

ExProfile™ Human Lung Cancer Gene qPCR Array

For focused group profiling of human lung cancer related genes expression

Cat. No. QG071-A (6 x 96-well plate, Format A)

Cat. No. QG071-B (6 x 96-well plate, Format B)

Cat. No. QG071-C (6 x 96-well plate, Format C)

Cat. No. QG071-D (6 x 96-well plate, Format D)

Cat. No. QG071-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 lung cancer gene qPCR array profiles 504 human genes to aberrantly expressed human genes involved in human lung cancer. 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 lung cancer gene array allows researchers to study the cancer-related genes to gain understanding of their roles of genes in head and lung cancer pathogenesis.

- QG071 plate 01: 84 unique gene PCR primer pairs
- QG071 plate 02: 84 unique gene PCR primer pairs
- QG071 plate 03: 84 unique gene PCR primer pairs
- QG071 plate 04: 84 unique gene PCR primer pairs
- QG071 plate 05: 84 unique gene PCR primer pairs
- QG071 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 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	EGFR	TP53	CYP1A1	KRAS	XRCC1	ERCC2	MPO	ERCC1	OGG1	EPHX1	NQO1	NAT2
B	MTHFR	CYP2E1	CHRNA5	XPC	MDM2	CHRNA3	TNF	APEX1	IL1B	MMP1	XRCC3	PTGS2
C	ABCB1	IL6	CYP1B1	SOD2	ERCC5	GSTM3	VEGFA	IL8	TGFB1	IL10	STK11	CCND1
D	FASLG	CYP3A5	CLPTM1L	MMP3	ERCC4	CYP2D6	CYP2C19	TP73	XRCC2	NOS3	NBN	MLH1
E	ERCC6	DRD2	TYMS	SULT1A1	BRCA2	PPARG	MMP9	MMP2	LIG1	IL13	IL1RN	IL1A
F	PARP1	COMT	EXO1	NAT1	BAT3	CASP8	XRCC4	BRAF	SLC19A1	PTEN	PSMA4	MTRR
G	MSH2	MMP12	MET	IL4	IGFBP3	HLA-A	ALDH2	AHR	CYP19A1	CYP17A1	CHRN4	CHEK2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG071 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	CDKN2A	TGFBR1	CCNH	MBD4	GGH	TNFRSF10A	CBS	AKR1C3	CASP3	VDR	RAD23B	SERPINA1
B	MTR	LIG3	IL12B	IL12A	IL8RA	IL6R	IL4R	IL2	IFNAR2	GSTA4	AKT1	ACE
C	CYP2C9	LOC123688	MYCL1	RRM1	HLA-DRB1	CDA	XRCC5	BCR	POLB	PCNA	LIG4	IFNGR2
D	HSD17B1	RGS17	FGFR4	ESR1	DRD4	CYP2A13	CSF2	RASSF1	SOD3	RET	CDH1	CDC25C
E	RGS6	TP53I3	TNKS	CASP7	TP53BP1	TERF2	TERF1	SFTPB	NOD2	BHMT	CCL2	RAD52
F	BARD1	ATR	PIM1	PIK3CA	REV1	NOS2A	MYC	ABCC1	LRMP	LEP	IL15RA	IL15
G	IL7R	IGF1R	IGF1	HSD3B1	APC	MSH6	GSTZ1	GSTA2	GPX3	GPX2	POT1	FHIT
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG071 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	ERCC3	ELA2	EGF	DHFR	CYP11A1	CTLA4	CSF3	SERPINA3	ADH1C	POLI	CHEK1	ALDH1L1
B	CDK5	ERBB4	DNMT3B	ABCC4	SLC2A4	SHBG	NR1H4	SLC23A1	CDC25B	CDC25A	RB1CC1	MATR3
C	CD81	CD40	MPDU1	GDF15	LITAF	SCARB1	CD86	CYP7B1	CD80	CER1	CD4	SLC28A1
D	PDCD5	MBD2	CBR3	CBR1	ABCC3	ABCB11	CAV1	RGS5	CASR	RAD54L	PLA2G6	CASP5
E	FZD7	AXIN2	NCOA3	COL18A1	CALCR	ZNF230	XBP1	WRN	VIL2	VCAM1	UCP3	TYR
F	TSG101	TNFRSF1B	TGM1	TFF1	TEP1	STAT1	SSTR3	SOD1	SLC4A2	SLAMF1	SFTPD	ABCG8
G	BLM	SEPP1	SELE	CCL5	RXRβ	RXRA	ROS1	RNASEL	BCL2	ZNF350	RAC1	PCTP
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG071 plate 03

Product Data Sheet

	1	2	3	4	5	6	7	8	9	10	11	12
A	BAK1	PTGS1	RAD18	MYNN	HIF1AN	FBXW7	CASC1	WDR79	UGT1A4	UGT1A7	POLD1	TLR9
B	PMS2	PMS1	PLK1	PLAUR	PLAU	PLA2G2A	SERPINE2	PHB	PGR	P2RX7	OPRM1	OCA2
C	ATP1B2	NINJ1	NFKBIE	NFKBIA	NFKB1	NCF2	NUBP1	MX1	MSH3	MIF	MEST	MBD1
D	MAOA	ARNT	LRP5	LRP6	LPL	LIPC	ARHGDI3	LEPR	LDLR	LCAT	KDR	RAB15
E	JAK3	ITGB3	IRS1	IRF3	IRF1	IREB2	INSR	IL18	IL12RB1	IL10RB	IL2RB	IL1R1
F	APOE	IGFBP6	IGFBP5	IGFBP2	IGFBP1	IGF2	IFNGR1	IFNAR1	APOA4	APOA2	HUS1	HTR1D
G	HTR1B	BIRC5	BIRC2	HSD17B4	HSD17B2	HSD3B2	APAF1	HMGCR	HIF1A	HFE	HADHA	SLC39A2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG071 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	RPA4	GSK3B	GRPR	GPX4	AMT	DHDH	GC	HSPB8	KRT23	LMOD1	XRCC6	ALOX15
B	ALOX5	ALOX12	AMACR	FPGS	SEZ6L	SEC14L2	FOS	ABCA5	ABCA6	FOXC1	ITGA11	GPC5
C	FANCA	ESR2	AKT2	EPHX2	DNAJC18	ABCA1	EFNB3	EDN1	DRD1	CYP27B1	CYP24A1	CYP2F1
D	CYBB	CX3CR1	CTSH	CTNBN1	CTH	CSTF1	PARP4	CSF1R	CCR5	SAT2	AKR1C4	CGA
E	MTHFD2	CETP	SLC01B1	AKR1A1	CDKN1C	CDK7	CDK4	AKAP9	NUBP2	PTH	PAK6	MASP1
F	OPRD1	LMO2	IGFALS	ALAD	FAM82A	TNIP1	LOC729230	SUMO1	KLF6	CDC2	KEAP1	HDAC9
G	GSTO1	41532	CD14	DDX18	NR112	CFLAR	EIF3S3	MFSD2	HAVCR2	CASP6	MAD1L1	TFPI2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG071 plate 05

	1	2	3	4	5	6	7	8	9	10	11	12
A	SUV39H2	NEIL1	BHLHB3	FTO	SEMA3B	PRDM2	SLC30A4	SLC30A2	SCGB1A1	UBE2I	TWIST1	C5AR1
B	BUB1B	TDG	STAT3	SLC22A3	SLC30A5	CYP3A43	MAP2K4	CXCL12	SAT1	S100A2	RGS2	RFC1
C	RB1	RASA1	RAGE	PXN	PTPRJ	PTPN13	CYP4F11	XPO5	PTGIS	PTGER2	MAP2K1	APOM
D	LAPTM4B	PPP2R1B	TRIT1	PPARD	PPARA	PON1	TERF2IP	POLE	HSPA14	SH3GLB1	GEMIN4	SERPINB2
E	SERPINE1	ATP2A3	NRAS	NME1	NFE2L2	PPP1R12A	MUC1	MSRA	MPG	MNAT1	MMP8	MLLT3
F	ARG2	ARG1	KIT	ITGB8	AR	IL8RB	IL5RA	IL5	IL1RAP	IDH2	IDH1	APOB
G	BIRC4	HSPB1	HSPA8	HRAS	HPRT1	HPGD	CARD10	GTF2H1	GSTM2	GSTA3	GSS	GSR
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure6. Illustration of QG071 plate 06

- **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
QG071-01	A01	HQP004605	NM_005228	EGFR
QG071-01	A02	HQP018175	NM_000546	TP53
QG071-01	A03	HQP003772	NM_000499	CYP1A1
QG071-01	A04	HQP010133	NM_004985	KRAS
QG071-01	A05	HQP018562	NM_006297	XRCC1

Product Data Sheet

QG071-01	A06	HQP004976	NM_000400	ERCC2
QG071-01	A07	HQP011309	NM_000250	MPO
QG071-01	A08	HQP004975	NM_202001	ERCC1
QG071-01	A09	HQP012021	NM_002542	OGG1
QG071-01	A10	HQP004948	NM_000120	EPHX1
QG071-01	A11	HQP004317	NM_000903	NQO1
QG071-01	A12	HQP001136	NM_000015	NAT2
QG071-01	B01	HQP011547	NM_005957	MTHFR
QG071-01	B02	HQP003817	NM_000773	CYP2E1
QG071-01	B03	HQP001643	NM_000745	CHRNA5
QG071-01	B04	HQP018556	NM_004628	XPC
QG071-01	B05	HQP011135	NM_002392	MDM2
QG071-01	B06	HQP001632	NM_000743	CHRNA3
QG071-01	B07	HQP018141	NM_000594	TNF
QG071-01	B08	HQP009061	NM_001641	APEX1
QG071-01	B09	HQP009641	NM_000576	IL1B
QG071-01	B10	HQP011255	NM_002421	MMP1
QG071-01	B11	HQP018564	NM_005432	XRCC3
QG071-01	B12	HQP015598	NM_000963	PTGS2
QG071-01	C01	HQP013100	NM_000927	ABCB1
QG071-01	C02	HQP009670	NM_000600	IL6
QG071-01	C03	HQP003775	NM_000104	CYP1B1
QG071-01	C04	HQP017616	NM_000636	SOD2
QG071-01	C05	HQP004985	NM_000123	ERCC5
QG071-01	C06	HQP008483	NM_000849	GSTM3
QG071-01	C07	HQP018475	NM_001025366	VEGFA
QG071-01	C08	HQP009678	NM_000584	IL8
QG071-01	C09	HQP018044	NM_000660	TGFB1
QG071-01	C10	HQP009685	NM_000572	IL10
QG071-01	C11	HQP017794	NM_000455	STK11
QG071-01	C12	HQP016204	NM_053056	CCND1
QG071-01	D01	HQP009671	NM_000639	FASLG
QG071-01	D02	HQP003841	NM_000777	CYP3A5
QG071-01	D03	HQP019829	NM_030782	CLPTM1L
QG071-01	D04	HQP011257	NM_002422	MMP3
QG071-01	D05	HQP004984	NM_005236	ERCC4
QG071-01	D06	HQP003814	NM_000106	CYP2D6
QG071-01	D07	HQP003809	NM_000769	CYP2C19
QG071-01	D08	HQP018180	NM_005427	TP73
QG071-01	D09	HQP018563	NM_005431	XRCC2
QG071-01	D10	HQP011868	NM_000603	NOS3
QG071-01	D11	HQP011687	NM_002485	NBN
QG071-01	D12	HQP011235	NM_000249	MLH1
QG071-01	E01	HQP004986	NM_000124	ERCC6
QG071-01	E02	HQP004446	NM_000795	DRD2
QG071-01	E03	HQP018342	NM_001071	TYMS
QG071-01	E04	HQP017811	NM_177536	SULT1A1
QG071-01	E05	HQP017753	NM_000059	BRCA2

Product Data Sheet

QG071-01	E06	HQP013633	NM_005037	PPARG
QG071-01	E07	HQP011263	NM_004994	MMP9
QG071-01	E08	HQP011256	NM_004530	MMP2
QG071-01	E09	HQP010609	NM_000234	LIG1
QG071-01	E10	HQP009697	NM_002188	IL13
QG071-01	E11	HQP009645	NM_000577	IL1RN
QG071-01	E12	HQP009640	NM_000575	IL1A
QG071-01	F01	HQP003120	NM_001618	PARP1
QG071-01	F02	HQP002671	NM_000754	COMT
QG071-01	F03	HQP054030	NM_130398	EXO1
QG071-01	F04	HQP023467	NM_000662	NAT1
QG071-01	F05	HQP018983	NM_004639	BAT3
QG071-01	F06	HQP018966	NM_001080124	CASP8
QG071-01	F07	HQP018565	NM_003401	XRCC4
QG071-01	F08	HQP017733	NM_004333	BRAF
QG071-01	F09	HQP017482	NM_194255	SLC19A1
QG071-01	F10	HQP015535	NM_000314	PTEN
QG071-01	F11	HQP015192	NM_002789	PSMA4
QG071-01	F12	HQP011555	NM_002454	MTRR
QG071-01	G01	HQP011491	NM_000251	MSH2
QG071-01	G02	HQP011266	NM_002426	MMP12
QG071-01	G03	HQP011181	NM_000245	MET
QG071-01	G04	HQP009662	NM_000589	IL4
QG071-01	G05	HQP009544	NM_000598	IGFBP3
QG071-01	G06	HQP008849	NM_002116	HLA-A
QG071-01	G07	HQP005087	NM_000690	ALDH2
QG071-01	G08	HQP004658	NM_001621	AHR
QG071-01	G09	HQP003904	NM_000103	CYP19A1
QG071-01	G10	HQP003888	NM_000102	CYP17A1
QG071-01	G11	HQP001667	NM_000750	CHRN4
QG071-01	G12	HQP001396	NM_001005735	CHEK2
QG071-01	H01	HGDC		
QG071-01	H02	HGDC		
QG071-01	H03	HQP006940	NM_002046	GAPDH
QG071-01	H04	HQP016381	NM_001101	ACTB
QG071-01	H05	HQP015171	NM_004048	B2M
QG071-01	H06	HQP006171	NM_012423	RPL13A
QG071-01	H07	HQP009026	NM_000194	HPRT1
QG071-01	H08	HQP054253	NR_003286	RN18S1
QG071-01	H09	RT		
QG071-01	H10	RT		
QG071-01	H11	PCR		
QG071-01	H12	PCR		
QG071-02	A01	HQP000370	NM_058195	CDKN2A
QG071-02	A02	HQP054057	BC071181	TGFB1
QG071-02	A03	HQP021906	NM_001239	CCNH
QG071-02	A04	HQP021736	NM_003925	MBD4
QG071-02	A05	HQP021603	NM_003878	GGH

Product Data Sheet

QG071-02	A06	HQP021557	NM_003844	TNFRSF10A
QG071-02	A07	HQP021518	NM_000071	CBS
QG071-02	A08	HQP021385	NM_003739	AKR1C3
QG071-02	A09	HQP020297	NM_004346	CASP3
QG071-02	A10	HQP018474	NM_000376	VDR
QG071-02	A11	HQP016076	NM_002874	RAD23B
QG071-02	A12	HQP013122	NM_000295	SERPINA1
QG071-02	B01	HQP011554	NM_000254	MTR
QG071-02	B02	HQP010612	NM_013975	LIG3
QG071-02	B03	HQP009693	NM_002187	IL12B
QG071-02	B04	HQP009692	NM_000882	IL12A
QG071-02	B05	HQP009679	NM_000634	IL8RA
QG071-02	B06	HQP009672	NM_000565	IL6R
QG071-02	B07	HQP009664	NM_000418	IL4R
QG071-02	B08	HQP009649	NM_000586	IL2
QG071-02	B09	HQP009460	NM_000874	IFNAR2
QG071-02	B10	HQP008479	NM_001512	GSTA4
QG071-02	B11	HQP004991	NM_001014431	AKT1
QG071-02	B12	HQP004081	NM_000789	ACE
QG071-02	C01	HQP003811	NM_000771	CYP2C9
QG071-02	C02	HQP002215	NM_001013619	LOC123688
QG071-02	C03	HQP011601	NM_005376	MYCL1
QG071-02	C04	HQP054684	NM_001033	RRM1
QG071-02	C05	HQP054047	BC008403	HLA-DRB1
QG071-02	C06	HQP023203	NM_001785	CDA
QG071-02	C07	HQP018568	NM_021141	XRCC5
QG071-02	C08	HQP016418	NM_004327	BCR
QG071-02	C09	HQP013422	NM_002690	POLB
QG071-02	C10	HQP012420	NM_002592	PCNA
QG071-02	C11	HQP010613	NM_002312	LIG4
QG071-02	C12	HQP009472	NM_005534	IFNGR2
QG071-02	D01	HQP009064	NM_000413	HSD17B1
QG071-02	D02	HQP007304	NM_012419	RGS17
QG071-02	D03	HQP005440	NM_022963	FGFR4
QG071-02	D04	HQP004998	NM_000125	ESR1
QG071-02	D05	HQP004450	NM_000797	DRD4
QG071-02	D06	HQP003804	NM_000766	CYP2A13
QG071-02	D07	HQP003159	NM_000758	CSF2
QG071-02	D08	HQP001374	NM_007182	RASSF1
QG071-02	D09	HQP017618	NM_003102	SOD3
QG071-02	D10	HQP054045	BC004257	RET
QG071-02	D11	HQP023466	NM_004360	CDH1
QG071-02	D12	HQP023409	NM_001790	CDC25C
QG071-02	E01	HQP023003	NM_004296	RGS6
QG071-02	E02	HQP022884	NM_004881	TP53I3
QG071-02	E03	HQP021405	NM_003747	TNKS
QG071-02	E04	HQP020481	NM_033338	CASP7
QG071-02	E05	HQP018176	NM_005657	TP53BP1

Product Data Sheet

QG071-02	E06	HQP018016	NM_005652	TERF2
QG071-02	E07	HQP018014	NM_003218	TERF1
QG071-02	E08	HQP016958	NM_000542	SFTPBP
QG071-02	E09	HQP016801	NM_022162	NOD2
QG071-02	E10	HQP016634	NM_001713	BHMT
QG071-02	E11	HQP016621	NM_002982	CCL2
QG071-02	E12	HQP016087	NM_134424	RAD52
QG071-02	F01	HQP015946	NM_000465	BARD1
QG071-02	F02	HQP013596	NM_001184	ATR
QG071-02	F03	HQP013152	NM_002648	PIM1
QG071-02	F04	HQP013150	NM_006218	PIK3CA
QG071-02	F05	HQP012755	NM_001037872	REV1
QG071-02	F06	HQP011866	NM_000625	NOS2A
QG071-02	F07	HQP011597	NM_002467	MYC
QG071-02	F08	HQP011325	NM_019899	ABCC1
QG071-02	F09	HQP010868	NM_006152	LRMP
QG071-02	F10	HQP010581	NM_000230	LEP
QG071-02	F11	HQP009710	NM_002189	IL15RA
QG071-02	F12	HQP009708	NM_000585	IL15
QG071-02	G01	HQP009677	NM_002185	IL7R
QG071-02	G02	HQP009523	NM_000875	IGF1R
QG071-02	G03	HQP009518	NM_000618	IGF1
QG071-02	G04	HQP009059	NM_000862	HSD3B1
QG071-02	G05	HQP009024	NM_000038	APC
QG071-02	G06	HQP008493	NM_000179	MSH6
QG071-02	G07	HQP008490	NM_001513	GSTZ1
QG071-02	G08	HQP008476	NM_000846	GSTA2
QG071-02	G09	HQP008282	NM_002084	GPX3
QG071-02	G10	HQP008281	NM_002083	GPX2
QG071-02	G11	HQP006855	NM_001042594	POT1
QG071-02	G12	HQP005447	NM_002012	FHIT
QG071-02	H01	HGDC		
QG071-02	H02	HGDC		
QG071-02	H03	HQP006940	NM_002046	GAPDH
QG071-02	H04	HQP016381	NM_001101	ACTB
QG071-02	H05	HQP015171	NM_004048	B2M
QG071-02	H06	HQP006171	NM_012423	RPL13A
QG071-02	H07	HQP009026	NM_000194	HPRT1
QG071-02	H08	HQP054253	NR_003286	RN18S1
QG071-02	H09	RT		
QG071-02	H10	RT		
QG071-02	H11	PCR		
QG071-02	H12	PCR		
QG071-03	A01	HQP004983	NM_000122	ERCC3
QG071-03	A02	HQP004689	NM_001972	ELA2
QG071-03	A03	HQP004599	NM_001963	EGF
QG071-03	A04	HQP004309	NM_000791	DHFR
QG071-03	A05	HQP003871	NM_000781	CYP11A1

Product Data Sheet

QG071-03	A06	HQP003499	NM_001037631	CTLA4
QG071-03	A07	HQP003173	NM_000759	CSF3
QG071-03	A08	HQP002590	NM_001085	SERPINA3
QG071-03	A09	HQP002413	NM_000669	ADH1C
QG071-03	A10	HQP001399	NM_007195	POLI
QG071-03	A11	HQP001282	NM_001274	CHEK1
QG071-03	A12	HQP000963	NM_012190	ALDH1L1
QG071-03	B01	HQP000261	NM_004935	CDK5
QG071-03	B02	HQP004972	NM_001042599	ERBB4
QG071-03	B03	HQP004414	NM_006892	DNMT3B
QG071-03	B04	HQP000320	NM_005845	ABCC4
QG071-03	B05	HQP053960	NM_001042	SLC2A4
QG071-03	B06	HQP053959	NM_001040	SHBG
QG071-03	B07	HQP023437	NM_005123	NR1H4
QG071-03	B08	HQP023414	NM_005847	SLC23A1
QG071-03	B09	HQP023399	NM_004358	CDC25B
QG071-03	B10	HQP023385	NM_001789	CDC25A
QG071-03	B11	HQP023240	NM_014781	RB1CC1
QG071-03	B12	HQP023195	NM_018834	MATR3
QG071-03	C01	HQP023168	NM_004356	CD81
QG071-03	C02	HQP022955	NM_001250	CD40
QG071-03	C03	HQP022863	NM_004870	MPDU1
QG071-03	C04	HQP022853	NM_004864	GDF15
QG071-03	C05	HQP022851	NM_004862	LITAF
QG071-03	C06	HQP022833	NM_005505	SCARB1
QG071-03	C07	HQP022747	NM_175862	CD86
QG071-03	C08	HQP022723	NM_004820	CYP7B1
QG071-03	C09	HQP022722	NM_005191	CD80
QG071-03	C10	HQP022580	NM_005454	CER1
QG071-03	C11	HQP022316	NM_000616	CD4
QG071-03	C12	HQP022202	NM_004213	SLC28A1
QG071-03	D01	HQP022170	NM_004708	PDCD5
QG071-03	D02	HQP021737	NM_003927	MBD2
QG071-03	D03	HQP021508	NM_001236	CBR3
QG071-03	D04	HQP021495	NM_001757	CBR1
QG071-03	D05	HQP021459	NM_003786	ABCC3
QG071-03	D06	HQP021388	NM_003742	ABCB11
QG071-03	D07	HQP021313	NM_001753	CAV1
QG071-03	D08	HQP021023	NM_003617	RGS5
QG071-03	D09	HQP020899	NM_000388	CASR
QG071-03	D10	HQP020707	NM_003579	RAD54L
QG071-03	D11	HQP020418	NM_001004426	PLA2G6
QG071-03	D12	HQP020384	NM_004347	CASP5
QG071-03	E01	HQP020136	NM_003507	FZD7
QG071-03	E02	HQP020116	NM_004655	AXIN2
QG071-03	E03	HQP020040	NM_006534	NCOA3
QG071-03	E04	HQP019768	NM_130445	COL18A1
QG071-03	E05	HQP019452	NM_001742	CALCR

Product Data Sheet

QG071-03	E06	HQP018739	NM_006300	ZNF230
QG071-03	E07	HQP018550	NM_001079539	XBP1
QG071-03	E08	HQP018545	NM_000553	WRN
QG071-03	E09	HQP018488	NM_003379	VIL2
QG071-03	E10	HQP018466	NM_080682	VCAM1
QG071-03	E11	HQP018404	NM_003356	UCP3
QG071-03	E12	HQP018343	NM_000372	TYR
QG071-03	F01	HQP018269	NM_006292	TSG101
QG071-03	F02	HQP018149	NM_001066	TNFRSF1B
QG071-03	F03	HQP018060	NM_000359	TGM1
QG071-03	F04	HQP018035	NM_003225	TFF1
QG071-03	F05	HQP018013	NM_007110	TEP1
QG071-03	F06	HQP017764	NM_007315	STAT1
QG071-03	F07	HQP017745	NM_001051	SSTR3
QG071-03	F08	HQP017615	NM_000454	SOD1
QG071-03	F09	HQP017367	NM_003040	SLC4A2
QG071-03	F10	HQP017303	NM_003037	SLAMF1
QG071-03	F11	HQP016980	NM_003019	SFTPD
QG071-03	F12	HQP016876	NM_022437	ABCG8
QG071-03	G01	HQP016844	NM_000057	BLM
QG071-03	G02	HQP016819	NM_005410	SEPP1
QG071-03	G03	HQP016744	NM_000450	SELE
QG071-03	G04	HQP016626	NM_002985	CCL5
QG071-03	G05	HQP016527	NM_021976	RXRБ
QG071-03	G06	HQP016526	NM_002957	RXRA
QG071-03	G07	HQP016380	NM_002944	ROS1
QG071-03	G08	HQP016303	NM_021133	RNASEL
QG071-03	G09	HQP016212	NM_000657	BCL2
QG071-03	G10	HQP016170	NM_021632	ZNF350
QG071-03	G11	HQP016064	NM_018890	RAC1
QG071-03	G12	HQP016003	NM_021213	PCTP
QG071-03	H01	HGDC		
QG071-03	H02	HGDC		
QG071-03	H03	HQP006940	NM_002046	GAPDH
QG071-03	H04	HQP016381	NM_001101	ACTB
QG071-03	H05	HQP015171	NM_004048	B2M
QG071-03	H06	HQP006171	NM_012423	RPL13A
QG071-03	H07	HQP009026	NM_000194	HPRT1
QG071-03	H08	HQP054253	NR_003286	RN18S1
QG071-03	H09	RT		
QG071-03	H10	RT		
QG071-03	H11	PCR		
QG071-03	H12	PCR		
QG071-04	A01	HQP015917	NM_001188	BAK1
QG071-04	A02	HQP015596	NM_000962	PTGS1
QG071-04	A03	HQP015190	NM_020165	RAD18
QG071-04	A04	HQP014808	NM_018657	MYNN
QG071-04	A05	HQP014564	NM_017902	HIF1AN

Product Data Sheet

QG071-04	A06	HQP014292	NM_018315	FBXW7
QG071-04	A07	HQP014243	NM_018272	CASC1
QG071-04	A08	HQP014089	NM_018081	WDR79
QG071-04	A09	HQP013614	NM_007120	UGT1A4
QG071-04	A10	HQP013581	NM_019077	UGT1A7
QG071-04	A11	HQP013423	NM_002691	POLD1
QG071-04	A12	HQP013388	NM_017442	TLR9
QG071-04	B01	HQP013352	NM_000535	PMS2
QG071-04	B02	HQP013312	NM_000534	PMS1
QG071-04	B03	HQP013262	NM_005030	PLK1
QG071-04	B04	HQP013205	NM_001005376	PLAUR
QG071-04	B05	HQP013204	NM_002658	PLAU
QG071-04	B06	HQP013194	NM_000300	PLA2G2A
QG071-04	B07	HQP013128	NM_006216	SERPINE2
QG071-04	B08	HQP013104	NM_002634	PHB
QG071-04	B09	HQP013099	NM_000926	PGR
QG071-04	B10	HQP012100	NM_002562	P2RX7
QG071-04	B11	HQP012062	NM_001008505	OPRM1
QG071-04	B12	HQP012008	NM_000275	OCA2
QG071-04	C01	HQP011852	NM_001678	ATP1B2
QG071-04	C02	HQP011836	NM_004148	NINJ1
QG071-04	C03	HQP011813	NM_004556	NFKBIE
QG071-04	C04	HQP011810	NM_020529	NFKBIA
QG071-04	C05	HQP011807	NM_003998	NFKB1
QG071-04	C06	HQP011693	NM_000433	NCF2
QG071-04	C07	HQP011685	NM_002484	NUBP1
QG071-04	C08	HQP011582	NM_002462	MX1
QG071-04	C09	HQP011492	NM_002439	MSH3
QG071-04	C10	HQP011219	NM_002415	MIF
QG071-04	C11	HQP011179	NM_002402	MEST
QG071-04	C12	HQP011073	NM_015844	MBD1
QG071-04	D01	HQP011007	NM_000240	MAOA
QG071-04	D02	HQP010924	NM_001668	ARNT
QG071-04	D03	HQP010878	NM_002335	LRP5
QG071-04	D04	HQP010877	NM_002336	LRP6
QG071-04	D05	HQP010847	NM_000237	LPL
QG071-04	D06	HQP010626	NM_000236	LIPC
QG071-04	D07	HQP010610	NM_001175	ARHGDI3
QG071-04	D08	HQP010584	NM_002303	LEPR
QG071-04	D09	HQP010577	NM_000527	LDLR
QG071-04	D10	HQP010564	NM_000229	LCAT
QG071-04	D11	HQP010070	NM_002253	KDR
QG071-04	D12	HQP010008	NM_198686	RAB15
QG071-04	E01	HQP009851	NM_000215	JAK3
QG071-04	E02	HQP009818	NM_000212	ITGB3
QG071-04	E03	HQP009788	NM_005544	IRS1
QG071-04	E04	HQP009780	NM_001571	IRF3
QG071-04	E05	HQP009778	NM_002198	IRF1

Product Data Sheet

QG071-04	E06	HQP009777	NM_004136	IREB2
QG071-04	E07	HQP009764	NM_000208	INSR
QG071-04	E08	HQP009718	NM_001562	IL18
QG071-04	E09	HQP009694	NM_005535	IL12RB1
QG071-04	E10	HQP009687	NM_000628	IL10RB
QG071-04	E11	HQP009658	NM_000878	IL2RB
QG071-04	E12	HQP009642	NM_000877	IL1R1
QG071-04	F01	HQP009556	NM_000041	APOE
QG071-04	F02	HQP009555	NM_002178	IGFBP6
QG071-04	F03	HQP009551	NM_000599	IGFBP5
QG071-04	F04	HQP009541	NM_000597	IGFBP2
QG071-04	F05	HQP009539	NM_000596	IGFBP1
QG071-04	F06	HQP009529	NM_000612	IGF2
QG071-04	F07	HQP009469	NM_000416	IFNGR1
QG071-04	F08	HQP009458	NM_000629	IFNAR1
QG071-04	F09	HQP009171	NM_000482	APOA4
QG071-04	F10	HQP009139	NM_001643	APOA2
QG071-04	F11	HQP009138	NM_004507	HUS1
QG071-04	F12	HQP009117	NM_000864	HTR1D
QG071-04	G01	HQP009116	NM_000863	HTR1B
QG071-04	G02	HQP009101	NM_001168	BIRC5
QG071-04	G03	HQP009072	NM_001166	BIRC2
QG071-04	G04	HQP009067	NM_000414	HSD17B4
QG071-04	G05	HQP009066	NM_002153	HSD17B2
QG071-04	G06	HQP009060	NM_000198	HSD3B2
QG071-04	G07	HQP008934	NM_001160	APAF1
QG071-04	G08	HQP008891	NM_000859	HMGCR
QG071-04	G09	HQP008832	NM_181054	HIF1A
QG071-04	G10	HQP008757	NM_000410	HFE
QG071-04	G11	HQP008718	NM_000182	HADHA
QG071-04	G12	HQP008659	NM_014579	SLC39A2
QG071-04	H01	HGDC		
QG071-04	H02	HGDC		
QG071-04	H03	HQP006940	NM_002046	GAPDH
QG071-04	H04	HQP016381	NM_001101	ACTB
QG071-04	H05	HQP015171	NM_004048	B2M
QG071-04	H06	HQP006171	NM_012423	RPL13A
QG071-04	H07	HQP009026	NM_000194	HPRT1
QG071-04	H08	HQP054253	NR_003286	RN18S1
QG071-04	H09	RT		
QG071-04	H10	RT		
QG071-04	H11	PCR		
QG071-04	H12	PCR		
QG071-05	A01	HQP008605	NM_013347	RPA4
QG071-05	A02	HQP008469	NM_002093	GSK3B
QG071-05	A03	HQP008464	NM_005314	GRPR
QG071-05	A04	HQP008285	NM_002085	GPX4
QG071-05	A05	HQP007737	NM_000481	AMT

Product Data Sheet

QG071-05	A06	HQP007643	NM_014475	DHDH
QG071-05	A07	HQP007230	NM_000583	GC
QG071-05	A08	HQP007224	NM_014365	HSPB8
QG071-05	A09	HQP006945	NM_015515	KRT23
QG071-05	A10	HQP006733	NM_012134	LMOD1
QG071-05	A11	HQP006547	NM_001469	XRCC6
QG071-05	A12	HQP006425	NM_001140	ALOX15
QG071-05	B01	HQP006359	NM_000698	ALOX5
QG071-05	B02	HQP006356	NM_000697	ALOX12
QG071-05	B03	HQP006261	NM_014324	AMACR
QG071-05	B04	HQP006228	NM_001018078	FPGS
QG071-05	B05	HQP006199	NM_021115	SEZ6L
QG071-05	B06	HQP006189	NM_012429	SEC14L2
QG071-05	B07	HQP006188	NM_005252	FOS
QG071-05	B08	HQP006120	NM_018672	ABCA5
QG071-05	B09	HQP006119	NM_080284	ABCA6
QG071-05	B10	HQP005629	NM_001453	FOXC1
QG071-05	B11	HQP005459	NM_001004439	ITGA11
QG071-05	B12	HQP005436	NM_004466	GPC5
QG071-05	C01	HQP005081	NM_000135	FANCA
QG071-05	C02	HQP005002	NM_001437	ESR2
QG071-05	C03	HQP004995	NM_001626	AKT2
QG071-05	C04	HQP004950	NM_001979	EPHX2
QG071-05	C05	HQP004849	NM_152686	DNAJC18
QG071-05	C06	HQP004727	NM_005502	ABCA1
QG071-05	C07	HQP004598	NM_001406	EFNB3
QG071-05	C08	HQP004557	NM_001955	EDN1
QG071-05	C09	HQP004445	NM_000794	DRD1
QG071-05	C10	HQP003923	NM_000785	CYP27B1
QG071-05	C11	HQP003917	NM_000782	CYP24A1
QG071-05	C12	HQP003818	NM_000774	CYP2F1
QG071-05	D01	HQP003741	NM_000397	CYBB
QG071-05	D02	HQP003691	NM_001337	CX3CR1
QG071-05	D03	HQP003620	NM_004390	CTSH
QG071-05	D04	HQP003539	NM_001904	CTNNB1
QG071-05	D05	HQP003493	NM_001902	CTH
QG071-05	D06	HQP003395	NM_001033521	CSTF1
QG071-05	D07	HQP003171	NM_006437	PARP4
QG071-05	D08	HQP003158	NM_005211	CSF1R
QG071-05	D09	HQP002210	NM_000579	CCR5
QG071-05	D10	HQP001449	NM_133491	SAT2
QG071-05	D11	HQP001262	NM_001818	AKR1C4
QG071-05	D12	HQP000956	NM_000735	CGA
QG071-05	E01	HQP000932	NM_006636	MTHFD2
QG071-05	E02	HQP000853	NM_000078	CETP
QG071-05	E03	HQP000720	NM_006446	SLCO1B1
QG071-05	E04	HQP000405	NM_006066	AKR1A1
QG071-05	E05	HQP000356	NM_000076	CDKN1C

Product Data Sheet

QG071-05	E06	HQP000285	NM_001799	CDK7
QG071-05	E07	HQP000245	NM_000075	CDK4
QG071-05	E08	HQP000179	NM_147185	AKAP9
QG071-05	E09	HQP000134	NM_012225	NUBP2
QG071-05	E10	HQP015595	NM_000315	PTH
QG071-05	E11	HQP015245	NM_020168	PAK6
QG071-05	E12	HQP015091	NM_001031849	MASP1
QG071-05	F01	HQP012054	NM_000911	OPRD1
QG071-05	F02	HQP010701	NM_005574	LMO2
QG071-05	F03	HQP009537	NM_004970	IGFALS
QG071-05	F04	HQP005011	NM_000031	ALAD
QG071-05	F05	HQP003625	NM_144713	FAM82A
QG071-05	F06	HQP000394	NM_006058	TNIP1
QG071-05	F07	HQP054686	NM_001123396	LOC729230
QG071-05	F08	HQP053990	NM_003352	SUMO1
QG071-05	F09	HQP053978	NM_001300	KLF6
QG071-05	F10	HQP023261	NM_001786	CDC2
QG071-05	F11	HQP023233	NM_012289	KEAP1
QG071-05	F12	HQP023134	NM_014707	HDAC9
QG071-05	G01	HQP022764	NM_004832	GSTO1
QG071-05	G02	HQP022686	NM_004261	SEP15
QG071-05	G03	HQP022490	NM_000591	CD14
QG071-05	G04	HQP021670	NM_006773	DDX18
QG071-05	G05	HQP021630	NM_003889	NR1I2
QG071-05	G06	HQP021604	NM_003879	CFLAR
QG071-05	G07	HQP021417	NM_003756	EIF3S3
QG071-05	G08	HQP020994	NM_032793	MFSD2
QG071-05	G09	HQP020984	NM_032782	HAVCR2
QG071-05	G10	HQP020427	NM_001226	CASP6
QG071-05	G11	HQP020331	NM_001013836	MAD1L1
QG071-05	G12	HQP019278	NM_006528	TFPI2
QG071-05	H01	HGDC		
QG071-05	H02	HGDC		
QG071-05	H03	HQP006940	NM_002046	GAPDH
QG071-05	H04	HQP016381	NM_001101	ACTB
QG071-05	H05	HQP015171	NM_004048	B2M
QG071-05	H06	HQP006171	NM_012423	RPL13A
QG071-05	H07	HQP009026	NM_000194	HPRT1
QG071-05	H08	HQP054253	NR_003286	RN18S1
QG071-05	H09	RT		
QG071-05	H10	RT		
QG071-05	H11	PCR		
QG071-05	H12	PCR		
QG071-06	A01	HQP019201	NM_024670	SUV39H2
QG071-06	A02	HQP019141	NM_024608	NEIL1
QG071-06	A03	HQP019015	NM_030762	BHLHB3
QG071-06	A04	HQP018895	NM_001080432	FTO
QG071-06	A05	HQP018813	NM_001005914	SEMA3B

Product Data Sheet

QG071-06	A06	HQP018755	NM_001007257	PRDM2
QG071-06	A07	HQP018747	NM_013309	SLC30A4
QG071-06	A08	HQP018744	NM_001004434	SLC30A2
QG071-06	A09	HQP018411	NM_003357	SCGB1A1
QG071-06	A10	HQP018379	NM_003345	UBE2I
QG071-06	A11	HQP018328	NM_000474	TWIST1
QG071-06	A12	HQP018326	NM_001736	C5AR1
QG071-06	B01	HQP018022	NM_001211	BUB1B
QG071-06	B02	HQP017996	NM_003211	TDG
QG071-06	B03	HQP017767	NM_003150	STAT3
QG071-06	B04	HQP017492	NM_021977	SLC22A3
QG071-06	B05	HQP017246	NM_022902	SLC30A5
QG071-06	B06	HQP017181	NM_022820	CYP3A43
QG071-06	B07	HQP016830	NM_003010	MAP2K4
QG071-06	B08	HQP016670	NM_001033886	CXCL12
QG071-06	B09	HQP016572	NM_002970	SAT1
QG071-06	B10	HQP016538	NM_005978	S100A2
QG071-06	B11	HQP016252	NM_002923	RGS2
QG071-06	B12	HQP016224	NM_002913	RFC1
QG071-06	C01	HQP016131	NM_000321	RB1
QG071-06	C02	HQP016125	NM_002890	RASA1
QG071-06	C03	HQP016084	NM_014226	RAGE
QG071-06	C04	HQP015978	NM_002859	PXN
QG071-06	C05	HQP015925	NM_002843	PTPRJ
QG071-06	C06	HQP015897	NM_006264	PTPN13
QG071-06	C07	HQP015892	NM_021187	CYP4F11
QG071-06	C08	HQP015667	NM_020750	XPO5
QG071-06	C09	HQP015586	NM_000961	PTGIS
QG071-06	C10	HQP015544	NM_000956	PTGER2
QG071-06	C11	HQP014907	NM_002755	MAP2K1
QG071-06	C12	HQP014846	NM_019101	APOM
QG071-06	D01	HQP014362	NM_018407	LAPTM4B
QG071-06	D02	HQP014161	NM_002716	PPP2R1B
QG071-06	D03	HQP013708	NM_017646	TRIT1
QG071-06	D04	HQP013627	NM_006238	PPARD
QG071-06	D05	HQP013617	NM_001001928	PPARA
QG071-06	D06	HQP013473	NM_000446	PON1
QG071-06	D07	HQP013450	NM_018975	TERF2IP
QG071-06	D08	HQP013425	NM_006231	POLE
QG071-06	D09	HQP012484	NM_016299	HSPA14
QG071-06	D10	HQP012393	NM_016009	SH3GLB1
QG071-06	D11	HQP012169	NM_015721	GEMIN4
QG071-06	D12	HQP012155	NM_002575	SERPINB2
QG071-06	E01	HQP012154	NM_000602	SERPINE1
QG071-06	E02	HQP011920	NM_005173	ATP2A3
QG071-06	E03	HQP011914	NM_002524	NRAS
QG071-06	E04	HQP011853	NM_000269	NME1
QG071-06	E05	HQP011800	NM_006164	NFE2L2

Product Data Sheet

QG071-06	E06	HQP011656	NM_002480	PPP1R12A
QG071-06	E07	HQP011559	NM_001018016	MUC1
QG071-06	E08	HQP011523	NM_012331	MSRA
QG071-06	E09	HQP011304	NM_001015052	MPG
QG071-06	E10	HQP011278	NM_002431	MNAT1
QG071-06	E11	HQP011262	NM_002424	MMP8
QG071-06	E12	HQP011245	NM_004529	MLLT3
QG071-06	F01	HQP010138	NM_001172	ARG2
QG071-06	F02	HQP010127	NM_000045	ARG1
QG071-06	F03	HQP010099	NM_000222	KIT
QG071-06	F04	HQP009828	NM_002214	ITGB8
QG071-06	F05	HQP009801	NM_000044	AR
QG071-06	F06	HQP009681	NM_001557	IL8RB
QG071-06	F07	HQP009667	NM_000564	IL5RA
QG071-06	F08	HQP009666	NM_000879	IL5
QG071-06	F09	HQP009643	NM_002182	IL1RAP
QG071-06	F10	HQP009353	NM_002168	IDH2
QG071-06	F11	HQP009351	NM_005896	IDH1
QG071-06	F12	HQP009218	NM_000384	APOB
QG071-06	G01	HQP009091	NM_001167	BIRC4
QG071-06	G02	HQP009089	NM_001540	HSPB1
QG071-06	G03	HQP009086	NM_006597	HSPA8
QG071-06	G04	HQP009036	NM_005343	HRAS
QG071-06	G05	HQP009026	NM_000194	HPRT1
QG071-06	G06	HQP009022	NM_000860	HPGD
QG071-06	G07	HQP008522	NM_014550	CARD10
QG071-06	G08	HQP008502	NM_005316	GTF2H1
QG071-06	G09	HQP008482	NM_000848	GSTM2
QG071-06	G10	HQP008478	NM_000847	GSTA3
QG071-06	G11	HQP008474	NM_000178	GSS
QG071-06	G12	HQP008473	NM_000637	GSR
QG071-06	H01	HGDC		
QG071-06	H02	HGDC		
QG071-06	H03	HQP006940	NM_002046	GAPDH
QG071-06	H04	HQP016381	NM_001101	ACTB
QG071-06	H05	HQP015171	NM_004048	B2M
QG071-06	H06	HQP006171	NM_012423	RPL13A
QG071-06	H07	HQP009026	NM_000194	HPRT1
QG071-06	H08	HQP054253	NR_003286	RN18S1
QG071-06	H09	RT		
QG071-06	H10	RT		
QG071-06	H11	PCR		
QG071-06	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Lung Cancer 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