

ExProfile™ Human Breast Cancer Gene qPCR Array

For focused group profiling of human breast cancer related genes expression

Cat. No. QG063-A (6 x 96-well plate, Format A)

Cat. No. QG063-B (6 x 96-well plate, Format B)

Cat. No. QG063-C (6 x 96-well plate, Format C)

Cat. No. QG063-D (6 x 96-well plate, Format D)

Cat. No. QG063-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 breast cancer gene qPCR array profiles 504 human genes to aberrantly expressed human genes involved in human breast 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 breast cancer gene array allows researchers to study the cancer-related genes to gain understanding of their roles in breast cancer pathogenesis.

- QG063 plate 01: 84 unique gene PCR primer pairs
- QG063 plate 02: 84 unique gene PCR primer pairs
- QG063 plate 03: 84 unique gene PCR primer pairs
- QG063 plate 04: 84 unique gene PCR primer pairs
- QG063 plate 05: 84 unique gene PCR primer pairs
- QG063 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

GeneCopia provides five qPCR array formats (A, B, C, D, and E) suitable for use with the following real-time cyclers.

Important note: Upon receiving, please check to make sure that the correct array format was ordered to ensure the compatibility with your qPCR instrument.

Plate format	Instrument provider	qPCR instrument model
A (96-well)	Applied Biosystems	5700, 7000, 7300, 7500, 7700, 7900HT (Standard 96-well block), ViiA™7 (Standard 96-well block)
B (96-well)	Applied Biosystems	7500 (Fast block), 7900HT (Fast block), StepOnePlus™, ViiA™7 (Fast block)
C (96-well)	Bio-Rad Laboratories	iCycler iQ®, MyiQ™, iQ™5
D (96-well)	Bio-Rad Laboratories	CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™
E (96-well)	Roche Applied Science	LightCycler® 480 (96-well block)

Quality control

1. Each pair of primers in the ExProfile gene qPCR array has been experimentally validated to yield a single dissociation curve peak and to generate a single amplicon of the correct size for the targeted gene.
2. The positive PCR controls (PCR) have been verified to amplify a single amplicon of the correct size with Ct values around **20±2**.
3. The Spike-in reverse transcription controls (RT) have been verified to amplify a single amplicon of the correct size with Ct values around **20±3**.
4. $R^2 > 0.99$ was observed for high inter/ intra-array reproducibility.

Materials required but not provided

All-in-One™ First-Strand cDNA Synthesis Kit
 All-in-One™ qPCR Mix
 Total RNA extraction kit (RNAzol® RT RNA extraction reagent is recommended)
 DNase/RNase free tips, PCR reaction tubes, 1.5 ml microcentrifuge tubes
 5 ml and 10 ml graduated pipettes, beakers, flasks, and cylinders
 10 µl to 1,000 µl adjustable single channel micropipettes with disposable tips
 5 µl to 20 µl adjustable multichannel micropipette, disposable tips, and reservoir
 qPCR instrument, compatible with gene qPCR arrays ordered

Array layout

	1	2	3	4	5	6	7	8	9	10	11	12
A	BRCA2	TP53	CHEK2	XRCC1	CYP17A1	CYP1B1	ESR1	XRCC3	COMT	SULT1A1	CYP1A1	CYP19A1
B	SOD2	TGFB1	VDR	CYP2D6	MTHFR	ERCC2	NAT2	FGFR2	NBN	TNF	IGF1	MDM2
C	AR	ESR2	PGR	CCND1	VEGFA	ABCB1	XRCC2	CASP8	IGFBP3	PIK3CA	ERCC4	MAP3K1
D	PTGS2	CYP3A5	HSD17B1	BARD1	NQO1	LIG4	NOS3	XPC	EGFR	IL6	TYMS	CDKN2A
E	ERCC5	IL10	LEPR	LSP1	NCOA3	CDH1	CYP2C19	AHR	MMP2	MPO	MTR	SERPINE1
F	RAD52	RAD50	ADH1B	CYP2C9	IRS1	SHBG	SRD5A2	TGFBR1	XRCC4	NAT1	ADH1C	CYP11A1
G	EPHX1	APEX1	IGFBP1	IL1B	ITGB3	MLH1	OGG1	PHB	ZNF350	SIPA1	NR1H2	CDK4
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG063 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	CCNE1	CCND2	XRCC5	TXNRD1	TXN	SOD1	CXCL12	RAD51C	BACH1	PPARG	ATR	MRE11A
B	MMP3	MMP1	LEP	FASLG	IGF1R	GH1	TXN2	XRCC6	ERCC6	EP300	CYP2B6	PARP1
C	TXNRD2	CDKN2D	IGFALS	HLA-DRB1	GSTO1	CBR3	WRN	UGT2B7	SULT1E1	BLM	BCL2	RB1
D	PPARBP	GHRL	NME1	MYC	MMP9	LIG1	KDR	IL8	IL1A	APOE	APC	MSH6
E	GSTM3	GSR	GPX4	FANCD2	ERCC1	CTLA4	CHEK1	CDKN2C	CDK2	DNMT3B	SST	UGT2B15
F	CLCA2	NCOR1	ADIPOQ	PERLD1	PTTG1	MBD2	CBS	CBR1	AKR1C3	PPP1R1B	BAP1	COL18A1
G	TTK	TOP2A	BUB1B	TERF2	TERF1	TEP1	SSTR2	SKP2	SHC1	RNASEL	RAD51L1	RAD23B
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG063 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	PTPRJ	PTGS1	PTGIS	PTGDS	PTEN	PRL	ECHDC1	FBXW7	KIAA1794	PON1	POLD1	ADIPOR1
B	NPAS2	NQO2	NFKBIA	MUC1	MTRR	STS	LIG3	ITGB4	IL1RN	KLK3	HSD17B2	HRAS
C	ACACA	HIF1A	HFE	HDAC2	GSTZ1	GSTM2	GSTA2	GNB3	GLO1	GHRH	GC	POT1
D	IKZF3	FGFR4	ALDH3A1	FANCF	FANCC	ALDH1A1	F5	F2	ERCC3	EGF	DNASE1	DMBT1
E	DHFR	ACE	ADRB2	CTNINB1	CREBBP	CCR5	AKAP13	RASSF1	PPARGC1A	CARM1	CDK6	TP73
F	NEUROD2	LUM	GHSR	GHRHR	ABCC4	FGFR1	EXO1	SLC6A4	DMTF1	CDC34	CDC25A	CDC20
G	SETDB1	EPM2AIP1	ELMO1	CDC2	SPOCK2	ESPL1	KLK4	SLC4A7	TBX4	ROCK2	C1orf38	CHST3
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG063 plate 03

Product Data Sheet

	1	2	3	4	5	6	7	8	9	10	11	12
A	NRXN2	KL	ITGB1BP1	CD4	BUB3	NMI	VNN2	PCAF	NRP1	NRP2	CES2	TNFRSF10A
B	TNFRSF10B	TNFSF10	RIPK1	TNKS	CAV1	CNTNAP4	LMO4	CNTNAP1	FBXL20	KLF11	RECK	CCDC98
C	AXIN2	AXIN1	NRIP1	CDT1	DPF3	CLPTM1L	ACAD10	TNKS2	COASY	CALCR	EHMT1	SHFM1
D	VTCN1	ADIPOR2	CXCR4	PRDM2	WEE1	VHL	USP1	UQCRRF31	UCP2	UBE2I	TWIST1	TSHR
E	TP53BP1	TNFRSF1B	TLR3	TGFB2	TFR2	BUB1	TBXAS1	TBX2	TBP	ADAM17	STAT3	STAT1
F	SSTR5	SSTR3	SREBF1	SRD5A1	SMARCA2	SLC20A1	SLC19A1	NOD2	BHMT	CCL1	RPS6KA1	MRPL23
G	RPA3	RPA2	RPA1	ACTB	REV3L	RELA	RBBP8	RAD17	RAD1	PTPRN2	LSM2	CALCOCO1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG063 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	MYH7B	MTUS1	MTA3	PTHLH	PTCH1	BCCIP	TEX14	MAPK9	MAPK8	PRKCA	IL17RB	PPP2R5E
B	PPP2R2A	PPP2R1B	ZWILCH	PPP1CB	ATRX	PPARA	UGT1A6	TERF2IP	POLB	41521	TLR9	PNMT
C	PLK1	PIK3CB	PGC	CRKRS	UIMC1	HSD17B7	REV1	TUBD1	SLC45A2	HSD17B12	PCNA	SH3GLB1
D	MLXIPL	GMNN	ODC1	NFKB1	NEK2	NCF4	NCF2	MYT1	MSH3	MSH2	MPG	MMP12
E	KITLG	MDM4	MBP	MAT2A	SMAD7	LRP1	LPL	RHOC	YPEL2	KPNB1	KLKB1	KISS1
F	JUN	ITGB5	ITGAV	ITGA6	INSR	INS	IL18	IL13	IL12RB2	IL4	IGHMMP2	IGFBP7
G	IGFBP5	IGFBP2	IGF2	HUS1	HTR3A	HSD11B1	HSD3B1	PRMT1	HP	FOXA1	HMMR	HLA-A
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG063 plate 05

	1	2	3	4	5	6	7	8	9	10	11	12
A	HGF	SLC40A1	GZMB	GTF2E1	GTF2A1	GSK3B	GRIK1	GPC1	GP1BA	GNRHR	GNAS	GLG1
B	PELP1	STEAP1	AATF	EHF	RAD54B	ALOX5AP	PLD3	FN1	RRP1B	CD93	PPM1E	RRAS2
C	FEN1	FCGR2B	FCGR2A	FCER1A	FBN1	FASN	ALDH2	FANCA	ESRRG	AKT1	ABCA1	E2F2
D	AGT	DRD2	ADAMTS18	HTR3C	DHX9	DCN	CYP24A1	CYP2E1	ADRB3	CYP2C8	BTLA	IL23R
E	CSTF1	CSNK1E	CSNK1D	CWF19L2	CRY2	JMY	GIPC3	ADH1A	CCR7	MMP21	AKAP10	IRAK3
F	YWHAQ	STIP1	EHMT2	MAGED2	PPP1R13L	ALDH1L1	CTCF	POSTN	CENPE	HOXB13	MAD2L2	TOMM40
G	EMG1	CDK7	AKAP9	MUC6	DIRAS3	SHARPIN	TPMT	SSTR4	SSTR1	KLK10	POU1F1	NOS1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure6. Illustration of QG063 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
QG063-01	A01	HQP017753	NM_000059	BRCA2
QG063-01	A02	HQP018175	NM_000546	TP53
QG063-01	A03	HQP001396	NM_001005735	CHEK2
QG063-01	A04	HQP018562	NM_006297	XRCC1
QG063-01	A05	HQP003888	NM_000102	CYP17A1

Product Data Sheet

QG063-01	A06	HQP003775	NM_000104	CYP1B1
QG063-01	A07	HQP004998	NM_000125	ESR1
QG063-01	A08	HQP018564	NM_005432	XRCC3
QG063-01	A09	HQP002671	NM_000754	COMT
QG063-01	A10	HQP017811	NM_177536	SULT1A1
QG063-01	A11	HQP003772	NM_000499	CYP1A1
QG063-01	A12	HQP003904	NM_000103	CYP19A1
QG063-01	B01	HQP017616	NM_000636	SOD2
QG063-01	B02	HQP018044	NM_000660	TGFB1
QG063-01	B03	HQP018474	NM_000376	VDR
QG063-01	B04	HQP003814	NM_000106	CYP2D6
QG063-01	B05	HQP011547	NM_005957	MTHFR
QG063-01	B06	HQP004976	NM_000400	ERCC2
QG063-01	B07	HQP001136	NM_000015	NAT2
QG063-01	B08	HQP054054	BC039243	FGFR2
QG063-01	B09	HQP011687	NM_002485	NBN
QG063-01	B10	HQP018141	NM_000594	TNF
QG063-01	B11	HQP009518	NM_000618	IGF1
QG063-01	B12	HQP011135	NM_002392	MDM2
QG063-01	C01	HQP009801	NM_000044	AR
QG063-01	C02	HQP005002	NM_001437	ESR2
QG063-01	C03	HQP013099	NM_000926	PGR
QG063-01	C04	HQP016204	NM_053056	CCND1
QG063-01	C05	HQP018475	NM_001025366	VEGFA
QG063-01	C06	HQP013100	NM_000927	ABCB1
QG063-01	C07	HQP018563	NM_005431	XRCC2
QG063-01	C08	HQP018966	NM_001080124	CASP8
QG063-01	C09	HQP009544	NM_000598	IGFBP3
QG063-01	C10	HQP013150	NM_006218	PIK3CA
QG063-01	C11	HQP004984	NM_005236	ERCC4
QG063-01	C12	HQP053966	NM_005921	MAP3K1
QG063-01	D01	HQP015598	NM_000963	PTGS2
QG063-01	D02	HQP003841	NM_000777	CYP3A5
QG063-01	D03	HQP009064	NM_000413	HSD17B1
QG063-01	D04	HQP015946	NM_000465	BARD1
QG063-01	D05	HQP004317	NM_000903	NQO1
QG063-01	D06	HQP010613	NM_002312	LIG4
QG063-01	D07	HQP011868	NM_000603	NOS3
QG063-01	D08	HQP018556	NM_004628	XPC
QG063-01	D09	HQP004605	NM_005228	EGFR
QG063-01	D10	HQP009670	NM_000600	IL6
QG063-01	D11	HQP018342	NM_001071	TYMS
QG063-01	D12	HQP000370	NM_058195	CDKN2A
QG063-01	E01	HQP004985	NM_000123	ERCC5
QG063-01	E02	HQP009685	NM_000572	IL10
QG063-01	E03	HQP010584	NM_002303	LEPR
QG063-01	E04	HQP010892	NM_001013253	LSP1
QG063-01	E05	HQP020040	NM_006534	NCOA3

Product Data Sheet

QG063-01	E06	HQP023466	NM_004360	CDH1
QG063-01	E07	HQP003809	NM_000769	CYP2C19
QG063-01	E08	HQP004658	NM_001621	AHR
QG063-01	E09	HQP011256	NM_004530	MMP2
QG063-01	E10	HQP011309	NM_000250	MPO
QG063-01	E11	HQP011554	NM_000254	MTR
QG063-01	E12	HQP012154	NM_000602	SERPINE1
QG063-01	F01	HQP016087	NM_134424	RAD52
QG063-01	F02	HQP000145	NM_005732	RAD50
QG063-01	F03	HQP002331	NM_000668	ADH1B
QG063-01	F04	HQP003811	NM_000771	CYP2C9
QG063-01	F05	HQP009788	NM_005544	IRS1
QG063-01	F06	HQP053959	NM_001040	SHBG
QG063-01	F07	HQP017698	NM_000348	SRD5A2
QG063-01	F08	HQP054057	BC071181	TGFB1
QG063-01	F09	HQP018565	NM_003401	XRCC4
QG063-01	F10	HQP023467	NM_000662	NAT1
QG063-01	F11	HQP002413	NM_000669	ADH1C
QG063-01	F12	HQP003871	NM_000781	CYP11A1
QG063-01	G01	HQP004948	NM_000120	EPHX1
QG063-01	G02	HQP009061	NM_001641	APEX1
QG063-01	G03	HQP009539	NM_000596	IGFBP1
QG063-01	G04	HQP009641	NM_000576	IL1B
QG063-01	G05	HQP009818	NM_000212	ITGB3
QG063-01	G06	HQP011235	NM_000249	MLH1
QG063-01	G07	HQP012021	NM_002542	OGG1
QG063-01	G08	HQP013104	NM_002634	PHB
QG063-01	G09	HQP016170	NM_021632	ZNF350
QG063-01	G10	HQP017259	NM_006747	SIPA1
QG063-01	G11	HQP021630	NM_003889	NR1H2
QG063-01	G12	HQP000245	NM_000075	CDK4
QG063-01	H01	HGDC		
QG063-01	H02	HGDC		
QG063-01	H03	HQP006940	NM_002046	GAPDH
QG063-01	H04	HQP016381	NM_001101	ACTB
QG063-01	H05	HQP015171	NM_004048	B2M
QG063-01	H06	HQP006171	NM_012423	RPL13A
QG063-01	H07	HQP009026	NM_000194	HPRT1
QG063-01	H08	HQP054253	NR_003286	RN18S1
QG063-01	H09	RT		
QG063-01	H10	RT		
QG063-01	H11	PCR		
QG063-01	H12	PCR		
QG063-02	A01	HQP021819	NM_001238	CCNE1
QG063-02	A02	HQP021754	NM_001759	CCND2
QG063-02	A03	HQP018568	NM_021141	XRCC5
QG063-02	A04	HQP018337	NM_182729	TXNRD1
QG063-02	A05	HQP018335	NM_003329	TXN

Product Data Sheet

QG063-02	A06	HQP017615	NM_000454	SOD1
QG063-02	A07	HQP016670	NM_001033886	CXCL12
QG063-02	A08	HQP016079	NM_002876	RAD51C
QG063-02	A09	HQP015496	NM_001186	BACH1
QG063-02	A10	HQP013633	NM_005037	PPARG
QG063-02	A11	HQP013596	NM_001184	ATR
QG063-02	A12	HQP011320	NM_005590	MRE11A
QG063-02	B01	HQP011257	NM_002422	MMP3
QG063-02	B02	HQP011255	NM_002421	MMP1
QG063-02	B03	HQP010581	NM_000230	LEP
QG063-02	B04	HQP009671	NM_000639	FASLG
QG063-02	B05	HQP009523	NM_000875	IGF1R
QG063-02	B06	HQP007385	NM_000515	GH1
QG063-02	B07	HQP006764	NM_012473	TXN2
QG063-02	B08	HQP006547	NM_001469	XRCC6
QG063-02	B09	HQP004986	NM_000124	ERCC6
QG063-02	B10	HQP004897	NM_001429	EP300
QG063-02	B11	HQP003808	NM_000767	CYP2B6
QG063-02	B12	HQP003120	NM_001618	PARP1
QG063-02	C01	HQP000708	NM_006440	TXNRD2
QG063-02	C02	HQP000408	NM_001800	CDKN2D
QG063-02	C03	HQP009537	NM_004970	IGFALS
QG063-02	C04	HQP054047	BC008403	HLA-DRB1
QG063-02	C05	HQP022764	NM_004832	GSTO1
QG063-02	C06	HQP021508	NM_001236	CBR3
QG063-02	C07	HQP018545	NM_000553	WRN
QG063-02	C08	HQP018418	NM_001074	UGT2B7
QG063-02	C09	HQP017784	NM_005420	SULT1E1
QG063-02	C10	HQP016844	NM_000057	BLM
QG063-02	C11	HQP016212	NM_000657	BCL2
QG063-02	C12	HQP016131	NM_000321	RB1
QG063-02	D01	HQP013635	NM_004774	PPARBP
QG063-02	D02	HQP012996	NM_016362	GHRL
QG063-02	D03	HQP011853	NM_000269	NME1
QG063-02	D04	HQP011597	NM_002467	MYC
QG063-02	D05	HQP011263	NM_004994	MMP9
QG063-02	D06	HQP010609	NM_000234	LIG1
QG063-02	D07	HQP010070	NM_002253	KDR
QG063-02	D08	HQP009678	NM_000584	IL8
QG063-02	D09	HQP009640	NM_000575	IL1A
QG063-02	D10	HQP009556	NM_000041	APOE
QG063-02	D11	HQP009024	NM_000038	APC
QG063-02	D12	HQP008493	NM_000179	MSH6
QG063-02	E01	HQP008483	NM_000849	GSTM3
QG063-02	E02	HQP008473	NM_000637	GSR
QG063-02	E03	HQP008285	NM_002085	GPX4
QG063-02	E04	HQP005084	NM_001018115	FANCD2
QG063-02	E05	HQP004975	NM_202001	ERCC1

Product Data Sheet

QG063-02	E06	HQP003499	NM_001037631	CTLA4
QG063-02	E07	HQP001282	NM_001274	CHEK1
QG063-02	E08	HQP000396	NM_001262	CDKN2C
QG063-02	E09	HQP000225	NM_001798	CDK2
QG063-02	E10	HQP004414	NM_006892	DNMT3B
QG063-02	E11	HQP053963	NM_001048	SST
QG063-02	E12	HQP053916	NM_001076	UGT2B15
QG063-02	F01	HQP023012	NM_006536	CLCA2
QG063-02	F02	HQP022978	NM_006311	NCOR1
QG063-02	F03	HQP022625	NM_004797	ADIPOQ
QG063-02	F04	HQP022529	NM_033419	PERLD1
QG063-02	F05	HQP022362	NM_004219	PTTG1
QG063-02	F06	HQP021737	NM_003927	MBD2
QG063-02	F07	HQP021518	NM_000071	CBS
QG063-02	F08	HQP021495	NM_001757	CBR1
QG063-02	F09	HQP021385	NM_003739	AKR1C3
QG063-02	F10	HQP020512	NM_181505	PPP1R1B
QG063-02	F11	HQP020117	NM_004656	BAP1
QG063-02	F12	HQP019768	NM_130445	COL18A1
QG063-02	G01	HQP018288	NM_003318	TTK
QG063-02	G02	HQP018172	NM_001067	TOP2A
QG063-02	G03	HQP018022	NM_001211	BUB1B
QG063-02	G04	HQP018016	NM_005652	TERF2
QG063-02	G05	HQP018014	NM_003218	TERF1
QG063-02	G06	HQP018013	NM_007110	TEP1
QG063-02	G07	HQP017744	NM_001050	SSTR2
QG063-02	G08	HQP017300	NM_005983	SKP2
QG063-02	G09	HQP017080	NM_003029	SHC1
QG063-02	G10	HQP016303	NM_021133	RNASEL
QG063-02	G11	HQP016082	NM_133509	RAD51L1
QG063-02	G12	HQP016076	NM_002874	RAD23B
QG063-02	H01	HGDC		
QG063-02	H02	HGDC		
QG063-02	H03	HQP006940	NM_002046	GAPDH
QG063-02	H04	HQP016381	NM_001101	ACTB
QG063-02	H05	HQP015171	NM_004048	B2M
QG063-02	H06	HQP006171	NM_012423	RPL13A
QG063-02	H07	HQP009026	NM_000194	HPRT1
QG063-02	H08	HQP054253	NR_003286	RN18S1
QG063-02	H09	RT		
QG063-02	H10	RT		
QG063-02	H11	PCR		
QG063-02	H12	PCR		
QG063-03	A01	HQP015925	NM_002843	PTPRJ
QG063-03	A02	HQP015596	NM_000962	PTGS1
QG063-03	A03	HQP015586	NM_000961	PTGIS
QG063-03	A04	HQP015539	NM_000954	PTGDS
QG063-03	A05	HQP015535	NM_000314	PTEN

Product Data Sheet

QG063-03	A06	HQP015024	NM_000948	PRL
QG063-03	A07	HQP014787	NM_018479	ECHDC1
QG063-03	A08	HQP014292	NM_018315	FBXW7
QG063-03	A09	HQP014181	NM_018193	KIAA1794
QG063-03	A10	HQP013473	NM_000446	PON1
QG063-03	A11	HQP013423	NM_002691	POLD1
QG063-03	A12	HQP012386	NM_015999	ADIPOR1
QG063-03	B01	HQP011885	NM_002518	NPAS2
QG063-03	B02	HQP011858	NM_000904	NQO2
QG063-03	B03	HQP011810	NM_020529	NFKBIA
QG063-03	B04	HQP011559	NM_001018016	MUC1
QG063-03	B05	HQP011555	NM_002454	MTRR
QG063-03	B06	HQP011009	NM_000351	STS
QG063-03	B07	HQP010612	NM_013975	LIG3
QG063-03	B08	HQP009819	NM_000213	ITGB4
QG063-03	B09	HQP009645	NM_000577	IL1RN
QG063-03	B10	HQP009633	NM_001030047	KLK3
QG063-03	B11	HQP009066	NM_002153	HSD17B2
QG063-03	B12	HQP009036	NM_005343	HRAS
QG063-03	C01	HQP008970	NM_198834	ACACA
QG063-03	C02	HQP008832	NM_181054	HIF1A
QG063-03	C03	HQP008757	NM_000410	HFE
QG063-03	C04	HQP008746	NM_001527	HDAC2
QG063-03	C05	HQP008490	NM_001513	GSTZ1
QG063-03	C06	HQP008482	NM_000848	GSTM2
QG063-03	C07	HQP008476	NM_000846	GSTA2
QG063-03	C08	HQP007767	NM_002075	GNB3
QG063-03	C09	HQP007705	NM_006708	GLO1
QG063-03	C10	HQP007396	NM_021081	GHRH
QG063-03	C11	HQP007230	NM_000583	GC
QG063-03	C12	HQP006855	NM_001042594	POT1
QG063-03	D01	HQP005462	NM_012481	IKZF3
QG063-03	D02	HQP005440	NM_022963	FGFR4
QG063-03	D03	HQP005100	NM_000691	ALDH3A1
QG063-03	D04	HQP005098	NM_022725	FANCF
QG063-03	D05	HQP005083	NM_000136	FANCC
QG063-03	D06	HQP005075	NM_000689	ALDH1A1
QG063-03	D07	HQP005058	NM_000130	F5
QG063-03	D08	HQP005052	NM_000506	F2
QG063-03	D09	HQP004983	NM_000122	ERCC3
QG063-03	D10	HQP004599	NM_001963	EGF
QG063-03	D11	HQP004392	NM_005223	DNASE1
QG063-03	D12	HQP004353	NM_004406	DMBT1
QG063-03	E01	HQP004309	NM_000791	DHFR
QG063-03	E02	HQP004081	NM_000789	ACE
QG063-03	E03	HQP003791	NM_000024	ADRB2
QG063-03	E04	HQP003539	NM_001904	CTNNB1
QG063-03	E05	HQP002920	NM_001079846	CREBBP

Product Data Sheet

QG063-03	E06	HQP002210	NM_000579	CCR5
QG063-03	E07	HQP001409	NM_006738	AKAP13
QG063-03	E08	HQP001374	NM_007182	RASSF1
QG063-03	E09	HQP001016	NM_013261	PPARGC1A
QG063-03	E10	HQP000602	NM_199141	CARM1
QG063-03	E11	HQP000274	NM_001259	CDK6
QG063-03	E12	HQP018180	NM_005427	TP73
QG063-03	F01	HQP011772	NM_006160	NEUROD2
QG063-03	F02	HQP010927	NM_002345	LUM
QG063-03	F03	HQP007399	NM_004122	GHSR
QG063-03	F04	HQP007397	NM_000823	GHRHR
QG063-03	F05	HQP000320	NM_005845	ABCC4
QG063-03	F06	HQP054052	BC015035	FGFR1
QG063-03	F07	HQP054030	NM_130398	EXO1
QG063-03	F08	HQP053962	NM_001045	SLC6A4
QG063-03	F09	HQP023451	NM_021145	DMTF1
QG063-03	F10	HQP023443	NM_004359	CDC34
QG063-03	F11	HQP023385	NM_001789	CDC25A
QG063-03	F12	HQP023365	NM_001255	CDC20
QG063-03	G01	HQP023308	NM_012432	SETDB1
QG063-03	G02	HQP023284	NM_014805	EPM2AIP1
QG063-03	G03	HQP023267	NM_001039459	ELMO1
QG063-03	G04	HQP023261	NM_001786	CDC2
QG063-03	G05	HQP023219	NM_014767	SPOCK2
QG063-03	G06	HQP023095	NM_012291	ESPL1
QG063-03	G07	HQP022998	NM_004917	KLK4
QG063-03	G08	HQP022828	NM_003615	SLC4A7
QG063-03	G09	HQP022827	NM_018488	TBX4
QG063-03	G10	HQP022805	NM_004850	ROCK2
QG063-03	G11	HQP022802	NM_001039477	C1orf38
QG063-03	G12	HQP022796	NM_004273	CHST3
QG063-03	H01	HGDC		
QG063-03	H02	HGDC		
QG063-03	H03	HQP006940	NM_002046	GAPDH
QG063-03	H04	HQP016381	NM_001101	ACTB
QG063-03	H05	HQP015171	NM_004048	B2M
QG063-03	H06	HQP006171	NM_012423	RPL13A
QG063-03	H07	HQP009026	NM_000194	HPRT1
QG063-03	H08	HQP054253	NR_003286	RN18S1
QG063-03	H09	RT		
QG063-03	H10	RT		
QG063-03	H11	PCR		
QG063-03	H12	PCR		
QG063-04	A01	HQP022640	NM_138732	NRXN2
QG063-04	A02	HQP022615	NM_004795	KL
QG063-04	A03	HQP022444	NM_004763	ITGB1BP1
QG063-04	A04	HQP022316	NM_000616	CD4
QG063-04	A05	HQP022268	NM_001007793	BUB3

Product Data Sheet

QG063-04	A06	HQP022093	NM_004688	NMI
QG063-04	A07	HQP021655	NM_004665	VNN2
QG063-04	A08	HQP021621	NM_003884	PCAF
QG063-04	A09	HQP021597	NM_003873	NRP1
QG063-04	A10	HQP021589	NM_003872	NRP2
QG063-04	A11	HQP021586	NM_198061	CES2
QG063-04	A12	HQP021557	NM_003844	TNFRSF10A
QG063-04	B01	HQP021553	NM_003842	TNFRSF10B
QG063-04	B02	HQP021502	NM_003810	TNFSF10
QG063-04	B03	HQP021492	NM_003804	RIPK1
QG063-04	B04	HQP021405	NM_003747	TNKS
QG063-04	B05	HQP021313	NM_001753	CAV1
QG063-04	B06	HQP021232	NM_033401	CNTNAP4
QG063-04	B07	HQP021225	NM_006769	LMO4
QG063-04	B08	HQP021130	NM_003632	CNTNAP1
QG063-04	B09	HQP021074	NM_032875	FBXL20
QG063-04	B10	HQP020830	NM_003597	KLF11
QG063-04	B11	HQP020699	NM_021111	RECK
QG063-04	B12	HQP020507	NM_139076	CCDC98
QG063-04	C01	HQP020116	NM_004655	AXIN2
QG063-04	C02	HQP020114	NM_003502	AXIN1
QG063-04	C03	HQP020042	NM_003489	NRIP1
QG063-04	C04	HQP019938	NM_030928	CDT1
QG063-04	C05	HQP019836	NM_012074	DPF3
QG063-04	C06	HQP019829	NM_030782	CLPTM1L
QG063-04	C07	HQP019718	NM_025247	ACAD10
QG063-04	C08	HQP019694	NM_025235	TNKS2
QG063-04	C09	HQP019689	NM_001042529	COASY
QG063-04	C10	HQP019452	NM_001742	CALCR
QG063-04	C11	HQP019284	NM_024757	EHMT1
QG063-04	C12	HQP019268	NM_006304	SHFM1
QG063-04	D01	HQP019159	NM_024626	VTCN1
QG063-04	D02	HQP019090	NM_024551	ADIPOR2
QG063-04	D03	HQP018802	NM_001008540	CXCR4
QG063-04	D04	HQP018755	NM_001007257	PRDM2
QG063-04	D05	HQP018519	NM_003390	WEE1
QG063-04	D06	HQP018485	NM_000551	VHL
QG063-04	D07	HQP018447	NM_001017415	USP1
QG063-04	D08	HQP018438	NM_006003	UQCRFS1
QG063-04	D09	HQP018403	NM_003355	UCP2
QG063-04	D10	HQP018379	NM_003345	UBE2I
QG063-04	D11	HQP018328	NM_000474	TWIST1
QG063-04	D12	HQP018271	NM_000369	TSHR
QG063-04	E01	HQP018176	NM_005657	TP53BP1
QG063-04	E02	HQP018149	NM_001066	TNFRSF1B
QG063-04	E03	HQP018115	NM_003265	TLR3
QG063-04	E04	HQP018047	NM_003238	TGFB2
QG063-04	E05	HQP018040	NM_003227	TFR2

Product Data Sheet

QG063-04	E06	HQP017999	NM_004336	BUB1
QG063-04	E07	HQP017938	NM_001061	TBXAS1
QG063-04	E08	HQP017929	NM_005994	TBX2
QG063-04	E09	HQP017928	NM_003194	TBP
QG063-04	E10	HQP017866	NM_003183	ADAM17
QG063-04	E11	HQP017767	NM_003150	STAT3
QG063-04	E12	HQP017764	NM_007315	STAT1
QG063-04	F01	HQP017747	NM_001053	SSTR5
QG063-04	F02	HQP017745	NM_001051	SSTR3
QG063-04	F03	HQP017703	NM_001005291	SREBF1
QG063-04	F04	HQP017697	NM_001047	SRD5A1
QG063-04	F05	HQP017505	NM_003070	SMARCA2
QG063-04	F06	HQP017483	NM_005415	SLC20A1
QG063-04	F07	HQP017482	NM_194255	SLC19A1
QG063-04	F08	HQP016801	NM_022162	NOD2
QG063-04	F09	HQP016634	NM_001713	BHMT
QG063-04	F10	HQP016620	NM_002981	CCL1
QG063-04	F11	HQP016466	NM_001006665	RPS6KA1
QG063-04	F12	HQP016426	NM_021134	MRPL23
QG063-04	G01	HQP016391	NM_002947	RPA3
QG063-04	G02	HQP016390	NM_002946	RPA2
QG063-04	G03	HQP016389	NM_002945	RPA1
QG063-04	G04	HQP016381	NM_001101	ACTB
QG063-04	G05	HQP016223	NM_002912	REV3L
QG063-04	G06	HQP016213	NM_021975	RELA
QG063-04	G07	HQP016155	NM_002894	RBBP8
QG063-04	G08	HQP016070	NM_002873	RAD17
QG063-04	G09	HQP015947	NM_002853	RAD1
QG063-04	G10	HQP015929	NM_002847	PTPRN2
QG063-04	G11	HQP015877	NM_021177	LSM2
QG063-04	G12	HQP015784	NM_020898	CALCOCO1
QG063-04	H01	HGDC		
QG063-04	H02	HGDC		
QG063-04	H03	HQP006940	NM_002046	GAPDH
QG063-04	H04	HQP016381	NM_001101	ACTB
QG063-04	H05	HQP015171	NM_004048	B2M
QG063-04	H06	HQP006171	NM_012423	RPL13A
QG063-04	H07	HQP009026	NM_000194	HPRT1
QG063-04	H08	HQP054253	NR_003286	RN18S1
QG063-04	H09	RT		
QG063-04	H10	RT		
QG063-04	H11	PCR		
QG063-04	H12	PCR		
QG063-05	A01	HQP015771	NM_020884	MYH7B
QG063-05	A02	HQP015662	NM_001001924	MTUS1
QG063-05	A03	HQP015657	NM_020744	MTA3
QG063-05	A04	HQP015610	NM_002820	PTHLH
QG063-05	A05	HQP015530	NM_000264	PTCH1

Product Data Sheet

QG063-05	A06	HQP015124	NM_016567	BCCIP
QG063-05	A07	HQP015009	NM_198393	TEX14
QG063-05	A08	HQP014896	NM_002752	MAPK9
QG063-05	A09	HQP014886	NM_002750	MAPK8
QG063-05	A10	HQP014706	NM_002737	PRKCA
QG063-05	A11	HQP014454	NM_018725	IL17RB
QG063-05	A12	HQP014301	NM_006246	PPP2R5E
QG063-05	B01	HQP014174	NM_002717	PPP2R2A
QG063-05	B02	HQP014161	NM_002716	PPP2R1B
QG063-05	B03	HQP014012	NM_017975	ZWILCH
QG063-05	B04	HQP013970	NM_002709	PPP1CB
QG063-05	B05	HQP013636	NM_000489	ATRX
QG063-05	B06	HQP013617	NM_001001928	PPARA
QG063-05	B07	HQP013583	NM_205862	UGT1A6
QG063-05	B08	HQP013450	NM_018975	TERF2IP
QG063-05	B09	HQP013422	NM_002690	POLB
QG063-05	B10	HQP013403	NM_004574	SEPT4
QG063-05	B11	HQP013388	NM_017442	TLR9
QG063-05	B12	HQP013382	NM_002686	PNMT
QG063-05	C01	HQP013262	NM_005030	PLK1
QG063-05	C02	HQP013151	NM_006219	PIK3CB
QG063-05	C03	HQP013087	NM_002630	PGC
QG063-05	C04	HQP013013	NM_016507	CRKRS
QG063-05	C05	HQP012986	NM_016290	UIMC1
QG063-05	C06	HQP012770	NM_016371	HSD17B7
QG063-05	C07	HQP012755	NM_001037872	REV1
QG063-05	C08	HQP012478	NM_016261	TUBD1
QG063-05	C09	HQP012457	NM_001012509	SLC45A2
QG063-05	C10	HQP012447	NM_016142	HSD17B12
QG063-05	C11	HQP012420	NM_002592	PCNA
QG063-05	C12	HQP012393	NM_016009	SH3GLB1
QG063-05	D01	HQP012371	NM_032951	MLXIPL
QG063-05	D02	HQP012335	NM_015895	GMNN
QG063-05	D03	HQP012013	NM_002539	ODC1
QG063-05	D04	HQP011807	NM_003998	NFKB1
QG063-05	D05	HQP011764	NM_002497	NEK2
QG063-05	D06	HQP011694	NM_000631	NCF4
QG063-05	D07	HQP011693	NM_000433	NCF2
QG063-05	D08	HQP011661	NM_004535	MYT1
QG063-05	D09	HQP011492	NM_002439	MSH3
QG063-05	D10	HQP011491	NM_000251	MSH2
QG063-05	D11	HQP011304	NM_001015052	MPG
QG063-05	D12	HQP011266	NM_002426	MMP12
QG063-05	E01	HQP011205	NM_003994	KITLG
QG063-05	E02	HQP011140	NM_002393	MDM4
QG063-05	E03	HQP011085	NM_001025081	MBP
QG063-05	E04	HQP011047	NM_005911	MAT2A
QG063-05	E05	HQP010966	NM_005904	SMAD7

Product Data Sheet

QG063-05	E06	HQP010870	NM_002332	LRP1
QG063-05	E07	HQP010847	NM_000237	LPL
QG063-05	E08	HQP010410	NM_001042678	RHOC
QG063-05	E09	HQP010260	NM_001005404	YPEL2
QG063-05	E10	HQP010124	NM_002265	KPNB1
QG063-05	E11	HQP010104	NM_000892	KLKB1
QG063-05	E12	HQP010098	NM_002256	KISS1
QG063-05	F01	HQP009853	NM_002228	JUN
QG063-05	F02	HQP009825	NM_002213	ITGB5
QG063-05	F03	HQP009808	NM_002210	ITGAV
QG063-05	F04	HQP009774	NM_000210	ITGA6
QG063-05	F05	HQP009764	NM_000208	INSR
QG063-05	F06	HQP009749	NM_000207	INS
QG063-05	F07	HQP009718	NM_001562	IL18
QG063-05	F08	HQP009697	NM_002188	IL13
QG063-05	F09	HQP009696	NM_001559	IL12RB2
QG063-05	F10	HQP009662	NM_000589	IL4
QG063-05	F11	HQP009571	NM_002180	IGHMBP2
QG063-05	F12	HQP009558	NM_001553	IGFBP7
QG063-05	G01	HQP009551	NM_000599	IGFBP5
QG063-05	G02	HQP009541	NM_000597	IGFBP2
QG063-05	G03	HQP009529	NM_000612	IGF2
QG063-05	G04	HQP009138	NM_004507	HUS1
QG063-05	G05	HQP009123	NM_000869	HTR3A
QG063-05	G06	HQP009062	NM_005525	HSD11B1
QG063-05	G07	HQP009059	NM_000862	HSD3B1
QG063-05	G08	HQP009053	NM_001536	PRMT1
QG063-05	G09	HQP009019	NM_005143	HP
QG063-05	G10	HQP008904	NM_004496	FOXA1
QG063-05	G11	HQP008896	NM_012484	HMMR
QG063-05	G12	HQP008849	NM_002116	HLA-A
QG063-05	H01	HGDC		
QG063-05	H02	HGDC		
QG063-05	H03	HQP006940	NM_002046	GAPDH
QG063-05	H04	HQP016381	NM_001101	ACTB
QG063-05	H05	HQP015171	NM_004048	B2M
QG063-05	H06	HQP006171	NM_012423	RPL13A
QG063-05	H07	HQP009026	NM_000194	HPRT1
QG063-05	H08	HQP054253	NR_003286	RN18S1
QG063-05	H09	RT		
QG063-05	H10	RT		
QG063-05	H11	PCR		
QG063-05	H12	PCR		
QG063-06	A01	HQP008801	NM_001010931	HGF
QG063-06	A02	HQP008694	NM_014585	SLC40A1
QG063-06	A03	HQP008690	NM_004131	GZMB
QG063-06	A04	HQP008498	NM_005513	GTF2E1
QG063-06	A05	HQP008494	NM_015859	GTF2A1

Product Data Sheet

QG063-06	A06	HQP008469	NM_002093	GSK3B
QG063-06	A07	HQP008340	NM_000830	GRIK1
QG063-06	A08	HQP007811	NM_002081	GPC1
QG063-06	A09	HQP007803	NM_000173	GP1BA
QG063-06	A10	HQP007781	NM_000406	GNRHR
QG063-06	A11	HQP007755	NM_000516	GNAS
QG063-06	A12	HQP007697	NM_012201	GLG1
QG063-06	B01	HQP007465	NM_014389	PELP1
QG063-06	B02	HQP007382	NM_012449	STEAP1
QG063-06	B03	HQP007303	NM_012138	AATF
QG063-06	B04	HQP007210	NM_012153	EHF
QG063-06	B05	HQP006715	NM_012415	RAD54B
QG063-06	B06	HQP006375	NM_001629	ALOX5AP
QG063-06	B07	HQP006299	NM_001031696	PLD3
QG063-06	B08	HQP006023	NM_054034	FN1
QG063-06	B09	HQP005730	NM_015056	RRP1B
QG063-06	B10	HQP005589	NM_012072	CD93
QG063-06	B11	HQP005496	NM_014906	PPM1E
QG063-06	B12	HQP005458	NM_012250	RRAS2
QG063-06	C01	HQP005391	NM_004111	FEN1
QG063-06	C02	HQP005292	NM_001002275	FCGR2B
QG063-06	C03	HQP005280	NM_021642	FCGR2A
QG063-06	C04	HQP005238	NM_002001	FCER1A
QG063-06	C05	HQP005210	NM_000138	FBN1
QG063-06	C06	HQP005134	NM_004104	FASN
QG063-06	C07	HQP005087	NM_000690	ALDH2
QG063-06	C08	HQP005081	NM_000135	FANCA
QG063-06	C09	HQP005005	NM_001438	ESRRG
QG063-06	C10	HQP004991	NM_001014431	AKT1
QG063-06	C11	HQP004727	NM_005502	ABCA1
QG063-06	C12	HQP004526	NM_004091	E2F2
QG063-06	D01	HQP004494	NM_000029	AGT
QG063-06	D02	HQP004446	NM_000795	DRD2
QG063-06	D03	HQP004275	NM_199355	ADAMTS18
QG063-06	D04	HQP004260	NM_130770	HTR3C
QG063-06	D05	HQP004156	NM_001357	DHX9
QG063-06	D06	HQP004072	NM_133504	DCN
QG063-06	D07	HQP003917	NM_000782	CYP24A1
QG063-06	D08	HQP003817	NM_000773	CYP2E1
QG063-06	D09	HQP003812	NM_000025	ADRB3
QG063-06	D10	HQP003810	NM_000770	CYP2C8
QG063-06	D11	HQP003654	NM_181780	BTLA
QG063-06	D12	HQP003495	NM_144701	IL23R
QG063-06	E01	HQP003395	NM_001033521	CSTF1
QG063-06	E02	HQP003249	NM_001894	CSNK1E
QG063-06	E03	HQP003240	NM_001893	CSNK1D
QG063-06	E04	HQP003161	NM_152434	CWF19L2
QG063-06	E05	HQP003084	NM_021117	CRY2

Product Data Sheet

QG063-06	E06	HQP002760	NM_152405	JMY
QG063-06	E07	HQP002370	NM_133261	GIPC3
QG063-06	E08	HQP002301	NM_000667	ADH1A
QG063-06	E09	HQP002217	NM_001838	CCR7
QG063-06	E10	HQP002019	NM_147191	MMP21
QG063-06	E11	HQP001414	NM_007202	AKAP10
QG063-06	E12	HQP001408	NM_007199	IRAK3
QG063-06	F01	HQP001109	NM_006826	YWHAQ
QG063-06	F02	HQP001103	NM_006819	STIP1
QG063-06	F03	HQP001051	NM_025256	EHMT2
QG063-06	F04	HQP001047	NM_014599	MAGED2
QG063-06	F05	HQP000970	NM_006663	PPP1R13L
QG063-06	F06	HQP000963	NM_012190	ALDH1L1
QG063-06	F07	HQP000809	NM_006565	CTCF
QG063-06	F08	HQP000764	NM_006475	POSTN
QG063-06	F09	HQP000759	NM_001813	CENPE
QG063-06	F10	HQP000583	NM_006361	HOXB13
QG063-06	F11	HQP000552	NM_006341	MAD2L2
QG063-06	F12	HQP000540	NM_006114	TOMM40
QG063-06	G01	HQP000519	NM_006331	EMG1
QG063-06	G02	HQP000285	NM_001799	CDK7
QG063-06	G03	HQP000179	NM_147185	AKAP9
QG063-06	G04	HQP054687	NM_005961	MUC6
QG063-06	G05	HQP022027	NM_004675	DIRAS3
QG063-06	G06	HQP020006	NM_030974	SHARPIN
QG063-06	G07	HQP018213	NM_000367	TPMT
QG063-06	G08	HQP017746	NM_001052	SSTR4
QG063-06	G09	HQP017743	NM_001049	SSTR1
QG063-06	G10	HQP015106	NM_001077500	KLK10
QG063-06	G11	HQP013519	NM_000306	POU1F1
QG063-06	G12	HQP011865	NM_000620	NOS1
QG063-06	H01	HGDC		
QG063-06	H02	HGDC		
QG063-06	H03	HQP006940	NM_002046	GAPDH
QG063-06	H04	HQP016381	NM_001101	ACTB
QG063-06	H05	HQP015171	NM_004048	B2M
QG063-06	H06	HQP006171	NM_012423	RPL13A
QG063-06	H07	HQP009026	NM_000194	HPRT1
QG063-06	H08	HQP054253	NR_003286	RN18S1
QG063-06	H09	RT		
QG063-06	H10	RT		
QG063-06	H11	PCR		
QG063-06	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Breast 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

GeneCopoeia Products are for Research Use Only

Trademarks: GeneCopoeia™, ExProfile™, All-in-One™, (GeneCopoeia Inc.); RNAsol® RT (Molecular Research Center, Inc.); SYBR® (Molecular Probes); iCycler iQ®, MyiQ™, iQ™5, CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™ (Bio-Rad); LightCycler® (Roche); Trizol™, ABI®, ROX®, ViiA™, StepOnePlus™ (Life Technologies).

Copyright © 2016 GeneCopoeia, Inc.

QG063-160202