

Proteinase K Antigen Retrieval Kit

Cat. No. VB-6034S

Introduction

The proteinase K based solution is designed to break the protein cross-links, therefore unmask the antigens and epitopes in formalin-fixed and paraffin embedded tissue sections, thus enhancing staining intensity of antibodies

Component

- Proteinase K Stock Solution (20X) 1ml
- Proteinase K Dilution Buffer 20ml

Storage

Proteinase K Stock Solution (20X) should be stored at –20 °C.

Proteinase K Stock Solution (20X) can be aliquoted and stored at –20 °C for 2-3 years.

Protocol

1. Deparaffinize sections in 2 changes of xylene, 5 minutes each.
2. Hydrate in 2 changes of 100% ethanol for 2 minutes each, 95% and 75% ethanol for 2 minute each. Then rinse in distilled water.
3. Proteinase K Stock Solution (20X) should be diluted 1:20 for working solution (1x) using Proteinase K Dilution Buffer.
4. Cover sections with Proteinase K working solution and incubate 10-20 minutes at 37 °C in humidified chamber (optimal incubation time may vary depending on tissue type and degree of fixation, and should be determined by user).
5. Allow sections to cool at room temperature for 10 minutes.
6. Rinse sections in PBS Tween 20 for 2x2 min.
7. Proceed to standard immunohistochemistry protocol.

Note: Microwave, pressure cooker and autoclave can be used as alternative heating source to replace steamer or water bath.

Note

Avoid contact with eyes, skin and clothing. Do not ingest. Wear gloves.