

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG013-01	A01	HQP017661	NM_199367	SPG7
QG013-01	A02	HQP010506	NM_198129	LAMA3
QG013-01	A03	HQP006023	NM_054034	FN1
QG013-01	A04	HQP022974	NM_001001390	CD44
QG013-01	A05	HQP009026	NM_000194	HPRT1
QG013-01	A06	HQP018502	NM_000638	VTN
QG013-01	A07	HQP053975	NM_001078	VCAM1
QG013-01	A08	HQP009140	NM_002160	TNC
QG013-01	A09	HQP018140	NM_003278	CLEC3B
QG013-01	A10	HQP018094	NM_000362	TIMP3
QG013-01	A11	HQP018093	NM_003255	TIMP2
QG013-01	A12	HQP018092	NM_003254	TIMP1
QG013-01	B01	HQP018070	NM_007112	THBS3
QG013-01	B02	HQP018069	NM_003247	THBS2
QG013-01	B03	HQP018068	NM_003246	THBS1
QG013-01	B04	HQP018050	NM_000358	TGFBI
QG013-01	B05	HQP017673	NM_000582	SPP1
QG013-01	B06	HQP017660	NM_003119	SPG7
QG013-01	B07	HQP017653	NM_003118	SPARC
QG013-01	B08	HQP021702	NM_003919	SGCE
QG013-01	B09	HQP016746	NM_003005	SELP
QG013-01	B10	HQP016745	NM_000655	SELL
QG013-01	B11	HQP016744	NM_000450	SELE
QG013-01	B12	HQP013015	NM_000442	PECAM1
QG013-01	C01	HQP011263	NM_004994	MMP9
QG013-01	C02	HQP011262	NM_002424	MMP8
QG013-01	C03	HQP011258	NM_002423	MMP7
QG013-01	C04	HQP011257	NM_002422	MMP3
QG013-01	C05	HQP011270	NM_005941	MMP16
QG013-01	C06	HQP011269	NM_002428	MMP15
QG013-01	C07	HQP011268	NM_004995	MMP14
QG013-01	C08	HQP011267	NM_002427	MMP13
QG013-01	C09	HQP011266	NM_002426	MMP12
QG013-01	C10	HQP011265	NM_005940	MMP11
QG013-01	C11	HQP011264	NM_002425	MMP10
QG013-01	C12	HQP011255	NM_002421	MMP1
QG013-01	D01	HQP010527	NM_002291	LAMB1
QG013-01	D02	HQP010505	NM_000227	LAMA3
QG013-01	D03	HQP010497	NM_000426	LAMA2
QG013-01	D04	HQP008000	NM_005559	LAMA1
QG013-01	D05	HQP009858	NM_000216	KAL1

QG013-01	D06	HQP009825	NM_002213	ITGB5
QG013-01	D07	HQP009819	NM_000213	ITGB4
QG013-01	D08	HQP009818	NM_000212	ITGB3
QG013-01	D09	HQP009815	NM_000211	ITGB2
QG013-01	D10	HQP009810	NM_002211	ITGB1
QG013-01	D11	HQP009808	NM_002210	ITGAV
QG013-01	D12	HQP009807	NM_000632	ITGAM
QG013-01	E01	HQP009806	NM_002209	ITGAL
QG013-01	E02	HQP021139	NM_003638	ITGA8
QG013-01	E03	HQP009800	NM_002206	ITGA7
QG013-01	E04	HQP009774	NM_000210	ITGA6
QG013-01	E05	HQP009799	NM_002205	ITGA5
QG013-01	E06	HQP009798	NM_000885	ITGA4
QG013-01	E07	HQP009794	NM_002203	ITGA2
QG013-01	E08	HQP009793	NM_181501	ITGA1
QG013-01	E09	HQP009184	NM_000201	ICAM1
QG013-01	E10	HQP006022	NM_002026	FN1
QG013-01	E11	HQP003544	NM_001331	CTNND1
QG013-01	E12	HQP003539	NM_001904	CTNNB1
QG013-01	F01	HQP003515	NM_001903	CTNNA1
QG013-01	F02	HQP003488	NM_001901	CTGF
QG013-01	F03	HQP003299	NM_004385	VCAN
QG013-01	F04	HQP002571	NM_001850	COL8A1
QG013-01	F05	HQP002551	NM_001849	COL6A2
QG013-01	F06	HQP002549	NM_001848	COL6A1
QG013-01	F07	HQP002541	NM_000093	COL5A1
QG013-01	F08	HQP002506	NM_001846	COL4A2
QG013-01	F09	HQP002462	NM_000088	COL1A1
QG013-01	F10	HQP002648	NM_001856	COL16A1
QG013-01	F11	HQP002644	NM_001855	COL15A1
QG013-01	F12	HQP018426	NM_021110	COL14A1
QG013-01	G01	HQP002612	NM_004370	COL12A1
QG013-01	G02	HQP002602	NM_080629	COL11A1
QG013-01	G03	HQP002435	NM_001843	CNTN1
QG013-01	G04	HQP022972	NM_000610	CD44
QG013-01	G05	HQP001257	NM_007037	ADAMTS8
QG013-01	G06	HQP022847	NM_006988	ADAMTS1
QG013-01	G07	HQP002566	NM_000094	COL7A1
QG013-01	G08	HQP002613	NM_080645	COL12A1
QG013-01	G09	HQP023466	NM_004360	CDH1
QG013-01	G10	HQP004541	NM_004425	ECM1
QG013-01	G11	HQP008723	NM_001523	HAS1
QG013-01	G12	HQP009796	NM_002204	ITGA3

QG013-01	H01	HGDC		
QG013-01	H02	HGDC		
QG013-01	H03	HQP006940	NM_002046	GAPDH
QG013-01	H04	HQP016381	NM_001101	ACTB
QG013-01	H05	HQP015171	NM_004048	B2M
QG013-01	H06	HQP006171	NM_012423	RPL13A
QG013-01	H07	HQP009026	NM_000194	HPRT1
QG013-01	H08	HQP054253	NR_003286	RN18S1
QG013-01	H09	RT		
QG013-01	H10	RT		
QG013-01	H11	PCR		
QG013-01	H12	PCR		