



# 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	Sci-70
SS-A (52kD & 60kD) <sup>2</sup>	Jo-1
SS-B	Centromere B
Vasculitis	
MPO	PR3
GBM	
Celiac IgA	
tTG	DGP
Celiac IgG	
tTG	DGP
Antiphospholipid Syndrome (APLS) IgG	
Cardiolipin	β2-GPI
Antiphospholipid Syndrome (APLS) IgM	
Cardiolipin	β2-GPI
Antiphospholipid Syndrome (APLS) IgA	
Cardiolipin	β2-GPI
Anti-CCP	
CCP	

## Special Chemistry Panels

25-OH Vitamin D	
25-OH Vitamin D	

Other Panels in Development – Lyme Borreliosis

## Infectious Disease Panels

EBV IgG	
VCA	EBNA
EA-D	
EBV IgM	
VCA	Heterophile
Syphilis Total & RPR	
Syphilis Total	RPR Titration
RPR	
ToRC IgG	
<i>Toxoplasma</i>	CMV
Rubella (Quantitative)	
ToRC IgM	
<i>Toxoplasma</i>	CMV
Rubella	
MMRV IgG	
Mumps	VZV
Measles	Rubella (Qualitative)
HSV-1 & HSV-2 IgG	
HSV-1	HSV-2
HIV Ag-Ab <sup>†</sup>	
HIV Ag-Ab <sup>3</sup>	HIV-1 p24 Antigen
HIV-1 Antibody <sup>2</sup> (Group M and O)	HIV-2 Antibody

<sup>1</sup> Composite roll up of all individual results

<sup>2</sup> Composite roll up of individual results. Individual results not available in the US

<sup>3</sup> Overall result of individual assays

<sup>†</sup> All individual assay results reported

