

# Acid Fast Bacteria Staining Kit

(Cat# K1424-30, -125; Store at RT)

## I. Introduction:

The **Acid Fast Bacteria (AFB) Staining Kit** is intended for use in the histological visualization of Acid Fast Bacteria and Tubercle Bacilli. This kit is a rapid 15-minute procedure. The lipoid capsule of the acid-fast organism takes up Carbol Fuchsin and resists decolorization. **Staining Interpretation: Acid Fast Organisms: Bright Red; Background: Light Green**

## II. Application:

- Histological applications, IHC-P (use at an assay dependent concentration)
- For *in vitro* Diagnostic Use

## III. Sample Type:

- Any well-fixed paraffin embedded or frozen tissue section cut at 3-5  $\mu\text{m}$
- Cell smear (air dried)

## IV. Kit Contents:

Components	K1424-30	K1424-125	Part Number	Storage Temperature
Carbol Fuchsin Solution	30 ml	125 ml	K1424-XX-1	RT
Acid Alcohol Solution (0.5%)	2 x 30 ml	500 ml	K1424-XX-2	RT
Light Green Solution	30 ml	125 ml	K1424-XX-3	RT

## V. User Supplied Reagents and Equipment:

- Distilled water
- Coplin jars
- Xylene/Xylene Substitute
- Synthetic resin

## VI. Shipment and Storage:

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

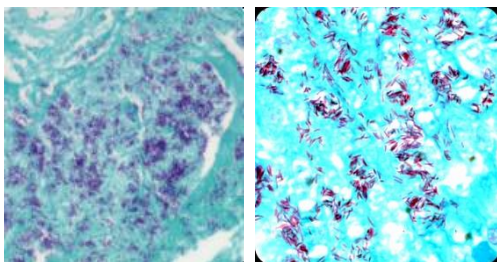
## VII. Reagent Preparation:

- Carbol Fuchsin Solution should be filtered when a thick sheen develops on top of solution. To avoid possible contamination, gloves should be worn when performing this procedure. Do not use tap water prior to application of Carbol Fuchsin Solution as it is reported that Acid Fast Bacteria can be found in some systems. Therefore, use of distilled water is recommended whenever possible.
- Do not use if reagents become cloudy.
- Do not use past expiration date.
- Use caution when handling reagents.
- Non-Sterile.

## VIII. Protocol:

1. Deparaffinize sections if necessary and hydrate to distilled water.
2. Incubate slide in Carbol Fuchsin Solution for 15 minutes.
3. Rinse for 2 minutes in running tap water followed by 2 changes of distilled water.
4. Decolorize in Acid Alcohol Solution (0.5%) until sections are a pale pink color.
5. Dip slide twice in distilled water to rinse.
6. Counterstain in Light Green Solution for 1-2 min.
7. Rinse in several changes of distilled water.
8. Dehydrate quickly through graded alcohols ending with 2 changes in absolute alcohol.
9. Clear, and mount in synthetic resin.

## IX. Data:



Staining using Acid Fast Bacteria Staining Kit

FOR RESEARCH USE ONLY! Not to be used on humans.