



List of EIA Microplate Assays

Blood Virus Assays

HIV	HAV	HBV	HCV
<ul style="list-style-type: none"> • HIV-1 • HIV-1/HIV-2 +O • HIV-2 	<ul style="list-style-type: none"> • anti-HAV • anti-HAV IgM 	<ul style="list-style-type: none"> • HBsAg • HBsAg confirmation • anti-HBs quantitative • anti-HBs qualitative • anti-HBc • anti-HBc IgM • HBeAg • Anti-HBe 	<ul style="list-style-type: none"> • HCV

Autoimmune Assays

Systemic	Anti-Phospholipid
<ul style="list-style-type: none"> • ANA Screen • ENA Plus Screen • ENA Profile • Anti-dsDNA • Anti-ssDNA • Anti-Jo-1 • Anti-SS-A • Anti-SS-B • Anti-Scl-70 • Anti-Sm • Anti-Sm/RNP 	<ul style="list-style-type: none"> • Anti-Centromere • Anti-CCP • Anti-GBM • Anti-MPO pANCA • Anti-PR3 cANCA • Anti-Tg • Anti-TPO • Anti-Mitochondrial • Anti-RO (SS-A) • Anti-La (SS-B)
	<ul style="list-style-type: none"> • aPS IgG • aPS IgM • aPS IgA • aCL IgG • aCL IgM • aCL IgA • aβ2GPI IgG • aβ2GPI IgM • aβ2GPI IgA

Infectious Disease Assays

Qualitative	Semi-Qualitative	Quantitative
<ul style="list-style-type: none"> • Rubella IgG • Rubella IgM • Toxoplasma IgG • Toxoplasma IgM • CMV IgG • CMV IgM • HSV IgG • VZV IgG • Measles IgG • Mumps IgG • EB NA IgG • EB VCA IgG • EB VCA IgM • Syphilis IgG • Pylori Detect IgG • Pylori Detect IgA • Chlamydia Ag • Anti-Borrelia (Lyme) • Aspergillus 	<ul style="list-style-type: none"> • Rubella IgG • Toxoplasma IgG • CMV IgG • HSV IgG • VZV IgG • Measles IgG • Mumps IgG • EB NA IgG • EB VCA IgG 	<ul style="list-style-type: none"> • Rubella IgG • Toxoplasma IgG • Toxoplasma IgM

**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. 1-800-2BIO-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 36-1-455-8800 India 91-124-4029300 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 65-6415-3188 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