

Restriction Endonuclease Nde II (Mbo I)

From *Neisseria denitrificans* NRCC 31009

Cat. No. 11 040 243 001 1000 units (5 U/μl)



Version 20

Content version: March 2016

Store at -15 to -25°C

Stability/Storage The undiluted enzyme solution is stable when stored at -15 to -25°C until the control date printed on the label. Do not store below -25°C to avoid freezing.

Sequence Specificity *Nde* II recognizes the sequence /GATC and generates fragments with 5'-cohesive termini (1).

Compatible ends *Nde* II generates compatible ends to *Bam* HI, *Bcl* I, *Bgl* II and *Xho* II.

Isoschizomers *Nde* II is an isoschizomer to *Mbo* I, *Sau* 3A and *Dpn* I.

Methylation sensitivity *Nde* II is inhibited by overlapping dam-methylation (*), in contrast to its isoschizomer *Sau* 3A which is not influenced by dam-methylation. See also *Dpn* I. *Nde* II is not inhibited by 5-methylcytosine.

Storage buffer 20 mM Tris-HCl, 50 mM NaCl, 0.1 mM EDTA, 1 mM dithioerythritol, 0.02% polydocanol, 0.01% gelatine, 50% glycerol (v/v), pH approx. 7.5 (at 4°C).

Supplied incubation buffer (2× conc.) 200 mM Tris-HCl, 300 mM NaCl, 20 mM MgCl₂, 2 mM dithioerythritol, pH 7.6 (at 37°C).

Activity in SuRE/Cut Buffer System

| | A | B | L | M | H |
|--|--------|--------|-------|-------|--------|
| | 10-25% | 10-25% | 0-10% | 0-10% | 10-25% |

Incubation temp. 37°C

Unit definition One unit is the enzyme activity that completely cleaves 1 μg λ dam⁻, dcm⁻ DNA in 1 h at 37°C in a total volume of 25 μl incubation buffer.

Typical experiment

| Component | Final concentration |
|-----------------------------|-------------------------------|
| DNA | 1 μg/l |
| 2× Suppl. incubation buffer | 12.5 μl |
| Repurified water | Up to a total volume of 25 μl |
| Restriction enzyme | 1 unit |

Incubate at 37°C for 1 h.

Heat inactivation There is no information about *Nde* II and heat-inactivation available.

Number of cleavage sites on different DNA's (2):

| | λ | Ad2 | SV40 | Φ X174 | M13mp7 | pBR322 | pBR328 | pUC18 |
|--|-----|-----|------|--------|--------|--------|--------|-------|
| | 116 | 87 | 8 | 0 | 8 | 22 | 27 | 15 |

Ligation and recutting assay *Nde* II fragments obtained by complete digestion of 1 μg λ dam⁻ DNA are ligated with 1 U T4-DNA ligase (Cat. No. 10 481 220 001) in a volume of 10 μl by incubation for 16 h at 4°C in 66 mM Tris-HCl, 5 mM MgCl₂, 5 mM dithioerythritol, 1 mM ATP, pH 7.5 (at 20°C) resulting in >95% recovery of 1 μg λdam⁻ DNA fragments. Subsequent re-cutting with *Nde* II yields >95% of the typical pattern of λdam⁻ DNA × *Nde* II fragments

Troubleshooting A critical component is the DNA substrate. Many compounds used in the isolation of DNA such as phenol, chloroform, EtOH, SDS, high levels of NaCl, metals (e.g. Hg²⁺, Mn²⁺) inhibit or alter recognition specificity of many restriction enzymes. Such compounds should be removed by EtOH precipitation followed by drying, before the DNA is added to the restriction digest reaction. Appropriate mixing of the enzyme is recommended.

Quality control

Lot-specific certificates of analysis are available at www.lifescience.roche.com/certificates

Absence of unspecific endonuclease activities 1 μg λ dam⁻, dcm⁻ or pBR322 DNA is incubated for 16 h in 50 μl incubation buffer with excess of *Nde* II. The number of enzyme units which do not change the enzyme-specific pattern is stated under "Endo" printed on the label.

Absence of exonuclease activity Approx. 5 μg [³H] labeled calf thymus DNA are incubated with 3 μl *Nde* II for 4 h at 37°C in a total volume of 100 μl 50 mM Tris-HCl, 10 mM MgCl₂, 1 mM dithioerythritol, pH approx. 7.5. Under these conditions, no release of radioactivity is detectable, as stated in the certificate of analysis

References

- 1 Watson, R. J. et al. (1982) *FEBS Lett.* **150**, 114-116.
- 2 Kessler, C. & Manta, V. (1990) *Gene* **92**, 1-250.
- 3 Rebase The Restriction Enzyme Database: <http://rebase.neb.com>

Ordering Information

| Product | Application | Packsizes | Cat. No. |
|--|--|---|--|
| Rapid DNA Ligation Kit | Ligation of sticky- or blunt-ended DNA fragments in just 5 min at +15 to +25 °C. | Kit (40 DNA ligations) | 11 635 379 001 |
| T4 DNA Ligase | Ligation of sticky- and blunt- ended DNA fragments. | 100 U 500 units (1 U/μl) | 10 481 220 001 10 716 359 001 |
| rAPid Phosphatase | Dephosphorylation of 5'-phosphate residues from nucleic acids | 1000 U 5000 U | 04 898 133 001 04 898 141 001 |
| rAPid Dephos and Ligation Kit | Dephosphorylation of nucleic acids. | 40 reactions 160 reactions | 04 898 117 001 04 898 125 001 |
| Alkaline Phosphatase (AP), special quality for molecular biology | Dephosphorylation of 5'-phosphate residues from nucleic acids. | 1000 U (20 U/μl) | 11 097 075 001 |
| Agarose MP | Multipurpose agarose for analytical and preparative electrophoresis of nucleic acids | 100 g 500 g | 11 388 983 001 11 388 991 001 |
| Agarose LE | Separation of nucleic acids in the range 0.2 - 1.5 kbp | 100 g 500 g | 11 685 660 001 11 685 678 001 |
| Agarose Gel DNA Extraction Kit | For the elution of DNA fragments from agarose gels. | 1 Kit (max. 100 reactions) | 11 696 505 001 |
| High Pure PCR Product Purification Kit | Purification of PCR or enzymatic modification reaction (e.g. restriction digest) | 50 purifications 250 purifications | 11 732 668 001 11 732 676 001 |
| SuRE/Cut Buffer Set for Restriction Enzymes | Incubation buffers A, B, L, M and H for restriction enzymes | 1 ml each (10× conc. solutions) | 11 082 035 001 |
| SuRE/Cut Buffer A | Restriction enzyme incubation | 5 × 1 ml (10× conc. solution) | 11 417 959 001 |
| SuRE/Cut Buffer B | Restriction enzyme incubation | 5 × 1 ml (10× conc. solution) | 11 417 967 001 |
| SuRE/Cut Buffer H | Restriction enzyme incubation | 5 × 1 ml (10× conc. solution) | 11 417 991 001 |
| SuRE/Cut Buffer L | Restriction enzyme incubation | 5 × 1 ml (10× conc. solution) | 11 417 975 001 |
| SuRE/Cut Buffer M | Restriction enzyme incubation | 5 × 1 ml (10× conc. solution) | 11 417 983 001 |
| Water, PCR Grade | Specially purified, double-distilled, deionized, and autoclaved | 100 ml (4 vials of 25 ml) 25 ml (25 vials of 1 ml) 25 ml (1 vial of 25 ml) | 03 315 843 001 03 315 932 001 03 315 959 001 |
| BSA, special quality for molecular biology | Maintaining enzyme stability | 20 mg (1 ml) | 10 711 454 001 |

Changes to previous version

- Update of quality control
- Editorial changes.

Trademarks

HIGH PURE and SURE/CUT are trademarks of Roche. All other product names and trademarks are the property of their respective owners.

Regulatory Disclaimer

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

Commonly used bacterial strains

| Strain | Genotype |
|-----------------------|--|
| BL21 | <i>E. coli</i> B F ⁻ <i>dcm ompT hsdS(r_B- m_B-) gal</i> (Studier, F.W. <i>et al</i> (1986) <i>J. Mol. Biol.</i> , 189 , 113.) |
| C600 ^e | <i>supE44 hsdR2 thi-1 thr-1 leuB6 lacY1 tonA21</i> ; (Hanahan, D. (1983) <i>J. Mol. Biol.</i> 166 , 557) |
| DH5α | <i>supE44 Δ(lacU169 (φ80d/lacZΔM15) hsdR17 recA1 endA1 gyrA96 thi-1 relA1</i> ; (Hanahan, D. (1983) <i>J. Mol. Biol.</i> 166 , 557) |
| HB101 | <i>supE44 hsdS20 recA13 ara-14 proA2 lacY1 galK2 rpsL20 xyl-5 mtl-1</i> ; (Hanahan, D., (1983) <i>J. Mol. Biol.</i> 166 , 557.) |
| JM108 | <i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB)</i> ; (Yanisch- Perron, C. <i>et al.</i> , (1985) <i>Gene</i> 33 , 103.) |
| JM109 | <i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB) F[traD36proAB⁺, lac^q lacZΔM15]</i> ; (Yanisch- Perron, C. <i>et al.</i> , (1985) <i>Gene</i> 33 , 103.) |
| JM110 | <i>rpsL (Str^r) thr leu thi-1 lacY galK galT ara tonA tsx dam dcm supE44 Δ(lac-proAB) F[traD36proAB⁺, lac^q lacZΔM15]</i> ; (Yanisch- Perron, C. <i>et al.</i> , (1985) <i>Gene</i> 33 , 103.) |
| K802 | <i>supE hsdR gal metB</i> ; (Raleigh, E. <i>et al.</i> , (1986) <i>Proc.Natl. Acad.Sci USA</i> , 83 , 9070.; Wood, W.B. (1966) <i>J. Mol. Biol.</i> , 16 , 118.) |
| SURE ^f | <i>recB recJ sbc C201 uvrC umuC::Tn5(kan^r) lac</i> , Δ(<i>hsdRMS</i>) <i>endA1 gyrA96 thi relA1 supE44 F[proAB⁺ lac^q lacZΔM15 Tn10 (tet^r)</i> ; (Greener, A. (1990) <i>Stratagies</i> , 3 , 5.) |
| TG1 | <i>supE hsd Δ5 thi Δ(lac-proAB) F[traD36proAB⁺, lac^q lacZΔM15]</i> ; (Gibson, T.J. (1984) <i>PhD Theses. Cambridge University, U.K.</i>) |
| XL1-Blue ^f | <i>supE44 hsdR17 recA1 endA1 gyrA46 thi relA1 lac F[proAB⁺, lac^q lacZΔM15 Tn10 (tet^r)</i> ; (Bullock <i>et al.</i> , (1987) <i>BioTechniques</i> , 5 , 376.) |

Contact and Support

To ask questions, solve problems, suggest enhancements and 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.



Roche Diagnostics GmbH
Sandhofer Strasse 116
68305 Mannheim
Germany