

Serology EIA Assays

Bio-Rad offers a continually expanding and comprehensive menu of Serology Infectious Disease products, including reagent kits, instrumentation, software, and accessories.

Convenient, Objective, Simple

- Easy-to-use kits with standardized procedure and common reagents
- Microplate strip format, with color-coded wells
- Designed for manual and automated use
- Rapid turnaround time
- Ready-to-use reagents

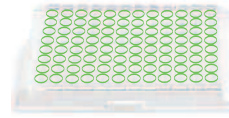
BIO-RAD

Serology EIA Kits

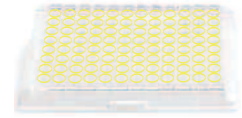
Ordering Information

Catalog No.	Description	Tests
Measles		
25181	Measles IgG EIA	96 tests
Mumps		
25182	Mumps IgG EIA	96 tests
Rubella		
25173	Rubella IgG EIA	96 tests
25174	Rubella IgM EIA	96 tests
Varicella-Zoster Virus (VZV)		
25180	VZV IgG EIA	96 tests
Epstein-Barr Virus		
25183	EB-NA IgG EIA	96 tests
25184	EB-VCA IgG EIA	96 tests
25185	EB-VCA IgM EIA	96 tests
Toxoplasmosis		
25175	Toxoplasma IgG EIA	96 tests
25176	Toxoplasma IgM EIA	96 tests
Cytomegalovirus		
25177	CMV IgG EIA	96 tests
25178	CMV IgM EIA	96 tests
Herpes Simplex Virus		
25179	HSV (1 and 2) IgG EIA	96 tests
Helicobacter Pylori*		
25225	Pylori Detect IgA Qualitative	96 tests
25226	Pylori Detect IgG Qualitative	96 tests
Borreliosis (Lyme Disease)*		
32507	Bio-Rad Anti-Borrelia (Lyme) EIA	96 tests

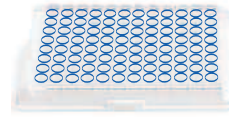
* No color-coded microplate
 * No common reagents
 * Different procedure



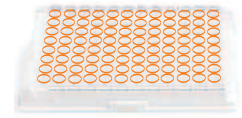
Rubella IgG EIA
Green



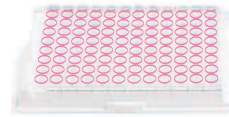
Rubella IgM EIA
Yellow



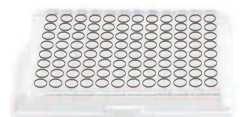
Toxoplasma IgG EIA
Blue



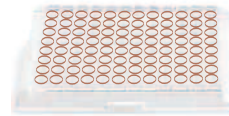
Toxoplasma IgM EIA
Orange



CMV IgG EIA
Pink



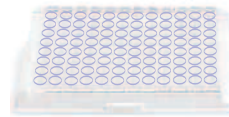
CMV IgM EIA
Grey



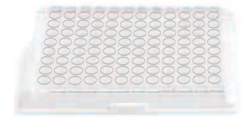
HSV IgG EIA
Brown



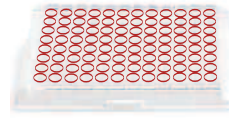
VZV IgG EIA
Clear



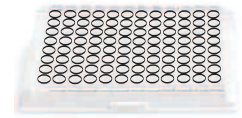
Measles IgG EIA
Lilac



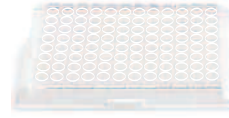
Mumps IgG EIA
Silver



EB NA IgG EIA
Red



EB VCA IgG EIA
Black



EB VCA IgM EIA
White



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-(55)-5200-0520 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