



Selected publications citing targets on OmicsArray™ antigen microarrays for autoimmune & other diseases

Systemic Lupus Erythematosus		
Citation	OmicsArray™ Antigen array	
• Martinez, J., et al. Noncanonical autophagy inhibits the autoinflammatory, lupus-like response to dying cells. <i>Nature</i> 2016; 533: 115. • Rajasinghe, L.D., et al. Omega-3 fatty acid intake suppresses induction of diverse autoantibody repertoire by crystalline silica in lupus-prone mice. <i>Autoimmunity</i> 2020; 53: 415. • Shiozawa, S., et al. DOCK8-expressing T follicular helper cells newly generated beyond self-organized criticality cause systemic lupus erythematosus. <i>iScience</i> 2021; 25: 103537.	Human autoantigens general survey (PA001)	
Sjögren's Syndrome		
Citation	OmicsArray™ Antigen array	
• Kramer, J.M., <i>et al.</i> Analysis of IgM antibody production and repertoire in a mouse model of Sjögren's syndrome. <i>J. Leukoc. Biol.</i> 2016; 99: 321.	Human autoantigens general survey (PA001)	
Rheumatoid Arthritis		
Citation	OmicsArray™ Antigen array	
• Huang, Q-q., <i>et al.</i> CD11c-mediated deletion of Flip promotes autoreactivity and inflammatory arthritis. <i>Nat. Commun.</i> 2015; 6: 7086.	Human autoantigens general survey (PA001)	
Chronic Obstructive Pulmonary Disease (COPD)		
Citation	OmicsArray™ Antigen array	
• Cass, S.P., et al. Differential expression of sputum and serum autoantibodies in patients with chronic obstructive pulmonary disease. <i>Am. J. Physiol. Lung Cell Mol. Physiol.</i> 2021; 320: L1169	Human autoantigens general survey (PA001)	
Allergy		
Citation	OmicsArray™ Antigen array	
• Turner, J.A., <i>et al.</i> Regulatory T Cell-Derived TGF-β1 Controls Multiple Checkpoints Governing Allergy and Autoimmunity. <i>Immunity</i> 2020; 53: 1202.	Human autoantigens general survey (PA001)	
Morphea		
Citation	OmicsArray™ Antigen array	
• Zhu, J.L., <i>et al.</i> Autoantigen microarrays reveal myelin basic protein autoantibodies in morphea. <i>J. Transl. Med.</i> 2022; 20: 41.	Human brain and central nervous system disorders (PA002)	

Hereditary alpha-tryptasemia		
Citation	OmicsArray™ Antigen array	
Konnikova, L., <i>et al.</i> Small intestinal immunopathology and GI-associated antibody formation in hereditary alpha-tryptasemia. <i>J. Allergy and Clinical Immunology</i> 2021; 148: 813.	Human autoantigens general survey (PA001)	
COVID-19		
Citation	OmicsArray™ Antigen array	
van Oers N.S.C., <i>et al.</i> SARS-CoV-2 infection associated with hepatitis in an infant with X-linked severe combined immunodeficiency. <i>Clin. Immunol.</i> 2021; 224: 108662.	SARS-CoV-2 coronavirus variant proteins (PA013)	
Muthukumar, A., <i>et al.</i> In-Depth Evaluation of a Case of Presumed Myocarditis After the Second Dose of COVID-19 mRNA Vaccine. <i>Circulation</i> 2021; 144: 487.		
Rojas, M., <i>et al.</i> Autoimmunity is a hallmark of post-COVID syndrome. <i>J. Transl. Med.</i> 2022; 20: 129.	Human Coronavirus-associated Autoimmunity (PA012)	
Cancer Immunotherapy		
Citation	OmicsArray™ Antigen array	
Khan, S., <i>et al.</i> Late-Onset Immunotherapy Toxicity and Delayed Autoantibody Changes: Checkpoint Inhibitor-Induced Raynaud's-Like Phenomenon. <i>Oncologist</i> 2020; 25: e753.	Human cancer and neoplasms (PA003)	
Ghosh, N., <i>et al.</i> Lower baseline autoantibody levels are associated with immune-related adverse events from immune checkpoint inhibition. <i>J. Immunother. Cancer</i> 2022; 10: e004008.		
Vaccine-related autoimmunity		
Citation	OmicsArray™ Antigen array	
Labombarde, J.G., <i>et al.</i> Induction of broadly reactive influenza antibodies increases susceptibility to autoimmunity. <i>Cell Rep.</i> 2022; 38: 110482.	Human autoantigens general survey (PA001)	
General autoimmunity		
Citation	OmicsArray™ Antigen array	
Mayer, C.T., <i>et al.</i> An apoptosis-dependent checkpoint for autoimmunity in memory B and plasma cells. <i>Proc. Natl. Acad. Sci. U.S.A.</i> 2020; 117: 24957.	Human autoantigens general survey (PA001)	
Rojas, M., <i>et al.</i> New insights into the taxonomy of autoimmune diseases based on polyautoimmunity. <i>J. Autoimm.</i> 2022; 126: 102780.		
Pichilingue-Reto, P., <i>et al.</i> Serum IgG Profiling of Toddlers Reveals a Subgroup with Elevated Seropositive Antibodies to Viruses Correlating with Increased Vaccine and Autoantigen Responses. <i>J. Clin. Immunol.</i> 2021; 41: 1031	Human Autoimmunity, Allergy, and Infection (PA010)	

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