

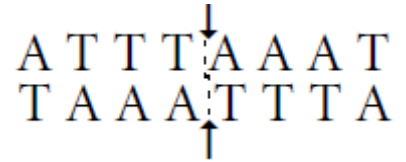
For life science research only.
Not for use in diagnostic procedures.



Restriction Endonuclease

Swa I

from *Staphylococcus warneri*



 **Version: 11**

Content Version: March 2020

Cat. No. 11 371 517 001 200 U
 10 U/μl

Store product at –15 to –25°C.

1.	General Information	3
1.1.	Contents	3
1.2.	Storage and Stability	3
	Storage Conditions (Product)	3
1.3.	Application	3
2.	How to Use this Product	4
2.1.	Protocols	4
2.2.	Parameters	4
	Activity in PCR Buffer	4
	Buffers	4
	Activity in SuRE/Cut Buffer System	4
	Cleavage Sites	4
	Number of cleavage sites on different DNAs	4
	Compatible Ends	5
	Inactivation	5
	Isoschizomers	5
	Methylation Sensitivity	5
	Recognition Sites	5
	Temperature Optimum	5
	Unit Definition	5
3.	Troubleshooting	6
4.	Additional Information on this Product	7
4.1.	Test Principle	7
	Genomic digest	7
	Commonly used bacterial strains	7
4.2.	References	8
4.3.	Quality Control	8
5.	Supplementary Information	9
5.1.	Conventions	9
5.2.	Changes to previous version	9
5.3.	Ordering Information	9
5.4.	Trademarks	10
5.5.	License Disclaimer	10
5.6.	Regulatory Disclaimer	10
5.7.	Safety Data Sheet	10
5.8.	Contact and Support	10

1. General Information

1.1. Contents

Vial / Bottle	Cap	Label	Function / Description	Content
Swa I	red	Swa I	Contains 20 mM Tris-HCl, 500 mM NaCl, 0.1 mM EDTA, 10 mM 2-mercaptoethanol, 50% glycerol (v/v), 0.01% polydocanol, pH approximately 8.0 (+4°C).	1 vial, 200 U (10 U/μl)
H	red	SuRE/Cut Buffer H for Restriction Enzymes, 10x conc.	Contains 500 mM Tris-HCl, 1 M NaCl, 100 mM MgCl ₂ , 10 mM dithioerythritol, pH 7.5 (+37°C).	1 vial, 1 ml

1.2. Storage and Stability

Storage Conditions (Product)

The product is shipped on dry ice.

When stored at –15 to –25°C, the product is stable through the expiration date printed on the label.

Vial / Bottle	Cap	Label	Storage
Swa I	red	Swa I	Store at –15 to –25°C. ⚠ Do not store below –25°C.
H	red	SuRE/Cut Buffer H, 10x conc.	Store at –15 to –25°C.

1.3. Application

Swa I recognizes the sequence ATTT/AAAT and generates fragments with blunt ends.

2. How to Use this Product

2.1. Protocols

The following steps describe a typical experiment.

- 1 Prepare the restriction digest according to the following table.

Reagent	Final conc.
DNA	1 µg
10x SuRE/Cut Buffer H	2.5 µl
Water, PCR Grade*	Up to total volume of 25 µl
Swa I	1 U

- 2 Incubate at +25°C for 1 hour.

2.2. Parameters

Activity in PCR Buffer

Not tested.

Buffers

Activity in SuRE/Cut Buffer System

A	H ⁽¹⁾	M
0 to 10%	100% ⁽²⁾	0 to 10%

⁽¹⁾ Supplied Buffer

⁽²⁾ Indicates recommended buffer for optimal activity.

Cleavage Sites

Number of cleavage sites on different DNAs

λ	Ad2	SV40	ΦX174	M13mp7	pBR322	pBR328	pUC18
0	1	1	0	1	0	0	-

Compatible Ends

Swa I generates compatible ends to Pme I.

Inactivation

Swa I cannot be heat inactivated by incubation at +65°C for 15 minutes.

Isoschizomers

The enzyme is not known to have isoschizomers.

Methylation Sensitivity

There is no information available.

Recognition Sites

ATTAAAT

Temperature Optimum

+25°C

 *Swa I has a special incubation temperature.*

Unit Definition

One unit is the enzyme activity that completely cleaves 1 µg Expand vector III DNA (15.2 kb) in 1 hour at +25°C in a total volume of 25 µl SuRE/Cut Buffer H. 1 µg M13 mp9 DNA is digested completely by 3 U of Swa I.

3. Troubleshooting

Observation	Possible cause	Recommendation
Inhibition or alteration of recognition specificity of restriction enzyme.	Compounds were used in the isolation of the DNA substrate, such as phenol, chloroform, ethanol, SDS, high levels of NaCl, and metal ions, such as Hg ²⁺ and Mn ²⁺ .	Remove compounds by ethanol precipitation followed by drying, before adding DNA to the restriction digest reaction. <hr/> Mix vial of restriction enzyme gently but completely prior to use.

4. Additional Information on this Product

4.1. Test Principle

Genomic digest

The enzyme is very useful for GC-rich bacterial genomes. 1 µg of *E. coli* genomic DNA embedded in agarose for PFGE analysis is completely digested with 20 U Swa I in 4 hours at +25°C in SuRE/Cut buffer H containing 100 µg/ml bovine serum albumin. The largest resulting fragment has approximately 230 kb. An excess of 100 U Swa I does not change the banding pattern during an incubation time of 16 hours. On *Staphylococcus aureus* DNA with a GC content of 34%, Swa I cleaves more frequently; the size range of the fragments obtained is in a range <50 kb. On GC-rich genomic DNA (*Rhodobacter sphaeroides* with a GC content of 66%), Swa I cleaves much less frequently: one fragment has a size of 160 kb, the residual cleavage products are in the range of 800 kb. When human DNA in agarose plugs was digested with Swa I, the fragments were found mainly in a size range of 30 to 100 kb.

Commonly used bacterial strains

Strain	Genotype
BL21	<i>E. coli</i> B F ⁻ dcm ompT hsdS(<i>r_B</i> - <i>m_B</i> -) gal (Studier FW, et al, 1986).
C600 ^o	<i>supE44 hsdR2 thi-1 thr-1 leuB6 lacY1 tonA21</i> (Hanahan D, 1983).
DH5α	<i>supE44 Δ(lacU169 (Φ80d)lacZΔM15) hsdR17 recA1 endA1 gyrA96 thi-1 relA1</i> (Hanahan D, 1983).
HB101	<i>supE44 hsdS20 recA13 ara-14 proA2 lacY1 galK2 rpsL20 xyl-5 mtl-1</i> (Hanahan D, 1983).
JM108	<i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB)</i> (Yanisch-Perron C, et al, 1985).
JM109	<i>recA1 supE44 endA1 hsdR17 gyrA96 relA1 thi Δ(lac-proAB) F'[traD36proAB⁺, lacI^q lacZΔM15]</i> (Yanisch-Perron C, et al, 1985).
JM110	<i>rpsL (Str^r) thr leu thi-I lacY galK galT ara tonA tsx dam dcm supE44 Δ(lac-proAB) F'[traD36proAB⁺, lacI^q lacZΔM15]</i> (Yanisch-Perron C, et al, 1985).
K802	<i>supE hsdR gal metB</i> (Raleigh E, et al, 1986; Wood WB, 1966).
SURE ^r	<i>recB recJ sbc C201 uvrC umuC::Tn5(kan^r) lac, Δ(hsdRMS) endA1 gyrA96 thi relA1 supE44 F'[proAB⁺ lacI^q lacZΔM15 Tn10 (tet^r)</i> (Greener A, 1990).
TG1	<i>supE hsd Δ5 thi Δ(lac-proAB) F'[traD36proAB⁺, lacI^q lacZΔM15]</i> (Gibson TJ, 1984).
XL1-Blue ^r	<i>supE44 hsdR17 recA1 endA1 gyrA46 thi relA1 lac F'[proAB⁺, lacI^q lacZΔM15 Tn10 (tet^r)</i> (Bullock WO, et al, 1987).

4. Additional Information on this Product

4.2. References

- Bullock WO, Fernandez JM, Short JM. XL1-Blue– a high-efficiency plasmid transforming recA Escherichia coli strain with β -galactosidase selection. BioTechniques. 1987;5:376-379.
- Gibson, TJ. PhD Theses. Cambridge University, U.K 1984.
- Greener, A. Strategies 1990;3:5.
- Hanahan D. Studies on transformation of Escherichia coli with plasmids. J Mol Biol.1983;166:557-580.
- Raleigh EA, Wilson G. Escherichia coli K-12 restricts DNA containing 5-methylcytosine. Proc Natl Acad Sci USA.1986;83:9070-9074.
- Studier FW, Moffatt BA. Use of bacteriophage T7 RNA polymerase to direct selective high-level expression of cloned genes. J Mol Biol.1986;189:113-130.
- Yanisch-Perron C, Vieira J, Messing J. Improved M13 phage cloning vectors and host strains: nucleotide sequences of the M13mp18 and pUC19 vectors. Gene.1985;33:103-19.
- Wood WB. Host specificity of DNA produced by Escherichia coli: bacterial mutations affecting the restriction and modification of DNA. J Mol Biol.1966;16:118-133.

4.3. Quality Control

For lot-specific certificates of analysis, see section **Contact and Support**.

5. Supplementary Information

5.1. Conventions

To make information consistent and easier to read, the following text conventions and symbols are used in this document to highlight important information:

Text convention and symbols

 *Information Note: Additional information about the current topic or procedure.*

 **Important Note: Information critical to the success of the current procedure or use of the product.**

① ② ③ etc. Stages in a process that usually occur in the order listed.

① ② ③ etc. Steps in a procedure that must be performed in the order listed.

* (Asterisk) The Asterisk denotes a product available from Roche Diagnostics.

5.2. Changes to previous version

Editorial changes.

5.3. Ordering Information

Product	Pack Size	Cat. No.
Reagents, kits		
T4 DNA Ligase	100 U, 1 U/μl	10 481 220 001
	500 U, 1 U/μl	10 716 359 001
	500 U, 5 U/μl	10 799 009 001
Water, PCR Grade	25 ml, 25 x 1 ml	03 315 932 001
	25 ml, 1 x 25 ml	03 315 959 001
	100 ml, 4 x 25 ml	03 315 843 001
SuRE/Cut Buffers	SuRE/Cut Buffer A, 5 x 1 ml	11 417 959 001
	SuRE/Cut Buffer M, 5 x 1 ml	11 417 983 001
	SuRE/Cut Buffer H, 5 x 1 ml	11 417 991 001
1,4-Dithiothreitol	2 g	10 197 777 001
	10 g	10 708 984 001
	25 g	11 583 786 001

5. Supplementary Information

5.4. Trademarks

EXPAND is a trademark of Roche.

All other product names and trademarks are the property of their respective owners.

5.5. License Disclaimer

For patent license limitations for individual products please refer to:

List of biochemical reagent products.

5.6. Regulatory Disclaimer

For life science research only. Not for use in diagnostic procedures.

5.7. Safety Data Sheet

Please follow the instructions in the Safety Data Sheet (SDS).

5.8. Contact and Support

To ask questions, solve problems, suggest enhancements or 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.

