

Version 2a Last updated 15 January 2021

ab245887

Picro-Sirius Red Stain Kit (Cardiac Muscle)

For the histological visualization of the thin septa and collagen fibers.

View Picro-Sirius Red Stain Kit datasheet:

<https://www.abcam.com/ab245887>

[use www.abcam.cn/ab245887 for China, or www.abcam.co.jp/ab245887 for Japan]

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Materials Supplied and Storage	2
3. General guidelines, precautions, and troubleshooting	3
4. Assay Procedure	4
5. Notes	5

1. Overview

The Picro-Sirius Red Stain Kit (For Cardiac Muscle) is intended for use in the histological visualization of thin septa and collagen fibers. This modification of the Picro-Sirius Red stain eliminates the yellow cytoplasmic staining that can obscure thin collagenous septa. Using this procedure easily allows viewing of collagenous septa as thin as 0.2-0.5 microns. The PSR stain may be viewed using standard light microscopy or polarized light resulting in birefringence of the collagen fibers.

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

Staining Interpretation:

Collagen:	Red
Septa:	Red
Cytoplasm:	Green/Blue

ΔNote: Collagen appears Yellow-Orange and Green Birefringence under Polarized Light Microscopy.

Control Tissue:

Cardiac muscle, uterus, lung and kidney tissues.

2. Materials Supplied and Storage

Store kit at Room temperature immediately on receipt and check below for storage for individual components. 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.

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
Phosphomolybdic Acid Solution (0.2%)	250 mL	RT	RT
Picro-Sirius Red Solution	250 mL	RT	RT
Acetic Acid Solution (0.5%)	250 mL	RT	RT

3. 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.

4. Assay Procedure

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

- 1.1 Deparaffinize sections if necessary and hydrate to distilled water.
- 1.2 Apply adequate Phosphomolybdic Acid Solution (0.2%) to completely cover tissue section and incubate for 1-5 minutes.
- 1.3 Dip slide one time in distilled water.
- 1.4 Apply adequate Picro-Sirius Red Solution to completely cover tissue section and incubate for 60-90 minutes.
- 1.5 Rinse slide quickly in two changes of Acetic Acid Solution (0.5%).
- 1.6 Rinse slide using absolute alcohol.
- 1.7 Dehydrate in 2 changes of absolute alcohol, clear, and mount in synthetic resin.

5. Notes

Technical Support

Copyright © 2021 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)