



PhD Ix Method List

Autoimmune Systemic IFA (Immunofluorescence Assay)

Catalog Number	Kit Name	Wells
32583	Kallestad™ HEP-2 Kit — 48 tests	8 x 6
30471	Kallestad HEP-2 Kit — 60 tests	5 x 12
30472	Kallestad HEP-2 Kit — 240 tests	20 x 12
26101	Kallestad HEP-2 Slides	20 x 6
26102	Kallestad HEP-2 Slides	5 x 12
26103	Kallestad HEP-2 Slides	20 x 12
26104	Kallestad HEP-2 Slides	100 x 12
30404	Kallestad <i>Crithidia luciliae</i> Kit — 24 tests	6 x 4
31069	Kallestad <i>Crithidia luciliae</i> Kit — 48 tests	6 x 8
26108	Kallestad <i>Crithidia luciliae</i> Slides	20 x 4
26109	Kallestad <i>Crithidia luciliae</i> Slides	6 x 8
26110	Kallestad <i>Crithidia luciliae</i> Slides	20 x 8

ANCA (Anti-Neutrophil Cytoplasmic Antibody) IFA

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

Organ-Specific IFA

Catalog Number	Kit Name	Wells
30443	Kallestad Mouse Stomach/Kidney (MSK) Kit — 48 tests	6 x 8
26111	Kallestad MSK Slides	6 x 8
26115	Kallestad MSK Slides	20 x 8
26112	Kallestad MSK Slides	100 x 8
29020*	Kallestad Rat Liver/Kidney with Medulla/Stomach Slides — 96 tests	12 x 8
29021*	Kallestad Rat Liver/Kidney with Medulla/Stomach Slides — 48 tests	12 x 4
29003*	Kallestad Monkey Endomysium — 48 tests	12 x 4
29060*	Kallestad Monkey Endomysium — 96 tests	12 x 8

*Not available in the U.S.

Autoimmune Systemic EIA (Enzyme Immunoassay)

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

Infectious Disease EIA (U.S. 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

Bio-Rad and Kallestad are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner.



Bio-Rad
Laboratories, Inc.

Clinical
Diagnostics Group

Website www.bio-rad.com/diagnostics **U.S.** 1 800 224 6723 **Australia** +61 (2) 9914 2800 **Austria** +43 (0) 1 877 89 01 9 **Belgium** +32 (0) 3 710 53 00 **Brazil** +55 11 3065 7550 **Canada** +1 514 334 4372 **China** +86 21 6169 8500 **Czech Republic** +420 241 431 660 **Denmark** +45 44 52 10 00 **Finland** +358 9 804 22 00 **France** +33 (0) 1 47 95 60 00 **Germany** +49 (0) 89 31884 393 **Greece** +30 210 7774396 **Hong Kong** +85 2 2789 3300 **Hungary** +36 1 459 6190 **India** +91 124 4029300 **Israel** +972 03 963 6025 **Italy** +39 024 94 86 600 **Japan** +81 3 6361 7070 **Mexico** +52 (55) 5488 7670 **The Netherlands** +31 (0) 318 540 666 **New Zealand** +64 (9)415 2280 **Norway** +47 23 38 41 30 **Poland** +48 22 331 99 99 **Portugal** +351 21 47 27 700 **Republic of Korea** +82 080 007 7373 **Russia** +7 (495) 721-14-04 **Singapore** +65 6415 3170 **South Africa** +27 11 442 8508 **Spain** +34 91 490 6580 **Sweden** +46 844 98053 **Switzerland** +41 (0) 61 717 9555 **Taiwan** +886 (2) 2578-7189 **Thailand** (662) 651 8311 **United Kingdom** +44 (0) 1923 471301