

Product datasheet

Pneumocystis Stain Kit (Microorganism Stain) ab150682

2 Images

Overview

Product name Pneumocystis Stain Kit (Microorganism Stain)
Product overview The Pneumocystis (Microorganism Stain) is intended for use in the histological visualization of *Pneumocystis carinii* in cytology smears, and paraffin or frozen tissue sections.

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

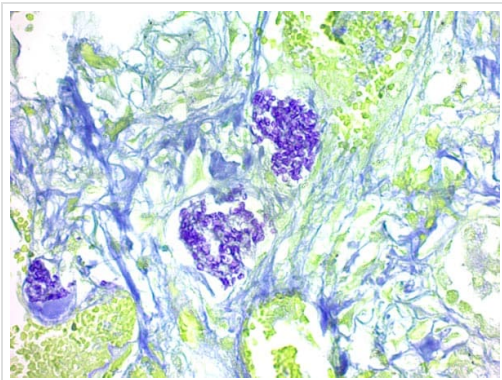
<i>Pneumocystis carinii</i>	Violet/Purple
Connective Tissue	Blue/Green
Erythrocytes	Yellow
Mucin	Rose/Purple
Cartilage	Rose/Purple

Control Tissue: Any prepared slide that contains *Pneumocystis carinii*.

Properties

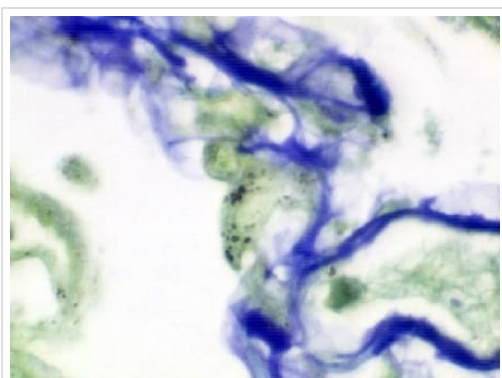
Components	100 tests
Cresyl Echt Violet Solution	1 x 125ml
Naphthol Yellow S Solution	1 x 125ml
Staining Jar	1 unit

Images



ab150682 - Pneumocystis Stain Kit (Microorganism Stain)

ab150682 Pneumocystis Stain Kit (Microorganism Stain) staining formalin-fixed-paraffin embedded *Pneumocystis carinii*.



ab150682 - Pneumocystis Stain Kit (Microorganism Stain)

Staining using ab150682 - Pneumocystis 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