



BioPlex™ 2200 Panel Menu

Autoimmune Disease Panels

ANA Screen with MDSS	
• ANA Screen***	• Sm
• dsDNA	• Sm/RNP
• Chromatin	• RNP (A and 68)***
• Ribosomal P	• SCL-70
• SS-A (52kD & 60kD)**	• Jo-1
• SS-B	• Centromere B

Vasculitis		
• MPO	• PR3	• GBM

Gastrointestinal IgG* and IgA*		
• Gliadin	• tTG	• ASCA

Anti-Phospholipid Syndrome IgG* and IgM*	
• Cardiolipin	• β2-GPI
• Phosphatidylserine	• Prothrombin

Rheumatoid Arthritis*	
• Rheumatoid Factor	• CCP

Other Panels in Development

• Cardiac	• Diabetes
• Toxicology	

* In development

** Individual results available outside US only

*** ANA Screen is a composite roll up of all individual results

Infectious Disease Panels

EBV IgG and IgM	
• VCA (IgG and IgM)	• EBNA (IgG)
• EA-D (IgG)	• Heterophile (IgM)

Syphilis IgG	
• <i>T. pallidum</i> 15kD**	• <i>T. pallidum</i> 47kD**
• <i>T. pallidum</i> 17kD**	

Syphilis IgM*	
• <i>T. pallidum</i> 17kD	• <i>T. pallidum</i> 47kD

ToRC IgG* and IgM*	
• <i>Toxoplasma gondii</i>	• Cytomegalovirus
• Rubella	

Immunity IgG*	
• Mumps	• Varicella-Zoster virus
• Measles	• Rubella

Lyme Borreliosis*	
• IgG	• IgM

HSV*	
• HSV-1 IgG	• HSV-1/2 IgM
• HSV-2 IgG	

