



PRODUCT INFORMATION

KpnI

#ER0522 5x4000 U

Lot: ____ **Expiry Date:** __

5'...**G G T A C↓C**...3'
3'...**C↑C A T G G**...5'

Concentration: 10 U/μL
Source: *Klebsiella pneumoniae* OK8
Supplied with: 4x1 mL of 10X Buffer KpnI
1 mL of 10X Buffer Tango

Store at -20°C



In total 10 vials.

BSA included

www.thermoscientific.com/onebio

RECOMMENDATIONS

1X Buffer KpnI (for 100% KpnI digestion)
10 mM Tris-HCl (pH 7.5), 10 mM MgCl₂,
0.02% Triton X-100, 0.1 mg/mL BSA.

Incubation temperature
37°C.

Unit Definition

One unit is defined as the amount of KpnI required to digest 1 μg of lambda DNA-BamHI fragments in 1 hour at 37°C in 50 μL of recommended reaction buffer.

Dilution

Dilute with Dilution Buffer (#B19): 10 mM Tris-HCl (pH 7.4 at 25°C), 100 mM KCl, 1 mM EDTA, 1 mM DTT, 0.2 mg/mL BSA and 50% glycerol.

Double Digests

Thermo Scientific Tango Buffer is provided to simplify buffer selection for double digests. 98% of Thermo Scientific restriction enzymes are active in a 1X or 2X concentration of Tango™ Buffer. Please refer to www.thermoscientific.com/doubledigest to choose the best buffer for your experiments.

1X Tango Buffer: 33 mM Tris-acetate (pH 7.9 at 37°C), 10 mM magnesium acetate, 66 mM potassium acetate, 0.1 mg/mL BSA.

Storage Buffer

KpnI is supplied in: 10 mM Tris-HCl (pH 7.5 at 25°C), 50 mM KCl, 1 mM DTT, 0.1 mM EDTA, 0.2 mg/mL BSA and 50% glycerol.

Recommended Protocol for Digestion

- Add:
 - nuclease-free water 16 µL
 - 10X Buffer KpnI 2 µL
 - DNA (0.5-1 µg/µL) 1 µL
 - KpnI 0.5-2 µL^{*,**}
 - Mix gently and spin down for a few seconds.
 - Incubate at 37°C for 1-16 hours^{**}.
- The digestion reaction may be scaled either up or down.

Recommended Protocol for Digestion of PCR Products Directly after Amplification

- Add:
 - PCR reaction mixture 10 µL (~0.1-0.5 µg of DNA)
 - nuclease-free water 18 µL
 - 10X Buffer KpnI 2 µL
 - KpnI 1-2 µL^{*,**}
- Mix gently and spin down for a few seconds.
- Incubate at 37°C for 1-16 hours^{**}.

* This volume of the enzyme is recommended for preparations of standard concentrations (10 U/µL), whereas HC enzymes (50 U/µL) should be diluted with Dilution Buffer to obtain 10 U/µL concentration.

** See Overdigestion Assay.

Thermal Inactivation

KpnI is inactivated by incubation at 80°C for 20 min.

ENZYME PROPERTIES

Enzyme Activity in Thermo Scientific REase Buffers, %

KpnI	B	G	O	R	Tango	2X Tango
100	20-50	0-20	0-20	0-20	20-50	0-20

Methylation Effects on Digestion

Dam: never overlaps – no effect.
Dcm: may overlap – no effect.
CpG: may overlap – no effect.
EcoKI: never overlaps – no effect.
EcoBI: never overlaps – no effect.

Stability during Prolonged Incubation

A minimum of 0.2 units of the enzyme is required for complete digestion of 1 µg of lambda DNA in 16 hours at 37°C.

Digestion of Agarose-embedded DNA

A minimum of 5 units of the enzyme is required for complete digestion of 1 µg of agarose-embedded lambda DNA in 16 hours.

Number of Recognition Sites in DNA

λ	ΦX174	pBR322	pUC57	pUC18/19	pTZ19R/U	M13mp18/19
2	0	0	1	1	1	1

For **CERTIFICATE OF ANALYSIS** see back page

CERTIFICATE OF ANALYSIS

Overdigestion Assay

No detectable change in the specific fragmentation pattern is observed after an 80-fold overdigestion with KpnI (5 U/μg lambda DNA × 16 hours).

Ligation and Recleavage (L/R) Assay

The ligation and recleavage assay was replaced with LO test after validating experiments showed LO test ability to trace nuclease and phosphatase activities with sensitivity that is higher than L/R by a factor of 100.

Labeled Oligonucleotide (LO) Assay

No detectable degradation of single-stranded or double-stranded labeled oligonucleotides occurred during incubation with 10 units of KpnI for 4 hours.

Blue/White (B/W) Cloning Assay

The B/W assay was replaced with LO test after validating experiments showed LO test ability to detect nuclease and phosphatase activities with sensitivity that equals to that of B/W test..

Quality authorized by:



Jurgita Zilinskiene

PRODUCT USE LIMITATION

This product is developed, designed and sold exclusively *for research purposes and in vitro use only*. The product was not tested for use in diagnostics or for drug development, nor is it suitable for administration to humans or animals.

Please refer to www.thermoscientific.com/onebio for Material Safety Data Sheet of the product.

© 2012 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.