015101	NM_001012964	KLK6
QG096-06	C07	HQP015107	NM_002777	PRTN3
QG096-06	C08	HQP015130	NM_000447	PSEN2
QG096-06	C09	HQP015182	NM_002787	PSMA2
QG096-06	C10	HQP015185	NM_002788	PSMA3

QG096-06	C11	HQP015193	NM_002790	PSMA5
QG096-06	C12	HQP015201	NM_002792	PSMA7
QG096-06	D01	HQP015285	NM_002799	PSMB7
QG096-06	D02	HQP017668	NM_001032367	SPINT1
QG096-06	D03	HQP019702	NM_003474	ADAM12
QG096-06	D04	HQP020030	NM_006012	CLPP
QG096-06	D05	HQP020063	NM_001039590	USP9X
QG096-06	D06	HQP021134	NM_006983	MMP23B
QG096-06	D07	HQP021182	NM_001014447	CPZ
QG096-06	D08	HQP021505	NM_003813	ADAM21
QG096-06	D09	HQP021117	NM_001080481	USP45
QG096-06	D10	HQP021515	NM_001005845	ADAM9
QG096-06	D11	HQP021634	NM_003891	PROZ
QG096-06	D12	HQP021964	NM_138355	SCRN2
QG096-06	E01	HQP022744	NM_004826	ECEL1
QG096-06	E02	HQP000094	NM_005700	DPP3
QG096-06	E03	HQP000259	NM_005800	USPL1
QG096-06	E04	HQP022016	NM_033280	SEC11C
QG096-06	E05	HQP006097	NM_012217	TPSD1
QG096-06	E06	HQP022077	NM_139159	DPP9
QG096-06	E07	HQP007029	NM_001042472	ABHD12
QG096-06	E08	HQP007428	NM_001014443	USP21
QG096-06	E09	HQP012238	NM_021044	DHH
QG096-06	E10	HQP012486	NM_016302	CRBN
QG096-06	E11	HQP012637	NM_001033569	AMZ2
QG096-06	E12	HQP012918	NM_016023	OTUD6B
QG096-06	F01	HQP022481	NM_052866	ADAMTSL1
QG096-06	F02	HQP013523	NM_019029	CPVL
QG096-06	F03	HQP013525	NM_019032	ADAMTSL4
QG096-06	F04	HQP013810	NM_017743	DPP8
QG096-06	F05	HQP014415	NM_001037639	PARL
QG096-06	F06	HQP015103	NM_021801	MMP26
QG096-06	F07	HQP015128	NM_019894	TMPRSS4
QG096-06	F08	HQP003527	NM_145652	WFDC5
QG096-06	F09	HQP015789	NM_020903	USP29
QG096-06	F10	HQP016751	NM_022119	PRSS22
QG096-06	F11	HQP016802	NM_022164	TINAGL1
QG096-06	F12	HQP016828	NM_022350	LRAP
QG096-06	G01	HQP016837	NM_022355	DPEP2
QG096-06	G02	HQP016887	NM_022450	RHBDF1
QG096-06	G03	HQP016947	NM_004142	MMP25
QG096-06	G04	HQP017095	NM_024022	TMPRSS3
QG096-06	G05	HQP018831	NM_023112	OTUB2

QG096-06	G06	HQP018987	NM_001018055	BRCC3
QG096-06	G07	HQP019193	NM_024663	NPEPL1
QG096-06	G08	HQP019307	NM_001031855	LONRF3
QG096-06	G09	HQP004025	NM_175882	IMP5
QG096-06	G10	HQP019876	NM_030789	HM13
QG096-06	G11	HQP019939	NM_030929	KAZALD1
QG096-06	G12	HQP020190	NM_031301	APH1B
QG096-06	H01	HGDC		
QG096-06	H02	HGDC		
QG096-06	H03	HQP006940	NM_002046	GAPDH
QG096-06	H04	HQP016381	NM_001101	ACTB
QG096-06	H05	HQP015171	NM_004048	B2M
QG096-06	H06	HQP006171	NM_012423	RPL13A
QG096-06	H07	HQP009026	NM_000194	HPRT1
QG096-06	H08	HQP054253	NR_003286	RN18S1
QG096-06	H09	RT		
QG096-06	H10	RT		
QG096-06	H11	PCR		
QG096-06	H12	PCR		