

PRODUCT INFORMATION

PacI

#ER2201 250 U

Lot: ____ **Expiry Date:** __

5'...**T T A A T↓T A A**...3'

3'...**A A T↑T A A T T**...5'

Concentration: 10 U/μL

Source: *E.coli* that carries the cloned *pacIR* gene from *Pseudomonas alcaligenes*

Supplied with: 1 mL of 10X Buffer PacI

Store at -20°C



BSA included

www.thermoscientific.com/onebio

RECOMMENDATIONS

1X Buffer PacI (for 100% PacI digestion)

10 mM Bis-Tris Propane-HCl (pH 6.5), 10 mM MgCl₂,
0.1 mg/mL BSA.

Incubation temperature

37°C.

Unit Definition

One unit is defined as the amount of PacI required to digest 1 μg of control DNA in 1 hour at 37°C in 50 μL of recommended reaction buffer. The control DNA is linearized pJET1 DNA with inserted PacI recognition site.

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.

Storage Buffer

PacI is supplied in: 10 mM potassium phosphate (pH 7.5 at 25°C), 200 mM NaCl, 1 mM DTT, 1 mM EDTA, 0.2 mg/mL BSA and 50% glycerol.

Rev.4



53

Recommended Protocol for Digestion

- Add:

nuclease-free water	16 μ L
10X Buffer Pacl	2 μ L
DNA (0.5-1 μ g/ μ L)	1 μ L
Pacl	0.5-2 μ L
- 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 μ L (~0.1-0.5 μ g of DNA)
nuclease-free water	18 μ L
10X Buffer Pacl	2 μ L
Pacl	1-2 μ L
- Mix gently and spin down for a few seconds.
- Incubate at 37°C for 1-16 hours.

Thermal Inactivation

Pacl is inactivated by incubation at 65°C for 20 min.

ENZYME PROPERTIES

Enzyme Activity in Thermo Scientific REase Buffers, %

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

Methylation Effects on Digestion

Dam: never overlaps – no effect.
Dcm: never overlaps – no effect.
CpG: never overlaps – no effect.
EcoKI: may overlap – effect not determined.
EcoBI: never overlaps – no effect.

Stability during Prolonged Incubation

A minimum of 0.3 unit of the enzyme is required for complete digestion of 1 μ g of 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 μ g of agarose-embedded pJET1 DNA with inserted Pacl recognition sequence in 16 hours.

Compatible Ends

Bsh1285I, BstKTI, SfaNI (AsiSI), PvuI.

Number of Recognition Sites in DNA

Ad2	λ	Φ X174	pBR322	pUC57	pUC18/19	pTZ19R/U	M13mp18/19
1	0	0	0	0	0	0	1

For **CERTIFICATE OF ANALYSIS** see back page

CERTIFICATE OF ANALYSIS

Overdigestion Assay

No detectable change in the specific fragmentation pattern is observed after a 160-fold overdigestion with PacI (10 U/ μ g DNA x 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 PacI 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

This product or portions thereof is manufactured and sold under license under U.S. Patent Nos. 5,874,259 and 6,472,177 and other foreign patents.

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.

