



# List of EIA Microplate Assays

## Blood Virus Assays

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

## Autoimmune Assays

Systemic	Anti-Phospholipid
<ul style="list-style-type: none"> <li>• ANA Screen</li> <li>• ENA Plus Screen</li> <li>• ENA Profile</li> <li>• Anti-dsDNA</li> <li>• Anti-ssDNA</li> <li>• Anti-Jo-1</li> <li>• Anti-SS-A</li> <li>• Anti-SS-B</li> <li>• Anti-Scl-70</li> <li>• Anti-Sm</li> <li>• Anti-Sm/RNP</li> </ul>	<ul style="list-style-type: none"> <li>• Anti-Centromere</li> <li>• Anti-CCP</li> <li>• Anti-GBM</li> <li>• Anti-MPO pANCA</li> <li>• Anti-PR3 cANCA</li> <li>• Anti-Tg</li> <li>• Anti-TPO</li> <li>• Anti-Mitochondrial</li> <li>• Anti-RO (SS-A)</li> <li>• Anti-La (SS-B)</li> </ul>
	<ul style="list-style-type: none"> <li>• aPS IgG</li> <li>• aPS IgM</li> <li>• aPS IgA</li> <li>• aCL IgG</li> <li>• aCL IgM</li> <li>• aCL IgA</li> <li>• a<math>\beta</math>2GPI IgG</li> <li>• a<math>\beta</math>2GPI IgM</li> <li>• a<math>\beta</math>2GPI IgA</li> </ul>

## Infectious Disease Assays

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

**Bio-Rad  
Laboratories**

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

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 36-1-455-8800 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