

## EpiQuik Red Taq DNA Polymerase

(Catalog No. R12011)

### Description

EpiQuik Red Taq DNA Polymerase has a special formulation with added inert red dye. This makes it very suitable for standard applications. Strong red color of the enzyme allows user to check the polymerase addition and verify adequate mixing. Reaction products are ready for direct gel loading and the dye serves as a marker for electrophoresis progress monitoring. EpiQuik Red Taq DNA Polymerase has 5'→3' DNA synthesis activity.

### Concentration

1 unit/μl (One unit of the enzyme catalyzes the incorporation of 10 nanomoles of deoxyribonucleotides into a polynucleotide fraction in 30 min at 70°C).

### Composition

- EpiQuik Red Taq DNA Polymerase in Storage Buffer: 20 mM Tris-HCl (pH 8.0), 1mM DTT, 0.1 mM EDTA, 100 mM KCl, 0.5% Nonidet P40, 0.5% Tween 20 and 50% glycerol
- 10x Reaction Buffer: 100 mM Tris-HCl (pH 8.8 at 25°C), 500 mM KCl, 0.8% Nonidet P40
- 10x Reaction Buffer with (NH<sub>4</sub>)<sub>2</sub>SO<sub>4</sub>: 750mM Tris-HCl (pH 8.8 at 25°C), 200 mM (NH<sub>4</sub>)<sub>2</sub>SO<sub>4</sub>, 0.1% Tween 20
- 25 mM MgCl<sub>2</sub> Solution

### Quality Data

Activity and stability tested at 20, 30 and 40 cycles of PCR reactions at 95°C. Tested for the absence of human DNA contamination by PCR with Alu-specific primers.

### Applications

- Suited for a wide range of PCR assays
- Easy visualization of enzyme addition
- Visualization of complete reaction mixing
- Direct loading of samples following amplification

### Recommended PCR Reaction Mix

| Component  | Quantity            |
|--|---------------------|
| EpiQuik Red Taq (1 U/μl)   | 1.25-2.5 U          |
| 10x Reaction Buffer (or with (NH <sub>4</sub> ) <sub>2</sub> SO <sub>4</sub> ) | 5 μl (1x)           |
| 25 mM MgCl <sub>2</sub>  | 3-5 μl (1.5-2.5 mM) |
| 10 mM dNTP mix   | 1 μl (200 μM)       |
| Primer Forward   | 0.3 -1 μM           |
| Primer Reverse   | 0.3 -1 μM           |
| DNA template   | 1-100 ng/μl         |
| H <sub>2</sub> O PCR grade   | Up to 50 μl         |
| <b>Total</b>   | <b>50 μl</b>        |

*This product is for research purposes only. Not intended for use in diagnostic procedures.*

### Recommended PCR Cycles

| <u>Cycle step</u>    | <u>Temp.</u> | <u>Time</u> | <u>Cycles</u> |
|----------------------|--------------|-------------|---------------|
| Initial denaturation | 95°C         | 3-5 min     | 1             |
| Denaturation         | 95°C         | 30-60 s     | 26-35         |
| Annealing            | 50-68°C      | 30-60 s     | 26-35         |
| Elongation           | 72°C         | 1-4 min     | 26-35         |
| Final elongation     | 72°C         | 5-10 min    | 1             |

IMPORTANT: Annealing temperature should be 2-6°C lower than the primer melting temperature.

### Safety Warnings and Precautions

This product is designed for research purposes and in vitro use only. According to common laboratory safety practice, it is recommended to wear protective clothing, gloves and safety glasses. Please refer to [www.epigentek.com](http://www.epigentek.com) for Material Safety Data Sheet of the product. Some applications this product is used in may require a license which is not provided by the purchase of this product. Users should obtain the license if required.

### Storage Conditions

Store at -20°C. Guaranteed stable for 12 months when properly stored.

### Ordering Information

| <b>Products</b>                | <b>Size</b> | <b>Cat. No.</b> |
|--------------------------------|-------------|-----------------|
| EpiQuik Red Taq DNA Polymerase | 500 Units   | R12011-1        |
| EpiQuik Red Taq DNA Polymerase | 1000 Units  | R12011-2        |
| EpiQuik Red Taq DNA Polymerase | 2500 Units  | R12011-3        |