



BioPlex® 2200 Panel Menu

Autoimmune Disease Panels

ANA Screen with MDSS	
ANA Screen ¹	Sm
dsDNA	Sm/RNP
Chromatin	RNP (A and 68) ²
Ribosomal P	Scl-70
SS-A (52kD & 60kD) ²	Jo-1
SS-B	Centromere B

Vasculitis		
MPO	PR3	GBM

Celiac IgA ³	
Deamidated Gliadin Peptide	tTG

Celiac IgG ³	
Deamidated Gliadin Peptide	tTG

Antiphospholipid Syndrome (APLS) IgG ³	
Cardiolipin	β2-GPI

Antiphospholipid Syndrome (APLS) IgM ³	
Cardiolipin	β2-GPI

Antiphospholipid Syndrome (APLS) IgA ³	
Cardiolipin	β2-GPI

Anti-CCP	
CCP	

Other Panels in Development

HAV	HBV	Vitamin D
HBSAg	HCV	Lyme Borreliosis

¹ ANA Screen is a composite roll up of all individual results

² Individual results available outside US only

³ Available outside US only

⁴ Available in US only

⁵ In development

Infectious Disease Panels

EBV IgG	
VCA	EBNA
EA-D	

EBV IgM	
VCA	Heterophile

Syphilis IgG	
<i>T. pallidum</i> 15kD ²	<i>T. pallidum</i> 47kD ²
<i>T. pallidum</i> 17kD ²	

Syphilis IgM ³	
<i>T. pallidum</i> 17kD ²	<i>T. pallidum</i> 47kD ²

ToRC IgG	
<i>Toxoplasma gondii</i>	Cytomegalovirus
Rubella	

ToRC IgM ³	
<i>Toxoplasma gondii</i>	Cytomegalovirus
Rubella	

Rubella & CMV IgM ⁴	
Rubella	Cytomegalovirus

MMRV IgG	
Mumps	Varicella-Zoster virus
Measles	Rubella (Qualitative)

HSV-1 & HSV-2 IgG	
HSV-1	HSV-2

HIV Ag-Ab ⁵		
HIV-1 p24 Ag	HIV-1 Ab (Group M & O)	HIV-2 Ab

