

PRODUCT INFORMATION

Ppu21I (BsaAI)

#ER1971 500 U

Lot: ____ **Expiry Date:** __

5'...**Y A C↓G T R**...3'

3'...**R T G↑C A Y**...5'

Concentration: 10 U/μL

Source: *E.coli* that carries the cloned *ppu21IR* gene from *Pseudomonas putida* RFL21

Supplied with: 1 mL of 10X Buffer Ppu21I

Store at -20°C



BSA included

www.thermoscientific.com/onebio

RECOMMENDATIONS

1X Buffer Ppu21I (for 100% Ppu21I digestion)
10 mM Tris-HCl (pH 7.2 at 37°C), 3 mM MgCl₂,
150 mM NaCl, 0.1 mg/mL BSA.

Incubation Temperature

30°C*.

Unit Definition

One unit is defined as the amount of Ppu21I required to digest 1 μg of lambda DNA in 1 hour at 30°C in 50 μL of 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

Please refer to www.thermoscientific.com/doubledigest to choose the best buffer for your experiments.

Storage Buffer

Ppu21I is supplied in: 10 mM potassium phosphate (pH 7.4 at 25°C), 100 mM NaCl, 1 mM EDTA, 1 mM DTT, 0.2 mg/mL BSA and 50% glycerol.

* Incubation at 37°C results in less than 30% activity.

Recommended Protocol for Digestion

- Add:

nuclease-free water	16 µL
10X Buffer Ppu21I	2 µL
DNA (0.5-1 µg/µL)	1 µL
Ppu21I	0.5-1 µL
- Mix gently and spin down for a few seconds.
- Incubate at 30°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)
Water, nuclease-free	18 µL
10X Buffer Ppu21I	2 µL
Ppu21I	1-2 µL
- Mix gently and spin down for a few seconds.
- Incubate at 30°C for 1-16 hours.

Thermal Inactivation

Ppu21I is inactivated by incubation at 65°C for 20 min.

ENZYME PROPERTIES

Enzyme Activity in Thermo Scientific REase Buffers, %

Ppu21I	B	G	O	R	Tango	2X Tango
100	50-100***	100***	20-50	NR	NR	NR

*** – Star activity appears at a greater than 5-fold overdigestion (5 u x 1h).

NR – Buffer is not recommended, because of high star activity.

Methylation Effects on Digestion

Dam: never overlaps – no effect.

Dcm: never overlaps – no effect.

CpG: completely overlaps – blocked.

EcoKI: may overlap – effect not determined.

EcoBI: may overlap – effect not determined.

Stability during Prolonged Incubation

A minimum of 0.5 units of the enzyme is required for complete digestion of 1 µg of lambda DNA in 16 hours at 30°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
14	2	1	0	0	1	5

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 Ppu21I (10 U/μg lambda 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 Ppu21I 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.

