



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 (β 2GPI) IgG
425-2080	Anti-Beta 2 Glycoprotein I (β 2GPI) IgM
425-2100	Anti-Beta 2 Glycoprotein I (β 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

*Clinical
Diagnostics Group*

Website 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

Printed on recycled paper with soy-based inks

© 2010 Bio-Rad Laboratories, Inc. Printed in USA 06/10 DG10-0889 Q-784