

## ab287884 – GMS Staining ELISA Kit

For use in the histologic visualization of fungi, basement membrane and some opportunistic organisms such as *Pneumocystis carinii*  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab287884>

### Storage and Stability

All the reagents are shipped at room temperature (RT) and stored at multiple temperatures.

### Materials Supplied

Item	Quantity	Storage Condition
Silver Nitrate Solution (0.2%)	125 mL	2-8°C
Methenamine Solution	125 mL	2-8°C
Gold Chloride Solution (0.2%)	125 mL	2-8°C
Borax Solution	15 mL	RT
Sodium Bisulfite Solution	125 mL	RT
Chromic Acid Solution	125 mL	RT
Sodium Thiosulfate Solution (5%)	125 mL	RT
Light Green Solution	125 mL	RT

### Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully utilize this assay:

- Distilled water
- Coplin jars
- Plastic forceps
- Absolute alcohol
- Synthetic resin

### Reagent Preparation

- Do not use if reagents become cloudy.
- Do not use past expiration date.
- Use caution when handling reagents.
- All glassware used in this procedure should be chemically cleaned and rinsed thoroughly in distilled water.
- Failure to adequately remove the alcohol used in deparaffination will result in reduction of the chromic acid solution.
- Reduction of the chromic acid solution will result in a change in color from orange to brown. Discard the reagent if color change is noted.
- Do not use metal forceps to remove slides from reagents. Use plastic forceps only.
- Pre-warm all reagents to room temperature prior to use.

### Assay Procedure - Standard

1. Deparaffinize sections if necessary and hydrate to distilled water.
2. Incubate slide in Chromic Acid Solution for 10 minutes.
3. Rinse in tap water followed by 2 changes of distilled water.
4. Incubate slide in Sodium Bisulfite Solution for 1 minute (to remove any residual chromic acid).

5. Rinse in tap water followed by 2 changes of distilled water.
6. Combine the following for a working GMS solution:

- a. 25 ml Silver Nitrate Solution (0.2%)
- b. 25 ml Methenamine Solution
- c. 2 ml Borax Solution

**ΔNote:** Mixed solution may not be stored for reuse later.

7. Place working GMS solution in 60° centigrade water bath and allow temperature to equilibrate.
8. Incubate slide in working GMS solution for 10-15 minutes. Using plastic forceps, dip slide in distilled water and check under a microscope for evaluation of silver impregnation. Fungi should be dark brown. If color is not sufficient, return the slide to working GMS solution for 2-3 minutes and check again.
9. Rinse in 4 changes of distilled water.
10. Incubate slide in Gold Chloride Solution for 15-30 seconds.
11. Rinse in 4 changes of distilled water.
12. Incubate slide in Sodium Thiosulfate Solution (5%) for 2 minutes.
13. Rinse in tap water followed by 2 changes of distilled water.
14. Incubate slide in Light Green Solution for 2 minutes.
15. Rinse slide using absolute alcohol.
16. Dehydrate in 2 changes of absolute alcohol, clear, and mount in synthetic resin.

### Assay Procedure – Microwave

**Note:** These instructions were developed using a standard 500-watt microwave oven. Heating times should be modified as needed depending on the microwave oven used.

1. Deparaffinize sections if necessary and hydrate to distilled water.
2. Place slide in plastic coplin jar filled with Chromic Acid solution. Cap jar loosely!
3. Place jar in microwave oven and heat on high power for 10 seconds. Allow slide to remain in warm solution for 3 minutes.
4. Rinse in tap water followed by 2 changes of distilled water.
5. Incubate slide in Sodium Bisulfite solution for 1 minute (to remove any residual chromic acid).
6. Rinse in tap water followed by 2 changes of distilled water.
7. Combine the following for a working GMS solution:

- a. 25 ml Silver Nitrate
- b. 25 ml Methenamine
- c. 2 ml Borax Solution

**ΔNote:** Mixed solution may not be stored for reuse later.

8. Place working GMS solution (loosely capped) in microwave oven for 40 seconds. Remove and pour several times between coplin jar and a clear graduated cylinder to mix thoroughly (use protective glove to avoid burning hand). Mixed solution remains in coplin jar.
9. Incubate slide in working GMS solution (heated) for 2-6 minutes until the tissue is medium brown in color. Using plastic forceps, dip slide in distilled water and check under a microscope for evaluation of silver impregnation. Fungi should be dark brown. If color is not sufficient, return the slide to working GMS solution for 1-2 minutes and check again. Reheat solution if needed.
10. Rinse in 4 changes of distilled water.

11. Incubate slide in Gold Chloride solution for 15-30 seconds.
12. Rinse in 4 changes of distilled water.
13. Incubate slide in Sodium Thiosulfate for 2 minutes.
14. Rinse in tap water followed by 2 changes of distilled water.
15. Incubate slide in Light Green Solution for 2 minutes.
16. Rinse slide using absolute alcohol.
17. Dehydrate in 2 changes of absolute alcohol, clear, and mount in synthetic resin.

#### **Technical Support**

Copyright © 2021 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)