

Version 3 Last updated 25 June 2018

ab150663 Congo Red Stain Kit (Amyloid Stain)

For the histological staining of Amyloid in tissue sections.

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

Table of Contents

1. Overview	3
2. Materials Supplied and Storage	4
3. Materials Required, Not Supplied	5
4. General guidelines, precautions, and troubleshooting	6
5. Assay Procedure	7
6. FAQs / Troubleshooting	8
7. Typical Data	9
8. Notes	10

1. Overview

The Amyloid Stain (Congo Red) is intended for use in the histological visualization of amyloid in tissue sections. Examination under a polarizing microscope results in green birefringence of amyloid.

Staining Interpretation:

Amyloid	Red to Pink
Erythrocytes	Light Orange
Eosinophil Granules	Orange to Red
Nuclei	Blue

Control Tissue:

Freshly cut sections containing amyloid.

Cut sections 6-12 microns to show smaller amyloid deposits.

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 Data Sheet.

Congo Red solution is flammable.

Item	Quantity	Storage temperature (before prep)	Storage temperature (after prep)
Congo Red Solution	500 mL	Room temperature	Room temperature
Hematoxylin	500 mL	Room temperature	Room temperature
Bluing Reagent	500 mL	Room temperature	Room temperature

3. Materials Required, Not Supplied

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

- 95% Ethanol
- 100% Ethanol
- Distilled water

4. 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.

5. Reagent Preparation

5.1 Congo Red Solution

Ready to use as supplied.

5.2 Hematoxylin

Ready to use as supplied.

5.3 Bluing Reagent

Ready to use as supplied.

6. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature just prior to use and gently agitate.
- 6.1 Deparaffinize sections if necessary and hydrate to distilled water.
- 6.2 Add 4-6 drops of Hematoxylin and incubate for 5 minutes.
- 6.3 Rinse slide in tap water.
- 6.4 Apply 4-6 drops of Bluing Reagent and incubate for 30 seconds.
- 6.5 Rinse slide in distilled water.
- 6.6 Dip slide in 95% alcohol for 5 seconds
- 6.7 Stain slide with Congo Red Solution for 20 minutes.
- 6.8 Dip twice (quickly) in 100% alcohol.
- 6.9 Dip repeatedly (4-5 dips) in clearing agent, and mount in synthetic resin.

7. FAQs / Troubleshooting

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

8. Notes

Technical Support

Copyright © 2012-2018 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | 400 921 0189 / +86 21 2070 0500

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abc.com | +64-(0)9-909-7829