

ExProfile™ Human Leukemia Gene qPCR Array

For focused group profiling of human leukemia related genes expression

Cat. No. QG070-A (6 x 96-well plate, Format A)

Cat. No. QG070-B (6 x 96-well plate, Format B)

Cat. No. QG070-C (6 x 96-well plate, Format C)

Cat. No. QG070-D (6 x 96-well plate, Format D)

Cat. No. QG070-E (6 x 96-well plate, Format E)

Available as 1 set or 6 sets. Each set contains 504 unique gene primers deposited in six 96-well plates.

Introduction

The ExProfile human leukemia gene qPCR array profiles 504 human genes to aberrantly expressed human genes involved in human leukemia. 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 leukemia cancer gene array allows researchers to study the cancer-related genes to gain understanding of their roles of genes in leukemia cancer pathogenesis.

- QG070 plate 01: 84 unique gene PCR primer pairs
- QG070 plate 02: 84 unique gene PCR primer pairs
- QG070 plate 03: 84 unique gene PCR primer pairs
- QG070 plate 04: 84 unique gene PCR primer pairs
- QG070 plate 05: 84 unique gene PCR primer pairs
- QG070 plate 06: 84 unique gene PCR primer pairs

Shipping and storage condition

Shipped at room temperature

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

Array format

GeneCopeia 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 gene 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	FLT3	MTHFR	HLA-A	HLA-DRB1	TP53	CYP1A1	CEBPA	NQO1	JAK2	ABCB1	XRCC1	WT1
B	CYP3A5	NRAS	MDM2	ERCC2	IDH1	KIT	TPMT	TYMS	TNF	MLL	CYP2E1	CYP2D6
C	KIAA1546	IDH2	RUNX1	XRCC3	BCL2	P2RX7	KRAS	IRF4	IL10	HFE	MLH1	NAT2
D	SLC22A1	NBN	MPO	KIR3DL1	KIR2DL1	ERCC5	VEGFA	NOD2	CCND1	MTRR	KIR2DL3	JAK3
E	IL1A	ERCC1	ASXL1	CYP2C9	CYP2C19	NAT1	CD5	CASP3	CASP1	CASP8	XRCC2	TNFRSF1B
F	KIR2DS3	KIR2DL4	IL4	IL1B	BIRC5	APEX1	FCGR2A	KIR3DL3	IKZF1	KIR3DL2	CD38	TNFRSF10A
G	ARID5B	CASP7	XRCC4	XPC	VDR	VCAM1	TYK2	BRCA2	SOD2	BCR	BCL2L2	MSH3
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG070 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	MSH2	JAK1	IL8	FASLG	IL1RN	APAF1	IL19	GNB3	FMO2	ERCC6	ERCC4	EPHX1
B	CYP2B6	SP140	CDKN2A	ULK3	CD28	SLC28A1	CCNH	TNFSF18	TNFRSF10B	FADD	CBL	CBFB
C	CASP5	TLR10	CXCR4	XRCC5	TLR1	TGFB1	STAT3	STAT1	SPI1	SELPLG	SELP	CXCL12
D	CCL2	RASA1	RAD23B	POLB	PIM1	PDGFRB	IRAK4	ORM1	NOS2A	NFKB1	NF1	MTR
E	MPL	LIG4	LIG1	KDR	IL18	IL13	IL12RB1	IL12A	IL8RB	IL6ST	IL6	IL5RA
F	IL1R1	APOE	IFNGR2	IFNGR1	SP110	HSPA6	GZMB	STRN4	GSTM3	NR3C1	ABO	GNAS
G	LILRA4	F5	ETV6	AKT1	CYP21A2	CYP1B1	CTLA4	CSF3	CSF2	PARP1	CCR8	CCR7
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG070 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	TOPBP1	NOD1	TLR6	RAD50	TNFSF13	CEBPE	RYR2	ZNF527	ICHTHYIN	PRF1	EXO1	CASP8AP2
B	CDC25C	CDC6	CDA	CD81	CD79B	CD79A	ESPL1	N4BP1	MDC1	CLCA2	KLK4	CD40
C	TP53I3	CD14	DLG5	AURKB	CD4	IL1RL1	SLC28A2	PDCD5	SOCS3	BCL10	STK19	CFLAR
D	SOCS2	IL1RL2	TNFRSF11A	RIPK2	TNFRSF14	TNFSF10	RIPK1	ADAM19	TRADD	ABCC3	SOCS1	AKR1C3
E	MADD	TSLP	CNTNAP1	FOXN1	RAD54L	ACAD11	CASP6	CASP4	CAPN3	CAPN2	DEFB126	COL18A1
F	RMI1	C7orf24	DDR1	ZNF230	ZBTB16	XDH	SCGB1A1	C8G	C8B	C7	TYR	C5AR1
G	C4BPB	C4BPA	TRAF6	TRAF5	TRAF3	TRAF2	C2	TP53BP1	TOP2A	C1QC	TLR3	C1QB
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG070 plate 03

Product Data Sheet

	1	2	3	4	5	6	7	8	9	10	11	12
A	BUB1B	C2orf3	TAL1	SUOX	SULT1A1	STAT4	STAT2	BPI	SOS1	SOD1	SLC2A11	SLC19A1
B	CPEB1	DCLRE1C	BLM	SLC28A3	SELE	SECTM1	CCL24	CCL20	CCL18	CCL13	CCL11	CCL5
C	CCL1	SERPINB3	CFB	SERPING1	RPA3	RPA1	BCL7A	REV3L	RELB	RELA	REL	RBL2
D	RAP1GAP	RAC1	RAD1	BAK1	SFRS15	AICDA	PTGDR	PTCH1	GPR108	DHX33	KLK6	KLK7
E	PARD3	ANKH	FOXJ2	KLK15	FBXW7	USP47	ATR	TOLLIP	HAO1	POLE	TLR9	SERPINI1
F	PFKFB2	PECAM1	REV1	TNFRSF12A	C1RL	SH3GLB1	GMNN	DEF6	SERPINB2	SERPINE1	ORM2	ORC4L
G	OPRM1	NTRK1	NOS3	NQO2	NFKBIE	NFKBIA	NFKB2	NEDD4	NDUFS2	NCF4	NCF2	NAB2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG070 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	SERPINC1	MYH11	MYD88	MYC	ABCC1	MRE11A	MMP14	MMP10	MMP9	MMP8	MIF	CD46
B	MCM5	MCM4	MC4R	MBP	MBD1	MAPT	MAN2A1	MAL	SMAD7	CD180	LMAN1	LIG3
C	LEPR	LBP	LAMA2	KLKB1	KLK1	ITGB4	ITGAE	IRF5	IRAK2	INHBC	TNFRSF9	IL15
D	IL12RB2	IL12B	IL10RB	IL8RA	IL6R	IL5	IL4R	IL2	IKKBK	IFNAR2	IFNAR1	IRF8
E	ICAM4	NLRP14	APOA2	BIRC4	HOXD4	ACAA1	CD209	HGF	CXXC1	HIST1H1T	ICOS	GTF2E1
F	GSTZ1	GSTA2	GPX4	GPX3	GPX2	ATG9B	SLCO1B3	GLS	POLL	GATA2	SENP3	XRCC6
G	ALOX15B	ALOX15	DEFB119	ALOX5AP	ALOX5	ALOX12	KCNE4	LY96	ORC3L	CASP14	TRAM1	FMO3
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG070 plate 05

	1	2	3	4	5	6	7	8	9	10	11	12
A	CD93	NLRP1	FGF4	FGF2	FES	FCER2	FCER1G	FCER1A	F2	ESR1	ERCC3	EPHX2
B	NLRP11	ENO2	EGFR	EGF	EDN1	E2F2	DOK1	DNMT3A	DHFR	ZNF169	DEFA5	DEFA4
C	DDC	ACE	DCK	CYP4B1	ESCO2	CYP2C18	CYP2C8	CTNNA1	TICAM1	NLRP4	HIPK4	CSF1
D	DEFB127	CPT2	CPT1B	NLRP8	CCR4	CCR1	SOCS4	DEFB118	ERCC8	CHUK	TWF2	GPR44
E	IRAK3	POLI	CHEK2	CHIT1	WDHD1	NUDT6	FAF1	DSCR1L2	CCL27	CAPN9	ARL6IP5	SEC23B
F	CCL26	CDK6	CDK4	TANK	MUC6	BAT5	TP73	C1QB	SOD3	XCL1	C12orf5	MASP1
G	CXXC5	NOS1	MUC7	LPO	KLK2	ICAM3	ZP1	DNMT3B	DEFA6	NLRP5	NLRP13	CLCA1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure6. Illustration of QG070 plate 06

- **Gene primer pairs:** 84 wells from A to G in each 96-well plate 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
QG070-01	A01	HQP005890	NM_004119	FLT3
QG070-01	A02	HQP011547	NM_005957	MTHFR
QG070-01	A03	HQP008849	NM_002116	HLA-A
QG070-01	A04	HQP054047	BC008403	HLA-DRB1
QG070-01	A05	HQP018175	NM_000546	TP53

Product Data Sheet

QG070-01	A06	HQP003772	NM_000499	CYP1A1
QG070-01	A07	HQP000615	NM_004364	CEBPA
QG070-01	A08	HQP004317	NM_000903	NQO1
QG070-01	A09	HQP009850	NM_004972	JAK2
QG070-01	A10	HQP013100	NM_000927	ABCB1
QG070-01	A11	HQP018562	NM_006297	XRCC1
QG070-01	A12	HQP018546	NM_000378	WT1
QG070-01	B01	HQP003841	NM_000777	CYP3A5
QG070-01	B02	HQP011914	NM_002524	NRAS
QG070-01	B03	HQP011135	NM_002392	MDM2
QG070-01	B04	HQP004976	NM_000400	ERCC2
QG070-01	B05	HQP009351	NM_005896	IDH1
QG070-01	B06	HQP010099	NM_000222	KIT
QG070-01	B07	HQP018213	NM_000367	TPMT
QG070-01	B08	HQP018342	NM_001071	TYMS
QG070-01	B09	HQP018141	NM_000594	TNF
QG070-01	B10	HQP011241	NM_005933	MLL
QG070-01	B11	HQP003817	NM_000773	CYP2E1
QG070-01	B12	HQP003814	NM_000106	CYP2D6
QG070-01	C01	HQP013696	NM_017628	KIAA1546
QG070-01	C02	HQP009353	NM_002168	IDH2
QG070-01	C03	HQP021347	NM_001754	RUNX1
QG070-01	C04	HQP018564	NM_005432	XRCC3
QG070-01	C05	HQP016212	NM_000657	BCL2
QG070-01	C06	HQP012100	NM_002562	P2RX7
QG070-01	C07	HQP010133	NM_004985	KRAS
QG070-01	C08	HQP009781	NM_002460	IRF4
QG070-01	C09	HQP009685	NM_000572	IL10
QG070-01	C10	HQP008757	NM_000410	HFE
QG070-01	C11	HQP011235	NM_000249	MLH1
QG070-01	C12	HQP001136	NM_000015	NAT2
QG070-01	D01	HQP017490	NM_003057	SLC22A1
QG070-01	D02	HQP011687	NM_002485	NBN
QG070-01	D03	HQP011309	NM_000250	MPO
QG070-01	D04	HQP010095	NM_013289	KIR3DL1
QG070-01	D05	HQP010083	NM_014218	KIR2DL1
QG070-01	D06	HQP004985	NM_000123	ERCC5
QG070-01	D07	HQP018475	NM_001025366	VEGFA
QG070-01	D08	HQP016801	NM_022162	NOD2
QG070-01	D09	HQP016204	NM_053056	CCND1
QG070-01	D10	HQP011555	NM_002454	MTRR
QG070-01	D11	HQP010086	NM_015868	KIR2DL3
QG070-01	D12	HQP009851	NM_000215	JAK3
QG070-01	E01	HQP009640	NM_000575	IL1A
QG070-01	E02	HQP004975	NM_202001	ERCC1
QG070-01	E03	HQP004287	NM_015338	ASXL1
QG070-01	E04	HQP003811	NM_000771	CYP2C9
QG070-01	E05	HQP003809	NM_000769	CYP2C19

Product Data Sheet

QG070-01	E06	HQP023467	NM_000662	NAT1
QG070-01	E07	HQP022337	NM_014207	CD5
QG070-01	E08	HQP020297	NM_004346	CASP3
QG070-01	E09	HQP020207	NM_001223	CASP1
QG070-01	E10	HQP018966	NM_001080124	CASP8
QG070-01	E11	HQP018563	NM_005431	XRCC2
QG070-01	E12	HQP018149	NM_001066	TNFRSF1B
QG070-01	F01	HQP010092	NM_012313	KIR2DS3
QG070-01	F02	HQP010089	NM_002255	KIR2DL4
QG070-01	F03	HQP009662	NM_000589	IL4
QG070-01	F04	HQP009641	NM_000576	IL1B
QG070-01	F05	HQP009101	NM_001168	BIRC5
QG070-01	F06	HQP009061	NM_001641	APEX1
QG070-01	F07	HQP005280	NM_021642	FCGR2A
QG070-01	F08	HQP001810	NM_153443	KIR3DL3
QG070-01	F09	HQP000397	NM_006060	IKZF1
QG070-01	F10	HQP010096	NM_006737	KIR3DL2
QG070-01	F11	HQP022870	NM_001775	CD38
QG070-01	F12	HQP021557	NM_003844	TNFRSF10A
QG070-01	G01	HQP020515	NM_032199	ARID5B
QG070-01	G02	HQP020481	NM_033338	CASP7
QG070-01	G03	HQP018565	NM_003401	XRCC4
QG070-01	G04	HQP018556	NM_004628	XPC
QG070-01	G05	HQP018474	NM_000376	VDR
QG070-01	G06	HQP018466	NM_080682	VCAM1
QG070-01	G07	HQP018340	NM_003331	TYK2
QG070-01	G08	HQP017753	NM_000059	BRCA2
QG070-01	G09	HQP017616	NM_000636	SOD2
QG070-01	G10	HQP016418	NM_004327	BCR
QG070-01	G11	HQP016260	NM_004050	BCL2L2
QG070-01	G12	HQP011492	NM_002439	MSH3
QG070-01	H01	HGDC		
QG070-01	H02	HGDC		
QG070-01	H03	HQP006940	NM_002046	GAPDH
QG070-01	H04	HQP016381	NM_001101	ACTB
QG070-01	H05	HQP015171	NM_004048	B2M
QG070-01	H06	HQP006171	NM_012423	RPL13A
QG070-01	H07	HQP009026	NM_000194	HPRT1
QG070-01	H08	HQP054253	NR_003286	RN18S1
QG070-01	H09	RT		
QG070-01	H10	RT		
QG070-01	H11	PCR		
QG070-01	H12	PCR		
QG070-02	A01	HQP011491	NM_000251	MSH2
QG070-02	A02	HQP009849	NM_002227	JAK1
QG070-02	A03	HQP009678	NM_000584	IL8
QG070-02	A04	HQP009671	NM_000639	FASLG
QG070-02	A05	HQP009645	NM_000577	IL1RN

Product Data Sheet

QG070-02	A06	HQP008934	NM_001160	APAF1
QG070-02	A07	HQP008620	NM_013371	IL19
QG070-02	A08	HQP007767	NM_002075	GNB3
QG070-02	A09	HQP005951	NM_001460	FMO2
QG070-02	A10	HQP004986	NM_000124	ERCC6
QG070-02	A11	HQP004984	NM_005236	ERCC4
QG070-02	A12	HQP004948	NM_000120	EPHX1
QG070-02	B01	HQP003808	NM_000767	CYP2B6
QG070-02	B02	HQP001475	NM_001005176	SP140
QG070-02	B03	HQP000370	NM_058195	CDKN2A
QG070-02	B04	HQP054690	NM_001099436	ULK3
QG070-02	B05	HQP022699	NM_006139	CD28
QG070-02	B06	HQP022202	NM_004213	SLC28A1
QG070-02	B07	HQP021906	NM_001239	CCNH
QG070-02	B08	HQP021834	NM_005092	TNFSF18
QG070-02	B09	HQP021553	NM_003842	TNFRSF10B
QG070-02	B10	HQP021526	NM_003824	FADD
QG070-02	B11	HQP021430	NM_005188	CBL
QG070-02	B12	HQP021408	NM_022845	CBFB
QG070-02	C01	HQP020384	NM_004347	CASP5
QG070-02	C02	HQP019982	NM_001017388	TLR10
QG070-02	C03	HQP018802	NM_001008540	CXCR4
QG070-02	C04	HQP018568	NM_021141	XRCC5
QG070-02	C05	HQP018113	NM_003263	TLR1
QG070-02	C06	HQP018044	NM_000660	TGFB1
QG070-02	C07	HQP017767	NM_003150	STAT3
QG070-02	C08	HQP017764	NM_007315	STAT1
QG070-02	C09	HQP017663	NM_003120	SPI1
QG070-02	C10	HQP016747	NM_003006	SELPLG
QG070-02	C11	HQP016746	NM_003005	SELP
QG070-02	C12	HQP016670	NM_001033886	CXCL12
QG070-02	D01	HQP016621	NM_002982	CCL2
QG070-02	D02	HQP016125	NM_002890	RASA1
QG070-02	D03	HQP016076	NM_002874	RAD23B
QG070-02	D04	HQP013422	NM_002690	POLB
QG070-02	D05	HQP013152	NM_002648	PIM1
QG070-02	D06	HQP012889	NM_002609	PDGFRB
QG070-02	D07	HQP012441	NM_016123	IRAK4
QG070-02	D08	HQP012078	NM_000607	ORM1
QG070-02	D09	HQP011866	NM_000625	NOS2A
QG070-02	D10	HQP011807	NM_003998	NFKB1
QG070-02	D11	HQP011774	NM_000267	NF1
QG070-02	D12	HQP011554	NM_000254	MTR
QG070-02	E01	HQP011308	NM_005373	MPL
QG070-02	E02	HQP010613	NM_002312	LIG4
QG070-02	E03	HQP010609	NM_000234	LIG1
QG070-02	E04	HQP010070	NM_002253	KDR
QG070-02	E05	HQP009718	NM_001562	IL18

Product Data Sheet

QG070-02	E06	HQP009697	NM_002188	IL13
QG070-02	E07	HQP009694	NM_005535	IL12RB1
QG070-02	E08	HQP009692	NM_000882	IL12A
QG070-02	E09	HQP009681	NM_001557	IL8RB
QG070-02	E10	HQP009675	NM_175767	IL6ST
QG070-02	E11	HQP009670	NM_000600	IL6
QG070-02	E12	HQP009667	NM_000564	IL5RA
QG070-02	F01	HQP009642	NM_000877	IL1R1
QG070-02	F02	HQP009556	NM_000041	APOE
QG070-02	F03	HQP009472	NM_005534	IFNGR2
QG070-02	F04	HQP009469	NM_000416	IFNGR1
QG070-02	F05	HQP009399	NM_004509	SP110
QG070-02	F06	HQP009085	NM_002155	HSPA6
QG070-02	F07	HQP008690	NM_004131	GZMB
QG070-02	F08	HQP008565	NM_001039877	STRN4
QG070-02	F09	HQP008483	NM_000849	GSTM3
QG070-02	F10	HQP008402	NM_001020825	NR3C1
QG070-02	F11	HQP008362	NM_020469	ABO
QG070-02	F12	HQP007755	NM_000516	GNAS
QG070-02	G01	HQP006202	NM_012276	LILRA4
QG070-02	G02	HQP005058	NM_000130	F5
QG070-02	G03	HQP005022	NM_001987	ETV6
QG070-02	G04	HQP004991	NM_001014431	AKT1
QG070-02	G05	HQP003908	NM_000500	CYP21A2
QG070-02	G06	HQP003775	NM_000104	CYP1B1
QG070-02	G07	HQP003499	NM_001037631	CTLA4
QG070-02	G08	HQP003173	NM_000759	CSF3
QG070-02	G09	HQP003159	NM_000758	CSF2
QG070-02	G10	HQP003120	NM_001618	PARP1
QG070-02	G11	HQP002222	NM_005201	CCR8
QG070-02	G12	HQP002217	NM_001838	CCR7
QG070-02	H01	HGDC		
QG070-02	H02	HGDC		
QG070-02	H03	HQP006940	NM_002046	GAPDH
QG070-02	H04	HQP016381	NM_001101	ACTB
QG070-02	H05	HQP015171	NM_004048	B2M
QG070-02	H06	HQP006171	NM_012423	RPL13A
QG070-02	H07	HQP009026	NM_000194	HPRT1
QG070-02	H08	HQP054253	NR_003286	RN18S1
QG070-02	H09	RT		
QG070-02	H10	RT		
QG070-02	H11	PCR		
QG070-02	H12	PCR		
QG070-03	A01	HQP001228	NM_007027	TOPBP1
QG070-03	A02	HQP000459	NM_006092	NOD1
QG070-03	A03	HQP000415	NM_006068	TLR6
QG070-03	A04	HQP000145	NM_005732	RAD50
QG070-03	A05	HQP021498	NM_003808	TNFSF13

Product Data Sheet

QG070-03	A06	HQP000648	NM_001805	CEBPE
QG070-03	A07	HQP054701	NM_001035	RYR2
QG070-03	A08	HQP054688	NM_032453	ZNF527
QG070-03	A09	HQP054326	NM_001099287	ICHTHYIN
QG070-03	A10	HQP054058	NM_005041	PRF1
QG070-03	A11	HQP054030	NM_130398	EXO1
QG070-03	A12	HQP023464	NM_012115	CASP8AP2
QG070-03	B01	HQP023409	NM_001790	CDC25C
QG070-03	B02	HQP023354	NM_001254	CDC6
QG070-03	B03	HQP023203	NM_001785	CDA
QG070-03	B04	HQP023168	NM_004356	CD81
QG070-03	B05	HQP023155	NM_000626	CD79B
QG070-03	B06	HQP023145	NM_021601	CD79A
QG070-03	B07	HQP023095	NM_012291	ESPL1
QG070-03	B08	HQP023074	NM_153029	N4BP1
QG070-03	B09	HQP023038	NM_014641	MDC1
QG070-03	B10	HQP023012	NM_006536	CLCA2
QG070-03	B11	HQP022998	NM_004917	KLK4
QG070-03	B12	HQP022955	NM_001250	CD40
QG070-03	C01	HQP022884	NM_004881	TP53I3
QG070-03	C02	HQP022490	NM_000591	CD14
QG070-03	C03	HQP022361	NM_004747	DLG5
QG070-03	C04	HQP022323	NM_004217	AURKB
QG070-03	C05	HQP022316	NM_000616	CD4
QG070-03	C06	HQP022241	NM_003856	IL1RL1
QG070-03	C07	HQP022197	NM_004212	SLC28A2
QG070-03	C08	HQP022170	NM_004708	PDCD5
QG070-03	C09	HQP021889	NM_003955	SOCS3
QG070-03	C10	HQP021725	NM_003921	BCL10
QG070-03	C11	HQP021636	NM_032454	STK19
QG070-03	C12	HQP021604	NM_003879	CFLAR
QG070-03	D01	HQP021602	NM_003877	SOCS2
QG070-03	D02	HQP021569	NM_003854	IL1RL2
QG070-03	D03	HQP021550	NM_003839	TNFRSF11A
QG070-03	D04	HQP021524	NM_003821	RIPK2
QG070-03	D05	HQP021522	NM_003820	TNFRSF14
QG070-03	D06	HQP021502	NM_003810	TNFSF10
QG070-03	D07	HQP021492	NM_003804	RIPK1
QG070-03	D08	HQP021484	NM_033274	ADAM19
QG070-03	D09	HQP021462	NM_003789	TRADD
QG070-03	D10	HQP021459	NM_003786	ABCC3
QG070-03	D11	HQP021399	NM_003745	SOCS1
QG070-03	D12	HQP021385	NM_003739	AKR1C3
QG070-03	E01	HQP021290	NM_003682	MADD
QG070-03	E02	HQP021260	NM_033035	TSLP
QG070-03	E03	HQP021130	NM_003632	CNTNAP1
QG070-03	E04	HQP020808	NM_003593	FOXN1
QG070-03	E05	HQP020707	NM_003579	RAD54L

Product Data Sheet

QG070-03	E06	HQP020499	NM_032169	ACAD11
QG070-03	E07	HQP020427	NM_001226	CASP6
QG070-03	E08	HQP020332	NM_001225	CASP4
QG070-03	E09	HQP020071	NM_000070	CAPN3
QG070-03	E10	HQP020070	NM_001748	CAPN2
QG070-03	E11	HQP019941	NM_030931	DEFB126
QG070-03	E12	HQP019768	NM_130445	COL18A1
QG070-03	F01	HQP019462	NM_024945	RMI1
QG070-03	F02	HQP018854	NM_024051	C7orf24
QG070-03	F03	HQP018768	NM_001954	DDR1
QG070-03	F04	HQP018739	NM_006300	ZNF230
QG070-03	F05	HQP018682	NM_001018011	ZBTB16
QG070-03	F06	HQP018552	NM_000379	XDH
QG070-03	F07	HQP018411	NM_003357	SCGB1A1
QG070-03	F08	HQP018391	NM_000606	C8G
QG070-03	F09	HQP018380	NM_000066	C8B
QG070-03	F10	HQP018353	NM_000587	C7
QG070-03	F11	HQP018343	NM_000372	TYR
QG070-03	F12	HQP018326	NM_001736	C5AR1
QG070-03	G01	HQP018276	NM_000716	C4BPB
QG070-03	G02	HQP018259	NM_000715	C4BPA
QG070-03	G03	HQP018237	NM_004620	TRAF6
QG070-03	G04	HQP018236	NM_001033910	TRAF5
QG070-03	G05	HQP018234	NM_003300	TRAF3
QG070-03	G06	HQP018233	NM_021138	TRAF2
QG070-03	G07	HQP018226	NM_000063	C2
QG070-03	G08	HQP018176	NM_005657	TP53BP1
QG070-03	G09	HQP018172	NM_001067	TOP2A
QG070-03	G10	HQP018170	NM_172369	C1QC
QG070-03	G11	HQP018115	NM_003265	TLR3
QG070-03	G12	HQP018106	NM_001212	C1QBP
QG070-03	H01	HGDC		
QG070-03	H02	HGDC		
QG070-03	H03	HQP006940	NM_002046	GAPDH
QG070-03	H04	HQP016381	NM_001101	ACTB
QG070-03	H05	HQP015171	NM_004048	B2M
QG070-03	H06	HQP006171	NM_012423	RPL13A
QG070-03	H07	HQP009026	NM_000194	HPRT1
QG070-03	H08	HQP054253	NR_003286	RN18S1
QG070-03	H09	RT		
QG070-03	H10	RT		
QG070-03	H11	PCR		
QG070-03	H12	PCR		
QG070-04	A01	HQP018022	NM_001211	BUB1B
QG070-04	A02	HQP017964	NM_003203	C2orf3
QG070-04	A03	HQP017895	NM_003189	TAL1
QG070-04	A04	HQP017816	NM_000456	SUOX
QG070-04	A05	HQP017811	NM_177536	SULT1A1

Product Data Sheet

QG070-04	A06	HQP017770	NM_003151	STAT4
QG070-04	A07	HQP017766	NM_005419	STAT2
QG070-04	A08	HQP017702	NM_001725	BPI
QG070-04	A09	HQP017625	NM_005633	SOS1
QG070-04	A10	HQP017615	NM_000454	SOD1
QG070-04	A11	HQP017549	NM_030807	SLC2A11
QG070-04	A12	HQP017482	NM_194255	SLC19A1
QG070-04	B01	HQP017022	NM_001079533	CPEB1
QG070-04	B02	HQP016982	NM_001033855	DCLRE1C
QG070-04	B03	HQP016844	NM_000057	BLM
QG070-04	B04	HQP016763	NM_022127	SLC28A3
QG070-04	B05	HQP016744	NM_000450	SELE
QG070-04	B06	HQP016738	NM_003004	SECTM1
QG070-04	B07	HQP016644	NM_002991	CCL24
QG070-04	B08	HQP016639	NM_004591	CCL20
QG070-04	B09	HQP016637	NM_002988	CCL18
QG070-04	B10	HQP016630	NM_005408	CCL13
QG070-04	B11	HQP016629	NM_002986	CCL11
QG070-04	B12	HQP016626	NM_002985	CCL5
QG070-04	C01	HQP016620	NM_002981	CCL1
QG070-04	C02	HQP016581	NM_006919	SERPINB3
QG070-04	C03	HQP016565	NM_001710	CFB
QG070-04	C04	HQP016500	NM_001032295	SERPING1
QG070-04	C05	HQP016391	NM_002947	RPA3
QG070-04	C06	HQP016389	NM_002945	RPA1
QG070-04	C07	HQP016344	NM_001024808	BCL7A
QG070-04	C08	HQP016223	NM_002912	REV3L
QG070-04	C09	HQP016214	NM_006509	RELB
QG070-04	C10	HQP016213	NM_021975	RELA
QG070-04	C11	HQP016208	NM_002908	REL
QG070-04	C12	HQP016172	NM_005611	RBL2
QG070-04	D01	HQP016107	NM_002885	RAP1GAP
QG070-04	D02	HQP016064	NM_018890	RAC1
QG070-04	D03	HQP015947	NM_002853	RAD1
QG070-04	D04	HQP015917	NM_001188	BAK1
QG070-04	D05	HQP015625	NM_020706	SFRS15
QG070-04	D06	HQP015565	NM_020661	AICDA
QG070-04	D07	HQP015537	NM_000953	PTGDR
QG070-04	D08	HQP015530	NM_000264	PTCH1
QG070-04	D09	HQP015248	NM_001080452	GPR108
QG070-04	D10	HQP015240	NM_020162	DHX33
QG070-04	D11	HQP015102	NM_001012965	KLK6
QG070-04	D12	HQP015096	NM_005046	KLK7
QG070-04	E01	HQP015061	NM_019619	PARD3
QG070-04	E02	HQP015023	NM_054027	ANKH
QG070-04	E03	HQP014732	NM_018416	FOXJ2
QG070-04	E04	HQP014462	NM_017509	KLK15
QG070-04	E05	HQP014292	NM_018315	FBXW7

Product Data Sheet

QG070-04	E06	HQP013991	NM_017944	USP47
QG070-04	E07	HQP013596	NM_001184	ATR
QG070-04	E08	HQP013496	NM_019009	TOLLIP
QG070-04	E09	HQP013447	NM_017545	HAO1
QG070-04	E10	HQP013425	NM_006231	POLE
QG070-04	E11	HQP013388	NM_017442	TLR9
QG070-04	E12	HQP013133	NM_005025	SERPINI1
QG070-04	F01	HQP013071	NM_006212	PFKFB2
QG070-04	F02	HQP013015	NM_000442	PECAM1
QG070-04	F03	HQP012755	NM_001037872	REV1
QG070-04	F04	HQP012651	NM_016639	TNFRSF12A
QG070-04	F05	HQP012584	NM_016546	C1RL
QG070-04	F06	HQP012393	NM_016009	SH3GLB1
QG070-04	F07	HQP012335	NM_015895	GMNN
QG070-04	F08	HQP012166	NM_022047	DEF6
QG070-04	F09	HQP012155	NM_002575	SERPINB2
QG070-04	F10	HQP012154	NM_000602	SERPINE1
QG070-04	F11	HQP012079	NM_000608	ORM2
QG070-04	F12	HQP012073	NM_002552	ORC4L
QG070-04	G01	HQP012062	NM_001008505	OPRM1
QG070-04	G02	HQP011938	NM_001007792	NTRK1
QG070-04	G03	HQP011868	NM_000603	NOS3
QG070-04	G04	HQP011858	NM_000904	NQO2
QG070-04	G05	HQP011813	NM_004556	NFKBIE
QG070-04	G06	HQP011810	NM_020529	NFKBIA
QG070-04	G07	HQP011808	NM_001077493	NFKB2
QG070-04	G08	HQP011741	NM_006154	NEDD4
QG070-04	G09	HQP011728	NM_004550	NDUFS2
QG070-04	G10	HQP011694	NM_000631	NCF4
QG070-04	G11	HQP011693	NM_000433	NCF2
QG070-04	G12	HQP011663	NM_005967	NAB2
QG070-04	H01	HGDC		
QG070-04	H02	HGDC		
QG070-04	H03	HQP006940	NM_002046	GAPDH
QG070-04	H04	HQP016381	NM_001101	ACTB
QG070-04	H05	HQP015171	NM_004048	B2M
QG070-04	H06	HQP006171	NM_012423	RPL13A
QG070-04	H07	HQP009026	NM_000194	HPRT1
QG070-04	H08	HQP054253	NR_003286	RN18S1
QG070-04	H09	RT		
QG070-04	H10	RT		
QG070-04	H11	PCR		
QG070-04	H12	PCR		
QG070-05	A01	HQP011620	NM_000488	SERPINC1
QG070-05	A02	HQP011616	NM_001040113	MYH11
QG070-05	A03	HQP011603	NM_002468	MYD88
QG070-05	A04	HQP011597	NM_002467	MYC
QG070-05	A05	HQP011325	NM_019899	ABCC1

Product Data Sheet

QG070-05	A06	HQP011320	NM_005590	MRE11A
QG070-05	A07	HQP011268	NM_004995	MMP14
QG070-05	A08	HQP011264	NM_002425	MMP10
QG070-05	A09	HQP011263	NM_004994	MMP9
QG070-05	A10	HQP011262	NM_002424	MMP8
QG070-05	A11	HQP011219	NM_002415	MIF
QG070-05	A12	HQP011113	NM_002389	CD46
QG070-05	B01	HQP011109	NM_006739	MCM5
QG070-05	B02	HQP011108	NM_005914	MCM4
QG070-05	B03	HQP011097	NM_005912	MC4R
QG070-05	B04	HQP011085	NM_001025081	MBP
QG070-05	B05	HQP011073	NM_015844	MBD1
QG070-05	B06	HQP011024	NM_016835	MAPT
QG070-05	B07	HQP011004	NM_002372	MAN2A1
QG070-05	B08	HQP010996	NM_022438	MAL
QG070-05	B09	HQP010966	NM_005904	SMAD7
QG070-05	B10	HQP010932	NM_005582	CD180
QG070-05	B11	HQP010660	NM_005570	LMAN1
QG070-05	B12	HQP010612	NM_013975	LIG3
QG070-05	C01	HQP010584	NM_002303	LEPR
QG070-05	C02	HQP010560	NM_004139	LBP
QG070-05	C03	HQP010497	NM_000426	LAMA2
QG070-05	C04	HQP010104	NM_000892	KLKB1
QG070-05	C05	HQP010100	NM_002257	KLK1
QG070-05	C06	HQP009819	NM_000213	ITGB4
QG070-05	C07	HQP009805	NM_002208	ITGAE
QG070-05	C08	HQP009782	NM_002200	IRF5
QG070-05	C09	HQP009776	NM_001570	IRAK2
QG070-05	C10	HQP009745	NM_005538	INHBC
QG070-05	C11	HQP009716	NM_001561	TNFRSF9
QG070-05	C12	HQP009708	NM_000585	IL15
QG070-05	D01	HQP009696	NM_001559	IL12RB2
QG070-05	D02	HQP009693	NM_002187	IL12B
QG070-05	D03	HQP009687	NM_000628	IL10RB
QG070-05	D04	HQP009679	NM_000634	IL8RA
QG070-05	D05	HQP009672	NM_000565	IL6R
QG070-05	D06	HQP009666	NM_000879	IL5
QG070-05	D07	HQP009664	NM_000418	IL4R
QG070-05	D08	HQP009649	NM_000586	IL2
QG070-05	D09	HQP009639	NM_001556	IKKB
QG070-05	D10	HQP009460	NM_000874	IFNAR2
QG070-05	D11	HQP009458	NM_000629	IFNAR1
QG070-05	D12	HQP009251	NM_002163	IRF8
QG070-05	E01	HQP009201	NM_001039132	ICAM4
QG070-05	E02	HQP009177	NM_176822	NLRP14
QG070-05	E03	HQP009139	NM_001643	APOA2
QG070-05	E04	HQP009091	NM_001167	BIRC4
QG070-05	E05	HQP009011	NM_014621	HOXD4

Product Data Sheet

QG070-05	E06	HQP008846	NM_001607	ACAA1
QG070-05	E07	HQP008808	NM_021155	CD209
QG070-05	E08	HQP008801	NM_001010931	HGF
QG070-05	E09	HQP008799	NM_014593	CXXC1
QG070-05	E10	HQP008700	NM_005323	HIST1H1T
QG070-05	E11	HQP008554	NM_012092	ICOS
QG070-05	E12	HQP008498	NM_005513	GTF2E1
QG070-05	F01	HQP008490	NM_001513	GSTZ1
QG070-05	F02	HQP008476	NM_000846	GSTA2
QG070-05	F03	HQP008285	NM_002085	GPX4
QG070-05	F04	HQP008282	NM_002084	GPX3
QG070-05	F05	HQP008281	NM_002083	GPX2
QG070-05	F06	HQP008198	NM_173681	ATG9B
QG070-05	F07	HQP007821	NM_019844	SLCO1B3
QG070-05	F08	HQP007722	NM_014905	GLS
QG070-05	F09	HQP007689	NM_013274	POLL
QG070-05	F10	HQP007152	NM_032638	GATA2
QG070-05	F11	HQP007100	NM_015670	SENP3
QG070-05	F12	HQP006547	NM_001469	XRCC6
QG070-05	G01	HQP006427	NM_001039130	ALOX15B
QG070-05	G02	HQP006425	NM_001140	ALOX15
QG070-05	G03	HQP006396	NM_153289	DEFB119
QG070-05	G04	HQP006375	NM_001629	ALOX5AP
QG070-05	G05	HQP006359	NM_000698	ALOX5
QG070-05	G06	HQP006356	NM_000697	ALOX12
QG070-05	G07	HQP006320	NM_080671	KCNE4
QG070-05	G08	HQP006296	NM_015364	LY96
QG070-05	G09	HQP006249	NM_012381	ORC3L
QG070-05	G10	HQP006233	NM_012114	CASP14
QG070-05	G11	HQP006131	NM_014294	TRAM1
QG070-05	G12	HQP005959	NM_001002294	FMO3
QG070-05	H01	HGDC		
QG070-05	H02	HGDC		
QG070-05	H03	HQP006940	NM_002046	GAPDH
QG070-05	H04	HQP016381	NM_001101	ACTB
QG070-05	H05	HQP015171	NM_004048	B2M
QG070-05	H06	HQP006171	NM_012423	RPL13A
QG070-05	H07	HQP009026	NM_000194	HPRT1
QG070-05	H08	HQP054253	NR_003286	RN18S1
QG070-05	H09	RT		
QG070-05	H10	RT		
QG070-05	H11	PCR		
QG070-05	H12	PCR		
QG070-06	A01	HQP005589	NM_012072	CD93
QG070-06	A02	HQP005511	NM_001033053	NLRP1
QG070-06	A03	HQP005405	NM_002007	FGF4
QG070-06	A04	HQP005403	NM_002006	FGF2
QG070-06	A05	HQP005395	NM_002005	FES

Product Data Sheet

QG070-06	A06	HQP005242	NM_002002	FCER2
QG070-06	A07	HQP005240	NM_004106	FCER1G
QG070-06	A08	HQP005238	NM_002001	FCER1A
QG070-06	A09	HQP005052	NM_000506	F2
QG070-06	A10	HQP004998	NM_000125	ESR1
QG070-06	A11	HQP004983	NM_000122	ERCC3
QG070-06	A12	HQP004950	NM_001979	EPHX2
QG070-06	B01	HQP004934	NM_145007	NLRP11
QG070-06	B02	HQP004864	NM_001975	ENO2
QG070-06	B03	HQP004605	NM_005228	EGFR
QG070-06	B04	HQP004599	NM_001963	EGF
QG070-06	B05	HQP004557	NM_001955	EDN1
QG070-06	B06	HQP004526	NM_004091	E2F2
QG070-06	B07	HQP004426	NM_001381	DOK1
QG070-06	B08	HQP004411	NM_022552	DNMT3A
QG070-06	B09	HQP004309	NM_000791	DHFR
QG070-06	B10	HQP004242	NM_194320	ZNF169
QG070-06	B11	HQP004186	NM_021010	DEFA5
QG070-06	B12	HQP004182	NM_001925	DEFA4
QG070-06	C01	HQP004115	NM_000790	DDC
QG070-06	C02	HQP004081	NM_000789	ACE
QG070-06	C03	HQP004066	NM_000788	DCK
QG070-06	C04	HQP003855	NM_000779	CYP4B1
QG070-06	C05	HQP003826	NM_001017420	ESCO2
QG070-06	C06	HQP003813	NM_000772	CYP2C18
QG070-06	C07	HQP003810	NM_000770	CYP2C8
QG070-06	C08	HQP003539	NM_001904	CTNNB1
QG070-06	C09	HQP003420	NM_182919	TICAM1
QG070-06	C10	HQP003408	NM_134444	NLRP4
QG070-06	C11	HQP003393	NM_144685	HIPK4
QG070-06	C12	HQP003151	NM_172211	CSF1
QG070-06	D01	HQP003069	NM_139074	DEFB127
QG070-06	D02	HQP002873	NM_000098	CPT2
QG070-06	D03	HQP002869	NM_004377	CPT1B
QG070-06	D04	HQP002353	NM_176811	NLRP8
QG070-06	D05	HQP002209	NM_005508	CCR4
QG070-06	D06	HQP002198	NM_001295	CCR1
QG070-06	D07	HQP002181	NM_080867	SOCS4
QG070-06	D08	HQP001954	NM_054112	DEFB118
QG070-06	D09	HQP001862	NM_000082	ERCC8
QG070-06	D10	HQP001708	NM_001278	CHUK
QG070-06	D11	HQP001608	NM_007284	TWF2
QG070-06	D12	HQP001457	NM_004778	GPR44
QG070-06	E01	HQP001408	NM_007199	IRAK3
QG070-06	E02	HQP001399	NM_007195	POLI
QG070-06	E03	HQP001396	NM_001005735	CHEK2
QG070-06	E04	HQP001384	NM_003465	CHIT1
QG070-06	E05	HQP001351	NM_001008396	WDHD1

Product Data Sheet

QG070-06	E06	HQP001341	NM_007083	NUDT6
QG070-06	E07	HQP001288	NM_007051	FAF1
QG070-06	E08	HQP001287	NM_013441	DSCR1L2
QG070-06	E09	HQP000973	NM_006664	CCL27
QG070-06	E10	HQP000888	NM_006615	CAPN9
QG070-06	E11	HQP000657	NM_006407	ARL6IP5
QG070-06	E12	HQP000586	NM_006363	SEC23B
QG070-06	F01	HQP000421	NM_006072	CCL26
QG070-06	F02	HQP000274	NM_001259	CDK6
QG070-06	F03	HQP000245	NM_000075	CDK4
QG070-06	F04	HQP000020	NM_133484	TANK
QG070-06	F05	HQP054687	NM_005961	MUC6
QG070-06	F06	HQP018996	NM_021160	BAT5
QG070-06	F07	HQP018180	NM_005427	TP73
QG070-06	F08	HQP018159	NM_000491	C1QB
QG070-06	F09	HQP017618	NM_003102	SOD3
QG070-06	F10	HQP016651	NM_002995	XCL1
QG070-06	F11	HQP015394	NM_020375	C12orf5
QG070-06	F12	HQP015091	NM_001031849	MASP1
QG070-06	G01	HQP012805	NM_016463	CXXC5
QG070-06	G02	HQP011865	NM_000620	NOS1
QG070-06	G03	HQP011571	NM_152291	MUC7
QG070-06	G04	HQP010851	NM_006151	LPO
QG070-06	G05	HQP010101	NM_001002231	KLK2
QG070-06	G06	HQP009192	NM_002162	ICAM3
QG070-06	G07	HQP005588	NM_207341	ZP1
QG070-06	G08	HQP004414	NM_006892	DNMT3B
QG070-06	G09	HQP004189	NM_001926	DEFA6
QG070-06	G10	HQP002354	NM_153447	NLRP5
QG070-06	G11	HQP002352	NM_176810	NLRP13
QG070-06	G12	HQP001979	NM_001285	CLCA1
QG070-06	H01	HGDC		
QG070-06	H02	HGDC		
QG070-06	H03	HQP006940	NM_002046	GAPDH
QG070-06	H04	HQP016381	NM_001101	ACTB
QG070-06	H05	HQP015171	NM_004048	B2M
QG070-06	H06	HQP006171	NM_012423	RPL13A
QG070-06	H07	HQP009026	NM_000194	HPRT1
QG070-06	H08	HQP054253	NR_003286	RN18S1
QG070-06	H09	RT		
QG070-06	H10	RT		
QG070-06	H11	PCR		
QG070-06	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Leukemia 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