

# Proteinase K, Recombinant

Catalog Numbers 25530015 and 25530031

Doc. Part No. 25530015.pps Pub. No. MAN0001455 Rev. C



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](https://www.thermofisher.com/support).

## Product description

Proteinase K is a non-specific serine protease. It is not activated by metal ions, chelating agents (for example, EDTA), sulfhydryl reagents, or by trypsin or chymotrypsin inhibitors. It is stable over a wide pH range (4–12.5), with optimal activity at pH 6.5–9.5. Activity can be stimulated by addition of denaturing agents (SDS and urea). The optimal temperature for the enzyme is 65°C (12 times more active at 65°C than at 25°C). Rapid denaturation of the enzyme occurs at temperatures above 65°C.

Autolysis of the enzyme occurs increasingly at alkaline pH. However, Proteinase K is not completely inactivated by autolysis. Some enzyme fragments continue to maintain their complete proteolytic activity, even after extensive autolysis.

## Contents and storage

Item	Cat. No. <a href="#">25530015</a>	Cat. No. <a href="#">25530031</a>	Storage
Proteinase K, Recombinant, ≥30 Units/mg <sup>[1]</sup>	100 mg	1 g	2°C–8°C

<sup>[1]</sup> One mAnson unit is described as that amount of enzyme that liberates 1 μmole of Folin-positive amino acid within one minute at 37°C, using hemoglobin as a substrate.

## Reconstitute Proteinase K

Reconstitute the Proteinase K, using one of the following options:

- Dissolve in 10 mM Tris HCl (pH 7.5), 20 mM calcium chloride, and 50% glycerol, then store at –20°C.
- Dissolve in 50 mM Tris HCl (pH 8.0), 1–5 mM calcium acetate, then store at 2°C–8°C.

**Note:** Storage at –20°C in the absence of glycerol can cause precipitation of the Proteinase K. Bacterial growth can occur in solutions stored at 2°C–8°C over prolonged periods of time. Ca<sup>2+</sup> can serve as a stabilizer to suppress autolysis.

## Applications

Proteinase K is used to rapidly inactivate endogenous nucleases, such as RNases and DNases when isolating RNA or DNA from tissues and cell lines. This enzyme can also be used to remove nucleases in the preparation of tissue sections for *in situ* hybridization.

## Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](https://www.thermofisher.com/support).

[thermofisher.com/support](https://www.thermofisher.com/support) | [thermofisher.com/askaquestion](https://www.thermofisher.com/askaquestion)



Life Technologies Corporation | 5781 Van Allen Way | Carlsbad, California 92008 USA

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2016–2024 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

**For Research Use Only. Not for use in diagnostic procedures.**