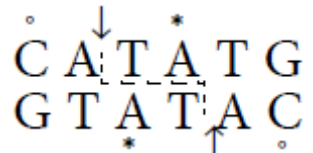


For life science research only.  
Not for use in diagnostic procedures.



# Restriction Endonuclease Nde I from *Neisseria denitrificans* NRCC 31009



 **Version: 17**

Content Version: February 2019

**Cat. No. 11 040 227 001**    1,000 U  
   10 U/μl

**Store product at –15 to –25°C.**

<b>1.</b>	<b>General Information .....</b>	<b>3</b>
1.1.	Contents .....	3
1.2.	Storage and Stability .....	3
	Storage Conditions (Product) .....	3
1.3.	Application .....	3
<b>2.</b>	<b>How to Use this Product .....</b>	<b>4</b>
2.1.	Protocols .....	4
2.2.	Parameters .....	4
	Activity in PCR Buffer .....	4
	Buffers .....	4
	Activity in SuRE/Cut Buffer System .....	4
	Cleavage Sites .....	4
	Number of cleavage sites on different DNAs .....	4
	Compatible Ends .....	4
	Inactivation .....	5
	Isoschizomers .....	5
	Methylation Sensitivity .....	5
	Recognition Sites .....	5
	Temperature Optimum .....	5
	Unit Definition .....	5
<b>3.</b>	<b>Troubleshooting .....</b>	<b>6</b>
<b>4.</b>	<b>Additional Information on this Product .....</b>	<b>7</b>
4.1.	Test Principle .....	7
	Absence of nonspecific endonuclease activities .....	7
	Absence of exonuclease activity .....	7
	Ligation and recutting assay .....	7
	Commonly used bacterial strains .....	7
4.2.	References .....	8
4.3.	Quality Control .....	8
<b>5.</b>	<b>Supplementary Information .....</b>	<b>9</b>
5.1.	Conventions .....	9
5.2.	Changes to previous version .....	9
5.3.	Ordering Information .....	9
5.4.	Trademarks .....	10
5.5.	License Disclaimer .....	10
5.6.	Regulatory Disclaimer .....	10
5.7.	Safety Data Sheet .....	10
5.8.	Contact and Support .....	10

# 1. General Information

## 1.1. Contents

Vial / Bottle	Cap	Label	Function / Description	Content
Nde I	red	Nde I	Contains 20 mM Tris-HCl, 100 mM NaCl, 0.1 mM EDTA, 1 mM dithioerythritol, 0.02% polydocanol, 50% glycerol (v/v), 0.01% gelatin, pH approximately 7.5 (+4°C).	1 vial, 1,000 U (10 U/μl)
H	red	SuRE/Cut Buffer H for Restriction Enzymes, 10x conc.	Contains 500 mM Tris-HCl, 1 M NaCl, 100 mM MgCl <sub>2</sub> , 10 mM dithiothreitol, pH 7.5 (+37°C).	1 vial, 1 ml

## 1.2. Storage and Stability

### Storage Conditions (Product)

The product is shipped on dry ice.

When stored at –15 to –25°C, the product is stable through the expiration date printed on the label.

Vial / Bottle	Cap	Label	Storage
Nde I	red	Nde I	Store at –15 to –25°C. <b>⚠ Do not store below –25°C.</b>
H	red	SuRE/Cut Buffer H, 10x conc.	Store at –15 to –25°C.

## 1.3. Application

Nde I recognizes the sequence CA/TATG and generates fragments with 5'-cohesive termini (Watson RJ, et al, 1982).

## 2. How to Use this Product

### 2.1. Protocols

The following steps describe a typical experiment.

- 1 Prepare the restriction digest according to the following table.

Reagent	Final conc.
DNA	1 µg
10x SuRE/Cut Buffer H	2.5 µl
Water, PCR Grade*	Up to total volume of 25 µl
Nde I	1 U

- 2 Incubate at +37°C for 1 hour.

### 2.2. Parameters

#### Activity in PCR Buffer

0%  
Relative activity in PCR mix (Taq DNA Polymerase buffer) is <5%. The PCR mix contained λ target DNA, primers, 10 mM Tris-HCl (pH 8.3, +20°C), 50 mM KCl, 1.5 mM MgCl<sub>2</sub>, 200 µM dNTPs, 2.5 U Taq DNA polymerase. The mix was subjected to 25 amplification cycles.

#### Buffers

#### Activity in SuRE/Cut Buffer System

A	H <sup>(1)</sup>	M
25 to 50%	100% <sup>(2)</sup>	50 to 75%

<sup>(1)</sup> Supplied Buffer

<sup>(2)</sup> Indicates recommended buffer for optimal activity.

#### Cleavage Sites

#### Number of cleavage sites on different DNAs

λ	Ad2	SV40	ΦX174	M13mp7	pBR322	pBR328	pUC18
7	2	2	0	3	1	0	1

## Compatible Ends

Nde I generates compatible ends to Asn I, Mae I, and Tru 9 I.

Enzyme with compatible ends	Recognition sequence	New sequence if Nde I is ligated to enzyme with compatible ends		Enzyme that can cut this new sequence
		Nde I – Enzyme	Enzyme – Nde I	
Asn I	AT/TAAT	CATAAT	ATTATG	–
Mae I	C/TAG	CATAG	CTATG	–
<b>Nde I</b>	<b>CA/TATG</b>	<b>CA/TATG</b>	<b>CA/TATG</b>	<b>Nde I</b>
Tru 9 I	T/TAA	CATAA	TTATG	–

## Inactivation

Nde I can be heat inactivated by incubation at +65°C for 15 minutes.

## Isoschizomers

The enzyme is not known to have isoschizomers.

## Methylation Sensitivity

Nde I is inhibited by internal 6-methyladenine as indicated (\*), but not by 5-methylcytosine (°).

## Recognition Sites

°CAT\*ATG

**i** \* indicates methylation sensitivity.

## Temperature Optimum

+37°C

## Unit Definition

One unit is the enzyme activity that completely cleaves 1 µg λDNA in 1 hour at +37°C in a total volume of 25 µl SuRE/Cut Buffer H.

## 3. Troubleshooting

Observation	Possible cause	Recommendation
Inhibition or alteration of recognition specificity of restriction enzyme.	Compounds were used in the isolation of the DNA substrate, such as phenol, chloroform, ethanol, SDS, high levels of NaCl, and metal ions, such as Hg <sup>2+</sup> and Mn <sup>2+</sup> .	Remove compounds by ethanol precipitation followed by drying, before adding DNA to the restriction digest reaction. <hr/> Mix vial of restriction enzyme gently but completely prior to use.

## 4. Additional Information on this Product

### 4.1. Test Principle

#### Absence of nonspecific endonuclease activities

1 µg λDNA is incubated for 16 hours in 50 µl SuRE/Cut buffer H with excess of Nde I. The number of enzyme units which do not change the enzyme-specific pattern is stated in the certificate of analysis.

#### Absence of exonuclease activity

Approximately 5 µg [<sup>3</sup>H]-labeled calf thymus DNA are incubated with 3 µl Nde I for 4 hours at +37°C in a total volume of 100 µl 50 mM Tris-HCl, 10 mM MgCl<sub>2</sub>, 1 mM dithioerythritol, pH approximately 7.5. Under these conditions, no release of radioactivity is detectable as stated in the certificate of analysis.

#### Ligation and recutting assay

Nde I fragments obtained by full digestion of 1 µg λDNA are ligated with 1 U T4 DNA Ligase\* in a volume of 10 µl by incubation for 16 hours at +4°C in 66 mM Tris-HCl, 5 mM MgCl<sub>2</sub>, 5 mM Dithiothreitol\*, 1 mM ATP, pH 7.5 (+20°C) resulting in >95 % recovery of 1 µg λDNA fragments. Subsequent recutting with Nde I yields >95% of the typical pattern of Nde I digested λDNA fragments.

#### Commonly used bacterial strains

Strain	Genotype
BL21	<i>E. coli</i> B F <sup>-</sup> <i>dcm ompT hsdS</i> (r <sub>B</sub> - m <sub>B</sub> -) <i>gal</i> (Studier FW, et al, 1986).
C600 <sup>e</sup>	<i>supE44 hsdR2 thi-1 thr-1 leuB6 lacY1 tonA21</i> (Hanahan D, 1983).
DH5α	<i>supE44 Δ(lacU169 (Φ80dlacZΔM15) hsdR17 recA1 endA1 gyrA96 thi-1 relA1</i> (Hanahan D, 1983).
HB101	<i>supE44 hsdS20 recA13 ara-14 proA2 lacY1 galk2 rpsL20 xyl-5 mtl-1</i> (Hanahan D, 1983).
JM108	<i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB)</i> (Yanisch-Perron C, et al, 1985).
JM109	<i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB) F'[traD36proAB<sup>+</sup>, lacI<sup>q</sup> lacZΔM15]</i> (Yanisch-Perron C, et al, 1985).
JM110	<i>rpsL (Str<sup>r</sup>) thr leu thi-1 lacY galk galT ara tonA tsx dam dcm supE44 Δ(lac-proAB) F'[traD36proAB<sup>+</sup>, lacI<sup>q</sup> lacZΔM15]</i> (Yanisch-Perron C, et al, 1985).
K802	<i>supE hsdR gal metB</i> (Raleigh E, et al, 1986; Wood WB, 1966).
SURE <sup>r</sup>	<i>recB recJ sbc C201 uvrC umuC::Tn5(kan<sup>r</sup>) lac, Δ(hsdRMS) endA1 gyrA96 thi relA1 supE44 F'[proAB<sup>+</sup> lacI<sup>q</sup> lacZΔM15 Tn10 (tet<sup>r</sup>)</i> (Greener A, 1990).
TG1	<i>supE hsd Δ5 thi Δ(lac-proAB) F'[traD36proAB<sup>+</sup>, lacI<sup>q</sup> lacZΔM15]</i> (Gibson TJ, 1984).
XL1-Blue <sup>r</sup>	<i>supE44 hsdR17 recA1 endA1 gyrA46 thi relA1 lac F'[proAB<sup>+</sup>, lacI<sup>q</sup> lacZΔM15 Tn10 (tet<sup>r</sup>)</i> (Bullock WO, et al, 1987).

## 4. Additional Information on this Product

### 4.2. References

- Bullock WO, Fernandez JM, Short JM. XL1-Blue– a high-efficiency plasmid transforming recA Escherichia coli strain with  $\beta$ -galactosidase selection. *BioTechniques*. 1987;5:376-379.
- Gibson, TJ. PhD Theses. Cambridge University, U.K 1984.
- Greener, A. *Strategies* 1990;3:5.
- Hanahan D. Studies on transformation of Escherichia coli with plasmids. *J Mol Biol.*1983;166:557-580.
- Raleigh EA, Wilson G. Escherichia coli K-12 restricts DNA containing 5-methylcytosine. *Proc Natl Acad Sci USA.*1986;83:9070-9074.
- Wood WB. Host specificity of DNA produced by Escherichia coli: bacterial mutations affecting the restriction and modification of DNA. *J Mol Biol.*1966;16:118-133.
- Studier FW, Moffatt BA. Use of bacteriophage T7 RNA polymerase to direct selective high-level expression of cloned genes. *J Mol Biol.*1986;189:113-130.
- Watson RJ, Schildraut I, Qiang BQ, Martin SM, Visentin LP. NdeI: a restriction endonuclease from *Neisseria denitrificans* which cleaves DNA at 5'-CATATG-3' sequences. *FEBS Lett.*1982;150:114-116.
- Yanisch-Perron C, Vieira J, Messing J. Improved M13 phage cloning vectors and host strains: nucleotide sequences of the M13mp18 and pUC19 vectors. *Gene.*1985;33:103-19.

### 4.3. Quality Control



For lot-specific certificates of analysis, see section **Contact and Support**.



## 5. Supplementary Information

### 5.1. Conventions

To make information consistent and easier to read, the following text conventions and symbols are used in this document to highlight important information:

Text convention and symbols	
 Information Note: Additional information about the current topic or procedure.	
 <b>Important Note: Information critical to the success of the current procedure or use of the product.</b>	
① ② ③ etc.	Stages in a process that usually occur in the order listed.
① ② ③ etc.	Steps in a procedure that must be performed in the order listed.
* (Asterisk)	The Asterisk denotes a product available from Roche Diagnostics.

### 5.2. Changes to previous version

Layout changes.  
Editorial changes.

### 5.3. Ordering Information

Product	Pack Size	Cat. No.
Reagents, kits		
1,4-Dithiothreitol	2 g	10 197 777 001
	10 g	10 708 984 001
	25 g	11 583 786 001
T4 DNA Ligase	100 U, 1 U/μl	10 481 220 001
	500 U, 1 U/μl	10 716 359 001
	500 U, 5 U/μl	10 799 009 001
Water, PCR Grade	25 ml, 25 x 1 ml	03 315 932 001
	25 ml, 1 x 25 ml	03 315 959 001
	100 ml, 4 x 25 ml	03 315 843 001
SuRE/Cut Buffer A	5 x 1 ml	11 417 959 001
SuRE/Cut Buffer M	5 x 1 ml	11 417 983 001
SuRE/Cut Buffer H	5 x 1 ml	11 417 991 001

## 5. Supplementary Information

### 5.4. Trademarks

SURE/CUT is a trademark of Roche.

All other product names and trademarks are the property of their respective owners.

### 5.5. License Disclaimer

For patent license limitations for individual products please refer to:

**List of biochemical reagent products.**

### 5.6. Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

### 5.7. Safety Data Sheet

Please follow the instructions in the Safety Data Sheet (SDS).

### 5.8. Contact and Support

To ask questions, solve problems, suggest enhancements or report new applications, please visit our

**Online Technical Support Site.**

To call, write, fax, or email us, visit **sigma-aldrich.com**, and select your home country. Country-specific contact information will be displayed.

