

Restriction Endonuclease *Mun* I (*Mfe* I)

From *Mycoplasma species*



Cat. No. 11 441 337 001

200 units (10 U/ μ l)

Version 17
Content version: July 2011

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: The product is shipped on dry ice.

Sequence specificity *Mun* I recognizes the sequence C/AATTG and generates fragments with 5'-cohesive termini (1).

Compatible ends The enzyme generates compatible ends to *Acs* I and *Eco* RI.

Enzyme with compatible ends	Recognition sequence	New sequence if <i>Mun</i> I is ligated to enzyme with compatible ends		Enzyme that can cut this new sequence
		<i>Mun</i> I-Enzyme	Enzyme- <i>Mun</i> I	
<i>Acs</i> I	(A,G)/AATT(C,T)	C/AATT(C,T)	(A,G)/AATTG	<i>Mun</i> I, <i>Tsp</i> EI
<i>Eco</i> RI	G/AATTC	C/AATTC	G/AATTG	<i>Tsp</i> EI
<i>Mun</i> I	C/AATTG	C/AATG	C/AATTG	<i>Mun</i> I + <i>Mfe</i> I

Isoschizomers *Mun* I is an isoschizomer to *Mfe* I.

Methylation sensitivity *Mun* I is inhibited by the presence of N⁵-methyladenine in the sequence C/A^mATTG.

Storage buffer 10 mM Tris-HCl, 100 mM KCl, 1 mM EDTA, 1 mM Dithioerythritol, 200 μ g/ml bovine serum albumin, 50% glycerol (v/v), pH approx. 8.0 (at 4 °C).

Incubation buffer, (10 \times , 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 Bold face printed buffer indicates the recommended buffer for optimal activity:

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

Incubation temperature **37 °C**

Unit definition One unit is the enzyme activity that completely cleaves 1 μ g λ DNA in 1 h at 37 °C in a total volume of 25 μ l SuRE/Cut buffer **M**.

Typical experiment

Component	Final concentration
DNA	1 μ g
10 \times SuRE/Cut Buffer M	2.5 μ l
Repurified water	Up to a total volume of 25 μ l
Restriction enzyme	1 unit

Incubate at **37 °C** for 1 h.

Heat inactivation *Mun* I can be heat-inactivated by 15 min incubation at 65 °C up to a concentration of 15 U/ μ g DNA, higher concentrations of enzyme can no more be completely inactivated under these conditions.

Number of cleavage sites on different DNAs (2):

λ	Ad2	SV40	Φ X174	M13mp7	pBR322	pBR328	pUC18
8	4	4	1	0	0	0	2

PFGE tested

Mun I has been tested in Pulsed-Field Gel Electrophoresis (test system bacterial chromosomes). For cleavage of genomic DNA (*E. coli* C600) embedded in agarose for PFGE analysis 10 units of enzyme/ μ g DNA and 4 h incubation time are recommended.

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

The recognition specificity of *Mun* I is relaxed in extended incubation with excess enzyme, or at low ionic strength.

Quality control

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

Absence of unspecific endonuclease activities

1 μ g λ DNA is incubated for 16 h in 50 μ l SuRE/Cut buffer **M** with excess of *Mun* 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 μ g [³H] labeled calf thymus DNA are incubated with 3 μ l *Mun* I for 4 h at 37 °C in a total volume of 100 μ l 50 mM Tris-HCl, 10 mM MgCl₂, 1 mM Dithioerythritol, pH approx. 7.5. The release of radioactivity is calculated as a percentage value of liberated to input radioactivity per unit of enzyme (stated in the certificate of analysis).

Ligation and recutting assay

Mun I fragments obtained by complete digestion of 1 μ g λ DNA are ligated with 1 U T4-DNA ligase (Cat. No. 10 481 220 001) in a volume of 10 μ l by incubation for 16 h at 4 °C in 66 mM Tris-HCl, 5 mM MgCl₂, 5 mM dithioerythritol, 1 mM ATP, pH 7.5 (at 20 °C).

The percentage of ligation and subsequent recutting with *Mun* I which yields the typical pattern of $\lambda \times$ *Mun* I fragments is determined and stated in the certificate of analysis.

References

- 1 Janulaitis, A. et al., unpublished observations.
- 2 Kessler, C. & Manta, V. (1990) *Gene* **92**, 1-248.
- 3 Rebase The Restriction Enzyme Database: <http://rebase.neb.com>
- 4 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 home page, 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. package insert) of the selected restriction enzyme.

Product	Application	Packsize	Cat. No.
Restriction Enzymes	DNA restriction digestion	Please refer to website or catalogue	
Rapid DNA Ligation Kit	Ligation of sticky- or blunt-ended DNA fragments in just 5 min at 15 - 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

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

Lot-specific information is no longer printed on the label of the product. Instead, the address for certificates of analysis is provided (www.roche-applied-science.com/certificates).

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 ⁻ <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⁺, lacI^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⁺, lacI^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⁺ 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 <i>et al.</i> , (1987) <i>BioTechniques</i> , 5 , 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

To call, write, fax, or email us, visit the Roche Applied Science home page, 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