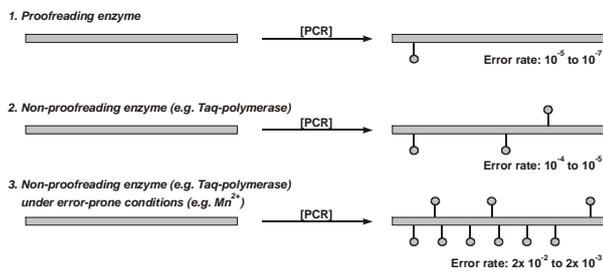




JBS Error-Prone Kit

Random Mutagenesis by Error-Prone PCR

Cat. No.	Amount
PP-102	15 reactions



Enhanced mutational rate by error-prone PCR compared to standard PCR reactions.

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at $-20\text{ }^{\circ}\text{C}$

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Description:

JBS Mutagenesis Series

Within three billion years of evolution, nature has produced a plethora of proteins simply by repeated cycles of random mutagenesis followed by *in vivo* selection for superior function of the encoded proteins. This example of natural evolution has guided researchers within the last two decades to develop strategies for *in vitro* permutation of proteins.

Among the variety of strategies applied, three major powerful techniques have emerged.

Random Mutagenesis by dNTP Analogs

This method is based on the incorporation of mutagenic dNTP analogs, such as 8-oxo-dGTP and dPTP, into an amplified DNA fragment by PCR. The mutagenic dNTPs are eliminated by a second PCR step in the presence of the four natural dNTPs only, resulting in a rate of mutagenesis of up to 20%.

→ **JBS dNTP-Mutagenesis Kit #PP-101**

Random Mutagenesis by Error-Prone PCR

Developed by Caldwell & Joyce (1992) this method introduces mutations in the gene of interest using a PCR reaction under conditions that induce an increased error-rate of the DNA-polymerase. The rate of mutagenesis achieved by error-prone PCR is in the range of 0.6-2.0%.

→ **JBS Error-prone Kit #PP-102**

Random Mutagenesis by DNA Shuffling

Developed by Stemmer (1994) DNA shuffling generates libraries by random fragmentation of one gene or a pool of related genes, followed by the reassembly of the fragments in a self-priming PCR reaction. This method allows the recombination of sequences from different, related genes. The overall rate of mutagenesis is approx. 0.7%.

→ **JBS DNA-Shuffling Kit #PP-103**

Jena Bioscience now offers all components necessary for each of these techniques 'ready-to-go' in a separate kit, accompanied by a streamlined documentation that maximizes success.

Content:

Taq Polymerase (red cap)

5 units/ μl , 20 μl

10x Reaction Buffer (blue cap)

10x concentration, 100 μl

10x Error-prone Solution (yellow cap)

10x concentration, 100 μl

dNTP Error-prone Mix (white cap)

unbalanced dNTP ratio (dATP, dCTP, dGTP, dTTP), 40 μl



JBS Error-Prone Kit

Random Mutagenesis by Error-Prone PCR

PCR-grade Water (white cap)
1 ml

Random Mutagenesis by Error-Prone PCR

The standard DNA polymerases used in conventional PCR reactions display error rates that are usually not suitable for directed mutagenesis experiments. For example, proofreading enzymes such as *Pfu* exhibit error rates in the range from 10^{-6} to 10^{-7} whereas non-proofreading enzymes like *Taq* Polymerase show error rates in the range from 10^{-4} to 10^{-5} . This rate however, can be significantly enhanced by modifying the following parameters of a PCR-reaction:

- Higher Mg²⁺-concentration of up to 7 mM
- Partial substitution of Mg²⁺ by Mn²⁺
- Optimized dNTP concentrations at unbalanced rates

Recommended assay preparation

- For a 50 µl reaction, take 5 µl of 10x Reaction Buffer in a sterile vial and refer to Tab. 1.
- Add 2 µl dNTP Error-prone Mix.
- Add template and appropriate primers. Note that depending on the template the required concentration can be up to 10 times higher compared to standard PCR.
- Add 0.4-1 µl *Taq* Polymerase as recommended.
- Add PCR-grade Water to a final volume of 45 µl.
- Add 5 µl of 10x Error-prone Solution. Note that the solution should be added last to the reaction mixture to prevent precipitation and must be protected from oxidation (oxidation of Mn²⁺ to Mn³⁺ may destroy the polymerase).

Tab. 1: Amounts of components for error-prone PCR conditions (50 µl PCR assay)

Component	Amount	Final conc.	Cap
10x Reaction Buffer	5 µl	1x	blue
dNTP Error-prone Mix	2 µl	unbalanced ratio	white
Primers		20-100 pmol	
Template		3-100 fmol / 2 - 50 ng	
<i>Taq</i> Polymerase	0.4-1 µl	2-5 units	red
PCR-grade Water	Fill up to 45 µl		white
10x Error-prone Solution	5 µl	1x	yellow

Recommended thermocycling conditions

Denaturation	94° C	30 sec
Annealing ¹⁾	approx. 45-68° C	30 sec
Extension ²⁾	72° C	1 min

Number of cycles: 30

¹⁾ The annealing temperature depends on the melting temperature of the primers.

²⁾ The elongation time depends on the length of the fragments to be amplified. A time of 1 min per kbp is recommended.

For optimal specificity and amplification an individual optimization of the recommended parameters may be necessary for each new template DNA and/or primer pair.

Selected References:

Kim *et al.* (2001) Improvement of tagatose conversion rate by genetic evolution of thermostable galactose isomerase. *Biotechnol. Appl. Biochem.* **34**:99.

Daugherty *et al.* (2000) Quantitative analysis of the effect of the mutation frequency on the affinity maturation of single chain Fv antibodies. *PNAS* **97**:2029.

Wan *et al.* (1998) In vitro evolution of horse heart myoglobin to increase peroxidase activity. *Proc. Natl. Acad. Sci. USA* **95**:12825.

Cline *et al.* (1996) PCR fidelity of *Pfu* DNA polymerase and other thermostable DNA polymerases. *Nucleic Acids Res.* **24**:3546.

Vartanian *et al.* (1996) Hypermutagenic PCR involving all four transitions and a sizeable proportion of transversions. *Nucleic Acids Res.* **24**:2627.

Cadwell *et al.* (1992) Randomization of genes by PCR mutagenesis. *PCR Meth. Appl.* **2**:28.

Kunkel (1992) DNA replication fidelity. *J. Biol. Chem.* **267**:18251.