

## ExProfile™ Human Bladder Cancer Gene qPCR Array

For focused group profiling of human bladder cancer related genes expression

Cat. No. QG064-A (5 x 96-well plate, Format A)

Cat. No. QG064-B (5 x 96-well plate, Format B)

Cat. No. QG064-C (5 x 96-well plate, Format C)

Cat. No. QG064-D (5 x 96-well plate, Format D)

Cat. No. QG064-E (5 x 96-well plate, Format E)

Available as 1 set or 6 sets. Each set contains 420 unique gene primers deposited in five 96-well plates.

### Introduction

The ExProfile human bladder cancer gene qPCR array profiles 420 human genes to aberrantly expressed human genes involved in human bladder 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 bladder cancer gene array allows researchers to study the cancer-related genes to gain understanding of their roles in bladder cancer pathogenesis.

- QG064 plate 01: 84 unique gene PCR primer pairs
- QG064 plate 02: 84 unique gene PCR primer pairs
- QG064 plate 03: 84 unique gene PCR primer pairs
- QG064 plate 04: 84 unique gene PCR primer pairs
- QG064 plate 05: 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
<b>A</b> (96-well)	Applied Biosystems	5700, 7000, 7300, 7500, 7700, 7900HT (Standard 96-well block), ViiA™7 (Standard 96-well block)
<b>B</b> (96-well)	Applied Biosystems	7500 (Fast block), 7900HT (Fast block), StepOnePlus™, ViiA™7 (Fast block)
<b>C</b> (96-well)	Bio-Rad Laboratories	iCycler iQ®, MyiQ™, iQ™5
<b>D</b> (96-well)	Bio-Rad Laboratories	CFX96™, DNA Engine Opticon™, DNA Engine Opticon 2™, Chromo4™
<b>E</b> (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	NAT2	XRCC1	TP53	ERCC2	XRCC3	NAT1	MTHFR	CYP1A1	XPC	FGFR3	OGG1	NQO1
B	APEX1	TNF	CCND1	ERCC5	SULT1A1	SOD2	NBN	MTR	MDM2	ERCC6	ERCC1	CYP2E1
C	CDH1	XRCC4	TGFB1	MPO	MMP1	IL6	IL1RN	IL1B	ERCC4	CYP2D6	CYP1B1	CASP8
D	XRCC5	XRCC2	RAD23B	PTGS2	NFKB1	MSH2	KRAS	IL4	HRAS	MSH6	GSTM3	COMT
E	VEGFA	UGT2B7	TYMS	BRCA2	RAD52	POLD1	POLB	PCNA	NOS2A	MTRR	MMP9	MMP3
F	MLH1	LIG1	IL10	IL8	APOB	EPHX1	AHR	EGFR	PARP1	CDKN2A	TGFB1	RGS6
G	CCNH	NFRSF10A	CBS	AKR1C3	VDR	BRAF	SLC19A1	ABCG8	BLM	CCL2	ZNF350	PPARG
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG064 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	UGT1A4	PMS2	PMS1	MYC	MSH3	MMP12	MMP8	MMP2	ARNT	LIG3	AR	IL12A
B	FASLG	IL4R	IL2	IL1A	APC	FANCA	ERCC3	EPHX2	EGF	DRD2	CYP2C9	CYP2C19
C	CTH	PARP4	ADH1C	CETP	ARL6IP5	CDK4	PARP3	CYP7A1	LOC729230	MUC5B	RET	EXO1
D	SLC2A4	SHBG	NR1H4	SLC23A1	CDC25C	CDC25B	CDC25A	RB1CC1	MATR3	CD81	CD40	VPS26A
E	TP53I3	MPDU1	GDF15	LITAF	SCARB1	GSTO1	CD86	CYP7B1	CD80	41532	RECQL4	CD14
F	CD4	CG018	CCDC97	CCNE1	MBD2	MBD4	CCKAR	GGH	CD80	41532	RECQL4	CD14
G	PTCH2	CAV1	RGS5	CASR	RAD54L	NICN1	PLA2G6	CASP3	FZD7	AXIN2	NCOA3	CLPTM1L
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG064 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	COL18A1	MGC4093	COASY	CALCR	SLC30A4	ZNF230	XBP1	WRN	VIL2	VCAM1	UCP3	TYR
B	TSG101	TSC2	TGM1	TFF1	TERF2	TERF1	TEP1	TDG	STK11	STAT1	SSTR3	BPI
C	SOD1	SMO	SLC4A2	SLAMF1	SFTPD	SEPP1	NOD2	SELE	CXCL12	BHMT	CCL5	RXRβ
D	RXRA	BCR	RPA3	RPA2	RPA1	ROS1	RNASEL	RHEB	RGS2	RFC5	RFC3	RFC2
E	RFC1	REV3L	GNB4	RAD51C	RAC1	PCTP	BARD1	BAK1	MYH7B	PTGS1	PTEN	MYNN
F	HIF1AN	FBXW7	WDR79	UGT1A7	PON1	POLRMT	POLR2K	POLR2J	POLR2F	POLR2E	POLR2D	POLR2C
G	POLR2B	POLG	POLE2	POLE	POLD2	PLK1	PLAU	PLA2G2A	PIM1	PHB	ABCB1	PGR
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG064 plate 03

## Product Data Sheet

	1	2	3	4	5	6	7	8	9	10	11	12
A	ENPP1	POLK	HAO2	P2RX7	OPRM1	OCA2	NRAS	NOS3	NQO2	ATP1B2	NINJ1	NFKBIE
B	NCF2	MX1	MRE11A	MPG	METTL1	MET	MEST	MAOA	LRPAP1	LRP5	LRP6	LPL
C	LIPC	LIG4	ARHGDI3	LEPR	LEP	LDLR	LCAT	KIR3DL1	RAB15	JAK3	IRS1	IRF3
D	IRF1	INSR	IL15RA	IL15	IL13	IL12B	IL8RA	IL7R	IL6R	IL1R1	APOE	IGFBP6
E	SLC6A18	IGFBP5	IGFBP3	IGFBP2	IGFBP1	IGF2	IGF1R	IGF1	IFNGR2	IFNGR1	IFNAR2	IFI27
F	APOA4	APOA2	HUS1	HTR1D	HTR1B	BIRC2	HSD17B4	HSD17B2	HSD17B1	HSD3B2	HSD3B1	APAF1
G	HMGCR	HIF1A	HFE	HADHA	SLC39A2	RPA4	GSTZ1	GSTA4	GSK3B	GRPR	GPX4	GPX3
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG064 plate 04

	1	2	3	4	5	6	7	8	9	10	11	12
A	GPX2	GNB3	GNAS	POLL	DHDH	RGS17	GC	HSPB8	KRT23	POT1	LMOD1	ANKK1
B	XRCC6	FUT7	ALOX15	ALOX5	ALOX12	POLA2	AMACR	SEC14L2	FOS	ABCA5	ABCA6	FOXC1
C	FGFR4	FCGR2A	ALDH2	FANCD2	FANCC	ESR2	ESR1	AKT1	DNAJC18	ABCA1	EFNB3	EDN1
D	DRD4	DRD1	DHFR	DDB2	ACE	C11orf65	CYP24A1	CYP19A1	CYP17A1	CYP3A5	CYP2A13	CYBB
E	CX3CR1	CTSH	CTNNA1	CTLA4	CSTF1	CSF3	CSF2	CSF1R	CR1	CCR5	ERCC8	CHRNA3
F	SAT2	DDX20	POLI	CHEK2	CHEK1	AKR1C4	PPP1R13L	ALDH1L1	CGA	MTHFD2	POLD3	TXNRD2
G	TACC3	AKR1A1	CDKN1C	CDK7	CDK6	CDK5	RAD50	NUBP2	AKT3	WDFY4	TCTA	SOD3
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure5. Illustration of QG064 plate 05

- **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
QG064-01	A01	HQP001136	NM_000015	NAT2
QG064-01	A02	HQP018562	NM_006297	XRCC1
QG064-01	A03	HQP018175	NM_000546	TP53
QG064-01	A04	HQP004976	NM_000400	ERCC2
QG064-01	A05	HQP018564	NM_005432	XRCC3
QG064-01	A06	HQP023467	NM_000662	NAT1
QG064-01	A07	HQP011547	NM_005957	MTHFR
QG064-01	A08	HQP003772	NM_000499	CYP1A1
QG064-01	A09	HQP018556	NM_004628	XPC
QG064-01	A10	HQP005434	NM_000142	FGFR3
QG064-01	A11	HQP012021	NM_002542	OGG1
QG064-01	A12	HQP004317	NM_000903	NQO1
QG064-01	B01	HQP009061	NM_001641	APEX1
QG064-01	B02	HQP018141	NM_000594	TNF

## Product Data Sheet

QG064-01	B03	HQP016204	NM_053056	CCND1
QG064-01	B04	HQP004985	NM_000123	ERCC5
QG064-01	B05	HQP017811	NM_177536	SULT1A1
QG064-01	B06	HQP017616	NM_000636	SOD2
QG064-01	B07	HQP011687	NM_002485	NBN
QG064-01	B08	HQP011554	NM_000254	MTR
QG064-01	B09	HQP011135	NM_002392	MDM2
QG064-01	B10	HQP004986	NM_000124	ERCC6
QG064-01	B11	HQP004975	NM_202001	ERCC1
QG064-01	B12	HQP003817	NM_000773	CYP2E1
QG064-01	C01	HQP023466	NM_004360	CDH1
QG064-01	C02	HQP018565	NM_003401	XRCC4
QG064-01	C03	HQP018044	NM_000660	TGFB1
QG064-01	C04	HQP011309	NM_000250	MPO
QG064-01	C05	HQP011255	NM_002421	MMP1
QG064-01	C06	HQP009670	NM_000600	IL6
QG064-01	C07	HQP009645	NM_000577	IL1RN
QG064-01	C08	HQP009641	NM_000576	IL1B
QG064-01	C09	HQP004984	NM_005236	ERCC4
QG064-01	C10	HQP003814	NM_000106	CYP2D6
QG064-01	C11	HQP003775	NM_000104	CYP1B1
QG064-01	C12	HQP018966	NM_001080124	CASP8
QG064-01	D01	HQP018568	NM_021141	XRCC5
QG064-01	D02	HQP018563	NM_005431	XRCC2
QG064-01	D03	HQP016076	NM_002874	RAD23B
QG064-01	D04	HQP015598	NM_000963	PTGS2
QG064-01	D05	HQP011807	NM_003998	NFKB1
QG064-01	D06	HQP011491	NM_000251	MSH2
QG064-01	D07	HQP010133	NM_004985	KRAS
QG064-01	D08	HQP009662	NM_000589	IL4
QG064-01	D09	HQP009036	NM_005343	HRAS
QG064-01	D10	HQP008493	NM_000179	MSH6
QG064-01	D11	HQP008483	NM_000849	GSTM3
QG064-01	D12	HQP002671	NM_000754	COMT
QG064-01	E01	HQP018475	NM_001025366	VEGFA
QG064-01	E02	HQP018418	NM_001074	UGT2B7
QG064-01	E03	HQP018342	NM_001071	TYMS
QG064-01	E04	HQP017753	NM_000059	BRCA2
QG064-01	E05	HQP016087	NM_134424	RAD52
QG064-01	E06	HQP013423	NM_002691	POLD1
QG064-01	E07	HQP013422	NM_002690	POLB
QG064-01	E08	HQP012420	NM_002592	PCNA
QG064-01	E09	HQP011866	NM_000625	NOS2A
QG064-01	E10	HQP011555	NM_002454	MTRR
QG064-01	E11	HQP011263	NM_004994	MMP9
QG064-01	E12	HQP011257	NM_002422	MMP3
QG064-01	F01	HQP011235	NM_000249	MLH1
QG064-01	F02	HQP010609	NM_000234	LIG1

Product Data Sheet

QG064-01	F03	HQP009685	NM_000572	IL10
QG064-01	F04	HQP009678	NM_000584	IL8
QG064-01	F05	HQP009218	NM_000384	APOB
QG064-01	F06	HQP004948	NM_000120	EPHX1
QG064-01	F07	HQP004658	NM_001621	AHR
QG064-01	F08	HQP004605	NM_005228	EGFR
QG064-01	F09	HQP003120	NM_001618	PARP1
QG064-01	F10	HQP000370	NM_058195	CDKN2A
QG064-01	F11	HQP054057	BC071181	TGFBR1
QG064-01	F12	HQP023003	NM_004296	RGS6
QG064-01	G01	HQP021906	NM_001239	CCNH
QG064-01	G02	HQP021557	NM_003844	TNFRSF10A
QG064-01	G03	HQP021518	NM_000071	CBS
QG064-01	G04	HQP021385	NM_003739	AKR1C3
QG064-01	G05	HQP018474	NM_000376	VDR
QG064-01	G06	HQP017733	NM_004333	BRAF
QG064-01	G07	HQP017482	NM_194255	SLC19A1
QG064-01	G08	HQP016876	NM_022437	ABCG8
QG064-01	G09	HQP016844	NM_000057	BLM
QG064-01	G10	HQP016621	NM_002982	CCL2
QG064-01	G11	HQP016170	NM_021632	ZNF350
QG064-01	G12	HQP013633	NM_005037	PPARG
QG064-01	H01	HGDC		
QG064-01	H02	HGDC		
QG064-01	H03	HQP006940	NM_002046	GAPDH
QG064-01	H04	HQP016381	NM_001101	ACTB
QG064-01	H05	HQP015171	NM_004048	B2M
QG064-01	H06	HQP006171	NM_012423	RPL13A
QG064-01	H07	HQP009026	NM_000194	HPRT1
QG064-01	H08	HQP054253	NR_003286	RN18S1
QG064-01	H09	RT		
QG064-01	H10	RT		
QG064-01	H11	PCR		
QG064-01	H12	PCR		
QG064-02	A01	HQP013614	NM_007120	UGT1A4
QG064-02	A02	HQP013352	NM_000535	PMS2
QG064-02	A03	HQP013312	NM_000534	PMS1
QG064-02	A04	HQP011597	NM_002467	MYC
QG064-02	A05	HQP011492	NM_002439	MSH3
QG064-02	A06	HQP011266	NM_002426	MMP12
QG064-02	A07	HQP011262	NM_002424	MMP8
QG064-02	A08	HQP011256	NM_004530	MMP2
QG064-02	A09	HQP010924	NM_001668	ARNT
QG064-02	A10	HQP010612	NM_013975	LIG3
QG064-02	A11	HQP009801	NM_000044	AR
QG064-02	A12	HQP009692	NM_000882	IL12A
QG064-02	B01	HQP009671	NM_000639	FASLG
QG064-02	B02	HQP009664	NM_000418	IL4R

## Product Data Sheet

QG064-02	B03	HQP009649	NM_000586	IL2
QG064-02	B04	HQP009640	NM_000575	IL1A
QG064-02	B05	HQP009024	NM_000038	APC
QG064-02	B06	HQP005081	NM_000135	FANCA
QG064-02	B07	HQP004983	NM_000122	ERCC3
QG064-02	B08	HQP004950	NM_001979	EPHX2
QG064-02	B09	HQP004599	NM_001963	EGF
QG064-02	B10	HQP004446	NM_000795	DRD2
QG064-02	B11	HQP003811	NM_000771	CYP2C9
QG064-02	B12	HQP003809	NM_000769	CYP2C19
QG064-02	C01	HQP003493	NM_001902	CTH
QG064-02	C02	HQP003171	NM_006437	PARP4
QG064-02	C03	HQP002413	NM_000669	ADH1C
QG064-02	C04	HQP000853	NM_000078	CETP
QG064-02	C05	HQP000657	NM_006407	ARL6IP5
QG064-02	C06	HQP000245	NM_000075	CDK4
QG064-02	C07	HQP000050	NM_001003931	PARP3
QG064-02	C08	HQP003859	NM_000780	CYP7A1
QG064-02	C09	HQP054686	NM_001123396	LOC729230
QG064-02	C10	HQP054195	NM_002458	MUC5B
QG064-02	C11	HQP054045	BC004257	RET
QG064-02	C12	HQP054030	NM_130398	EXO1
QG064-02	D01	HQP053960	NM_001042	SLC2A4
QG064-02	D02	HQP053959	NM_001040	SHBG
QG064-02	D03	HQP023437	NM_005123	NR1H4
QG064-02	D04	HQP023414	NM_005847	SLC23A1
QG064-02	D05	HQP023409	NM_001790	CDC25C
QG064-02	D06	HQP023399	NM_004358	CDC25B
QG064-02	D07	HQP023385	NM_001789	CDC25A
QG064-02	D08	HQP023240	NM_014781	RB1CC1
QG064-02	D09	HQP023195	NM_018834	MATR3
QG064-02	D10	HQP023168	NM_004356	CD81
QG064-02	D11	HQP022955	NM_001250	CD40
QG064-02	D12	HQP022910	NM_001035260	VPS26A
QG064-02	E01	HQP022884	NM_004881	TP53I3
QG064-02	E02	HQP022863	NM_004870	MPDU1
QG064-02	E03	HQP022853	NM_004864	GDF15
QG064-02	E04	HQP022851	NM_004862	LITAF
QG064-02	E05	HQP022833	NM_005505	SCARB1
QG064-02	E06	HQP022764	NM_004832	GSTO1
QG064-02	E07	HQP022747	NM_175862	CD86
QG064-02	E08	HQP022723	NM_004820	CYP7B1
QG064-02	E09	HQP022722	NM_005191	CD80
QG064-02	E10	HQP022686	NM_004261	15-Sep
QG064-02	E11	HQP022676	NM_004260	RECQL4
QG064-02	E12	HQP022490	NM_000591	CD14
QG064-02	F01	HQP022316	NM_000616	CD4
QG064-02	F02	HQP022001	NM_001079691	CG018

Product Data Sheet

QG064-02	F03	HQP021913	NM_052848	CCDC97
QG064-02	F04	HQP021819	NM_001238	CCNE1
QG064-02	F05	HQP021737	NM_003927	MBD2
QG064-02	F06	HQP021736	NM_003925	MBD4
QG064-02	F07	HQP021647	NM_000730	CCKAR
QG064-02	F08	HQP021603	NM_003878	GGH
QG064-02	F09	HQP021508	NM_001236	CBR3
QG064-02	F10	HQP021495	NM_001757	CBR1
QG064-02	F11	HQP021405	NM_003747	TNKS
QG064-02	F12	HQP021388	NM_003742	ABCB11
QG064-02	G01	HQP021384	NM_003738	PTCH2
QG064-02	G02	HQP021313	NM_001753	CAV1
QG064-02	G03	HQP021023	NM_003617	RGS5
QG064-02	G04	HQP020899	NM_000388	CASR
QG064-02	G05	HQP020707	NM_003579	RAD54L
QG064-02	G06	HQP020618	NM_032316	NICN1
QG064-02	G07	HQP020418	NM_001004426	PLA2G6
QG064-02	G08	HQP020297	NM_004346	CASP3
QG064-02	G09	HQP020136	NM_003507	FZD7
QG064-02	G10	HQP020116	NM_004655	AXIN2
QG064-02	G11	HQP020040	NM_006534	NCOA3
QG064-02	G12	HQP019829	NM_030782	CLPTM1L
QG064-02	H01	HGDC		
QG064-02	H02	HGDC		
QG064-02	H03	HQP006940	NM_002046	GAPDH
QG064-02	H04	HQP016381	NM_001101	ACTB
QG064-02	H05	HQP015171	NM_004048	B2M
QG064-02	H06	HQP006171	NM_012423	RPL13A
QG064-02	H07	HQP009026	NM_000194	HPRT1
QG064-02	H08	HQP054253	NR_003286	RN18S1
QG064-02	H09	RT		
QG064-02	H10	RT		
QG064-02	H11	PCR		
QG064-02	H12	PCR		
QG064-03	A01	HQP019768	NM_130445	COL18A1
QG064-03	A02	HQP019764	NM_030578	MGC4093
QG064-03	A03	HQP019689	NM_001042529	COASY
QG064-03	A04	HQP019452	NM_001742	CALCR
QG064-03	A05	HQP018747	NM_013309	SLC30A4
QG064-03	A06	HQP018739	NM_006300	ZNF230
QG064-03	A07	HQP018550	NM_001079539	XBP1
QG064-03	A08	HQP018545	NM_000553	WRN
QG064-03	A09	HQP018488	NM_003379	VIL2
QG064-03	A10	HQP018466	NM_080682	VCAM1
QG064-03	A11	HQP018404	NM_003356	UCP3
QG064-03	A12	HQP018343	NM_000372	TYR
QG064-03	B01	HQP018269	NM_006292	TSG101
QG064-03	B02	HQP018265	NM_000548	TSC2

Product Data Sheet

QG064-03	B03	HQP018060	NM_000359	TGM1
QG064-03	B04	HQP018035	NM_003225	TFF1
QG064-03	B05	HQP018016	NM_005652	TERF2
QG064-03	B06	HQP018014	NM_003218	TERF1
QG064-03	B07	HQP018013	NM_007110	TEP1
QG064-03	B08	HQP017996	NM_003211	TDG
QG064-03	B09	HQP017794	NM_000455	STK11
QG064-03	B10	HQP017764	NM_007315	STAT1
QG064-03	B11	HQP017745	NM_001051	SSTR3
QG064-03	B12	HQP017702	NM_001725	BPI
QG064-03	C01	HQP017615	NM_000454	SOD1
QG064-03	C02	HQP017563	NM_005631	SMO
QG064-03	C03	HQP017367	NM_003040	SLC4A2
QG064-03	C04	HQP017303	NM_003037	SLAMF1
QG064-03	C05	HQP016980	NM_003019	SFTPD
QG064-03	C06	HQP016819	NM_005410	SEPP1
QG064-03	C07	HQP016801	NM_022162	NOD2
QG064-03	C08	HQP016744	NM_000450	SELE
QG064-03	C09	HQP016670	NM_001033886	CXCL12
QG064-03	C10	HQP016634	NM_001713	BHMT
QG064-03	C11	HQP016626	NM_002985	CCL5
QG064-03	C12	HQP016527	NM_021976	RXR
QG064-03	D01	HQP016526	NM_002957	RXRA
QG064-03	D02	HQP016418	NM_004327	BCR
QG064-03	D03	HQP016391	NM_002947	RPA3
QG064-03	D04	HQP016390	NM_002946	RPA2
QG064-03	D05	HQP016389	NM_002945	RPA1
QG064-03	D06	HQP016380	NM_002944	ROS1
QG064-03	D07	HQP016303	NM_021133	RNASEL
QG064-03	D08	HQP016276	NM_005614	RHEB
QG064-03	D09	HQP016252	NM_002923	RGS2
QG064-03	D10	HQP016231	NM_181578	RFC5
QG064-03	D11	HQP016227	NM_002915	RFC3
QG064-03	D12	HQP016225	NM_002914	RFC2
QG064-03	E01	HQP016224	NM_002913	RFC1
QG064-03	E02	HQP016223	NM_002912	REV3L
QG064-03	E03	HQP016168	NM_021629	GNB4
QG064-03	E04	HQP016079	NM_002876	RAD51C
QG064-03	E05	HQP016064	NM_018890	RAC1
QG064-03	E06	HQP016003	NM_021213	PCTP
QG064-03	E07	HQP015946	NM_000465	BARD1
QG064-03	E08	HQP015917	NM_001188	BAK1
QG064-03	E09	HQP015771	NM_020884	MYH7B
QG064-03	E10	HQP015596	NM_000962	PTGS1
QG064-03	E11	HQP015535	NM_000314	PTEN
QG064-03	E12	HQP014808	NM_018657	MYNN
QG064-03	F01	HQP014564	NM_017902	HIF1AN
QG064-03	F02	HQP014292	NM_018315	FBXW7



Product Data Sheet

QG064-03	F03	HQP014089	NM_018081	WDR79
QG064-03	F04	HQP013581	NM_019077	UGT1A7
QG064-03	F05	HQP013473	NM_000446	PON1
QG064-03	F06	HQP013459	NM_005035	POLRMT
QG064-03	F07	HQP013454	NM_005034	POLR2K
QG064-03	F08	HQP013452	NM_006234	POLR2J
QG064-03	F09	HQP013444	NM_021974	POLR2F
QG064-03	F10	HQP013443	NM_002695	POLR2E
QG064-03	F11	HQP013438	NM_004805	POLR2D
QG064-03	F12	HQP013434	NM_032940	POLR2C
QG064-03	G01	HQP013431	NM_000938	POLR2B
QG064-03	G02	HQP013428	NM_002693	POLG
QG064-03	G03	HQP013427	NM_002692	POLE2
QG064-03	G04	HQP013425	NM_006231	POLE
QG064-03	G05	HQP013424	NM_006230	POLD2
QG064-03	G06	HQP013262	NM_005030	PLK1
QG064-03	G07	HQP013204	NM_002658	PLAU
QG064-03	G08	HQP013194	NM_000300	PLA2G2A
QG064-03	G09	HQP013152	NM_002648	PIM1
QG064-03	G10	HQP013104	NM_002634	PHB
QG064-03	G11	HQP013100	NM_000927	ABCB1
QG064-03	G12	HQP013099	NM_000926	PGR
QG064-03	H01	HGDC		
QG064-03	H02	HGDC		
QG064-03	H03	HQP006940	NM_002046	GAPDH
QG064-03	H04	HQP016381	NM_001101	ACTB
QG064-03	H05	HQP015171	NM_004048	B2M
QG064-03	H06	HQP006171	NM_012423	RPL13A
QG064-03	H07	HQP009026	NM_000194	HPRT1
QG064-03	H08	HQP054253	NR_003286	RN18S1
QG064-03	H09	RT		
QG064-03	H10	RT		
QG064-03	H11	PCR		
QG064-03	H12	PCR		
QG064-04	A01	HQP012957	NM_006208	ENPP1
QG064-04	A02	HQP012725	NM_016218	POLK
QG064-04	A03	HQP012482	NM_001005783	HAO2
QG064-04	A04	HQP012100	NM_002562	P2RX7
QG064-04	A05	HQP012062	NM_001008505	OPRM1
QG064-04	A06	HQP012008	NM_000275	OCA2
QG064-04	A07	HQP011914	NM_002524	NRAS
QG064-04	A08	HQP011868	NM_000603	NOS3
QG064-04	A09	HQP011858	NM_000904	NQO2
QG064-04	A10	HQP011852	NM_001678	ATP1B2
QG064-04	A11	HQP011836	NM_004148	NINJ1
QG064-04	A12	HQP011813	NM_004556	NFKBIE
QG064-04	B01	HQP011693	NM_000433	NCF2
QG064-04	B02	HQP011582	NM_002462	MX1

Product Data Sheet

QG064-04	B03	HQP011320	NM_005590	MRE11A
QG064-04	B04	HQP011304	NM_001015052	MPG
QG064-04	B05	HQP011182	NM_005371	METTL1
QG064-04	B06	HQP011181	NM_000245	MET
QG064-04	B07	HQP011179	NM_002402	MEST
QG064-04	B08	HQP011007	NM_000240	MAOA
QG064-04	B09	HQP010885	NM_002337	LRPAP1
QG064-04	B10	HQP010878	NM_002335	LRP5
QG064-04	B11	HQP010877	NM_002336	LRP6
QG064-04	B12	HQP010847	NM_000237	LPL
QG064-04	C01	HQP010626	NM_000236	LIPC
QG064-04	C02	HQP010613	NM_002312	LIG4
QG064-04	C03	HQP010610	NM_001175	ARHGDI1B
QG064-04	C04	HQP010584	NM_002303	LEPR
QG064-04	C05	HQP010581	NM_000230	LEP
QG064-04	C06	HQP010577	NM_000527	LDLR
QG064-04	C07	HQP010564	NM_000229	LCAT
QG064-04	C08	HQP010095	NM_013289	KIR3DL1
QG064-04	C09	HQP010008	NM_198686	RAB15
QG064-04	C10	HQP009851	NM_000215	JAK3
QG064-04	C11	HQP009788	NM_005544	IRS1
QG064-04	C12	HQP009780	NM_001571	IRF3
QG064-04	D01	HQP009778	NM_002198	IRF1
QG064-04	D02	HQP009764	NM_000208	INSR
QG064-04	D03	HQP009710	NM_002189	IL15RA
QG064-04	D04	HQP009708	NM_000585	IL15
QG064-04	D05	HQP009697	NM_002188	IL13
QG064-04	D06	HQP009693	NM_002187	IL12B
QG064-04	D07	HQP009679	NM_000634	IL8RA
QG064-04	D08	HQP009677	NM_002185	IL7R
QG064-04	D09	HQP009672	NM_000565	IL6R
QG064-04	D10	HQP009642	NM_000877	IL1R1
QG064-04	D11	HQP009556	NM_000041	APOE
QG064-04	D12	HQP009555	NM_002178	IGFBP6
QG064-04	E01	HQP009552	NM_182632	SLC6A18
QG064-04	E02	HQP009551	NM_000599	IGFBP5
QG064-04	E03	HQP009544	NM_000598	IGFBP3
QG064-04	E04	HQP009541	NM_000597	IGFBP2
QG064-04	E05	HQP009539	NM_000596	IGFBP1
QG064-04	E06	HQP009529	NM_000612	IGF2
QG064-04	E07	HQP009523	NM_000875	IGF1R
QG064-04	E08	HQP009518	NM_000618	IGF1
QG064-04	E09	HQP009472	NM_005534	IFNGR2
QG064-04	E10	HQP009469	NM_000416	IFNGR1
QG064-04	E11	HQP009460	NM_000874	IFNAR2
QG064-04	E12	HQP009385	NM_005532	IFI27
QG064-04	F01	HQP009171	NM_000482	APOA4
QG064-04	F02	HQP009139	NM_001643	APOA2

Product Data Sheet

QG064-04	F03	HQP009138	NM_004507	HUS1
QG064-04	F04	HQP009117	NM_000864	HTR1D
QG064-04	F05	HQP009116	NM_000863	HTR1B
QG064-04	F06	HQP009072	NM_001166	BIRC2
QG064-04	F07	HQP009067	NM_000414	HSD17B4
QG064-04	F08	HQP009066	NM_002153	HSD17B2
QG064-04	F09	HQP009064	NM_000413	HSD17B1
QG064-04	F10	HQP009060	NM_000198	HSD3B2
QG064-04	F11	HQP009059	NM_000862	HSD3B1
QG064-04	F12	HQP008934	NM_001160	APAF1
QG064-04	G01	HQP008891	NM_000859	HMGCR
QG064-04	G02	HQP008832	NM_181054	HIF1A
QG064-04	G03	HQP008757	NM_000410	HFE
QG064-04	G04	HQP008718	NM_000182	HADHA
QG064-04	G05	HQP008659	NM_014579	SLC39A2
QG064-04	G06	HQP008605	NM_013347	RPA4
QG064-04	G07	HQP008490	NM_001513	GSTZ1
QG064-04	G08	HQP008479	NM_001512	GSTA4
QG064-04	G09	HQP008469	NM_002093	GSK3B
QG064-04	G10	HQP008464	NM_005314	GRPR
QG064-04	G11	HQP008285	NM_002085	GPX4
QG064-04	G12	HQP008282	NM_002084	GPX3
QG064-04	H01	HGDC		
QG064-04	H02	HGDC		
QG064-04	H03	HQP006940	NM_002046	GAPDH
QG064-04	H04	HQP016381	NM_001101	ACTB
QG064-04	H05	HQP015171	NM_004048	B2M
QG064-04	H06	HQP006171	NM_012423	RPL13A
QG064-04	H07	HQP009026	NM_000194	HPRT1
QG064-04	H08	HQP054253	NR_003286	RN18S1
QG064-04	H09	RT		
QG064-04	H10	RT		
QG064-04	H11	PCR		
QG064-04	H12	PCR		
QG064-05	A01	HQP008281	NM_002083	GPX2
QG064-05	A02	HQP007767	NM_002075	GNB3
QG064-05	A03	HQP007755	NM_000516	GNAS
QG064-05	A04	HQP007689	NM_013274	POLL
QG064-05	A05	HQP007643	NM_014475	DHDH
QG064-05	A06	HQP007304	NM_012419	RGS17
QG064-05	A07	HQP007230	NM_000583	GC
QG064-05	A08	HQP007224	NM_014365	HSPB8
QG064-05	A09	HQP006945	NM_015515	KRT23
QG064-05	A10	HQP006855	NM_001042594	POT1
QG064-05	A11	HQP006733	NM_012134	LMOD1
QG064-05	A12	HQP006583	NM_178510	ANKK1
QG064-05	B01	HQP006547	NM_001469	XRCC6
QG064-05	B02	HQP006463	NM_004479	FUT7

Product Data Sheet

QG064-05	B03	HQP006425	NM_001140	ALOX15
QG064-05	B04	HQP006359	NM_000698	ALOX5
QG064-05	B05	HQP006356	NM_000697	ALOX12
QG064-05	B06	HQP006305	NM_002689	POLA2
QG064-05	B07	HQP006261	NM_014324	AMACR
QG064-05	B08	HQP006189	NM_012429	SEC14L2
QG064-05	B09	HQP006188	NM_005252	FOS
QG064-05	B10	HQP006120	NM_018672	ABCA5
QG064-05	B11	HQP006119	NM_080284	ABCA6
QG064-05	B12	HQP005629	NM_001453	FOXC1
QG064-05	C01	HQP005440	NM_022963	FGFR4
QG064-05	C02	HQP005280	NM_021642	FCGR2A
QG064-05	C03	HQP005087	NM_000690	ALDH2
QG064-05	C04	HQP005084	NM_001018115	FANCD2
QG064-05	C05	HQP005083	NM_000136	FANCC
QG064-05	C06	HQP005002	NM_001437	ESR2
QG064-05	C07	HQP004998	NM_000125	ESR1
QG064-05	C08	HQP004991	NM_001014431	AKT1
QG064-05	C09	HQP004849	NM_152686	DNAJC18
QG064-05	C10	HQP004727	NM_005502	ABCA1
QG064-05	C11	HQP004598	NM_001406	EFNB3
QG064-05	C12	HQP004557	NM_001955	EDN1
QG064-05	D01	HQP004450	NM_000797	DRD4
QG064-05	D02	HQP004445	NM_000794	DRD1
QG064-05	D03	HQP004309	NM_000791	DHFR
QG064-05	D04	HQP004114	NM_000107	DDB2
QG064-05	D05	HQP004081	NM_000789	ACE
QG064-05	D06	HQP003932	NM_152587	C11orf65
QG064-05	D07	HQP003917	NM_000782	CYP24A1
QG064-05	D08	HQP003904	NM_000103	CYP19A1
QG064-05	D09	HQP003888	NM_000102	CYP17A1
QG064-05	D10	HQP003841	NM_000777	CYP3A5
QG064-05	D11	HQP003804	NM_000766	CYP2A13
QG064-05	D12	HQP003741	NM_000397	CYBB
QG064-05	E01	HQP003691	NM_001337	CX3CR1
QG064-05	E02	HQP003620	NM_004390	CTSH
QG064-05	E03	HQP003539	NM_001904	CTNNB1
QG064-05	E04	HQP003499	NM_001037631	CTLA4
QG064-05	E05	HQP003395	NM_001033521	CSTF1
QG064-05	E06	HQP003173	NM_000759	CSF3
QG064-05	E07	HQP003159	NM_000758	CSF2
QG064-05	E08	HQP003158	NM_005211	CSF1R
QG064-05	E09	HQP002881	NM_000651	CR1
QG064-05	E10	HQP002210	NM_000579	CCR5
QG064-05	E11	HQP001862	NM_000082	ERCC8
QG064-05	E12	HQP001632	NM_000743	CHRNA3
QG064-05	F01	HQP001449	NM_133491	SAT2
QG064-05	F02	HQP001416	NM_007204	DDX20

Product Data Sheet

QG064-05	F03	HQP001399	NM_007195	POLI
QG064-05	F04	HQP001396	NM_001005735	CHEK2
QG064-05	F05	HQP001282	NM_001274	CHEK1
QG064-05	F06	HQP001262	NM_001818	AKR1C4
QG064-05	F07	HQP000970	NM_006663	PPP1R13L
QG064-05	F08	HQP000963	NM_012190	ALDH1L1
QG064-05	F09	HQP000956	NM_000735	CGA
QG064-05	F10	HQP000932	NM_006636	MTHFD2
QG064-05	F11	HQP000847	NM_006591	POLD3
QG064-05	F12	HQP000708	NM_006440	TXNRD2
QG064-05	G01	HQP000554	NM_006342	TACC3
QG064-05	G02	HQP000405	NM_006066	AKR1A1
QG064-05	G03	HQP000356	NM_000076	CDKN1C
QG064-05	G04	HQP000285	NM_001799	CDK7
QG064-05	G05	HQP000274	NM_001259	CDK6
QG064-05	G06	HQP000261	NM_004935	CDK5
QG064-05	G07	HQP000145	NM_005732	RAD50
QG064-05	G08	HQP000134	NM_012225	NUBP2
QG064-05	G09	HQP000001	NM_005465	AKT3
QG064-05	G10	HQP054694	NM_020945	WDFY4
QG064-05	G11	HQP017991	NM_022171	TCTA
QG064-05	G12	HQP017618	NM_003102	SOD3
QG064-05	H01	HGDC		
QG064-05	H02	HGDC		
QG064-05	H03	HQP006940	NM_002046	GAPDH
QG064-05	H04	HQP016381	NM_001101	ACTB
QG064-05	H05	HQP015171	NM_004048	B2M
QG064-05	H06	HQP006171	NM_012423	RPL13A
QG064-05	H07	HQP009026	NM_000194	HPRT1
QG064-05	H08	HQP054253	NR_003286	RN18S1
QG064-05	H09	RT		
QG064-05	H10	RT		
QG064-05	H11	PCR		
QG064-05	H12	PCR		

### Limited Use License

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