

GS HIV-2 EIA

The only licensed HIV assay specifically for the detection of HIV-2 in blood screening and diagnostic testing of human serum and plasma

The GS HIV-2 EIA assay can be used as part of the CDC recommended algorithm for HIV-1/HIV-2 EIA testing.

Confidence in Results

- Sensitivity 100%*
- Specificity 99.8%
- Reproducible results

Convenience

- Microplate strip format
- Results in 3-4 hours

* In a study of known positives, determination of sensitivity for the HIV-2 was limited due to the difficulty of obtaining an adequate number of samples



GS HIV-2 EIA

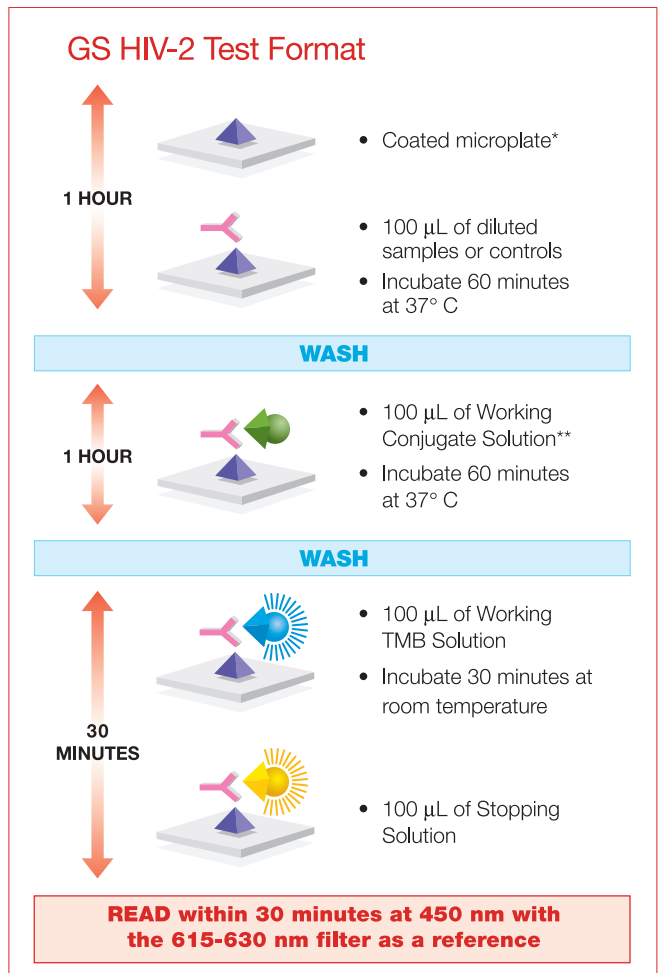
Table 1: Reproducibility of the HIV-2 EIA

Inter-Assay Reproducibility				
Specimen	N	Mean OD/CO	SD	% CD
1	36	4.927	1.010	20.5%
2	36	3.530	0.789	22.4%
3	36	2.799	0.897	32.0%
4	36	2.366	0.449	19.0%
5	36	1.821	0.401	22.0%
6	18	0.838	0.279	33.3%
7	16	0.205	0.056	27.3%
HIV-2 EIA Positive Control	276	1.228*	0.226	18.4%
HIV-2 EIA Negative Control	412	0.063*	0.018	28.6%

* Controls are expressed as mean OD values.

Intra-Assay Reproducibility				
Specimen	N	Mean OD	SD	% CD
HIV-2 EIA Positive Control	288	0.900	0.040	4.4%
HIV-2 EIA Negative Control	288	0.060	0.008	13.3%
Positive Sample	288	0.585	0.028	4.8%

Reproducibility was determined by assaying seven specimens, along with the HIV-2 EIA Positive and HIV-2 EIA Negative Controls, at six different sites. Three lots of materials were used. Intra-assay reproducibility was determined by assaying replicates of the HIV-2 EIA Positive and HIV-2 EIA Negative Control as well as one positive sample. Three lots of materials were also used in this testing, and the resultant reproducibility was analyzed by applying the pooled estimate of standard deviation. Results of reproducibility testing are shown in Table 1.



- * Microplate coated with: HIV-2 gp 36.
- ** Conjugate: Goat anti-human IgM and IgG horseradish peroxidase.

Ordering Information

Catalog No.	Description	
32536	GS HIV-2 EIA.....	480 tests



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 +36-1-459-6100 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