

ExProfile™ Human Targets of Wnt/ β -catenin Signaling Related Gene qPCR Array

For focused group profiling of human targets of Wnt/ β -catenin signaling genes expression

Cat. No. QG060-A (1 x 96-well plate, Format A)

Cat. No. QG060-B (1 x 96-well plate, Format B)

Cat. No. QG060-C (1 x 96-well plate, Format C)

Cat. No. QG060-D (1 x 96-well plate, Format D)

Cat. No. QG060-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 targets of Wnt/ β -catenin signaling related gene qPCR array profiles the expression of 84 human genes related to Wnt/ β -catenin signaling pathway. These genes are carefully chosen for their close correlation based on a thorough literature search of peer-reviewed publications, mainly including Wnt/ β -catenin signaling target genes. This array allows researchers to study the pathway-related genes to gain understanding of their roles in Wnt/ β -catenin signaling pathway.

- QG060 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

GeneCopeia 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	WISP2	WISP1	VEGFC	VEGFA	VCAN	TWIST1	TLE1	TIAM1	TGFB2	TCF7L2	TCF7L1	TCF7
B	TCF4	TCF3	TCF2	TCF1	TBL1XR1	TBL1X	T	SOX9	SALL4	RUNX2	RHOA	PTTG1
C	PTGS2	PPARD	POSTN	PLAUR	NRCAM	NOS2A	NANOG	MYCBP	MYC	MMP9	MMP7	MMP3
D	MMP2	MET	LRP6	LEF1	L1CAM	JUN	JAG1	ISL1	IL6	IGF2	IGF1	ID2
E	GJA1	GCG	FZD8	FZD7	FOXP1	FOSL1	FGF9	FGF4	FGF20	EPHB3	EPHB2	EGFR
F	EFNB1	EDN1	EDA	DLL1	DKK1	CTNNA1	CLDN1	CD44	BTRC	BMP4	BIRC5	BGLAP
G	AXIN2	AES	ATOH1	CDX1	CDX4	FST	NEUROG1	CDH1	ISLR	TNFSF11	GSN	IRX3
H	HGDC	HGDC	GAPDH	ACTB	B2M	RPL13A	HPRT1	RN18S1	RT	RT	PCR	PCR

Figure1. Illustration of QG060 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
QG060-01	A01	HQP021608	NM_003881	WISP2
QG060-01	A02	HQP021610	NM_003882	WISP1
QG060-01	A03	HQP018483	NM_005429	VEGFC
QG060-01	A04	HQP018475	NM_001025366	VEGFA
QG060-01	A05	HQP003299	NM_004385	VCAN
QG060-01	A06	HQP018328	NM_000474	TWIST1
QG060-01	A07	HQP018104	NM_005077	TLE1
QG060-01	A08	HQP018090	NM_003253	TIAM1
QG060-01	A09	HQP018047	NM_003238	TGFB2
QG060-01	A10	HQP017962	NM_030756	TCF7L2
QG060-01	A11	HQP020164	NM_031283	TCF7L1
QG060-01	A12	HQP017958	NM_003202	TCF7
QG060-01	B01	HQP017951	NM_003199	TCF4
QG060-01	B02	HQP017957	NM_003200	TCF3
QG060-01	B03	HQP017955	NM_000458	TCF2
QG060-01	B04	HQP017954	NM_000545	TCF1
QG060-01	B05	HQP019196	NM_024665	TBL1XR1
QG060-01	B06	HQP017927	NM_005647	TBL1X
QG060-01	B07	HQP017857	NM_003181	T
QG060-01	B08	HQP017635	NM_000346	SOX9
QG060-01	B09	HQP015461	NM_020436	SALL4
QG060-01	B10	HQP016478	NM_001015051	RUNX2
QG060-01	B11	HQP015997	NM_021205	RHOA
QG060-01	B12	HQP022362	NM_004219	PTTG1
QG060-01	C01	HQP015598	NM_000963	PTGS2
QG060-01	C02	HQP013627	NM_006238	PPARD
QG060-01	C03	HQP000764	NM_006475	POSTN
QG060-01	C04	HQP013205	NM_001005376	PLAUR
QG060-01	C05	HQP011915	NM_001037132	NRCAM
QG060-01	C06	HQP011866	NM_000625	NOS2A
QG060-01	C07	HQP019390	NM_024865	NANOG
QG060-01	C08	HQP007208	NM_012333	MYCBP
QG060-01	C09	HQP011597	NM_002467	MYC
QG060-01	C10	HQP011263	NM_004994	MMP9
QG060-01	C11	HQP011258	NM_002423	MMP7
QG060-01	C12	HQP011257	NM_002422	MMP3
QG060-01	D01	HQP011256	NM_004530	MMP2
QG060-01	D02	HQP011181	NM_000245	MET
QG060-01	D03	HQP010877	NM_002336	LRP6
QG060-01	D04	HQP012480	NM_016269	LEF1
QG060-01	D05	HQP010390	NM_000425	L1CAM
QG060-01	D06	HQP009853	NM_002228	JUN

QG060-01	D07	HQP004470	NM_000214	JAG1
QG060-01	D08	HQP009791	NM_002202	ISL1
QG060-01	D09	HQP009670	NM_000600	IL6
QG060-01	D10	HQP009529	NM_000612	IGF2
QG060-01	D11	HQP009518	NM_000618	IGF1
QG060-01	D12	HQP009273	NM_002166	ID2
QG060-01	E01	HQP007411	NM_000165	GJA1
QG060-01	E02	HQP007233	NM_002054	GCG
QG060-01	E03	HQP020140	NM_031866	FZD8
QG060-01	E04	HQP020136	NM_003507	FZD7
QG060-01	E05	HQP020808	NM_003593	FOXN1
QG060-01	E06	HQP019708	NM_005438	FOSL1
QG060-01	E07	HQP005416	NM_002010	FGF9
QG060-01	E08	HQP005405	NM_002007	FGF4
QG060-01	E09	HQP007198	NM_019851	FGF20
QG060-01	E10	HQP004942	NM_004443	EPHB3
QG060-01	E11	HQP004939	NM_004442	EPHB2
QG060-01	E12	HQP004605	NM_005228	EGFR
QG060-01	F01	HQP004596	NM_004429	EFNB1
QG060-01	F02	HQP004557	NM_001955	EDN1
QG060-01	F03	HQP004544	NM_001005609	EDA
QG060-01	F04	HQP008111	NM_005618	DLL1
QG060-01	F05	HQP005612	NM_012242	DKK1
QG060-01	F06	HQP003539	NM_001904	CTNNB1
QG060-01	F07	HQP022026	NM_021101	CLDN1
QG060-01	F08	HQP022972	NM_000610	CD44
QG060-01	F09	HQP021752	NM_003939	BTRC
QG060-01	F10	HQP017396	NM_001202	BMP4
QG060-01	F11	HQP009099	NM_001012270	BIRC5
QG060-01	F12	HQP016599	NM_199173	BGLAP
QG060-01	G01	HQP020116	NM_004655	AXIN2
QG060-01	G02	HQP004183	NM_001130	AES
QG060-01	G03	HQP011762	NM_005172	ATOH1
QG060-01	G04	HQP000534	NM_001804	CDX1
QG060-01	G05	HQP000567	NM_005193	CDX4
QG060-01	G06	HQP000564	NM_006350	FST
QG060-01	G07	HQP011773	NM_006161	NEUROG1
QG060-01	G08	HQP023466	NM_004360	CDH1
QG060-01	G09	HQP009792	NM_005545	ISLR
QG060-01	G10	HQP021321	NM_003701	TNFSF11
QG060-01	G11	HQP008470	NM_000177	GSN
QG060-01	G12	HQP018993	NM_024336	IRX3
QG060-01	H01	HGDC		
QG060-01	H02	HGDC		
QG060-01	H03	HQP006940	NM_002046	GAPDH
QG060-01	H04	HQP016381	NM_001101	ACTB

QG060-01	H05	HQP015171	NM_004048	B2M
QG060-01	H06	HQP006171	NM_012423	RPL13A
QG060-01	H07	HQP009026	NM_000194	HPRT1
QG060-01	H08	HQP054253	NR_003286	RN18S1
QG060-01	H09	RT		
QG060-01	H10	RT		
QG060-01	H11	PCR		
QG060-01	H12	PCR		

Limited Use License

Following terms and conditions apply to use of ExProfile™ Human Targets of Wnt/β-catenin 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