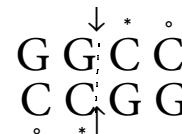


Restriction Endonuclease Hae III

From *Haemophilus aegyptius*

Cat. No. 10 693 944 001 5000 units (10 U/μl)



Version 20

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 *Hae* III recognizes the sequence GG/CC and generates fragments with blunt ends (1).

Compatible ends *Hae* III generates compatible ends to any blunt end.

Isoschizomers The enzyme is an isoschizomer to BsuR I and Pal I.

Methylation sensitivity *Hae* III is inhibited by 5-methyl (*) and 5-hydroxymethylcytosine at the internal C residue. At the external C residue the enzyme is only inhibited by the presence of hydroxymethylcytosine, but not by 5-methylcytosine (°).

Storage buffer 20 mM Tris-HCl, 400 mM NaCl, 0.1 mM EDTA, 10 mM 2-mercaptoethanol, 50% glycerol (v/v), 0.01 polydocanol, pH approx. 7.7 (at 4°C).

Suppl. Incubation buffer, (10x) 100 mM Tris-HCl, 500 mM NaCl, 100 mM MgCl₂, 10 mM Dithioerythritol, pH 7.5 (at 37°C), (= SuRE/Cut Buffer **M**)

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

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

Incubation temperatur **37°C**

Unit definition One unit is the enzyme activity that completely cleaves 1 μg λDNA in 1 h at **37°C** in SuRE/Cut buffer **M** in a total volume of 25 μl. 1 μg pBR322 DNA is digested completely by ca. 2 units of *Hae* III on account of the larger number of cleavage sites per μg pBR322 DNA as compared to λDNA.

Typical experiment

Component	Final concentration
DNA	1 μg
10 × SuRE/Cut Buffer M	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 The enzyme can not 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
149	216	18	11	15	22	30	11

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 MgCl₂, 200 μM dNTPs, 2.5 U Taq DNA polymerase. The mix was subjected to 25 amplification cycles.

Ligation and recutting assay

Hae III fragments obtained by complete digestion of 1 μg λDNA 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 MgCl₂, 5 mM dithiothreitol, 1 mM ATP, pH 7.5 (at 20°C) resulting in >50% recovery of 1 μg λDNA × *Hae* III fragments. Subsequent re-cutting with *Hae* III yields >95% of the typical pattern of λDNA × *Hae* III 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.

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 M with excess of *Hae* 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 *Hae* III 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 Bron, S. & Murray, K. (1975) *Mol. Gen. Genet.* **143**, 25.
- 2 Kessler, C. & Manta, V. (1990) *Gene* **92**, 1-248.
- 3 Rebase The Restriction Enzyme Database: <http://rebase.neb.com>

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) 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

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(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, Δ(hsdRMS) 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