

Antigen Protein Microarrays

Full length antigen proteins

GeneCopoeia's OmicsArray™ antigen protein microarrays are engineered for several powerful applications, including autoantibody detection, biomarker profiling, and characterization of protein-molecule interactions. Antigen protein microarrays enable parallel detection of antibodies or other biomarkers present in patient samples for autoimmune diseases including rheumatoid arthritis, muscular dystrophy, systemic lupus erythematosus, and type 1 diabetes as well as pathogen antibodies.

Each array contains superior-quality purified proteins spotted onto nitrocellulose filters, which are adhered to glass slides. Proteins known to be associated with specific diseases are chosen based on a thorough review of peer-reviewed publications.

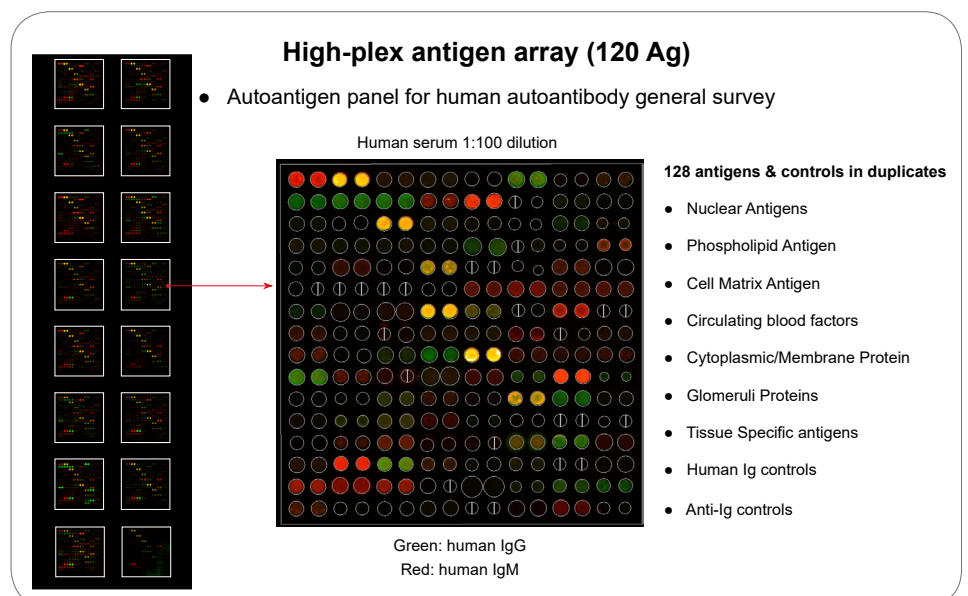
In addition to predesigned arrays, arrays containing customized sets of proteins are available, as well as array profiling services and data analysis.

Advantages

- **Multiplexing.** OmicsArray™ antigen protein microarrays can assay up to 120 antigens per well, compared with 1 protein per well for ELISA.
- **High throughput.** Each slide can process up to 15 samples in parallel.
- **High sensitivity.** Each array can detect as little as 1 pg/ml of antibody, which is 100-fold more sensitive than ELISA.
- **Full length antigen proteins.**

Antibody Profiling:

- Autoimmune diseases
- Neurological disorders
- Cancer immunotherapy
- Pathogens & allergens
- Coronavirus variants



Antigen Protein Microarrays

Pre-designed antigen protein microarrays

Catalog Number	Number of Antigens	Description
PA001	120	Human autoantigens general survey
PA002	120	Human brain and central nervous system disorders
PA003	120	Human cancer and neoplasms
PA006	75	Human common allergens
PA010	120	Human autoimmunity, allergy, and infection
PA012	120	Human coronavirus-associated autoimmunity (CAA)
PA013	58	NEW! SARS-CoV-2 coronavirus variant proteins

Custom antigen protein microarray services

- **Sample processing.** Send us your blood, plasma, tissue, or other biological sample and we will prepare it for processing and incubation with any of our predesigned antigen microarrays or custom-built antigen microarrays
- **Data analysis.** Once samples are processed and incubated with an antigen microarray, we will analyze the raw data. The standard analysis service includes: 1) An Excel file of the Net Signal Intensity (NSI) for each antigen on the array, normalized to internal controls; and 2) a heat map.

Additional analysis services, including proteomic analysis, pathway analysis, and more, are also available.

To inquire about custom antigen microarray services, please contact us at inquiry@genecopoeia.com.

Contact us:

GeneCopoeia, Inc.

9620 Medical Center Dr. Suite 101,
Rockville, MD 20850 USA



Email

Tel

Fax

Website

inquiry@genecopoeia.com

+1 (301) 762-0888

+1 (301) 762-3888

www.genecopoeia.com