

Version 4b Last updated 21 April 2026

ab150675

Luxol Fast Blue Stain

(Myelin Stain)

For the histological identification of basic neuronal structures in brain or spinal cord sections.

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

Table of Contents

1. Overview	5
2. General guidelines, precautions, and troubleshooting	7
3. Materials Supplied and Storage	3
4. Materials required, not supplied	3
5. Staining Protocol	4
6. FAQs / Troubleshooting	4
8. Notes	5

1. Overview

The Luxol Fast Blue Stain (Myelin Stain)(ab150675) is designed for staining myelin/myelinated axons and Nissil substance on formalin fixed, paraffin-embedded tissue as well as frozen tissue. This product is used for identifying the basic neuronal structure in brain or spinal cord sections.

Staining Interpretation:

Myelinated Fibers	Blue
Nissil Substance	Violet
Nerve cells	Violet

Control Tissue:

Cerebral cortex, Spinal cord.

2. 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.

If applicable, please refer to the current Safety Data Sheet (SDS) provided with this product for safety, handling, and disposal information. The most up to date and current versions are available on our website <https://www.abcam.com/en-us>.

3. Materials Supplied and Storage

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

Keep away from open flame and refer to the safety datasheet.

Item	Quantity (30 mL pack size)	Quantity (100 Test/125 mL pack size)	Storage temperature (before prep)
Cresyl Echt Violet Solution	30 mL	125 mL	4°C
Luxol Fast Blue Solution	60 mL	125 mL	RT
Lithium Carbonate Solution (0.05%)	60 mL	500 mL	RT
Alcohol Reagent (70%)	60 mL	500 mL	RT

4. Materials Required, Not Supplied

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

- Absolute alcohol

5. Staining Protocol

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

- 5.1** Deparaffinize sections if necessary and hydrate in distilled water.
- 5.2** Incubate slide in Luxol Fast Blue Solution for 24 hours at room temperature or 2 hours at 60°C.
- 5.3** Rinse slide thoroughly in distilled water.
- 5.4** Differentiate the section by dipping in Lithium Carbonate Solution several times (up to 20 seconds).
- 5.5** Continue differentiation by repeatedly dipping in Alcohol Reagent until gray-matter is colorless and the white-matter remains blue.
- 5.6** Rinse slide in distilled water.
- 5.7** Incubate slide in Cresyl Echt Violet for 2-5 minutes.
- 5.8** Rinse quickly in distilled water.
- 5.9** Dehydrate quickly in 3 changes of absolute alcohol.
- 5.10** Clear and mount in synthetic resin.

6. FAQs / Troubleshooting

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

7. Notes

Technical Support

Copyright © 2026 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:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)