

ExProfile[™] Human Ubiquitination (Ubiquitylation) Related Gene qPCR Array

For focused group profiling of human ubiquitination (ubiquitylation) genes expression

Cat. No. QG059-A (1 x 96-well plate, Format A) Cat. No. QG059-B (1 x 96-well plate, Format B) Cat. No. QG059-C (1 x 96-well plate, Format C) Cat. No. QG059-D (1 x 96-well plate, Format D) Cat. No. QG059-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 ubiquitination (ubiquitylation) related gene qPCR array profiles the expression of 84 human genes related to regulated degradation of cellular proteins or foreign proteins by the ubiquitinproteasome system. These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including genes that encode ubiquitin-activating enzymes, ubiquitin-conjugating enzymes and ubiquitin-protein ligases. This array allows researchers to study the pathway-related genes to gain understanding of their roles in ubiquitin degradation.

• QG059 plate 01: 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

GeneCopoeia provides five qPCR array formats (A, B, C, D, and E) suitable for use with the following realtime 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
Α	UBE2C	UBE2C	FBXO4	PARK2	HPRT1	WWP1	VHL	UBR2	UBR1	UBE4B	UBE2Z	UBE2W
В	UBE2T	UBE2R2	UBE2Q1	UBE2N	UBE2M	UBE2L3	HIP2	UBE2J2	UBE2J1	UBE2I	UBE2G2	UBE2G1
С	UBE2E3	UBE2E2	UBE2D3	UBE2D2	UBE2D1	UBE2C	UBE1L2	UBE1C	SAE2	UBE1	TP53	Kua
D	SMURF2	SMURF1	SKP2	SKP1A	SAE1	RNF123	PARK2	PARC	APPBP1	MIB1	MDM2	41703
Е	HUWE1	HECW2	HECW1	FBXW10	FBXO4	FBX03	DZIP3	DDB1	CUL7	CUL5	CUL4B	CUL4A
F	CUL3	CUL2	CUL1	CDC34	CBL	BTRC	BRCC3	BRCA1	BARD1	ATG7	ARIH1	ANAPC2
G	UBE2E1	UBE2H	ANAPC11	FBX031	UBE2S	NEDD8	RFWD2	RNF148	SYVN1	UBE1DC1	UBE2A	UBE2B
Н	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG059 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 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 predeposited DNA template with its specific pre-deposited primer pairs.



Gene primer list

Plate	Position	Catalog No. of Primer	Catalog No. Accession No. of Primer of Gene	
QG059-01	A01	HQP001215	NM_181799	UBE2C
QG059-01	A02	HQP001217	NM_181801	UBE2C
QG059-01	A03	HQP007187	NM_033484	FBXO4
QG059-01	A04	HQP012195	NM_013988	PARK2
QG059-01	A05	HQP009026	NM_000194	HPRT1
QG059-01	A06	HQP001203	NM_007013	WWP1
QG059-01	A07	HQP018485	NM_000551	VHL
QG059-01	A08	HQP005973	NM_015255	UBR2
QG059-01	A09	HQP004660	NM_174916	UBR1
QG059-01	A10	HQP000338	NM 006048	UBE4B
QG059-01	A11	HQP017382	NM_023079	UBE2Z
QG059-01	A12	HQP014272	NM 018299	UBE2W
QG059-01	B01	HQP008400	NM 014176	UBE2T
QG059-01	B02	HQP013865	NM 017811	UBE2R2
QG059-01	B03	HQP014486	NM 017582	UBE2Q1
QG059-01	B04	HQP018383	NM 003348	UBE2N
QG059-01	B05	HQP021936	NM 003969	UBE2M
QG059-01	B06	HQP018381	NM 003347	UBE2L3
QG059-01	B07	HQP008834	NM 005339	HIP2
QG059-01	B08	HQP054042	NM 194458	UBE2J2
QG059-01	B09	HQP012761	NM 016021	UBE2J1
QG059-01	B10	HQP018379	NM 003345	UBE2I
QG059-01	B11	HQP018376	NM 182688	UBE2G2
QG059-01	B12	HQP018373	NM 003342	UBE2G1
QG059-01	C01	HQP000577	NM 006357	UBE2E3
QG059-01	C02	HQP018372	NM 152653	UBE2E2
QG059-01	C03	HQP018369	NM 181893	UBE2D3
QG059-01	C04	HQP018366	NM 181838	UBE2D2
QG059-01	C05	HQP018364	NM 003338	UBE2D1
QG059-01	C06	HQP001218	NM 181803	UBE2C
QG059-01	C07	HQP014208	NM 018227	UBE1L2
QG059-01	C08	HQP021932	NM 003968	UBE1C
QG059-01	C09	HQP000071	NM 005499	SAE2
QG059-01	C10	HQP018357	NM 003334	UBE1
QG059-01	C11	HQP018175	NM_000546	TP53
QG059-01	C12	HQP010193	NM 199129	Kua
QG059-01	D01	HQP017122	NM 022739	SMURF2
QG059-01	D02	HQP015449	NM 020429	SMURF1
QG059-01	D03	HQP017300	NM 005983	SKP2
QG059-01	D04	HQP017289	NM 006930	SKP1A
QG059-01	D05	HQP000072	NM 005500	SAE1
QG059-01	D06	HQP016673	NM_022064	RNF123



QG059-01	D07	HQP012193	NM 004562	PARK2
QG059-01	D08	HQP005770	NM_015089	PARC
QG059-01	D09	HQP021668	NM_003905	APPBP1
QG059-01	D10	HQP015685	NM_020774	MIB1
QG059-01	D11	HQP011135	NM_002392	MDM2
QG059-01	D12	HQP013642	NM_017824	5-Mar
QG059-01	E01	HQP000096	NM_031407	HUWE1
QG059-01	E02	HQP015673	NM_020760	HECW2
QG059-01	E03	HQP005726	NM_015052	HECW1
QG059-01	E04	HQP000620	NM_031456	FBXW10
QG059-01	E05	HQP007186	NM_012176	FBXO4
QG059-01	E06	HQP007188	NM_012175	FBXO3
QG059-01	E07	HQP023052	NM_014648	DZIP3
QG059-01	E08	HQP004111	NM_001923	DDB1
QG059-01	E09	HQP023238	NM_014780	CUL7
QG059-01	E10	HQP019709	NM_003478	CUL5
QG059-01	E11	HQP020764	NM_003588	CUL4B
QG059-01	E12	HQP020774	NM_003589	CUL4A
QG059-01	F01	HQP020782	NM_003590	CUL3
QG059-01	F02	HQP020788	NM_003591	CUL2
QG059-01	F03	HQP020798	NM_003592	CUL1
QG059-01	F04	HQP023443	NM_004359	CDC34
QG059-01	F05	HQP021430	NM_005188	CBL
QG059-01	F06	HQP021753	NM_033637	BTRC
QG059-01	F07	HQP018988	NM_024332	BRCC3
QG059-01	F08	HQP017713	NM_007294	BRCA1
QG059-01	F09	HQP015946	NM_000465	BARD1
QG059-01	F10	HQP000641	NM_006395	ATG7
QG059-01	F11	HQP006752	NM_005744	ARIH1
QG059-01	F12	HQP008560	NM_013366	ANAPC2
QG059-01	G01	HQP018371	NM_182666	UBE2E1
QG059-01	G02	HQP018378	NM_182697	UBE2H
QG059-01	G03	HQP012810	NM_001002244	ANAPC11
QG059-01	G04	HQP019258	NM_024735	FBXO31
QG059-01	G05	HQP007682	NM_014501	UBE2S
QG059-01	G06	HQP011745	NM_006156	NEDD8
QG059-01	G07	HQP016911	NM_022457	RFWD2
QG059-01	G08	HQP010064	NM_198085	RNF148
QG059-01	G09	HQP020731	NM_172230	SYVN1
QG059-01	G10	HQP019348	NM_198329	UBE1DC1
QG059-01	G11	HQP018359	NM_003336	UBE2A
QG059-01	G12	HQP018363	NM_003337	UBE2B
QG059-01	H01	HGDC		
QG059-01	H02	HGDC		
QG059-01	H03	HQP006940	NM_002046	GAPDH
QG059-01	H04	HQP016381	NM_001101	ACTB



QG059-01	H05	HQP015171	NM_004048	B2M
QG059-01	H06	HQP006171	NM_012423	RPL13A
QG059-01	H07	HQP009026	NM_000194	HPRT1
QG059-01	H08	HQP054253	NR_003286	RN18S1
QG059-01	H09	RT		
QG059-01	H10	RT		
QG059-01	H11	PCR		
QG059-01	H12	PCR		



Limited Use License

Following terms and conditions apply to use of ExProfile[™] Human Ubiquitination (Ubiquitylation) 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). QG059-160202