

miProfile[™] Mouse Lung Cancer miRNA qPCR Array

For focused group profiling of mouse lung cancer related miRNA expression

Cat. No. QM023-A (1 x 96-well plate, Format A) Cat. No. QM023-B (1 x 96-well plate, Format B)

Cat. No. QM023-C (1 x 96-well plate, Format C)

Cat. No. QM023-D (1 x 96-well plate, Format D)

Cat. No. QM023-E (1 x 96-well plate, Format E)

Available as 1 set or 6 sets. Each set contains 84 unique miRNA primers deposited in one 96-well plate.

Introduction

The miProfile mouse lung cancer miRNA qPCR array profiles 84 mouse miRNAs homologous to aberrantly expressed human miRNAs involved in lung cancer. These miRNAs are carefully chosen for their close cancer correlation based on a thorough literature search of peer-reviewed publications. miRNAs are known for their extensive involvement in cancer development and progression and their altered pattern of expression may affect the cell cycle and survival program. The miPofile mouse lung cancer array allows researchers to study the lung cancer-related miRNAs to gain understanding of the role of miRNAs in lung cancer pathogenesis.

QM023 plate 01: 84 unique miRNA PCR primer pairs

Shipping and storage condition

Shipped at room temperate Stable for at least 6 months when stored at -20 ℃

Array format

GeneCopoeia 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

- Each miRNA-specific primer in the miProfile miRNA 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 miRNA.
- 2. The positive PCR controls (PCR) have been verified to amplify a single amplicon of the correct size with Ct values around 20±2.
- 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. R₂ > 0.99 was observed for high inter/ intra-array reproducibility.

Materials required but not provided

All-in-OneTM miRNA First-Strand cDNA Synthesis Kit
All-in-OneTM 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	1	2	3	4	5	6	7	8	9	10	11	12
В	13	14	15	16	17	18	19	20	21	22	23	24
С	25	26	27	28	29	30	31	32	33	34	35	36
D	37	38	39	40	41	42	43	44	45	46	47	48
Е	49	50	51	52	53	54	55	56	57	58	59	60
F	61	62	63	64	65	66	67	68	69	70	71	72
G	73	74	75	76	77	78	79	80	81	82	83	84
Н	NC	NC	HK1	HK2	HK3	HK4	HK5	HK6	RT	RT	PCR	PCR

Figure 1. Illustration of miProfile miRNA qPCR array (96-well plate)

- miRNA primer pairs: Wells 1-84 are designated wells for pre-deposited miRNA primer pairs.
- NC: Negative controls, which only have the pre-deposited reverse universal primers.
- **HK1-6:** Six pre-deposited housekeeping snRNAs primer pairs, which can be used as endogenous positive controls as well as for array normalization.
- RT: Three replicates of spike-in reverse transcription controls, which can be used to monitor the
 efficiency of the RT reaction. These pre-deposited primer pairs specifically amplify the cDNA
 template reversed transcribed from the spike-in exogenous RNA in the sample.
- **PCR:** Three replicates of 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.

miRNA primer list

The primer list can be downloaded from http://www.genecopoeia.com/product/qpcr-arrays/mirna/cancer.php.

Limited Use License

Following terms and conditions apply to use of miProfileTM Mouse Lung Cancer miRNA 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