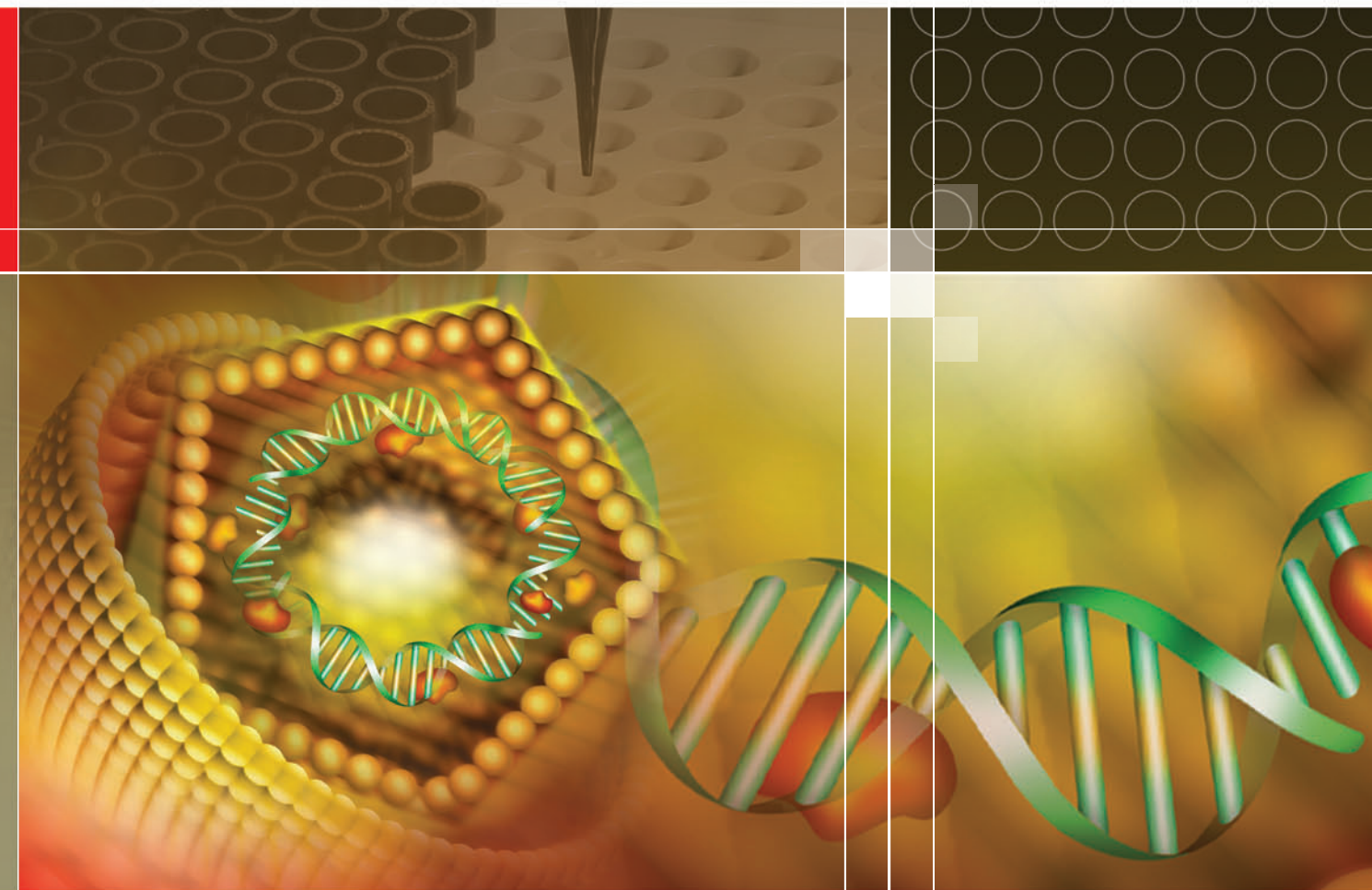


# GS HBsAg EIA 3.0

**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 ≥ 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).

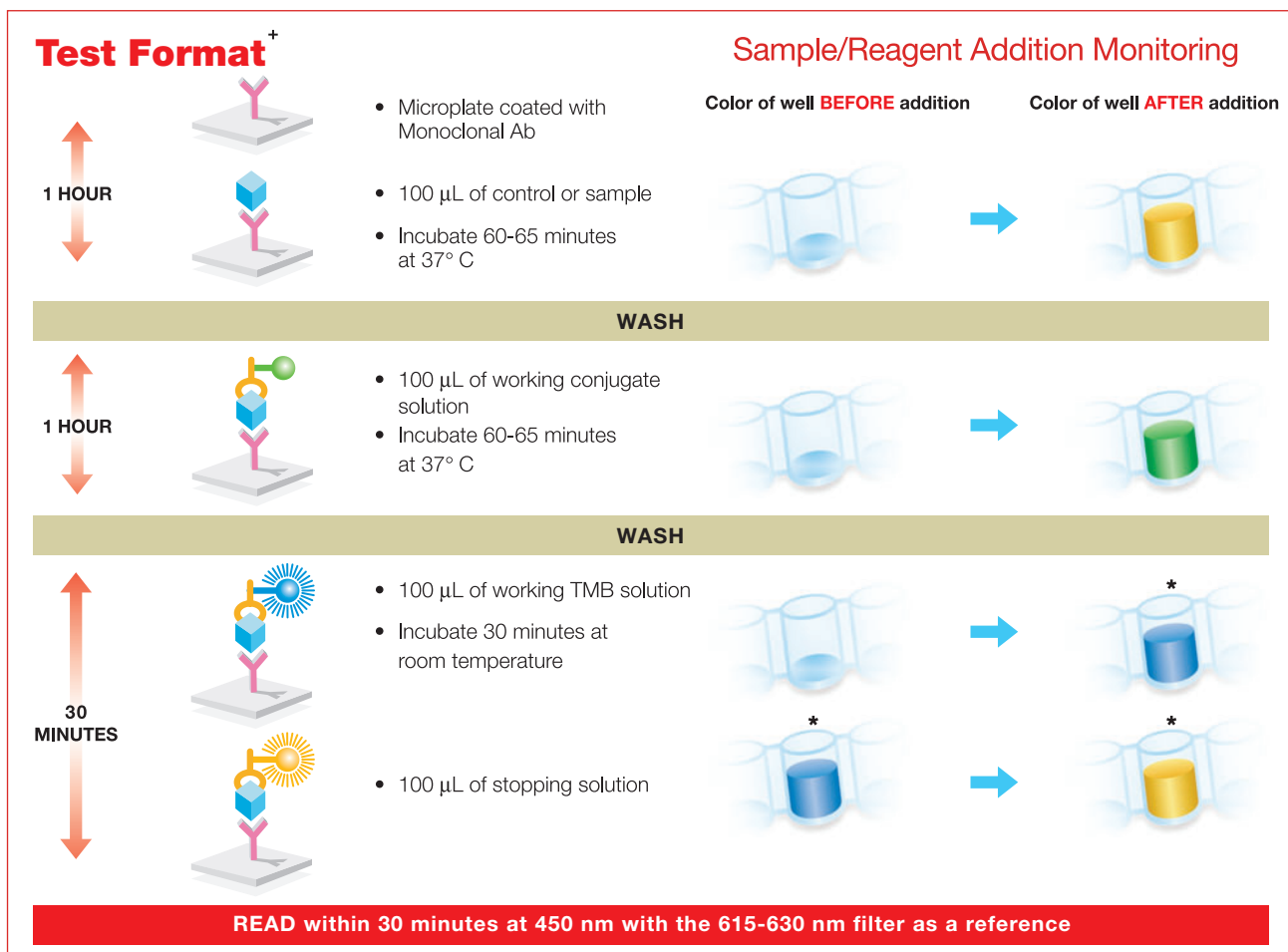
### GS HBsAg 3.0 Static Procedure

- 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](http://www.bio-rad.com/diagnostics).*

Clinical Diagnostics Group

**Website** [www.bio-rad.com/diagnostics](http://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