

Product datasheet

Phosphotungstic Acid Hematoxylin (PTAH) Stain Kit (Connective Tissue Stain) ab150683

3 Images

Overview

Product name	Phosphotungstic Acid Hematoxylin (PTAH) Stain Kit (Connective Tissue Stain)
Sample type	Tissue sections
Product overview	The PTAH Stain Kit (Connective Tissue Stain) is intended for use in the histological visualization of collagen, striated muscle, glial fibers and collagen without using Zenker's Fixative containing mercuric chloride.

This kit may be used on formalin-fixed, paraffin-embedded or frozen 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

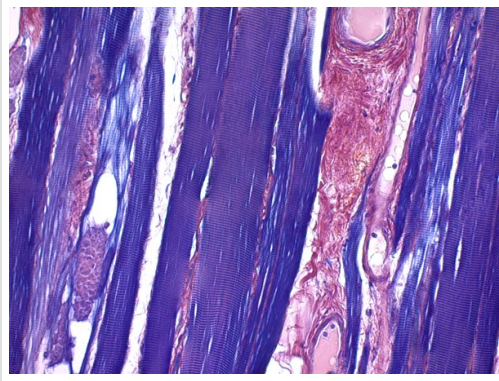
Fibrin, striated muscle, glial fibers	Blue
Collagen	Brown/Red
Nuclei	Blue

Control Tissue: Striated muscle.

Properties

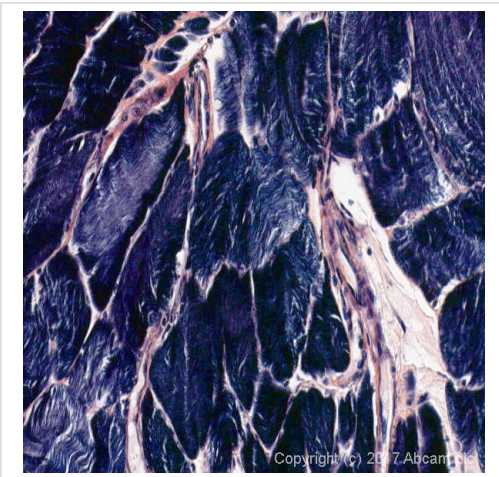
Storage instructions Store at room temperature. Please refer to protocols.

Components	100 tests
Ferric Ammonium Sulfate Solution	1 x 125ml
PTAH Solution	1 x 125ml
Zinc chloride solution	1 x 500ml



ab150683 - Phosphotungstic Acid Hematoxylin (PTAH) Stain Kit (Connective Tissue Stain)

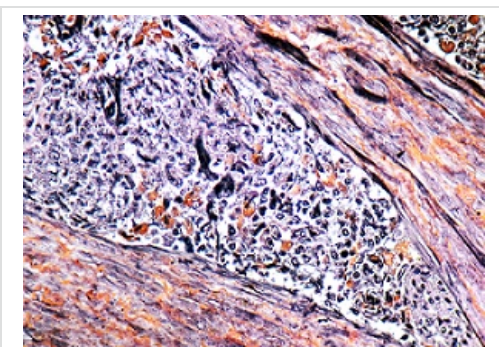
ab150683 Phosphotungstic Acid Hematoxylin (PTAH) Stain Kit (Connective Tissue Stain) staining formalin-fixed-paraffin embedded human skeletal muscle.



ab150683 - Phosphotungstic Acid Hematoxylin (PTAH) Stain Kit (Connective Tissue Stain)

ab150683 (Phosphotungstic Acid Hematoxylin (PTAH) Stain) staining striated muscle (blue) and collagen (red) in formalin fixed paraffin embedded normal human muscle

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



ab150683 - Phosphotungstic Acid Hematoxylin (PTAH) Stain Kit (Connective Tissue Stain)

Staining using ab150683 - Phosphotungstic acid Hematoxylin (PTAH) Stain Kit.

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