

Reticulum Stain Kit (Connective Tissue Stain) ab150684

4 References 1 Image

Overview

Product name	Reticulum Stain Kit (Connective Tissue Stain)
Sample type	Tissue sections
Product overview	The Reticulum Stain Kit (Connective Tissue Stain) is intended for use in histological demonstration of reticular fibers. The main function of reticular fibers is to provide support. They are normally found throughout the body, particularly in liver, lymph node, spleen and kidney. Ammoniacal silver stains are the most commonly used methods for demonstration of reticular fibers.

Other products for staining tissue sections

Find more kits and reagents in the [special stains guide](#), or products for antigen retrieval, blocking, signal amplification, visualization, counterstaining, and mounting in the [IHC kits and reagents guide](#).

Notes

Staining Interpretation

Reticulum	Black
Nuclei	Red

Control Tissue: Liver, Kidney, Lymph Node, Spleen

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	100 tests
Ferric Ammonium Sulfate Solution (3%)	1 x 125ml
Formalin Solution (20%)	1 x 125ml

Components	100 tests
Gold Chloride Solution (0.1%)	1 x 125ml
Nuclear Fast Red Solution (Enhanced)	1 x 125ml
Potassium Metabisulfite Solution (3%)	1 x 125ml
Potassium Permanganate Solution (1%)	1 x 125ml
Silver Nitrate Solution (10%)	5 x 10ml
Sodium Hydroxide Solution (3%)	1 x 125ml
Sodium Thiosulfate Solution (5%)	1 x 125ml

Images



Staining in Liver using ab150684 - Reticulum Stain Kit.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

-
- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors