

ExProfile™ Human Stem Cell Related Gene qPCR Array

For focused group profiling of human stem cell genes expression

Cat. No. QG049-A (5 x 96-well plate, Format A)

Cat. No. QG049-B (5 x 96-well plate, Format B)

Cat. No. QG049-C (5 x 96-well plate, Format C)

Cat. No. QG049-D (5 x 96-well plate, Format D)

Cat. No. QG049-E (5 x 96-well plate, Format E)

Plates available individually or as a set of 6. Each set contains 420 unique gene primer pairs deposited in one 96-well plate.

Introduction

The ExProfile human stem cell related gene qPCR array profiles the expression of 420 human genes related to osteogenic differentiation. These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including genes involved in identification, growth and differentiation of stem cells. This array allows researchers to study the related genes to gain understanding of their roles in the functioning and characterization of stem cell.

- QG049 plate 01: 84 unique gene PCR primer pairs
- QG049 plate 02: 84 unique gene PCR primer pairs
- QG049 plate 03: 84 unique gene PCR primer pairs
- QG049 plate 04: 84 unique gene PCR primer pairs
- QG049 plate 05: 84 unique gene PCR primer pairs

Shipping and storage condition

Shipped at room temperate

Stable for at least 6 months when stored at -20 °C

Array format

GeneCopoeia provides five qPCR array formats (A, B, C, D, and E) suitable for use with the following real-time cyclers.

Important note: Upon receiving, please check to make sure that the correct array format was ordered to ensure the compatibility with your qPCR instrument.

Plate format	Instrument provider	qPCR instrument model
A (96-well)	Applied Biosystems	5700, 7000, 7300, 7500, 7700, 7900HT (Standard 96-well block), ViiA™7 (Standard 96-well block)
B (96-well)	Applied Biosystems	7500 (Fast block), 7900HT (Fast block), StepOnePlus™, ViiA™7 (Fast block)
C (96-well)	Bio-Rad Laboratories	iCycler iQ®, MyiQ™, iQ™5
D (96-well)	Bio-Rad Laboratories	CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™
E (96-well)	Roche Applied Science	LightCycler® 480 (96-well block)

Quality control

1. Each pair of primers in the ExProfile gene qPCR array has been experimentally validated to yield a single dissociation curve peak and to generate a single amplicon of the correct size for the targeted gene.
2. The positive PCR controls (PCR) have been verified to amplify a single amplicon of the correct size with Ct values around **20±2**.
3. The Spike-in reverse transcription controls (RT) have been verified to amplify a single amplicon of the correct size with Ct values around **20±3**.
4. $R^2 > 0.99$ was observed for high inter/ intra-array reproducibility.

Materials required but not provided

All-in-One™ First-Strand cDNA Synthesis Kit

All-in-One™ qPCR Mix

Total RNA extraction kit (RNAzol® RT RNA extraction reagent is recommended)

DNase/RNase free tips, PCR reaction tubes, 1.5 ml microcentrifuge tubes

5 ml and 10 ml graduated pipettes, beakers, flasks, and cylinders

10 µl to 1,000 µl adjustable single channel micropipettes with disposable tips

5 µl to 20 µl adjustable multichannel micropipette, disposable tips, and reservoir

qPCR instrument, compatible with gene qPCR arrays ordered

Array layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	FGFR1	DVL1	COL9A1	CD44	HPRT1	ADAR	WNT1	SOX1	PPARG	PPARD	OPRS1	NOTCH2
B	MYST1	KRT15	HSPA9	HDAC2	GJB2	GDF2	GCN5L2	FZD1	FRAT1	EP300	DVL1	DTX2
C	DTX1	DLL3	DLL1	CTNNA1	CDH2	CDC42	CDC2	CD8B	CD3D	CCNE1	CCNA2	BTRC
D	BMP3	BMP1	BGLAP	AXIN1	ALPI	ALDH2	ALDH1A1	ADAR	ACTC1	WEE1	VIM	VEGFA
E	VDR	VCAM1	UBE3A	TUBB3	TRIM14	TP73L	TP53	TNFSF10	TNFRSF1A	TNF	TNC	TLR4
F	TIMP1	THY1	THPO	TGFB2	TGFB1	TFRC	TFEB	TFAP2C	TFAP2A	TEK	TEAD2	TDGF1
G	TCL1A	TCF7L1	TCF3	TCEA3	TAL1	TAGLN	TACC3	T	SYP	SULT1A1	STAT5B	STAT5A
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG049 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	STAT3	SST	SREBF1	SPP1	SPN	SPI1	SOX9	SOX6	SOX2	SOX17	SOCS3	SMARCA4
B	SLC2A3	SLC22A18	SLAMF1	SLAIN1	SIPA1	SHMT1	SHH	SHC1	SH2B2	SFRP2	SFRP1	SELE
C	SCD5	SALL4	SALL1	S100B	S100A6	S100A4	RUNX2	RPS8KA5	ROCK1	RNF2	RB1	RARG
D	RAC1	RAB3B	PUM2	PTPRF	PTPRC	PTPN11	PTEN	PTCH1	PSEN1	PROM2	PROM1	PPIC
E	PPARG	POU5F1	PODXL	PLAT	PIWIL2	PHF21A	PECAM1	PDGFB	PCGF2	PAX6	PARP4	PAFAH1B1
F	P2RX7	OTX2	NUMB	NR4A2	NOTCH1	NODAL	NOD2	NKD2	NFKB1	NES	NANOG	MYF5
G	MYC	MTRR	MTHFR	MSH2	MS4A1	MPL	MMP9	MME	MITF	MEST	MEN1	MELK
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG049 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	MCFD2	MBL2	MAT2B	MAT2A	MAPK14	MAP3K1	MAP2K1	MAML3	LY6E	LXN	LRIG1	LIFR
B	LIF	LHX2	LEP	LEFTY2	LAMA1	KRT19	KRT17	KITLG	KIT	KDR	JAM2	JAK2
C	JAG2	JAG1	ITGB2	ITGB1	ITGAM	ITGA6	ITGA2B	ITGA2	ITGA1	ISL1	INHBB	IL7
D	IL6ST	IL6	IL5	IL4	IL2	IL15	IL12B	IL12A	IL10	IGF2R	IGF2	IGF1
E	IFNG	ICAM1	HSPG2	HSPB1	HSPA1A	HSP90B1	HOXB4	HOXA9	HOXA10	HMGA1	HIF1A	HGF
F	HES2	HES1	HDAC9	HDAC1	GSK3B	GNMT	GNAT1	GLI1	GJB1	GJA1	GH1	GDNF
G	GDF11	GCM1	GATA6	GATA4	GATA3	GATA2	GABRB3	GABPA	G6PD	FZD7	FUT4	FRAP1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG049 plate 03

	1	2	3	4	5	6	7	8	9	10	11	12
A	FPR1	FOXP3	FOXO3A	FOXL1	FOXK1	FOXA2	FOS	FOLR1	FN1	FLT3	FLT1	FLI1
B	FLCN	FGFR3	FGF7	FGF5	FGF4	FGF2	FGF1	FBLN2	FBLN1	FASLG	FAS	F2
C	EZH2	ETV5	ESR1	EPHB2	EPHA2	EPAS1	EOMES	EMP1	ELOVL6	EGR3	EGFR	E2F1
D	DPPA4	DPPA2	DNMT3L	DNMT3A	DLG7	DKK3	DKC1	DES	CYP7A1	CYP26A1	CXCR4	CXCL12
E	CTNNB1	CSF3R	CSF3	CSF2	CSF1R	CSF1	CPT1A	COL3A1	COL2A1	COL1A1	CFD	CEBPA
F	CDX2	CDKN2C	CDKN2A	CDKN1B	CDKN1A	CDH5	CDC42	CD96	CD9	CD68	CD44	CD4
G	CD34	CD19	CCND2	CCND1	CCL5	CBS	CASP3	CADM1	BYSL	BRCA1	BMPR2	BMP4
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG049 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	BMP2	BMI1	BHMT2	BCR	BCL2	ARNT	APP	APC	ANXA5	ANPEP	ALPL	AKT1
B	AGT	AFP	ADIPOQ	ACE	ACAN	ABCG2	ABCB1	CDX4	EPO	FST	GBX2	IL9
C	MEPE	MYOD1	NR0B1	NR2E1	COL9A1	MSX1	NEUROG2	ADCYAP1	ANXA2	CAT	LTB4R	CLDN7
D	DAB2	DNMT1	DNMT3B	DPPA3	DPPA5	FGF13	FGF3	FGFR1	FLT3LG	FOXO3	GDF3	GJA3
E	GYPB	H2AFX	HAND1	HOXB6	IL11	IL3	KLF2	NCAM1	NEUROD1	NFYA	NR6A1	OCIAD1
F	PDX1	PHB2	PHC1	PITX2	PIWIL1	PSCA	PUM1	RUNX1	SC4MOL	SMN1	STMN2	SUMO1
G	SUZ12	TERT	USP9X	ASCL2	CDH1	DHH	MYST2	CD8A	PARD6A	BMP1	NCAM1	CBX4
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG049 plate 05

- **Gene primer pairs:** 84 wells (A row to G row) are designated for a real-time PCR assay for genes (see the primer list).
- **HK1-6:** Six pre-deposited housekeeping gene (HK1-6) primer pairs, which can be used as endogenous positive controls as well as for array normalization.
- **GDC:** Genomic DNA controls, which can be used to specifically detect genomic DNA contamination with a high level of sensitivity.
- **RT:** Spike-in reverse transcription controls, which can be used to monitor the efficiency of the RT reactions. These pre-deposited primer pairs specifically amplify the cDNA template reversed transcribed from the spike-in control RNA in the sample.
- **PCR:** Positive PCR controls, which are used to verify the PCR efficiency by amplifying the pre-deposited DNA template with its specific pre-deposited primer pairs.

Gene primer list

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	SYT
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		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Human Stem Cell Related Gene qPCR Array (the Product). If the terms and conditions are not acceptable, the Product in its entirety must be returned to GeneCopoeia within 5 calendar days. A limited End-User license is granted to the purchaser of the Product. The Product shall be used by the purchaser for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products or deliver information obtained in service without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund. This limited warranty shall not extend to anyone other than the original purchaser of the Product. Notice of nonconforming products must be made to GeneCopoeia within 30 days of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product or a refund limited to the actual purchase price.

GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2016 GeneCopoeia, Inc.

GeneCopoeia, Inc.
9620 Medical Center Drive, Suite 101
Rockville, MD 20850
+1 (301) 762-0888
+1 (866) 360-9531
inquiry@genecopoeia.com