

Restriction Endonuclease *Bcl* I

From *Bacillus caldolyticus*



Cat. No. 10 693 952 001

500 units (10 U/μl)

Cat. No. 11 097 059 001

2500 units, high concentration (40 U/μl)

Version 20

Content version: March 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.
Note: Product is shipped on dry ice.

Sequence specificity *Bcl* I recognizes the sequence T/GATCA and generates fragments with 5'-cohesive termini (1).

Compatible ends *Bcl* I generates compatible ends to *Bam* HI, *Bgl* II, *Nde* II, *Sau* 3A and *Xho* II.

Isoschizomers *Bcl* I is an isoschizomer to *Bsi* QI.

Methylation sensitivity *Bcl* I digestion of DNA is inhibited by the *dam* gene product of *E. coli*, which methylates the 6N position of adenine within the sequence GATC (2). Therefore *Bcl* I digestion of DNA isolated from *E. coli dam* cells is inhibited in methylated TG*ATCA sequences (overlapping *Bcl* I and *dam* recognition sequence) as indicated. *Bcl* I is further inhibited by presence of 5-hydroxymethylcytosine. *Bcl* I is not inhibited by 5-methylcytosine.

Storage buffer 20 mM Tris-HCl, 200 mM NaCl, 0.1 mM EDTA, 10 mM 2-mercaptoethanol, 0.2% Triton X-100, 50% glycerol (v/v), pH approx. 8.0 (at 4 °C).

Incubation buffer (10×, included) 100 mM Tris-HCl, 500 mM NaCl, 100 mM MgCl₂, 10 mM dithioerythritol, pH 7.5 (at 37 °C), (Δ SuRE/Cut Buffer **M**).

Activity in SuRE/Cut Buffer System

A	B	L	M	H
100%	100%	25–50%	100%	100%

Incubation temperature 50 °C

Unit definition One unit is the enzyme activity that completely cleaves 1 μg λ *dam*⁻, *dcm*⁻ DNA in 1 h at 50 °C in SuRE/Cut buffer **M** in a total volume of 25 μl.

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 50 °C for 1 h.

Heat Inactivation The enzyme cannot be heat-inactivated by heating to 65 °C for 15 min.

Number of cleavage sites on different DNA's (2):

λ	Ad2	SV40	Φ X174	M13mp7	pBR322	pBR328	pUC18
8	5	1	0	0	0	1	0

Ligation and recutting assay

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

Troubleshooting

A critical component is the DNA substrate. Many compounds used in the isolation of DNA e.g., phenol, chloroform, EtOH, SDS, high levels of NaCl, metals (e.g., Hg²⁺, Mn²⁺), 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.lifescience.roche.com/certificates

Absence of unspecific endonuclease activities

1 μg λ *dam*⁻, *dcm*⁻ or pBR322 DNA is incubated for 16 h in 50 μl SuRE/Cut buffer **M** with excess of *Bcl* 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 [³H] labeled calf thymus DNA are incubated with 3 μl *Bcl* I 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

- Bingham, A. H. A. et al. (1978) *Nucleic Acids Res.* **5**, 3457.
- McClelland, M. (1981) *Nucleic Acids Res.* **9**, 5859.
- Kessler, C. & Manta, V. (1990) *Gene* **92**, 1–250.
- Rebase The Restriction Enzyme Database: <http://rebase.neb.com>

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

- Update of quality control
- 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. et al (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. et al., (1985) <i>Gene</i> 33 , 103.)
JM109	<i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB) F[traD36proAB⁺, lacI^q lacZΔM15]</i> ; (Yanisch-Perron, C. et al., (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⁺, lacI^q lacZΔM15]</i> ; (Yanisch-Perron, C. et al., (1985) <i>Gene</i> 33 , 103.)
K802	<i>supE hsdR gal metB</i> ; (Raleigh, E. et al., (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⁺ lacI^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⁺, lacI^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⁺, lacI^q lacZΔM15 Tn10 (tet^r)</i> ; (Bullock et al., (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