

Oil Red O Stain Kit (Lipid Stain) ab150678

★★★★★ [1 Abreviews](#) [79 References](#) [画像数 4](#)

製品の概要

製品名 Oil Red O Stain Kit (Lipid Stain)

製品の概要 Oil Red O (Lipid Stain) kit 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.

If you prefer to prepare your own Oil Red O solution, we recommend our solid [Oil Red O stain ab146295](#).

- Oil Red O staining protocol summary:
- prepare fresh or frozen tissue sections
 - incubate slide in propylene glycol for 2 min
 - incubate slide in oil red o solution for 6 min
 - differentiate section in 85% propylene glycol for 1 min
 - rinse slide twice in water
 - incubate in hematoxylin for 1-2 min
 - rinse slide three times in water
 - coverslip with an aqueous mounting medium

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](#).

特記事項

Oil Red O is a fat-soluble dye that stains neutral triglycerides and lipids. It cannot be used with formaldehyde-fixed paraffin embedded sections as the alcohols used remove most lipids.

Staining Interpretation

Fat Cells	Red
Neutral Fat	Red
Nuclei	Blue

Control Tissue: Any frozen section containing fat.

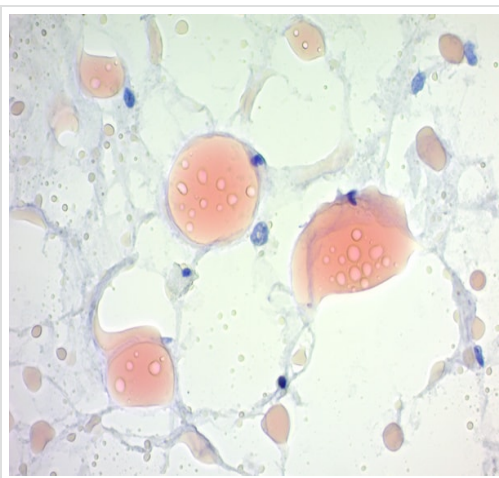
製品の特性

保存方法

Store at room temperature. Please refer to protocols.

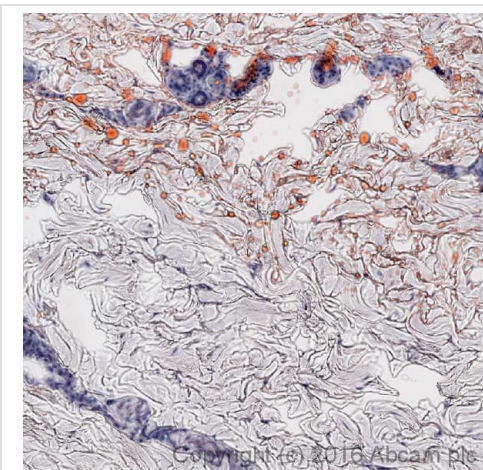
内容	100 tests
Hematoxylin (Modified Mayer's Solution)	1 x 125ml
Oil Red O Solution	1 x 125ml
Propylene Glycol	1 x 500ml

画像



ab150678 - Oil Red O Stain Kit (Lipid Stain)

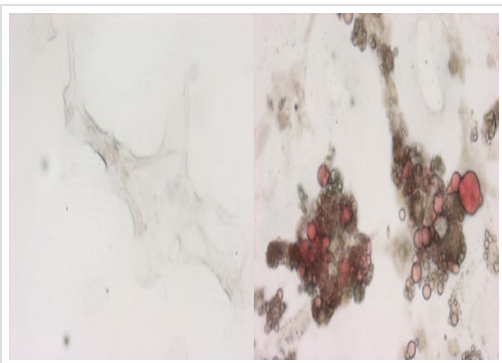
ab150678 Oil Red O Stain Kit (Lipid Stain) staining frozen human human adipose.



ab150678 - Oil Red O Stain Kit (Lipid Stain)

ab150678 (Oil Red O stain) staining lipid deposits (red) in frozen normal human skin. The dermis is unstained, whereas the hypodermis is stained, demonstrating dye specificity. Hematoxylin (blue) is used as a counterstain.

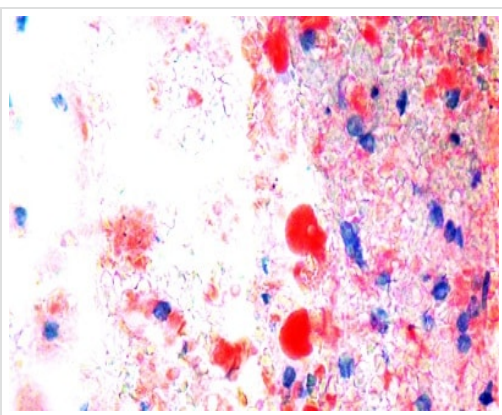
*Tissue obtained from the Human Research Tissue Bank, supported by the NIHR Cambridge Biomedical Research Centre.



ab150678 - Oil Red O Stain Kit (Lipid Stain)

Oil Red O staining of 3T3-L1 and adipocyte-like cells

During adipocyte differentiation, cells accumulate lipid. Oil Red O staining kit (ab150678) was used to visualize lipid in 3T3-L1 cells (left) and adipocyte-like cells (right). As expected, lipid accumulates in 3T3-L1 cells after differentiation into adipocyte-like cells.



ab150678 - Oil Red O Stain Kit (Lipid Stain)

Staining of Frozen Breast Tissue using ab150678 - Oil Red O 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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors