

# Pneumocystis Staining Kit

(Cat# K1434-30, -125; Store at Multiple Temperatures)

## I. Introduction:

The Pneumocystis Staining Kit is intended for use in the histological visualization of *Pneumocystis carinii* in cytology smears, and paraffin or frozen tissue sections.

***Pneumocystis carinii*: Violet/Purple**

**Connective Tissue: Blue / Green**

**Erythrocytes: Yellow / Green**

**Mucin: Rose/Purple**

**Cartilage: Rose/Purple**

## II. Application:

- Histological applications
- For *in vitro* diagnostic use

## III. Sample Type:

- Paraffin or frozen tissue sections.
- Cytology smears
- Control Tissue: Any prepared slide that contains *Pneumocystis carinii*

## IV. Kit Contents:

Components	K1434-30	K1434-125	Part Number	Storage Temperature
Cresyl Echt Violet Solution (0.1%)	30 ml	125 ml	K1434-XX-1	2-8°C
Naphthol Yellow S Solution	30 ml	125 ml	K1434-XX-2	RT
Staining Jar	30 ml	60 ml	K1434-XX-3	RT

## V. User Supplied Reagents and Equipment:

- Distilled water
- Forceps
- Absolute alcohol
- Synthetic resin
- Glacial Acetic Acid
- Sulfuric Acid

## VI. Shipment and Storage:

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

## VII. Reagent Preparation:

- Do not use if reagents become cloudy.
- Do not use past expiration date.
- Use caution when handling reagents.
- Non-Sterile

## VIII. Procedure (Standard):

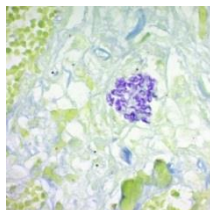
**Preparation of Sulfation Reagent Prior to Beginning (Not Included in Kit):** NOTE: Wear protective clothing, gloves, and eyewear when mixing and handling this reagent. Make fresh for each use.

1. Pour 15 ml of Glacial Acetic Acid into the Staining Jar provided with this kit.
2. Slowly add 5 ml of Sulfuric Acid to the Staining Jar.
3. Screw cap tightly on staining jar and invert several times to thoroughly mix acids.
4. Wait 5-10 minutes before proceeding with stain procedure to allow mixed acids to cool.
5. Use Staining Jar only for Sulfation procedure.

### Protocol:

1. Deparaffinize sections if necessary and hydrate to distilled water.
2. Place slide in freshly mixed Sulfation Reagent for 10 min. Note: Agitate staining Jar every few min to keep acids mixed.
3. Rinse in 3 changes of distilled water.
4. Incubate slide in Cresyl Echt Violet Solution (0.1%) for 10-15 min. Note: Agitate slide several times during incubation step.
5. Rinse quickly in distilled water.
6. Apply Naphthol Yellow S Solution to tissue section for no longer than 4 sec. Note: Excessive Naphthol Yellow S Solution decolorizes *Pneumocystis* and connective tissue.
7. Rinse very quickly using absolute alcohol.
8. Dehydrate very quickly in 2 changes of fresh Absolute Alcohol. Alternative Method: Dip slide twice in Absolute Alcohol and air-dry slide.
9. Clear, and mount in synthetic resin.

## IX. Data:



Pneumocystis Staining Kit

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