

## ExProfile™ Human Proteases Related Gene qPCR Array

For focused group profiling of human proteases genes expression

Cat. No. QG096-A (6 x 96-well plate, Format A)

Cat. No. QG096-B (6 x 96-well plate, Format B)

Cat. No. QG096-C (6 x 96-well plate, Format C)

Cat. No. QG096-D (6 x 96-well plate, Format D)

Cat. No. QG096-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 proteases related gene qPCR array profiles the expression of 504 human genes related to proteases. These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including genes that encode various proteases. This array allows researchers to study the related genes to gain understanding of their roles in the functioning and characterization of proteases.

- QG096 plate 01: 84 unique gene PCR primer pairs
- QG096 plate 02: 84 unique gene PCR primer pairs
- QG096 plate 03: 84 unique gene PCR primer pairs
- QG096 plate 04: 84 unique gene PCR primer pairs
- QG096 plate 05: 84 unique gene PCR primer pairs
- QG096 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
<b>A</b>	TYSDN1	ASTL	DDI1	OVCH1	OVCH2	MPN2	C9orf43	MAP1D	ABHD7	USP12	SENP5	ADAM32
<b>B</b>	HTRA4	UNQ9391	IMMP1L	ADAMTS19	ADAMTS17	ADAMTS16	ADAMTS15	PRSS35	RHBDL3	OTUD7A	USP54	USP51
<b>C</b>	SASP	FLJ32569	PRSS36	PSMA8	TRY1	ACY1L2	CD109	AGBL1	UNQ1887	CPXM2	MMP21	OMA1
<b>D</b>	BIRC8	HTRA3	CAPN13	LONRF1	XRCC6BP1	LMLN	ATG4D	ABHD13	ATG4C	SPPL2A	AGBL4	CNDP1
<b>E</b>	USP32	RHBDL1	USP48	USP42	USP44	IMMP2L	PRSS27	LONP2	NRIP2	ADAMTS10	ADAMTS12	VCPIP1
<b>F</b>	THSD4	AGBL2	SCRN3	MMP28	USP46	TPSB2	DPEP3	OSGEPL1	MMP27	XPNPEP3	ABHD4	ELA2A
<b>G</b>	PAPPA2	SPCS3	AGBL5	SENP2	SCPEP1	ACE2	USP37	USP28	DPP10	USP36	STAMBPL1	NLN
<b>H</b>	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG096 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
<b>A</b>	TMEM27	SENP7	RNPEPL1	CPA6	ADAMTS9	OTUD7B	DPYSL5	NRIP3	CNDP2	OSGEP	TASP1	OTUB1
<b>B</b>	OTUD5	KLK15	YOD1	ABHD10	C4orf20	USP40	USP47	RHBDL2	PGPEP1	ZRANB1	PCSK4	USP53
<b>C</b>	ARTS-1	UCHL5	MBTPS2	C1RL	CPA4	ABHD5	LAP3	KLK14	DPP7	SENP1	USP25	TSP50
<b>D</b>	PYCARD	TMPRSS11E	SPCS1	HTRA2	ADAMDEC1	TINAG	SENP3	KLK13	SENP6	DKFZP586H2123	USP49	BACE2
<b>E</b>	KLK5	QPCT	TMEFF2	BACE1	CASP14	DNPEP	SEC11A	CAPN7	AGTPBP1	PMPCA	ATG4B	METAP1
<b>F</b>	PEG10	USP33	CTRC	USP18	KLK8	SUPT16H	ADAMTS7	CAPN10	CAPN11	PRSS23	ADAMTS5	ADAMTS8
<b>G</b>	ADAM29	ADAM30	METAP2	COPB5	AFG3L2	MMP24	MALT1	USP19	USP20	ADAM28	CLPX	CAPN9
<b>H</b>	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG096 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	MASP2	YME1L1	USP39	CORIN	STAMPB	USP16	PITRM1	WFDC2	PGCP	PRSS16	PSMD14	ELA3A
B	PIGK	NAALADL1	NAALAD2	RCE1	USP3	USP15	USP52	SPCS2	USP34	ADAMTSL2	ECE2	ESPL1
C	KLK4	PREPL	NPEPP8	ADAMT81	ADAMT82	ADAMT83	ADAMT84	NAPSA	LONP1	MMP20	PCSK7	USP8
D	USP10	USP6	USP14	USP13	P11	CFLAR	GGH	ADAM7	ADAM15	ADAM18	ADAM23	CRADD
E	ADAM19	CTSF	MBTPS1	PRSS12	BAP1	USP9Y	USP11	USP5	USP7	XPNPEP2	XPNPEP1	USP1
F	UQCRC2	UQCRC1	USP4	UCHL3	UCHL1	TPSAB1	TPP2	TNFAIP3	TMPRSS2	TLL2	TLL1	THOP1
G	TFRC	TAF2	ADAM17	ST14	SPG7	SOLH	SHH	RNPEP	REN	PZP	PSMD7	PSMB10
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG096 plate 03

	1	2	3	4	5	6	7	8	9	10	11	12
A	PSMB9	PSMB8	PSMB6	PSMB5	PSMB4	PSMB3	PSMB2	PSMB1	PSMA6	PSMA4	PSMA1	PSEN1
B	HTRA1	PRSS8	PRSS7	KLK7	PRSS2	PRSS1	LGMN	PROC	PREP	PLG	PLAU	PLAT
C	SERPINB5	SERPINA1	PHOX	PGC	PEPD	PCSK2	PCSK5	PCSK1	SERPINA5	PAPPA	SERPINB2	SERPINE1
D	FURIN	NSF	NRD1	MMP17	MMP16	MMP15	MMP14	MMP13	MMP12	MMP11	MMP10	MMP9
E	MMP8	MMP7	MMP3	MMP1	MME	MIPEP	MEP1A	ADAM11	LTF	LTA4H	LNPEP	KLKB1
F	KLK1	KEL	IHH	HP	HGF	HABP2	GZMK	GZMB	GZMA	GZMH	PDIA3	FOLH1
G	F12	F11	F10	F2	EPHX1	ENPEP	ENO1	ELA2	ELA1	ECE1	DPP6	DPP4
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG096 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	CFD	ACE	CYLD	CTSS	CTSO	CTSL1	CTSK	CTSH	CTSG	CTSE	CTSD	CTSB
B	CTRB1	CPN1	CPM	CPE	CPD	CPB1	CPA3	CPA1	CMA1	TPP1	CTSC	CASP10
C	CASP9	CASP8	CASP7	CASP6	CASP5	CASP4	CASP3	CASP2	CASP1	CAPN6	CAPN3	CAPN2
D	CAPN1	CAPN5	C3	C2	C1S	C1R	BLMH	CFB	KLK3	BIRC4	BIRC2	APEH
E	ANPEP	AGA	AEBP1	ADAM10	ACY1	ACR	AZU1	COL7A1	F7	FAP	GZMM	KLK2
F	MASP1	RBP3	RHBDL1	TMPRSS11D	KLK11	ADAMT86	CPAMD8	KLK12	RHBDF2	ADAMT20	TMPRSS5	USP26
G	PAPLN	TMPRSS11B	ADAMT14	AMZ1	FREM1	TMPRSS6	PCSK9	KLK9	SERPINA9	TESSP2	TMPRSS7	TMPRSS9
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG096 plate 05

	1	2	3	4	5	6	7	8	9	10	11	12
A	ADAM8	AFG3L1	BMP1	CAPNS1	CPA2	CPB2	CPN2	CTRL	CTSL2	CTSW	CTSZ	DPEP1
B	TMPRSS13	GGT1	HGFAC	HPN	IDE	CFI	MEP1B	MMP2	USP38	MST1	PA2G4	PCSK6
C	CHMP1A	PGA5	CTSA	PRCP	C9orf3	KLK6	PRTN3	PSEN2	PSMA2	PSMA3	PSMA5	PSMA7
D	PSMB7	SPINT1	ADAM12	CLPP	USP9X	MMP23B	CPZ	ADAM21	USP45	ADAM9	PROZ	SCRN2
E	ECEL1	DPP3	USPL1	SEC11C	TPSD1	DPP9	ABHD12	USP21	DHH	CRBN	AMZ2	OTUD6B
F	ADAMTSL1	CPVL	ADAMTSL4	DPP8	PARL	MMP26	TMPRSS4	WFDC5	USP29	PRSS22	TINAGL1	LRAP
G	DPEP2	RHBDF1	MMP25	TMPRSS3	OTUB2	BRCC3	NPEPL1	LONRF3	IMP5	HM13	KAZALD1	APH1B
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure6. Illustration of QG096 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
QG096-01	A01	HQP005152	NM_173555	TYSND1
QG096-01	A02	HQP011260	NM_001002036	ASTL
QG096-01	A03	HQP011040	NM_001001711	DDI1
QG096-01	A04	HQP009339	NM_183378	OVCH1
QG096-01	A05	HQP009338	NM_198185	OVCH2
QG096-01	A06	HQP009253	NM_183062	MPN2
QG096-01	A07	HQP006680	NM_152786	C9orf43
QG096-01	A08	HQP006518	NM_199227	MAP1D
QG096-01	A09	HQP006473	NM_173567	ABHD7
QG096-01	A10	HQP005110	NM_182488	USP12
QG096-01	A11	HQP004954	NM_152699	SENP5
QG096-01	A12	HQP004885	NM_145004	ADAM32
QG096-01	B01	HQP004884	NM_153692	HTRA4
QG096-01	B02	HQP004881	NM_198464	UNQ9391
QG096-01	B03	HQP004619	NM_144981	IMMP1L
QG096-01	B04	HQP004286	NM_133638	ADAMTS19
QG096-01	B05	HQP004273	NM_139057	ADAMTS17
QG096-01	B06	HQP004272	NM_139056	ADAMTS16
QG096-01	B07	HQP004271	NM_139055	ADAMTS15
QG096-01	B08	HQP004199	NM_153362	PRSS35
QG096-01	B09	HQP004021	NM_138328	RHBDL3
QG096-01	B10	HQP003993	NM_130901	OTUD7A
QG096-01	B11	HQP003916	NM_152586	USP54
QG096-01	B12	HQP003903	NM_201286	USP51
QG096-01	C01	HQP003631	NM_152792	SASP
QG096-01	C02	HQP003467	NM_152491	FLJ32569
QG096-01	C03	HQP003313	NM_173502	PRSS36
QG096-01	C04	HQP003141	NM_001025096	PSMA8
QG096-01	C05	HQP002849	NM_001001317	TRY1
QG096-01	C06	HQP002810	NM_001010853	ACY1L2
QG096-01	C07	HQP002808	NM_133493	CD109
QG096-01	C08	HQP002214	NM_152336	AGBL1
QG096-01	C09	HQP002146	NM_139015	UNQ1887
QG096-01	C10	HQP002050	NM_198148	CPXM2
QG096-01	C11	HQP002019	NM_147191	MMP21
QG096-01	C12	HQP001781	NM_145243	OMA1
QG096-01	D01	HQP001440	NM_033341	BIRC8
QG096-01	D02	HQP022682	NM_053044	HTRA3
QG096-01	D03	HQP022353	NM_144575	CAPN13
QG096-01	D04	HQP022234	NM_152271	LONRF1

QG096-01	D05	HQP022169	NM_033276	XRCC6BP1
QG096-01	D06	HQP021775	NM_033029	LMLN
QG096-01	D07	HQP021086	NM_032885	ATG4D
QG096-01	D08	HQP021058	NM_032859	ABHD13
QG096-01	D09	HQP021051	NM_032852	ATG4C
QG096-01	D10	HQP021002	NM_032802	SPPL2A
QG096-01	D11	HQP020987	NM_032785	AGBL4
QG096-01	D12	HQP020923	NM_032649	CNDP1
QG096-01	E01	HQP020865	NM_032582	USP32
QG096-01	E02	HQP020573	NM_032276	RHBDD1
QG096-01	E03	HQP020542	NM_001032730	USP48
QG096-01	E04	HQP020501	NM_032172	USP42
QG096-01	E05	HQP020484	NM_001042403	USP44
QG096-01	E06	HQP020402	NM_032549	IMMP2L
QG096-01	E07	HQP020371	NM_031948	PRSS27
QG096-01	E08	HQP020323	NM_031490	LONP2
QG096-01	E09	HQP020303	NM_031474	NRIP2
QG096-01	E10	HQP019983	NM_030957	ADAMTS10
QG096-01	E11	HQP019981	NM_030955	ADAMTS12
QG096-01	E12	HQP019526	NM_025054	VCPIP1
QG096-01	F01	HQP019346	NM_024817	THSD4
QG096-01	F02	HQP019313	NM_024783	AGBL2
QG096-01	F03	HQP019117	NM_024583	SCRN3
QG096-01	F04	HQP018955	NM_001032278	MMP28
QG096-01	F05	HQP017206	NM_022832	USP46
QG096-01	F06	HQP017018	NM_024164	TPSB2
QG096-01	F07	HQP016839	NM_022357	DPEP3
QG096-01	F08	HQP016833	NM_022353	OSGEPL1
QG096-01	F09	HQP016754	NM_022122	MMP27
QG096-01	F10	HQP016705	NM_022098	XPNPEP3
QG096-01	F11	HQP016665	NM_022060	ABHD4
QG096-01	F12	HQP016571	NM_033440	ELA2A
QG096-01	G01	HQP016354	NM_020318	PAPPA2
QG096-01	G02	HQP016338	NM_021928	SPCS3
QG096-01	G03	HQP016331	NM_001035507	AGBL5
QG096-01	G04	HQP016166	NM_021627	SEN2
QG096-01	G05	HQP016165	NM_021626	SCPEP1
QG096-01	G06	HQP016137	NM_021804	ACE2
QG096-01	G07	HQP015809	NM_020935	USP37
QG096-01	G08	HQP015773	NM_020886	USP28
QG096-01	G09	HQP015762	NM_001004360	DPP10
QG096-01	G10	HQP015746	NM_025090	USP36
QG096-01	G11	HQP015710	NM_020799	STAMBPL1

QG096-01	G12	HQP015643	NM_020726	NLN
QG096-01	H01	HGDC		
QG096-01	H02	HGDC		
QG096-01	H03	HQP006940	NM_002046	GAPDH
QG096-01	H04	HQP016381	NM_001101	ACTB
QG096-01	H05	HQP015171	NM_004048	B2M
QG096-01	H06	HQP006171	NM_012423	RPL13A
QG096-01	H07	HQP009026	NM_000194	HPRT1
QG096-01	H08	HQP054253	NR_003286	RN18S1
QG096-01	H09	RT		
QG096-01	H10	RT		
QG096-01	H11	PCR		
QG096-01	H12	PCR		
QG096-02	A01	HQP015571	NM_020665	TMEM27
QG096-02	A02	HQP015549	NM_001077203	SEN7
QG096-02	A03	HQP015431	NM_018226	RNPEPL1
QG096-02	A04	HQP015385	NM_020361	CPA6
QG096-02	A05	HQP015322	NM_182920	ADAMTS9
QG096-02	A06	HQP015282	NM_020205	OTUD7B
QG096-02	A07	HQP015212	NM_020134	DPYSL5
QG096-02	A08	HQP015151	NM_020645	NRIP3
QG096-02	A09	HQP014659	NM_018235	CNDP2
QG096-02	A10	HQP014544	NM_017807	OSGEP
QG096-02	A11	HQP014516	NM_017714	TASP1
QG096-02	A12	HQP014508	NM_017670	OTUB1
QG096-02	B01	HQP014494	NM_017602	OTUD5
QG096-02	B02	HQP014462	NM_017509	KLK15
QG096-02	B03	HQP014401	NM_018566	YOD1
QG096-02	B04	HQP014354	NM_018394	ABHD10
QG096-02	B05	HQP014328	NM_018359	C4orf20
QG096-02	B06	HQP014204	NM_018218	USP40
QG096-02	B07	HQP013991	NM_017944	USP47
QG096-02	B08	HQP013872	NM_017821	RHBDL2
QG096-02	B09	HQP013781	NM_017712	PGPEP1
QG096-02	B10	HQP013680	NM_017580	ZRANB1
QG096-02	B11	HQP013677	NM_017573	PCSK4
QG096-02	B12	HQP013547	NM_019050	USP53
QG096-02	C01	HQP013009	NM_001040458	ARTS-1
QG096-02	C02	HQP012692	NM_015984	UCHL5
QG096-02	C03	HQP012673	NM_015884	MBTPS2
QG096-02	C04	HQP012584	NM_016546	C1RL
QG096-02	C05	HQP012509	NM_016352	CPA4
QG096-02	C06	HQP012392	NM_016006	ABHD5

QG096-02	C07	HQP012337	NM_015907	LAP3
QG096-02	C08	HQP011329	NM_022046	KLK14
QG096-02	C09	HQP008625	NM_013379	DPP7
QG096-02	C10	HQP008550	NM_014554	SEN1
QG096-02	C11	HQP008516	NM_013396	USP25
QG096-02	C12	HQP008440	NM_013270	TSP50
QG096-02	D01	HQP008424	NM_013258	PYCARD
QG096-02	D02	HQP008345	NM_014058	TMPRSS11E
QG096-02	D03	HQP008333	NM_014041	SPCS1
QG096-02	D04	HQP007710	NM_013247	HTRA2
QG096-02	D05	HQP007649	NM_014479	ADAMDEC1
QG096-02	D06	HQP007631	NM_014464	TINAG
QG096-02	D07	HQP007100	NM_015670	SEN3
QG096-02	D08	HQP007024	NM_015596	KLK13
QG096-02	D09	HQP007008	NM_015571	SEN6
QG096-02	D10	HQP006828	NM_001001991	DKFZP586H2123
QG096-02	D11	HQP006799	NM_018561	USP49
QG096-02	D12	HQP006760	NM_012105	BACE2
QG096-02	E01	HQP006748	NM_001077491	KLK5
QG096-02	E02	HQP006726	NM_012413	QPCT
QG096-02	E03	HQP006312	NM_016192	TMEFF2
QG096-02	E04	HQP006277	NM_012104	BACE1
QG096-02	E05	HQP006233	NM_012114	CASP14
QG096-02	E06	HQP006204	NM_012100	DNPEP
QG096-02	E07	HQP006137	NM_014300	SEC11A
QG096-02	E08	HQP006132	NM_014296	CAPN7
QG096-02	E09	HQP005957	NM_015239	AGTPBP1
QG096-02	E10	HQP005864	NM_015160	PMPCA
QG096-02	E11	HQP005850	NM_013325	ATG4B
QG096-02	E12	HQP005828	NM_015143	METAP1
QG096-02	F01	HQP005745	NM_001040152	PEG10
QG096-02	F02	HQP005686	NM_015017	USP33
QG096-02	F03	HQP001582	NM_007272	CTRC
QG096-02	F04	HQP001496	NM_017414	USP18
QG096-02	F05	HQP001400	NM_007196	KLK8
QG096-02	F06	HQP001391	NM_007192	SUPT16H
QG096-02	F07	HQP001357	NM_014272	ADAMTS7
QG096-02	F08	HQP001299	NM_021251	CAPN10
QG096-02	F09	HQP001298	NM_007058	CAPN11
QG096-02	F10	HQP001260	NM_007173	PRSS23
QG096-02	F11	HQP001258	NM_007038	ADAMTS5
QG096-02	F12	HQP001257	NM_007037	ADAMTS8
QG096-02	G01	HQP001248	NM_014269	ADAM29

QG096-02	G02	HQP001247	NM_021794	ADAM30
QG096-02	G03	HQP001123	NM_006838	METAP2
QG096-02	G04	HQP001122	NM_006837	COPS5
QG096-02	G05	HQP001071	NM_006796	AFG3L2
QG096-02	G06	HQP001019	NM_006690	MMP24
QG096-02	G07	HQP001017	NM_006785	MALT1
QG096-02	G08	HQP000993	NM_006677	USP19
QG096-02	G09	HQP000992	NM_001008563	USP20
QG096-02	G10	HQP000984	NM_014265	ADAM28
QG096-02	G11	HQP000967	NM_006660	CLPX
QG096-02	G12	HQP000888	NM_006615	CAPN9
QG096-02	H01	HGDC		
QG096-02	H02	HGDC		
QG096-02	H03	HQP006940	NM_002046	GAPDH
QG096-02	H04	HQP016381	NM_001101	ACTB
QG096-02	H05	HQP015171	NM_004048	B2M
QG096-02	H06	HQP006171	NM_012423	RPL13A
QG096-02	H07	HQP009026	NM_000194	HPRT1
QG096-02	H08	HQP054253	NR_003286	RN18S1
QG096-02	H09	RT		
QG096-02	H10	RT		
QG096-02	H11	PCR		
QG096-02	H12	PCR		
QG096-03	A01	HQP000882	NM_006610	MASP2
QG096-03	A02	HQP000864	NM_014263	YME1L1
QG096-03	A03	HQP000846	NM_006590	USP39
QG096-03	A04	HQP000841	NM_006587	CORIN
QG096-03	A05	HQP000747	NM_006463	STAMPB
QG096-03	A06	HQP000723	NM_001001992	USP16
QG096-03	A07	HQP000640	NM_014889	PITRM1
QG096-03	A08	HQP000481	NM_006103	WFDC2
QG096-03	A09	HQP000480	NM_016134	PGCP
QG096-03	A10	HQP000341	NM_005865	PRSS16
QG096-03	A11	HQP000265	NM_005805	PSMD14
QG096-03	A12	HQP000168	NM_005747	ELA3A
QG096-03	B01	HQP000045	NM_005482	PIGK
QG096-03	B02	HQP000007	NM_005468	NAALADL1
QG096-03	B03	HQP000006	NM_005467	NAALAD2
QG096-03	B04	HQP023448	NM_001032279	RCE1
QG096-03	B05	HQP023411	NM_006537	USP3
QG096-03	B06	HQP023408	NM_006313	USP15
QG096-03	B07	HQP023370	NM_014871	USP52
QG096-03	B08	HQP023202	NM_014752	SPCS2



QG096-03	B09	HQP023140	NM_014709	USP34
QG096-03	B10	HQP023116	NM_014694	ADAMTSL2
QG096-03	B11	HQP023114	NM_001037324	ECE2
QG096-03	B12	HQP023095	NM_012291	ESPL1
QG096-03	C01	HQP022998	NM_004917	KLK4
QG096-03	C02	HQP022936	NM_001042385	PREPL
QG096-03	C03	HQP022857	NM_006310	NPEPPS
QG096-03	C04	HQP022847	NM_006988	ADAMTS1
QG096-03	C05	HQP022844	NM_014244	ADAMTS2
QG096-03	C06	HQP022843	NM_014243	ADAMTS3
QG096-03	C07	HQP022842	NM_005099	ADAMTS4
QG096-03	C08	HQP022806	NM_004851	NAPSA
QG096-03	C09	HQP022602	NM_004793	LONP1
QG096-03	C10	HQP022517	NM_004771	MMP20
QG096-03	C11	HQP022211	NM_004716	PCSK7
QG096-03	C12	HQP022076	NM_005154	USP8
QG096-03	D01	HQP022072	NM_005153	USP10
QG096-03	D02	HQP022065	NM_004505	USP6
QG096-03	D03	HQP022063	NM_001037334	USP14
QG096-03	D04	HQP021764	NM_003940	USP13
QG096-03	D05	HQP021700	NM_006025	P11
QG096-03	D06	HQP021604	NM_003879	CFLAR
QG096-03	D07	HQP021603	NM_003878	GGH
QG096-03	D08	HQP021517	NM_003817	ADAM7
QG096-03	D09	HQP021509	NM_003815	ADAM15
QG096-03	D10	HQP021507	NM_014237	ADAM18
QG096-03	D11	HQP021504	NM_003812	ADAM23
QG096-03	D12	HQP021493	NM_003805	CRADD
QG096-03	E01	HQP021483	NM_023038	ADAM19
QG096-03	E02	HQP021475	NM_003793	CTSF
QG096-03	E03	HQP021471	NM_003791	MBTPS1
QG096-03	E04	HQP021042	NM_003619	PRSS12
QG096-03	E05	HQP020117	NM_004656	BAP1
QG096-03	E06	HQP020088	NM_004654	USP9Y
QG096-03	E07	HQP020062	NM_004651	USP11
QG096-03	E08	HQP019770	NM_003481	USP5
QG096-03	E09	HQP018818	NM_003470	USP7
QG096-03	E10	HQP018560	NM_003399	XPNPEP2
QG096-03	E11	HQP018559	NM_020383	XPNPEP1
QG096-03	E12	HQP018447	NM_001017415	USP1
QG096-03	F01	HQP018437	NM_003366	UQCRC2
QG096-03	F02	HQP018436	NM_003365	UQCRC1
QG096-03	F03	HQP018429	NM_003363	USP4

QG096-03	F04	HQP018398	NM_006002	UCHL3
QG096-03	F05	HQP018397	NM_004181	UCHL1
QG096-03	F06	HQP018221	NM_003294	TPSAB1
QG096-03	F07	HQP018219	NM_003291	TPP2
QG096-03	F08	HQP018145	NM_006290	TNFAIP3
QG096-03	F09	HQP018136	NM_005656	TMPRSS2
QG096-03	F10	HQP018110	NM_012465	TLL2
QG096-03	F11	HQP018109	NM_012464	TLL1
QG096-03	F12	HQP018073	NM_003249	THOP1
QG096-03	G01	HQP018041	NM_003234	TFRC
QG096-03	G02	HQP017875	NM_003184	TAF2
QG096-03	G03	HQP017866	NM_003183	ADAM17
QG096-03	G04	HQP017759	NM_021978	ST14
QG096-03	G05	HQP017660	NM_003119	SPG7
QG096-03	G06	HQP017620	NM_005632	SOLH
QG096-03	G07	HQP017098	NM_000193	SHH
QG096-03	G08	HQP016333	NM_020216	RNPEP
QG096-03	G09	HQP016215	NM_000537	REN
QG096-03	G10	HQP016039	NM_002864	PZP
QG096-03	G11	HQP015430	NM_002811	PSMD7
QG096-03	G12	HQP015323	NM_002801	PSMB10
QG096-03	H01	HGDC		
QG096-03	H02	HGDC		
QG096-03	H03	HQP006940	NM_002046	GAPDH
QG096-03	H04	HQP016381	NM_001101	ACTB
QG096-03	H05	HQP015171	NM_004048	B2M
QG096-03	H06	HQP006171	NM_012423	RPL13A
QG096-03	H07	HQP009026	NM_000194	HPRT1
QG096-03	H08	HQP054253	NR_003286	RN18S1
QG096-03	H09	RT		
QG096-03	H10	RT		
QG096-03	H11	PCR		
QG096-03	H12	PCR		
QG096-04	A01	HQP015311	NM_002800	PSMB9
QG096-04	A02	HQP015291	NM_004159	PSMB8
QG096-04	A03	HQP015273	NM_002798	PSMB6
QG096-04	A04	HQP015262	NM_002797	PSMB5
QG096-04	A05	HQP015253	NM_002796	PSMB4
QG096-04	A06	HQP015241	NM_002795	PSMB3
QG096-04	A07	HQP015228	NM_002794	PSMB2
QG096-04	A08	HQP015219	NM_002793	PSMB1
QG096-04	A09	HQP015194	NM_002791	PSMA6
QG096-04	A10	HQP015192	NM_002789	PSMA4

QG096-04	A11	HQP015177	NM_002786	PSMA1
QG096-04	A12	HQP015123	NM_000021	PSEN1
QG096-04	B01	HQP015105	NM_002775	HTRA1
QG096-04	B02	HQP015100	NM_002773	PRSS8
QG096-04	B03	HQP015097	NM_002772	PRSS7
QG096-04	B04	HQP015096	NM_005046	KLK7
QG096-04	B05	HQP015084	NM_002770	PRSS2
QG096-04	B06	HQP015083	NM_002769	PRSS1
QG096-04	B07	HQP015082	NM_001008530	LGMN
QG096-04	B08	HQP015041	NM_000312	PROC
QG096-04	B09	HQP014431	NM_002726	PREP
QG096-04	B10	HQP013257	NM_000301	PLG
QG096-04	B11	HQP013204	NM_002658	PLAU
QG096-04	B12	HQP013201	NM_000930	PLAT
QG096-04	C01	HQP013125	NM_002639	SERPINB5
QG096-04	C02	HQP013122	NM_000295	SERPINA1
QG096-04	C03	HQP013108	NM_000444	PHEX
QG096-04	C04	HQP013087	NM_002630	PGC
QG096-04	C05	HQP013045	NM_000285	PEPD
QG096-04	C06	HQP012576	NM_002594	PCSK2
QG096-04	C07	HQP012566	NM_006200	PCSK5
QG096-04	C08	HQP012539	NM_000439	PCSK1
QG096-04	C09	HQP012332	NM_000624	SERPINA5
QG096-04	C10	HQP012189	NM_002581	PAPPA
QG096-04	C11	HQP012155	NM_002575	SERPINB2
QG096-04	C12	HQP012154	NM_000602	SERPINE1
QG096-04	D01	HQP012121	NM_002569	FURIN
QG096-04	D02	HQP011931	NM_006178	NSF
QG096-04	D03	HQP011918	NM_002525	NRD1
QG096-04	D04	HQP011272	NM_016155	MMP17
QG096-04	D05	HQP011270	NM_005941	MMP16
QG096-04	D06	HQP011269	NM_002428	MMP15
QG096-04	D07	HQP011268	NM_004995	MMP14
QG096-04	D08	HQP011267	NM_002427	MMP13
QG096-04	D09	HQP011266	NM_002426	MMP12
QG096-04	D10	HQP011265	NM_005940	MMP11
QG096-04	D11	HQP011264	NM_002425	MMP10
QG096-04	D12	HQP011263	NM_004994	MMP9
QG096-04	E01	HQP011262	NM_002424	MMP8
QG096-04	E02	HQP011258	NM_002423	MMP7
QG096-04	E03	HQP011257	NM_002422	MMP3
QG096-04	E04	HQP011255	NM_002421	MMP1
QG096-04	E05	HQP011254	NM_000902	MME

QG096-04	E06	HQP011222	NM_005932	MIPEP
QG096-04	E07	HQP011177	NM_005588	MEP1A
QG096-04	E08	HQP011129	NM_002390	ADAM11
QG096-04	E09	HQP010919	NM_002343	LTF
QG096-04	E10	HQP010906	NM_000895	LTA4H
QG096-04	E11	HQP010792	NM_005575	LNPEP
QG096-04	E12	HQP010104	NM_000892	KLKB1
QG096-04	F01	HQP010100	NM_002257	KLK1
QG096-04	F02	HQP010071	NM_000420	KEL
QG096-04	F03	HQP009632	NM_002181	IHH
QG096-04	F04	HQP009019	NM_005143	HP
QG096-04	F05	HQP008800	NM_000601	HGF
QG096-04	F06	HQP008711	NM_004132	HABP2
QG096-04	F07	HQP008691	NM_002104	GZMK
QG096-04	F08	HQP008690	NM_004131	GZMB
QG096-04	F09	HQP008689	NM_006144	GZMA
QG096-04	F10	HQP008675	NM_033423	GZMH
QG096-04	F11	HQP008463	NM_005313	PDIA3
QG096-04	F12	HQP006129	NM_001014986	FOLH1
QG096-04	G01	HQP005068	NM_000505	F12
QG096-04	G02	HQP005066	NM_000128	F11
QG096-04	G03	HQP005064	NM_000504	F10
QG096-04	G04	HQP005052	NM_000506	F2
QG096-04	G05	HQP004948	NM_000120	EPHX1
QG096-04	G06	HQP004867	NM_001977	ENPEP
QG096-04	G07	HQP004860	NM_001428	ENO1
QG096-04	G08	HQP004689	NM_001972	ELA2
QG096-04	G09	HQP004688	NM_001971	ELA1
QG096-04	G10	HQP004537	NM_001397	ECE1
QG096-04	G11	HQP004435	NM_001039350	DPP6
QG096-04	G12	HQP004434	NM_001935	DPP4
QG096-04	H01	HGDC		
QG096-04	H02	HGDC		
QG096-04	H03	HQP006940	NM_002046	GAPDH
QG096-04	H04	HQP016381	NM_001101	ACTB
QG096-04	H05	HQP015171	NM_004048	B2M
QG096-04	H06	HQP006171	NM_012423	RPL13A
QG096-04	H07	HQP009026	NM_000194	HPRT1
QG096-04	H08	HQP054253	NR_003286	RN18S1
QG096-04	H09	RT		
QG096-04	H10	RT		
QG096-04	H11	PCR		
QG096-04	H12	PCR		

QG096-05	A01	HQP004198	NM_001928	CFD
QG096-05	A02	HQP004081	NM_000789	ACE
QG096-05	A03	HQP003761	NM_001042355	CYLD
QG096-05	A04	HQP003668	NM_004079	CTSS
QG096-05	A05	HQP003658	NM_001334	CTSO
QG096-05	A06	HQP003630	NM_001912	CTSL1
QG096-05	A07	HQP003626	NM_000396	CTSK
QG096-05	A08	HQP003620	NM_004390	CTSH
QG096-05	A09	HQP003612	NM_001911	CTSG
QG096-05	A10	HQP003605	NM_001910	CTSE
QG096-05	A11	HQP003599	NM_001909	CTSD
QG096-05	A12	HQP003596	NM_001908	CTSB
QG096-05	B01	HQP003577	NM_001906	CTRB1
QG096-05	B02	HQP002857	NM_001308	CPN1
QG096-05	B03	HQP002855	NM_001005502	CPM
QG096-05	B04	HQP002847	NM_001873	CPE
QG096-05	B05	HQP002842	NM_001304	CPD
QG096-05	B06	HQP002835	NM_001871	CPB1
QG096-05	B07	HQP002832	NM_001870	CPA3
QG096-05	B08	HQP002821	NM_001868	CPA1
QG096-05	B09	HQP002142	NM_001836	CMA1
QG096-05	B10	HQP002078	NM_000391	TPP1
QG096-05	B11	HQP000894	NM_001814	CTSC
QG096-05	B12	HQP020709	NM_001230	CASP10
QG096-05	C01	HQP020648	NM_001229	CASP9
QG096-05	C02	HQP018966	NM_001080124	CASP8
QG096-05	C03	HQP020480	NM_001227	CASP7
QG096-05	C04	HQP020427	NM_001226	CASP6
QG096-05	C05	HQP020384	NM_004347	CASP5
QG096-05	C06	HQP020332	NM_001225	CASP4
QG096-05	C07	HQP020297	NM_004346	CASP3
QG096-05	C08	HQP020249	NM_032982	CASP2
QG096-05	C09	HQP020207	NM_001223	CASP1
QG096-05	C10	HQP020086	NM_014289	CAPN6
QG096-05	C11	HQP020071	NM_000070	CAPN3
QG096-05	C12	HQP020070	NM_001748	CAPN2
QG096-05	D01	HQP020065	NM_005186	CAPN1
QG096-05	D02	HQP018286	NM_004055	CAPN5
QG096-05	D03	HQP018238	NM_000064	C3
QG096-05	D04	HQP018226	NM_000063	C2
QG096-05	D05	HQP018206	NM_001734	C1S
QG096-05	D06	HQP018179	NM_001733	C1R
QG096-05	D07	HQP016897	NM_000386	BLMH

QG096-05	D08	HQP016565	NM_001710	CFB
QG096-05	D09	HQP009633	NM_001030047	KLK3
QG096-05	D10	HQP009091	NM_001167	BIRC4
QG096-05	D11	HQP009072	NM_001166	BIRC2
QG096-05	D12	HQP009056	NM_001640	APEH
QG096-05	E01	HQP008414	NM_001150	ANPEP
QG096-05	E02	HQP004378	NM_000027	AGA
QG096-05	E03	HQP004154	NM_001129	AEBP1
QG096-05	E04	HQP000372	NM_001110	ADAM10
QG096-05	E05	HQP022963	NM_000666	ACY1
QG096-05	E06	HQP012072	NM_001097	ACR
QG096-05	E07	HQP015156	NM_001700	AZU1
QG096-05	E08	HQP002566	NM_000094	COL7A1
QG096-05	E09	HQP005059	NM_000131	F7
QG096-05	E10	HQP005101	NM_004460	FAP
QG096-05	E11	HQP008692	NM_005317	GZMM
QG096-05	E12	HQP010101	NM_001002231	KLK2
QG096-05	F01	HQP015091	NM_001031849	MASP1
QG096-05	F02	HQP016196	NM_002900	RBP3
QG096-05	F03	HQP021904	NM_003961	RHBDL1
QG096-05	F04	HQP022693	NM_004262	TMPRSS11D
QG096-05	F05	HQP001148	NM_006853	KLK11
QG096-05	F06	HQP001358	NM_197941	ADAMTS6
QG096-05	F07	HQP007550	NM_015692	CPAMD8
QG096-05	F08	HQP011330	NM_019598	KLK12
QG096-05	F09	HQP019131	NM_001005498	RHBDF2
QG096-05	F10	HQP019504	NM_025003	ADAMTS20
QG096-05	F11	HQP019815	NM_030770	TMPRSS5
QG096-05	F12	HQP020337	NM_031907	USP26
QG096-05	G01	HQP021825	NM_173462	PAPLN
QG096-05	G02	HQP002727	NM_182502	TMPRSS11B
QG096-05	G03	HQP003053	NM_080722	ADAMTS14
QG096-05	G04	HQP003800	NM_133463	AMZ1
QG096-05	G05	HQP003867	NM_144966	FREM1
QG096-05	G06	HQP004119	NM_153609	TMPRSS6
QG096-05	G07	HQP006606	NM_174936	PCSK9
QG096-05	G08	HQP008030	NM_012315	KLK9
QG096-05	G09	HQP009051	NM_001042518	SERPINA9
QG096-05	G10	HQP009274	NM_182702	TESSP2
QG096-05	G11	HQP009439	NM_001042575	TMPRSS7
QG096-05	G12	HQP009712	NM_182973	TMPRSS9
QG096-05	H01	HGDC		
QG096-05	H02	HGDC		

QG096-05	H03	HQP006940	NM_002046	GAPDH
QG096-05	H04	HQP016381	NM_001101	ACTB
QG096-05	H05	HQP015171	NM_004048	B2M
QG096-05	H06	HQP006171	NM_012423	RPL13A
QG096-05	H07	HQP009026	NM_000194	HPRT1
QG096-05	H08	HQP054253	NR_003286	RN18S1
QG096-05	H09	RT		
QG096-05	H10	RT		
QG096-05	H11	PCR		
QG096-05	H12	PCR		
QG096-06	A01	HQP000246	NM_001109	ADAM8
QG096-06	A02	HQP004322	NM_001031805	AFG3L1
QG096-06	A03	HQP017280	NM_001199	BMP1
QG096-06	A04	HQP020082	NM_001003962	CAPNS1
QG096-06	A05	HQP002824	NM_001869	CPA2
QG096-06	A06	HQP002836	NM_001872	CPB2
QG096-06	A07	HQP002860	NM_001080513	CPN2
QG096-06	A08	HQP003587	NM_001907	CTRL
QG096-06	A09	HQP003635	NM_001333	CTSL2
QG096-06	A10	HQP003677	NM_001335	CTSW
QG096-06	A11	HQP003680	NM_001336	CTSZ
QG096-06	A12	HQP004430	NM_004413	DPEP1
QG096-06	B01	HQP020429	NM_001077263	TMPRSS13
QG096-06	B02	HQP007378	NM_001032364	GGT1
QG096-06	B03	HQP008811	NM_001528	HGFAC
QG096-06	B04	HQP009023	NM_002151	HPN
QG096-06	B05	HQP009349	NM_004969	IDE
QG096-06	B06	HQP009376	NM_000204	CFI
QG096-06	B07	HQP011178	NM_005925	MEP1B
QG096-06	B08	HQP054072	NM_004530	MMP2
QG096-06	B09	HQP020839	NM_032557	USP38
QG096-06	B10	HQP011524	NM_020998	MST1
QG096-06	B11	HQP012112	NM_006191	PA2G4
QG096-06	B12	HQP012122	NM_002570	PCSK6
QG096-06	C01	HQP012507	NM_001083314	CHMP1A
QG096-06	C02	HQP013084	NM_014224	PGA5
QG096-06	C03	HQP013686	NM_000308	CTSA
QG096-06	C04	HQP014413	NM_005040	PRCP
QG096-06	C05	HQP021022	NM_032823	C9orf3
QG096-06	C06	HQP015101	NM_001012964	KLK6
QG096-06	C07	HQP015107	NM_002777	PRTN3
QG096-06	C08	HQP015130	NM_000447	PSEN2
QG096-06	C09	HQP015182	NM_002787	PSMA2

QG096-06	C10	HQP015185	NM_002788	PSMA3
QG096-06	C11	HQP015193	NM_002790	PSMA5
QG096-06	C12	HQP015201	NM_002792	PSMA7
QG096-06	D01	HQP015285	NM_002799	PSMB7
QG096-06	D02	HQP017668	NM_001032367	SPINT1
QG096-06	D03	HQP019702	NM_003474	ADAM12
QG096-06	D04	HQP020030	NM_006012	CLPP
QG096-06	D05	HQP020063	NM_001039590	USP9X
QG096-06	D06	HQP021134	NM_006983	MMP23B
QG096-06	D07	HQP021182	NM_001014447	CPZ
QG096-06	D08	HQP021505	NM_003813	ADAM21
QG096-06	D09	HQP021117	NM_001080481	USP45
QG096-06	D10	HQP021515	NM_001005845	ADAM9
QG096-06	D11	HQP021634	NM_003891	PROZ
QG096-06	D12	HQP021964	NM_138355	SCRN2
QG096-06	E01	HQP022744	NM_004826	ECEL1
QG096-06	E02	HQP000094	NM_005700	DPP3
QG096-06	E03	HQP000259	NM_005800	USPL1
QG096-06	E04	HQP022016	NM_033280	SEC11C
QG096-06	E05	HQP006097	NM_012217	TPSD1
QG096-06	E06	HQP022077	NM_139159	DPP9
QG096-06	E07	HQP007029	NM_001042472	ABHD12
QG096-06	E08	HQP007428	NM_001014443	USP21
QG096-06	E09	HQP012238	NM_021044	DHH
QG096-06	E10	HQP012486	NM_016302	CRBN
QG096-06	E11	HQP012637	NM_001033569	AMZ2
QG096-06	E12	HQP012918	NM_016023	OTUD6B
QG096-06	F01	HQP022481	NM_052866	ADAMTSL1
QG096-06	F02	HQP013523	NM_019029	CPVL
QG096-06	F03	HQP013525	NM_019032	ADAMTSL4
QG096-06	F04	HQP013810	NM_017743	DPP8
QG096-06	F05	HQP014415	NM_001037639	PARL
QG096-06	F06	HQP015103	NM_021801	MMP26
QG096-06	F07	HQP015128	NM_019894	TMPRSS4
QG096-06	F08	HQP003527	NM_145652	WFDC5
QG096-06	F09	HQP015789	NM_020903	USP29
QG096-06	F10	HQP016751	NM_022119	PRSS22
QG096-06	F11	HQP016802	NM_022164	TINAGL1
QG096-06	F12	HQP016828	NM_022350	LRAP
QG096-06	G01	HQP016837	NM_022355	DPEP2
QG096-06	G02	HQP016887	NM_022450	RHBDF1
QG096-06	G03	HQP016947	NM_004142	MMP25
QG096-06	G04	HQP017095	NM_024022	TMPRSS3



QG096-06	G05	HQP018831	NM_023112	OTUB2
QG096-06	G06	HQP018987	NM_001018055	BRCC3
QG096-06	G07	HQP019193	NM_024663	NPEPL1
QG096-06	G08	HQP019307	NM_001031855	LONRF3
QG096-06	G09	HQP004025	NM_175882	IMP5
QG096-06	G10	HQP019876	NM_030789	HM13
QG096-06	G11	HQP019939	NM_030929	KAZALD1
QG096-06	G12	HQP020190	NM_031301	APH1B
QG096-06	H01	HGDC		
QG096-06	H02	HGDC		
QG096-06	H03	HQP006940	NM_002046	GAPDH
QG096-06	H04	HQP016381	NM_001101	ACTB
QG096-06	H05	HQP015171	NM_004048	B2M
QG096-06	H06	HQP006171	NM_012423	RPL13A
QG096-06	H07	HQP009026	NM_000194	HPRT1
QG096-06	H08	HQP054253	NR_003286	RN18S1
QG096-06	H09	RT		
QG096-06	H10	RT		
QG096-06	H11	PCR		
QG096-06	H12	PCR		

### Limited Use License

Following terms and conditions apply to use of ExProfile™ Human Proteases 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](mailto:inquiry@genecopoeia.com)