

ExProfile™ Human Hypoxia Signaling Related Gene qPCR Array

For focused group profiling of human hypoxia signaling genes expression

Cat. No. QG026-A (1 x 96-well plate, Format A)

Cat. No. QG026-B (1 x 96-well plate, Format B)

Cat. No. QG026-C (1 x 96-well plate, Format C)

Cat. No. QG026-D (1 x 96-well plate, Format D)

Cat. No. QG026-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 hypoxia signaling related gene qPCR array profiles the expression of 84 human genes related to hypoxia signaling. 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 the response to hypoxia and oxidative stress, as well as physiological and pathophysiological processes induced by hypoxia. This array allows researchers to study the related genes to gain understanding of their roles in the functioning and characterization of hypoxia signaling pathway.

- QG026 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

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 | SPTBN1 | PPARA | IL6ST | HPRT1 | CASP1 | VEGFA | UCP2 | TUBA4A | TST | TH | SPTBN1 | SNRP70 |
| B | SLC2A4 | SLC2A1 | SAE1 | RPS7 | RPS2 | RPL32 | RPL28 | RARA | PSMB3 | PRKAA1 | PPARA | PLOD3 |
| C | PLAU | PIP3-E | NUDT2 | NPY | NOTCH1 | NOS2A | MYBL2 | MT3 | MAN2B1 | LEP | LCT | KIT |
| D | KHSRP | IQGAP1 | IL6ST | IL6 | IL1A | IGFBP1 | IGF2 | HYOU1 | HTATIP | HMOX1 | HK2 | HIF1AN |
| E | HIF1A | HBB | GPX1 | GPI | GNA11 | EPAS1 | EP300 | ENO1 | EEF1A1 | ECE1 | DCTN2 | DAPK3 |
| F | CYGB | CSTB | CREBBP | COL1A1 | CHGA | CDC42 | CASP1 | CA1 | BIRC5 | BHLHB2 | BAX | ARNT2 |
| G | ANGPTL4 | AGTPBP1 | ADM | AGPAT2 | DR1 | EPO | MOCS3 | ARD1A | CAT | SSSCA1 | HIF3A | PDIA2 |
| H | HGDC | HGDC | GAPDH | ACTB | B2M | RPL13A | HPRT1 | RN18S1 | RT | RT | PCR | PCR |

Figure1. Illustration of QG026 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 |
|----------|----------|-----------------------|-----------------------|--------|
| QG026-01 | A01 | HQP017693 | NM_178313 | SPTBN1 |
| QG026-01 | A02 | HQP013618 | NM_032644 | PPARA |
| QG026-01 | A03 | HQP009675 | NM_175767 | IL6ST |
| QG026-01 | A04 | HQP009026 | NM_000194 | HPRT1 |
| QG026-01 | A05 | HQP020210 | NM_033294 | CASP1 |
| QG026-01 | A06 | HQP018481 | NM_003376 | VEGFA |
| QG026-01 | A07 | HQP018403 | NM_003355 | UCP2 |
| QG026-01 | A08 | HQP018298 | NM_006000 | TUBA4A |
| QG026-01 | A09 | HQP018280 | NM_003312 | TST |
| QG026-01 | A10 | HQP018064 | NM_000360 | TH |
| QG026-01 | A11 | HQP017692 | NM_003128 | SPTBN1 |
| QG026-01 | A12 | HQP017587 | NM_003089 | SNRP70 |
| QG026-01 | B01 | HQP053960 | NM_001042 | SLC2A4 |
| QG026-01 | B02 | HQP017350 | NM_006516 | SLC2A1 |
| QG026-01 | B03 | HQP000072 | NM_005500 | SAE1 |
| QG026-01 | B04 | HQP053973 | NM_001011 | RPS7 |
| QG026-01 | B05 | HQP016454 | NM_002952 | RPS2 |
| QG026-01 | B06 | HQP016435 | NM_000994 | RPL32 |
| QG026-01 | B07 | HQP016432 | NM_000991 | RPL28 |
| QG026-01 | B08 | HQP016114 | NM_000964 | RARA |
| QG026-01 | B09 | HQP015241 | NM_002795 | PSMB3 |
| QG026-01 | B10 | HQP014530 | NM_006251 | PRKAA1 |
| QG026-01 | B11 | HQP054001 | NM_005036 | PPARA |
| QG026-01 | B12 | HQP021796 | NM_001084 | PLOD3 |
| QG026-01 | C01 | HQP013204 | NM_002658 | PLAU |
| QG026-01 | C02 | HQP006987 | NM_015553 | PIP3-E |
| QG026-01 | C03 | HQP008954 | NM_001161 | NUDT2 |
| QG026-01 | C04 | HQP011874 | NM_000905 | NPY |
| QG026-01 | C05 | HQP011873 | NM_017617 | NOTCH1 |
| QG026-01 | C06 | HQP011866 | NM_000625 | NOS2A |
| QG026-01 | C07 | HQP011593 | NM_002466 | MYBL2 |
| QG026-01 | C08 | HQP011539 | NM_005954 | MT3 |
| QG026-01 | C09 | HQP011005 | NM_000528 | MAN2B1 |
| QG026-01 | C10 | HQP010581 | NM_000230 | LEP |
| QG026-01 | C11 | HQP010570 | NM_002299 | LCT |
| QG026-01 | C12 | HQP010099 | NM_000222 | KIT |
| QG026-01 | D01 | HQP021301 | NM_003685 | KHSRP |
| QG026-01 | D02 | HQP021588 | NM_003870 | IQGAP1 |
| QG026-01 | D03 | HQP009674 | NM_002184 | IL6ST |
| QG026-01 | D04 | HQP009670 | NM_000600 | IL6 |
| QG026-01 | D05 | HQP009640 | NM_000575 | IL1A |
| QG026-01 | D06 | HQP009539 | NM_000596 | IGFBP1 |

| | | | | |
|----------|-----|-----------|--------------|---------|
| QG026-01 | D07 | HQP009529 | NM_000612 | IGF2 |
| QG026-01 | D08 | HQP000633 | NM_006389 | HYOU1 |
| QG026-01 | D09 | HQP000630 | NM_006388 | HTATIP |
| QG026-01 | D10 | HQP008898 | NM_002133 | HMOX1 |
| QG026-01 | D11 | HQP008843 | NM_000189 | HK2 |
| QG026-01 | D12 | HQP014564 | NM_017902 | HIF1AN |
| QG026-01 | E01 | HQP008831 | NM_001530 | HIF1A |
| QG026-01 | E02 | HQP008729 | NM_000518 | HBB |
| QG026-01 | E03 | HQP008279 | NM_000581 | GPX1 |
| QG026-01 | E04 | HQP007814 | NM_000175 | GPI |
| QG026-01 | E05 | HQP007743 | NM_002067 | GNA11 |
| QG026-01 | E06 | HQP004903 | NM_001430 | EPAS1 |
| QG026-01 | E07 | HQP004897 | NM_001429 | EP300 |
| QG026-01 | E08 | HQP004860 | NM_001428 | ENO1 |
| QG026-01 | E09 | HQP004570 | NM_001402 | EEF1A1 |
| QG026-01 | E10 | HQP004537 | NM_001397 | ECE1 |
| QG026-01 | E11 | HQP000649 | NM_006400 | DCTN2 |
| QG026-01 | E12 | HQP003982 | NM_001348 | DAPK3 |
| QG026-01 | F01 | HQP001679 | NM_134268 | CYGB |
| QG026-01 | F02 | HQP003388 | NM_000100 | CSTB |
| QG026-01 | F03 | HQP002921 | NM_004380 | CREBBP |
| QG026-01 | F04 | HQP002462 | NM_000088 | COL1A1 |
| QG026-01 | F05 | HQP001309 | NM_001275 | CHGA |
| QG026-01 | F06 | HQP053981 | NM_001791 | CDC42 |
| QG026-01 | F07 | HQP020208 | NM_033292 | CASP1 |
| QG026-01 | F08 | HQP018641 | NM_001738 | CA1 |
| QG026-01 | F09 | HQP009101 | NM_001168 | BIRC5 |
| QG026-01 | F10 | HQP021270 | NM_003670 | BHLHB2 |
| QG026-01 | F11 | HQP015964 | NM_004324 | BAX |
| QG026-01 | F12 | HQP023361 | NM_014862 | ARNT2 |
| QG026-01 | G01 | HQP012432 | NM_001039667 | ANGPTL4 |
| QG026-01 | G02 | HQP005957 | NM_015239 | AGTPBP1 |
| QG026-01 | G03 | HQP002765 | NM_001124 | ADM |
| QG026-01 | G04 | HQP000663 | NM_006412 | AGPAT2 |
| QG026-01 | G05 | HQP004443 | NM_001938 | DR1 |
| QG026-01 | G06 | HQP004956 | NM_000799 | EPO |
| QG026-01 | G07 | HQP007660 | NM_014484 | MOCS3 |
| QG026-01 | G08 | HQP020079 | NM_003491 | ARD1A |
| QG026-01 | G09 | HQP020946 | NM_001752 | CAT |
| QG026-01 | G10 | HQP000642 | NM_006396 | SSSCA1 |
| QG026-01 | G11 | HQP016928 | NM_152794 | HIF3A |
| QG026-01 | G12 | HQP017106 | NM_006849 | PDIA2 |
| QG026-01 | H01 | HGDC | | |
| QG026-01 | H02 | HGDC | | |
| QG026-01 | H03 | HQP006940 | NM_002046 | GAPDH |
| QG026-01 | H04 | HQP016381 | NM_001101 | ACTB |

| | | | | |
|----------|-----|-----------|-----------|--------|
| QG026-01 | H05 | HQP015171 | NM_004048 | B2M |
| QG026-01 | H06 | HQP006171 | NM_012423 | RPL13A |
| QG026-01 | H07 | HQP009026 | NM_000194 | HPRT1 |
| QG026-01 | H08 | HQP054253 | NR_003286 | RN18S1 |
| QG026-01 | H09 | RT | | |
| QG026-01 | H10 | RT | | |
| QG026-01 | H11 | PCR | | |
| QG026-01 | H12 | PCR | | |

Limited Use License

Following terms and conditions apply to use of ExProfile™ Human Hypoxia 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

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). QG026-160202

Copyright © 2016 GeneCopoeia, Inc.