

GS HIV-1 Western Blot

A supplemental and more specific test for the detection of antibodies to HIV-1

Confidence in Results

- 100% sensitivity in a study of known positives
- Clear band identification for a precise interpretation of results

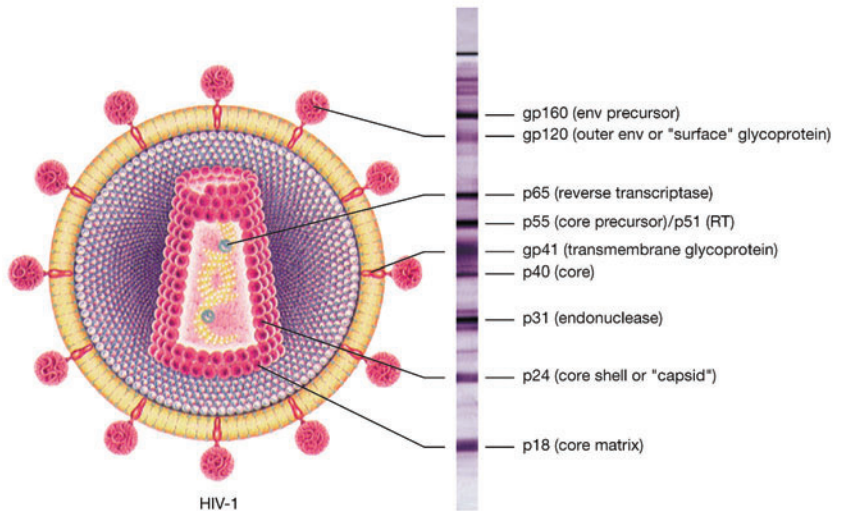
Convenience

- Fast turnaround for a same day result
- Easy storage and retrieval with pre-wetted strips in a resealable pouch
- User-friendly procedure with ready-to-use reagents



GS HIV-1 Western Blot

Detection and identification of specific HIV-1 antibodies in human serum, plasma, or dried blood spots.



100% Sensitivity in a Study of Known Positives

Competitive Performance (Serum/Plasma)						
Clinical Group	EIA Status	Study Results	GS HIV-1 Western Blot	Competitor A	Competitor B	Competitor C
AIDS/ARC	RR	Number Studied	172	36	184	52
		Positive	172 (100%)	29 (80.5%)	178 (96.7%)	52 (100%)
		Indeterminate	0 (0%)	7 (19.5%)	6 (3.3%)	0 (0%)
		Negative	0 (0%)	0 (0%)	0 (0%)	0 (0%)

10.7% Indeterminate Rate in Low Risk, EIA NR Populations

Competitive Performance (Serum/Plasma)						
Clinical Group	EIA Status	Study Results	GS HIV-1 Western Blot	Competitor A	Competitor B	Competitor C
Low Risk	NR	Number Studied	298	709	254	86
		Positive	0 (0%)	0 (0%)	0 (0%)	0 (0%)
		Indeterminate	32 (10.7%)	92 (13.0%)	41 (16.1%)	13 (15.1%)
		Negative	266 (89.3%)	617 (87.0%)	213 (83.9%)	73 (84.9%)

Ordering Information

Catalog No.	Description	
32508	GS HIV-1 Western Blot.....	40 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 36-1-455-8800 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