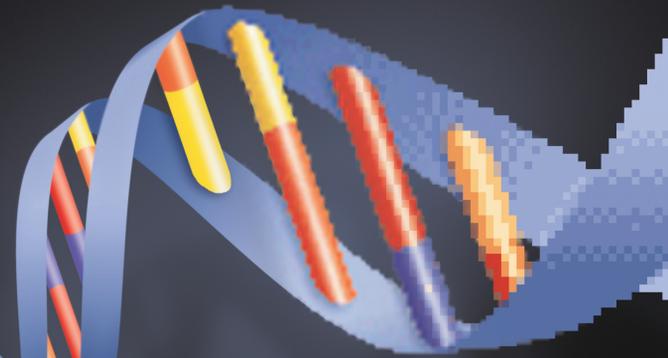


PCR

POLYMERASE CHAIN REACTION



The Cell is where it all starts. The library of genetic information is contained here.

Chromosomes are bundles of DNA that carry the individual sequences that we're interested in.

The PCR Tube contains the DNA and all of the chemicals necessary for the reaction to occur.

A DNA Sequence of interest is what we're looking for, and it's just a tiny fraction of the entire chromosome.

The PCR Revolution

“Beginning with a single molecule of the genetic material DNA, PCR can generate 100 billion similar molecules in an afternoon. The reaction is easy to execute. It requires no more than a test tube, a few simple reagents, and a source of heat.”

DNA is the molecule that guides every aspect of life. It contains the genetic code, or blueprint, needed to make precisely one of you, a blade of grass, or a worm, and the exact same DNA is found in almost every cell in your body. DNA is composed of four basic repeating units called nucleotides or bases. The nucleotides are represented by the first letters of their names: A (adenine), G (guanine), T (thymine), and C (cytosine). In total, each cell in your body contains about 3 billion base pairs of DNA.

PCR produces exponentially large amounts of a target piece of DNA from trace amounts of starting material (template). The template can be any form of DNA. A researcher can take trace amounts of DNA from a drop of blood, a single hair follicle, a flower petal, or a virus particle and use PCR to generate billions of copies of a desired DNA fragment. In theory, only a single intact strand of template DNA is needed to generate millions of new DNA molecules via PCR. Before PCR was invented, it was extremely difficult and often impossible to do forensic or genetic studies with such a small amount of DNA. PCR has become one of the most important and most commonly used tools in molecular biology.

PCR relies on thermal cycling, during which cycles of heating and cooling cause the double-stranded DNA to separate (denature), followed by binding (annealing) of the primers and elongation (extension) of the DNA from the primer position. These three steps make up one PCR cycle. The DNA copies that are created at each cycle become templates for successive cycles, and a chain reaction resulting in exponential DNA amplification is set in motion.

Denaturation + Annealing + Extension = Complete Cycle of PCR

Denaturation Step / The sample is heated to 94°C, which causes the hydrogen bonds between base pairs to separate (or melt).

Annealing Step / The temperature is lowered to 50–65°C to allow the binding of primers, short fragments of DNA (usually 15–30 bases in length). The primers are complementary to the DNA sequences flanking the target sequence. The precise temperature for annealing depends on the length and sequence of the primer.

Extension Step / The temperature is increased to 72°C to speed up the polymerase activity. The DNA polymerase (for example, *Taq* polymerase) adds dNTPs (deoxynucleotide triphosphates — dATP, dTTP, dCTP, and dGTP) onto the ends of primers. Extension follows the base pairing rule and proceeds in the 5' to 3' direction.

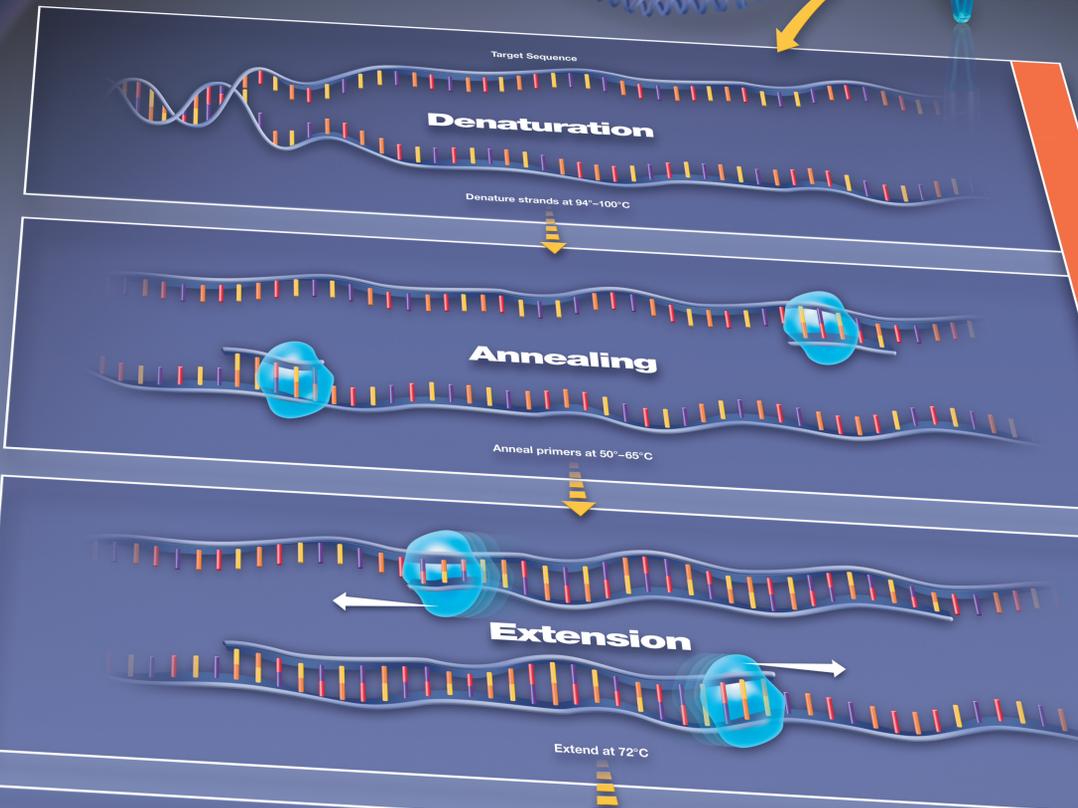
PCR generates DNA of a precise length and sequence. On the first cycle, the two primers anneal to the original genomic template DNA strands. After the first complete temperature cycle, two new strands are generated that are shorter than the original template strands but still longer than the target DNA. It isn't until the third thermal cycle that fragments of the precise length are generated.

Cycle 40 / It is the template strands of the precise length that are amplified exponentially (X^n , where X = the number of original template strands and n = the number of cycles). After just 30 cycles, there would be over 1 billion (1,073,741,824) copies of every original target sequence.

The DNA polymerase used in PCR must be thermally stable because the polymerase chain reaction cycles between temperatures in the range of 50°C to 100°C. The first and most commonly used thermostable DNA polymerase was isolated from a thermophilic bacterium, *Thermus aquaticus* (*Taq*), which lives in high-temperature steam vents such as those found in Yellowstone National Park.



In 1983, Kary Mullis, then working at Cetus Corporation, developed the molecular biology technique that revolutionized genetic research and earned him the Nobel Prize in 1993. PCR transformed molecular biology into a multidisciplinary research tool. PCR has had a tremendous impact on these areas of biotechnology: gene mapping, cloning, DNA sequencing, gene expression, gene detection, and DNA profiling. In fact, PCR is now routinely used in diverse applications such as genomic studies, forensics, medical research and diagnostics, and evolutionary and ecological studies.



Three Step Cycle Repeats

Cycle 2

Cycle 5

Cycle 40

Four Copies

Thirty-two Copies

Billions of Copies of the Target Sequence

Applications

Genomics / The Human Genome Project (HGP) was one of the greatest scientific endeavors in history. Completed in 2003, it took a coordinated international research effort to sequence and map the complete human genome. One surprising outcome of the HGP was just how few genes are encoded in the human genome, and how much of the genome is made up of noncoding sequences. Knowing the sequence is just the beginning.



Forensics / Forensic sciences describe the boundary between science and the law. Science can as easily convict someone of a crime as free someone wrongly convicted. Today, molecular genetic methods are used to determine the exact genotype of a DNA sample. Using PCR, it is possible to analyze DNA from the tiniest of biological samples and to distinguish any two people on the planet (with the exception of identical twins), living or dead.



Medicine / With increased understanding of molecular genetics, it has become clear that many diseases have a genetic component. Certain genes are linked to diseases such as sickle cell anemia or breast cancer (BRCA1 gene). In other cases, environmental damage (UV light, cigarette smoke) can cause mutations in DNA that lead to cancer. In the future, our unique genetic makeup will determine which treatments will be most effective in treating our diseases.



Evolution / Why is there a duck-billed platypus (*Ornithorhynchus anatinus*)? This unusual animal — which both lays eggs and has mammary glands — appears to be an evolutionary bridge between mammals and reptiles. Genomic studies provide molecular evidence of the genetic divergence between mammals and their reptilian relatives. The duck-billed platypus data have been used to deepen our understanding of developmental biology, comparative physiology, and neurobiology.



Ecology / Research groups like NOAA's National Marine Fisheries Service use variation in molecular genetic markers to understand the migratory and dispersal patterns of species like the beluga whale. The genetic data provide tools to understand the breeding behavior and social organization of the species. By combining genetic data with traditional ecological information, NOAA researchers are better able to design species management plans.



explorer.bio-rad.com

BIO-RAD

