



# BioPlex® 2200 Panel Menu

## Autoimmune Disease Panels

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

Vasculitis		
MPO	PR3	GBM

Celiac IgA <sup>3</sup>	
Deamidated Gliadin Peptide	tTG

Celiac IgG <sup>3</sup>	
Deamidated Gliadin Peptide	tTG

Antiphospholipid Syndrome (APLS) IgG <sup>3</sup>	
Cardiolipin	β2-GPI

Antiphospholipid Syndrome (APLS) IgM <sup>3</sup>	
Cardiolipin	β2-GPI

Antiphospholipid Syndrome (APLS) IgA <sup>3</sup>	
Cardiolipin	β2-GPI

Anti-CCP	
CCP	

## Other Panels in Development

HAV	HBV	Vitamin D
HBSAg	HCV	Lyme Borreliosis

<sup>1</sup> ANA Screen is a composite roll up of all individual results

<sup>2</sup> Individual results available outside US only

<sup>3</sup> Available outside US only

<sup>4</sup> Available in US only

<sup>5</sup> In development

## Infectious Disease Panels

EBV IgG	
VCA	EBNA
EA-D	

EBV IgM	
VCA	Heterophile

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

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

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

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

Rubella & CMV IgM <sup>4</sup>	
Rubella	Cytomegalovirus

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

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

HIV Ag-Ab <sup>5</sup>		
HIV-1 p24 Ag	HIV-1 Ab (Group M & O)	HIV-2 Ab

