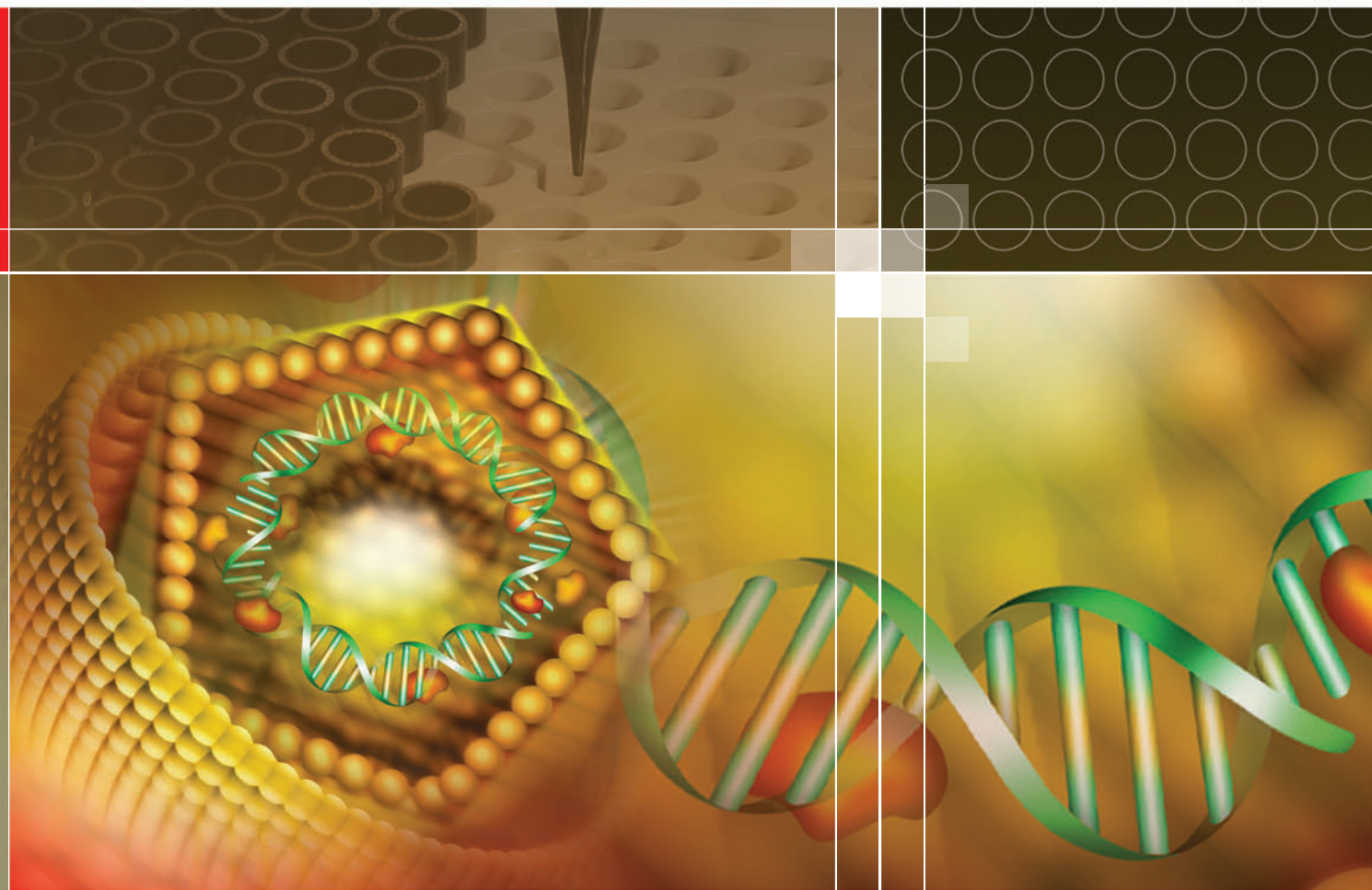


GS HBsAg EIA 3.0

Superior Hepatitis Testing from Bio-Rad





GS HBsAg EIA 3.0

Outstanding sensitivity of less than 0.1 ng/mL of ad/ay HBsAg*

The GS HBsAg EIA 3.0 is an EIA for the detection of Hepatitis B Surface Antigen (HBsAg) in human serum, plasma, and cadaveric serum samples.

GS HBsAg Confirmatory Assay 3.0

For Confirmatory Neutralization of GS HBsAg EIA 3.0 Reactive Specimens

* Calibrated versus PHA 805 commercial panel

Ease-of-Use

- No sample predilution
- Total incubation time = 2.5 hours
- Common reagents can be used with other Bio-Rad EIA assays
- Manual and automated use

Advanced Safety Features

- Bar-coded reagents
- Color-coded reagent labels
- Microplate strip identification
- Color changes during protocol steps

Earliest Hepatitis B Infection Detection Demonstrated on 21 Commercial Seroconversion Panels*

GS HBsAg EIA 3.0 Reactivity on Commercial HBsAg Seroconversion Panels Versus Competitors				
Panel I.D.#	GS HBsAg EIA 3.0		HBsAg EIA Competitor #1 (overnight)	HBsAg EIA Competitor #2
	Static	Shaker		
902	7	7	9	9
903	3	2	5	5
904	2	1	3	3
905	1	1	1	1
906	2	2	2	3
907	6	5	6	6
908	6	4	7	7
909	4	3	4	5
910	3	3	3	4
911	20	20	21	22
912	7	6	8	8
914	2	2	5	5
915	1	1	9	11
917	2	2	3	3
918	2	2	3	3
919	5	4	6	7
920	3	3	3	3
924	3	2	3	4
931	5	3	NT	7
932	10	8	NT	10
40565L	2	1	3	NT

* The sensitivity of GS HBsAg EIA 3.0, procedures A and B, was assessed using 20 seroconversion panels purchased from Boston Biomedica, Inc. (West Bridgewater, MA), and 1 panel from Serologicals, Inc. (Clarkston, GA). The number of the bleed at which HBsAg was detected (S/C \geq 1.0) by the GS HBsAg EIA 3.0, procedures A and B, is compared to other licensed HBsAg assays (as recorded in the certificates of analysis).

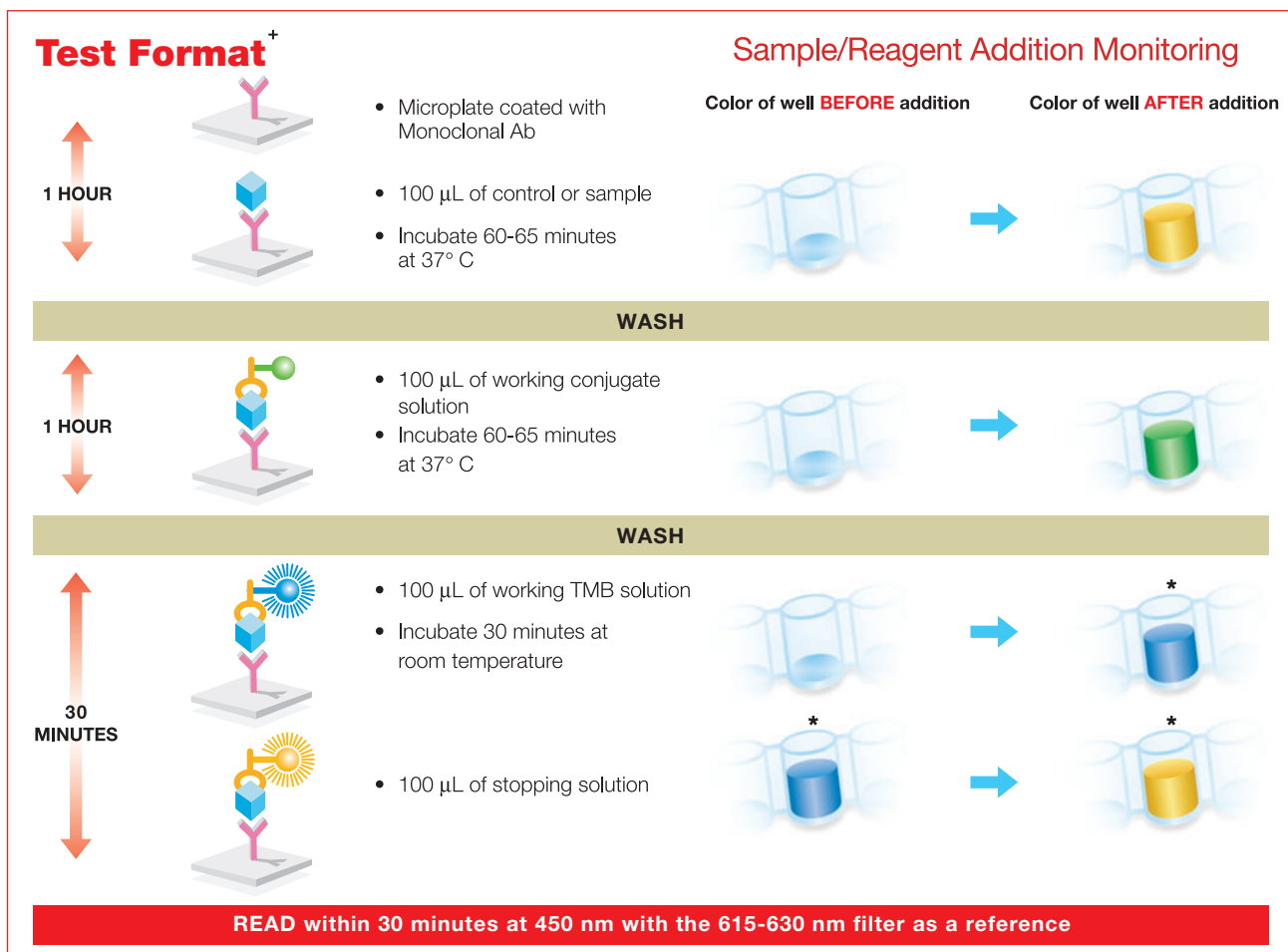
GS HBsAg 3.0 Static Procedure

- Earlier detection on 12 out of 19 seroconversions for HBsAg 3.0 versus competitor assay number 1
- Earlier detection on 16 out of 20 seroconversions for HBsAg 3.0 versus competitor assay number 2

GS HBsAg 3.0 Shaking Procedure

- Earlier detection on 15 out of 19 seroconversions versus competitor assay number 1
- Earlier detection on 18 out of 20 seroconversions versus competitor assay number 2

GS HBsAg EIA 3.0 Testing Procedure



Anti-HBsAg Monoclonal Ab
 HBsAg
 Anti HBsAg Monoclonal Ab Conjugate

+ Static or Shaker. * Reactive samples after 30 minute incubation.

Ordering Information

Enzyme immunoassay (EIA) for the detection of Hepatitis B Surface Antigen (HBsAg) in Human Serum, Plasma, and Cadaveric Serum Specimens:

Catalog No.	Description	Tests
32591	GS HBsAg EIA 3.0	480 tests
32592	GS HBsAg EIA 3.0	960 tests
25258	GS HBsAg EIA 3.0	4800 tests

For Confirmatory Neutralization of GS HBsAg EIA 3.0 Reactive Specimens:

Catalog No.	Description	Tests
32594	GS HBsAg Confirmatory Assay 3.0.....	25 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