



Technical notes provide unique applications, innovative methods, and clear protocols designed specifically for Agilent reagents and instruments.

# High-Fidelity PCR Enzymes: Properties and Error Rate Determinations

The polymerase chain reaction (PCR) has revolutionized biological sciences, in particular genetics and proteomics. Since the introduction of Taq DNA polymerase in the late 1980s, significant progress has been made in developing PCR enzyme formulations with improved fidelity, PCR performance, and speed. This Technical Note surveys commercial PCR enzymes developed for high-fidelity PCR applications, such as cloning, mutation detection, and site-directed mutagenesis. We provide detailed information regarding the composition, PCR characteristics, and applications of proofreading DNA polymerases and DNA polymerase blends. We discuss methods for determining DNA polymerase error rates, and provide an in-depth description of the procedure and results obtained using the *lacl*-based phenotypic mutation assay.

### Introduction

High-fidelity PCR enzymes are valuable for minimizing the introduction of amplification errors in products that will be cloned, sequenced, and expressed. Significant time and effort can be saved by employing high-fidelity amplification procedures that eliminate the need for downstream error-correction steps and minimize the number of clones that must be sequenced in order to obtain error free constructs or accurate consensus sequences. Moreover, the use of high-fidelity amplification conditions is essential when analyzing very small amounts of template DNA or rare molecules in heterogeneous populations<sup>1</sup>. Amplifications employing small amounts of template DNA are especially prone to high mutant frequencies due to PCR-generated errors in early cycles ("jackpot" artifacts) and high target doublings<sup>1</sup>. When analyzing rare sequences, such as allelic polymorphisms in individual mRNA transcripts<sup>2</sup>, allelic stages of single cells<sup>3</sup>, or rare mutations in human cells<sup>4</sup>, it is essential that polymerase-generated errors ("PCR-induced noise") are minimized to prevent masking of rare DNA sequences.

PCR fidelity is largely determined by the intrinsic error rate of a DNA polymerase under the reaction conditions employed. Parameters contributing to DNA polymerase fidelity have been reviewed<sup>5-8</sup> and include the tendency of a polymerase to incorporate incorrect nucleotides, the rate at which the enzyme can extend from mispaired 3' primer termini, and the presence of an integral 3'-5' exonuclease (proofreading) activity, which can remove mispaired bases. The importance of proofreading is evident in comparisons of base substitution error rates between non-proofreading ( $10^{-6}$  to  $10^{-7}$ ) DNA polymerase<sup>5.9</sup>. DNA polymerase error rates are influenced by PCR reaction conditions, and can be minimized by optimizing pH, Mg<sup>2+</sup> concentration, and nucleotide concentrations <sup>9-12</sup>.

*Taq* DNA polymerase is suitable for a number of PCR applications, and is still considered by many to be the industry standard. However, the performance of *Taq* is limited in more challenging applications, such as those requiring high fidelity, synthesis of long (> 2 kb) amplicons, and amplification of GC-rich sequences. *Taq* DNA polymerase lacks proofreading activity, and as a result, exhibits relatively poor fidelity.

## Background

#### **Proofreading Archaeal DNA Polymerases**

High-fidelity PCR enzymes include proofreading archaeal DNA polymerases (Table 1) and DNA polymerase blends (Table 2). Commercial proofreading DNA polymerases have been obtained from Thermococcus and Pyrococcus species of hyperthermophilic archaea and are classified as Family B-type DNA polymerases<sup>13</sup>. Unlike thermophilic eubacterial DNA polymerases (e.g., *Taq*), which may or may not possess 3'-5' exonuclease activity, all archaeal B-type DNA polymerases possess proofreading activity and lack an associated 5'-3' exonuclease activity.

The kinetic properties of several thermostable DNA polymerases have been reported<sup>13-15</sup>. Comparisons of steady-state kinetic parameters indicate that archaeal proofreading DNA polymerases exhibit lower K<sub>m</sub> [DNA] values (0.01–0.7 nM) and similar K<sub>m</sub> [dNTPs] values (16–57  $\mu$ M) compared to those reported for *Taq* (1–4 nM, K<sub>m</sub> [DNA]; 16–24  $\mu$ M, K<sub>m</sub> [dNTPs]). Most archaeal proofreading DNA polymerases (*Pfu*, Deep Vent) exhibit limited processivity (< 20 bases) in vitro (Table 1). The only known exceptions are KOD DNA polymerase, which is reported to be 10- to 15-fold more processive than *Pfu* and Deep Vent DNA polymerases<sup>14</sup>, and archaeal DNA polymerases that have been engineered for

increased processivity by fusion to DNA-binding proteins (see Archaeal DNA Polymerase Fusions section). Polymerization rates determined for thermostable DNA polymerases range from 9–25 nucleotides/second (*Pfu*) up to 47–61 nucleotides/second (*Taq*) and 106–138 nucleotides/ second (KOD)<sup>14, 15</sup>.

Unlike *Taq*, which possesses a structure-specific 5'-3' endonuclease activity that cleaves 5' flap structures<sup>16</sup>, archaeal DNA polymerases exhibit temperature-dependent strand displacement activity (e.g., detectable at  $\geq$  70°C for *Pfu*<sup>15,17</sup>). *Taq* DNA polymerase also adds extra non-template directed nucleotide(s) to the 3' ends of PCR fragments, and as a result, *Taq*-generated PCR products can be directly cloned into vectors containing 3'-T overhangs<sup>18,19</sup>. In contrast, archaeal DNA polymerases lack terminal extendase activity, and hence, produce blunt fragments that can be cloned directly into blunt-ended vectors<sup>18,20</sup>.

#### **Uracil Poisoning of Archaeal DNA Polymerases**

Unlike *Taq*, archaeal DNA polymerases possess a "read-ahead" function that detects uracil (dU) residues in the template strand and stalls synthesis<sup>21</sup>. Uracil detection is unique to archaeal DNA polymerases (e.g., *Pfu*), and is thought to represent the first step in a pathway to repair DNA cytosine deamination (dCMP  $\rightarrow$  dUMP) in archaea<sup>21</sup>. Stalling of DNA synthesis opposite uracil has significant implications for high-fidelity amplification with archaeal DNA polymerases. Techniques requiring dUTP (e.g., dUTP/UDG decontamination methods<sup>22</sup>) or uracil-containing oligonucleotides cannot be performed with proofreading DNA polymerases<sup>23,24</sup>. Even more importantly, uracil stalling has been shown to compromise the performance of archaeal DNA polymerases under standard PCR conditions<sup>25</sup>.

We found that during PCR amplification, a small amount of dCTP undergoes deamination to dUTP (%dUTP varies with cycling time), and is subsequently incorporated by archaeal DNA polymerases. Once incorporated, uracil-containing DNA inhibits archaeal DNA polymerases, limiting their efficiency. We found that adding a thermostable dUTPase (dUTP  $\rightarrow$  dUMP + PPi) to amplification reactions carried out with *Pfu* and Deep Vent DNA polymerases significantly increases PCR product yields by preventing dUTP incorporation<sup>25</sup>. Moreover, the target-length capability of *Pfu* DNA polymerase is dramatically improved in the presence of dUTPase (e.g., increased from < 2 kb to 14 kb<sup>25</sup>). Long-range PCR is particularly susceptible to dUTP poisoning due to the use of prolonged extension times (1–2 minutes per kb at 72°C) that promote dUTP formation.

#### Archaeal DNA Polymerase Fusions

In an effort to increase processivity, various DNA-binding proteins have been fused to the termini of DNA polymerases to increase template binding affinity. For example, fusing the small basic chromatin-like Sulfolobus solfataricus 7d (Sso7d) protein to the C-terminus of Pfu was shown to increase processivity by 8.6-fold<sup>26</sup>. When tested in PCR, the resulting Pfu-Sso7d fusion amplified longer targets in less time compared to native (unfused) Pfu. Several archaeal DNA polymerase fusions have been commercialized that differ with respect to DNA polymerase and/or DNA-binding domain employed, and the inclusion of various PCR-enhancing supplements. For example, the PfuUltra II Fusion HS DNA Polymerase is formulated with a Pfu-based DNA polymerase fused to a proprietary double-stranded DNA binding protein (and supplemented with P. furiosus dUTPase and hotstart antibody; see paragraph below), while Phusion DNA Polymerase consists of a chimeric Deep Vent/Pfu (Pyrococcus sp. GB-D/furiosus) DNA polymerase fused to Sso7d<sup>27</sup>. Fusion DNA polymerases also differ with respect to target-length capability (Table 1); however, all fusions support the use of shorter extension times (15-30 seconds/kb), and thereby provide shorter time-to-results and increased throughput.

#### PCR Characteristics of Proofreading DNA Polymerases

The source, composition, and PCR characteristics of commercial proofreading enzymes are provided in Table 1. PfuUltra and PfuUltra II (fusion) DNA polymerases are formulated with a proprietary Pfu mutant that provides 3-fold higher fidelity than Pfu. In addition, the PfuTurbo and PfuUltra enzymes contain P. furiosus dUTPase (ArchaeMaxx Polymerase Enhancing Factor) to minimize uracil poisoning. As a result, both yield and target-length capability are vastly improved, and genomic targets up to 19 kb in length have been amplified<sup>28,29</sup>. With PfuUltra II fusion HS DNA polymerase, the use of shorter extension times (15 seconds/ kb for < 10 kb targets) means that a 19 kb genomic fragment can be amplified in 5 hours (same-day analysis), instead of > 19 hours (next-day analysis) which is required for non-fusion archaeal DNA polymerases. Other archaeal DNA polymerase formulations that lack dUTPase exhibit comparatively shorter length-capability.

Several proofreading DNA polymerases are available as hotstart formulations. Heat-reversible inactivation is achieved by adding monoclonal antibodies that neutralize polymerase and 3´-5´ exonuclease activities (*PfuUltra* II fusion HS DNA polymerase, Platinum Superfi; no pre-activation required). With proofreading DNA polymerases, high background and/or low product yield may result from extension of non-specifically annealed primers at ambient temperature (common with *Taq*;<sup>30</sup>) or from degradation of primers and DNA template during room-temperature reaction assembly (unique to proofreading enzymes). In our experience, hotstart formulations provide improved yield and/or specificity when amplifying low-copy-number targets in complex backgrounds<sup>31</sup> or longer targets with KOD DNA polymerase (B.Arezi and W. Xing, personal communication).

Each manufacturer recommends somewhat different PCR conditions for optimal performance (Table 1). All manufacturers of proofreading enzymes recommend taking measures to minimize non-specific degradation of PCR primers or products, including using relatively high nucleotide concentrations (200-300 µM each), adding proofreading enzymes last to PCR reactions (after dNTPs), titrating the amount of enzyme, and using sufficient PCR primer concentrations. When testing different proofreading PCR enzymes, researchers are strongly encouraged to follow each manufacturer's recommendation for enzyme amount and extension time. With all proofreading enzymes, synthesizing longer targets or amplifying GC-rich (> 70 %) sequences typically requires additional optimization. In general, amplification of longer targets requires more enzyme units, higher nucleotide concentrations, and/or longer extension times. To enhance amplification of problematic or GC-rich templates, researchers can add DMSO to Pfu formulations (e.g., Herculase II fusion DNA polymerase plus 3–10 % DMSO; titrated in 1 % increments) or use the proprietary PCR additives that are provided with Phusion (GC buffer plus DMSO), Platinum Superfi (PCR, Solution), and DNA polymerases (Table 1).

#### Table 1. Characteristics of High-Fidelity PCR Enzymes<sup>+</sup>

DNA Polymerase	e Exonuclease Activity		Processivity	Polymerization	Uracil	Product Name	Notes and	Recommended	HotStart
(Fusion Domain)	3'-5'	3'-5'	(bases)	Rate (sec-1)	Stalling	(Manufacturer)	Recommendations for Use	Target Length	. lototal t
P. furiosus	Yes	No	10 <sup>15</sup> < 20 <sup>14</sup> 6.4 <sup>26</sup> , 15 <sup>47</sup>	9.3 <sup>15</sup> , 25 <sup>14</sup>	Yes <sup>25</sup> (dU-DNA formuation minimized by ArchaeMaxx factor)	<i>PfuTurbo</i> DNA Polymerase	Pfu PCR buffer optimized for fidelity; Formulated with ArchaeMaxx factor; Genomic < 10 kb: use 2.5 U/50 μl, 200 μM dNTPs and either 1 min/kb (≤ 6 kb) or 2 min/kb (> 6 kb) at 72°C extensions; Genomic > 10 kb: use 5 U/50 μl rxn, 500 μM dNTPs and 2 min/kb at 68°C extensions	Up to 19 kb genomic <sup>29</sup>	Yes*
						<i>PfuUltra</i> DNA Polymerase	Formulated with ArchaeMaxx factor and <i>Pfu</i> mutant that improves fidelity; See <i>PfuTurbo</i> recommendations	Up to 17 kb genomic	Yes*
P. furiosus fusion (double	Yes			ND	Yes	<i>PfuUltra</i> II Fusion HS DNA Polymerase	Formulated with ArchaeMaxx factor, hotstart antibody, and <i>Pfu</i> mutant that improves fidelity; Unique 10X buffer required for optimal activity of fusion; Targets < 10 kb: use 1 µl/50 µl, 250 µM dNTPs, and 15 sec (< 1 kb) or 15 sec/kb (> 1 kb) at 72°C extensions; Targets > 10 kb: use 1 µl/50 µl rxn, 500 µM dNTPs and 30 sec/kb at 68°C extensions	Up to 19 kb genomic	Yes*
(double- stranded DNA binding protein)		No	18547			Herculase II Fusion DNA Polymerase	Formulated with ArchaeMaxx factor; Includes unique 5X buffer and DMSO to enhance PCR of difficult targets; Targets < 12 kb: use 0.5 µl (< 1 kb) or 1 µl (< 1 kb)/50 µl, 250 µM dNTPs and 30 sec (< 1 kb) or 30 sec /kb (> 1 kb) at 72°C extensions; GC-rich targets: add DMSO (0-8 % in 1 % increments) and increase denaturation from 95°C to 98°C	Up to 12 kb genomic	No
P. sp. GB-D	Yes	No	< 2014	2314	Yes <sup>25</sup>	Deep Vent DNA Polymerase (New England BioLabs)	See manufacturer's recommendations	NR	No
P. sp. GB-D/ furiosus chimera fusion (Sso7d)	Yes	No	30-35 <sup>47</sup> (relative processivity: 10X Pfu, 1.6X Taq) <sup>48</sup>	ND	Yes	Phusion DNA Polymerase (ThermoFisher, New England BioLabs); iProof DNA Polymerase (BioRad)	See manufacturer's recommendations	NR	No
T. kodakaraensis KOD1	Yes	No	> 30014	106-13814	Yes <sup>25</sup>	KOD HiFi (Millipore Sigma)	See manufacturer's recommendations	Up to 6 kb	Yes*
Thermus aquaticus	No	Yes	10 <sup>15</sup> , 42 <sup>17</sup>	46.7 <sup>15</sup> , 61 <sup>14</sup>	No	Numerous	See manufacturer's recommendations	Up to 5 kb	Yes

 $\dagger$  Information from product manuals, unless otherwise specified; low-fidelity Taq included for comparative purposes

\* Hotstart formulation contains polymerase- and exonuclease-neutralizing monoclonal antibodies

# Source identified by manufacturer as Pyrococcus sp. strain KOD, but reclassified as T. kodakaraensis  $KOD1^{(44)}$ NR = no recommendations provided by manufacturer; ND = no data; *P. = Pyrococcus, T. = Thermococcus* 

#### **High-Fidelity DNA Polymerase Blends**

In addition to proofreading DNA polymerases, several DNA polymerase blends have been introduced for high-fidelity PCR (Table 2). Commercial DNA polymerase blends consist predominantly of *Taq* plus a lesser amount of a proofreading DNA polymerase (e.g., *Pfu*, Deep Vent) to enhance PCR product yields, amplification of long targets, and fidelity<sup>32</sup>. The fidelity of *Taq*-based blends is typically improved by increasing the proportion of proofreading to non-proofreading DNA polymerase and by modifying the PCR reaction buffer to optimize yield. Since product yield and target-length capability decrease with increasing proofreading: non-proofreading polymerase ratios<sup>32</sup>, higher fidelity *Taq*-based blends typically exhibit reduced performance compared to blends optimized for yield and length (i.e., blends with lower proofreading: non-proofreading polymerase ratios). In general, high-fidelity *Taq*-based blends provide superior performance compared to *Taq* alone with respect to fidelity, yield, and target-length capability (Table 2).

DNA Polymerase		Blend Composition	UstOtestt	Recommended Target		
(Manufacturer)	Major Polymerase	Minor Polymerase	Additives	HOUSTAIL*	Length	
TaqPlus Precision PCR System	Taq	Pfu	None	No	Up to 10 kb genomic and 15 kb vector	
Expand High Fidelity PCR System (Millipore Sigma)	Taq	Tgo	None	No	Up to 5 kb genomic	
Platinum <i>Taq</i> High Fidelity (Invitrogen/ThermoFisher)	Taq	Deep Vent	Taq- neutralizing mAb	Yes (only version available)	Up to 12 kb; up to 20 kb with optimization	
Advantage HF 2 PCR Kit (Takara)	Titanium Taq	Proofreading DNA Polymerase	Taq- neutralizing mAb	Yes (only version available)	Up to 5 kb	

Table 2. Characteristics of High-Fidelity DNA Polymerase Blends<sup>+</sup>

+ Information from manufacturers' catalog or product manual, unless otherwise specified; mAb, monoclonal antibody

\* HotStart formulation contains Pfu- and Taq- neutralizing monoclonal antibodies

# **Protocol/Experimental Methods**

#### **Error Rate Measurements**

DNA polymerase fidelity is expressed in terms of error rate, which corresponds to the number of misincorporated nucleotides per base synthesized. In PCR-based fidelity assays, error rate (E.R.) is calculated as:

> E.R.= number of mutations per bp number of amplicon doublings

where number of amplicon doublings (d) is quantified from the amount of input target DNA and amplicon yield, as:

$$2^{d} = \frac{\text{amplicon yield}}{\text{input target DNA}}$$

The error rates of Pfu and Tag DNA polymerases have been measured using several different methods, including DNA sequencing, denaturing gradient gel electrophoresis (DGGE), and phenotypic forward and reversion mutation assays<sup>1</sup>. Analyses employing direct sequencing or DGGE methods may provide more accurate estimates since all mutations, including silent and lethal mutations are taken into account. However, DNA sequencing is generally impractical for determining error rates of high-fidelity PCR enzymes due to the large number of clones that must be sequenced in order to obtain statistically significant results (e.g., > 23,000 clones must be sequenced to determine the error rate of the PfuUltra enzyme, assuming a mutation rate of 1 per 2.3 x 10<sup>6</sup> bases, 500 bases sequenced per clone, and 5X overage). Moreover, to minimize sequence bias, error rate measurements should employ multiple templates with varying sequence contexts (e.g., GC content, homopolymeric runs, etc.), which further increases cost and labor associated with direct methods. Indirect phenotypic methods are routinely employed by enzyme manufacturers for obvious reasons of simplicity and cost, and underestimates of mutation frequency can be avoided by choosing a well-characterized target gene, such as lacl.

The error rate of *Pfu* DNA polymerase has been estimated at 1.3 x 10<sup>-6</sup> mutations per bp per doubling using a PCR-based phenotypic assay<sup>9</sup> (see *lacl* Phenotypic Mutation Assay below). This is consistent with estimates obtained from DGGE (0.7 x 10<sup>-6</sup> for a 96 bp human mitochondrial sequence<sup>4</sup>; 1.8 x 10<sup>-6</sup> for a 121 bp human APC cDNA sequence<sup>35</sup>) and from DNA sequencing (< 3 x 10<sup>-6</sup>)<sup>36</sup>. At this rate, the probability of a base being mutated in a single round of replication is ~1-3 per 1,500,000 nucleotides, and after 20 doublings

(10<sup>6</sup>-fold amplification), ~1-2.5 % of 1 kb amplification products will contain mutations. In comparison, published error rates for *Taq* range from 0.5-21 x 10<sup>-5</sup> mutations per bp per doubling, and include: 7.2-21 x 10<sup>-5</sup> using DGGE<sup>12, 37</sup> 0.8-1.0 x 10<sup>-5</sup> (*lacl*) and 1.8 x 10<sup>-5</sup> (p53) using PCR-based phenotypic assays<sup>9,38,39</sup>, 2 x 10<sup>-5</sup> using a gap-filling *lacZ* assay<sup>10</sup>, and 0.5-2.7 x 10<sup>-5</sup> by DNA sequencing of PCR products<sup>36,40</sup>. At these rates, anywhere from 10 % to 100 % of 1 kb products amplified with *Taq* will contain one or more mutations (doublings = 20; mutation-containing products = 10-420 %).

Variation in published error rates reflects differences in the reaction conditions (e.g., pH, [dNTPs], [Mg<sup>2+</sup>], DNA template sequence) and types of fidelity assays employed<sup>1,11,12</sup>. Because different assays are likely to measure different parameters, error rates should only be compared among PCR enzymes tested in the same assay<sup>13</sup>, and preferably, according to manufacturers' recommendations.

#### lacl Phenotypic Mutation Assay

Our laboratory routinely employs a PCR-based forward mutation assay that utilizes the well-characterized lacl target gene<sup>9,38</sup>. In this assay, a 1.9 kb sequence encoding *lac*IOZ $\alpha$  is amplified and cloned, and the percentage of clones containing a mutation in lacl (% blue) is determined in a color-screening assay (Figure 1). To accurately determine mutation rates with a phenotypic assay, it is essential that the number of base changes producing a scorable mutant phenotype is known. Otherwise, mutation rates can be greatly underestimated by not taking into account silent mutations that alter DNA sequence without producing a change in protein sequence or function. The sensitivity of lacl to mutation is well known. More than 30,000 lacl mutants have been sequenced, and the results indicate that 349 single-base substitutions occurring at 179 amino acid positions in the 1080 bp lacl-coding region can be identified by color screening<sup>41</sup>.

Therefore, in the *lacl* assay, error rates are calculated as mutation frequency per 349 bp per duplication:

E.R.= <u>lacl-mutant frequency</u> (349 bases) (d)

where (d) = the number of amplicon doublings

We have measured the error rates of several DNA polymerases using the *lacl* assay (Table 3). Error rates were measured in each enzyme's recommended PCR buffer, and whenever possible, identical PCR conditions were used,

including DNA template concentration, PCR cycling parameters, and number of PCR cycles performed. The only exceptions were that each manufacturer's recommendations were followed with respect to number of enzyme units, nucleotide concentration, primer concentration, extension temperature, and extension time (shorter times were employed with fusion enzymes) (Table 3). To allow assay-toassay comparisons, *Pfu* DNA polymerase was run in every assay, and error rates were normalized relative to the mean value of  $1.3 \times 10^{-6}$  mutations per bp per doubling as determined for *Pfu* in study #1<sup>9</sup>. *Taq* DNA polymerase, serving as a second internal control, exhibited mean error rates of  $8.0 \times 10^{-6}$  (study #1; 11 PCRs) and 9.1 x 10<sup>-6</sup> (mean of studies #2-5; 14 PCRs) mutations per bp per doubling.

#### Results

As expected, proofreading DNA polymerases exhibited significantly lower error rates (1-3 errors per 10<sup>6</sup> bases) compared to *Taq* DNA polymerase (8-9 errors per 10<sup>6</sup> bases). The *PfuUltra* mutant DNA polymerase (non-fusion and fusion) formulations exhibited error rates (4 x 10<sup>-7</sup> mutations per bp per duplication) that were 3-fold lower than the error rates of *Pfu* and Phusion DNA polymerases. Relative differences in error rate observed with the *lacl* assay (Table 3) are consistent with those obtained using a p53-based forward mutation assay (e.g., *Pfu* < Taq<sup>39</sup>) and DGGE (e.g., *Pfu* < *Taq*<sup>1</sup>). In general, the error rates of high-fidelity DNA polymerase blends (3-6 errors per 10<sup>6</sup> bases) are intermediate between proofreading DNA polymerases and *Taq* (Table 3).

The use of high-fidelity DNA polymerases, especially those that support fast cycling, becomes increasingly important as amplicon size increases (Table 3). With Tag, the percentage of clones expected to contain mutations in a 106-fold amplification reaction increases from 4 % (for 250 bp amplicon) to 16 % (1 kb amplicon) to 80 % (5 kb amplicon), while the number of clones that should be sequenced to obtain an error-free clone (95 % confidence) increases from 1 to 2 to 14, respectively  $(0.95=1-(1-f)^n)$ , where f = frequency of error-free clones and  $n = number of clones sequenced^{42}$ ). When amplifying a broader range of targets (0.25 to 10 kb) with high-fidelity blends (E.R.=2.8-5.8 x 10<sup>-6</sup>), the percentage of clones likely to contain mutations increases from 1-3 % (250 bp amplicon) to 5–11 % (1 kb amplicon) to 28–58 % (5 kb amplicon) to 56-100 % (10 kb amplicon), and the number of clones that should be sequenced increases from 1-2 (up to 1 kb amplicon) to 3-5 (5 kb amplicon) to > 6 (10 kb amplicon). When amplifying similarly sized targets with the *PfuUltra* enzyme (E.R.= $4 \times 10^{-7}$ ), the frequency of error-containing clones is: < 1 % (up to 1 kb amplicon), 4 % (5 kb amplicon), and 8 % (10 kb amplicon), and sequencing 1 (up to 6 kb amplicon) or 2 (6–10 kb amplicon) clones should be sufficient for identifying an error-free clone. In addition, with the faster *PfuUltra* II fusion HS DNA polymerase, long fragments can be amplified with the same degree of accuracy in a fraction of the time; for example, a 5 or 10 kb fragment can be amplified with PfuUltra II enzyme in 1 or 3 hours respectively, compared to 3 or 10 hours required for amplification by a non-fusion high-fidelity PCR enzyme.

#### Table 3. Error Rates of High-Fidelity PCR Enzymes.

DNA Delementer	Number	Number	Error Rate# (x 10 <sup>-6</sup> ± S.D.)	Accuracy	Percentage of Clones with Mutations (10 <sup>6</sup> -fold Amplification)			
DNA Polymerase	of Studies	of PCRs		in Bases)	1 kb Amplicon	5 kb Amplicon	10 kb Amplicon	
Proofreading DNA Polymerases								
PfuUltra II Fusion HS DNA Polymerase	6	8	0.4 ± 0.06	2,500,000	0.8	4	8	
PfuUltra DNA Polymerase	5	12	0.4 ± 0.04	2,500,000	0.8	4	8	
Pfu DNA Polymerase	1	10	1.3 ± 0.2 <sup>9, 28</sup>	770,000	2.6	13	26	
Herculase II Fusion DNA Polymerase	6	6	1.3 ± 0.2	770,000	2.6	13	26	
Phusion DNA Polymerase/ iProof DNA Polymerase	6	5	1.3 + 0.4	770,000	2.6	13	26	
Deep Vent DNA Polymerase	1	4	2.7 ± 0.2°	370,000	5.4	NR	NR	
High Fidelity Blends								
TaqPlus Precision PCR System	2-3	13	4.0 ± 1.3 <sup>33</sup>	250,000	8	40	80	
Platinum Taq High Fidelity	3	2	5.8 ± 0.3 <sup>33</sup>	170,000	11.6	58	100	
Advantage-HF	3	2	6.1 ± 0.0 <sup>33</sup>	160,000	12.2	NR	NR	
	1	11	8.0 ± 3.9 <sup>9</sup>	125,000	16	80	NR	
rad DNA Polymerase	2-5	14	9.1 ± 2.4	110,000	18.2	91	NR	

NR = not recommended for 5 to 10 kb target sizes

<sup>#</sup>Error rates were measured in each enzyme's recommended PCR buffer. Cycling conditions were described in<sup>9</sup>, or were as follows: (*Taq*, *PfuUltra* DNA polymerases): 95°C 1 min. (1 cycle); 95°C 30 sec, 58°C 30 sec, 72°C 6 min. (30 cycles); 72°C 10 min. (1 cycle); (*PfuUltra* II, Herculase II, Phusion DNA polymerases): 95°C 1 min. (1 cycle); 95°C 30 sec, 58°C 30 sec, 72°C 45 sec (30 cycles); 72°C 5 min. (1 cycle); PCR reactions (50 µl) contained 0.2 µM each primer, 200 µM each nucleotide, 2.5 ng target DNA, and 2.5 U DNA polymerase, with the following manufacturer-recommended exceptions: Platinum Pfx- 300 µM each nucleotide, 1.25 U enzyme, and 68°C extension temperature; Deep Vent 1 U enzyme; *PfuUltra* II and Herculase II DNA polymerases-1 µl; Phusion-1 U  $\label{eq:constraint} \textbf{Table 4.} \ \text{Polymerase ordering guide.} \ \text{Volume can be customized to your needs.}$ 

PCR Enzyme	Fidelity	Speed	Yield	Target Length (genomic DNA)	Sensitivity			
High-Fidelity & Difficult/GC Rich PCR								
PfuUltra II Fusion HotStart DNA Polymerase Engineered to be the highest fidelity and fastest polymerase available	1 error/2.5 million bp	15 sec/kb		0-19 kb				
Herculase II Fusion DNA Polymerase High-fidelity polymerase for difficult targets. Provides superior yields over a broad range of targets. Economical enough for routine use	1 error/770,000 bp	15 sec/kb		"0-12kb 12-23 kb (optimized)"				
PfuUltra High-Fidelity DNA Polymerase AD Engineered for high-fidelity	1 error/2.25 million bp	1 min/kb		19 kb (optimized)				
<b>PfuTurbo DNA Polymerase AD</b> First high-fidelity polymerase to include the ArchaeMaxx Polymerase-Enhancing factor	1 error/770,000 bp	1 min/kb		19 kb (optimized)				
Herculase Enhanced DNA Polymerase Designed for difficult targets	1 error/375,000 bp	1 min/kb		12 kb				
Cloned Pfu DNA Polymerase AD Cloned to ensure ultrapure manufacturing of Pfu	1 error/770,000 bp	2 min/kb		"1 kb 5 kb (optimized)"				
<i>Pfu</i> DNA Polymerase Stratagene introduced the first thermophilic proofreading polymerase	1 error/770,000 bp	2 min/kb		(up to 1 kb)				
Specialty Enzymes	Specialty Enzymes							
<b>PfuTurbo Cx HotStart DNA Polymerase</b> The only high-fidelity polymerase that can read through dUTP in the template and extending strand	1 error/770,000 bp	1 min/kb		0-10 kb				
PicoMaxx High-Fidelity PCR System Most sensitive polymerase offered	2x Taq	1 min/kb		0-10 kb				
Easy-A High-Fidelity PCR Cloning Enzyme Proofreading DNA polymerase that adds 3'A overhangs to PCR amplicons	1 error/770,000 bp	1 min/kb		0-6 kb				
Routine Enzymes	Routine Enzymes							
Paq5000 DNA Polymerase Fast and economical alternative to Taq		30 sec/kb		0-6 kb				
Taq2000 DNA Polymerase Ultrapure cloned Taq that eliminates unwanted background artifacts		1 min/kb		"1 kb 4 kb (optimized)"				
Taq DNA Polymerase First thermophilic PCR enzyme.		1 min/kb		"1 kb 4 kb (optimized)"				

Blunt or 3'-A Ends	ArchaeMaxx Advantage	Enzyme Only		HotStart		Master Mix	PCR Enzyme	
		100 U 1000 U	500 U 5000 U	100 U 1000 U	500 U 5000 U	100 rxn 400 rxn		
High-Fidelity & Difficult/GC Rich PCR								
Blunt	ArchaeMaxx Advantage			(40 rxn) 600670 (400 rxn) 600674	(200 rxn) 600672 –	600850 600852	PfuUltra II Fusion HotStart DNA Polymerase	
Blunt	ArchaeMaxx Advantage	(40 rxn) 600675 (400 rxn) 600679	(200 rxn) 600677 –				Herculase II Fusion DNA Polymerase	
Blunt	ArchaeMaxx Advantage	600385 600389	600387	600390 600394	600392	600630 _	PfuUltra High-Fidelity DNA Polymerase AD	
Blunt	ArchaeMaxx Advantage	600255 600259	600257	600320 600324	600322		PfuTurbo DNA Polymerase AD	
Mixed	ArchaeMaxx Advantage	600260 600264	600262 600266	600310 600314	600312		Herculase Enhanced DNA Polymerase	
Blunt		600353 600357	600355				Cloned Pfu DNA Polymerase AD	
Blunt		600135 600140	600136				Native Pfu DNA Polymerase	
Specialty Enzymes								
Blunt	Alternative uracil resistance (Pfu mutation)			600410 600414	600412		PfuTurbo Cx HotStart DNA Polymerase	
Mixed	ArchaeMaxx Advantage			600420 600424	600422	600650	PicoMaxx High-Fidelity PCR System	
3'-A	ArchaeMaxx Advantage			600400 600404	600402	600640 600642	Easy-A High-Fidelity PCR Cloning Enzyme	
Routine Enzymes								
Mixed	ArchaeMaxx Advantage	600682	600680 600684			600870 600872	Paq5000 DNA Polymerase	
3'-A		600195 600197	600196	600280 600284	600282		Taq2000 DNA Polymerase	
3'-A							Taq DNA Polymerase	

 Table 5. Polymerase ordering guide. Reagents can be customized to your needs.

#### **Optimizing PCR Fidelity**

PCR error rate can be minimized by employing the highest fidelity PCR enzyme available for the desired application. As discussed above, commercial high-fidelity DNA polymerases show considerable variation in error rates, ranging from 0.4-3.5 x 10<sup>-6</sup> for proofreading DNA polymerases, up to 2.8-6.1 x 10<sup>-6</sup> for DNA polymerase blends (Table 3). However, when selecting a PCR enzyme, parameters other than fidelity may have to be considered. Current high-fidelity PCR enzymes are incompatible with dUTP/UNG decontamination<sup>22,23</sup> and direct TA cloning methods<sup>19</sup>. However, post-amplification addition of 3' A overhangs with Tag improves the TA cloning efficiency of bluntended fragments amplified with proofreading enzymes. (Post-amplification A-addition requires incubation for 8-10 minutes at 72°C; see the StrataClone PCR Cloning Kit instruction manual for details.) Alternatively, researchers can generate amplicons with 3' A overhangs using the Easy-A DNA Polymerase, a proprietary PCR enzyme with Pfu-like fidelity. Thus, suitable high-fidelity enzyme formulations are available for nearly every PCR application.

In addition to enzyme choice, researchers should also consider optimizing reaction conditions to further reduce PCR mutation frequency. While error rate is an intrinsic property of DNA polymerases (under defined reaction conditions), observed mutation frequencies can vary from PCR to PCR, depending on the number of amplicon doublings. For example, assuming we amplify a 1 kb fragment using Tag (E.R., 8 x 10<sup>-6</sup> mutations per bp per doubling), a PCR generating 5 µg of amplicon from 5 pg of target DNA has undergone 20 target doublings and produced 1.6 mutations per 10,000 bases (~3/20 clones with mutations). In comparison, a PCR generating 5 µg of amplicon from 75 ng target DNA has undergone 6 target doublings (67-fold amplification) and introduced 0.5 mutations per 10,000 bases (~1/20 clones with mutations). Therefore, researchers can minimize mutation frequency by limiting the number of target duplications, for example, by increasing the amount of input DNA template or reducing the number of PCR cycles.

Additional reductions in mutation frequency may be achieved by optimizing buffer composition, nucleotide concentration, or polymerase amount. As discussed above, the error rates shown in Table 3 were obtained using the PCR buffer and nucleotide concentration recommended by each manufacturer, which may or may not be optimal with respect to fidelity. High-fidelity PCR reaction conditions have been developed for Tag, Deep Vent, and Pfu DNA polymerases<sup>9-12,43</sup>. For example, the error rate of Pfu decreases from 2.6- to 1.1x 10<sup>-6</sup> as the nucleotide concentration is lowered from 1 mM to 100 µM each<sup>9</sup>. Even greater changes in *Pfu's* error rate were observed as the Mg<sup>2+</sup> concentration was increased from 1 mM ( $4.9 \times 10^{-6}$ ) to 2 mM MgSO<sub>4</sub> ( $1.3 \times 10^{-6}$ ) (at 200  $\mu$ M each dNTP, pH 8.8) and the pH was increased from pH 7.5 (8.2 x 10-6) to pH 8.8 (1.3 x 10-6) (at 200 µM each dNTP, 2 mM MgSO<sub>4</sub>)<sup>9</sup>. For enzymes whose pH and Mg<sup>2+</sup> optima are unknown, researchers can expect to achieve lower mutation frequencies by using the lowest balanced nucleotide concentration compatible with yield (e.g., 25–150 µM each). In addition, using lower enzyme concentrations is also likely to minimize polymerase extension from mispaired or misaligned primer termini<sup>11</sup>.

# Conclusion

Since the introduction of *Taq* DNA polymerase in the late 1980s, significant progress has been made in developing PCR enzyme formulations with improved fidelity, PCR performance, and speed. Proofreading DNA polymerases offer significantly higher fidelity compared to *Taq*, and initial problems associated with their use (low yield, unreliability, speed) have been largely overcome by reducing uracil poisoning (the *Pfu* formulations), preparing blends with *Taq* DNA polymerase, and developing faster, more processive proofreading DNA polymerase fusions. In fact, Agilent's new high-fidelity enzyme formulations provide significantly improved yield, throughput, and target-length capability compared to *Taq*.

### References

- Cha, R.S. and Thilly, W.G. (1995) PCR Primer: A Laboratory Manual (eds, Dieffenbach, C.W. and G.S. Dveksler) Cold Spring Harbor Laboratory Press
- 2. Frohman, M.A., Dush, M.K., and Martin, G. (1988) *Proc. Natl. Acad. Sci.* 85:8998-9002
- 3. Li, H., Cui, X., and Arnheim, N. (1990) Proc. Natl. Acad. Sci. 87:4580-4584
- 4. Andre, P., et al. (1997) *Genome Res.* 7:843-852
- 5. Kunkel, T.A. (1992) J. Biol. Chem. 267:18251-18254
- 6. Goodman, M.F., et al. (1993) Crit. Rev. Biochem. Mol. Biol. 28:83-126
- 7. Goodman, M.F., and Fygenson, K.D. (1998) *Genetics* 148:1475-1482
- 8. Kunkel, T.A., and Bebenek, K. (2000) Annu. Rev. Biochem. 69:497-529
- 9. Cline, J., Braman, J.C., and Hogrefe, H.H. (1996) *Nucl. Acids Res.* 24:3546-3551
- 10. Eckert, K.A. and Kunkel, T.A. (1990) Nucleic Acids Res. 18:3739-3744
- 11. Eckert, K.A. and Kunkel, T.A. (1991) PCR Methods Appl. 1:17-24
- 12. Ling, L.L., et al. (1991) PCR Methods and Applications 1:63-69
- 13. Perler, F.B., Kumar, S., and Kong, H. (1996) Adv. Protein Chem. 48:377-435
- 14. Takagi, M., M., et al. (1997) Appl. Environ. Microbiol. 63:4504-4510
- 15. Hogrefe, H.H., et al. (2001) *Methods of Enzymology* (eds. M.W.W. Adams and R.M. Kelly) Academic Press, N.Y., vol. 334, pg. 91-116
- 16. Lyamichev, V., Brow, M.A., and Dahlberg, J.E. (1993) Science 260:778-783
- 17. Kong, H., Kucera, R.B., and Jack, W.E. (1993) J. Biol. Chem. 268:1965-1975
- 18. Hu, G. (1993) DNA Cell Biol. 12: 763-770
- 19. Zhou, M.Y. and Gomez-Sanchez, C.E. (2000) Universal TA cloning. Curr. Issues Mol. Biol. 2:1-7
- 20. Costa, G. L.,and Weiner, M. P. (1994) PCR Methods Appl. 3:95-106
- 21. Greagg, M.A., et al. (1999) Proc. Natl. Acad. Sci. 96:9045-9050
- 22. Longo, M.C., Berninger, M.S., and Hartley, J.L. (1990) Gene 93:125-128
- 23. Slupphaug, G., et al. (1993) Anal. Biochem. 211:164-169
- 24. Sakaguchi, A.Y., et al. (1996) BioTechnniques 21:368-369
- 25. Hogrefe, H. H., et al. (2002) Proc. Natl. Acad. Sci. 99:596-601
- 26. Wang, Y., et al (2004) Nucl. Acids Res. 32:1197-1207
- Finney, M.J. (2004) "Thermostable Enzymes from a Randomized Hybrid Library". Presented at Extremophiles 2004: 5th International Conference on Extremophiles. Cambridge, MA. (September 19-23, 2004)
- 28. Borns, M., Cline, J., and Hogrefe, H. (2000) Strategies 13(1):27-30
- 29. Hogrefe, H., et al. (1997) Strategies 10(3):93-96
- 30. Kellogg, D.E., et al. (1994) BioTechniques 16:1134-1137

- 31. Borns, M., Scott, B., and Hogrefe, H. (2001) Strategies 14(1):5-8
- 32. Barnes, W. M. (1994) Proc. Natl. Acad. Sci. 91:2216-2220
- 33. Borns, M., and Hogrefe, H. (2000) Strategies 13(1):1-3
- 34. Borns, M., and Hogrefe, H. (2000) Strategies 13(2):76-79
- 35. Li-Sucholeiki, X.C. and Thilly, W.G. (2000) Nucl. Acids. Res. 28:E44
- 36. Bracho, M.A., Moya, A. and Barrio, E. (1998) J. Gen. Virol. 79:2921-2928
- 37. Keohavong, P., and Thilly, W. G. (1989) Proc. Natl. Acad. Sci. 23:9253-7
- 38. Lundberg, K.S., et al. (1991) Gene 180:1-6
- 39. Flaman, J.-M., et al. (1994) Nucleic Acids Res. 22:3259-3260
- 40. Martell, M., et al. (1992) J. Virol. 66:3225-3229
- 41. Provost, G.S., et al. (1993) Mut. Res. 288:133-149
- 42. Jeltsch, A. and Lanio, T. (2002) *In Vitro Mutagenesis Protocols* (ed, Braman, J.) Humana Press, second edition
- 43. Mattila, P., et al. (1991) Nucl. Acids Res. 19:4967-4973
- 44. Nishioka, M., H., et al. (2001) J. Biotechnology 88:141-149
- 45. Expressions 8.5:13 (Invitrogen)
- 46. Qiagen Technical Services
- 47. Strategies (2006) 19(1)
- 48. Phusion<sup>™</sup> High Fidelity DNA Polymerase brochure (FinnZymes) Reproduced with permission from Cold Spring Harbor Laboratory Press.<sup>©</sup> 2003

### Legal

- Deep Vent is a registered trademark of New England Biolabs
- Platinum is a registered trademark of ThermoFisher/Invitrogen
- Phusion is a trademark of Finnzymes Oy
- iProof is a trademark of BioRad Laboratories

#### www.agilent.com/genomics/lifescience

For Research Use Only. Not for use in diagnostic procedures.

This information is subject to change without notice.

PR7000-2235 © Agilent Technologies, Inc. 2019 Printed in the USA, May, 2019 5994-0929EN

