

# Restriction Endonuclease *Sau3A* I

From *Staphylococcus aureus* 3A

Cat. No. 10 709 751 001

500 units (3–5 U/μl)



**Version 20**

Content version: November 2012

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** *Sau3A* I recognizes the sequence /G°AT°C generating fragments with 5'-cohesive termini (3) which contain the same tetranucleotide sequence GATC as the cohesive termini of the *Bam* HI-, *Bcl* I-, *Bgl* II-, or *Xho* II-fragments (1, 2).

**Compatible ends** The enzyme generates compatible ends to *Bam* HI, *Bcl* I, *Bgl* II, and *Xho* II.

Enzyme with compatible ends	Recognition sequence	New sequence if <i>Sau3A</i> I is ligated to enzyme with compatible ends		Enzyme that can cut this new sequence
		<i>Sau3A</i> I - Enzyme	Enzyme - <i>Sau3A</i> I	
<i>Bam</i> HI	G/GATCC	/GATCC	G/G°AT°C	<i>Dpn</i> I, <i>Nde</i> II, <i>Sau3A</i> I
<i>Bcl</i> I	T/GATCA	/GATCA	T/G°AT°C	<i>Dpn</i> I, <i>Nde</i> II, <i>Sau3A</i> I
<i>Bgl</i> II	A/GATCT	/GATCT	A/G°AT°C	<i>Dpn</i> I, <i>Nde</i> II, <i>Sau3A</i> I, <i>Xho</i> II
<i>Nde</i> II	/G°AT°C	/G°AT°C	/G°AT°C	<i>Dpn</i> I, <i>Nde</i> II, <i>Sau3A</i> I
<b><i>Sau3A</i> I</b>	<b>/G°AT°C</b>	<b>/G°AT°C</b>	<b>/G°AT°C</b>	<b><i>Sau3A</i> I+</b> <b>Isoschizomers</b>
<i>Xho</i> II	Pu/GATCPy	/GATCPy	Pu/G°AT°C	<i>Dpn</i> I, <i>Nde</i> II, <i>Sau3A</i> I

**Isoschizomers** *Sau3A* I is an isoschizomer to *Bsp* 143I, *Dpn* I, *Dpn* II, *Mbo* I, and *Nde* II (1,2).

**Methylation sensitivity** In contrast to the *Mbo* I and *Nde* II-isoschizomer, *Sau3A* I digestion of DNA is not inhibited by the dam gene product of *E. coli*, which methylates the 6N-position of adenine(ε) within the sequence GATC (1, 2). 5-methylcytosine, 4-methylcytosine, and 5-hydroxymethylcytosine at the C-position are inhibiting(\*)

**Storage buffer** 20 mM Tris-HCl, 250 mM NaCl, 0.1 mM EDTA, 5 mM 2-Mercaptoethanol, 0.01% Polydocanol (v/v), 50% Glycerol (v/v), pH approx. 7.5 (at 4°C).

**Suppl. Incubation buffer (10x)** 330 mM Tris-acetate, 100 mM Mg-acetate, 660 mM K-acetate, 5 mM Dithiothreitol, pH 7.9 (at 37°C), (Δ SuRE/Cut Buffer **A**).

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

<b>A</b>	B	L	M	H
<b>100%</b>	25–50%	25–50%	75–100%	0–10%

**Incubation temp.** **37°C**

**Unit definition** One unit is the enzyme activity that completely cleaves 1 μg λDNA in 1 h at **37°C** in the SuRE/Cut Buffer **A** in a total volume of 25 μl. 1 μg pBR322 DNA is digested completely by approx. 5 units of *Sau3A* I because 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 <b>A</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** There is no information about *Sau3A* I and heat inactivation available.

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

λ	Ad2	SV40	Φ X174	M13mp7	pBR322	pBR328	pUC18
116	87	8	0	8	22	27	15

## 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 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

*Sau3A* I fragments obtained by complete digestion of 1 μg λDNA fragments 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<sub>2</sub>, 5 mM Dithiothreitol, 1 mM ATP, pH 7.5 (at 20°C) resulting in >95 % recovery of 1 μg λDNA × *Sau3A* I fragments.

Subsequent re-cutting with *Sau3A* I yields > 95% of the typical pattern of λDNA × *Sau3A* I fragments.

## Troubleshooting

A critical component is the DNA substrate. Many compounds used in the isolation of DNA, for example, phenol, chloroform, EtOH, SDS, high levels of NaCl, metals (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 EtOH precipitation followed by drying, before the DNA is added to the restriction digest reaction. Appropriate mixing of the enzyme is recommended.

☎ Check out the Restrictions Enzymes Frequently Asked Questions at <http://www.roche-applied-science.com/support>.

## 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 A with excess of *Sau3A* 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 *Sau3A* 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

- 1 Roberts, R. J. (1983) *Nucleic Acids Res.* **11**, r135.
- 2 Kessler, C. & Manta, V. (1990) *Gene* **92**, 1–248.
- 3 Sussenbach, J. S. et al. (1976) *Nucleic Acids Res.* **3**, 3193.
- 4 Rexer, B. et al., Roche Diagnostics GmbH, unpublished.
- 5 Rebase The Restriction Enzyme Database: <http://rebase.neb.com>
- 6 Benchmate: <http://www.roche-applied-science.com/benchmate>

## Ordering Information

Roche Applied Science offers a large selection of reagents and systems for life science research. For a complete overview of related products and manuals, please visit and bookmark our homepage, [www.roche-applied-science.com](http://www.roche-applied-science.com), and our Special Interest Sites, including "Mapping & Cloning": <http://www.restriction-enzymes.com>.

The convenient RE Finder Program located on our Bench Mate website, <http://www.roche-applied-science.com/benchmate> helps you identify the enzymes that will cut your DNA sequence, and displays the names and recognition sequences of enzymes, and isoschizomers as well as links to detailed information (e.g., Instructions for Use) of the selected restriction enzyme.

Product	Application	Pack Size	Cat. No.
Restriction Enzymes	DNA restriction digestion.	Please refer to website or catalog	
T4 DNA Ligase	Ligation of sticky- and blunt-ended DNA fragments.	100 U 500 units (1 U/μl) 500 units (5 U/μl)	10 481 220 001 10 716 359 001 10 799 009 001
Rapid DNA Dephos & Ligation Kit	Upgrade from the Rapid DNA Ligation Kit for fast and efficient DNA dephosphorylation and ligation of sticky- or blunt-ended DNA fragments.	40 reactions 160 reactions	04 898 117 001 04 898 125 001
rAPid Alkaline Phosphatase	Dephosphorylation of 5'-phosphate residues from nucleic acids.	1,000 U 5,000 U	04 898 133 001 04 898 141 001
Rapid DNA Ligation Kit	Ligation of nucleic acids.	Kit (40 DNA ligations)	11 635 379 001
Alkaline Phosphatase (AP), special quality for molecular biology	Dephosphorylation of 5'-phosphate residues from nucleic acids.	1,000 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	Specialty 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

**Printed Materials** You can view the following manuals on our website:

Lab FAQs "Find a Quick Solution"
Restriction Enzyme Ordering Guide
Molecular Weight Markers for Nucleic Acids

## Changes to previous version

Star activity information removed.

## 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.

## 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. et al (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/lacΔM15) 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. et al., (1985) <i>Gene</i> <b>33</b> , 103.)
JM109	<i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB) F[traD36proAB<sup>+</sup>, lacI<sup>q</sup> lacZΔM15]</i> ; (Yanisch-Perron, C. et al., (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>, lacI<sup>q</sup> lacZΔM15]</i> ; (Yanisch-Perron, C. et al., (1985) <i>Gene</i> <b>33</b> , 103.)
K802	<i>supE hsdR gal metB</i> ; (Raleigh, E. et al., (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<sup>-</sup>, Δ(hsdRMS) endA1 gyrA96 thi relA1 supE44 F[proAB<sup>+</sup> lacI<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>, lacI<sup>q</sup> lacZΔM15]</i> ; (Gibson, T.J. (1984) PhD Theses. Cambridge University, U.K.)
XL1-Blue <sup>f</sup>	<i>supE44 hsdR17 recA1 endA1 gyrA46 thi relA1 lac F[proAB<sup>+</sup>, lacI<sup>q</sup> lacZΔM15 Tn10 (tet<sup>r</sup>)]</i> ; (Bullock et al., (1987) <i>BioTechniques</i> , <b>5</b> , 376.)

## Contact and Support

To ask questions, solve problems, suggest enhancements or report new applications, please visit our **Online Technical Support Site** at:

**[www.roche-applied-science.com/support](http://www.roche-applied-science.com/support)**

To call, write, fax, or email us, visit the Roche Applied Science home page, [www.roche-applied-science.com](http://www.roche-applied-science.com), and select your home country. Country-specific contact information will be displayed. Use the Product Search function to find Pack Inserts and Material Safety Data Sheets.



Roche Diagnostics GmbH  
Roche Applied Science  
68298 Mannheim  
Germany