



PhD™ Ix Method File 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 II

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

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- <i>Saccharomyces cerevisiae</i> (ASCA) IgA
425-2340	Anti- <i>Saccharomyces cerevisiae</i> (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
25180	VZV IgG

Vasculitis EIA

Catalog Number	Kit Name
425-2420	Anti-PR3 (cANCA)
425-2380	Anti-MPO (pANCA)
425-2400	Anti-GBM*

*Not available in the US

Autoimmune Systemic IFA

Catalog Number	Kit Name	
32583	Kallestad® Hep-2 Kit (48 tests)	8 x 6 wells
30471	Kallestad® Hep-2 Kit (60 tests)	5 x 12 wells
30472	Kallestad® Hep-2 Kit (240 tests)	20 x 12 wells
26100	Kallestad® Hep-2 Slides	8 x 6 wells
26101	Kallestad® Hep-2 Slides	20 x 6 wells
26102	Kallestad® Hep-2 Slides	5 x 12 wells
26103	Kallestad® Hep-2 Slides	20 x 12 wells
26104	Kallestad® Hep-2 Slides	100 x 12 wells
30404	Kallestad® <i>Crithidia luciliae</i> Kit (24 tests)	6 x 4 wells
31069	Kallestad® <i>Crithidia luciliae</i> Kit (48 tests)	6 x 8 wells
26107	Kallestad® <i>Crithidia luciliae</i> Slides	6 x 4 wells
26108	Kallestad® <i>Crithidia luciliae</i> Slides	20 x 4 wells
26109	Kallestad® <i>Crithidia luciliae</i> Slides	6 x 8 wells
26110	Kallestad® <i>Crithidia luciliae</i> Slides	20 x 8 wells

ANCA (Anti-Neutrophil Cytoplasmic Antibody) IFA

Catalog Number	Kit Name	
29403	ANCA Ethanol Kit (60 tests)	10 x 6 wells
29416	ANCA Slide Ethanol Fixation	5 x 6 wells
29417	ANCA Slide Ethanol Fixation	5 x 12 wells
29418	ANCA Slide Formalin Fixation	5 x 6 wells
29419	ANCA Combi Slide (6 ethanol + 6 formalin wells)	5 x 12 wells

Organ-Specific IFA

Catalog Number	Kit Name	
30443	Kallestad® MSK Kit (48 tests)	6 x 8 wells
26111	Kallestad® MSK Slides	6 x 8 wells
26115	Kallestad® MSK Slides	20 x 8 wells
26112	Kallestad® MSK Slides	100 x 8 wells



**Bio-Rad
Laboratories**

For further information, please contact the Bio-Rad office nearest you or visit our website at www.bio-rad.com/diagnostics

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics U.S. 1-800-224-6723 Australia 61-2-9914-2800 Austria 43-1-877-8901 Belgium +32 (3)710-53-00 Brazil +55 (31)3689-6600 Canada 1-514-334-4372 China 86-21-61698500 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 1800-180-1224 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 +7-495-721-1404 Singapore 65-6415-3170 South Africa 27-11-442-85-08 Spain 34-91-590-5200 Sweden +46-8-555-127-00 Switzerland +41 (0)26-674-55-05/06 Taiwan 886-2-2578-7189 Thailand 662-651-8311 United Kingdom +44 (0)20-8328-2000