



List of EIA Assays

Blood Virus Menu for HIV and Hepatitis

HIV	HAV	HBV	HCV
<ul style="list-style-type: none"> • HIV-1 • HIV-1/HIV-2 • HIV-2 	<ul style="list-style-type: none"> • HAV Total • HAV IgM 	<ul style="list-style-type: none"> • HBsAg • HBsAg confirmation • HBsAb quantitative • HBsAb qualitative • HBc Total • HBc IgM • HBeAg • HBeAb 	<ul style="list-style-type: none"> • HCV

List of EIA Assays

Autoimmune Menu

Systemic Tests	Anti-Phospholipid Tests
<ul style="list-style-type: none"> • ANA Screen • ENA Plus Screen • Anti-dsDNA • Anti-Jo-1 • Anti-SS-A • Anti-SS-B • Anti-Scl-70 • Anti-Sm • Anti-Sm/RNP • Anti-Centromere • Anti-GBM • Anti-mPOp ANCA • Anti-PR3c ANCA • Anti-Tg • Anti-TPO • Anti-Mitochondrial 	<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 Menu

Qualitative	Semi-Quantitative	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 	<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** 81-3-5811-6290 **Korea** 82-2-3473-4460 **Mexico** +52-(555)-200-0520 **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