

VitroView™ Congo Red Amyloid Stain Kit

Cat. No. VB-3011

Introduction

This kit is used for the detection of amyloid in formalin-fixed, paraffin-embedded tissue sections with amyloidosis, and may be used for frozen sections as well. Amyloid deposits will stain red, while nuclei will stain blue. The thickness of sections is usually 5 μ m. However, in cases containing low amyloid deposits, 10 μ m thick sections will give better results.

Kit Components

1. Congo Red Solution - 250 ml
2. Alkaline Alcohol Solution - 250 ml
3. Mayer's Hematoxylin Solution - 250 ml

Storage

Room temperature.

Protocol

1. For paraffin-embedded tissue sections, deparaffinize and hydrate to 70% alcohol. Rinse in distilled water. For frozen sections, dry sections for 5 minutes. Fix sections in neutral formalin solution for 20 minutes. Rinse in distilled water.
2. Stain in congo red solution for 15-20 minutes.
3. Rinse in distilled water.
4. Differentiate quickly (5-10 dips) in alkaline alcohol solution.
5. Rinse in tap water for 1 minute.
6. Counterstain with Mayer's hematoxylin for 30 seconds.
7. Rinse in tap water for 2 minutes.
8. Dehydrate through 95% alcohol, 2 changes of 100% alcohol, 3 minutes each.
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

- Amyloid, elastic fibers, eosinophil granules ----- red
- Nuclei ----- blue

Positive Controls

Know amyloidosis containing tissue (i.e. Alzheimer's disease).

Precautions

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