

VitroView™ Sudan Black B Stain Kit

Cat. No. VB-3012

Introduction

Sudan Black is a slightly basic dye that will combine with acidic groups in compound lipids. Sudan Black will also stain phospholipids. This kit is used for lipid and fat staining in formalin fixed frozen sections. It is not suitable for paraffin embedded tissue sections. Lipid and fat will stain blue-black, while cell nuclei will stain red.

Kit Components

1. Sudan Black B Solution - 100 ml
2. Pre-Stain Solution - 100 ml
3. Differentiation Solution - 100 ml
4. Nuclear Fast Red Solution - 100 ml

Storage

Room temperature.

Protocol

1. Fix frozen sections in 10% formalin for 20 minutes.
2. Wash well in tap water, rinse in distilled water, drain off excess water.
3. Wash with Pre-Stain Solution, two changes, 5 minutes each.
4. Treat with Sudan Black, 7 minutes, agitate.
5. Treat with Differentiation Solution, 3 minutes.
6. Rinse in distilled water.
7. Treat with Nuclear Fast Red, 3 minutes.
8. Wash in water.
9. Wash in tap water, rinse in distilled.
10. Mount with aqueous mounting media, Glycerin Jelly.

Results

- Fat blue-----black
- Nuclei----- red

Precautions

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