

**Dako
Proteinase K
Ready-to-use**

Code S3020

Intended use

For In Vitro Diagnostic Use.

Proteinase K is used for the proteolytic digestion of formalin-fixed, paraffin-embedded tissue sections prior to immunohistochemical (IHC) and in situ hybridization (ISH) procedures.

Refer to Dako's *General Instructions for Immunohistochemical Staining* or the detection system instructions of IHC procedures for: 1) Principle of Procedure, 2) Materials Required, Not Supplied, 3) Storage, 4) Specimen Preparation, 5) Staining Procedure, 6) Quality Control, 7) Troubleshooting, 8) Interpretation of Staining, 9) General Limitations.

Summary and principle

Proteolytic digestion of formalin-fixed, paraffin-embedded tissue sections improves the accessibility of antibodies and probes to target sites within the tissue. Proteolytic digestion exposes certain epitopes/sites which have been masked during fixation.

Reagent provided

Proteinase K: Proteolytic enzyme solution diluted in 0.05 mol/L Tris-HCl, 0.015 mol/L sodium azide, pH 7.5.

Ready-to-Use Proteinase K solution is available in manual and automated sizes.

Precautions

1. For professional users.
2. Wear appropriate Personal Protective Equipment to avoid contact with eyes and skin.
3. Unused solution should be disposed of according to local, State and Federal regulations.
4. Safety Data Sheet available for professional users on request.

Storage

Store at 2–8 °C.



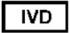






Procedure

The time required for optimal digestion of formalin-fixed, paraffin-embedded tissue will vary with the extent of fixation. Generally, digestion of tissue for 3 to 6 minutes at room temperature is sufficient. Digestion time may be extended up to 15 minutes if necessary.

Limitations

Over-digestion of tissue sections may result in the loss of tissue morphology or cause sections to detach from the slides. For IHC and ISH, either Silanized Slides or poly-L-lysine coated slides are recommended.

Explanation of symbols

 Catalogue number	 Temperature limitation	 In vitro diagnostic medical device
 Manufacturer	 Batch code	 Contains sufficient for <n> tests
 Use by	 Consult instructions for use	 Authorized representative in the European Community



Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051
United States

Tel. +44 161 492 7050
www.agilent.com

TX02438/01

Revision 2020.07