

## ExProfile™ Human Colorectal Cancer Gene qPCR Array

For focused group profiling of human colorectal cancer related genes expression

Cat. No. QG065-A (4 x 96-well plate, Format A)

Cat. No. QG065-B (4 x 96-well plate, Format B)

Cat. No. QG065-C (4 x 96-well plate, Format C)

Cat. No. QG065-D (4 x 96-well plate, Format D)

Cat. No. QG065-E (4 x 96-well plate, Format E)

Available as 1 set or 6 sets. Each set contains 336 unique gene primers deposited in four 96-well plates.

### Introduction

The ExProfile human colorectal cancer gene qPCR array profiles 336 human genes to aberrantly expressed human genes involved in human colorectal 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 colorectal cancer gene array allows researchers to study the cancer-related genes to gain understanding of their roles in colorectal cancer pathogenesis.

- QG065 plate 01: 84 unique gene PCR primer pairs
- QG065 plate 02: 84 unique gene PCR primer pairs
- QG065 plate 03: 84 unique gene PCR primer pairs
- QG065 plate 04: 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
<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	KRAS	MTHFR	MLH1	MSH2	TP53	BRAF	APC	TYMS	MSH6	NAT2	XRCC1	VDR
B	CYP1A1	ERCC2	NAT1	CCND1	PTGS2	PIK3CA	MTR	ABCB1	ERCC1	CYP2E1	CDH1	EPHX1
C	NQO1	IL6	CYP2C9	OGG1	VEGFA	IGF1	EGFR	TGFB1	PPARG	MTRR	SMAD7	TNF
D	SULT1A1	ALDH2	DPYD	CYP1B1	CHEK2	TGFBR1	XRCC3	NOD2	MMP1	APOE	UGT1A7	PMS2
E	IGFBP3	MDM2	IL10	HFE	RHPN2	UGT1A6	IRS1	IL8	IL1B	GSTM3	CYP2D6	ADH1C
F	BCR	PTGS1	NFKB1	MMP9	MMP3	EGF	CTNNB1	CDKN2A	ERCC5	CASR	MMP2	ESR1
G	ADH1B	TP73	UGT1A9	XPC	BRCA2	IL4	IL1RN	APEX1	HRAS	HIF1A	GH1	FCGR2A
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG065 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	CTLA4	PARP1	NUDT1	DNMT3B	EXO1	BMP4	ADIPOQ	CBS	EIF3S3	CASP8	SOD2	PTEN
B	ODC1	NOS3	MSH3	LEPR	LEP	LCT	ALOX12	ESR2	ACE	CYP19A1	CYP17A1	CYP3A5
C	CYP2C19	COMT	CES2	AXIN1	XRCC4	TCN2	SLC19A1	SLC10A2	SEPP1	CXCL12	RETN	PLAU
D	SERPINE1	NRAS	NOS2A	NBN	SMAD4	AR	IL4R	IGF1R	ALOX5	FOLH1	F5	AKT1
E	ERCC6	ERCC4	AHR	DRD2	CD36	CD14	CCNH	AXIN2	XRCC5	XRCC2	UMPS	SFRP2
F	BLM	SELE	RXRA	RAD23B	PTPRJ	PTPN13	PPARD	POU2AF1	POLB	PMS1	PLA2G2A	PGR
G	GHRL	NFE2L2	MYC	ABCC1	MC4R	ITGB3	INSR	INS	FASLG	IL1A	HSD17B2	HPGD
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG065 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	GPX4	FRZB	FRAP1	FMO3	EPHB2	DHFR	DCC	CYP27B1	CYP24A1	CYP2C8	CHEK1	PPP1R13L
B	CDK7	ADA	MYCL1	CYP7A1	SELS	SLC6A4	CASP8AP2	SLC23A1	CDA	HDAC9	GDF15	ATG5
C	CD86	CD28	ACVR2A	DLG5	ATG12	BTRC	MTMR3	NR1I2	WISP1	GGH	TNFRSF10A	TNFRSF10B
D	RIPK1	ABCC3	CAV1	CDC14A	SSH2	PTPN5	LTBP4	CASP7	CASP6	CASP3	CLPTM1L	RLM1
E	EHMT1	GALNT12	CALB2	CXCR4	PRDM2	CA6	WRN	WNT2B	VHL	UGT2B7	TXNRD1	TTK
F	TSC2	TSC1	TP53BP1	TLE1	TCF7	TCF1	BMPR2	SLC22A5	SLC22A4	MAP2K4	SELP	CX3CL1
G	BHMT	CCL2	CCL1	RPS6KB1	RPS6KA2	RPS6KA1	RNASEL	RFC1	GNB4	RB1	RAD52	BARD1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG065 plate 03

## Product Data Sheet

	1	2	3	4	5	6	7	8	9	10	11	12
A	PTPRS	PTGER4	PTGER2	PTCH1	PSMB9	RAD18	RPRM	PRNP	TDP1	PRKAA1	FAM46A	LAPTM4B
B	C14orf103	PPP1R3A	TRPM7	ATR	PON1	POMC	SSH1	PIK3R1	SERPINA1	SERPINF1	UPB1	NLK
C	PRKAG2	CDC40	PCNA	SH3GLB1	ADIPOR1	IL22	RRM2B	FURIN	NT5E	ATP2A3	ATP2A2	NME1
D	NFKBIA	ATIC	MTM1	MPO	MMP12	CIITA	MGST1	MAT2A	MANBA	SMAD3	SMAD2	ARNT
E	LTBP1	LIPC	LIG4	LIG3	LIG1	KDR	IL18	IL12B	IL12A	IKBKB	IGFBP1	C11orf53
F	BIRC5	HSD17B3	HMGCR	GSTM2	GSTA4	GSTA2	DDX25	NR3C1	GPX3	GPX2	ATG9B	GNAS
G	GLI3	DKK3	AMPD1	AMD1	GART	PYGO1	BAMBI	RAD54B	AMACR	FPGS	FOLR1	FLT1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG065 plate 04

- **Gene primer pairs:** 84 wells (A row to G row) are designated for a real-time PCR assay for genes (see the primer list).
- **HK1-6:** Six pre-deposited housekeeping gene (HK1-6) primer pairs, which can be used as endogenous positive controls as well as for array normalization.
- **GDC:** Genomic DNA controls, which can be used to specifically detect genomic DNA contamination with a high level of sensitivity.
- **RT:** Spike-in reverse transcription controls, which can be used to monitor the efficiency of the RT reactions. These pre-deposited primer pairs specifically amplify the cDNA template reversed transcribed from the spike-in control RNA in the sample.
- **PCR:** Positive PCR controls, which are used to verify the PCR efficiency by amplifying the pre-deposited DNA template with its specific pre-deposited primer pairs.

### Gene primer list

Plate	Position	Catalog No. of Primer	Accession No. of Gene	Symbol
QG065-01	A01	HQP010133	NM_004985	KRAS
QG065-01	A02	HQP011547	NM_005957	MTHFR
QG065-01	A03	HQP011235	NM_000249	MLH1
QG065-01	A04	HQP011491	NM_000251	MSH2
QG065-01	A05	HQP018175	NM_000546	TP53
QG065-01	A06	HQP017733	NM_004333	BRAF
QG065-01	A07	HQP009024	NM_000038	APC
QG065-01	A08	HQP018342	NM_001071	TYMS
QG065-01	A09	HQP008493	NM_000179	MSH6
QG065-01	A10	HQP001136	NM_000015	NAT2
QG065-01	A11	HQP018562	NM_006297	XRCC1
QG065-01	A12	HQP018474	NM_000376	VDR
QG065-01	B01	HQP003772	NM_000499	CYP1A1
QG065-01	B02	HQP004976	NM_000400	ERCC2
QG065-01	B03	HQP023467	NM_000662	NAT1
QG065-01	B04	HQP016204	NM_053056	CCND1
QG065-01	B05	HQP015598	NM_000963	PTGS2
QG065-01	B06	HQP013150	NM_006218	PIK3CA
QG065-01	B07	HQP011554	NM_000254	MTR
QG065-01	B08	HQP013100	NM_000927	ABCB1
QG065-01	B09	HQP004975	NM_202001	ERCC1
QG065-01	B10	HQP003817	NM_000773	CYP2E1

## Product Data Sheet

QG065-01	B11	HQP023466	NM_004360	CDH1
QG065-01	B12	HQP004948	NM_000120	EPHX1
QG065-01	C01	HQP004317	NM_000903	NQO1
QG065-01	C02	HQP009670	NM_000600	IL6
QG065-01	C03	HQP003811	NM_000771	CYP2C9
QG065-01	C04	HQP012021	NM_002542	OGG1
QG065-01	C05	HQP018475	NM_001025366	VEGFA
QG065-01	C06	HQP009518	NM_000618	IGF1
QG065-01	C07	HQP004605	NM_005228	EGFR
QG065-01	C08	HQP018044	NM_000660	TGFB1
QG065-01	C09	HQP013633	NM_005037	PPARG
QG065-01	C10	HQP011555	NM_002454	MTRR
QG065-01	C11	HQP010966	NM_005904	SMAD7
QG065-01	C12	HQP018141	NM_000594	TNF
QG065-01	D01	HQP017811	NM_177536	SULT1A1
QG065-01	D02	HQP005087	NM_000690	ALDH2
QG065-01	D03	HQP004439	NM_000110	DPYD
QG065-01	D04	HQP003775	NM_000104	CYP1B1
QG065-01	D05	HQP001396	NM_001005735	CHEK2
QG065-01	D06	HQP054057	BC071181	TGFBR1
QG065-01	D07	HQP018564	NM_005432	XRCC3
QG065-01	D08	HQP016801	NM_022162	NOD2
QG065-01	D09	HQP011255	NM_002421	MMP1
QG065-01	D10	HQP009556	NM_000041	APOE
QG065-01	D11	HQP013581	NM_019077	UGT1A7
QG065-01	D12	HQP013352	NM_000535	PMS2
QG065-01	E01	HQP009544	NM_000598	IGFBP3
QG065-01	E02	HQP011135	NM_002392	MDM2
QG065-01	E03	HQP009685	NM_000572	IL10
QG065-01	E04	HQP008757	NM_000410	HFE
QG065-01	E05	HQP021215	NM_033103	RHPN2
QG065-01	E06	HQP013583	NM_205862	UGT1A6
QG065-01	E07	HQP009788	NM_005544	IRS1
QG065-01	E08	HQP009678	NM_000584	IL8
QG065-01	E09	HQP009641	NM_000576	IL1B
QG065-01	E10	HQP008483	NM_000849	GSTM3
QG065-01	E11	HQP003814	NM_000106	CYP2D6
QG065-01	E12	HQP002413	NM_000669	ADH1C
QG065-01	F01	HQP016418	NM_004327	BCR
QG065-01	F02	HQP015596	NM_000962	PTGS1
QG065-01	F03	HQP011807	NM_003998	NFKB1
QG065-01	F04	HQP011263	NM_004994	MMP9
QG065-01	F05	HQP011257	NM_002422	MMP3
QG065-01	F06	HQP004599	NM_001963	EGF
QG065-01	F07	HQP003539	NM_001904	CTNNB1
QG065-01	F08	HQP000370	NM_058195	CDKN2A
QG065-01	F09	HQP004985	NM_000123	ERCC5
QG065-01	F10	HQP020899	NM_000388	CASR

## Product Data Sheet

QG065-01	F11	HQP011256	NM_004530	MMP2
QG065-01	F12	HQP004998	NM_000125	ESR1
QG065-01	G01	HQP002331	NM_000668	ADH1B
QG065-01	G02	HQP018180	NM_005427	TP73
QG065-01	G03	HQP013597	NM_021027	UGT1A9
QG065-01	G04	HQP018556	NM_004628	XPC
QG065-01	G05	HQP017753	NM_000059	BRCA2
QG065-01	G06	HQP009662	NM_000589	IL4
QG065-01	G07	HQP009645	NM_000577	IL1RN
QG065-01	G08	HQP009061	NM_001641	APEX1
QG065-01	G09	HQP009036	NM_005343	HRAS
QG065-01	G10	HQP008832	NM_181054	HIF1A
QG065-01	G11	HQP007385	NM_000515	GH1
QG065-01	G12	HQP005280	NM_021642	FCGR2A
QG065-01	H01	HGDC		
QG065-01	H02	HGDC		
QG065-01	H03	HQP006940	NM_002046	GAPDH
QG065-01	H04	HQP016381	NM_001101	ACTB
QG065-01	H05	HQP015171	NM_004048	B2M
QG065-01	H06	HQP006171	NM_012423	RPL13A
QG065-01	H07	HQP009026	NM_000194	HPRT1
QG065-01	H08	HQP054253	NR_003286	RN18S1
QG065-01	H09	RT		
QG065-01	H10	RT		
QG065-01	H11	PCR		
QG065-01	H12	PCR		
QG065-02	A01	HQP003499	NM_001037631	CTLA4
QG065-02	A02	HQP003120	NM_001618	PARP1
QG065-02	A03	HQP011544	NM_002452	NUDT1
QG065-02	A04	HQP004414	NM_006892	DNMT3B
QG065-02	A05	HQP054030	NM_130398	EXO1
QG065-02	A06	HQP053910	NM_130851	BMP4
QG065-02	A07	HQP022625	NM_004797	ADIPOQ
QG065-02	A08	HQP021518	NM_000071	CBS
QG065-02	A09	HQP021417	NM_003756	EIF3S3
QG065-02	A10	HQP018966	NM_001080124	CASP8
QG065-02	A11	HQP017616	NM_000636	SOD2
QG065-02	A12	HQP015535	NM_000314	PTEN
QG065-02	B01	HQP012013	NM_002539	ODC1
QG065-02	B02	HQP011868	NM_000603	NOS3
QG065-02	B03	HQP011492	NM_002439	MSH3
QG065-02	B04	HQP010584	NM_002303	LEPR
QG065-02	B05	HQP010581	NM_000230	LEP
QG065-02	B06	HQP010570	NM_002299	LCT
QG065-02	B07	HQP006356	NM_000697	ALOX12
QG065-02	B08	HQP005002	NM_001437	ESR2
QG065-02	B09	HQP004081	NM_000789	ACE
QG065-02	B10	HQP003904	NM_000103	CYP19A1

## Product Data Sheet

QG065-02	B11	HQP003888	NM_000102	CYP17A1
QG065-02	B12	HQP003841	NM_000777	CYP3A5
QG065-02	C01	HQP003809	NM_000769	CYP2C19
QG065-02	C02	HQP002671	NM_000754	COMT
QG065-02	C03	HQP021586	NM_198061	CES2
QG065-02	C04	HQP020114	NM_003502	AXIN1
QG065-02	C05	HQP018565	NM_003401	XRCC4
QG065-02	C06	HQP017979	NM_000355	TCN2
QG065-02	C07	HQP017482	NM_194255	SLC19A1
QG065-02	C08	HQP017462	NM_000452	SLC10A2
QG065-02	C09	HQP016819	NM_005410	SEPP1
QG065-02	C10	HQP016670	NM_001033886	CXCL12
QG065-02	C11	HQP015160	NM_020415	RETN
QG065-02	C12	HQP013204	NM_002658	PLAU
QG065-02	D01	HQP012154	NM_000602	SERPINE1
QG065-02	D02	HQP011914	NM_002524	NRAS
QG065-02	D03	HQP011866	NM_000625	NOS2A
QG065-02	D04	HQP011687	NM_002485	NBN
QG065-02	D05	HQP010961	NM_005359	SMAD4
QG065-02	D06	HQP009801	NM_000044	AR
QG065-02	D07	HQP009664	NM_000418	IL4R
QG065-02	D08	HQP009523	NM_000875	IGF1R
QG065-02	D09	HQP006359	NM_000698	ALOX5
QG065-02	D10	HQP006129	NM_001014986	FOLH1
QG065-02	D11	HQP005058	NM_000130	F5
QG065-02	D12	HQP004991	NM_001014431	AKT1
QG065-02	E01	HQP004986	NM_000124	ERCC6
QG065-02	E02	HQP004984	NM_005236	ERCC4
QG065-02	E03	HQP004658	NM_001621	AHR
QG065-02	E04	HQP004446	NM_000795	DRD2
QG065-02	E05	HQP022821	NM_000072	CD36
QG065-02	E06	HQP022490	NM_000591	CD14
QG065-02	E07	HQP021906	NM_001239	CCNH
QG065-02	E08	HQP020116	NM_004655	AXIN2
QG065-02	E09	HQP018568	NM_021141	XRCC5
QG065-02	E10	HQP018563	NM_005431	XRCC2
QG065-02	E11	HQP018425	NM_000373	UMPS
QG065-02	E12	HQP016874	NM_003013	SFRP2
QG065-02	F01	HQP016844	NM_000057	BLM
QG065-02	F02	HQP016744	NM_000450	SELE
QG065-02	F03	HQP016526	NM_002957	RXRA
QG065-02	F04	HQP016076	NM_002874	RAD23B
QG065-02	F05	HQP015925	NM_002843	PTPRJ
QG065-02	F06	HQP015897	NM_006264	PTPN13
QG065-02	F07	HQP013627	NM_006238	PPARD
QG065-02	F08	HQP013528	NM_006235	POU2AF1
QG065-02	F09	HQP013422	NM_002690	POLB
QG065-02	F10	HQP013312	NM_000534	PMS1

## Product Data Sheet

QG065-02	F11	HQP013194	NM_000300	PLA2G2A
QG065-02	F12	HQP013099	NM_000926	PGR
QG065-02	G01	HQP012996	NM_016362	GHRL
QG065-02	G02	HQP011800	NM_006164	NFE2L2
QG065-02	G03	HQP011597	NM_002467	MYC
QG065-02	G04	HQP011325	NM_019899	ABCC1
QG065-02	G05	HQP011097	NM_005912	MC4R
QG065-02	G06	HQP009818	NM_000212	ITGB3
QG065-02	G07	HQP009764	NM_000208	INSR
QG065-02	G08	HQP009749	NM_000207	INS
QG065-02	G09	HQP009671	NM_000639	FASLG
QG065-02	G10	HQP009640	NM_000575	IL1A
QG065-02	G11	HQP009066	NM_002153	HSD17B2
QG065-02	G12	HQP009022	NM_000860	HPGD
QG065-02	H01	HGDC		
QG065-02	H02	HGDC		
QG065-02	H03	HQP006940	NM_002046	GAPDH
QG065-02	H04	HQP016381	NM_001101	ACTB
QG065-02	H05	HQP015171	NM_004048	B2M
QG065-02	H06	HQP006171	NM_012423	RPL13A
QG065-02	H07	HQP009026	NM_000194	HPRT1
QG065-02	H08	HQP054253	NR_003286	RN18S1
QG065-02	H09	RT		
QG065-02	H10	RT		
QG065-02	H11	PCR		
QG065-02	H12	PCR		
QG065-03	A01	HQP008285	NM_002085	GPX4
QG065-03	A02	HQP006431	NM_001463	FRZB
QG065-03	A03	HQP006426	NM_004958	FRAP1
QG065-03	A04	HQP005959	NM_001002294	FMO3
QG065-03	A05	HQP004940	NM_017449	EPHB2
QG065-03	A06	HQP004309	NM_000791	DHFR
QG065-03	A07	HQP004053	NM_005215	DCC
QG065-03	A08	HQP003923	NM_000785	CYP27B1
QG065-03	A09	HQP003917	NM_000782	CYP24A1
QG065-03	A10	HQP003810	NM_000770	CYP2C8
QG065-03	A11	HQP001282	NM_001274	CHEK1
QG065-03	A12	HQP000970	NM_006663	PPP1R13L
QG065-03	B01	HQP000285	NM_001799	CDK7
QG065-03	B02	HQP000132	NM_000022	ADA
QG065-03	B03	HQP011601	NM_005376	MYCL1
QG065-03	B04	HQP003859	NM_000780	CYP7A1
QG065-03	B05	HQP054043	NM_203472	SELS
QG065-03	B06	HQP053962	NM_001045	SLC6A4
QG065-03	B07	HQP023464	NM_012115	CASP8AP2
QG065-03	B08	HQP023414	NM_005847	SLC23A1
QG065-03	B09	HQP023203	NM_001785	CDA
QG065-03	B10	HQP023134	NM_014707	HDAC9

## Product Data Sheet

QG065-03	B11	HQP022853	NM_004864	GDF15
QG065-03	B12	HQP022804	NM_004849	ATG5
QG065-03	C01	HQP022747	NM_175862	CD86
QG065-03	C02	HQP022699	NM_006139	CD28
QG065-03	C03	HQP022491	NM_001616	ACVR2A
QG065-03	C04	HQP022361	NM_004747	DLG5
QG065-03	C05	HQP022168	NM_004707	ATG12
QG065-03	C06	HQP021752	NM_003939	BTRC
QG065-03	C07	HQP021683	NM_021090	MTMR3
QG065-03	C08	HQP021630	NM_003889	NR1I2
QG065-03	C09	HQP021610	NM_003882	WISP1
QG065-03	C10	HQP021603	NM_003878	GGH
QG065-03	C11	HQP021557	NM_003844	TNFRSF10A
QG065-03	C12	HQP021553	NM_003842	TNFRSF10B
QG065-03	D01	HQP021492	NM_003804	RIPK1
QG065-03	D02	HQP021459	NM_003786	ABCC3
QG065-03	D03	HQP021313	NM_001753	CAV1
QG065-03	D04	HQP021276	NM_003672	CDC14A
QG065-03	D05	HQP021250	NM_033389	SSH2
QG065-03	D06	HQP020982	NM_001039970	PTPN5
QG065-03	D07	HQP020596	NM_001042544	LTBP4
QG065-03	D08	HQP020481	NM_033338	CASP7
QG065-03	D09	HQP020427	NM_001226	CASP6
QG065-03	D10	HQP020297	NM_004346	CASP3
QG065-03	D11	HQP019829	NM_030782	CLPTM1L
QG065-03	D12	HQP019462	NM_024945	RMI1
QG065-03	E01	HQP019284	NM_024757	EHMT1
QG065-03	E02	HQP019172	NM_024642	GALNT12
QG065-03	E03	HQP019051	NM_007088	CALB2
QG065-03	E04	HQP018802	NM_001008540	CXCR4
QG065-03	E05	HQP018755	NM_001007257	PRDM2
QG065-03	E06	HQP018659	NM_001215	CA6
QG065-03	E07	HQP018545	NM_000553	WRN
QG065-03	E08	HQP018540	NM_004185	WNT2B
QG065-03	E09	HQP018485	NM_000551	VHL
QG065-03	E10	HQP018418	NM_001074	UGT2B7
QG065-03	E11	HQP018337	NM_182729	TXNRD1
QG065-03	E12	HQP018288	NM_003318	TTK
QG065-03	F01	HQP018265	NM_000548	TSC2
QG065-03	F02	HQP018263	NM_000368	TSC1
QG065-03	F03	HQP018176	NM_005657	TP53BP1
QG065-03	F04	HQP018104	NM_005077	TLE1
QG065-03	F05	HQP017958	NM_003202	TCF7
QG065-03	F06	HQP017954	NM_000545	TCF1
QG065-03	F07	HQP017535	NM_001204	BMPR2
QG065-03	F08	HQP017496	NM_003060	SLC22A5
QG065-03	F09	HQP017495	NM_003059	SLC22A4
QG065-03	F10	HQP016830	NM_003010	MAP2K4



Product Data Sheet

QG065-03	F11	HQP016746	NM_003005	SELP
QG065-03	F12	HQP016652	NM_002996	CX3CL1
QG065-03	G01	HQP016634	NM_001713	BHMT
QG065-03	G02	HQP016621	NM_002982	CCL2
QG065-03	G03	HQP016620	NM_002981	CCL1
QG065-03	G04	HQP016471	NM_003161	RPS6KB1
QG065-03	G05	HQP016468	NM_001006932	RPS6KA2
QG065-03	G06	HQP016466	NM_001006665	RPS6KA1
QG065-03	G07	HQP016303	NM_021133	RNASEL
QG065-03	G08	HQP016224	NM_002913	RFC1
QG065-03	G09	HQP016168	NM_021629	GNB4
QG065-03	G10	HQP016131	NM_000321	RB1
QG065-03	G11	HQP016087	NM_134424	RAD52
QG065-03	G12	HQP015946	NM_000465	BARD1
QG065-03	H01	HGDC		
QG065-03	H02	HGDC		
QG065-03	H03	HQP006940	NM_002046	GAPDH
QG065-03	H04	HQP016381	NM_001101	ACTB
QG065-03	H05	HQP015171	NM_004048	B2M
QG065-03	H06	HQP006171	NM_012423	RPL13A
QG065-03	H07	HQP009026	NM_000194	HPRT1
QG065-03	H08	HQP054253	NR_003286	RN18S1
QG065-03	H09	RT		
QG065-03	H10	RT		
QG065-03	H11	PCR		
QG065-03	H12	PCR		
QG065-04	A01	HQP015939	NM_002850	PTPRS
QG065-04	A02	HQP015563	NM_000958	PTGER4
QG065-04	A03	HQP015544	NM_000956	PTGER2
QG065-04	A04	HQP015530	NM_000264	PTCH1
QG065-04	A05	HQP015311	NM_002800	PSMB9
QG065-04	A06	HQP015190	NM_020165	RAD18
QG065-04	A07	HQP015087	NM_019845	RPRM
QG065-04	A08	HQP015032	NM_000311	PRNP
QG065-04	A09	HQP014691	NM_001008744	TDP1
QG065-04	A10	HQP014530	NM_006251	PRKAA1
QG065-04	A11	HQP014500	NM_017633	FAM46A
QG065-04	A12	HQP014362	NM_018407	LAPTM4B
QG065-04	B01	HQP014052	NM_018036	C14orf103
QG065-04	B02	HQP014024	NM_002711	PPP1R3A
QG065-04	B03	HQP013734	NM_017672	TRPM7
QG065-04	B04	HQP013596	NM_001184	ATR
QG065-04	B05	HQP013473	NM_000446	PON1
QG065-04	B06	HQP013467	NM_000939	POMC
QG065-04	B07	HQP013463	NM_018984	SSH1
QG065-04	B08	HQP013155	NM_181504	PIK3R1
QG065-04	B09	HQP013122	NM_000295	SERPINA1
QG065-04	B10	HQP013026	NM_002615	SERPINF1

## Product Data Sheet

QG065-04	B11	HQP012993	NM_016327	UPB1
QG065-04	B12	HQP012972	NM_016231	NLK
QG065-04	C01	HQP012723	NM_016203	PRKAG2
QG065-04	C02	HQP012675	NM_015891	CDC40
QG065-04	C03	HQP012420	NM_002592	PCNA
QG065-04	C04	HQP012393	NM_016009	SH3GLB1
QG065-04	C05	HQP012386	NM_015999	ADIPOR1
QG065-04	C06	HQP012161	NM_020525	IL22
QG065-04	C07	HQP012131	NM_015713	RRM2B
QG065-04	C08	HQP012121	NM_002569	FURIN
QG065-04	C09	HQP011932	NM_002526	NT5E
QG065-04	C10	HQP011920	NM_005173	ATP2A3
QG065-04	C11	HQP011909	NM_001681	ATP2A2
QG065-04	C12	HQP011853	NM_000269	NME1
QG065-04	D01	HQP011810	NM_020529	NFKBIA
QG065-04	D02	HQP011727	NM_004044	ATIC
QG065-04	D03	HQP011549	NM_000252	MTM1
QG065-04	D04	HQP011309	NM_000250	MPO
QG065-04	D05	HQP011266	NM_002426	MMP12
QG065-04	D06	HQP011211	NM_000246	CIITA
QG065-04	D07	HQP011208	NM_020300	MGST1
QG065-04	D08	HQP011047	NM_005911	MAT2A
QG065-04	D09	HQP011006	NM_005908	MANBA
QG065-04	D10	HQP010960	NM_005902	SMAD3
QG065-04	D11	HQP010959	NM_001003652	SMAD2
QG065-04	D12	HQP010924	NM_001668	ARNT
QG065-04	E01	HQP010912	NM_206943	LTBP1
QG065-04	E02	HQP010626	NM_000236	LIPC
QG065-04	E03	HQP010613	NM_002312	LIG4
QG065-04	E04	HQP010612	NM_013975	LIG3
QG065-04	E05	HQP010609	NM_000234	LIG1
QG065-04	E06	HQP010070	NM_002253	KDR
QG065-04	E07	HQP009718	NM_001562	IL18
QG065-04	E08	HQP009693	NM_002187	IL12B
QG065-04	E09	HQP009692	NM_000882	IL12A
QG065-04	E10	HQP009639	NM_001556	IKBKB
QG065-04	E11	HQP009539	NM_000596	IGFBP1
QG065-04	E12	HQP009334	NM_198498	C11orf53
QG065-04	F01	HQP009101	NM_001168	BIRC5
QG065-04	F02	HQP009065	NM_000197	HSD17B3
QG065-04	F03	HQP008891	NM_000859	HMGCR
QG065-04	F04	HQP008482	NM_000848	GSTM2
QG065-04	F05	HQP008479	NM_001512	GSTA4
QG065-04	F06	HQP008476	NM_000846	GSTA2
QG065-04	F07	HQP008434	NM_013264	DDX25
QG065-04	F08	HQP008402	NM_001020825	NR3C1
QG065-04	F09	HQP008282	NM_002084	GPX3
QG065-04	F10	HQP008281	NM_002083	GPX2

Product Data Sheet

QG065-04	F11	HQP008198	NM_173681	ATG9B
QG065-04	F12	HQP007755	NM_000516	GNAS
QG065-04	G01	HQP007703	NM_000168	GLI3
QG065-04	G02	HQP007526	NM_001018057	DKK3
QG065-04	G03	HQP007507	NM_000036	AMPD1
QG065-04	G04	HQP007213	NM_001634	AMD1
QG065-04	G05	HQP007110	NM_000819	GART
QG065-04	G06	HQP007058	NM_015617	PYGO1
QG065-04	G07	HQP006736	NM_012342	BAMBI
QG065-04	G08	HQP006715	NM_012415	RAD54B
QG065-04	G09	HQP006261	NM_014324	AMACR
QG065-04	G10	HQP006228	NM_001018078	FPGS
QG065-04	G11	HQP006144	NM_000802	FOLR1
QG065-04	G12	HQP005879	NM_002019	FLT1
QG065-04	H01	HGDC		
QG065-04	H02	HGDC		
QG065-04	H03	HQP006940	NM_002046	GAPDH
QG065-04	H04	HQP016381	NM_001101	ACTB
QG065-04	H05	HQP015171	NM_004048	B2M
QG065-04	H06	HQP006171	NM_012423	RPL13A
QG065-04	H07	HQP009026	NM_000194	HPRT1
QG065-04	H08	HQP054253	NR_003286	RN18S1
QG065-04	H09	RT		
QG065-04	H10	RT		
QG065-04	H11	PCR		
QG065-04	H12	PCR		

### Limited Use License

Following terms and conditions apply to use of ExProfile™ Colorectal 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)

GeneCopoeia Products are for Research Use Only

Trademarks: GeneCopoeia™, ExProfile™, All-in-One™, (GeneCopoeia Inc.); RNAzol® 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.

QG065-160202