

ExProfile™ Human Protein Kinases Related Gene qPCR Array

For focused group profiling of human protein kinases genes expression

Cat. No. QG095-A (6 x 96-well plate, Format A)

Cat. No. QG095-B (6 x 96-well plate, Format B)

Cat. No. QG095-C (6 x 96-well plate, Format C)

Cat. No. QG095-D (6 x 96-well plate, Format D)

Cat. No. QG095-E (6 x 96-well plate, Format E)

Plates available individually or as a set of 6. Each set contains 504 unique gene primer pairs deposited in one 96-well plate.

Introduction

The ExProfile human protein kinases related gene qPCR array profiles the expression of 504 human genes related to protein kinase pathway. These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including genes involved in protein kinase pathway which has the ability to reversibly phosphorylate proteins. This array allows researchers to study the related genes to gain understanding of their roles in the functioning and characterization of protein kinase pathway.

- QG095 plate 01: 84 unique gene PCR primer pairs
- QG095 plate 02: 84 unique gene PCR primer pairs
- QG095 plate 03: 84 unique gene PCR primer pairs
- QG095 plate 04: 84 unique gene PCR primer pairs
- QG095 plate 05: 84 unique gene PCR primer pairs
- QG095 plate 06: 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

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	ACVR1C	UHMK1	CSNK1A1L	LRRK2	PIK3AP1	STK11IP	TP53RK	MGC16169	LYK5	MLCK	NEK9	XRCC6BP1
B	ADCK2	MYLK2	MASTL	TTBK1	KIAA1804	BRSK1	FLJ23356	TSSK6	TSSK1B	STK40	NUAK2	TSSK3
C	CDADC1	LRRRC27	ALPK1	ADCK4	LRRK1	FASTKD3	CAMKV	STK33	WNK2	WNK3	WNK4	WNK1
D	ALS2CR7	IQCH	C17orf75	MARK4	TRIB3	TAOK1	ALPK3	SCYL1	CLK4	CAMK1G	SCYL3	CAMK1D
E	CABC1	BCCIP	STK31	PBK	DOK4	SCYL2	BMP2K	CDC42BPG	ALS2CR2	STYK1	STK32B	PANK4
F	UBE2R2	PXK	SNRK	TRPM7	TOLLIP	RIPK4	CSNK1G1	ZAK	RP6-213H1	CRKR8	NLK	PRKAG2
G	TAOK3	CDKL3	CRIM1	VRK3	IRAK4	TNNI3K	TRAT1	PIK3R4	HUNK	SH3KBP1	PACSIN1	NRBP1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG095 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	PKN3	EEF2K	TBK1	HIPK2	STK39	RPS6KA6	STK36	EIF2AK1	RPS6KC1	SRPK3	LATS2	HSPB8
B	RIPK5	PRKD3	NUP62	TSSK2	DAPK2	BRD4	KIAA0999	SNF1LK2	PASK	MAST2	MAP3K7IP2	SMG1
C	TNIK	STK38L	FASTKD2	ICK	LMTK2	AAK1	TWF2	STK38	AKAP13	IRAK3	CHEK2	WIF1
D	MAP4K5	PIM2	RIPK3	FASTK	NEK6	PLK2	MAP3K2	PLK4	CAMKK2	TUSC4	ERN2	STK25
E	MERTK	TESK2	TFG	PAK4	BCKDK	TRIB1	TNK2	HIPK3	COL4A3BP	AKT3	GOLGA5	OXSR1
F	NUAK1	TLK1	MELK	SLK	ULK2	IKBKE	CDC42BPB	ROCK2	EIF2AK3	MAP4K4	TAOK2	STK17A
G	STK17B	RPS6KA5	TBRG4	AURKB	DCLK1	MAP3K13	LATS1	MAP3K14	CDKL2	KALRN	SKAP2	PRPF4B
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG095 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	CPNE3	CDK5R1	KSR1	NRP2	CDKL1	RIOK3	RIPK2	RIPK1	SKAP1	CDC2L5	STK16	PRKRA
B	CASK	MKNK1	CAMK1	IKBK	MAP4K3	CDC42BPA	DOC2B	DYRK2	DYRK3	LTBP4	ULK1	BRD3
C	MAPKAPK3	MAP3K12	YES1	WEE1	VRK2	TYRO3	TYK2	TXK	TTN	TTK	TRIO	TIE1
D	TGFBR3	TGFBR1	TGFB2	TGFA	TESK1	TEK	TEC	MAP3K7	TAF1	SYK	AURKC	STK11
E	STK10	AURKA	STK4	NEK4	SRC	SGK	MAP2K4	CCL4	CCL2	RYK	RPS6KB2	RPS6KB1
F	RPS6KA3	RPS6KA2	RPS6KA1	ROS1	ROCK1	BRD2	RNASEL	TRIM27	RAF1	RAGE	PTK7	PTK6
G	PTK2	PSKH1	LGMN	PRKX	EIF2AK2	MAP2K7	MAP2K6	MAP2K5	MAP2K3	MAP2K2	MAP2K1	MAPK13
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG095 plate 03

	1	2	3	4	5	6	7	8	9	10	11	12
A	MAPK10	MAPK9	MAPK11	MAPK8	MAPK6	MAPK4	MAPK3	MAPK1	PRKG2	PRKG1	PRKCZ	PRKCSH
B	PRKCQ	PRKD1	PKN2	PRKCI	PRKCH	PRKCG	PRKCD	PRKCB1	PRKCA	PRKAR2A	PRKAG1	PRKACG
C	PRKACA	PRKAA2	PRKAA1	PPP4C	EXOSC10	PLK1	PIM1	PHKG2	PHKG1	PHKB	PHKA2	PHKA1
D	PFTK1	PDPK1	PDK4	PDK3	PDK2	PDK1	PDGFRB	PDGFRL	PDGFRA	PCTK2	PCTK1	PAK3
E	PAK2	PAK1	DDR2	ROR2	NTRK3	NTRK1	NPR2	NEK3	NEK2	NEK1	NELL1	MUSK
F	MST1R	MOS	MAP3K11	MAP3K10	MAP3K9	MET	MAP3K4	MAP3K3	MAP3K1	MATK	MAK	LYN
G	LTK	LTBP1	LIMK2	LIMK1	LCK	KIT	KDR	JAK3	JAK2	JAK1	ITK	IRAK2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG095 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	IRAK1	INSRR	INSR	IKBKB	IGF1R	HUS1	HCK	GUCY2D	GUCY2C	GTF2H1	GSK3B	GSK3A
B	MKNK2	GRK5	GRK4	GAK	FYN	FRK	FLT4	FLT3	FLT1	FGR	FGFR4	FGFR2
C	FGFR3	FES	FER	PTK2B	ERN1	ERBB3	ERBB2	EPHB6	EPHB4	EPHB3	EPHB2	EPHA7
D	EPHA4	EPHA3	EPHA1	EPHA2	EGFR	EFNB3	EFNA4	EFNA3	DAPK3	DAPK1	CSNK2A2	CSNK2A1
E	CSNK1G2	CSNK1E	CSNK1D	CSNK1A1	CSK	CSF1R	MAPK14	CRKL	MAP3K8	PLK3	CKS2	CKS1B
F	CHUK	CHEK1	CDKN1A	CDK9	CDK7	CDK6	CDK5	CDK4	CDK2	CDC2	CCNH	CCND3
G	CAMK2D	CAMK2B	CAMK2A	CAMK4	CALM2	CAD	DDR1	BUB1B	BUB1	BTK	BRAF	BMX
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG095 plate 05

	1	2	3	4	5	6	7	8	9	10	11	12
A	BMPR2	BMPR1B	BMPR1A	BLK	BCR	CCND1	AXL	ATR	ATM	ARAF	AMHR2	ALK
B	AKT2	AKT1	ADRBK2	ADRBK1	ACVRL1	ACVR2B	ACVR2A	ACVR1B	ACVR1	ABL1	EPHB1	PRKACB
C	SRPK2	STK24	BRSK2	LOC91481	TPD52L3	OBSCN	GSG2	RIOK2	ABL2	TEX14	CAMK2G	CDC2L1
D	CDK3	CDK8	CDKN2B	CKMT2	CLK1	CLK3	CSNK1G3	CSNK2B	DMPK	FASTKD5	MAP3K6	EPHA5
E	EPHA8	DUSP21	FKBP1A	TLK2	MAP4K1	MARK3	MAP3K5	MAPKAP1	NTRK2	CCRK	PCTK3	PRKCE
F	PKN1	PRKDC	MAPK7	NEK11	SGK3	MAP4K2	RET	YSK4	PRKD2	RIOK1	TRIB2	STK3
G	CDKL5	TGFBR2	VRK1	ZAP70	MAPKAP5	MINK1	TNK1	DYRK4	NRP1	MYO3A	RPS6KA4	ULK4
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure6. Illustration of QG095 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 reversed transcribed from the spike-in control RNA in the sample.
- **PCR:** Positive PCR controls, which are used to verify the PCR efficiency by amplifying the pre-deposited DNA template with its specific pre-deposited primer pairs.

Gene primer list

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG095-01	A01	HQP002611	NM_145259	ACVR1C

QG095-01	A02	HQP002467	NM_175866	UHMK1
QG095-01	A03	HQP002148	NM_145203	CSNK1A1L
QG095-01	A04	HQP002101	NM_198578	LRRK2
QG095-01	A05	HQP002012	NM_152309	PIK3AP1
QG095-01	A06	HQP001700	NM_052902	STK11IP
QG095-01	A07	HQP001522	NM_033550	TP53RK
QG095-01	A08	HQP022604	NM_033115	MGC16169
QG095-01	A09	HQP022363	NM_001003786	LYK5
QG095-01	A10	HQP022258	NM_182493	MLCK
QG095-01	A11	HQP022249	NM_033116	NEK9
QG095-01	A12	HQP022169	NM_033276	XRCC6BP1
QG095-01	B01	HQP022059	NM_052853	ADCK2
QG095-01	B02	HQP021196	NM_033118	MYLK2
QG095-01	B03	HQP021043	NM_032844	MASTL
QG095-01	B04	HQP020831	NM_032538	TTBK1
QG095-01	B05	HQP020737	NM_032435	KIAA1804
QG095-01	B06	HQP020729	NM_032430	BRSK1
QG095-01	B07	HQP020544	NM_032237	FLJ23356
QG095-01	B08	HQP020412	NM_032037	TSSK6
QG095-01	B09	HQP020401	NM_032028	TSSK1B
QG095-01	B10	HQP020391	NM_032017	STK40
QG095-01	B11	HQP019977	NM_030952	NUAK2
QG095-01	B12	HQP019947	NM_052841	TSSK3
QG095-01	C01	HQP019919	NM_030911	CDADC1
QG095-01	C02	HQP019655	NM_030626	LRRC27
QG095-01	C03	HQP019596	NM_025144	ALPK1
QG095-01	C04	HQP019401	NM_024876	ADCK4
QG095-01	C05	HQP019183	NM_024652	LRRK1
QG095-01	C06	HQP018898	NM_024091	FASTKD3
QG095-01	C07	HQP018852	NM_024046	CAMKV
QG095-01	C08	HQP017508	NM_030906	STK33
QG095-01	C09	HQP017387	NM_006648	WNK2
QG095-01	C10	HQP017385	NM_001002838	WNK3
QG095-01	C11	HQP017384	NM_032387	WNK4
QG095-01	C12	HQP017348	NM_018979	WNK1
QG095-01	D01	HQP017313	NM_139158	ALS2CR7
QG095-01	D02	HQP017170	NM_001031715	IQCH
QG095-01	D03	HQP016818	NM_022344	C17orf75
QG095-01	D04	HQP015862	NM_031417	MARK4
QG095-01	D05	HQP015857	NM_021158	TRIB3
QG095-01	D06	HQP015705	NM_020791	TAOK1
QG095-01	D07	HQP015689	NM_020778	ALPK3
QG095-01	D08	HQP015587	NM_001048218	SCYL1

QG095-01	D09	HQP015572	NM_020666	CLK4
QG095-01	D10	HQP015467	NM_020439	CAMK1G
QG095-01	D11	HQP015440	NM_020423	SCYL3
QG095-01	D12	HQP015408	NM_020397	CAMK1D
QG095-01	E01	HQP015320	NM_020247	CABC1
QG095-01	E02	HQP015124	NM_016567	BCCIP
QG095-01	E03	HQP015017	NM_031414	STK31
QG095-01	E04	HQP014794	NM_018492	PBK
QG095-01	E05	HQP014622	NM_018110	DOK4
QG095-01	E06	HQP014588	NM_017988	SCYL2
QG095-01	E07	HQP014489	NM_017593	BMP2K
QG095-01	E08	HQP014471	NM_017525	CDC42BPG
QG095-01	E09	HQP014403	NM_018571	ALS2CR2
QG095-01	E10	HQP014366	NM_018423	STYK1
QG095-01	E11	HQP014360	NM_018401	STK32B
QG095-01	E12	HQP014201	NM_018216	PANK4
QG095-01	F01	HQP013865	NM_017811	UBE2R2
QG095-01	F02	HQP013842	NM_017771	PXK
QG095-01	F03	HQP013788	NM_017719	SNRK
QG095-01	F04	HQP013734	NM_017672	TRPM7
QG095-01	F05	HQP013496	NM_019009	TOLLIP
QG095-01	F06	HQP013385	NM_020639	RIPK4
QG095-01	F07	HQP013348	NM_022048	CSNK1G1
QG095-01	F08	HQP013028	NM_016653	ZAK
QG095-01	F09	HQP013022	NM_001042452	RP6-213H19.1
QG095-01	F10	HQP013013	NM_016507	CRKRS
QG095-01	F11	HQP012972	NM_016231	NLK
QG095-01	F12	HQP012722	NM_001040633	PRKAG2
QG095-01	G01	HQP012666	NM_016281	TAOK3
QG095-01	G02	HQP012572	NM_016508	CDKL3
QG095-01	G03	HQP012543	NM_016441	CRIM1
QG095-01	G04	HQP012541	NM_001025778	VRK3
QG095-01	G05	HQP012441	NM_016123	IRAK4
QG095-01	G06	HQP012375	NM_015978	TNNI3K
QG095-01	G07	HQP012241	NM_016388	TRAT1
QG095-01	G08	HQP008817	NM_014602	PIK3R4
QG095-01	G09	HQP008771	NM_014586	HUNK
QG095-01	G10	HQP008685	NM_001024666	SH3KBP1
QG095-01	G11	HQP008669	NM_020804	PACSIN1
QG095-01	G12	HQP008633	NM_013392	NRBP1
QG095-01	H01	HGDC		
QG095-01	H02	HGDC		
QG095-01	H03	HQP006940	NM_002046	GAPDH

QG095-01	H04	HQP016381	NM_001101	ACTB
QG095-01	H05	HQP015171	NM_004048	B2M
QG095-01	H06	HQP006171	NM_012423	RPL13A
QG095-01	H07	HQP009026	NM_000194	HPRT1
QG095-01	H08	HQP054253	NR_003286	RN18S1
QG095-01	H09	RT		
QG095-01	H10	RT		
QG095-01	H11	PCR		
QG095-01	H12	PCR		
QG095-02	A01	HQP008609	NM_013355	PKN3
QG095-02	A02	HQP008578	NM_013302	EEF2K
QG095-02	A03	HQP008428	NM_013254	TBK1
QG095-02	A04	HQP008358	NM_022740	HIPK2
QG095-02	A05	HQP007693	NM_013233	STK39
QG095-02	A06	HQP007676	NM_014496	RPS6KA6
QG095-02	A07	HQP007549	NM_015690	STK36
QG095-02	A08	HQP007510	NM_014413	EIF2AK1
QG095-02	A09	HQP007372	NM_012424	RPS6KC1
QG095-02	A10	HQP007305	NM_014370	SRPK3
QG095-02	A11	HQP007288	NM_014572	LATS2
QG095-02	A12	HQP007224	NM_014365	HSPB8
QG095-02	B01	HQP006710	NM_015375	RIPK5
QG095-02	B02	HQP006319	NM_005813	PRKD3
QG095-02	B03	HQP006291	NM_012346	NUP62
QG095-02	B04	HQP006274	NM_053006	TSSK2
QG095-02	B05	HQP006265	NM_014326	DAPK2
QG095-02	B06	HQP006135	NM_014299	BRD4
QG095-02	B07	HQP006056	NM_025164	KIAA0999
QG095-02	B08	HQP005897	NM_015191	SNF1LK2
QG095-02	B09	HQP005832	NM_015148	PASK
QG095-02	B10	HQP005791	NM_015112	MAST2
QG095-02	B11	HQP005773	NM_015093	MAP3K7IP2
QG095-02	B12	HQP005703	NM_015092	SMG1
QG095-02	C01	HQP005699	NM_015028	TNIK
QG095-02	C02	HQP005665	NM_015000	STK38L
QG095-02	C03	HQP005522	NM_014929	FASTKD2
QG095-02	C04	HQP005508	NM_014920	ICK
QG095-02	C05	HQP005505	NM_014916	LMTK2
QG095-02	C06	HQP005501	NM_014911	AAK1
QG095-02	C07	HQP001608	NM_007284	TWF2
QG095-02	C08	HQP001580	NM_007271	STK38
QG095-02	C09	HQP001409	NM_006738	AKAP13
QG095-02	C10	HQP001408	NM_007199	IRAK3

QG095-02	C11	HQP001396	NM_001005735	CHEK2
QG095-02	C12	HQP001390	NM_007191	WIF1
QG095-02	D01	HQP001370	NM_006575	MAP4K5
QG095-02	D02	HQP001184	NM_006875	PIM2
QG095-02	D03	HQP001177	NM_006871	RIPK3
QG095-02	D04	HQP001055	NM_006712	FASTK
QG095-02	D05	HQP000920	NM_014397	NEK6
QG095-02	D06	HQP000904	NM_006622	PLK2
QG095-02	D07	HQP000881	NM_006609	MAP3K2
QG095-02	D08	HQP000868	NM_014264	PLK4
QG095-02	D09	HQP000781	NM_006549	CAMKK2
QG095-02	D10	HQP000776	NM_006545	TUSC4
QG095-02	D11	HQP000718	NM_033266	ERN2
QG095-02	D12	HQP000598	NM_006374	STK25
QG095-02	E01	HQP000555	NM_006343	MERTK
QG095-02	E02	HQP000503	NM_007170	TESK2
QG095-02	E03	HQP000419	NM_001007565	TFG
QG095-02	E04	HQP000365	NM_001014831	PAK4
QG095-02	E05	HQP000361	NM_005881	BCKDK
QG095-02	E06	HQP000276	NM_025195	TRIB1
QG095-02	E07	HQP000231	NM_001010938	TNK2
QG095-02	E08	HQP000149	NM_001048200	HIPK3
QG095-02	E09	HQP000113	NM_005713	COL4A3BP
QG095-02	E10	HQP000001	NM_005465	AKT3
QG095-02	E11	HQP023402	NM_005113	GOLGA5
QG095-02	E12	HQP023391	NM_005109	OXSRI
QG095-02	F01	HQP023331	NM_014840	NUAK1
QG095-02	F02	HQP023312	NM_012290	TLK1
QG095-02	F03	HQP023256	NM_014791	MELK
QG095-02	F04	HQP023153	NM_014720	SLK
QG095-02	F05	HQP023102	NM_014683	ULK2
QG095-02	F06	HQP023021	NM_014002	IKBKE
QG095-02	F07	HQP022933	NM_006035	CDC42BPB
QG095-02	F08	HQP022805	NM_004850	ROCK2
QG095-02	F09	HQP022771	NM_004836	EIF2AK3
QG095-02	F10	HQP022766	NM_004834	MAP4K4
QG095-02	F11	HQP022570	NM_004783	TAOK2
QG095-02	F12	HQP022428	NM_004760	STK17A
QG095-02	G01	HQP022427	NM_004226	STK17B
QG095-02	G02	HQP022404	NM_004755	RPS6KA5
QG095-02	G03	HQP022380	NM_004749	TBRG4
QG095-02	G04	HQP022323	NM_004217	AURKB
QG095-02	G05	HQP022304	NM_004734	DCLK1

QG095-02	G06	HQP022250	NM_004721	MAP3K13
QG095-02	G07	HQP022099	NM_004690	LATS1
QG095-02	G08	HQP021888	NM_003954	MAP3K14
QG095-02	G09	HQP021841	NM_003948	CDKL2
QG095-02	G10	HQP021838	NM_001024660	KALRN
QG095-02	G11	HQP021742	NM_003930	SKAP2
QG095-02	G12	HQP021688	NM_003913	PRPF4B
QG095-02	H01	HGDC		
QG095-02	H02	HGDC		
QG095-02	H03	HQP006940	NM_002046	GAPDH
QG095-02	H04	HQP016381	NM_001101	ACTB
QG095-02	H05	HQP015171	NM_004048	B2M
QG095-02	H06	HQP006171	NM_012423	RPL13A
QG095-02	H07	HQP009026	NM_000194	HPRT1
QG095-02	H08	HQP054253	NR_003286	RN18S1
QG095-02	H09	RT		
QG095-02	H10	RT		
QG095-02	H11	PCR		
QG095-02	H12	PCR		
QG095-03	A01	HQP021681	NM_003909	CPNE3
QG095-03	A02	HQP021622	NM_003885	CDK5R1
QG095-03	A03	HQP021615	NM_014238	KSR1
QG095-03	A04	HQP021589	NM_003872	NRP2
QG095-03	A05	HQP021574	NM_004196	CDKL1
QG095-03	A06	HQP021534	NM_003831	RIOK3
QG095-03	A07	HQP021524	NM_003821	RIPK2
QG095-03	A08	HQP021492	NM_003804	RIPK1
QG095-03	A09	HQP021368	NM_001075099	SKAP1
QG095-03	A10	HQP021349	NM_003718	CDC2L5
QG095-03	A11	HQP021306	NM_001008910	STK16
QG095-03	A12	HQP021305	NM_003690	PRKRA
QG095-03	B01	HQP021303	NM_003688	CASK
QG095-03	B02	HQP021299	NM_003684	MKNK1
QG095-03	B03	HQP021198	NM_003656	CAMK1
QG095-03	B04	HQP021140	NM_003639	IKBKG
QG095-03	B05	HQP021033	NM_003618	MAP4K3
QG095-03	B06	HQP020935	NM_003607	CDC42BPA
QG095-03	B07	HQP020755	NM_003585	DOC2B
QG095-03	B08	HQP020743	NM_003583	DYRK2
QG095-03	B09	HQP020734	NM_001004023	DYRK3
QG095-03	B10	HQP020596	NM_001042544	LTBP4
QG095-03	B11	HQP020476	NM_003565	ULK1
QG095-03	B12	HQP019579	NM_007371	BRD3

QG095-03	C01	HQP018812	NM_004635	MAPKAPK3
QG095-03	C02	HQP018750	NM_006301	MAP3K12
QG095-03	C03	HQP018569	NM_005433	YES1
QG095-03	C04	HQP018519	NM_003390	WEE1
QG095-03	C05	HQP018500	NM_006296	VRK2
QG095-03	C06	HQP018346	NM_006293	TYRO3
QG095-03	C07	HQP018340	NM_003331	TYK2
QG095-03	C08	HQP018334	NM_003328	TXK
QG095-03	C09	HQP018289	NM_003319	TTN
QG095-03	C10	HQP018288	NM_003318	TTK
QG095-03	C11	HQP018246	NM_007118	TRIO
QG095-03	C12	HQP018091	NM_005424	TIE1
QG095-03	D01	HQP018055	NM_003243	TGFBR3
QG095-03	D02	HQP018051	NM_004612	TGFBR1
QG095-03	D03	HQP018047	NM_003238	TGFB2
QG095-03	D04	HQP018043	NM_003236	TGFA
QG095-03	D05	HQP018019	NM_006285	TESK1
QG095-03	D06	HQP018012	NM_000459	TEK
QG095-03	D07	HQP018008	NM_003215	TEC
QG095-03	D08	HQP017891	NM_003188	MAP3K7
QG095-03	D09	HQP017873	NM_004606	TAF1
QG095-03	D10	HQP017845	NM_003177	SYK
QG095-03	D11	HQP017795	NM_001015878	AURKC
QG095-03	D12	HQP017794	NM_000455	STK11
QG095-03	E01	HQP017793	NM_005990	STK10
QG095-03	E02	HQP017791	NM_003600	AURKA
QG095-03	E03	HQP017789	NM_006282	STK4
QG095-03	E04	HQP017787	NM_003157	NEK4
QG095-03	E05	HQP017696	NM_005417	SRC
QG095-03	E06	HQP017005	NM_005627	SGK
QG095-03	E07	HQP016830	NM_003010	MAP2K4
QG095-03	E08	HQP016625	NM_002984	CCL4
QG095-03	E09	HQP016621	NM_002982	CCL2
QG095-03	E10	HQP016530	NM_001005861	RYK
QG095-03	E11	HQP016472	NM_001007071	RPS6KB2
QG095-03	E12	HQP016471	NM_003161	RPS6KB1
QG095-03	F01	HQP016470	NM_004586	RPS6KA3
QG095-03	F02	HQP016468	NM_001006932	RPS6KA2
QG095-03	F03	HQP016466	NM_001006665	RPS6KA1
QG095-03	F04	HQP016380	NM_002944	ROS1
QG095-03	F05	HQP016371	NM_005406	ROCK1
QG095-03	F06	HQP016309	NM_005104	BRD2
QG095-03	F07	HQP016303	NM_021133	RNASEL

QG095-03	F08	HQP016233	NM_006510	TRIM27
QG095-03	F09	HQP016088	NM_002880	RAF1
QG095-03	F10	HQP016084	NM_014226	RAGE
QG095-03	F11	HQP015701	NM_002821	PTK7
QG095-03	F12	HQP015692	NM_005975	PTK6
QG095-03	G01	HQP015639	NM_005607	PTK2
QG095-03	G02	HQP015174	NM_006742	PSKH1
QG095-03	G03	HQP015082	NM_001008530	LGMN
QG095-03	G04	HQP014987	NM_005044	PRKX
QG095-03	G05	HQP014948	NM_002759	EIF2AK2
QG095-03	G06	HQP014926	NM_145185	MAP2K7
QG095-03	G07	HQP014919	NM_002758	MAP2K6
QG095-03	G08	HQP014917	NM_002757	MAP2K5
QG095-03	G09	HQP014914	NM_002756	MAP2K3
QG095-03	G10	HQP014909	NM_030662	MAP2K2
QG095-03	G11	HQP014907	NM_002755	MAP2K1
QG095-03	G12	HQP014906	NM_002754	MAPK13
QG095-03	H01	HGDC		
QG095-03	H02	HGDC		
QG095-03	H03	HQP006940	NM_002046	GAPDH
QG095-03	H04	HQP016381	NM_001101	ACTB
QG095-03	H05	HQP015171	NM_004048	B2M
QG095-03	H06	HQP006171	NM_012423	RPL13A
QG095-03	H07	HQP009026	NM_000194	HPRT1
QG095-03	H08	HQP054253	NR_003286	RN18S1
QG095-03	H09	RT		
QG095-03	H10	RT		
QG095-03	H11	PCR		
QG095-03	H12	PCR		
QG095-04	A01	HQP014900	NM_002753	MAPK10
QG095-04	A02	HQP014896	NM_002752	MAPK9
QG095-04	A03	HQP014895	NM_002751	MAPK11
QG095-04	A04	HQP014886	NM_002750	MAPK8
QG095-04	A05	HQP014876	NM_002748	MAPK6
QG095-04	A06	HQP014865	NM_002747	MAPK4
QG095-04	A07	HQP014854	NM_001040056	MAPK3
QG095-04	A08	HQP014848	NM_002745	MAPK1
QG095-04	A09	HQP014847	NM_006259	PRKG2
QG095-04	A10	HQP014844	NM_006258	PRKG1
QG095-04	A11	HQP014829	NM_001033581	PRKCZ
QG095-04	A12	HQP014813	NM_001001329	PRKCSH
QG095-04	B01	HQP014804	NM_006257	PRKCQ
QG095-04	B02	HQP014798	NM_002742	PRKD1

QG095-04	B03	HQP014791	NM_006256	PKN2
QG095-04	B04	HQP014771	NM_002740	PRKCI
QG095-04	B05	HQP014760	NM_006255	PRKCH
QG095-04	B06	HQP014751	NM_002739	PRKCG
QG095-04	B07	HQP014731	NM_006254	PRKCD
QG095-04	B08	HQP014718	NM_002738	PRKCB1
QG095-04	B09	HQP014706	NM_002737	PRKCA
QG095-04	B10	HQP014687	NM_004157	PRKAR2A
QG095-04	B11	HQP014627	NM_002733	PRKAG1
QG095-04	B12	HQP014594	NM_002732	PRKACG
QG095-04	C01	HQP014575	NM_002730	PRKACA
QG095-04	C02	HQP014541	NM_006252	PRKAA2
QG095-04	C03	HQP014530	NM_006251	PRKAA1
QG095-04	C04	HQP014321	NM_002720	PPP4C
QG095-04	C05	HQP013350	NM_001001998	EXOSC10
QG095-04	C06	HQP013262	NM_005030	PLK1
QG095-04	C07	HQP013152	NM_002648	PIM1
QG095-04	C08	HQP013119	NM_000294	PHKG2
QG095-04	C09	HQP013118	NM_006213	PHKG1
QG095-04	C10	HQP013115	NM_000293	PHKB
QG095-04	C11	HQP013114	NM_000292	PHKA2
QG095-04	C12	HQP013113	NM_002637	PHKA1
QG095-04	D01	HQP013082	NM_012395	PFTK1
QG095-04	D02	HQP012979	NM_002613	PDPK1
QG095-04	D03	HQP012953	NM_002612	PDK4
QG095-04	D04	HQP012942	NM_005391	PDK3
QG095-04	D05	HQP012931	NM_002611	PDK2
QG095-04	D06	HQP012923	NM_002610	PDK1
QG095-04	D07	HQP012889	NM_002609	PDGFRB
QG095-04	D08	HQP012871	NM_006207	PDGFRL
QG095-04	D09	HQP012866	NM_006206	PDGFRA
QG095-04	D10	HQP012596	NM_002595	PCTK2
QG095-04	D11	HQP012585	NM_006201	PCTK1
QG095-04	D12	HQP012174	NM_002578	PAK3
QG095-04	E01	HQP012170	NM_002577	PAK2
QG095-04	E02	HQP012156	NM_002576	PAK1
QG095-04	E03	HQP011955	NM_001014796	DDR2
QG095-04	E04	HQP011954	NM_004560	ROR2
QG095-04	E05	HQP011946	NM_001007156	NTRK3
QG095-04	E06	HQP011938	NM_001007792	NTRK1
QG095-04	E07	HQP011902	NM_003995	NPR2
QG095-04	E08	HQP011765	NM_002498	NEK3
QG095-04	E09	HQP011764	NM_002497	NEK2

QG095-04	E10	HQP011763	NM_012224	NEK1
QG095-04	E11	HQP011760	NM_006157	NELL1
QG095-04	E12	HQP011574	NM_005592	MUSK
QG095-04	F01	HQP011525	NM_002447	MST1R
QG095-04	F02	HQP011299	NM_005372	MOS
QG095-04	F03	HQP011240	NM_002419	MAP3K11
QG095-04	F04	HQP011237	NM_002446	MAP3K10
QG095-04	F05	HQP011236	NM_033141	MAP3K9
QG095-04	F06	HQP011181	NM_000245	MET
QG095-04	F07	HQP011166	NM_005922	MAP3K4
QG095-04	F08	HQP011164	NM_002401	MAP3K3
QG095-04	F09	HQP053966	NM_005921	MAP3K1
QG095-04	F10	HQP011048	NM_002378	MATK
QG095-04	F11	HQP010994	NM_005906	MAK
QG095-04	F12	HQP010935	NM_002350	LYN
QG095-04	G01	HQP010920	NM_002344	LTK
QG095-04	G02	HQP010911	NM_000627	LTBP1
QG095-04	G03	HQP010620	NM_001031801	LIMK2
QG095-04	G04	HQP010619	NM_002314	LIMK1
QG095-04	G05	HQP010565	NM_001042771	LCK
QG095-04	G06	HQP010099	NM_000222	KIT
QG095-04	G07	HQP010070	NM_002253	KDR
QG095-04	G08	HQP009851	NM_000215	JAK3
QG095-04	G09	HQP009850	NM_004972	JAK2
QG095-04	G10	HQP009849	NM_002227	JAK1
QG095-04	G11	HQP009835	NM_005546	ITK
QG095-04	G12	HQP009776	NM_001570	IRAK2
QG095-04	H01	HGDC		
QG095-04	H02	HGDC		
QG095-04	H03	HQP006940	NM_002046	GAPDH
QG095-04	H04	HQP016381	NM_001101	ACTB
QG095-04	H05	HQP015171	NM_004048	B2M
QG095-04	H06	HQP006171	NM_012423	RPL13A
QG095-04	H07	HQP009026	NM_000194	HPRT1
QG095-04	H08	HQP054253	NR_003286	RN18S1
QG095-04	H09	RT		
QG095-04	H10	RT		
QG095-04	H11	PCR		
QG095-04	H12	PCR		
QG095-05	A01	HQP009771	NM_001025242	IRAK1
QG095-05	A02	HQP009766	NM_014215	INSRR
QG095-05	A03	HQP009764	NM_000208	INSR
QG095-05	A04	HQP009639	NM_001556	IKBKB

QG095-05	A05	HQP009523	NM_000875	IGF1R
QG095-05	A06	HQP009138	NM_004507	HUS1
QG095-05	A07	HQP008739	NM_002110	HCK
QG095-05	A08	HQP008683	NM_000180	GUCY2D
QG095-05	A09	HQP008552	NM_004963	GUCY2C
QG095-05	A10	HQP008502	NM_005316	GTF2H1
QG095-05	A11	HQP054075	NM_002093	GSK3B
QG095-05	A12	HQP008468	NM_019884	GSK3A
QG095-05	B01	HQP008273	NM_017572	MKNK2
QG095-05	B02	HQP008260	NM_005308	GRK5
QG095-05	B03	HQP008257	NM_001004056	GRK4
QG095-05	B04	HQP006741	NM_005255	GAK
QG095-05	B05	HQP006484	NM_002037	FYN
QG095-05	B06	HQP006377	NM_002031	FRK
QG095-05	B07	HQP005915	NM_002020	FLT4
QG095-05	B08	HQP005890	NM_004119	FLT3
QG095-05	B09	HQP005879	NM_002019	FLT1
QG095-05	B10	HQP005444	NM_001042729	FGR
QG095-05	B11	HQP005439	NM_002011	FGFR4
QG095-05	B12	HQP005437	NM_000141	FGFR2
QG095-05	C01	HQP005434	NM_000142	FGFR3
QG095-05	C02	HQP005395	NM_002005	FES
QG095-05	C03	HQP005394	NM_005246	FER
QG095-05	C04	HQP005093	NM_004103	PTK2B
QG095-05	C05	HQP004992	NM_001433	ERN1
QG095-05	C06	HQP004970	NM_001005915	ERBB3
QG095-05	C07	HQP004968	NM_001005862	ERBB2
QG095-05	C08	HQP004947	NM_004445	EPHB6
QG095-05	C09	HQP004945	NM_004444	EPHB4
QG095-05	C10	HQP004942	NM_004443	EPHB3
QG095-05	C11	HQP004939	NM_004442	EPHB2
QG095-05	C12	HQP004930	NM_004440	EPHA7
QG095-05	D01	HQP004926	NM_004438	EPHA4
QG095-05	D02	HQP004924	NM_005233	EPHA3
QG095-05	D03	HQP004922	NM_005232	EPHA1
QG095-05	D04	HQP004657	NM_004431	EPHA2
QG095-05	D05	HQP004605	NM_005228	EGFR
QG095-05	D06	HQP004598	NM_001406	EFNB3
QG095-05	D07	HQP004592	NM_005227	EFNA4
QG095-05	D08	HQP004591	NM_004952	EFNA3
QG095-05	D09	HQP003982	NM_001348	DAPK3
QG095-05	D10	HQP003979	NM_004938	DAPK1
QG095-05	D11	HQP003285	NM_001896	CSNK2A2

QG095-05	D12	HQP003277	NM_001895	CSNK2A1
QG095-05	E01	HQP003257	NM_001319	CSNK1G2
QG095-05	E02	HQP003249	NM_001894	CSNK1E
QG095-05	E03	HQP003240	NM_001893	CSNK1D
QG095-05	E04	HQP003236	NM_001025105	CSNK1A1
QG095-05	E05	HQP003218	NM_004383	CSK
QG095-05	E06	HQP003158	NM_005211	CSF1R
QG095-05	E07	HQP003133	NM_001315	MAPK14
QG095-05	E08	HQP002989	NM_005207	CRKL
QG095-05	E09	HQP002725	NM_005204	MAP3K8
QG095-05	E10	HQP002379	NM_004073	PLK3
QG095-05	E11	HQP001894	NM_001827	CKS2
QG095-05	E12	HQP001884	NM_001826	CKS1B
QG095-05	F01	HQP001708	NM_001278	CHUK
QG095-05	F02	HQP001282	NM_001274	CHEK1
QG095-05	F03	HQP000331	NM_000389	CDKN1A
QG095-05	F04	HQP000321	NM_001261	CDK9
QG095-05	F05	HQP000285	NM_001799	CDK7
QG095-05	F06	HQP000274	NM_001259	CDK6
QG095-05	F07	HQP000261	NM_004935	CDK5
QG095-05	F08	HQP000245	NM_000075	CDK4
QG095-05	F09	HQP000225	NM_001798	CDK2
QG095-05	F10	HQP023261	NM_001786	CDC2
QG095-05	F11	HQP021906	NM_001239	CCNH
QG095-05	F12	HQP021757	NM_001760	CCND3
QG095-05	G01	HQP019986	NM_001221	CAMK2D
QG095-05	G02	HQP019959	NM_001220	CAMK2B
QG095-05	G03	HQP019917	NM_015981	CAMK2A
QG095-05	G04	HQP019874	NM_001744	CAMK4
QG095-05	G05	HQP019707	NM_001743	CALM2
QG095-05	G06	HQP018928	NM_004341	CAD
QG095-05	G07	HQP018768	NM_001954	DDR1
QG095-05	G08	HQP018022	NM_001211	BUB1B
QG095-05	G09	HQP017999	NM_004336	BUB1
QG095-05	G10	HQP017988	NM_000061	BTK
QG095-05	G11	HQP017733	NM_004333	BRAF
QG095-05	G12	HQP017566	NM_001721	BMX
QG095-05	H01	HGDC		
QG095-05	H02	HGDC		
QG095-05	H03	HQP006940	NM_002046	GAPDH
QG095-05	H04	HQP016381	NM_001101	ACTB
QG095-05	H05	HQP015171	NM_004048	B2M
QG095-05	H06	HQP006171	NM_012423	RPL13A

QG095-05	H07	HQP009026	NM_000194	HPRT1
QG095-05	H08	HQP054253	NR_003286	RN18S1
QG095-05	H09	RT		
QG095-05	H10	RT		
QG095-05	H11	PCR		
QG095-05	H12	PCR		
QG095-06	A01	HQP017535	NM_001204	BMPR2
QG095-06	A02	HQP017500	NM_001203	BMPR1B
QG095-06	A03	HQP017489	NM_004329	BMPR1A
QG095-06	A04	HQP016783	NM_001715	BLK
QG095-06	A05	HQP016418	NM_004327	BCR
QG095-06	A06	HQP016204	NM_053056	CCND1
QG095-06	A07	HQP014815	NM_001699	AXL
QG095-06	A08	HQP013596	NM_001184	ATR
QG095-06	A09	HQP011736	NM_000051	ATM
QG095-06	A10	HQP009832	NM_001654	ARAF
QG095-06	A11	HQP007421	NM_020547	AMHR2
QG095-06	A12	HQP006353	NM_004304	ALK
QG095-06	B01	HQP004995	NM_001626	AKT2
QG095-06	B02	HQP004991	NM_001014431	AKT1
QG095-06	B03	HQP003848	NM_005160	ADRBK2
QG095-06	B04	HQP003816	NM_001619	ADRBK1
QG095-06	B05	HQP022834	NM_000020	ACVRL1
QG095-06	B06	HQP022668	NM_001106	ACVR2B
QG095-06	B07	HQP022491	NM_001616	ACVR2A
QG095-06	B08	HQP022297	NM_004302	ACVR1B
QG095-06	B09	HQP022071	NM_001105	ACVR1
QG095-06	B10	HQP006954	NM_005157	ABL1
QG095-06	B11	HQP004933	NM_004441	EPHB1
QG095-06	B12	HQP014583	NM_002731	PRKACB
QG095-06	C01	HQP017725	NM_182691	SRPK2
QG095-06	C02	HQP020636	NM_001032296	STK24
QG095-06	C03	HQP021896	NM_003957	BRSK2
QG095-06	C04	HQP022182	NM_138370	LOC91461
QG095-06	C05	HQP021807	NM_001001874	TPD52L3
QG095-06	C06	HQP020433	NM_052843	OBSCN
QG095-06	C07	HQP020387	NM_031965	GSG2
QG095-06	C08	HQP014698	NM_018343	RIOK2
QG095-06	C09	HQP007785	NM_005158	ABL2
QG095-06	C10	HQP015008	NM_031272	TEX14
QG095-06	C11	HQP020019	NM_001222	CAMK2G
QG095-06	C12	HQP023275	NM_001787	CDC2L1
QG095-06	D01	HQP000234	NM_001258	CDK3

QG095-06	D02	HQP000312	NM_001260	CDK8
QG095-06	D03	HQP000382	NM_004936	CDKN2B
QG095-06	D04	HQP001845	NM_001825	CKMT2
QG095-06	D05	HQP002051	NM_001024646	CLK1
QG095-06	D06	HQP002068	NM_001292	CLK3
QG095-06	D07	HQP003259	NM_001031812	CSNK1G3
QG095-06	D08	HQP003289	NM_001320	CSNK2B
QG095-06	D09	HQP004379	NM_001081560	DMPK
QG095-06	D10	HQP016322	NM_021826	FASTKD5
QG095-06	D11	HQP022007	NM_004672	MAP3K6
QG095-06	D12	HQP004928	NM_004439	EPHA5
QG095-06	E01	HQP004931	NM_001006943	EPHA8
QG095-06	E02	HQP016685	NM_022076	DUSP21
QG095-06	E03	HQP005472	NM_000801	FKBP1A
QG095-06	E04	HQP001147	NM_006852	TLK2
QG095-06	E05	HQP001371	NM_001042600	MAP4K1
QG095-06	E06	HQP011031	NM_002376	MARK3
QG095-06	E07	HQP011168	NM_005923	MAP3K5
QG095-06	E08	HQP018931	NM_001006617	MAPKAP1
QG095-06	E09	HQP011941	NM_001007097	NTRK2
QG095-06	E10	HQP006208	NM_001039803	CCRK
QG095-06	E11	HQP012609	NM_002596	PCTK3
QG095-06	E12	HQP014742	NM_005400	PRKCE
QG095-06	F01	HQP014782	NM_002741	PKN1
QG095-06	F02	HQP014837	NM_001081640	PRKDC
QG095-06	F03	HQP014877	NM_002749	MAPK7
QG095-06	F04	HQP019328	NM_024800	NEK11
QG095-06	F05	HQP006316	NM_001033578	SGK3
QG095-06	F06	HQP016055	NM_004579	MAP4K2
QG095-06	F07	HQP016220	NM_020630	RET
QG095-06	F08	HQP019524	NM_001018046	YSK4
QG095-06	F09	HQP006801	NM_001079880	PRKD2
QG095-06	F10	HQP020310	NM_031480	RIOK1
QG095-06	F11	HQP008313	NM_021643	TRIB2
QG095-06	F12	HQP017788	NM_006281	STK3
QG095-06	G01	HQP017792	NM_001037343	CDKL5
QG095-06	G02	HQP018053	NM_001024847	TGFBR2
QG095-06	G03	HQP018499	NM_003384	VRK1
QG095-06	G04	HQP018578	NM_207519	ZAP70
QG095-06	G05	HQP021267	NM_003668	MAPKAPK5
QG095-06	G06	HQP012135	NM_001024937	MINK1
QG095-06	G07	HQP021457	NM_003985	TNK1
QG095-06	G08	HQP021558	NM_003845	DYRK4

QG095-06	G09	HQP021595	NM_001024628	NRP1
QG095-06	G10	HQP013335	NM_017433	MYO3A
QG095-06	G11	HQP021799	NM_001006944	RPS6KA4
QG095-06	G12	HQP053901	NM_017886	ULK4
QG095-06	H01	HGDC		
QG095-06	H02	HGDC		
QG095-06	H03	HQP006940	NM_002046	GAPDH
QG095-06	H04	HQP016381	NM_001101	ACTB
QG095-06	H05	HQP015171	NM_004048	B2M
QG095-06	H06	HQP006171	NM_012423	RPL13A
QG095-06	H07	HQP009026	NM_000194	HPRT1
QG095-06	H08	HQP054253	NR_003286	RN18S1
QG095-06	H09	RT		
QG095-06	H10	RT		
QG095-06	H11	PCR		
QG095-06	H12	PCR		

Limited Use License

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

Limited Warranty

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

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

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

© 2016 GeneCopoeia, Inc.

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