

Product datasheet

Colloidal Iron Stain Kit (Mucin Stain) ab150664

2 Images

Overview

Product name Colloidal Iron Stain Kit (Mucin Stain)
Product overview The Colloidal Iron (Mucin Stain) is designed for the histological visualization of acid mucopolysaccharides.

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

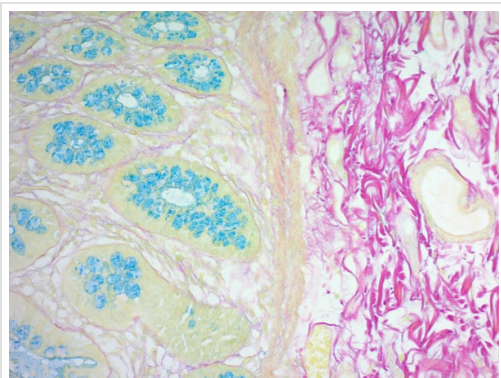
Acid Mucopolysaccharides	Blue
Collagen	Red-Purple

Control Tissue: Colon, Small Intestine

Properties

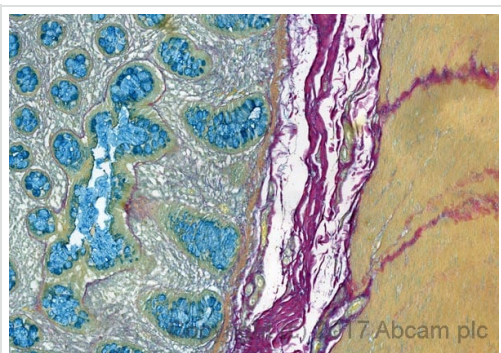
Storage instructions Store at room temperature. Please refer to protocols.

Components	100 tests
Acetic acid Solution	1 x 500ml
Colloidal Iron Stock Solution	1 x 125ml
Hydrochloric acid Solution	1 x 125ml
Potassium Ferrocyanide Solution	1 x 125ml
Van Gieson's Solution	1 x 125ml



ab150664 - Colloidal Iron Stain Kit (Mucin Stain)

ab150664 Colloidal Iron Stain Kit (Mucin Stain) staining formalin-fixed-paraffin embedded human GI tract.



ab150664 - Colloidal Iron Stain Kit (Mucin Stain)

ab150664 (Colloidal Iron stain) staining acid mucopolysaccharides (blue) and collagen (red-purple) in formalin fixed paraffin embedded human colon.

*Tissue obtained from the Human Research Tissue Bank, supported by the NIHR Cambridge Biomedical Research Centre.

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