

Catalog Number: NB900-66727

Product Name: Proteinase K

Format: Stabilized and clear 25X Concentrate

Volume/UOM: 2 mL

Intended Use: For In Vitro Diagnostic Use

Product Description

Proteinase K is specifically designed to achieve proteolytic digestion of formalin fixed, paraffin embedded (FFPE) tissue sections prior to performing in situ hybridization procedures.

Proteolytic digestion of formalin fixed tissues improves the accessibility of target DNA to biotinylated probes, which results in better in situ staining.

Storage and Handling

Store at 2-8C. Proteinase K is stable for 6 months. For long term storage, aliquot Proteinase K into smaller sizes and store at -20C.

Preparation of Working Solutions

Place 100 uL Proteinase K solution in a glass tube, add 2.4 mL distilled water, and mix.

Protocol Recommendations

Optimal digestion conditions will vary depending on fixation conditions and should be optimized by the end user. Incubation of tissue sections for 10-30 minutes at 37C is sufficient for in situ.

Performance Characteristics

The protocols for a specific application can vary. These include, but are not limited to: fixation, heat-retrieval method, incubation times, tissue section thickness and detection kit used. Due to the superior sensitivity of these unique reagents, the recommended incubation times and titers listed are not applicable to other detection systems, as results may vary. Ultimately, it is the responsibility of the investigator to determine optimal conditions.

Precautions

1. Wear disposable gloves when handling reagents.
2. Microbial contamination of reagents may result in an increase in nonspecific staining.
3. Incubation times or temperatures other than those specified may give erroneous results. The user must validate any such change.