

## ExProfile™ Human Mood Disorder Related Gene qPCR Array

For focused group profiling of human mood disorder disease genes expression

Cat. No. QG083-A (4 x 96-well plate, Format A)

Cat. No. QG083-B (4 x 96-well plate, Format B)

Cat. No. QG083-C (4 x 96-well plate, Format C)

Cat. No. QG083-D (4 x 96-well plate, Format D)

Cat. No. QG083-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 mood disorder related gene qPCR array profiles the expression of 336 human genes related to mood disorder. These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including key genes associated with bipolar affective disorder, schizophrenia and other mood disorders. This array allows researchers to study the related genes to gain understanding of their roles in the functioning and characterization of mood disorder.

- QG083 plate 01: 84 unique gene PCR primer pairs
- QG083 plate 02: 84 unique gene PCR primer pairs
- QG083 plate 03: 84 unique gene PCR primer pairs
- QG083 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

GeneCopoela 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™

### 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<sup>TM</sup> First-Strand cDNA Synthesis Kit

All-in-One<sup>TM</sup> 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	TRPV6	TRPM2	TRIP12	TRAF3IP1	TRAF1	TPH2	TPH1	TNF	TLE1	THRA	TGIF1	TGFB1
B	TFAP2B	TFAP2A	TF	TCF7	TCF4	SYN3	SYBL1	SULT1A1	STAR	ST8SIA2	SST	SPTBN4
C	SPTAN1	SPR	SOX10	SOD1	SNAP29	SMARCE1	SLC6A4	SLC6A2	SLC5A3	SLC18A1	SLC15A1	SFMBT2
D	SCG2	SAT1	S100A9	RUNX2	RTN4	RNF2	RIPK1	RGS4	RELN	RELA	RAP1GAP	RANBP9
E	RAB7A	PURA	PTGS2	PSMD8	PSMC6	PRKD3	PRKCZ	PRKCQ	PRKCI	PRKCG	PRKCD	PRKCB1
F	PRKCA	PRKACA	PREP	PPP2R5C	PPP2R2C	PPIE	PPFIA4	POLG	PLSCR4	PLP1	PLLP	PLCG1
G	PLA2G4C	PLA2G4A	PLA2G3	PLA2G2A	PLA2G1B	PITPNB	PIP5K2A	PIK4CB	PIK4CA	PIK3C3	PGK1	PFN2
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG083 plate 01

	1	2	3	4	5	6	7	8	9	10	11	12
A	PER3	PDE4D	PCNT	PCGF2	PAX5	PAWR	PARP1	PAFAH1B1	P2RX7	OTX2	NTRK1	NTF3
B	NRG1	NR3C1	NR1D1	NPY	NPAS2	NOS3	NOS1AP	NFKBIB	NFKB2	NFKB1	NDUFV2	NDUFV1
C	NDUF88	NDUF87	NDEL1	NAPG	MYC	MTRR	MTR	MTHFR	MOBP	MLC1	MERTK	ME2
D	MC5R	MBP	MAPK3	MAPK1	MAP1A	MAOA	MAL	LTA	LIN7A	LEF1	LARS2	KCNQ2
E	KCNN3	KCNK1	KATNB1	KATNAL1	KALRN	JUN	ITSN1	ITPR1	ISYNA1	INPP1	IMPA2	IMPA1
F	IL9R	IL6	IL2RB	IL1RN	IL1B	IL18	IGFBP2	IGF1	HTR3A	HTR2C	HTR2A	HTR1D
G	HTR1B	HSPA5	HSP90B1	HLA-DRA	HIP1R	HDAC1	GZMB	GSK3B	GRM4	GRM3	GRIN2D	GRIN2A
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure2. Illustration of QG083 plate 02

	1	2	3	4	5	6	7	8	9	10	11	12
<b>A</b>	GRIN1	GRIK4	GRIA1	GPX2	GPX1	GPC5	GOLGA2	GNB1L	GNAS	GFAP	GCH1	GALC
<b>B</b>	GAD1	GABRG2	GABRA5	GABRA1	G6PD	FZD3	FOSL1	FOS	FKBP5	FEZ1	FDFT1	FAT
<b>C</b>	FASLG	FAS	EZH2	ESR2	ESR1	ERBB3	EPB49	EP300	ENPP2	EMX2	EIF3S3	EIF2S2
<b>D</b>	EIF2S1	EIF2B1	EGFR	EED	EDG2	DUSP6	DTNBP1	DRD5	DRD4	DRD2	DRD1	DPYSL2
<b>E</b>	DOCK9	DNAJB1	DLG4	DISC1	DID01	DGKH	DBP	DAO	CXCL2	CX3CL1	CUTL2	CTBP1
<b>F</b>	CRTC1	CRHBP	CREBBP	CREB1	CPT2	COX6C	COX5A	COMT	CNP	CLOCK	CLDN11	CKM
<b>G</b>	CHUK	CHRNA7	CHRNA1	CHRM2	CEP290	CCND3	CCND2	CCND1	CCL3L1	CCL13	CCKAR	CBX2
<b>H</b>	<b>HGDC</b>	<b>HGDC</b>	<b>GAPDH</b>	<b>ACTB</b>	<b>B2M</b>	<b>RPL13A</b>	<b>HPRT1</b>	<b>RN18S1</b>	<b>RT</b>	<b>RT</b>	<b>PCR</b>	<b>PCR</b>

Figure3. Illustration of QG083 plate 03

	1	2	3	4	5	6	7	8	9	10	11	12
<b>A</b>	CASP8	CASP6	CASP2	CAMKK2	CALR	BPNT1	BMI1	BID	BCR	BCL2	BAX	BAK1
<b>B</b>	AVPR1B	ATXN1	ATP5J	ATP2A2	ATP1A3	ATM	ATF7IP	ATF5	ASPA	AR	APOL2	APOE
<b>C</b>	APAF1	ANKHD1	ALOX5	ALOX12	ALG9	AKT1	AGXT2L1	ADSS	ADRBK2	ADCY9	ACTN2	ACE
<b>D</b>	REG3A	SYNGR1	TBX1	TIMELESS	MOG	OLIG2	PENK	PRKACB	SYNJ1	TNNT1	ABCG1	ADARB1
<b>E</b>	ADCYAP1	ADRA1B	PLA2G4B	ATP5C1	ATP5G3	BDNF	CHMP1B	PLA2G4D	CLCF1	CRH	DIO2	DNMT1
<b>F</b>	DRD3	PMP22	GABRA3	GRIK3	HPR	HSPA12A	HTR1A	HTR4	IMMT	KCNQ5	KLK6	PRF1
<b>G</b>	MGAT3	MTHFD1	NCAM1	NDE1	NEDD4L	NTRK2	NUDT4	OLIG1	PAFAH1B3	PDLIM5	PER1	PER2
<b>H</b>	<b>HGDC</b>	<b>HGDC</b>	<b>GAPDH</b>	<b>ACTB</b>	<b>B2M</b>	<b>RPL13A</b>	<b>HPRT1</b>	<b>RN18S1</b>	<b>RT</b>	<b>RT</b>	<b>PCR</b>	<b>PCR</b>

Figure4. Illustration of QG083 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
QG083-01	A01	HQP014423	NM_018646	TRPV6
QG083-01	A02	HQP018256	NM_001001188	TRPM2
QG083-01	A03	HQP022528	NM_004238	TRIP12
QG083-01	A04	HQP007081	NM_015650	TRAF3IP1
QG083-01	A05	HQP018232	NM_005658	TRAF1
QG083-01	A06	HQP002121	NM_173353	TPH2
QG083-01	A07	HQP018195	NM_004179	TPH1

QG083-01	A08	HQP018141	NM_000594	TNF
QG083-01	A09	HQP018104	NM_005077	TLE1
QG083-01	A10	HQP018077	NM_003250	THRA
QG083-01	A11	HQP018056	NM_003244	TGIF1
QG083-01	A12	HQP018044	NM_000660	TGFB1
QG083-01	B01	HQP018026	NM_003221	TFAP2B
QG083-01	B02	HQP018023	NM_001032280	TFAP2A
QG083-01	B03	HQP018020	NM_001063	TF
QG083-01	B04	HQP017958	NM_003202	TCF7
QG083-01	B05	HQP017951	NM_003199	TCF4
QG083-01	B06	HQP020054	NM_003490	SYN3
QG083-01	B07	HQP017841	NM_005638	SYBL1
QG083-01	B08	HQP017810	NM_001055	SULT1A1
QG083-01	B09	HQP017762	NM_000349	STAR
QG083-01	B10	HQP019846	NM_006011	ST8SIA2
QG083-01	B11	HQP053963	NM_001048	SST
QG083-01	B12	HQP015850	NM_020971	SPTBN4
QG083-01	C01	HQP017688	NM_003127	SPTAN1
QG083-01	C02	HQP017676	NM_003124	SPR
QG083-01	C03	HQP017636	NM_006941	SOX10
QG083-01	C04	HQP017615	NM_000454	SOD1
QG083-01	C05	HQP022567	NM_004782	SNAP29
QG083-01	C06	HQP017556	NM_003079	SMARCE1
QG083-01	C07	HQP053962	NM_001045	SLC6A4
QG083-01	C08	HQP017397	NM_001043	SLC6A2
QG083-01	C09	HQP017388	NM_006933	SLC5A3
QG083-01	C10	HQP017478	NM_003053	SLC18A1
QG083-01	C11	HQP017471	NM_005073	SLC15A1
QG083-01	C12	HQP015830	NM_001029880	SFMBT2
QG083-01	D01	HQP018805	NM_003469	SCG2
QG083-01	D02	HQP016572	NM_002970	SAT1
QG083-01	D03	HQP016548	NM_002965	S100A9
QG083-01	D04	HQP016478	NM_001015051	RUNX2
QG083-01	D05	HQP015432	NM_007008	RTN4
QG083-01	D06	HQP016307	NM_007212	RNF2
QG083-01	D07	HQP021492	NM_003804	RIPK1
QG083-01	D08	HQP016259	NM_005613	RGS4
QG083-01	D09	HQP015094	NM_005045	RELN
QG083-01	D10	HQP016213	NM_021975	RELA
QG083-01	D11	HQP016107	NM_002885	RAP1GAP
QG083-01	D12	HQP000062	NM_005493	RANBP9

QG083-01	E01	HQP018819	NM_004637	RAB7A
QG083-01	E02	HQP015948	NM_005859	PURA
QG083-01	E03	HQP015598	NM_000963	PTGS2
QG083-01	E04	HQP015444	NM_002812	PSMD8
QG083-01	E05	HQP015369	NM_002806	PSMC6
QG083-01	E06	HQP006319	NM_005813	PRKD3
QG083-01	E07	HQP014829	NM_001033581	PRKCZ
QG083-01	E08	HQP014804	NM_006257	PRKCQ
QG083-01	E09	HQP014771	NM_002740	PRKCI
QG083-01	E10	HQP014751	NM_002739	PRKCG
QG083-01	E11	HQP014731	NM_006254	PRKCD
QG083-01	E12	HQP014718	NM_002738	PRKCB1
QG083-01	F01	HQP014706	NM_002737	PRKCA
QG083-01	F02	HQP014575	NM_002730	PRKACA
QG083-01	F03	HQP014431	NM_002726	PREP
QG083-01	F04	HQP014262	NM_002719	PPP2R5C
QG083-01	F05	HQP014202	NM_020416	PPP2R2C
QG083-01	F06	HQP000535	NM_006112	PPIE
QG083-01	F07	HQP021090	NM_015053	PPFIA4
QG083-01	F08	HQP013428	NM_002693	POLG
QG083-01	F09	HQP015379	NM_020353	PLSCR4
QG083-01	F10	HQP013272	NM_000533	PLP1
QG083-01	F11	HQP012381	NM_015993	PLLP
QG083-01	F12	HQP013238	NM_002660	PLCG1
QG083-01	G01	HQP021328	NM_003706	PLA2G4C
QG083-01	G02	HQP013195	NM_024420	PLA2G4A
QG083-01	G03	HQP012134	NM_015715	PLA2G3
QG083-01	G04	HQP013194	NM_000300	PLA2G2A
QG083-01	G05	HQP013193	NM_000928	PLA2G1B
QG083-01	G06	HQP006336	NM_012399	PITPNB
QG083-01	G07	HQP013171	NM_005028	PIP5K2A
QG083-01	G08	HQP013161	NM_002651	PIK4CB
QG083-01	G09	HQP013159	NM_002650	PIK4CA
QG083-01	G10	HQP013147	NM_002647	PIK3C3
QG083-01	G11	HQP013093	NM_000291	PGK1
QG083-01	G12	HQP013080	NM_002628	PFN2
QG083-01	H01	HGDC		
QG083-01	H02	HGDC		
QG083-01	H03	HQP006940	NM_002046	GAPDH
QG083-01	H04	HQP016381	NM_001101	ACTB
QG083-01	H05	HQP015171	NM_004048	B2M

QG083-01	H06	HQP006171	NM_012423	RPL13A
QG083-01	H07	HQP009026	NM_000194	HPRT1
QG083-01	H08	HQP054253	NR_003286	RN18S1
QG083-01	H09	RT		
QG083-01	H10	RT		
QG083-01	H11	PCR		
QG083-01	H12	PCR		
QG083-02	A01	HQP021640	NM_016831	PER3
QG083-02	A02	HQP012750	NM_006203	PDE4D
QG083-02	A03	HQP012474	NM_006031	PCNT
QG083-02	A04	HQP018681	NM_007144	PCGF2
QG083-02	A05	HQP012212	NM_016734	PAX5
QG083-02	A06	HQP012197	NM_002583	PAWR
QG083-02	A07	HQP003120	NM_001618	PARP1
QG083-02	A08	HQP012140	NM_000430	PAFAH1B1
QG083-02	A09	HQP012100	NM_002562	P2RX7
QG083-02	A10	HQP012085	NM_021728	OTX2
QG083-02	A11	HQP011938	NM_001007792	NTRK1
QG083-02	A12	HQP011933	NM_002527	NTF3
QG083-02	B01	HQP008818	NM_004495	NRG1
QG083-02	B02	HQP008401	NM_000176	NR3C1
QG083-02	B03	HQP022925	NM_021724	NR1D1
QG083-02	B04	HQP011874	NM_000905	NPY
QG083-02	B05	HQP011885	NM_002518	NPAS2
QG083-02	B06	HQP011868	NM_000603	NOS3
QG083-02	B07	HQP023119	NM_014697	NOS1AP
QG083-02	B08	HQP011811	NM_001001716	NFKBIB
QG083-02	B09	HQP011808	NM_001077493	NFKB2
QG083-02	B10	HQP011807	NM_003998	NFKB1
QG083-02	B11	HQP011735	NM_021074	NDUFV2
QG083-02	B12	HQP011730	NM_007103	NDUFV1
QG083-02	C01	HQP011734	NM_002496	NDUFS8
QG083-02	C02	HQP009875	NM_024407	NDUFS7
QG083-02	C03	HQP019903	NM_001025579	NDEL1
QG083-02	C04	HQP021529	NM_003826	NAPG
QG083-02	C05	HQP011597	NM_002467	MYC
QG083-02	C06	HQP011555	NM_002454	MTRR
QG083-02	C07	HQP011554	NM_000254	MTR
QG083-02	C08	HQP011547	NM_005957	MTHFR
QG083-02	C09	HQP011281	NM_182935	MOBP
QG083-02	C10	HQP005868	NM_015166	MLC1

QG083-02	C11	HQP000555	NM_006343	MERTK
QG083-02	C12	HQP011146	NM_002396	ME2
QG083-02	D01	HQP011098	NM_005913	MC5R
QG083-02	D02	HQP011085	NM_001025081	MBP
QG083-02	D03	HQP014854	NM_001040056	MAPK3
QG083-02	D04	HQP014848	NM_002745	MAPK1
QG083-02	D05	HQP011010	NM_002373	MAP1A
QG083-02	D06	HQP011007	NM_000240	MAOA
QG083-02	D07	HQP010995	NM_002371	MAL
QG083-02	D08	HQP010907	NM_000595	LTA
QG083-02	D09	HQP021587	NM_004664	LIN7A
QG083-02	D10	HQP012480	NM_016269	LEF1
QG083-02	D11	HQP006061	NM_015340	LARS2
QG083-02	D12	HQP010049	NM_004518	KCNQ2
QG083-02	E01	HQP010041	NM_002249	KCNN3
QG083-02	E02	HQP010024	NM_002245	KCNK1
QG083-02	E03	HQP000373	NM_005886	KATNB1
QG083-02	E04	HQP020437	NM_001014380	KATNAL1
QG083-02	E05	HQP021838	NM_001024660	KALRN
QG083-02	E06	HQP009853	NM_002228	JUN
QG083-02	E07	HQP017035	NM_001001132	ITSN1
QG083-02	E08	HQP009842	NM_002222	ITPR1
QG083-02	E09	HQP012769	NM_016368	ISYNA1
QG083-02	E10	HQP009747	NM_002194	INPP1
QG083-02	E11	HQP009727	NM_014214	IMPA2
QG083-02	E12	HQP009726	NM_005536	IMPA1
QG083-02	F01	HQP009683	NM_002186	IL9R
QG083-02	F02	HQP009670	NM_000600	IL6
QG083-02	F03	HQP009658	NM_000878	IL2RB
QG083-02	F04	HQP009645	NM_000577	IL1RN
QG083-02	F05	HQP009641	NM_000576	IL1B
QG083-02	F06	HQP009718	NM_001562	IL18
QG083-02	F07	HQP009541	NM_000597	IGFBP2
QG083-02	F08	HQP009518	NM_000618	IGF1
QG083-02	F09	HQP009123	NM_000869	HTR3A
QG083-02	F10	HQP009122	NM_000868	HTR2C
QG083-02	F11	HQP009120	NM_000621	HTR2A
QG083-02	F12	HQP009117	NM_000864	HTR1D
QG083-02	G01	HQP009116	NM_000863	HTR1B
QG083-02	G02	HQP009083	NM_005347	HSPA5
QG083-02	G03	HQP018231	NM_003299	HSP90B1

QG083-02	G04	HQP008866	NM_019111	HLA-DRA
QG083-02	G05	HQP021900	NM_003959	HIP1R
QG083-02	G06	HQP008745	NM_004964	HDAC1
QG083-02	G07	HQP008690	NM_004131	GZMB
QG083-02	G08	HQP008469	NM_002093	GSK3B
QG083-02	G09	HQP008449	NM_000841	GRM4
QG083-02	G10	HQP008448	NM_000840	GRM3
QG083-02	G11	HQP008380	NM_000836	GRIN2D
QG083-02	G12	HQP008371	NM_000833	GRIN2A
QG083-02	H01	HGDC		
QG083-02	H02	HGDC		
QG083-02	H03	HQP006940	NM_002046	GAPDH
QG083-02	H04	HQP016381	NM_001101	ACTB
QG083-02	H05	HQP015171	NM_004048	B2M
QG083-02	H06	HQP006171	NM_012423	RPL13A
QG083-02	H07	HQP009026	NM_000194	HPRT1
QG083-02	H08	HQP054253	NR_003286	RN18S1
QG083-02	H09	RT		
QG083-02	H10	RT		
QG083-02	H11	PCR		
QG083-02	H12	PCR		
QG083-03	A01	HQP008367	NM_000832	GRIN1
QG083-03	A02	HQP008363	NM_014619	GRIK4
QG083-03	A03	HQP008302	NM_000827	GRIA1
QG083-03	A04	HQP008281	NM_002083	GPX2
QG083-03	A05	HQP008279	NM_000581	GPX1
QG083-03	A06	HQP005436	NM_004466	GPC5
QG083-03	A07	HQP007788	NM_004486	GOLGA2
QG083-03	A08	HQP013589	NM_053004	GNB1L
QG083-03	A09	HQP007755	NM_000516	GNAS
QG083-03	A10	HQP007360	NM_002055	GFAP
QG083-03	A11	HQP007235	NM_000161	GCH1
QG083-03	A12	HQP006750	NM_000153	GALC
QG083-03	B01	HQP006683	NM_000817	GAD1
QG083-03	B02	HQP006651	NM_000816	GABRG2
QG083-03	B03	HQP006613	NM_000810	GABRA5
QG083-03	B04	HQP006599	NM_000806	GABRA1
QG083-03	B05	HQP006515	NM_000402	G6PD
QG083-03	B06	HQP019238	NM_017412	FZD3
QG083-03	B07	HQP019708	NM_005438	FOSL1
QG083-03	B08	HQP006188	NM_005252	FOS



QG083-03	B09	HQP005563	NM_004117	FKBP5
QG083-03	B10	HQP023016	NM_005103	FEZ1
QG083-03	B11	HQP005357	NM_004462	FDFT1
QG083-03	B12	HQP005141	NM_005245	FAT
QG083-03	C01	HQP009671	NM_000639	FASLG
QG083-03	C02	HQP009651	NM_000043	FAS
QG083-03	C03	HQP005050	NM_004456	EZH2
QG083-03	C04	HQP005001	NM_001040275	ESR2
QG083-03	C05	HQP004998	NM_000125	ESR1
QG083-03	C06	HQP004970	NM_001005915	ERBB3
QG083-03	C07	HQP004918	NM_001978	EPB49
QG083-03	C08	HQP004897	NM_001429	EP300
QG083-03	C09	HQP012960	NM_001040092	ENPP2
QG083-03	C10	HQP004842	NM_004098	EMX2
QG083-03	C11	HQP021417	NM_003756	EIF3S3
QG083-03	C12	HQP021680	NM_003908	EIF2S2
QG083-03	D01	HQP004643	NM_004094	EIF2S1
QG083-03	D02	HQP004650	NM_001414	EIF2B1
QG083-03	D03	HQP004605	NM_005228	EGFR
QG083-03	D04	HQP021480	NM_003797	EED
QG083-03	D05	HQP004555	NM_001401	EDG2
QG083-03	D06	HQP004504	NM_001946	DUSP6
QG083-03	D07	HQP020444	NM_032122	DTNBP1
QG083-03	D08	HQP004451	NM_000798	DRD5
QG083-03	D09	HQP004450	NM_000797	DRD4
QG083-03	D10	HQP004446	NM_000795	DRD2
QG083-03	D11	HQP004445	NM_000794	DRD1
QG083-03	D12	HQP004441	NM_001386	DPYSL2
QG083-03	E01	HQP006013	NM_015296	DOCK9
QG083-03	E02	HQP009103	NM_006145	DNAJB1
QG083-03	E03	HQP004342	NM_001365	DLG4
QG083-03	E04	HQP007578	NM_001012957	DISC1
QG083-03	E05	HQP001244	NM_022105	DIDO1
QG083-03	E06	HQP003959	NM_152910	DGKH
QG083-03	E07	HQP004032	NM_001352	DBP
QG083-03	E08	HQP003970	NM_001917	DAO
QG083-03	E09	HQP008458	NM_002089	CXCL2
QG083-03	E10	HQP016652	NM_002996	CX3CL1
QG083-03	E11	HQP005986	NM_015267	CUTL2
QG083-03	E12	HQP003464	NM_001012614	CTBP1
QG083-03	F01	HQP006044	NM_015321	CRTC1

QG083-03	F02	HQP002967	NM_001882	CRHBP
QG083-03	F03	HQP002920	NM_001079846	CREBBP
QG083-03	F04	HQP002907	NM_004379	CREB1
QG083-03	F05	HQP002873	NM_000098	CPT2
QG083-03	F06	HQP002788	NM_004374	COX6C
QG083-03	F07	HQP022636	NM_004255	COX5A
QG083-03	F08	HQP002671	NM_000754	COMT
QG083-03	F09	HQP002400	NM_033133	CNP
QG083-03	F10	HQP022927	NM_004898	CLOCK
QG083-03	F11	HQP012083	NM_005602	CLDN11
QG083-03	F12	HQP001829	NM_001824	CKM
QG083-03	G01	HQP001708	NM_001278	CHUK
QG083-03	G02	HQP001644	NM_000746	CHRNA7
QG083-03	G03	HQP001614	NM_000079	CHRNA1
QG083-03	G04	HQP001535	NM_000739	CHRM2
QG083-03	G05	HQP019569	NM_025114	CEP290
QG083-03	G06	HQP021757	NM_001760	CCND3
QG083-03	G07	HQP021754	NM_001759	CCND2
QG083-03	G08	HQP016204	NM_053056	CCND1
QG083-03	G09	HQP053971	NM_021006	CCL3L1
QG083-03	G10	HQP016630	NM_005408	CCL13
QG083-03	G11	HQP021647	NM_000730	CCKAR
QG083-03	G12	HQP020920	NM_005189	CBX2
QG083-03	H01	HGDC		
QG083-03	H02	HGDC		
QG083-03	H03	HQP006940	NM_002046	GAPDH
QG083-03	H04	HQP016381	NM_001101	ACTB
QG083-03	H05	HQP015171	NM_004048	B2M
QG083-03	H06	HQP006171	NM_012423	RPL13A
QG083-03	H07	HQP009026	NM_000194	HPRT1
QG083-03	H08	HQP054253	NR_003286	RN18S1
QG083-03	H09	RT		
QG083-03	H10	RT		
QG083-03	H11	PCR		
QG083-03	H12	PCR		
QG083-04	A01	HQP018966	NM_001080124	CASP8
QG083-04	A02	HQP020427	NM_001226	CASP6
QG083-04	A03	HQP020249	NM_032982	CASP2
QG083-04	A04	HQP000781	NM_006549	CAMKK2
QG083-04	A05	HQP019841	NM_004343	CALR
QG083-04	A06	HQP000445	NM_006085	BPNT1

QG083-04	A07	HQP017234	NM_005180	BMI1
QG083-04	A08	HQP016653	NM_001196	BID
QG083-04	A09	HQP016418	NM_004327	BCR
QG083-04	A10	HQP016211	NM_000633	BCL2
QG083-04	A11	HQP015964	NM_004324	BAX
QG083-04	A12	HQP015917	NM_001188	BAK1
QG083-04	B01	HQP014386	NM_000707	AVPR1B
QG083-04	B02	HQP016578	NM_000332	ATXN1
QG083-04	B03	HQP013091	NM_001003696	ATP5J
QG083-04	B04	HQP011909	NM_001681	ATP2A2
QG083-04	B05	HQP011806	NM_152296	ATP1A3
QG083-04	B06	HQP011736	NM_000051	ATM
QG083-04	B07	HQP014637	NM_018179	ATF7IP
QG083-04	B08	HQP005471	NM_012068	ATF5
QG083-04	B09	HQP011497	NM_000049	ASPA
QG083-04	B10	HQP009801	NM_000044	AR
QG083-04	B11	HQP006348	NM_030882	APOL2
QG083-04	B12	HQP009556	NM_000041	APOE
QG083-04	C01	HQP008934	NM_001160	APAF1
QG083-04	C02	HQP013821	NM_017747	ANKHD1
QG083-04	C03	HQP006359	NM_000698	ALOX5
QG083-04	C04	HQP006356	NM_000697	ALOX12
QG083-04	C05	HQP019261	NM_001077690	ALG9
QG083-04	C06	HQP004991	NM_001014431	AKT1
QG083-04	C07	HQP017203	NM_031279	AGXT2L1
QG083-04	C08	HQP003928	NM_001126	ADSS
QG083-04	C09	HQP003848	NM_005160	ADRBK2
QG083-04	C10	HQP001836	NM_001116	ADCY9
QG083-04	C11	HQP021691	NM_001103	ACTN2
QG083-04	C12	HQP004081	NM_000789	ACE
QG083-04	D01	HQP012188	NM_002580	REG3A
QG083-04	D02	HQP022179	NM_004711	SYNGR1
QG083-04	D03	HQP017910	NM_005992	TBX1
QG083-04	D04	HQP021724	NM_003920	TIMELESS
QG083-04	D05	HQP011291	NM_001008228	MOG
QG083-04	D06	HQP000268	NM_005806	OLIG2
QG083-04	D07	HQP013033	NM_006211	PENK
QG083-04	D08	HQP014583	NM_002731	PRKACB
QG083-04	D09	HQP021643	NM_003895	SYNJ1
QG083-04	D10	HQP018154	NM_003283	TNNT1
QG083-04	D11	HQP022987	NM_004915	ABCG1

QG083-04	D12	HQP000604	NM_001033049	ADARB1
QG083-04	E01	HQP001925	NM_001117	ADCYAP1
QG083-04	E02	HQP003417	NM_000679	ADRA1B
QG083-04	E03	HQP021431	NM_005090	PLA2G4B
QG083-04	E04	HQP012286	NM_001001973	ATP5C1
QG083-04	E05	HQP013049	NM_001002258	ATP5G3
QG083-04	E06	HQP016545	NM_001709	BDNF
QG083-04	E07	HQP015425	NM_020412	CHMP1B
QG083-04	E08	HQP007944	NM_178034	PLA2G4D
QG083-04	E09	HQP006178	NM_013246	CLCF1
QG083-04	E10	HQP002962	NM_000756	CRH
QG083-04	E11	HQP004331	NM_000793	DIO2
QG083-04	E12	HQP004407	NM_001379	DNMT1
QG083-04	F01	HQP004448	NM_000796	DRD3
QG083-04	F02	HQP013311	NM_000304	PMP22
QG083-04	F03	HQP006604	NM_000808	GABRA3
QG083-04	F04	HQP008361	NM_000831	GRIK3
QG083-04	F05	HQP009025	NM_020995	HPR
QG083-04	F06	HQP006867	NM_025015	HSPA12A
QG083-04	F07	HQP009115	NM_000524	HTR1A
QG083-04	F08	HQP009126	NM_000870	HTR4
QG083-04	F09	HQP001124	NM_006839	IMMT
QG083-04	F10	HQP015090	NM_019842	KCNQ5
QG083-04	H05	HQP009026	NM_000194	HPRT0
QG083-04	H04	HQP054253	NR_003285	RN18S0
QG083-04	H03	RT		
QG083-04	H02	RT		
QG083-04	H01	PCR		
QG083-04	H00	PCR		
QG083-04	G05	HQP005996	NM_015277	NEDD4L
QG083-04	G06	HQP011941	NM_001007097	NTRK2
QG083-04	G07	HQP001343	NM_019094	NUDT4
QG083-04	G08	HQP001890	NM_138983	OLIG1
QG083-04	G09	HQP012146	NM_002573	PAFAH1B3
QG083-04	G10	HQP000737	NM_001011513	PDLIM5
QG083-04	G11	HQP013046	NM_002616	PER1
QG083-04	G12	HQP021641	NM_003894	PER2
QG083-04	H01	HGDC		
QG083-04	H02	HGDC		
QG083-04	H03	HQP006940	NM_002046	GAPDH
QG083-04	H04	HQP016381	NM_001101	ACTB

QG083-04	H05	HQP015171	NM_004048	B2M
QG083-04	H06	HQP006171	NM_012423	RPL13A
QG083-04	H07	HQP009026	NM_000194	HPRT1
QG083-04	H08	HQP054253	NR_003286	RN18S1
QG083-04	H09	RT		
QG083-04	H10	RT		
QG083-04	H11	PCR		
QG083-04	H12	PCR		

---

### Limited Use License

Following terms and conditions apply to use of ExProfile™ Human Mood Disorder Related Gene qPCR Array (the Product). If the terms and conditions are not acceptable, the Product in its entirety must be returned to GeneCopoeia within 5 calendar days. A limited End-User license is granted to the purchaser of the Product. The Product shall be used by the purchaser for internal research purposes only. The Product is expressly not designed, intended, or warranted for use in humans or for therapeutic or diagnostic use. The Product must not be resold, repackaged or modified for resale, or used to manufacture commercial products or deliver information obtained in service without prior written consent from GeneCopoeia. This Product should be used in accordance with the NIH guidelines developed for recombinant DNA and genetic research. Use of any part of the Product constitutes acceptance of the above terms.

### Limited Warranty

GeneCopoeia warrants that the Product meets the specifications described in the accompanying Product Datasheet. If it is proven to the satisfaction of GeneCopoeia that the Product fails to meet these specifications, GeneCopoeia will replace the Product. In the event a replacement cannot be provided, GeneCopoeia will provide the purchaser with a refund. This limited warranty shall not extend to anyone other than the original purchaser of the Product. Notice of nonconforming products must be made to GeneCopoeia within 30 days of receipt of the Product. GeneCopoeia's liability is expressly limited to replacement of Product or a refund limited to the actual purchase price.

GeneCopoeia's liability does not extend to any damages arising from use or improper use of the Product, or losses associated with the use of additional materials or reagents. This limited warranty is the sole and exclusive warranty. GeneCopoeia does not provide any other warranties of any kind, expressed or implied, including the merchantability or fitness of the Product for a particular purpose.

GeneCopoeia is committed to providing our customers with high-quality products. If you should have any questions or concerns about any GeneCopoeia products, please contact us at 301-762-0888.

© 2016 GeneCopoeia, Inc.

GeneCopoeia, Inc.  
9620 Medical Center Drive, Suite 101  
Rockville, MD 20850  
+1 (301) 762-0888  
+1 (866) 360-9531  
[inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)