

ab150681 – Picro Sirius Red Stain Kit (Connective Tissue Stain)

For the histological visualization of collagen I and III fibers in addition to muscle in tissue sections.

For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

www.abcam.com/ab150681 (use www.abcam.cn/ab150681 for China, or www.abcam.co.jp/ab150681 for Japan)

Introduction:

Picro Sirius Red Stain Kit (Connective Tissue Stain) ab150681 is intended for use in the histological visualization of collagen I and III fibers in addition to muscle in tissue sections.

Picro Sirius Red Stain Kit may not be suitable for frozen sections.

The Picro Sirius red staining may be viewed using standard light microscopy. Picro Sirius red staining may also be viewed using polarized light to more easily differentiate the collagen fibers from the background. Studies have used the differences between the appearance of different collagen fibers under polarized light to differentiate between different types of collagen fibers; it is unclear if the differences in appearance of the fibers is due to orientation, thickness and packing of the fibers rather than reflecting different types of collagen fibers.

Staining Interpretation:

Light Microscopy

Collagen	Red
Muscle Fibers	Yellow
Cytoplasm	Yellow

Polarized Light Microscopy

Type I (Thick fibers)	Yellow/Orange Birefringence
Type III (Thin fibers)	Green Birefringence

Control Tissue:

Lung, Uterus, Muscle, Kidney.

Storage and Stability: Store kit at room temperature immediately on receipt. Kit can be stored for 1 year from receipt, if components have not been reconstituted. Keep away from open flame and refer to the safety datasheet.

Materials Supplied:

Item	Quantity	Storage temperature (before prep)
Picro Sirius Red Solution	250 mL	Room Temperature
Acetic Acid Solution (0.5%)	2 x 250 mL	Room Temperature

Materials Required, Not Supplied:

These materials are not included in the kit, but will be required to successfully perform this assay:

Absolute alcohol

Distilled water

General guidelines, precautions, and troubleshooting:

Please observe safe laboratory practice and consult the safety datasheet.

For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide: www.abcam.com/assaykitguidelines

For typical data produced using the assay, please see the assay kit datasheet on our website.

Staining Protocol:

Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.

1. Deparaffinize sections if necessary and hydrate in distilled water.
2. Apply adequate Picro Sirius Red Solution to completely cover the tissue section and incubate for 60 minutes.
3. Rinse slide quickly in 2 changes of Acetic Acid Solution.
4. Rinse slide in absolute alcohol.
5. Dehydrate in 2 changes of absolute alcohol.
6. Clear slide and mount in synthetic resin.

FAQs / Troubleshooting:

General troubleshooting points are found at www.abcam.com/assaykitguidelines.

Technical Support

Copyright © 2020 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

www.abcam.com/contactus

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)