



## OmicsArray™ Antigen Microarray Processing Kit

Catalog No. PA100, PA101, PA102, PA103

### User Manual

GeneCopoeia, Inc.  
9620 Medical Center Drive, #101  
Rockville, MD 20850  
USA

301-762-0888

[inquiry@genecopoeia.com](mailto:inquiry@genecopoeia.com)

[genecopoeia.com](http://genecopoeia.com)

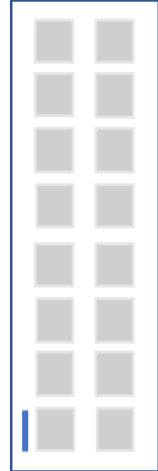
## I. Materials provided by GeneCopoeia

### 1. OmicsArray™ Antigen Microarray slide (sold separately)

The slides should be stored at 4°C upon receipt and are stable for up to 6 months.

### 2. OmicsArray™ Slide Processing kit

Front view of an  
OmicsArray™ slide



Product name	Contents	Cat. No.	Quantity	Storage temperature/ conditions
<b>OmicsArray™ antigen microarray processing kit (Human/Mouse) (PA100/PA101)</b>	<b>Blocking buffer I</b>	PA100-01	10 ml	4°C (Stable for 12months)
	<b>Blocking buffer II</b>	PA100-02	10 ml	4°C (Stable for 12months)
	<b>Wash buffer I (10x PBST)</b>	PA100-03	20 ml	4°C (Stable for 12months)
	<b>Wash buffer II (10x PBS)</b>	PA100-04	20 ml	4°C (Stable for 12months)
	<b>DNase I</b>	PA100-05	35 µl	-20°C (Stable for 12months)
	<b>DNase I buffer (10x)</b>	PA100-06	150 µl	-20°C (Stable for 12months)
	<b>0.1M DTT</b>	PA100-07	150 ul	-20°C (Stable for 12months)
	<b>anti-hlgG Cy3 anti-hlgM Cy5</b>	PA100-08 <sup>*1</sup> PA100-09 <sup>*1</sup>	10 µl	-20°C (Stable for 12months)
	<b>anti-mlgG Cy3 anti-mlgM Cy5</b>	PA101-08 <sup>2</sup> PA101-09 <sup>2</sup>	10 µl	-20°C (Stable for 12months)
<b>OmicsArray™ 4-slide Tray and gasket set (PA102)</b>	<b>4-slide tray and gasket set</b>	PA102	1 set	Room temperature
<b>OmicsArray™ 16-well slide gasket (PA103)</b>	<b>OmicsArray™ 16-Well Slide gasket</b>	PA103	2	Room temperature

<sup>1</sup>Only supplied in PA100; <sup>2</sup>Only supplied in PA101;

## II. User-Supplied Equipment

Equipment	Supplier
Orbitron Rotator II, Model 260250	Boekel Scientific
Centrifuge with swinging bucket rotor	Beckman Coulter or other vendor
Vacuum filtering flask connected with vacuum pump or vacuum system	Various

## III. Array Processing

### 1. Slide Preparation

- Take the slides out from 4°C and warm up to room temperature (RT) for 15 min.
- Mark slide number on the back of the slide (optional).
- Assemble the slide with the 16-well array gasket module by following the instructions (16-well array gasket module assembly). Make sure the array side is facing up (the line mark should be on the bottom left side as show in the figure).
- Add 100 µl wash buffer I to each well
- Wash with shaking (300rpm) for 5 min
- Remove the washing buffer with a pipette tip connected to a vacuum system.
- Blocking slides: Add 100 µl blocking buffer I into each well, incubate at RT for 1 hour using the orbital shaker (300 rpm).

### 2. Sample preparation

- Pre-treat serum or plasma samples with DNase I for 30min at RT. For samples that are not serum or plasma, DNase I treatment is not necessary. You can dilute those samples directly in 1x wash buffer I and proceed to Step 3 (Sample hybridization with arrays).
- Prepare DNase I treatment mixture:

	1 slide (15 samples)	2 slides (30 samples)	3 slides (45 samples)	4 slides (60 samples)
Nuclease free water	112.5 µl	225 µl	337.5 µl	450 µl
10x DNase I buffer	15 µl	30 µl	45 µl	60 µl
0.1M DTT	15 µl	30 µl	45 µl	60 µl
DNase I Enzyme	7.5 µl	15 µl	22.5 µl	30 µl

- Add 8ul of the above mixture into 2 µl each serum or plasma sample in a tube. Vortex briefly, and then spin down briefly.
- Incubate at RT for 30 min with slow shaking (120 rpm).
- Add 190 µl 1x wash buffer I into the tube of each serum or plasma mixture (total volume = 200 µl) and mix well. Now the samples are ready to be added onto the arrays.

### 3. Sample hybridization with arrays

- a) Remove blocking buffer I from each well of the array slide using a pipette tip connected to the vacuum filtering flask.
- b) Add 100  $\mu$ l wash buffer I to each well using a multi-channel pipette.
- c) Shake for 5 min.
- d) Repeat the washing step two times.
- e) Remove the washing buffer from each well of the slide using a vacuum
- f) Transfer 100  $\mu$ l diluted or treated sample into each well.
- g) Cover the array plates and incubate the arrays at RT for 1 hour with shaking (300rpm).

### 4. Washing

- a) Remove the samples from the wells of the array slide using a vacuum.
- b) Add 100  $\mu$ l wash buffer I to each well. Incubate 5 min. at RT with shaking (300rpm).
- c) Remove wash buffer I. Add 100  $\mu$ l blocking buffer I and shake for 5 min at RT.
- d) Remove blocking buffer II, then add 100ul wash buffer I to each well and shake for 5 min.
- e) Repeat the washing steps two times.

### 5. Secondary antibody processing

- a) Prepare the dilution mixture of secondary antibodies (1:1,000 dilutions of Cy3 anti-human or anti-mouse IgG and Cy5 anti-human or anti-mouse) IgM using the following table.

	1 slide	2 slides	3 slides	4 slides
1x wash buffer I	2 ml	4 ml	6 ml	8 ml
Cy3 anti-human IgG	2 $\mu$ l	4 $\mu$ l	6 $\mu$ l	8 $\mu$ l
Cy5 anti-human IgM	2 $\mu$ l	4 $\mu$ l	6 $\mu$ l	8 $\mu$ l

- b) Wrap the tube with foil to protect from light. Then mix and spin.
- c) Remove wash buffer from each well of the slide. Add 100  $\mu$ l freshly prepared secondary antibody dilution to each well.
- d) Incubate the slide with secondary antibodies for 1 hour at RT with shaking (300rpm). The plate should be incubated in the dark to prevent photobleaching.

### 5. Slide Washing

- a) Remove the secondary antibody mixture from each well. Add 100  $\mu$ l wash buffer I to each well and wash for 5 min. at RT with shaking (300rpm).
- b) Repeat the washing step two more times, for a total of three washes.

- c) Remove wash buffer I.
- d) Disassemble the gasket frame from the slide.
- e) Place the slide into a 50 ml conical tube containing 45 ml 1x wash buffer II (PBS buffer).
- f) Wash for 5 min. at RT with slow agitation.
- g) Transfer the slide to a new conical tube containing 45 ml of ddH<sub>2</sub>O.
- h) Rinse slide up and down at RT for 5 min.

Note: After removing buffer from each well, add new buffer immediately to prevent drying of the array between each step. Drying of the array during processing may generate high background.

## 6. Slide Scanning

- a) Transfer the slide to an empty conical tube with no cap. Centrifuge using a swinging bucket rotor at 500 rpm for 5 min. The slide should be dry after centrifugation.
- b) Place the slide into a GenePix 4000B or equivalent two-channel laser scanner. Make sure the slide is facing down (the line mark should be located in the bottom right in the slide holder of the scanner).
- c) Scan Settings: Please follow the manual of the manufacturer for scanner settings.

The following are the suggested settings used on the Genepix 4000B scanner

- For the Cy3 channel, use 532nm, PMT 280-400, and power 33.
  - For the Cy-5 channel, use 635nm, PMT 400-460 and power 33.
- d) Save the scanned images at both 532nm and 635nm.
  - e) Use GenePix (or equivalent) software to analyze the images. Apply the \*.gal file to the image and generate a GPR (GenePix Report) file.

## **IV. Limited Use License and Warranty**

### **Limited Use License**

The following terms and conditions apply to use of all OmicsArray™ antigen microarray processing kits (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 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.