

## Product Information

### REDTaq® Genomic DNA Polymerase with MgCl<sub>2</sub>

Catalog Number **D8312**  
Storage Temperature –20 °C

## TECHNICAL BULLETIN

### Product Description

REDTaq Genomic DNA Polymerase is Sigma's high quality Taq DNA Polymerase mixed with an inert red dye designed to provide an enhanced amplification of more complex or genomic templates. The dye allows quick recognition of reactions to which enzyme has been added, as well as visual confirmation of complete mixing. The enzyme is provided at 1 unit/μL for more accurate volume measurement and less waste.

Reactions using REDTaq Genomic DNA polymerase and 10× PCR reaction buffer are formulated as any PCR mixtures when optimizing individual components. There are no additional preparation steps or protocol changes required. The formulation allows aliquots (5–10 μL) from the PCR to be directly loaded onto an agarose gel without addition of electrophoresis loading buffers. The dye migrates at the same rate as a 125 bp fragment in a 1% agarose gel. Because a gel loading buffer is not added to the reaction mix, a sample can be re-amplified, such as in nested PCR.

If desired, the dye can be removed from the amplicon by any standard purification method. The presence of the dye has no effect on manual or automated DNA sequencing, ligation, and transformations. Though exceptions may exist, the dye is generally inert in restriction enzyme digestions.

**Unit Definition:** One unit incorporates 10 nmol of total deoxyribonucleoside-triphosphates into acid precipitable DNA in 30 min at 74 °C.

### Reagents provided

- REDTaq Genomic DNA Polymerase, Catalog Number D0688.  
1 unit/μL in 20 mM Tris-HCl, pH 8.0, 100 mM KCl, 0.1mM EDTA, 1 mM DTT, 0.5% TWEEN® 20, inert dye, 50% glycerol. Provided as 50, 250, 1,000 or 2,500 units
- 10× PCR Reaction Buffer, Catalog Number P2192, 100 mM Tris-HCl, pH 8.3, 500 mM KCl, 15 mM MgCl<sub>2</sub> and 0.01% gelatin. Provided as 1.5 ml vial, 1 vial per 50, or 250 units of REDTaq, 4 vials per 1,000 units of REDTaq, or 10 vials per 2,500 units of REDTaq.

### Equipment and reagents required but not provided

- Deoxynucleotide Mix, Catalog Number D7295  
10 mM each dATP, dCTP, dGTP, TTP
- Water, PCR Reagent, Catalog Number W1754
- Mineral Oil, Catalog Number M8662 (optional)
- Primers
- DNA template
- Dedicated pipettes
- PCR pipette tips
- 0.5 ml or 0.2 ml thin-walled PCR tubes, Catalog Numbers P3114 and P3364
- Thermal cycler

### Storage

Store all components at –20 °C.

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

### Procedure

Optimal concentrations of template DNA, MgCl<sub>2</sub>, KCl, and PCR adjuncts as well as pH are often target specific. It may be necessary to determine the optimal concentration of each component.

1. Add the following reagents to a 0.2 ml or 0.5 ml PCR tube. A master mix is highly recommended when performing multiple PCR reactions.

Volume	Reagent	Final Concentration
5 µL	10× PCR Buffer	1 ×
1 µL	Deoxynucleotide Mix	200 µM each dNTP
- µL	Primer	0.1-0.5 µM
- µL	Primer	0.1-0.5 µM
2.5 µL	REDTaq Genomic DNA Polymerase	0.05 unit/µL
- µL	Template DNA (typically 10 ng)	200 pg/µL
q.s.	Water	
50 µL	Total reaction	

2. Mix gently by vortex and briefly centrifuge to collect all components to the bottom of the tube.
3. Add 50 µL of mineral oil to the top of each tube to prevent evaporation if not using a thermal cycler with a heated lid.
4. The amplification parameters should be optimized for individual primers, template, and thermal cycler.

#### Typical cycling parameters for 0.2–2 kb fragments:

Initial denaturation	94 °C for 2 min
<b>25-30 cycles:</b>	
Denaturation	94 °C for 30 sec
Annealing	55 °C to 68 °C for 30 sec
Extension	72 °C for 2 min
Final extension:	72 °C for 5 min
Hold	4 °C

5. The amplified DNA can be evaluated by loading 5-10 µL of the PCR reaction directly onto agarose gel. It is not necessary to add a separate loading buffer/tracking dye. Amplification products can be visualized by standard methodologies such as ethidium bromide staining.

**Note:** A minimum of 1.5 units of REDTaq Genomic DNA polymerase must be added per 50 µL reaction to ensure enough glycerol is present for direct gel loading. The red dye co-migrates with a 125 bp fragment in a 1% agarose gel.

REDTaq is a registered trademark of Sigma Biotechnology LP. TWEEN is a registered trademark of Croda International PLC.

#### **NOTICE TO PURCHASER: DISCLAIMER OF LICENSE**

No license is conveyed with the purchase of this product under any of US Patents Nos. 5,804,375, 5,994,056, 6,171,785, 6,214,979, 5,538,848, 5,723,591, 5,876,930, 6,030,787, and 6,258,569, and corresponding patents outside the United States, or any other patents or patent applications, relating to the 5' Nuclease and dsDNA-Binding Dye Processes. For further information contact the Director of Licensing, Applied Biosystems, 850 Lincoln Centre Drive, Foster City, California 94404, USA.

AH,RS,PHC 08/10-1