## **AMPIGENE<sup>®</sup> Taq DNA Polymerase**

Optimized standard PCR for enhanced speed, yield, and specificity

AMPIGENE® Taq DNA Polymerase uses the latest developments in polymerase technology and buffer chemistry to optimize PCR. The AMPIGENE® Taq DNA Polymerase and 5x AMPIGENE® reaction buffer are separate for flexibility and end-user customization. It can be used for routine PCR applications including genotyping, screening, and library construction.

## **Ordering Information**

Order Online »

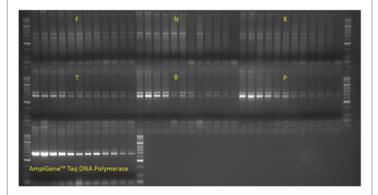
ENZ-PRT100-0500

500U

Manuals, SDS & CofA

**View Online** »

- For everyday PCR applications
- Consistent DNA polymerase
- Flexible format with separate buffer



Higher yield and sensitivity. 450bp fragment of the human myc gene was amplified with AmpiGene™ Taq and compared with Taq Polymerases from other suppliers. A serial dilution of human genomic DNA was incubated under standard PCR conditions. AmpiGene™ Taq delivers higher yield and sensitivity as compared with all six competing products.

## **Handling & Storage**

**Long Term Storage** -20°C

**Shipping** Blue Ice

Regulatory Status RUO - Research Use Only

**Product Details** 

**Application PCR** 

**Application Notes** For standard PCR applications

**Contents** 

AMPIGENE R Taq DNA polymerase (5u/ $\mu$ l) 5x AMPIGENE reaction buffer (15mM MgCl $_2$ , 5mM

dNTPs, enhancers and stabilizers)

Last modified: May 29, 2024



info-