

# Restriction Endonuclease Nco I

From *Rhodococcus species* (formerly *Nocardia corallina*)

**Cat. No. 10 835 315 001** 200 units (10 U/μl)  
**Cat. No. 10 835 323 001** 1000 units (10 U/μl)  
**Cat. No. 11 047 698 001** 1000 units, high concentration (40 U/μl)



**Version 20**

Content version: April 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** *Nco* I recognizes the sequence C/CATGG and generates fragments with 5'-cohesive termini (1).

**Compatible ends** *Nco* I generates compatible ends to *Afl* III, *Bsp* LU 111, *Dsa* I, *Rca* I.

Enzyme with compatible ends	Recognition sequence	New sequence if <i>Nco</i> I is ligated to enzyme with compatible ends		Enzyme that can cut this new sequence
		<i>Nco</i> I - Enzyme	Enzyme - <i>Nco</i> I	
<i>Afl</i> III	A/C(A,G)(T,C)GT	C/C(A,G)(T,C)GT	A/CATGG	<i>Mae</i> II, <i>Mvu</i> I, <i>Nla</i> III
<i>Bsp</i> LU 111	A/CATGT	C/CATGT	A/CATGG	<i>Nla</i> III
<i>Dsa</i> I	C/C(A,G)(C,T)GG	C/C(A,G)(C,T)GG	C/CATGG	<i>Dsa</i> I, <i>Ksp</i> I, <i>Mvn</i> I, <i>Nco</i> I, <i>Nla</i> III, <i>Sty</i> I
<b><i>Nco</i> I</b>	<b>C/CATGG</b>	<b>C/CATGG</b>	<b>C/CATGG</b>	<b><i>Nco</i> I</b> , <i>Dsa</i> I, <i>Sty</i> I
<i>Rca</i> I	T/CATGA	C/CATGA	T/CATGG	<i>Nla</i> III

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

**Methylation sensitivity** *Nco* I is inhibited by the presence of 5-methylcytosine as indicated (\*). Presence of 6-methyladenine does not influence the cleavage (°).

**Storage buffer** 20 mM Tris-HCl, 100 mM NaCl, 0.1 mM EDTA, 10 mM 2-Mercaptoethanol, 0.2% Triton<sup>®</sup> X-100 (v/v), 50% Glycerol (v/v), pH approx. 8.0 (at 4°C).

**Incubation buffer, 10x** 500 mM Tris-HCl, 1 M NaCl, 100 mM MgCl<sub>2</sub>, 10 mM Dithioerythritol, pH 7.5 (at 37°C), (= SuRE/Cut Buffer **H**)

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

A	B	L	M	<b>H</b>
50-75%	50-75%	50-75%	50-75%	<b>100%</b>

**Incubation temperature** **37°C**

**Unit definition** One unit is the enzyme activity that completely cleaves 1 μg λDNA in 1 h at **37°C** in a total volume of 25 μl in SuRE/Cut buffer **H**

**Typical experiment**

Component	Final concentration
DNA	1 μg
10 × SuRE/Cut Buffer <b>H</b>	2.5 μl
Repurified water	Up to a total volume of 25 μl
Restriction enzyme	1 unit

Incubate at **37°C** for 1 h.

**Heat inactivation** Inactivated by 15 min incubation at 65°C (tested up to 100 U/μg DNA).

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

λ	Ad2	SV40	Φ X174	M13mp7	pBR322	pBR328	pUC18
4	20	3	0	0	0	1	0

**Activity in PCR buffer** Relative activity in PCR mix (Taq DNA Polymerase buffer) is 50%. 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** *Nco* I fragments obtained by complete digestion of 1 μg λDNA are ligated with 0.1 units 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>, 1 mM Dithioerythritol, 1 mM ATP, pH 7.5 (at 20°C) resulting in >95% recovery of 1 μg λDNA fragments. Subsequent re-cutting with *Nco* I yields >95% of the typical pattern of λDNA × *Nco* I fragments.

**Troubleshooting** A critical component is the DNA substrate. Many compounds used in the isolation of DNA such as phenol, chloroform, ethanol, SDS, high levels of NaCl, metal ions (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 ethanol 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 λDNA is incubated for 16 h in 50 μl SuRE/Cut buffer H with excess of *Nco* 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 *Nco* 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**

- Langdale, J. A., Myers, P. A., & Roberts, R. J., unpublished observations.
- Kessler, C. & Manta, V. (1990) *Gene* **92**, 1-248.
- Rebase The Restriction Enzyme Database: <http://rebase.neb.com>
- Benchmark: <http://www.roche-applied-science.com/benchmark/>

## 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

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.

### Disclaimer of License

For patent license limitations for individual products please refer to: [List of biochemical reagent products](#)

## 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, 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>, lac<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>, lac<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</i> , Δ( <i>hsdRMS</i> ) <i>endA1 gyrA96 thi relA1 supE44 F[proAB<sup>+</sup> lac<sup>q</sup> lacZΔM15 Tn10 (tet<sup>r</sup>)</i> ; (Greener, A. (1990) <i>Stratagies</i> , <b>3</b> , 5.)
TG1	<i>supE hsd Δ5 thi Δ(lac-proAB) F[traD36proAB<sup>+</sup>, lac<sup>q</sup> lacZΔM15]</i> ; (Gibson, T.J. (1984) <i>PhD Theses</i> . Cambridge University, U.K.)
XL1-Blue <sup>f</sup>	<i>supE44 hsdR17 recA1 endA1 gyrA46 thi relA1 lac F[proAB<sup>+</sup>, lac<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 and report new applications, please visit our [Online Technical Support Site](#).

To call, write, fax, or email us, visit [sigma-aldrich.com](http://sigma-aldrich.com), and select your home country. Country-specific contact information will be displayed.



Roche Diagnostics GmbH  
Sandhofer Strasse 116  
68305 Mannheim  
Germany