

VitroView™ Iron Prussian Blue Stain Kit

Cat. No. VB-3009

Introduction

Small amounts of iron are found normally in spleen and bone marrow. Excessive amounts are present in hemochromatosis, with deposits found in the liver and pancreas, as well as in hemosiderosis, with deposits in the liver, spleen, and lymph nodes. This kit is designed for staining ferric iron in tissue sections.

Kit Components

Potassium Ferrocyanide Solution - 125 ml
Hydrochloric Acid Solution - 125 ml
Nuclear Fast Red Solution - 250 ml

Storage

Room temperature. Stable for 6 months.

Protocol

1. Deparaffinize and hydrate to 70% alcohol.
2. Rinse in distilled water.
3. Prepare Working Solution: Mix potassium ferrocyanide solution and hydrochloric acid solution in a 1:1 ratio, and let stand for 30 minutes at room temperature.
4. Note: Make Working Solution fresh. Discard after use.
5. Allow slides to stand in Working Solution for 5 minutes in the fume hood.
6. Rinse in distilled water.
7. Let stand in Nuclear-fast red solution for 5 minutes. Wash in tap water.
8. Dehydrate with 2 changes of 95% Ethanol and 2 changes of 100% Ethanol (2 minute per change).
9. Clear with 3 changes of xylene (5 minute per change).
10. Mount coverslip onto glass slide with Permount or some other suitable organic mounting medium.

Results

- Iron (hemosiderin) blue
- Nuclei red
- Background pink

Precautions

Handle with care. Avoid contact with eyes, skin and clothing. Do not ingest. Wear gloves.