

Restriction Endonuclease *Eco* 47 III

From *Escherichia coli* RFL 47

Cat. No. 11 167 103 001 100 units (5 U/ μ l)



Version 16

Content version: July 2017

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	<i>Eco</i> 47 III recognizes the sequence AGC/GCT and generates fragments blunt ends (1).										
Compatible ends	The enzyme generates compatible ends to any blunt end.										
Isoschizomers	The enzyme is not known to have isoschizomers.										
Methylation sensitivity	<i>Eco</i> 47 III is inhibited by the presence of 5-methylcytosine as indicated (*). 6-methyladenine is not inhibiting, as indicated (°).										
Storage buffer	10 mM Tris-HCl, 100 mM KCl, 0.1 mM EDTA, 10 mM dithiothreitol, 200 $\mu\text{g}/\text{ml}$ bovine serum albumin, 50% glycerol (v/v), pH approx. 8.0 (at 4°C).										
Incubation buffer, (10x)	500 mM Tris-HCl, 1 M NaCl, 100 mM MgCl_2 , 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:										
	<table border="1"> <thead> <tr> <th>A</th> <th>B</th> <th>L</th> <th>M</th> <th>H</th> </tr> </thead> <tbody> <tr> <td>25-50%</td> <td>50-75%</td> <td>0-10%</td> <td>25-50%</td> <td>100%</td> </tr> </tbody> </table>	A	B	L	M	H	25-50%	50-75%	0-10%	25-50%	100%
A	B	L	M	H							
25-50%	50-75%	0-10%	25-50%	100%							
Incubation temperature	37°C										
Unit definition	One unit is the enzyme activity that completely cleaves 1 μg λ DNA in 1 h at 37°C in the SuRE/Cut buffer H in a total volume of 25 μl . 1 μg pBR322 is digested completely by 4 units of <i>Eco</i> 47 III.										
Typical experiment	<table border="1"> <thead> <tr> <th>Component</th> <th>Final concentration</th> </tr> </thead> <tbody> <tr> <td>DNA</td> <td>1 μg</td> </tr> <tr> <td>10 \times SuRE/Cut Buffer H</td> <td>2.5 μl</td> </tr> <tr> <td>Repurified water</td> <td>Up to a total volume of 25 μl</td> </tr> <tr> <td>Restriction enzyme</td> <td>1 unit</td> </tr> </tbody> </table> <p>Incubate at 37°C for 1 h.</p>	Component	Final concentration	DNA	1 μg	10 \times SuRE/Cut Buffer H	2.5 μl	Repurified water	Up to a total volume of 25 μl	Restriction enzyme	1 unit
Component	Final concentration										
DNA	1 μg										
10 \times SuRE/Cut Buffer H	2.5 μl										
Repurified water	Up to a total volume of 25 μl										
Restriction enzyme	1 unit										
Heat inactivation	The enzyme can be heat inactivated by 15 min incubation at 65°C (tested up to 100 units/ μg DNA).										

Number of cleavage sites on different DNAs (2):

λ	Ad2	SV40	Φ X174	M13mp7	pBR322	pBR328	pUC18
2	13	1	0	2	4	3	0

PFGE tested

Eco 47 III has been tested in Pulsed-Field-Gel Electrophoresis (test system bacterial chromosomes). For cleavage of genomic DNA (*E.coli* C600) embedded in agarose for PFGE analysis 10 units of enzyme/ μg DNA and 4 h incubation time are recommended.

Activity in PCR buffer

Relative activity in PCR mix (Taq DNA Polymerase buffer) is **0%**. The PCR mix contained λ target DNA, primers, 10 mM Tris-HCl (pH 8.3, 20°C), 50 mM KCl, 1.5 mM MgCl_2 , 200 μM dNTPs, 2.5 U Taq DNA polymerase. The mix was subjected to 25 amplification cycles.

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^{2+} , Mn^{2+}) 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.lifescience.roche.com/certificates.

Absence of unspecific endonuclease activities

1 μg λ DNA is incubated for 16 h in 50 μl SuRE/Cut buffer H with excess of *Eco* 47 III. 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 [^3H] labeled calf thymus DNA are incubated with 3 μl *Eco* 47 III for 4 h at 37°C in a total volume of 100 μl 50 mM Tris-HCl, 10 mM MgCl_2 , 1 mM Dithioerythritol, pH approx. 7.5. The release of radioactivity is calculated as a percentage value of liberated to input radioactivity per unit of enzyme (stated in the certificate of analysis).

Ligation and recutting assay

Eco 47 III fragments obtained by complete digestion of 1 μg pBR322 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 MgCl_2 , 5 mM Dithiothreitol, 1 mM ATP, pH 7.5 (at 20°C).

The percentage of ligation and subsequent recutting with *Eco* 47 III which yields the typical pattern of pBR322 \times *Eco* 47 III fragments are determined and stated in the certificate of analysis.

References

- Janulaitis, A., Petrusytė, M. & Butkus, V. (1983) *FEBS Lett.* **161**, 213-216.
- Kessler, C. & Manta, V. (1980) *Gene* **92**, 1-248.
- Sagawa, H. et al. (1992) *Nucl. Acids Res.* **20**, 365;
- Rebase The Restriction Enzyme Database: <http://rebase.neb.com>

Ordering Information

Product	Application	Packsizes	Cat. No.
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
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)	03 315 843 001
		25 ml (25 vials of 1 ml)	03 315 932 001
		25 ml (1 vial of 25 ml)	03 315 959 001

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 ⁻ <i>dcm ompT hsdS</i> (r _B - m _B -) <i>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