

ExProfile™ Human General Toxicology Related Gene qPCR Array

For focused group profiling of human general toxicology genes expression

Cat. No. QG020-A (4 x 96-well plate, Format A)

Cat. No. QG020-B (4 x 96-well plate, Format B)

Cat. No. QG020-C (4 x 96-well plate, Format C)

Cat. No. QG020-D (4 x 96-well plate, Format D)

Cat. No. QG020-E (4 x 96-well plate, Format E)

Plates available individually or as a set of 6. Each set contains 336 unique gene primer pairs deposited in one 96-well plate.

Introduction

The ExProfile human general toxicology related gene qPCR array profiles the expression of 336 human genes related to the metabolic responses against drugs and other toxins. These genes are carefully chosen for their close pathway correlation based on a thorough literature search of peer-reviewed publications, mainly including genes involved in DNA damage and repair processes, apoptosis, cell cycle, electron transport, growth and proliferation, metabolite transporters, as well as stress response genes. This array allows researchers to study the pathway-related genes to gain understanding of their roles in the metabolic responses against drugs and other toxins.

- QG020 plate 01: 84 unique gene PCR primer pairs
- QG020 plate 02: 84 unique gene PCR primer pairs
- QG020 plate 03: 84 unique gene PCR primer pairs
- QG020 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
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	TNFRSF1A	TNFRSF11B	TNFRSF11A	TNFRSF10A	TNFAIP3	TNF	TH	TGFBR1	TGFB2	TGFB1	TGFA	TFDP1
B	TDP1	TDG	TCP1	TAP1	SULT1E1	SULT1A1	STAT5A	SRD5A2	SOD2	SOD1	SLC01C1	SLC2A1
C	SLC27A4	SLC27A1	SLC10A2	SLC10A1	SIRT1	SHH	SERPINE1	SDHB	SCP2	SCD	SAA1	RXRβ
D	RXRA	RELB	RB1	RARA	RAG1	RAF1	RAD52	RAD51	RAD50	RAD23A	RAD21	RAD18
E	RAD1	RAC1	PTGS2	PTGS1	PTEN	PTCH1	PRDX1	PPARGC1A	PPARG	PPARD	POR	PON3
F	POLL	POLK	POLI	POLB	PMS2	PLTP	PLA2G2A	PKLR	PGR	PERP	PCNA	ORM1
G	NRAS	NR1H2	NQO1	NOS3	NOS2A	NNMT	NFKB1	NF1	NDUFB1	NDUFB4	NDUFB5	NAT2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG020 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	NAT1	MYC	MVP	MTTP	MT3	MT2A	MSH6	MSH2	MPO	MPG	MMP9	MMP2
B	MLL	MLH1	MIF	MGST1	MET	MDM2	MCL1	MBL2	MBD4	MB	MAPK8	MAPK1
C	MAOB	MAOA	LTBR	LTA	LIG4	LIG1	LDLR	LBP	JUNB	JUN	IVD	ITGB2
D	ITGAL	INS	IL6	IL1B	IL1A	IL18	IGFBP6	IGFBP3	IGFBP2	IGFBP1	IGF2R	IGF1R
E	IGF1	HYOU1	HUS1	HSPD1	HSPB2	HSPB1	HSPA9	HSPA5	HSPA2	HSPA1L	HSPA1B	HSPA1A
F	HSP90AA1	HSF1	HPX	HPRT1	HMOX1	HIF1A	HGF	HBEGF	HAO1	GTF2H1	GSTA4	GSR
G	GSK3B	GPX5	GPX2	GPX1	GPD2	GAP43	GADD45A	FXYD2	FOXO3A	FOXO1	FOS	FMO4
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG020 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
A	FMO3	FMO1	FGFR2	FGF2	FGF1	FEN1	FDXR	FASLG	FAS	FANCC	FANCA	FADD
B	FABP7	FABP1	ESR2	ESR1	ERCC3	ERCC2	ERCC1	ERBB3	ERBB2	EPX	EPHX1	ELK1
C	EGR1	EGFR	EGF	EDNRB	EDN1	E2F4	E2F1	DPYD	DNAJB1	DNAJA1	DIABLO	DHFR
D	DDIT3	DDC	DDB2	DCLRE1C	DAP	CYP8B1	CYP7B1	CYP7A1	CYP51A1	CYP4B1	CYP4A11	CYP46A1
E	CYP3A5	CYP3A4	CYP2S1	CYP2F1	CYP2E1	CYP2D6	CYP2C9	CYP2B6	CYP2A13	CYP27B1	CYP27A1	CYP26A1
F	CYP24A1	CYP1B1	CYP1A2	CYP1A1	CYP19A1	CYP17A1	CYP11A1	CYB5R3	CXCL10	CTPS	CTNNB1	CSK
G	CSF2	CPT1A	CP	COMT	COL1A1	CLU	CHEK2	CHAT	CES1	CDKN2A	CDKN1B	CDKN1A
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure3. Illustration of QG020 plate 03

	1	2	3	4	5	6	7	8	9	10	11	12
A	CDK4	CDK2	CDC6	CDC37	CDC25C	CDC25B	CDC25A	CDC20	CD40	CD36	CCNG2	CCNE1
B	CCND2	CCND1	CCNC	CCNB1	CCNA2	CCL4	CASP9	CASP8	CASP3	BRCA2	BRCA1	BOK
C	BNIP3L	BNIP3	BLM	BIRC5	BID	BCR	BCL6	BCL2L1	BCL2	BCHE	BAX	BAK1
D	BAG1	BAD	ATP2A1	ATM	ARNT	AR	APP	APOE	APOC2	APOA1	APEX1	APC
E	APAF1	AOC3	ANXA5	ALOX5	ALDH1A1	AKT1	AHR	ADFP	ACAT1	ABP1	ABL1	ABCG8
F	ABCG5	ABCG2	ABCD3	ABCD2	ABCD1	ABCC6	ABCC5	ABCC3	ABCC2	ABCC1	ABCB4	ABCB11
G	ABCB1	ABCA1	GAB1	GADD45B	GCDH	NDUFB7	NTHL1	PKMYT1	PMAIP1	APRT	ABCG1	ACADSB
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure4. Illustration of QG020 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
QG020-01	A01	HQP018148	NM_001065	TNFRSF1A
QG020-01	A02	HQP012049	NM_002546	TNFRSF11B
QG020-01	A03	HQP021550	NM_003839	TNFRSF11A
QG020-01	A04	HQP021557	NM_003844	TNFRSF10A
QG020-01	A05	HQP018145	NM_006290	TNFAIP3
QG020-01	A06	HQP018141	NM_000594	TNF
QG020-01	A07	HQP018064	NM_000360	TH
QG020-01	A08	HQP018051	NM_004612	TGFBR1
QG020-01	A09	HQP018047	NM_003238	TGFB2
QG020-01	A10	HQP018044	NM_000660	TGFB1

QG020-01	A11	HQP018043	NM_003236	TGFA
QG020-01	A12	HQP018032	NM_007111	TFDP1
QG020-01	B01	HQP014691	NM_001008744	TDP1
QG020-01	B02	HQP017996	NM_003211	TDG
QG020-01	B03	HQP017984	NM_001008897	TCP1
QG020-01	B04	HQP017899	NM_000593	TAP1
QG020-01	B05	HQP017784	NM_005420	SULT1E1
QG020-01	B06	HQP017810	NM_001055	SULT1A1
QG020-01	B07	HQP017771	NM_003152	STAT5A
QG020-01	B08	HQP017698	NM_000348	SRD5A2
QG020-01	B09	HQP017616	NM_000636	SOD2
QG020-01	B10	HQP017615	NM_000454	SOD1
QG020-01	B11	HQP013340	NM_017435	SLCO1C1
QG020-01	B12	HQP017350	NM_006516	SLC2A1
QG020-01	C01	HQP001134	NM_005094	SLC27A4
QG020-01	C02	HQP010011	NM_198580	SLC27A1
QG020-01	C03	HQP017462	NM_000452	SLC10A2
QG020-01	C04	HQP017461	NM_003049	SLC10A1
QG020-01	C05	HQP006080	NM_012238	SIRT1
QG020-01	C06	HQP017098	NM_000193	SHH
QG020-01	C07	HQP012154	NM_000602	SERPINE1
QG020-01	C08	HQP016689	NM_003000	SDHB
QG020-01	C09	HQP016613	NM_001007098	SCP2
QG020-01	C10	HQP016583	NM_005063	SCD
QG020-01	C11	HQP016555	NM_000331	SAA1
QG020-01	C12	HQP016527	NM_021976	RXRβ
QG020-01	D01	HQP016526	NM_002957	RXRA
QG020-01	D02	HQP016214	NM_006509	RELB
QG020-01	D03	HQP016131	NM_000321	RB1
QG020-01	D04	HQP016114	NM_000964	RARA
QG020-01	D05	HQP016089	NM_000448	RAG1
QG020-01	D06	HQP016088	NM_002880	RAF1
QG020-01	D07	HQP016087	NM_134424	RAD52
QG020-01	D08	HQP016077	NM_002875	RAD51
QG020-01	D09	HQP000145	NM_005732	RAD50
QG020-01	D10	HQP016075	NM_005053	RAD23A
QG020-01	D11	HQP016074	NM_006265	RAD21
QG020-01	D12	HQP015190	NM_020165	RAD18
QG020-01	E01	HQP015947	NM_002853	RAD1
QG020-01	E02	HQP016063	NM_006908	RAC1
QG020-01	E03	HQP015598	NM_000963	PTGS2
QG020-01	E04	HQP015596	NM_000962	PTGS1
QG020-01	E05	HQP015535	NM_000314	PTEN
QG020-01	E06	HQP015530	NM_000264	PTCH1
QG020-01	E07	HQP012152	NM_002574	PRDX1
QG020-01	E08	HQP001016	NM_013261	PPARGC1A

QG020-01	E09	HQP013633	NM_005037	PPARG
QG020-01	E10	HQP013627	NM_006238	PPARD
QG020-01	E11	HQP013504	NM_000941	POR
QG020-01	E12	HQP013493	NM_000940	PON3
QG020-01	F01	HQP007689	NM_013274	POLL
QG020-01	F02	HQP012725	NM_016218	POLK
QG020-01	F03	HQP001399	NM_007195	POLI
QG020-01	F04	HQP013422	NM_002690	POLB
QG020-01	F05	HQP013352	NM_000535	PMS2
QG020-01	F06	HQP013280	NM_006227	PLTP
QG020-01	F07	HQP013194	NM_000300	PLA2G2A
QG020-01	F08	HQP013181	NM_000298	PKLR
QG020-01	F09	HQP013099	NM_000926	PGR
QG020-01	F10	HQP016753	NM_022121	PERP
QG020-01	F11	HQP012420	NM_002592	PCNA
QG020-01	F12	HQP012078	NM_000607	ORM1
QG020-01	G01	HQP011914	NM_002524	NRAS
QG020-01	G02	HQP021630	NM_003889	NR1I2
QG020-01	G03	HQP004317	NM_000903	NQO1
QG020-01	G04	HQP011868	NM_000603	NOS3
QG020-01	G05	HQP011866	NM_000625	NOS2A
QG020-01	G06	HQP011860	NM_006169	NNMT
QG020-01	G07	HQP011807	NM_003998	NFKB1
QG020-01	G08	HQP011774	NM_000267	NF1
QG020-01	G09	HQP011730	NM_007103	NDUFV1
QG020-01	G10	HQP011731	NM_002495	NDUFS4
QG020-01	G11	HQP011726	NM_005006	NDUFS1
QG020-01	G12	HQP001136	NM_000015	NAT2
QG020-01	H01	HGDC		
QG020-01	H02	HGDC		
QG020-01	H03	HQP006940	NM_002046	GAPDH
QG020-01	H04	HQP016381	NM_001101	ACTB
QG020-01	H05	HQP015171	NM_004048	B2M
QG020-01	H06	HQP006171	NM_012423	RPL13A
QG020-01	H07	HQP009026	NM_000194	HPRT1
QG020-01	H08	HQP054253	NR_003286	RN18S1
QG020-01	H09	RT		
QG020-01	H10	RT		
QG020-01	H11	PCR		
QG020-01	H12	PCR		
QG020-02	A01	HQP023467	NM_000662	NAT1
QG020-02	A02	HQP011597	NM_002467	MYC
QG020-02	A03	HQP023412	NM_005115	MVP
QG020-02	A04	HQP011553	NM_000253	MTTP
QG020-02	A05	HQP011539	NM_005954	MT3
QG020-02	A06	HQP011538	NM_005953	MT2A

QG020-02	A07	HQP008493	NM_000179	MSH6
QG020-02	A08	HQP011491	NM_000251	MSH2
QG020-02	A09	HQP011309	NM_000250	MPO
QG020-02	A10	HQP011304	NM_001015052	MPG
QG020-02	A11	HQP011263	NM_004994	MMP9
QG020-02	A12	HQP011256	NM_004530	MMP2
QG020-02	B01	HQP011241	NM_005933	MLL
QG020-02	B02	HQP011235	NM_000249	MLH1
QG020-02	B03	HQP011219	NM_002415	MIF
QG020-02	B04	HQP011208	NM_020300	MGST1
QG020-02	B05	HQP011181	NM_000245	MET
QG020-02	B06	HQP011135	NM_002392	MDM2
QG020-02	B07	HQP011104	NM_021960	MCL1
QG020-02	B08	HQP011077	NM_000242	MBL2
QG020-02	B09	HQP021736	NM_003925	MBD4
QG020-02	B10	HQP011071	NM_005368	MB
QG020-02	B11	HQP014886	NM_002750	MAPK8
QG020-02	B12	HQP014848	NM_002745	MAPK1
QG020-02	C01	HQP011008	NM_000898	MAOB
QG020-02	C02	HQP011007	NM_000240	MAOA
QG020-02	C03	HQP010915	NM_002342	LTBR
QG020-02	C04	HQP010907	NM_000595	LTA
QG020-02	C05	HQP010613	NM_002312	LIG4
QG020-02	C06	HQP010609	NM_000234	LIG1
QG020-02	C07	HQP010577	NM_000527	LDLR
QG020-02	C08	HQP010560	NM_004139	LBP
QG020-02	C09	HQP009854	NM_002229	JUNB
QG020-02	C10	HQP009853	NM_002228	JUN
QG020-02	C11	HQP009845	NM_002225	IVD
QG020-02	C12	HQP009815	NM_000211	ITGB2
QG020-02	D01	HQP009806	NM_002209	ITGAL
QG020-02	D02	HQP009749	NM_000207	INS
QG020-02	D03	HQP009670	NM_000600	IL6
QG020-02	D04	HQP009641	NM_000576	IL1B
QG020-02	D05	HQP009640	NM_000575	IL1A
QG020-02	D06	HQP009718	NM_001562	IL18
QG020-02	D07	HQP009555	NM_002178	IGFBP6
QG020-02	D08	HQP009544	NM_000598	IGFBP3
QG020-02	D09	HQP009541	NM_000597	IGFBP2
QG020-02	D10	HQP009539	NM_000596	IGFBP1
QG020-02	D11	HQP009532	NM_000876	IGF2R
QG020-02	D12	HQP009523	NM_000875	IGF1R
QG020-02	E01	HQP009518	NM_000618	IGF1
QG020-02	E02	HQP000633	NM_006389	HYOU1
QG020-02	E03	HQP009138	NM_004507	HUS1
QG020-02	E04	HQP009098	NM_002156	HSPD1

QG020-02	E05	HQP009090	NM_001541	HSPB2
QG020-02	E06	HQP009089	NM_001540	HSPB1
QG020-02	E07	HQP009088	NM_004134	HSPA9
QG020-02	E08	HQP009083	NM_005347	HSPA5
QG020-02	E09	HQP009080	NM_021979	HSPA2
QG020-02	E10	HQP009079	NM_005527	HSPA1L
QG020-02	E11	HQP009078	NM_005346	HSPA1B
QG020-02	E12	HQP009077	NM_005345	HSPA1A
QG020-02	F01	HQP009092	NM_001017963	HSP90AA1
QG020-02	F02	HQP009068	NM_005526	HSF1
QG020-02	F03	HQP009035	NM_000613	HPX
QG020-02	F04	HQP009026	NM_000194	HPRT1
QG020-02	F05	HQP008898	NM_002133	HMOX1
QG020-02	F06	HQP008831	NM_001530	HIF1A
QG020-02	F07	HQP008800	NM_000601	HGF
QG020-02	F08	HQP004493	NM_001945	HBEGF
QG020-02	F09	HQP013447	NM_017545	HAO1
QG020-02	F10	HQP008502	NM_005316	GTF2H1
QG020-02	F11	HQP008479	NM_001512	GSTA4
QG020-02	F12	HQP008473	NM_000637	GSR
QG020-02	G01	HQP008469	NM_002093	GSK3B
QG020-02	G02	HQP008288	NM_001509	GPX5
QG020-02	G03	HQP008281	NM_002083	GPX2
QG020-02	G04	HQP008279	NM_000581	GPX1
QG020-02	G05	HQP007813	NM_000408	GPD2
QG020-02	G06	HQP006930	NM_002045	GAP43
QG020-02	G07	HQP004125	NM_001924	GADD45A
QG020-02	G08	HQP011894	NM_001680	FXYD2
QG020-02	G09	HQP005759	NM_001455	FOXO3A
QG020-02	G10	HQP005712	NM_021953	FOXM1
QG020-02	G11	HQP006188	NM_005252	FOS
QG020-02	G12	HQP005968	NM_002022	FMO4
QG020-02	H01	HGDC		
QG020-02	H02	HGDC		
QG020-02	H03	HQP006940	NM_002046	GAPDH
QG020-02	H04	HQP016381	NM_001101	ACTB
QG020-02	H05	HQP015171	NM_004048	B2M
QG020-02	H06	HQP006171	NM_012423	RPL13A
QG020-02	H07	HQP009026	NM_000194	HPRT1
QG020-02	H08	HQP054253	NR_003286	RN18S1
QG020-02	H09	RT		
QG020-02	H10	RT		
QG020-02	H11	PCR		
QG020-02	H12	PCR		
QG020-03	A01	HQP005959	NM_001002294	FMO3
QG020-03	A02	HQP005939	NM_002021	FMO1

QG020-03	A03	HQP005437	NM_000141	FGFR2
QG020-03	A04	HQP005403	NM_002006	FGF2
QG020-03	A05	HQP005400	NM_000800	FGF1
QG020-03	A06	HQP005391	NM_004111	FEN1
QG020-03	A07	HQP005387	NM_004110	FDXR
QG020-03	A08	HQP009671	NM_000639	FASLG
QG020-03	A09	HQP009651	NM_000043	FAS
QG020-03	A10	HQP005083	NM_000136	FANCC
QG020-03	A11	HQP005081	NM_000135	FANCA
QG020-03	A12	HQP021526	NM_003824	FADD
QG020-03	B01	HQP005080	NM_001446	FABP7
QG020-03	B02	HQP005073	NM_001443	FABP1
QG020-03	B03	HQP005001	NM_001040275	ESR2
QG020-03	B04	HQP004998	NM_000125	ESR1
QG020-03	B05	HQP004983	NM_000122	ERCC3
QG020-03	B06	HQP004976	NM_000400	ERCC2
QG020-03	B07	HQP004974	NM_001983	ERCC1
QG020-03	B08	HQP004970	NM_001005915	ERBB3
QG020-03	B09	HQP004968	NM_001005862	ERBB2
QG020-03	B10	HQP020089	NM_000502	EPX
QG020-03	B11	HQP004948	NM_000120	EPHX1
QG020-03	B12	HQP004749	NM_005229	ELK1
QG020-03	C01	HQP004612	NM_001964	EGR1
QG020-03	C02	HQP004605	NM_005228	EGFR
QG020-03	C03	HQP004599	NM_001963	EGF
QG020-03	C04	HQP004564	NM_000115	EDNRB
QG020-03	C05	HQP004557	NM_001955	EDN1
QG020-03	C06	HQP004528	NM_001950	E2F4
QG020-03	C07	HQP004524	NM_005225	E2F1
QG020-03	C08	HQP004439	NM_000110	DPYD
QG020-03	C09	HQP009103	NM_006145	DNAJB1
QG020-03	C10	HQP009076	NM_001539	DNAJA1
QG020-03	C11	HQP015116	NM_019887	DIABLO
QG020-03	C12	HQP004309	NM_000791	DHFR
QG020-03	D01	HQP004127	NM_004083	DDIT3
QG020-03	D02	HQP004115	NM_000790	DDC
QG020-03	D03	HQP004114	NM_000107	DDB2
QG020-03	D04	HQP016982	NM_001033855	DCLRE1C
QG020-03	D05	HQP003975	NM_004394	DAP
QG020-03	D06	HQP003865	NM_004391	CYP8B1
QG020-03	D07	HQP022723	NM_004820	CYP7B1
QG020-03	D08	HQP003859	NM_000780	CYP7A1
QG020-03	D09	HQP003924	NM_000786	CYP51A1
QG020-03	D10	HQP003855	NM_000779	CYP4B1
QG020-03	D11	HQP003847	NM_000778	CYP4A11
QG020-03	D12	HQP000977	NM_006668	CYP46A1

QG020-03	E01	HQP003841	NM_000777	CYP3A5
QG020-03	E02	HQP003836	NM_017460	CYP3A4
QG020-03	E03	HQP008529	NM_030622	CYP2S1
QG020-03	E04	HQP003818	NM_000774	CYP2F1
QG020-03	E05	HQP003817	NM_000773	CYP2E1
QG020-03	E06	HQP003814	NM_000106	CYP2D6
QG020-03	E07	HQP003811	NM_000771	CYP2C9
QG020-03	E08	HQP003808	NM_000767	CYP2B6
QG020-03	E09	HQP003804	NM_000766	CYP2A13
QG020-03	E10	HQP003923	NM_000785	CYP27B1
QG020-03	E11	HQP003922	NM_000784	CYP27A1
QG020-03	E12	HQP003919	NM_000783	CYP26A1
QG020-03	F01	HQP003917	NM_000782	CYP24A1
QG020-03	F02	HQP003775	NM_000104	CYP1B1
QG020-03	F03	HQP003774	NM_000761	CYP1A2
QG020-03	F04	HQP003772	NM_000499	CYP1A1
QG020-03	F05	HQP003904	NM_000103	CYP19A1
QG020-03	F06	HQP003888	NM_000102	CYP17A1
QG020-03	F07	HQP003871	NM_000781	CYP11A1
QG020-03	F08	HQP004315	NM_000398	CYB5R3
QG020-03	F09	HQP009746	NM_001565	CXCL10
QG020-03	F10	HQP003572	NM_001905	CTPS
QG020-03	F11	HQP003539	NM_001904	CTNNB1
QG020-03	F12	HQP003218	NM_004383	CSK
QG020-03	G01	HQP003159	NM_000758	CSF2
QG020-03	G02	HQP002867	NM_001031847	CPT1A
QG020-03	G03	HQP002820	NM_000096	CP
QG020-03	G04	HQP002671	NM_000754	COMT
QG020-03	G05	HQP002462	NM_000088	COL1A1
QG020-03	G06	HQP002037	NM_001831	CLU
QG020-03	G07	HQP001396	NM_001005735	CHEK2
QG020-03	G08	HQP001182	NM_020549	CHAT
QG020-03	G09	HQP000815	NM_001025194	CES1
QG020-03	G10	HQP000369	NM_000077	CDKN2A
QG020-03	G11	HQP000342	NM_004064	CDKN1B
QG020-03	G12	HQP000331	NM_000389	CDKN1A
QG020-03	H01	HGDC		
QG020-03	H02	HGDC		
QG020-03	H03	HQP006940	NM_002046	GAPDH
QG020-03	H04	HQP016381	NM_001101	ACTB
QG020-03	H05	HQP015171	NM_004048	B2M
QG020-03	H06	HQP006171	NM_012423	RPL13A
QG020-03	H07	HQP009026	NM_000194	HPRT1
QG020-03	H08	HQP054253	NR_003286	RN18S1
QG020-03	H09	RT		
QG020-03	H10	RT		

QG020-03	H11	PCR		
QG020-03	H12	PCR		
QG020-04	A01	HQP000245	NM_000075	CDK4
QG020-04	A02	HQP000225	NM_001798	CDK2
QG020-04	A03	HQP023354	NM_001254	CDC6
QG020-04	A04	HQP001310	NM_007065	CDC37
QG020-04	A05	HQP023409	NM_001790	CDC25C
QG020-04	A06	HQP023399	NM_004358	CDC25B
QG020-04	A07	HQP023385	NM_001789	CDC25A
QG020-04	A08	HQP023365	NM_001255	CDC20
QG020-04	A09	HQP022955	NM_001250	CD40
QG020-04	A10	HQP022821	NM_000072	CD36
QG020-04	A11	HQP021882	NM_004354	CCNG2
QG020-04	A12	HQP021819	NM_001238	CCNE1
QG020-04	B01	HQP021754	NM_001759	CCND2
QG020-04	B02	HQP016204	NM_053056	CCND1
QG020-04	B03	HQP021734	NM_001013399	CCNC
QG020-04	B04	HQP021727	NM_031966	CCNB1
QG020-04	B05	HQP021701	NM_001237	CCNA2
QG020-04	B06	HQP016625	NM_002984	CCL4
QG020-04	B07	HQP020648	NM_001229	CASP9
QG020-04	B08	HQP018966	NM_001080124	CASP8
QG020-04	B09	HQP020297	NM_004346	CASP3
QG020-04	B10	HQP017753	NM_000059	BRCA2
QG020-04	B11	HQP017713	NM_007294	BRCA1
QG020-04	B12	HQP017642	NM_032515	BOK
QG020-04	C01	HQP017631	NM_004331	BNIP3L
QG020-04	C02	HQP017619	NM_004052	BNIP3
QG020-04	C03	HQP016844	NM_000057	BLM
QG020-04	C04	HQP009099	NM_001012270	BIRC5
QG020-04	C05	HQP016653	NM_001196	BID
QG020-04	C06	HQP016418	NM_004327	BCR
QG020-04	C07	HQP016328	NM_001706	BCL6
QG020-04	C08	HQP016237	NM_001191	BCL2L1
QG020-04	C09	HQP016211	NM_000633	BCL2
QG020-04	C10	HQP016108	NM_000055	BCHE
QG020-04	C11	HQP015964	NM_004324	BAX
QG020-04	C12	HQP015917	NM_001188	BAK1
QG020-04	D01	HQP015574	NM_004323	BAG1
QG020-04	D02	HQP015538	NM_004322	BAD
QG020-04	D03	HQP011898	NM_004320	ATP2A1
QG020-04	D04	HQP011736	NM_000051	ATM
QG020-04	D05	HQP010924	NM_001668	ARNT
QG020-04	D06	HQP009801	NM_000044	AR
QG020-04	D07	HQP009578	NM_000484	APP
QG020-04	D08	HQP009556	NM_000041	APOE

QG020-04	D09	HQP009447	NM_000483	APOC2
QG020-04	D10	HQP009125	NM_000039	APOA1
QG020-04	D11	HQP009061	NM_001641	APEX1
QG020-04	D12	HQP009024	NM_000038	APC
QG020-04	E01	HQP008934	NM_001160	APAF1
QG020-04	E02	HQP021378	NM_003734	AOC3
QG020-04	E03	HQP008829	NM_001154	ANXA5
QG020-04	E04	HQP006359	NM_000698	ALOX5
QG020-04	E05	HQP005075	NM_000689	ALDH1A1
QG020-04	E06	HQP004991	NM_001014431	AKT1
QG020-04	E07	HQP004658	NM_001621	AHR
QG020-04	E08	HQP002234	NM_001122	ADFP
QG020-04	E09	HQP010411	NM_000019	ACAT1
QG020-04	E10	HQP007422	NM_001091	ABP1
QG020-04	E11	HQP006954	NM_005157	ABL1
QG020-04	E12	HQP016876	NM_022437	ABCG8
QG020-04	F01	HQP016875	NM_022436	ABCG5
QG020-04	F02	HQP022745	NM_004827	ABCG2
QG020-04	F03	HQP015971	NM_002858	ABCD3
QG020-04	F04	HQP005426	NM_005164	ABCD2
QG020-04	F05	HQP005065	NM_000033	ABCD1
QG020-04	F06	HQP009816	NM_001079528	ABCC6
QG020-04	F07	HQP000074	NM_001023587	ABCC5
QG020-04	F08	HQP021459	NM_003786	ABCC3
QG020-04	F09	HQP002260	NM_000392	ABCC2
QG020-04	F10	HQP011322	NM_004996	ABCC1
QG020-04	F11	HQP013101	NM_000443	ABCB4
QG020-04	F12	HQP021388	NM_003742	ABCB11
QG020-04	G01	HQP013100	NM_000927	ABCB1
QG020-04	G02	HQP004727	NM_005502	ABCA1
QG020-04	G03	HQP006558	NM_002039	GAB1
QG020-04	G04	HQP011604	NM_015675	GADD45B
QG020-04	G05	HQP007231	NM_000159	GCDH
QG020-04	G06	HQP011720	NM_004146	NDUFB7
QG020-04	G07	HQP011937	NM_002528	NTHL1
QG020-04	G08	HQP022046	NM_004203	PKMYT1
QG020-04	G09	HQP013296	NM_021127	PMAIP1
QG020-04	G10	HQP009626	NM_000485	APRT
QG020-04	G11	HQP022987	NM_004915	ABCG1
QG020-04	G12	HQP009833	NM_001609	ACADSB
QG020-04	H01	HGDC		
QG020-04	H02	HGDC		
QG020-04	H03	HQP006940	NM_002046	GAPDH
QG020-04	H04	HQP016381	NM_001101	ACTB
QG020-04	H05	HQP015171	NM_004048	B2M
QG020-04	H06	HQP006171	NM_012423	RPL13A

QG020-04	H07	HQP009026	NM_000194	HPRT1
QG020-04	H08	HQP054253	NR_003286	RN18S1
QG020-04	H09	RT		
QG020-04	H10	RT		
QG020-04	H11	PCR		
QG020-04	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Human General Toxicology 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

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.

QG020-160202