



# PhD™ Validated Assay List

## Autoimmune Systemic EIA

Catalog Number	Kit Name
96AN	ANA Screen
96DS	Anti-dsDNA
96EP	ENA Plus Screen
96SA	Anti-SSA
96SB	Anti-SSB
96SM	Anti-Sm
96SR	Anti-SmRNP
96CE	Anti-Centromere
96JO	Anti-Jo-1
96SC	Anti-Scl-70
96SS	Anti-ssDNA
96CP	Anti-CCP

## Antiphospholipid Syndrome (APS) EIA

Catalog Number	Kit Name
425-2000	Anti-Cardiolipin IgG
425-2020	Anti-Cardiolipin IgM
425-2040	Anti-Cardiolipin IgA
425-2060	Anti-Beta 2 Glycoprotein I ( $\beta$ 2GPI) IgG
425-2080	Anti-Beta 2 Glycoprotein I ( $\beta$ 2GPI) IgM
425-2100	Anti-Beta 2 Glycoprotein I ( $\beta$ 2GPI) IgA
425-2120	Anti-Phosphatidylserine (aPS) IgG
425-2140	Anti-Phosphatidylserine (aPS) IgM
425-2160	Anti-Phosphatidylserine (aPS) IgA
425-2180	Anti-Prothrombin IgG
425-2200	Anti-Prothrombin IgM

## Gastrointestinal (GI) EIA

Catalog Number	Kit Name
425-2260	Anti-tissue Transglutaminase (tTG) IgA
425-2280	Anti-tissue Transglutaminase (tTG) IgG
425-2220	Anti-Gliadin IgA
425-2240	Anti-Gliadin IgG
425-2320	Anti-Saccharomyces cerevisiae (ASCA) IgA
425-2340	Anti-Saccharomyces cerevisiae (ASCA) IgG

## Infectious Disease EIA (US Only)

Catalog Number	Kit Name
32507	Anti-Borrelia (Lyme)
25177	CMV IgG
25178	CMV IgM
25184	EB VCA IgG
25185	EB VCA IgM
25183	EB NA IgG
25225	<i>Helicobacter pylori</i> ( <i>H. pylori</i> ) IgA
25226	<i>Helicobacter pylori</i> ( <i>H. pylori</i> ) IgG
25179	HSV IgG
25181	Measles IgG
25182	Mumps IgG
25173	Rubella IgG
25174	Rubella IgM
25175	Toxoplasma IgG
25176	Toxoplasma IgM
25180	VZV IgG

**PhD™ Validated Assay List**

**Autoimmune Systemic IFA**

Catalog Number	Kit Name
32583	Kallestad™ HEp-2 Kit (8x6w slides)
30471	Kallestad™ HEp-2 Kit (5x12w slides)
30472	Kallestad™ HEp-2 Kit (20x12w slides)
30404	Kallestad™ <i>Crithidia luciliae</i> (6x4w slides)
31069	Kallestad™ <i>Crithidia luciliae</i> (6x8w slides)

**ANCA (Anti-Neutrophil Cytoplasmic Antibody) IFA**

Catalog Number	Kit Name
29403	Kallestad™ ANCA IFA Kit (10x6w ethanol fixed slides), 60 Tests
29417	Kallestad™ ANCA (5x12w ethanol fixed slides)
29418	Kallestad™ ANCA (5x6w formalin fixed slides)
29419	Kallestad™ ANCA COMBI Slides (5x12w ethanol/formalin fixed slides, 6 wells each)

**Organ-Specific IFA**

Catalog Number	Kit Name
30443	Kallestad™ Mouse Stomach/Kidney (6x8w slides)



**Bio-Rad  
Laboratories**

*For further information, please contact the Bio-Rad office nearest you or visit our website at [www.bio-rad.com/autoimmune](http://www.bio-rad.com/autoimmune)*

Clinical  
Diagnostics Group

**Website** [www.bio-rad.com/diagnostics](http://www.bio-rad.com/diagnostics) **U.S.** 1-800-2BIO-RAD **Australia** 61-2-9914-2800 **Austria** 43-1-877-8901 **Belgium** 32-9-385-5511 **Brazil** 5521-3237-9400 **Canada** 1-514-334-4372 **China** 86-21-64260808 **Czech Republic** 420-241-430-532 **Denmark** +45-4452-1000 **Finland** 358-9-804-22-00 **France** 33-1-47-95-60-00 **Germany** +49-(0)89-318-840 **Greece** 30-210-7774396 **Hong Kong** 852-2789-3300 **Hungary** +36-1-459-6100 **India** 1-800-180-1224 **Israel** 972-3-9636050 **Italy** +39-02-216091 **Japan** 81-3-6361-7070 **Korea** 82-2-3473-4460 **Mexico** +52 (55) 5488-7670 **The Netherlands** +31-318-540666 **New Zealand** 64-9-415-2280 **Norway** 47-23-38-41-30 **Poland** 48-22-3319999 **Portugal** 351-21-472-7700 **Russia** 7-495-721-14-04 **Singapore** 65-6415-3170 **South Africa** 27-11-442-85-08 **Spain** 34-91-590-5200 **Sweden** 46-8-555-127-00 **Switzerland** 41-61-717-95-55 **Thailand** 662-651-8311 **United Kingdom** +44-(0)20-8328-2000