

ExProfile™ Human DNA Damage and Repair Related Gene qPCR Array

For focused group profiling of human DNA damage and repair genes expression

Cat. No. QG012-A (1 x 96-well plate, Format A)

Cat. No. QG012-B (1 x 96-well plate, Format B)

Cat. No. QG012-C (1 x 96-well plate, Format C)

Cat. No. QG012-D (1 x 96-well plate, Format D)

Cat. No. QG012-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 DNA damage and repair related gene qPCR array profiles the expression of 84 human genes related to DNA damage and repair. These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including genes involved in apoptosis, cell cycle, various DNA repair processes. This array allows researchers to study the related genes to gain understanding of their roles in the process of DNA damage and repair.

- QG012 plate 01: 84 unique gene PCR primer pairs

Shipping and storage condition

Shipped at room temperature

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
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	XRCC4	XRCC3	XRCC2	XRCC1	XPC	XAB2	UHRF1	TP53BP1	TP53	TDP1	TDG	SMUG1
B	REV3L	REV1	RAD52	RAD51C	RAD51	RAD23B	RAD23A	RAD21	RAD18	RAD17	RAD1	POLL
C	POLK	POLI	POLE	POLD1	POLB	PMS2	PCNA	NEIL3	NEIL2	NEIL1	MSH6	MSH3
D	MSH2	MRE11A	MPG	MLL	MLH1	MBD4	LIG4	LIG1	HUS1	GTF2H4	GTF2H3	GTF2H1
E	GADD45G	GADD45A	FEN1	FANCC	FANCA	ERCC4	ERCC3	ERCC2	ERCC1	DDB2	DDB1	DCLRE1C
F	DCLRE1B	DCLRE1A	BRCA2	BRCA1	BLM	ATRX	ATM	APEX2	APEX1	NTHL1	FANCG	H2AFX
G	UNG	XPA	MSH5	MUTYH	NBN	PARP2	PNKP	POLH	PRKDC	MSH4	RFC5	TREX1
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG012 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 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
QG012-01	A01	HQP018565	NM_003401	XRCC4
QG012-01	A02	HQP018564	NM_005432	XRCC3
QG012-01	A03	HQP018563	NM_005431	XRCC2
QG012-01	A04	HQP018562	NM_006297	XRCC1
QG012-01	A05	HQP018556	NM_004628	XPC
QG012-01	A06	HQP015272	NM_020196	XAB2
QG012-01	A07	HQP008445	NM_001048201	UHRF1
QG012-01	A08	HQP018176	NM_005657	TP53BP1
QG012-01	A09	HQP018175	NM_000546	TP53
QG012-01	A10	HQP014691	NM_001008744	TDP1
QG012-01	A11	HQP017996	NM_003211	TDG
QG012-01	A12	HQP006236	NM_014311	SMUG1
QG012-01	B01	HQP016223	NM_002912	REV3L
QG012-01	B02	HQP012755	NM_001037872	REV1
QG012-01	B03	HQP016087	NM_134424	RAD52
QG012-01	B04	HQP016079	NM_002876	RAD51C
QG012-01	B05	HQP016077	NM_002875	RAD51
QG012-01	B06	HQP016076	NM_002874	RAD23B
QG012-01	B07	HQP016075	NM_005053	RAD23A
QG012-01	B08	HQP016074	NM_006265	RAD21
QG012-01	B09	HQP015190	NM_020165	RAD18
QG012-01	B10	HQP016070	NM_002873	RAD17
QG012-01	B11	HQP015947	NM_002853	RAD1
QG012-01	B12	HQP007689	NM_013274	POLL
QG012-01	C01	HQP012725	NM_016218	POLK
QG012-01	C02	HQP001399	NM_007195	POLI
QG012-01	C03	HQP013425	NM_006231	POLE
QG012-01	C04	HQP013423	NM_002691	POLD1
QG012-01	C05	HQP013422	NM_002690	POLB
QG012-01	C06	HQP013352	NM_000535	PMS2
QG012-01	C07	HQP012420	NM_002592	PCNA
QG012-01	C08	HQP014222	NM_018248	NEIL3
QG012-01	C09	HQP006460	NM_145043	NEIL2
QG012-01	C10	HQP019141	NM_024608	NEIL1
QG012-01	C11	HQP008493	NM_000179	MSH6
QG012-01	C12	HQP011492	NM_002439	MSH3
QG012-01	D01	HQP011491	NM_000251	MSH2

QG012-01	D02	HQP011320	NM_005590	MRE11A
QG012-01	D03	HQP011304	NM_001015052	MPG
QG012-01	D04	HQP011241	NM_005933	MLL
QG012-01	D05	HQP011235	NM_000249	MLH1
QG012-01	D06	HQP021736	NM_003925	MBD4
QG012-01	D07	HQP010613	NM_002312	LIG4
QG012-01	D08	HQP010609	NM_000234	LIG1
QG012-01	D09	HQP009138	NM_004507	HUS1
QG012-01	D10	HQP008505	NM_001517	GTF2H4
QG012-01	D11	HQP008504	NM_001516	GTF2H3
QG012-01	D12	HQP008502	NM_005316	GTF2H1
QG012-01	E01	HQP001042	NM_006705	GADD45G
QG012-01	E02	HQP004125	NM_001924	GADD45A
QG012-01	E03	HQP005391	NM_004111	FEN1
QG012-01	E04	HQP005083	NM_000136	FANCC
QG012-01	E05	HQP005081	NM_000135	FANCA
QG012-01	E06	HQP004984	NM_005236	ERCC4
QG012-01	E07	HQP004983	NM_000122	ERCC3
QG012-01	E08	HQP004976	NM_000400	ERCC2
QG012-01	E09	HQP004974	NM_001983	ERCC1
QG012-01	E10	HQP004114	NM_000107	DDB2
QG012-01	E11	HQP004111	NM_001923	DDB1
QG012-01	E12	HQP016982	NM_001033855	DCLRE1C
QG012-01	F01	HQP017213	NM_022836	DCLRE1B
QG012-01	F02	HQP023381	NM_014881	DCLRE1A
QG012-01	F03	HQP017753	NM_000059	BRCA2
QG012-01	F04	HQP017713	NM_007294	BRCA1
QG012-01	F05	HQP016844	NM_000057	BLM
QG012-01	F06	HQP013636	NM_000489	ATRX
QG012-01	F07	HQP011736	NM_000051	ATM
QG012-01	F08	HQP007655	NM_014481	APEX2
QG012-01	F09	HQP009061	NM_001641	APEX1
QG012-01	F10	HQP011937	NM_002528	NTHL1
QG012-01	F11	HQP005099	NM_004629	FANCG
QG012-01	F12	HQP008703	NM_002105	H2AFX
QG012-01	G01	HQP018427	NM_003362	UNG
QG012-01	G02	HQP018555	NM_000380	XPA
QG012-01	G03	HQP011494	NM_002441	MSH5
QG012-01	G04	HQP011576	NM_001048171	MUTYH
QG012-01	G05	HQP011686	NM_001024688	NBN
QG012-01	G06	HQP000048	NM_001042618	PARP2

QG012-01	G07	HQP001521	NM_007254	PNKP
QG012-01	G08	HQP013429	NM_006502	POLH
QG012-01	G09	HQP014837	NM_001081640	PRKDC
QG012-01	G10	HQP011493	NM_002440	MSH4
QG012-01	G11	HQP016230	NM_007370	RFC5
QG012-01	G12	HQP001503	NM_016381	TREX1
QG012-01	H01	HGDC		
QG012-01	H02	HGDC		
QG012-01	H03	HQP006940	NM_002046	GAPDH
QG012-01	H04	HQP016381	NM_001101	ACTB
QG012-01	H05	HQP015171	NM_004048	B2M
QG012-01	H06	HQP006171	NM_012423	RPL13A
QG012-01	H07	HQP009026	NM_000194	HPRT1
QG012-01	H08	HQP054253	NR_003286	RN18S1
QG012-01	H09	RT		
QG012-01	H10	RT		
QG012-01	H11	PCR		
QG012-01	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Human DNA Damage and Repair 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