

# Restriction Endonuclease *Nhe* I

From *Neisseria mucosa*

**Cat. No. 10 885 843 001** 200 units (10 U/μl)  
**Cat. No. 10 885 851 001** 1,000 units (10 U/μl)  
**Cat. No. 10 885 860 001** 1,500 units, high concentration (40 U/μl)



**Version 20**  
 Content version: November 2012  
 Store at -15 to -25°C

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

**Sequence specificity** *Nhe* I recognizes the sequence G/CTAGC and generates fragments with 5'-cohesive termini (1, 2).

**Compatible ends** *Nhe* I generates compatible ends to *Avr* II, *Spe* I, and *Xba* I.

| Enzyme with compatible ends | Recognition sequence | New sequence if <i>Nhe</i> I is ligated to enzyme with compatible ends |                       | Enzyme that can cut this new sequence |
|-----------------------------|----------------------|--|-----------------------|---------------------------------------|
|                             |                      | <i>Nhe</i> I - Enzyme  | Enzyme - <i>Nhe</i> I |                                       |
| <b><i>Avr</i> II</b>        | C/CTAGG              | GC/TAGG  | CC/TAGC               | <i>Bfa</i> I, <i>Mae</i> I            |
| <b><i>Nhe</i> I</b>         | <b>G/CTAGC</b>       | <b>G/CTAGC</b>   | <b>G/CTAGC</b>        | <b><i>Nhe</i> I</b> + Iso-schizomers  |
| <b><i>Spe</i> I</b>         | A/CTAGT              | GC/TAGT  | AC/TAGC               | <i>Bfa</i> I, <i>Mae</i> I            |
| <b><i>Xba</i> I</b>         | T/CTAGA              | GC/TAGA  | TC/TAGC               | <i>Bfa</i> I, <i>Mae</i> I            |

**Isoschizomers** The enzyme is not known to have isoschizomers.

**Methylation sensitivity** *Nhe* I is inhibited by the presence of 5'-methylcytosine in the sequence G/CTAGC.

**Storage buffer** 20 mM Tris-HCl, 100 mM NaCl, 0.1 mM EDTA, 10 mM 2-Mercaptoethanol, 0.2% Polydocanol, 0.01% Gelatine, 50% Glycerol (v/v), pH approx. 7.8 (at 4°C).

**Suppl. Incubation buffer (10x)** 100 mM Tris-HCl, 500 mM NaCl, 100 mM MgCl<sub>2</sub>, 10 mM Dithioerythritol, pH 7.5 (at 37°C), (Δ SuRE/Cut Buffer **M**).

**Activity in SuRE/Cut Buffer System** Bold face printed buffer indicates the recommended buffer for optimal activity:

| A    | B      | L    | <b>M</b>    | H      |
|------|--------|------|-------------|--------|
| 100% | 25-50% | 100% | <b>100%</b> | 10-25% |

**Incubation temperature** **37°C**

**Unit definition** One unit is the enzyme activity that completely cleaves 1 μg λ x *Eco* RI DNA in 1 h at 37°C in a total volume of 25 μl SuRE/Cut buffer **M**. 1 μg pBR322 DNA is digested completely by approx. 8 units of *Nhe* I because of the larger number of cleavage sites per μg of pBR322 DNA as compared to λDNA.

**Typical experiment**

| Component                    | Final concentration           |
|------------------------------|-------------------------------|
| DNA                          | 1 μg                          |
| 10x SuRE/Cut Buffer <b>M</b> | 2.5 μl                        |
| Sterile double-dist. water   | Up to a total volume of 25 μl |
| Restriction enzyme           | 1 unit                        |

Incubate at 37°C for 1 h.

**Heat Inactivation** *Nhe* I can be heat inactivated by heating to 65°C for 15 min.

**Number of cleavage sites on different DNAs (2):**

| λ | Ad2 | SV40 | Φ X174 | M13mp7 | pBR322 | pBR328 | pUC18 |
|---|-----|------|--------|--------|--------|--------|-------|
| 1 | 4   | 0    | 0      | 0      | 1      | 1      | 0     |

**Activity in PCR buffer** Relative activity in PCR mix (Taq DNA Polymerase buffer) is **100%**. 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.

**Ligation and recutting assay** λ x *Eco* RI/*Nhe* I fragments obtained by complete digestion of 1 μg λ x *Eco* RI/*Nhe* I DNA are ligated with 1 U T4 DNA Ligase in a volume of 10 μl by incubation for 16 h at 4°C in 66 mM Tris-HCl, 5 mM MgCl<sub>2</sub>, 5 mM dithiothreitol, 1 mM ATP, pH 7.5 (at 20°C) resulting in >90% recovery of 1 μg λDNA x *Eco* RI/*Nhe* I fragments. Subsequent re-cutting with *Eco* RI/*Nhe* I yields > 80% of the typical pattern of λDNA x *Eco* RI/*Nhe* I 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., Hg<sup>2+</sup>, Mn<sup>2+</sup>), 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.roche-applied-science.com/certificates](http://www.roche-applied-science.com/certificates).

**Absence of unspecific endonuclease activities** 1 μg λ x *Eco* RI DNA is incubated for 16 h in 50 μl SuRE/Cut buffer **M** with excess of *Nhe* 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** Approx. 5 μg [<sup>3</sup>H] labeled calf thymus DNA are incubated with 3 μl *Nhe* I for 4 h at 37°C in a total volume of 100 μl 50 mM Tris-HCl, 10 mM MgCl<sub>2</sub>, 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 Comb, D. G. & Schildkraut, I., unpublished observations
- 2 Kessler, C. & Manta, V. (1990) *Gene* **92**, 1-250.
- 3 Rebase The Restriction Enzyme Database: <http://rebase.neb.com>
- 4 Benchmate: <http://www.roche-applied-science.com/benchmate>

## Ordering Information

Roche Applied Science offers a large selection of reagents and systems for life science research. For a complete overview of related products and manuals, please visit and bookmark our homepage, [www.rocche-applied-science.com](http://www.rocche-applied-science.com), and our Special Interest Sites, including "Mapping & Cloning": <http://www.restriction-enzymes.com>.

The convenient RE Finder Program located on our Bench Mate website, <http://www.rocche-applied-science.com/benchmate> helps you identify the enzymes that will cut your DNA sequence, and displays the names and recognition sequences of enzymes and isoschizomers as well as links to detailed information (e.g. Instructions for Use) of the selected restriction enzyme.

| Product  | Application  | Pack Size   | Cat. No.   |
|--|--|---|--|
| Restriction Enzymes  | DNA restriction digestion.   | Please refer to website or catalog  |  |
| T4 DNA Ligase  | Ligation of sticky- and blunt-ended DNA fragments.   | 100 U<br>500 U (1 U/μl)<br>500 U (5 U/μl)   | 10 481 220 001<br>10 716 359 001<br>10 799 009 001 |
| Rapid DNA Dephos & Ligation Kit                                  | Upgrade from the Rapid DNA Ligation Kit for fast and efficient DNA dephosphorylation and ligation of sticky- or blunt-ended DNA fragments. | 40 reactions<br>160 reactions   | 04 898 117 001<br>04 898 125 001                   |
| rAPid Alkaline Phosphatase                                       | Dephosphorylation of 5'-phosphate residues from nucleic acids.   | 1,000 U<br>5,000 U  | 04 898 133 001<br>04 898 141 001                   |
| Rapid DNA Ligation Kit   | Ligation of nucleic acids.   | Kit<br>(40 DNA ligations)   | 11 635 379 001                                     |
| Alkaline Phosphatase (AP), special quality for molecular biology | Dephosphorylation of 5'-phosphate residues from nucleic acids.   | 1,000 U<br>(20 U/μl)  | 11 097 075 001                                     |
| Agarose MP   | Multipurpose agarose for analytical and preparative electrophoresis of nucleic acids.  | 100 g<br>500 g  | 11 388 983 001<br>11 388 991 001                   |
| Agarose LE   | Separation of nucleic acids in the range 0.2 - 1.5 kbp.  | 100 g<br>500 g  | 11 685 660 001<br>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<br>250 purifications   | 11 732 668 001<br>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<br>(4 vials of 25 ml)<br>25 ml<br>(25 vials of 1 ml)<br>25 ml<br>(1 vial of 25 ml) | 03 315 843 001<br>03 315 932 001<br>03 315 959 001 |
| BSA, special quality for molecular biology                       | Maintaining enzyme stability.  | 20 mg (1 ml)  | 10 711 454 001                                     |

## Printed Materials

You can view the following manuals on our website:

|  |
|--|
| Lab FAQs "Find a Quick Solution"           |
| Restriction Enzyme Ordering Guide          |
| Molecular Weight Markers for Nucleic Acids |

## Changes to previous version

Star activity information removed.

## 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 <sup>-</sup> <i>dcm ompT hsdS(r<sub>B</sub><sup>-</sup> m<sub>B</sub><sup>-</sup>) gal</i> (Studier, F.W. <i>et al.</i> (1986) <i>J. Mol. Biol.</i> , <b>189</b> , 113.)  |
| C600 <sup>e</sup>     | <i>supE44 hsdR2 thi-1 thr-1 leuB6 lacY1 tonA21</i> ; (Hanahan, D. (1983) <i>J. Mol. Biol.</i> <b>166</b> , 557.)   |
| DH5α                  | <i>supE44 Δ(lacU169 (φ80d/lacZΔM15) hsdR17 recA1 endA1 gyrA96 thi-1 relA1</i> ; (Hanahan, D. (1983) <i>J. Mol. Biol.</i> <b>166</b> , 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> <b>166</b> , 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> <b>33</b> , 103.)  |
| 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. <i>et al.</i> , (1985) <i>Gene</i> <b>33</b> , 103.)  |
| 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. <i>et al.</i> , (1985) <i>Gene</i> <b>33</b> , 103.)             |
| K802                  | <i>supE hsdR gal metB</i> ; (Raleigh, E. <i>et al.</i> , (1986) <i>Proc. Natl. Acad. Sci. USA</i> , <b>83</b> , 9070.; Wood, W.B. (1966) <i>J. Mol. Biol.</i> , <b>16</b> , 118.)  |
| SURE <sup>f</sup>     | <i>recB recJ sbc C201 uvrC umuC::Tn5(kan<sup>r</sup>) lac<sup>+</sup>, 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) <i>Strategies</i> , <b>3</b> , 5.) |
| TG1                   | <i>supE hsd Δ5 thi Δ(lac-proAB) F[traD36proAB<sup>+</sup>, lacI<sup>q</sup> lacZΔM15]</i> ; (Gibson, T.J. (1984) PhD Theses. Cambridge University, U.K.)   |
| XL1-Blue <sup>f</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 <i>et al.</i> , (1987) <i>BioTechniques</i> , <b>5</b> , 376.)  |

## Contact and Support

To ask questions, solve problems, suggest enhancements or report new applications, please visit our **Online Technical Support Site** at:

**[www.rocche-applied-science.com/support](http://www.rocche-applied-science.com/support)**

To call, write, fax, or email us, visit the Roche Applied Science home page, [www.rocche-applied-science.com](http://www.rocche-applied-science.com), and select your home country. Country-specific contact information will be displayed. Use the Product Search function to find Pack Inserts and Material Safety Data Sheets.



Roche Diagnostics GmbH  
Roche Applied Science  
68298 Mannheim  
Germany