

Proteinase K

Promega Corporation
2800 Woods Hollow Rd., Madison, WI 53711



**This product uses less packaging
and generates less waste to support
sustainable practices in your lab.**

Materials Affected

- Proteinase K (Lyophilized) (Cat.# V3021)
- Proteinase K (PK) Solution, 4ml (Cat.# MC5005)
- Proteinase K (PK) Solution, 16ml (Cat.# MC5008)
- Proteinase K (PK) Solution, 23ml (Cat.# A5051)
- Maxwell® HT 96 gDNA Blood, Isolation System, 1 × 96 preps (Cat.# A2670)
- Maxwell® HT 96 gDNA Blood Isolation System, 4 × 96 preps (Cat.# A2671)
- ReliaPrep™ Large Volume HT gDNA Isolation System (Cat.# A2751)

Introduction

We are committed to reducing the environmental effects from our operations globally, and with the products that we provide. We are working to reduce greenhouse gas emissions, energy, water, waste and effects from product distribution as part of our Corporate Responsibility Program. To learn more, visit: promega.com/responsibility

This green sheet discloses the basis of our environmental claims and highlights how the Proteinase K formulation and packaging reduces its environmental effects.

Product Description

Proteinase K is a serine protease that exhibits broad cleavage activity. This protease cleaves peptide bonds adjacent to the carboxylic group of aliphatic and aromatic amino acids and is useful for general digestion of protein in biological samples. In addition to its ability to digest native proteins, Proteinase K is stable in urea and SDS, making it useful for a variety of applications, including chromosomal DNA preparation for pulsed-field gel electrophoresis, protein fingerprinting and nuclease removal from purified DNA and RNA.

Sustainability Features

Proteinase K is stable across a variety of temperature ranges and retains >80% activity at temperatures up to 60°C. This stability has the following environmental benefits:

- **Fewer Packaging Materials and Less Waste:** No dry ice, gel ice or expanded polystyrene (EPS) foam cooler is required for transportation.
- **Reduced Greenhouse Gas Emissions:** Shipping smaller and lighter boxes decreases fuel consumption during transport.
- **Energy Conservation:** Not needing ultra-low temperature storage at Promega and customer sites saves energy.

Note: Proteinase K (Lyophilized) does not require cold storage as shipped, but does when reconstituted.