

ExProfile™ Human Surface Antigens (CD) Related Gene qPCR Array

For focused group profiling of human surface antigens (CD) genes expression

Cat. No. QG097-A (3 x 96-well plate, Format A)

Cat. No. QG097-B (3 x 96-well plate, Format B)

Cat. No. QG097-C (3 x 96-well plate, Format C)

Cat. No. QG097-D (3 x 96-well plate, Format D)

Cat. No. QG097-E (3 x 96-well plate, Format E)

Plates available individually or as a set of 6. Each set contains 252 unique gene primer pairs deposited in one 96-well plate.

Introduction

The ExProfile human surface antigens (CD) related gene qPCR array profiles the expression of 252 human genes related to surface antigens (CD). These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including genes that encode various surface antigens (CD) and related molecules involved in a series of important physiological and pathological processes, such as cell growth and differentiation, immune response, inflammation and so on. This array allows researchers to study the related genes to gain understanding of their roles in the different physiological and pathological processes.

- QG097 plate 01: 84 unique gene PCR primer pairs
- QG097 plate 02: 84 unique gene PCR primer pairs
- QG097 plate 03: 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	BTN3A1	CD300C	CD3EAP	TNFSF13B	CD226	CXCR6	PROCR	SEMA4D	TLR6	CLEC4M	CD96	MRC2
B	NCR1	ABCG2	IGSF2	CD163	CD83	PROM1	CD84	TNFRSF10A	TNFRSF10B	TNFRSF10C	TNFRSF10D	TNFRSF11A
C	SIGLEC5	TNFSF10	TNFSF14	IFITM1	SEMA7A	FZD9	FZD4	CXCR4	VCAM1	TNFSF4	SPAN7	TLR4
D	TLR3	TLR2	TLR1	THY1	THBD	TFRC	SPN	SLC4A1	SLAMF1	ST6GAL1	SELPLG	SELP
E	SELL	SDC1	RHAG	PVRL2	PVRL1	PVR	PTPRJ	PTPRC	PTGFRN	PRNP	PLAUR	ABCB1
F	PECAM1	PDCD1	NT5E	NGFR	MUC1	MST1R	MRC1	MPL	CD200	MME	CD99	MF12
G	CD46	MCAM	TACSTD1	LY75	CD180	LY9	BCAM	LTBR	LRP1	LIFR	LEPR	LAIR2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG097 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	LAIR1	L1CAM	KLRD1	KLRB1	KIT	KIR3DL1	KIR2DL1	KEL	KDR	ITGB4	ITGB3	ITGB2
B	ITGB1	ITGAV	ITGAE	ITGAD	ITGA2B	ISG20	INSR	TNFRSF9	IL12RB1	IL9R	IL8RB	IL8RA
C	IL7R	IL6ST	IL6R	IL5RA	IL4R	IL3RA	IL2RG	IL2RB	IL2RA	IGF2R	IGF1R	IFNGR1
D	ICAM4	ICAM2	ICAM1	HMMR	GYPC	CXCR3	GP1BA	DARC	FUT4	FLT3	FGFR4	FGFR2
E	FGFR3	FCGR3A	FCGR2B	FCGR2A	FCGR1A	FCER2	F3	ERBB2	ENPEP	ENG	DPP4	ACE
F	CD55	CTLA4	CSF3R	CSF2RB	CSF1R	CR2	CCR7	CCR6	CCR5	CCR1	CEACAM8	CEACAM5
G	CD52	CDH5	CDH2	CD151	CD97	CD81	CD79B	CD72	CD70	CD69	CD68	CD63
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG097 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	CD59	CD58	CD48	CD44	CD40LG	CD40	ENTPD5	ENTPD3	ENTPD6	ENTPD2	ENTPD1	CD38
B	CD37	SCARB2	CD36	CD34	TNFSF8	TNFRSF8	CD86	CD80	CD28	CD27	CD22	CD19
C	CD14	CD9	CD8B	CD7	CD6	CD5L	CD5	CD4	CD247	CD3G	CD3E	CD3D
D	CD2	CD1E	CD1D	CD1C	CD1B	CD1A	DDR1	C5AR1	BST2	BST1	BSG	BMPR1A
E	BLR1	CEACAM1	TNFRSF17	ATP1B3	ART4	FASLG	FAS	ANPEP	ALK	JAG1	ART1	CCR3
F	ICAM3	KIR2DL4	MSR1	ADAM8	LAG3	MS4A3	CD33	SIGLEC6	TNFSF11	CD53	CD74	CD79A
G	CDH1	CEACAM3	CR1	CSF2RA	FCAR	FGFR1	FUT3	GGT1	GP5	CD302	CD82	LILRB2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG097 plate 03

- **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
QG097-01	A01	HQP001281	NM_007048	BTN3A1
QG097-01	A02	HQP000996	NM_006678	CD300C
QG097-01	A03	HQP000971	NM_012099	CD3EAP
QG097-01	A04	HQP000821	NM_006573	TNFSF13B
QG097-01	A05	HQP000811	NM_006566	CD226
QG097-01	A06	HQP000808	NM_006564	CXCR6
QG097-01	A07	HQP000652	NM_006404	PROCR
QG097-01	A08	HQP000613	NM_006378	SEMA4D
QG097-01	A09	HQP000415	NM_006068	TLR6
QG097-01	A10	HQP000411	NM_014257	CLEC4M
QG097-01	A11	HQP000279	NM_005816	CD96
QG097-01	A12	HQP023346	NM_006039	MRC2
QG097-01	B01	HQP022750	NM_004829	NCR1
QG097-01	B02	HQP022745	NM_004827	ABCG2
QG097-01	B03	HQP022665	NM_004258	IGSF2
QG097-01	B04	HQP022548	NM_004244	CD163
QG097-01	B05	HQP022500	NM_001040280	CD83
QG097-01	B06	HQP021613	NM_006017	PROM1
QG097-01	B07	HQP021599	NM_003874	CD84
QG097-01	B08	HQP021557	NM_003844	TNFRSF10A

QG097-01	B09	HQP021553	NM_003842	TNFRSF10B
QG097-01	B10	HQP021552	NM_003841	TNFRSF10C
QG097-01	B11	HQP021551	NM_003840	TNFRSF10D
QG097-01	B12	HQP021550	NM_003839	TNFRSF11A
QG097-01	C01	HQP021533	NM_003830	SIGLEC5
QG097-01	C02	HQP021502	NM_003810	TNFSF10
QG097-01	C03	HQP021496	NM_003807	TNFSF14
QG097-01	C04	HQP021142	NM_003641	IFITM1
QG097-01	C05	HQP020964	NM_003612	SEMA7A
QG097-01	C06	HQP020141	NM_003508	FZD9
QG097-01	C07	HQP020134	NM_012193	FZD4
QG097-01	C08	HQP018802	NM_001008540	CXCR4
QG097-01	C09	HQP018466	NM_080682	VCAM1
QG097-01	C10	HQP018329	NM_003326	TNFSF4
QG097-01	C11	HQP018120	NM_004615	TSPAN7
QG097-01	C12	HQP018116	NM_138554	TLR4
QG097-01	D01	HQP018115	NM_003265	TLR3
QG097-01	D02	HQP018114	NM_003264	TLR2
QG097-01	D03	HQP018113	NM_003263	TLR1
QG097-01	D04	HQP018083	NM_006288	THY1
QG097-01	D05	HQP018067	NM_000361	THBD
QG097-01	D06	HQP018041	NM_003234	TFRC
QG097-01	D07	HQP017670	NM_001030288	SPN
QG097-01	D08	HQP017365	NM_000342	SLC4A1
QG097-01	D09	HQP017303	NM_003037	SLAMF1
QG097-01	D10	HQP017179	NM_003032	ST6GAL1
QG097-01	D11	HQP016747	NM_003006	SELPLG
QG097-01	D12	HQP016746	NM_003005	SELP
QG097-01	E01	HQP016745	NM_000655	SELL
QG097-01	E02	HQP016659	NM_001006946	SDC1
QG097-01	E03	HQP016270	NM_000324	RHAG
QG097-01	E04	HQP015962	NM_001042724	PVRL2
QG097-01	E05	HQP015956	NM_002855	PVRL1
QG097-01	E06	HQP015954	NM_006505	PVR
QG097-01	E07	HQP015925	NM_002843	PTPRJ
QG097-01	E08	HQP015908	NM_002838	PTPRC
QG097-01	E09	HQP015570	NM_020440	PTGFRN
QG097-01	E10	HQP015032	NM_000311	PRNP
QG097-01	E11	HQP013205	NM_001005376	PLAUR
QG097-01	E12	HQP013100	NM_000927	ABCB1
QG097-01	F01	HQP013015	NM_000442	PECAM1
QG097-01	F02	HQP012662	NM_005018	PDCD1
QG097-01	F03	HQP011932	NM_002526	NT5E

QG097-01	F04	HQP011828	NM_002507	NGFR
QG097-01	F05	HQP011559	NM_001018016	MUC1
QG097-01	F06	HQP011525	NM_002447	MST1R
QG097-01	F07	HQP011319	NM_002438	MRC1
QG097-01	F08	HQP011308	NM_005373	MPL
QG097-01	F09	HQP011301	NM_001004196	CD200
QG097-01	F10	HQP011254	NM_000902	MME
QG097-01	F11	HQP011212	NM_002414	CD99
QG097-01	F12	HQP011189	NM_005929	MFI2
QG097-01	G01	HQP011113	NM_002389	CD46
QG097-01	G02	HQP011099	NM_006500	MCAM
QG097-01	G03	HQP010942	NM_002354	TACSTD1
QG097-01	G04	HQP010933	NM_002349	LY75
QG097-01	G05	HQP010932	NM_005582	CD180
QG097-01	G06	HQP010930	NM_001033667	LY9
QG097-01	G07	HQP010922	NM_001013257	BCAM
QG097-01	G08	HQP010915	NM_002342	LTBR
QG097-01	G09	HQP010870	NM_002332	LRP1
QG097-01	G10	HQP010608	NM_002310	LIFR
QG097-01	G11	HQP010582	NM_001003679	LEPR
QG097-01	G12	HQP010478	NM_002288	LAIR2
QG097-01	H01	HGDC		
QG097-01	H02	HGDC		
QG097-01	H03	HQP006940	NM_002046	GAPDH
QG097-01	H04	HQP016381	NM_001101	ACTB
QG097-01	H05	HQP015171	NM_004048	B2M
QG097-01	H06	HQP006171	NM_012423	RPL13A
QG097-01	H07	HQP009026	NM_000194	HPRT1
QG097-01	H08	HQP054253	NR_003286	RN18S1
QG097-01	H09	RT		
QG097-01	H10	RT		
QG097-01	H11	PCR		
QG097-01	H12	PCR		
QG097-02	A01	HQP010467	NM_002287	LAIR1
QG097-02	A02	HQP010390	NM_000425	L1CAM
QG097-02	A03	HQP010112	NM_002262	KLRD1
QG097-02	A04	HQP010106	NM_002258	KLRB1
QG097-02	A05	HQP010099	NM_000222	KIT
QG097-02	A06	HQP010095	NM_013289	KIR3DL1
QG097-02	A07	HQP010083	NM_014218	KIR2DL1
QG097-02	A08	HQP010071	NM_000420	KEL
QG097-02	A09	HQP010070	NM_002253	KDR
QG097-02	A10	HQP009819	NM_000213	ITGB4

QG097-02	A11	HQP009818	NM_000212	ITGB3
QG097-02	A12	HQP009815	NM_000211	ITGB2
QG097-02	B01	HQP009810	NM_002211	ITGB1
QG097-02	B02	HQP009808	NM_002210	ITGAV
QG097-02	B03	HQP009805	NM_002208	ITGAE
QG097-02	B04	HQP009804	NM_005353	ITGAD
QG097-02	B05	HQP009795	NM_000419	ITGA2B
QG097-02	B06	HQP009789	NM_002201	ISG20
QG097-02	B07	HQP009764	NM_000208	INSR
QG097-02	B08	HQP009716	NM_001561	TNFRSF9
QG097-02	B09	HQP009694	NM_005535	IL12RB1
QG097-02	B10	HQP009683	NM_002186	IL9R
QG097-02	B11	HQP009681	NM_001557	IL8RB
QG097-02	B12	HQP009679	NM_000634	IL8RA
QG097-02	C01	HQP009677	NM_002185	IL7R
QG097-02	C02	HQP009674	NM_002184	IL6ST
QG097-02	C03	HQP009672	NM_000565	IL6R
QG097-02	C04	HQP009667	NM_000564	IL5RA
QG097-02	C05	HQP009664	NM_000418	IL4R
QG097-02	C06	HQP009661	NM_002183	IL3RA
QG097-02	C07	HQP009659	NM_000206	IL2RG
QG097-02	C08	HQP009658	NM_000878	IL2RB
QG097-02	C09	HQP009650	NM_000417	IL2RA
QG097-02	C10	HQP009532	NM_000876	IGF2R
QG097-02	C11	HQP009523	NM_000875	IGF1R
QG097-02	C12	HQP009469	NM_000416	IFNGR1
QG097-02	D01	HQP009201	NM_001039132	ICAM4
QG097-02	D02	HQP009187	NM_000873	ICAM2
QG097-02	D03	HQP009184	NM_000201	ICAM1
QG097-02	D04	HQP008896	NM_012484	HMMR
QG097-02	D05	HQP008634	NM_002101	GYPC
QG097-02	D06	HQP007900	NM_001504	CXCR3
QG097-02	D07	HQP007803	NM_000173	GP1BA
QG097-02	D08	HQP006479	NM_002036	DARC
QG097-02	D09	HQP006455	NM_002033	FUT4
QG097-02	D10	HQP005890	NM_004119	FLT3
QG097-02	D11	HQP005439	NM_002011	FGFR4
QG097-02	D12	HQP005437	NM_000141	FGFR2
QG097-02	E01	HQP005434	NM_000142	FGFR3
QG097-02	E02	HQP005308	NM_000569	FCGR3A
QG097-02	E03	HQP005290	NM_001002273	FCGR2B
QG097-02	E04	HQP005280	NM_021642	FCGR2A
QG097-02	E05	HQP005251	NM_000566	FCGR1A

QG097-02	E06	HQP005242	NM_002002	FCER2
QG097-02	E07	HQP005057	NM_001993	F3
QG097-02	E08	HQP004968	NM_001005862	ERBB2
QG097-02	E09	HQP004867	NM_001977	ENPEP
QG097-02	E10	HQP004856	NM_000118	ENG
QG097-02	E11	HQP004434	NM_001935	DPP4
QG097-02	E12	HQP004081	NM_000789	ACE
QG097-02	F01	HQP003948	NM_000574	CD55
QG097-02	F02	HQP003499	NM_001037631	CTLA4
QG097-02	F03	HQP003185	NM_000760	CSF3R
QG097-02	F04	HQP003170	NM_000395	CSF2RB
QG097-02	F05	HQP003158	NM_005211	CSF1R
QG097-02	F06	HQP002889	NM_001006658	CR2
QG097-02	F07	HQP002217	NM_001838	CCR7
QG097-02	F08	HQP002212	NM_004367	CCR6
QG097-02	F09	HQP002210	NM_000579	CCR5
QG097-02	F10	HQP002198	NM_001295	CCR1
QG097-02	F11	HQP001014	NM_001816	CEACAM8
QG097-02	F12	HQP000593	NM_004363	CEACAM5
QG097-02	G01	HQP000525	NM_001803	CD52
QG097-02	G02	HQP000052	NM_001795	CDH5
QG097-02	G03	HQP000018	NM_001792	CDH2
QG097-02	G04	HQP023192	NM_001039490	CD151
QG097-02	G05	HQP023178	NM_001025160	CD97
QG097-02	G06	HQP023168	NM_004356	CD81
QG097-02	G07	HQP023155	NM_000626	CD79B
QG097-02	G08	HQP023117	NM_001782	CD72
QG097-02	G09	HQP023108	NM_001252	CD70
QG097-02	G10	HQP023094	NM_001781	CD69
QG097-02	G11	HQP023083	NM_001040059	CD68
QG097-02	G12	HQP023069	NM_001040034	CD63
QG097-02	H01	HGDC		
QG097-02	H02	HGDC		
QG097-02	H03	HQP006940	NM_002046	GAPDH
QG097-02	H04	HQP016381	NM_001101	ACTB
QG097-02	H05	HQP015171	NM_004048	B2M
QG097-02	H06	HQP006171	NM_012423	RPL13A
QG097-02	H07	HQP009026	NM_000194	HPRT1
QG097-02	H08	HQP054253	NR_003286	RN18S1
QG097-02	H09	RT		
QG097-02	H10	RT		
QG097-02	H11	PCR		
QG097-02	H12	PCR		

QG097-03	A01	HQP023056	NM_000611	CD59
QG097-03	A02	HQP023046	NM_001779	CD58
QG097-03	A03	HQP023005	NM_001778	CD48
QG097-03	A04	HQP022972	NM_000610	CD44
QG097-03	A05	HQP022962	NM_000074	CD40LG
QG097-03	A06	HQP022955	NM_001250	CD40
QG097-03	A07	HQP022934	NM_001249	ENTPD5
QG097-03	A08	HQP022921	NM_001248	ENTPD3
QG097-03	A09	HQP022912	NM_001247	ENTPD6
QG097-03	A10	HQP022894	NM_001246	ENTPD2
QG097-03	A11	HQP022883	NM_001776	ENTPD1
QG097-03	A12	HQP022870	NM_001775	CD38
QG097-03	B01	HQP022855	NM_001040031	CD37
QG097-03	B02	HQP022846	NM_005506	SCARB2
QG097-03	B03	HQP022821	NM_000072	CD36
QG097-03	B04	HQP022812	NM_001025109	CD34
QG097-03	B05	HQP022769	NM_001244	TNFSF8
QG097-03	B06	HQP022753	NM_001243	TNFRSF8
QG097-03	B07	HQP022746	NM_006889	CD86
QG097-03	B08	HQP022722	NM_005191	CD80
QG097-03	B09	HQP022699	NM_006139	CD28
QG097-03	B10	HQP022667	NM_001242	CD27
QG097-03	B11	HQP022561	NM_001771	CD22
QG097-03	B12	HQP022504	NM_001770	CD19
QG097-03	C01	HQP022490	NM_000591	CD14
QG097-03	C02	HQP022474	NM_001769	CD9
QG097-03	C03	HQP022438	NM_004931	CD8B
QG097-03	C04	HQP022399	NM_006137	CD7
QG097-03	C05	HQP022384	NM_006725	CD6
QG097-03	C06	HQP022357	NM_005894	CD5L
QG097-03	C07	HQP022337	NM_014207	CD5
QG097-03	C08	HQP022316	NM_000616	CD4
QG097-03	C09	HQP022295	NM_000734	CD247
QG097-03	C10	HQP022256	NM_000073	CD3G
QG097-03	C11	HQP022236	NM_000733	CD3E
QG097-03	C12	HQP022212	NM_000732	CD3D
QG097-03	D01	HQP022190	NM_001767	CD2
QG097-03	D02	HQP022159	NM_001042583	CD1E
QG097-03	D03	HQP022129	NM_001766	CD1D
QG097-03	D04	HQP022111	NM_001765	CD1C
QG097-03	D05	HQP022090	NM_001764	CD1B
QG097-03	D06	HQP022070	NM_001763	CD1A
QG097-03	D07	HQP018768	NM_001954	DDR1

QG097-03	D08	HQP018326	NM_001736	C5AR1
QG097-03	D09	HQP017844	NM_004335	BST2
QG097-03	D10	HQP017834	NM_004334	BST1
QG097-03	D11	HQP017820	NM_001728	BSG
QG097-03	D12	HQP017489	NM_004329	BMPR1A
QG097-03	E01	HQP016959	NM_001716	BLR1
QG097-03	E02	HQP016623	NM_001024912	CEACAM1
QG097-03	E03	HQP016367	NM_001192	TNFRSF17
QG097-03	E04	HQP011863	NM_001679	ATP1B3
QG097-03	E05	HQP011153	NM_021071	ART4
QG097-03	E06	HQP009671	NM_000639	FASLG
QG097-03	E07	HQP009651	NM_000043	FAS
QG097-03	E08	HQP008414	NM_001150	ANPEP
QG097-03	E09	HQP006353	NM_004304	ALK
QG097-03	E10	HQP004470	NM_000214	JAG1
QG097-03	E11	HQP011127	NM_004314	ART1
QG097-03	E12	HQP002207	NM_001837	CCR3
QG097-03	F01	HQP009192	NM_002162	ICAM3
QG097-03	F02	HQP010087	NM_001080770	KIR2DL4
QG097-03	F03	HQP011520	NM_002445	MSR1
QG097-03	F04	HQP000246	NM_001109	ADAM8
QG097-03	F05	HQP010462	NM_002286	LAG3
QG097-03	F06	HQP022541	NM_001031666	MS4A3
QG097-03	F07	HQP021167	NM_001082618	CD33
QG097-03	F08	HQP022797	NM_001245	SIGLEC6
QG097-03	F09	HQP021321	NM_003701	TNFSF11
QG097-03	F10	HQP023019	NM_000560	CD53
QG097-03	F11	HQP023127	NM_001025158	CD74
QG097-03	F12	HQP023144	NM_001783	CD79A
QG097-03	G01	HQP023466	NM_004360	CDH1
QG097-03	G02	HQP000972	NM_001815	CEACAM3
QG097-03	G03	HQP002880	NM_000573	CR1
QG097-03	G04	HQP003163	NM_006140	CSF2RA
QG097-03	G05	HQP005227	NM_002000	FCAR
QG097-03	G06	HQP005427	NM_015850	FGFR1
QG097-03	G07	HQP006454	NM_000149	FUT3
QG097-03	G08	HQP007378	NM_001032364	GGT1
QG097-03	G09	HQP007809	NM_004488	GP5
QG097-03	G10	HQP023380	NM_014880	CD302
QG097-03	G11	HQP009860	NM_001024844	CD82
QG097-03	G12	HQP000353	NM_001080978	LILRB2
QG097-03	H01	HGDC		
QG097-03	H02	HGDC		

QG097-03	H03	HQP006940	NM_002046	GAPDH
QG097-03	H04	HQP016381	NM_001101	ACTB
QG097-03	H05	HQP015171	NM_004048	B2M
QG097-03	H06	HQP006171	NM_012423	RPL13A
QG097-03	H07	HQP009026	NM_000194	HPRT1
QG097-03	H08	HQP054253	NR_003286	RN18S1
QG097-03	H09	RT		
QG097-03	H10	RT		
QG097-03	H11	PCR		
QG097-03	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Human Surface Antigens (CD) 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