

VitroView™ Modified Gomori's Trichrome Stain Kit for Frozen Section

Cat. No. VB-3014

Introduction

Trichrome stain is one of the most highly utilized special stains in the Histopathology Laboratory. Most of common uses for trichrome staining are liver biopsies, renal biopsies, dermatopathology, cardiac biopsies and muscle and nerve biopsies. Gomori's trichrome is a one-step staining procedure that combines plasma stain (chromotrope 2R) and connective fiber stain (fast green FCF) in a phosphotungstic acid solution to which glacial acetic acid has been added. This kit is designed for frozen tissue section.

Kit Components

1. Gomori's Trichrome solution -----250 ml
2. Harris Hematoxylin solution-----250 ml

Storage

Room temperature.

Protocol

1. Sample preparation: Cut 10 - 16 micron (12 µm) sections in a cryostat from snap frozen tissue. Fixation is not needed.
2. Immerse sections in Harris Hematoxylin for 5 minutes.
3. Wash with tap water until the water is clear.
4. Immerse sections in Gomori trichrome stain for 10 minutes.
5. Differentiate using 0.2% acetic acid. A few dips should be sufficient.
6. Immerse rack with sections directly into 95 % alcohol
7. Continue to dehydrate in ascending alcohol solutions (95% x 2, 100% x 2).
8. Clear with xylene (3 - 4 x) in a columbia staining dish (jar)s - Thomas Scientific #8542-E30.
9. Mount coverslip onto a labeled glass slide with Permount or some other suitable organic mounting medium.

Results

1. Nuclei: Dark blue
2. muscle myofibrils: Green-blue
3. Mitochondria and endoplasmic reticulum stain -red
4. Connective tissue stains-pale green-blue
5. Myelin stains-purple red
6. Type 1 fibers satin darker blue green as compared to type 2 fibers.

Precautions

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