

Version 3a Last updated 24 September 2025

ab150687 Von Kossa (Calcium Stain)

For the histological visualization of calcium deposits in tissue sections.

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. Materials Required, Not Supplied	3
4. General guidelines, precautions, and troubleshooting	4
5. Staining Protocol	5
6. FAQs / Troubleshooting	6
7. Typical Data	7
8. Notes	8

1. Overview

The Von Kossa (Calcium Stain)(150687) is intended for use in the histological visualization of calcium deposits in paraffin sections.

Staining Interpretation:

Calcium in mass deposits	Black
Calcium in dispersed deposits	Gray
Nuclei	Red
Cytoplasm	Pink

Control Tissue:

Any paraffin-embedded tissue containing calcium 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 datasheet.

Item	Quantity (30 mL pack size)	Quantity (100 test/ 125 mL pack size)	Storage temperatur e (before prep)
Silver Nitrate Solution (5%)	30 mL	125 mL	4°C
Sodium Thiosulfate Solution (5%)	30 mL	125 mL	RT
Nuclear Fast Red Solution	30 mL	125 mL	RT

3. Materials Required, Not Supplied

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

- Absolute alcohol

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. 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 Silver Nitrate Solution (5%) for 30-60 minutes with exposure to ultraviolet light or 100 watt incandescent.
- 5.3** Rinse in several changes of distilled water.
- 5.4** Incubate slide in Sodium Thiosulfate Solution (5%) for 2-3 minutes.
- 5.5** Rinse for 2 minutes in running tap water followed by 2 changes of distilled water.
- 5.6** Incubate slide in Nuclear Fast Red Solution for 5 minutes.
- 5.7** Rinse for 2 minutes in running tap water followed by 2 changes of distilled water.
- 5.8** Dehydrate very quickly in 3 changes of fresh Absolute Alcohol.
- 5.9** Clear and mount in organic mounting medium or synthetic resin.

6. FAQs / Troubleshooting

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

7. Notes

Technical Support

Copyright © 2025 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)