

## Product datasheet

# Reticulum Stain Kit (Modified Gomori's) ab236473

1 Image

### Overview

<b>Product name</b>	Reticulum Stain Kit (Modified Gomori's)
<b>Sample type</b>	Tissue sections
<b>Product overview</b>	The Reticulum Stain Kit (Modified Gomori's) 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.

### Notes

#### Staining Interpretation

Reticulum	Grey/Black
Nuclei	Pink/Red

**Control Tissue:** Liver, Kidney, Lymph Node, Spleen

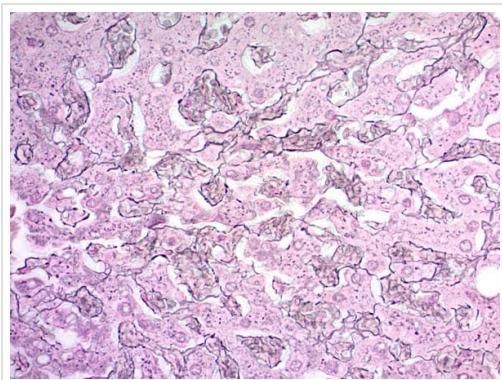
### Properties

**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
Gold Chloride Solution (0.2%)	1 x 125ml
Nuclear Fast Red Solution (Enhanced)	1 x 125ml
Potassium Hydroxide Solution (10%)	1 x 15ml
Potassium Metabisulfite Solution (3%)	1 x 125ml
Potassium Permanganate Solution	1 x 250ml
Silver Nitrate Solution (10%)	1 x 65ml

Components	100 tests
Sodium Thiosulfate Solution (5%)	1 x 125ml
Sulfuric Acid Solution (1N)	1 x 15ml

## Images



ab236473 - Reticulum Stain Kit (Modified Gomori's)

ab236473 Reticulum Stain Kit (Modified Gomori's) staining formalin-fixed-paraffin embedded human liver.

**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