

ab150678 – Oil Red O Stain Kit (Lipid Stain)

For the histological visualization of fat cells and neutral fat.
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit:

www.abcam.com/ab150678 (use www.abcam.cn/ab150678 for China, or www.abcam.co.jp/ab150678 for Japan)

Overview:

The Oil Red O (Lipid Stain) (ab150678) is intended for use in the histological visualization of fat cells and neutral fat. This kit may be used only on frozen tissue sections, fresh smears, or touch preps.

Staining Interpretation:

Fat Cells	Red
Neutral Fat	Red
Nuclei	Blue

Control Tissue: Any frozen section containing fat.

Storage and Stability: 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.

Materials Supplied:

Item	Quantity	Storage temperature (before prep)
Hematoxylin (Modified Mayer's)	125 mL	RT
Propylene Glycol	500 mL	RT
Oil Red O Solution	125 mL	RT

Materials Required, Not Supplied:

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

- Aqueous Mounting Medium.

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 prior to use.

ΔNote: Heat Oil Red O Solution to 60°C prior to use.

1. Prepare fresh or frozen tissue sections as usual.
2. Place slide in Propylene Glycol for 5 minutes.
3. Incubate slide in heated (60°C) Oil Red O Solution for 6-10 minutes or overnight at room temperature.

ΔNote: To improve staining, incubation time may be increased.

4. Prepare a mixture of 85% Propylene Glycol in distilled water and differentiate the tissue section in this mixture for 1 minute.

ΔNote: To improve staining, differentiation time may be reduced.

5. Rinse slide in 2 changes of distilled water.
6. Incubate slide in Hematoxylin for 1-2 minutes.
7. Rinse slide thoroughly in tap water.
8. Rinse slide in 2 changes of distilled water.
9. Coverslip using an aqueous mounting medium.

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)