

## ExProfile™ Human JAK/STAT Signaling Related Gene qPCR Array

For focused group profiling of human JAK/STAT signaling related gene expression

Cat. No. QG034-A (1 x 96-well plate, Format A)

Cat. No. QG034-B (1 x 96-well plate, Format B)

Cat. No. QG034-C (1 x 96-well plate, Format C)

Cat. No. QG034-D (1 x 96-well plate, Format D)

Cat. No. QG034-E (1 x 96-well plate, Format E)

Plates available individually or as a set of 6. Each set contains 84 unique gene primer pairs deposited in one 96-well plate.

### Introduction

The ExProfile human JAK/STAT signaling related gene qPCR array profiles the expression of 84 human genes related to JAK/STAT-mediated signal transduction. These genes are carefully chosen for their close pathway correlation based on a thorough literature search of peer-reviewed publications. This array allows researchers to study pathway-related genes to gain understanding of their roles in the JAK/STAT signaling pathway.

- QG034 plate 01: 84 unique gene PCR primer pairs

### Shipping and storage conditions

Shipped at room temperature

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 receipt, please check to make sure that the correct array format was ordered to ensure 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	A2M	SH2B2	BCL2L1	CCND1	CDKN1A	CRK	CRP	CSF1R	CSF2RB	CXCL9	EGFR	EPOR
B	F2	F2R	FAS	FCER1A	FCGR1A	ISG15	GATA3	GBP1	GHR	IFNAR1	IFNG	IFNGR1
C	IL10RA	IL20	IL2RA	IL2RG	IL4	IL4R	IL6ST	INSR	IRF1	JAK1	JAK2	JAK3
D	JUN	JUNB	MMP3	MPL	MYC	NFKB1	NOS2A	NR3C1	OAS1	PDGFRA	PIAS1	PIAS2
E	PPP2R1A	PRLR	PTPN1	PTPRC	SLA2	SMAD1	SMAD2	SMAD3	SMAD4	SMAD5	SOCS1	SOCS2
F	SOCS3	SOCS4	SOCS5	SP1	SPI1	SRC	STAM	STAT1	STAT2	STAT3	STAT4	STAT5A
G	STAT5B	STAT6	TYK2	YY1	HPRT1	FAS	IL6ST	PTPRC	STAT3	CEBPB	OSM	STUB1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG034 plate 01

- **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
QG034-01	A01	HQP008678	NM_000014	A2M
QG034-01	A02	HQP000726	NM_020979	SH2B2
QG034-01	A03	HQP016238	NM_138578	BCL2L1
QG034-01	A04	HQP016204	NM_053056	CCND1
QG034-01	A05	HQP000331	NM_000389	CDKN1A
QG034-01	A06	HQP002988	NM_016823	CRK
QG034-01	A07	HQP002994	NM_000567	CRP
QG034-01	A08	HQP003158	NM_005211	CSF1R
QG034-01	A09	HQP003170	NM_000395	CSF2RB
QG034-01	A10	HQP011220	NM_002416	CXCL9
QG034-01	A11	HQP004605	NM_005228	EGFR
QG034-01	A12	HQP004958	NM_000121	EPOR
QG034-01	B01	HQP005052	NM_000506	F2
QG034-01	B02	HQP005053	NM_001992	F2R
QG034-01	B03	HQP009651	NM_000043	FAS
QG034-01	B04	HQP005238	NM_002001	FCER1A
QG034-01	B05	HQP005251	NM_000566	FCGR1A
QG034-01	B06	HQP023013	NM_005101	ISG15
QG034-01	B07	HQP007166	NM_002051	GATA3
QG034-01	B08	HQP007221	NM_002053	GBP1
QG034-01	B09	HQP007395	NM_000163	GHR
QG034-01	B10	HQP009458	NM_000629	IFNAR1
QG034-01	B11	HQP009467	NM_000619	IFNG
QG034-01	B12	HQP009469	NM_000416	IFNGR1
QG034-01	C01	HQP009686	NM_001558	IL10RA
QG034-01	C02	HQP012157	NM_018724	IL20
QG034-01	C03	HQP009650	NM_000417	IL2RA
QG034-01	C04	HQP009659	NM_000206	IL2RG
QG034-01	C05	HQP009662	NM_000589	IL4
QG034-01	C06	HQP009664	NM_000418	IL4R
QG034-01	C07	HQP009674	NM_002184	IL6ST
QG034-01	C08	HQP009764	NM_000208	INSR
QG034-01	C09	HQP009778	NM_002198	IRF1
QG034-01	C10	HQP009849	NM_002227	JAK1
QG034-01	C11	HQP009850	NM_004972	JAK2
QG034-01	C12	HQP009851	NM_000215	JAK3
QG034-01	D01	HQP009853	NM_002228	JUN
QG034-01	D02	HQP009854	NM_002229	JUNB
QG034-01	D03	HQP011257	NM_002422	MMP3
QG034-01	D04	HQP011308	NM_005373	MPL
QG034-01	D05	HQP011597	NM_002467	MYC

Product Data Sheet

QG034-01	D06	HQP011807	NM_003998	NFKB1
QG034-01	D07	HQP011866	NM_000625	NOS2A
QG034-01	D08	HQP008401	NM_000176	NR3C1
QG034-01	D09	HQP011984	NM_002534	OAS1
QG034-01	D10	HQP012866	NM_006206	PDGFRA
QG034-01	D11	HQP021271	NM_016166	PIAS1
QG034-01	D12	HQP022005	NM_004671	PIAS2
QG034-01	E01	HQP014147	NM_014225	PPP2R1A
QG034-01	E02	HQP015027	NM_000949	PRLR
QG034-01	E03	HQP015828	NM_002827	PTPN1
QG034-01	E04	HQP015908	NM_002838	PTPRC
QG034-01	E05	HQP020528	NM_175077	SLA2
QG034-01	E06	HQP054006	NM_005900	SMAD1
QG034-01	E07	HQP054007	NM_005901	SMAD2
QG034-01	E08	HQP010960	NM_005902	SMAD3
QG034-01	E09	HQP010961	NM_005359	SMAD4
QG034-01	E10	HQP054008	NM_005903	SMAD5
QG034-01	E11	HQP021399	NM_003745	SOCS1
QG034-01	E12	HQP021602	NM_003877	SOCS2
QG034-01	F01	HQP021889	NM_003955	SOCS3
QG034-01	F02	HQP002181	NM_080867	SOCS4
QG034-01	F03	HQP054032	NM_144949	SOCS5
QG034-01	F04	HQP017640	NM_138473	SP1
QG034-01	F05	HQP017663	NM_003120	SPI1
QG034-01	F06	HQP017696	NM_005417	SRC
QG034-01	F07	HQP019637	NM_003473	STAM
QG034-01	F08	HQP017764	NM_007315	STAT1
QG034-01	F09	HQP017766	NM_005419	STAT2
QG034-01	F10	HQP017767	NM_003150	STAT3
QG034-01	F11	HQP017770	NM_003151	STAT4
QG034-01	F12	HQP017771	NM_003152	STAT5A
QG034-01	G01	HQP017774	NM_012448	STAT5B
QG034-01	G02	HQP017775	NM_003153	STAT6
QG034-01	G03	HQP018340	NM_003331	TYK2
QG034-01	G04	HQP018570	NM_003403	YY1
QG034-01	G05	HQP009026	NM_000194	HPRT1
QG034-01	G06	HQP009653	NM_152872	FAS
QG034-01	G07	HQP009675	NM_175767	IL6ST
QG034-01	G08	HQP015909	NM_080921	PTPRC
QG034-01	G09	HQP017769	NM_213662	STAT3
QG034-01	G10	HQP000623	NM_005194	CEBPB
QG034-01	G11	HQP012081	NM_020530	OSM
QG034-01	G12	HQP000334	NM_005861	STUB1
QG034-01	H01	HGDC		
QG034-01	H02	HGDC		
QG034-01	H03	HQP006940	NM_002046	GAPDH
QG034-01	H04	HQP016381	NM_001101	ACTB

Product Data Sheet

QG034-01	H05	HQP015171	NM_004048	B2M
QG034-01	H06	HQP006171	NM_012423	RPL13A
QG034-01	H07	HQP009026	NM_000194	HPRT1
QG034-01	H08	HQP054253	NR_003286	RN18S1
QG034-01	H09	RT		
QG034-01	H10	RT		
QG034-01	H11	PCR		
QG034-01	H12	PCR		

### Limited Use License

Following terms and conditions apply to use of ExProfile™ Human JAK/STAT Signaling 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)