

Ordering Information

Product	Application	Packsizes	Cat. No.
T4 DNA Ligase	Ligation of sticky- and blunt- ended DNA fragments.	100 U 500 units (1 U/ μ l)	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 \times conc. solutions)	11 082 035 001
SuRE/Cut Buffer A	Restriction enzyme incubation	5 \times 1 ml (10 \times conc. solution)	11 417 959 001
SuRE/Cut Buffer B	Restriction enzyme incubation	5 \times 1 ml (10 \times conc. solution)	11 417 967 001
SuRE/Cut Buffer H	Restriction enzyme incubation	5 \times 1 ml (10 \times conc. solution)	11 417 991 001
SuRE/Cut Buffer L	Restriction enzyme incubation	5 \times 1 ml (10 \times conc. solution)	11 417 975 001
SuRE/Cut Buffer M	Restriction enzyme incubation	5 \times 1 ml (10 \times conc. solution)	11 417 983 001
Water, PCR Grade	Specially purified, double-distilled, deionized, and autoclaved	100 ml (4 vials of 25 ml)	03 315 843 001
		25 ml (25 vials of 1 ml)	03 315 932 001
		25 ml (1 vial of 25 ml)	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</i> (r _B - m _B -) <i>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.)
DH5 α	<i>supE44</i> Δ (<i>lac</i> U169 (ϕ 80d/ <i>lacZ</i> Δ M15) <i>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</i> Δ (<i>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</i> Δ (<i>lac-proAB</i>) F[<i>traD36proAB</i> ⁺ , <i>lacI</i> ^q <i>lacZ</i> Δ M15]; (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</i> Δ (<i>lac-proAB</i>) F[<i>traD36proAB</i> ⁺ , <i>lacI</i> ^q <i>lacZ</i> Δ M15]; (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</i> F[<i>proAB</i> ⁺ <i>lacI</i> ^q <i>lacZ</i> Δ M15 Tn10 (<i>tet</i> ^r); (Greener, A. (1990) <i>Stratagies</i> , 3 , 5.)
TG1	<i>supE hsd</i> Δ 5 <i>thi</i> Δ (<i>lac-proAB</i>) F[<i>traD36proAB</i> ⁺ , <i>lacI</i> ^q <i>lacZ</i> Δ M15]; (Gibson, T.J. (1984) <i>PhD Theses. Cambridge University, U.K.</i>)
XL1-Blue ^f	<i>supE44 hsdR17 recA1 endA1 gyrA46 thi relA1 lac</i> F[<i>proAB</i> ⁺ , <i>lacI</i> ^q <i>lacZ</i> Δ M15 Tn10 (<i>tet</i> ^r); (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