

ExProfile™ Human Lymphoma Gene qPCR Array

For focused group profiling of human lymphoma related genes expression

Cat. No. QG073-A (5 x 96-well plate, Format A)

Cat. No. QG073-B (5 x 96-well plate, Format B)

Cat. No. QG073-C (5 x 96-well plate, Format C)

Cat. No. QG073-D (5 x 96-well plate, Format D)

Cat. No. QG073-E (5 x 96-well plate, Format E)

Available as 1 set or 6 sets. Each set contains 420 unique gene primers deposited in five 96-well plates.

Introduction

The ExProfile human lymphoma gene qPCR array profiles 420 human genes to aberrantly expressed human genes involved in human lymphoma. These genes are carefully chosen for their close cancer correlation based on a thorough literature search of peer-reviewed publications. Abnormal gene expression is often observed in cancer development and progression. The ExProfile human lymphoma gene array allows researchers to study the cancer-related genes to gain understanding of their roles in lymphoma pathogenesis.

- QG073 plate 01: 84 unique gene PCR primer pairs
- QG073 plate 02: 84 unique gene PCR primer pairs
- QG073 plate 03: 84 unique gene PCR primer pairs
- QG073 plate 04: 84 unique gene PCR primer pairs
- QG073 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

GeneCopia 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	MTHFR	IL10	TNF	CYP1A1	NQO1	HLA-DRB1	MTR	CYP2E1	XRCC1	TP53	TYMS	HLA-A
B	ERCC2	TPMT	FCGR2A	NAT2	MPO	IL4	NAT1	ABCB1	MTRR	NOD2	IL6	EPHX1
C	CTLA4	VDR	SOD2	SLC19A1	RFC1	BCL2	NBN	MYC	MDM2	IL4R	IL1RN	IL1B
D	CYP2D6	CYP2C9	CBS	BRCA2	FBXW7	NOS3	GSTM3	CYP1B1	CDKN2A	OGG1	ERCC5	CD40
E	XRCC3	VEGFA	TGFB1	CCL5	CCND1	NOS2A	NFKB1	LEP	IL12B	IL12A	IL8	IL2
F	IL1A	CD14	GGH	AKR1C3	ARID5B	CASP3	CASP8	XRCC2	WRN	VCAM1	TLR1	STAT3
G	SOD1	CXCL12	PTGS2	PTEN	TLR9	MSH2	MRE11A	LIG4	LEPR	JAK2	IRF4	IL13
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG073 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	IL5	APOE	FPGS	FLT3	F5	AHR	DHFR	CYP21A2	CYP17A1	CYP3A5	CX3CR1	CCR5
B	ALDH1L1	MTHFS	AKR1A1	RAD50	LOC729230	PRF1	CD38	CD28	MS4A1	SOCS3	RIPK1	CASP7
C	CASP6	CASP5	CASP4	CASP1	TLR10	XRCC4	XDTH	C8B	C7	TYR	C5AR1	C4BPB
D	C4BPA	C2	C1QC	TNFRSF1B	TCN2	TCN1	STAT4	STAT1	BLM	SELE	BHMT	CCL2
E	CFB	SERPING1	BCL2L2	RAC1	BAK1	BAD	PRL	PON1	GHRL	C1RL	SERPINE1	NPY
F	NFKBIB	NFKBIA	NCF4	NCF2	MYD88	MMP2	MIF	CD46	KRAS	KIR3DL1	KIR2DS3	KIR2DL3
G	KIR2DL1	JAK3	IL15	IL8RB	IL7R	IFNGR2	APOA2	NR3C1	GPX4	ATG9B	CD93	FCER2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG073 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	F2	ESR1	ERCC6	ERCC1	EDN1	CYP2B6	ADR2	PARP1	COMT	TLR6	C1QB	SOD3
B	NOS1	BCL2L10	ZNF527	ICHTHYIN	TGFBR1	BAG5	SHBG	CASP8AP2	SLC23A1	CDC6	CDA	BCLAF1
C	ESPL1	CLCA2	KLK4	TP53I3	BAG3	GSTO1	CD80	ADIPOQ	DLG5	CD5	TNFSF18	MBD2
D	BCL10	STK19	CFLAR	SOCS2	TNFRSF10A	RIPK2	TNFRSF14	ADAM19	TRADD	SERPINA6	SOCS1	RUNX1
E	TSLP	CNTNAP1	PARG	FOXN1	RAD54L	ACAD11	STARD3NL	BCL2L12	NCOA3	DEFB126	CFHR5	BCL2L14
F	BAT3	C7orf24	CXCR4	DDR1	PTP4A1	XPC	NR1H2	SCGB1A1	C8G	TYK2	TRAF6	TRAF5
G	TRAF2	TOP2A	TLR5	TLR3	C1QBP	ADAM17	STAT2	STAR	BPI	SRD5A1	BNIP3	SLC2A11
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG073 plate 03

Product Data Sheet

	1	2	3	4	5	6	7	8	9	10	11	12
A	SLC13A1	ACBD3	SELPLG	SELP	SECTM1	CCL24	CCL20	CCL18	CCL13	CCL11	CCL1	SERPINB3
B	BCL7A	REL	RB1	RASA1	RAD23B	PTPN1	SFRS15	PTGS1	AICDA	PTGDR	GPR108	DHX33
C	KLK6	KLK7	PARD3	FOXJ2	KLK15	USP47	ACP5	KIAA1546	PPARG	TOLLIP	POMC	POLB
D	PIM1	PIK3CA	SERPINI1	PGR	PFKFB2	IRAK4	PHF11	PAX5	PAX4	DEF6	SERPINB2	OCA2
E	NRAS	NPAS2	NNMT	NDUFS2	NAB2	SERPINC1	MTAP	ABCC1	MMP10	MMP9	MMP3	MMP1
F	MLH1	MITF	MC4R	MBP	MAL	CD180	LRP5	LPL	LMAN1	LIPC	LIG3	LIG1
G	LBP	LAMA2	KLKB1	KLK1	JAK1	ITPA	AR	IRAK2	INHBC	IL18	IL15RA	IL12RB2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG073 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	IL12RB1	IL8RA	IL2RA	IKBKB	IGFBP3	IGFBP1	IGF2	IGF1R	IGF1	IFNGR1	IDH1	APOB
B	ICAM4	APOA4	HSPA6	HSD17B2	HSD17B1	APEX1	HSD3B2	HSD3B1	HP	HMGCS2	HMGCS1	ACAA1
C	CD209	HGF	IL19	ICOS	GTF2E1	GRIA1	GPX3	GPX2	SLCO1B3	GLS	GH1	SENP3
D	ALOX15B	ALOX15	DEFB119	ALOX5AP	ALOX5	ALOX12	BCL2L13	LY96	PDSS1	CASP14	LILRA4	TRAM1
E	SIRT1	FLT1	FGF2	BZRPL1	FDFT1	FCGR2B	FCER1G	MS4A2	FCER1A	ETV6	ESR2	AKT1
F	ERCC4	ERCC3	EPHX2	AIF1	DRD5	DRD2	ZNF169	CFD	DEFA5	DEFA4	ACE	DCK
G	DAXX	CYP19A1	ADRB3	CYP2C19	CTNNB1	TICAM1	HIPK4	CSF3	CRY2	DEFB127	CR1	CPT2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG073 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 reverse 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
QG073-01	A01	HQP011547	NM_005957	MTHFR
QG073-01	A02	HQP009685	NM_000572	IL10
QG073-01	A03	HQP018141	NM_000594	TNF
QG073-01	A04	HQP003772	NM_000499	CYP1A1
QG073-01	A05	HQP004317	NM_000903	NQO1
QG073-01	A06	HQP054047	BC008403	HLA-DRB1
QG073-01	A07	HQP011554	NM_000254	MTR
QG073-01	A08	HQP003817	NM_000773	CYP2E1
QG073-01	A09	HQP018562	NM_006297	XRCC1
QG073-01	A10	HQP018175	NM_000546	TP53
QG073-01	A11	HQP018342	NM_001071	TYMS
QG073-01	A12	HQP008849	NM_002116	HLA-A
QG073-01	B01	HQP004976	NM_000400	ERCC2
QG073-01	B02	HQP018213	NM_000367	TPMT

Product Data Sheet

QG073-01	B03	HQP005280	NM_021642	FCGR2A
QG073-01	B04	HQP001136	NM_000015	NAT2
QG073-01	B05	HQP011309	NM_000250	MPO
QG073-01	B06	HQP009662	NM_000589	IL4
QG073-01	B07	HQP023467	NM_000662	NAT1
QG073-01	B08	HQP013100	NM_000927	ABCB1
QG073-01	B09	HQP011555	NM_002454	MTRR
QG073-01	B10	HQP016801	NM_022162	NOD2
QG073-01	B11	HQP009670	NM_000600	IL6
QG073-01	B12	HQP004948	NM_000120	EPHX1
QG073-01	C01	HQP003499	NM_001037631	CTLA4
QG073-01	C02	HQP018474	NM_000376	VDR
QG073-01	C03	HQP017616	NM_000636	SOD2
QG073-01	C04	HQP017482	NM_194255	SLC19A1
QG073-01	C05	HQP016224	NM_002913	RFC1
QG073-01	C06	HQP016212	NM_000657	BCL2
QG073-01	C07	HQP011687	NM_002485	NBN
QG073-01	C08	HQP011597	NM_002467	MYC
QG073-01	C09	HQP011135	NM_002392	MDM2
QG073-01	C10	HQP009664	NM_000418	IL4R
QG073-01	C11	HQP009645	NM_000577	IL1RN
QG073-01	C12	HQP009641	NM_000576	IL1B
QG073-01	D01	HQP003814	NM_000106	CYP2D6
QG073-01	D02	HQP003811	NM_000771	CYP2C9
QG073-01	D03	HQP021518	NM_000071	CBS
QG073-01	D04	HQP017753	NM_000059	BRCA2
QG073-01	D05	HQP014292	NM_018315	FBXW7
QG073-01	D06	HQP011868	NM_000603	NOS3
QG073-01	D07	HQP008483	NM_000849	GSTM3
QG073-01	D08	HQP003775	NM_000104	CYP1B1
QG073-01	D09	HQP000370	NM_058195	CDKN2A
QG073-01	D10	HQP012021	NM_002542	OGG1
QG073-01	D11	HQP004985	NM_000123	ERCC5
QG073-01	D12	HQP022955	NM_001250	CD40
QG073-01	E01	HQP018564	NM_005432	XRCC3
QG073-01	E02	HQP018475	NM_001025366	VEGFA
QG073-01	E03	HQP018044	NM_000660	TGFB1
QG073-01	E04	HQP016626	NM_002985	CCL5
QG073-01	E05	HQP016204	NM_053056	CCND1
QG073-01	E06	HQP011866	NM_000625	NOS2A
QG073-01	E07	HQP011807	NM_003998	NFKB1
QG073-01	E08	HQP010581	NM_000230	LEP
QG073-01	E09	HQP009693	NM_002187	IL12B
QG073-01	E10	HQP009692	NM_000882	IL12A
QG073-01	E11	HQP009678	NM_000584	IL8
QG073-01	E12	HQP009649	NM_000586	IL2
QG073-01	F01	HQP009640	NM_000575	IL1A
QG073-01	F02	HQP022490	NM_000591	CD14

Product Data Sheet

QG073-01	F03	HQP021603	NM_003878	GGH
QG073-01	F04	HQP021385	NM_003739	AKR1C3
QG073-01	F05	HQP020515	NM_032199	ARID5B
QG073-01	F06	HQP020297	NM_004346	CASP3
QG073-01	F07	HQP018966	NM_001080124	CASP8
QG073-01	F08	HQP018563	NM_005431	XRCC2
QG073-01	F09	HQP018545	NM_000553	WRN
QG073-01	F10	HQP018466	NM_080682	VCAM1
QG073-01	F11	HQP018113	NM_003263	TLR1
QG073-01	F12	HQP017767	NM_003150	STAT3
QG073-01	G01	HQP017615	NM_000454	SOD1
QG073-01	G02	HQP016670	NM_001033886	CXCL12
QG073-01	G03	HQP015598	NM_000963	PTGS2
QG073-01	G04	HQP015535	NM_000314	PTEN
QG073-01	G05	HQP013388	NM_017442	TLR9
QG073-01	G06	HQP011491	NM_000251	MSH2
QG073-01	G07	HQP011320	NM_005590	MRE11A
QG073-01	G08	HQP010613	NM_002312	LIG4
QG073-01	G09	HQP010584	NM_002303	LEPR
QG073-01	G10	HQP009850	NM_004972	JAK2
QG073-01	G11	HQP009781	NM_002460	IRF4
QG073-01	G12	HQP009697	NM_002188	IL13
QG073-01	H01	HGDC		
QG073-01	H02	HGDC		
QG073-01	H03	HQP006940	NM_002046	GAPDH
QG073-01	H04	HQP016381	NM_001101	ACTB
QG073-01	H05	HQP015171	NM_004048	B2M
QG073-01	H06	HQP006171	NM_012423	RPL13A
QG073-01	H07	HQP009026	NM_000194	HPRT1
QG073-01	H08	HQP054253	NR_003286	RN18S1
QG073-01	H09	RT		
QG073-01	H10	RT		
QG073-01	H11	PCR		
QG073-01	H12	PCR		
QG073-02	A01	HQP009666	NM_000879	IL5
QG073-02	A02	HQP009556	NM_000041	APOE
QG073-02	A03	HQP006228	NM_001018078	FPGS
QG073-02	A04	HQP005890	NM_004119	FLT3
QG073-02	A05	HQP005058	NM_000130	F5
QG073-02	A06	HQP004658	NM_001621	AHR
QG073-02	A07	HQP004309	NM_000791	DHFR
QG073-02	A08	HQP003908	NM_000500	CYP21A2
QG073-02	A09	HQP003888	NM_000102	CYP17A1
QG073-02	A10	HQP003841	NM_000777	CYP3A5
QG073-02	A11	HQP003691	NM_001337	CX3CR1
QG073-02	A12	HQP002210	NM_000579	CCR5
QG073-02	B01	HQP000963	NM_012190	ALDH1L1
QG073-02	B02	HQP000709	NM_006441	MTHFS

Product Data Sheet

QG073-02	B03	HQP000405	NM_006066	AKR1A1
QG073-02	B04	HQP000145	NM_005732	RAD50
QG073-02	B05	HQP054686	NM_001123396	LOC729230
QG073-02	B06	HQP054058	NM_005041	PRF1
QG073-02	B07	HQP022870	NM_001775	CD38
QG073-02	B08	HQP022699	NM_006139	CD28
QG073-02	B09	HQP022527	NM_021950	MS4A1
QG073-02	B10	HQP021889	NM_003955	SOCS3
QG073-02	B11	HQP021492	NM_003804	RIPK1
QG073-02	B12	HQP020481	NM_033338	CASP7
QG073-02	C01	HQP020427	NM_001226	CASP6
QG073-02	C02	HQP020384	NM_004347	CASP5
QG073-02	C03	HQP020332	NM_001225	CASP4
QG073-02	C04	HQP020207	NM_001223	CASP1
QG073-02	C05	HQP019982	NM_001017388	TLR10
QG073-02	C06	HQP018565	NM_003401	XRCC4
QG073-02	C07	HQP018552	NM_000379	XDH
QG073-02	C08	HQP018380	NM_000066	C8B
QG073-02	C09	HQP018353	NM_000587	C7
QG073-02	C10	HQP018343	NM_000372	TYR
QG073-02	C11	HQP018326	NM_001736	C5AR1
QG073-02	C12	HQP018276	NM_000716	C4BPB
QG073-02	D01	HQP018259	NM_000715	C4BPA
QG073-02	D02	HQP018226	NM_000063	C2
QG073-02	D03	HQP018170	NM_172369	C1QC
QG073-02	D04	HQP018149	NM_001066	TNFRSF1B
QG073-02	D05	HQP017979	NM_000355	TCN2
QG073-02	D06	HQP017978	NM_001062	TCN1
QG073-02	D07	HQP017770	NM_003151	STAT4
QG073-02	D08	HQP017764	NM_007315	STAT1
QG073-02	D09	HQP016844	NM_000057	BLM
QG073-02	D10	HQP016744	NM_000450	SELE
QG073-02	D11	HQP016634	NM_001713	BHMT
QG073-02	D12	HQP016621	NM_002982	CCL2
QG073-02	E01	HQP016565	NM_001710	CFB
QG073-02	E02	HQP016500	NM_001032295	SERPING1
QG073-02	E03	HQP016260	NM_004050	BCL2L2
QG073-02	E04	HQP016064	NM_018890	RAC1
QG073-02	E05	HQP015917	NM_001188	BAK1
QG073-02	E06	HQP015538	NM_004322	BAD
QG073-02	E07	HQP015024	NM_000948	PRL
QG073-02	E08	HQP013473	NM_000446	PON1
QG073-02	E09	HQP012996	NM_016362	GHRL
QG073-02	E10	HQP012584	NM_016546	C1RL
QG073-02	E11	HQP012154	NM_000602	SERPINE1
QG073-02	E12	HQP011874	NM_000905	NPY
QG073-02	F01	HQP011812	NM_002503	NFKBIB
QG073-02	F02	HQP011810	NM_020529	NFKBIA

Product Data Sheet

QG073-02	F03	HQP011694	NM_000631	NCF4
QG073-02	F04	HQP011693	NM_000433	NCF2
QG073-02	F05	HQP011603	NM_002468	MYD88
QG073-02	F06	HQP011256	NM_004530	MMP2
QG073-02	F07	HQP011219	NM_002415	MIF
QG073-02	F08	HQP011113	NM_002389	CD46
QG073-02	F09	HQP010133	NM_004985	KRAS
QG073-02	F10	HQP010095	NM_013289	KIR3DL1
QG073-02	F11	HQP010092	NM_012313	KIR2DS3
QG073-02	F12	HQP010086	NM_015868	KIR2DL3
QG073-02	G01	HQP010083	NM_014218	KIR2DL1
QG073-02	G02	HQP009851	NM_000215	JAK3
QG073-02	G03	HQP009708	NM_000585	IL15
QG073-02	G04	HQP009681	NM_001557	IL8RB
QG073-02	G05	HQP009677	NM_002185	IL7R
QG073-02	G06	HQP009472	NM_005534	IFNGR2
QG073-02	G07	HQP009139	NM_001643	APOA2
QG073-02	G08	HQP008402	NM_001020825	NR3C1
QG073-02	G09	HQP008285	NM_002085	GPX4
QG073-02	G10	HQP008198	NM_173681	ATG9B
QG073-02	G11	HQP005589	NM_012072	CD93
QG073-02	G12	HQP005242	NM_002002	FCER2
QG073-02	H01	HGDC		
QG073-02	H02	HGDC		
QG073-02	H03	HQP006940	NM_002046	GAPDH
QG073-02	H04	HQP016381	NM_001101	ACTB
QG073-02	H05	HQP015171	NM_004048	B2M
QG073-02	H06	HQP006171	NM_012423	RPL13A
QG073-02	H07	HQP009026	NM_000194	HPRT1
QG073-02	H08	HQP054253	NR_003286	RN18S1
QG073-02	H09	RT		
QG073-02	H10	RT		
QG073-02	H11	PCR		
QG073-02	H12	PCR		
QG073-03	A01	HQP005052	NM_000506	F2
QG073-03	A02	HQP004998	NM_000125	ESR1
QG073-03	A03	HQP004986	NM_000124	ERCC6
QG073-03	A04	HQP004975	NM_202001	ERCC1
QG073-03	A05	HQP004557	NM_001955	EDN1
QG073-03	A06	HQP003808	NM_000767	CYP2B6
QG073-03	A07	HQP003791	NM_000024	ADRB2
QG073-03	A08	HQP003120	NM_001618	PARP1
QG073-03	A09	HQP002671	NM_000754	COMT
QG073-03	A10	HQP000415	NM_006068	TLR6
QG073-03	A11	HQP018159	NM_000491	C1QB
QG073-03	A12	HQP017618	NM_003102	SOD3
QG073-03	B01	HQP011865	NM_000620	NOS1
QG073-03	B02	HQP000027	NM_020396	BCL2L10

Product Data Sheet

QG073-03	B03	HQP054688	NM_032453	ZNF527
QG073-03	B04	HQP054326	NM_001099287	ICHTHYIN
QG073-03	B05	HQP054057	BC071181	TGFBR1
QG073-03	B06	HQP053998	NM_004873	BAG5
QG073-03	B07	HQP053959	NM_001040	SHBG
QG073-03	B08	HQP023464	NM_012115	CASP8AP2
QG073-03	B09	HQP023414	NM_005847	SLC23A1
QG073-03	B10	HQP023354	NM_001254	CDC6
QG073-03	B11	HQP023203	NM_001785	CDA
QG073-03	B12	HQP023186	NM_014739	BCLAF1
QG073-03	C01	HQP023095	NM_012291	ESPL1
QG073-03	C02	HQP023012	NM_006536	CLCA2
QG073-03	C03	HQP022998	NM_004917	KLK4
QG073-03	C04	HQP022884	NM_004881	TP53I3
QG073-03	C05	HQP022872	NM_004281	BAG3
QG073-03	C06	HQP022764	NM_004832	GSTO1
QG073-03	C07	HQP022722	NM_005191	CD80
QG073-03	C08	HQP022625	NM_004797	ADIPOQ
QG073-03	C09	HQP022361	NM_004747	DLG5
QG073-03	C10	HQP022337	NM_014207	CD5
QG073-03	C11	HQP021834	NM_005092	TNFSF18
QG073-03	C12	HQP021737	NM_003927	MBD2
QG073-03	D01	HQP021725	NM_003921	BCL10
QG073-03	D02	HQP021636	NM_032454	STK19
QG073-03	D03	HQP021604	NM_003879	CFLAR
QG073-03	D04	HQP021602	NM_003877	SOCS2
QG073-03	D05	HQP021557	NM_003844	TNFRSF10A
QG073-03	D06	HQP021524	NM_003821	RIPK2
QG073-03	D07	HQP021522	NM_003820	TNFRSF14
QG073-03	D08	HQP021484	NM_033274	ADAM19
QG073-03	D09	HQP021462	NM_003789	TRADD
QG073-03	D10	HQP021420	NM_001756	SERPINA6
QG073-03	D11	HQP021399	NM_003745	SOCS1
QG073-03	D12	HQP021347	NM_001754	RUNX1
QG073-03	E01	HQP021260	NM_033035	TSLP
QG073-03	E02	HQP021130	NM_003632	CNTNAP1
QG073-03	E03	HQP021129	NM_003631	PARG
QG073-03	E04	HQP020808	NM_003593	FOXN1
QG073-03	E05	HQP020707	NM_003579	RAD54L
QG073-03	E06	HQP020499	NM_032169	ACAD11
QG073-03	E07	HQP020390	NM_032016	STARD3NL
QG073-03	E08	HQP020245	NM_001040668	BCL2L12
QG073-03	E09	HQP020040	NM_006534	NCOA3
QG073-03	E10	HQP019941	NM_030931	DEFB126
QG073-03	E11	HQP019873	NM_030787	CFHR5
QG073-03	E12	HQP019021	NM_030766	BCL2L14
QG073-03	F01	HQP018983	NM_004639	BAT3
QG073-03	F02	HQP018854	NM_024051	C7orf24

Product Data Sheet

QG073-03	F03	HQP018802	NM_001008540	CXCR4
QG073-03	F04	HQP018768	NM_001954	DDR1
QG073-03	F05	HQP018760	NM_003463	PTP4A1
QG073-03	F06	HQP018556	NM_004628	XPC
QG073-03	F07	HQP018431	NM_007121	NR1H2
QG073-03	F08	HQP018411	NM_003357	SCGB1A1
QG073-03	F09	HQP018391	NM_000606	C8G
QG073-03	F10	HQP018340	NM_003331	TYK2
QG073-03	F11	HQP018237	NM_004620	TRAF6
QG073-03	F12	HQP018236	NM_001033910	TRAF5
QG073-03	G01	HQP018233	NM_021138	TRAF2
QG073-03	G02	HQP018172	NM_001067	TOP2A
QG073-03	G03	HQP018118	NM_003268	TLR5
QG073-03	G04	HQP018115	NM_003265	TLR3
QG073-03	G05	HQP018106	NM_001212	C1QBP
QG073-03	G06	HQP017866	NM_003183	ADAM17
QG073-03	G07	HQP017766	NM_005419	STAT2
QG073-03	G08	HQP017762	NM_000349	STAR
QG073-03	G09	HQP017702	NM_001725	BPI
QG073-03	G10	HQP017697	NM_001047	SRD5A1
QG073-03	G11	HQP017619	NM_004052	BNIP3
QG073-03	G12	HQP017549	NM_030807	SLC2A11
QG073-03	H01	HGDC		
QG073-03	H02	HGDC		
QG073-03	H03	HQP006940	NM_002046	GAPDH
QG073-03	H04	HQP016381	NM_001101	ACTB
QG073-03	H05	HQP015171	NM_004048	B2M
QG073-03	H06	HQP006171	NM_012423	RPL13A
QG073-03	H07	HQP009026	NM_000194	HPRT1
QG073-03	H08	HQP054253	NR_003286	RN18S1
QG073-03	H09	RT		
QG073-03	H10	RT		
QG073-03	H11	PCR		
QG073-03	H12	PCR		
QG073-04	A01	HQP017469	NM_022444	SLC13A1
QG073-04	A02	HQP017117	NM_022735	ACBD3
QG073-04	A03	HQP016747	NM_003006	SELPLG
QG073-04	A04	HQP016746	NM_003005	SELP
QG073-04	A05	HQP016738	NM_003004	SECTM1
QG073-04	A06	HQP016644	NM_002991	CCL24
QG073-04	A07	HQP016639	NM_004591	CCL20
QG073-04	A08	HQP016637	NM_002988	CCL18
QG073-04	A09	HQP016630	NM_005408	CCL13
QG073-04	A10	HQP016629	NM_002986	CCL11
QG073-04	A11	HQP016620	NM_002981	CCL1
QG073-04	A12	HQP016581	NM_006919	SERPINB3
QG073-04	B01	HQP016344	NM_001024808	BCL7A
QG073-04	B02	HQP016208	NM_002908	REL

Product Data Sheet

QG073-04	B03	HQP016131	NM_000321	RB1
QG073-04	B04	HQP016125	NM_002890	RASA1
QG073-04	B05	HQP016076	NM_002874	RAD23B
QG073-04	B06	HQP015828	NM_002827	PTPN1
QG073-04	B07	HQP015625	NM_020706	SFRS15
QG073-04	B08	HQP015596	NM_000962	PTGS1
QG073-04	B09	HQP015565	NM_020661	AICDA
QG073-04	B10	HQP015537	NM_000953	PTGDR
QG073-04	B11	HQP015248	NM_001080452	GPR108
QG073-04	B12	HQP015240	NM_020162	DHX33
QG073-04	C01	HQP015102	NM_001012965	KLK6
QG073-04	C02	HQP015096	NM_005046	KLK7
QG073-04	C03	HQP015061	NM_019619	PARD3
QG073-04	C04	HQP014732	NM_018416	FOXJ2
QG073-04	C05	HQP014462	NM_017509	KLK15
QG073-04	C06	HQP013991	NM_017944	USP47
QG073-04	C07	HQP013957	NM_001611	ACP5
QG073-04	C08	HQP013696	NM_017628	KIAA1546
QG073-04	C09	HQP013633	NM_005037	PPARG
QG073-04	C10	HQP013496	NM_019009	TOLLIP
QG073-04	C11	HQP013467	NM_000939	POMC
QG073-04	C12	HQP013422	NM_002690	POLB
QG073-04	D01	HQP013152	NM_002648	PIM1
QG073-04	D02	HQP013150	NM_006218	PIK3CA
QG073-04	D03	HQP013133	NM_005025	SERPINI1
QG073-04	D04	HQP013099	NM_000926	PGR
QG073-04	D05	HQP013071	NM_006212	PFKFB2
QG073-04	D06	HQP012441	NM_016123	IRAK4
QG073-04	D07	HQP012436	NM_001040443	PHF11
QG073-04	D08	HQP012212	NM_016734	PAX5
QG073-04	D09	HQP012211	NM_006193	PAX4
QG073-04	D10	HQP012166	NM_022047	DEF6
QG073-04	D11	HQP012155	NM_002575	SERPINB2
QG073-04	D12	HQP012008	NM_000275	OCA2
QG073-04	E01	HQP011914	NM_002524	NRAS
QG073-04	E02	HQP011885	NM_002518	NPAS2
QG073-04	E03	HQP011860	NM_006169	NNMT
QG073-04	E04	HQP011728	NM_004550	NDUFS2
QG073-04	E05	HQP011663	NM_005967	NAB2
QG073-04	E06	HQP011620	NM_000488	SERPINC1
QG073-04	E07	HQP011540	NM_002451	MTAP
QG073-04	E08	HQP011325	NM_019899	ABCC1
QG073-04	E09	HQP011264	NM_002425	MMP10
QG073-04	E10	HQP011263	NM_004994	MMP9
QG073-04	E11	HQP011257	NM_002422	MMP3
QG073-04	E12	HQP011255	NM_002421	MMP1
QG073-04	F01	HQP011235	NM_000249	MLH1
QG073-04	F02	HQP011223	NM_000248	MITF

Product Data Sheet

QG073-04	F03	HQP011097	NM_005912	MC4R
QG073-04	F04	HQP011085	NM_001025081	MBP
QG073-04	F05	HQP010996	NM_022438	MAL
QG073-04	F06	HQP010932	NM_005582	CD180
QG073-04	F07	HQP010878	NM_002335	LRP5
QG073-04	F08	HQP010847	NM_000237	LPL
QG073-04	F09	HQP010660	NM_005570	LMAN1
QG073-04	F10	HQP010626	NM_000236	LIPC
QG073-04	F11	HQP010612	NM_013975	LIG3
QG073-04	F12	HQP010609	NM_000234	LIG1
QG073-04	G01	HQP010560	NM_004139	LBP
QG073-04	G02	HQP010497	NM_000426	LAMA2
QG073-04	G03	HQP010104	NM_000892	KLKB1
QG073-04	G04	HQP010100	NM_002257	KLK1
QG073-04	G05	HQP009849	NM_002227	JAK1
QG073-04	G06	HQP009837	NM_033453	ITPA
QG073-04	G07	HQP009801	NM_000044	AR
QG073-04	G08	HQP009776	NM_001570	IRAK2
QG073-04	G09	HQP009745	NM_005538	INHBC
QG073-04	G10	HQP009718	NM_001562	IL18
QG073-04	G11	HQP009710	NM_002189	IL15RA
QG073-04	G12	HQP009696	NM_001559	IL12RB2
QG073-04	H01	HGDC		
QG073-04	H02	HGDC		
QG073-04	H03	HQP006940	NM_002046	GAPDH
QG073-04	H04	HQP016381	NM_001101	ACTB
QG073-04	H05	HQP015171	NM_004048	B2M
QG073-04	H06	HQP006171	NM_012423	RPL13A
QG073-04	H07	HQP009026	NM_000194	HPRT1
QG073-04	H08	HQP054253	NR_003286	RN18S1
QG073-04	H09	RT		
QG073-04	H10	RT		
QG073-04	H11	PCR		
QG073-04	H12	PCR		
QG073-05	A01	HQP009694	NM_005535	IL12RB1
QG073-05	A02	HQP009679	NM_000634	IL8RA
QG073-05	A03	HQP009650	NM_000417	IL2RA
QG073-05	A04	HQP009639	NM_001556	IKBKB
QG073-05	A05	HQP009544	NM_000598	IGFBP3
QG073-05	A06	HQP009539	NM_000596	IGFBP1
QG073-05	A07	HQP009529	NM_000612	IGF2
QG073-05	A08	HQP009523	NM_000875	IGF1R
QG073-05	A09	HQP009518	NM_000618	IGF1
QG073-05	A10	HQP009469	NM_000416	IFNGR1
QG073-05	A11	HQP009351	NM_005896	IDH1
QG073-05	A12	HQP009218	NM_000384	APOB
QG073-05	B01	HQP009201	NM_001039132	ICAM4
QG073-05	B02	HQP009171	NM_000482	APOA4

Product Data Sheet

QG073-05	B03	HQP009085	NM_002155	HSPA6
QG073-05	B04	HQP009066	NM_002153	HSD17B2
QG073-05	B05	HQP009064	NM_000413	HSD17B1
QG073-05	B06	HQP009061	NM_001641	APEX1
QG073-05	B07	HQP009060	NM_000198	HSD3B2
QG073-05	B08	HQP009059	NM_000862	HSD3B1
QG073-05	B09	HQP009019	NM_005143	HP
QG073-05	B10	HQP008893	NM_005518	HMGCS2
QG073-05	B11	HQP008892	NM_002130	HMGCS1
QG073-05	B12	HQP008846	NM_001607	ACAA1
QG073-05	C01	HQP008808	NM_021155	CD209
QG073-05	C02	HQP008801	NM_001010931	HGF
QG073-05	C03	HQP008620	NM_013371	IL19
QG073-05	C04	HQP008554	NM_012092	ICOS
QG073-05	C05	HQP008498	NM_005513	GTF2E1
QG073-05	C06	HQP008302	NM_000827	GRIA1
QG073-05	C07	HQP008282	NM_002084	GPX3
QG073-05	C08	HQP008281	NM_002083	GPX2
QG073-05	C09	HQP007821	NM_019844	SLCO1B3
QG073-05	C10	HQP007722	NM_014905	GLS
QG073-05	C11	HQP007385	NM_000515	GH1
QG073-05	C12	HQP007100	NM_015670	SENP3
QG073-05	D01	HQP006427	NM_001039130	ALOX15B
QG073-05	D02	HQP006425	NM_001140	ALOX15
QG073-05	D03	HQP006396	NM_153289	DEFB119
QG073-05	D04	HQP006375	NM_001629	ALOX5AP
QG073-05	D05	HQP006359	NM_000698	ALOX5
QG073-05	D06	HQP006356	NM_000697	ALOX12
QG073-05	D07	HQP006350	NM_015367	BCL2L13
QG073-05	D08	HQP006296	NM_015364	LY96
QG073-05	D09	HQP006245	NM_014317	PDSS1
QG073-05	D10	HQP006233	NM_012114	CASP14
QG073-05	D11	HQP006202	NM_012276	LILRA4
QG073-05	D12	HQP006131	NM_014294	TRAM1
QG073-05	E01	HQP006080	NM_012238	SIRT1
QG073-05	E02	HQP005879	NM_002019	FLT1
QG073-05	E03	HQP005403	NM_002006	FGF2
QG073-05	E04	HQP005369	NM_001010873	BZRPL1
QG073-05	E05	HQP005357	NM_004462	FDFT1
QG073-05	E06	HQP005292	NM_001002275	FCGR2B
QG073-05	E07	HQP005240	NM_004106	FCER1G
QG073-05	E08	HQP005239	NM_000139	MS4A2
QG073-05	E09	HQP005238	NM_002001	FCER1A
QG073-05	E10	HQP005022	NM_001987	ETV6
QG073-05	E11	HQP005002	NM_001437	ESR2
QG073-05	E12	HQP004991	NM_001014431	AKT1
QG073-05	F01	HQP004984	NM_005236	ERCC4
QG073-05	F02	HQP004983	NM_000122	ERCC3

Product Data Sheet

QG073-05	F03	HQP004950	NM_001979	EPHX2
QG073-05	F04	HQP004724	NM_001623	AIF1
QG073-05	F05	HQP004451	NM_000798	DRD5
QG073-05	F06	HQP004446	NM_000795	DRD2
QG073-05	F07	HQP004242	NM_194320	ZNF169
QG073-05	F08	HQP004198	NM_001928	CFD
QG073-05	F09	HQP004186	NM_021010	DEFA5
QG073-05	F10	HQP004182	NM_001925	DEFA4
QG073-05	F11	HQP004081	NM_000789	ACE
QG073-05	F12	HQP004066	NM_000788	DCK
QG073-05	G01	HQP003992	NM_001350	DAXX
QG073-05	G02	HQP003904	NM_000103	CYP19A1
QG073-05	G03	HQP003812	NM_000025	ADRB3
QG073-05	G04	HQP003809	NM_000769	CYP2C19
QG073-05	G05	HQP003539	NM_001904	CTNNB1
QG073-05	G06	HQP003420	NM_182919	TICAM1
QG073-05	G07	HQP003393	NM_144685	HIPK4
QG073-05	G08	HQP003173	NM_000759	CSF3
QG073-05	G09	HQP003084	NM_021117	CRY2
QG073-05	G10	HQP003069	NM_139074	DEFB127
QG073-05	G11	HQP002881	NM_000651	CR1
QG073-05	G12	HQP002873	NM_000098	CPT2
QG073-05	H01	HGDC		
QG073-05	H02	HGDC		
QG073-05	H03	HQP006940	NM_002046	GAPDH
QG073-05	H04	HQP016381	NM_001101	ACTB
QG073-05	H05	HQP015171	NM_004048	B2M
QG073-05	H06	HQP006171	NM_012423	RPL13A
QG073-05	H07	HQP009026	NM_000194	HPRT1
QG073-05	H08	HQP054253	NR_003286	RN18S1
QG073-05	H09	RT		
QG073-05	H10	RT		
QG073-05	H11	PCR		
QG073-05	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Lymphoma 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