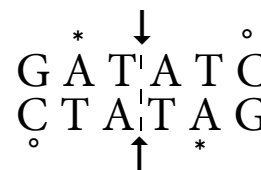


Restriction Endonuclease *EcoR V*

From *Escherichia coli* J62 pLG74

Cat. No. 10 667 145 001 2000 units (10 U/μl)
Cat. No. 10 667 153 001 10 000 units (10 U/μl)
Cat. No. 11 040 197 001 10 000 units, high concentration (40 U/μl)



Version 19

Content version: June 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 *EcoR V* recognizes the sequence GAT/ATC and generates fragments with blunt ends (1).

Compatible ends *EcoR V* generates fragments which are compatible to any blunt end.

Isoschizomers *EcoR V* is an isoschizomer of *Eco 32I*.

Methylation sensitivity *EcoR V* is inhibited by the presence of N6-Methyladenine, as indicated (*). *EcoR V* is not inhibited by the presence of 5-Methylcytosine (°) or 5-Hydroxymethylcytosine.

Storage buffer 20 mM Tris-HCl, 100 mM NaCl, 1 mM EDTA, 10 mM 2-Mercaptoethanol, 0.02% Polydocanol, 50% Glycerol (v/v), pH approx. 7.5 (at 4°C)

Suppl Incubation buffer (10x) 100 mM Tris-HCl, 1 M NaCl, 50 mM MgCl₂, 10 mM 2-Mercaptoethanol, pH 8.0 (at 37°C), (Δ SuRE/Cut Buffer **B**)

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

A	B	L	M	H
25-50%	100%	0-10%	25-50%	50-75%

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 SuRE/Cut buffer **B**.

Typical experiment

Component	Final concentration
DNA	1 μg
10 × SuRE/Cut Buffer 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 *EcoR V* is not heat-inactivated by 15 min incubation at 65°C.

Number of cleavage sites on different DNAs (2):

λ	Ad2	SV40	Φ X174	M13mp7	pBR322	pBR328	pUC18
21	9	1	0	0	1	1	0

Activity in PCR buffer Relative activity in PCR mix (Taq DNA Polymerase buffer) is **10%**. The PCR mix contained λ target DNA, primers, 10 mM Tris-HCl (pH 8.3, 20°C), 50 mM KCl, 1.5 mM MgCl₂, 200 μM dNTPs, 2.5 U Taq DNA polymerase. The mix was subjected to 25 amplification cycles.

Ligation and recutting assay

EcoR V fragments obtained by complete digestion of 1 μg λDNA are ligated with 1 U T4-DNA ligase (Cat. No. 10 481 220 001) in a volume of 10 μl in incubation for 16 h at 25°C in 66 mM Tris-HCl, 5 mM MgCl₂, 5 mM Dithiothreitol, 1 mM ATP, pH 7.5 (at 20°C) resulting in >85% recovery of 1 μg λDNA × *EcoR V* fragments. Subsequent re-cutting with *Xho I* yields >95% of the typical pattern of λDNA × *EcoR V* 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., Hg²⁺, Mn²⁺) 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.

Star activity

EcoR V exhibits star activity under non-optimal conditions.

Quality control

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

Absence of unspecific endonuclease activities

1 μg λDNA is incubated for 16 h in 50 μl SuRE/Cut buffer B with excess of *EcoR V*. 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 *EcoR V* 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

- Schildkraut, J. et al. (1984) *Gene* **27**, 327.
- Kessler, C. & Manta, V. (1990) *Gene* **92**, 1-248.
- Rebase The Restriction Enzyme Database: <http://rebase.neb.com>
- Bougueleret, L. et al. (1985) "Overproduction of the *EcoRV* endonuclease and methylase" *Nucl. Acids Res.* **13**, 3823-3839.
- Cholmin, G.V. et al. (1980) *Biochimica* **253**, 495-497.
- Cosstick, R. et al. (1990) "Molecular recognition in the minor groove of the DNA helix. Studies on the synthesis of oligonucleotides and polynucleotides containing 3-deaza-2'-deoxyadenosine. Interaction of the oligonucleotides with the restriction endonuclease *EcoRV*" *Nucl. Acids Res.* **18**, 4771-4778.
- Gubler, U. and Hoffman, B.J. (1983) "A simple and very efficient method for generating cDNA libraries" *Gene* **25**, 263-269.
- Kholmina, G.V. et al. (1980) "Isolation and characterization of a new site-specific endonuclease *EcoRV*" *Dokl. Akad. Nauk SSSR* **253**, 495-497.
- Nwosu, V.U. et al. (1988) "The cloning, purification and characterization of the *EcoRV* modification methylase" *Nucl. Acids Res.* **16**, 3705-3720.
- Taylor, J.D. and Halford, S. E. (1989) "Discrimination between DNA sequences by the *EcoRV* restriction endonuclease" *Biochem.* **28**, 6198-6207
- Winkler, F.K. et al. (1993) "The crystal structure of *EcoRV* endonuclease and of its complexes with cognate and non-cognate DNA fragments" *EMBO J.* **12**, 1781-1795.

Ordering Information

Product	Application	Packsizes	Cat. No.
Restriction Enzymes	DNA restriction digestion	Please refer to website	
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(lacU)169 (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