

# AMPIGENE<sup>®</sup> Hot Start Taq DNA Polymerase

Enhanced standard Taq DNA Polymerase  
with increased sensitivity for a broader range  
of samples

AMPIGENE<sup>®</sup> Hot Start *Taq* DNA Polymerase uses the latest developments in polymerase technology and buffer chemistry to enhance PCR speed, yield and specificity. The enzyme uses advanced hot-start technology for superior sensitivity. The enzyme and buffer system allow for superior PCR performance on complex templates such as mammalian genomic DNA. Due to enhanced efficiency and specificity the enzyme is perfectly suited to difficult PCR.

AMPIGENE<sup>®</sup> Hot Start *Taq* DNA Polymerase is a robust enzyme for all your everyday PCR applications. AMPIGENE<sup>®</sup> Hot Start *Taq* DNA Polymerase can perform consistently well on a broad range of templates (including both GC and AT rich).

AMPIGENE<sup>®</sup> Hot Start *Taq* DNA Polymerase has an error rate of approximately 1 error per  $2.0 \times 10^5$  nucleotides incorporated. PCR products generated with AMPIGENE<sup>®</sup> Hot Start *Taq* DNA Polymerase are A-tailed and may be cloned into TA cloning vectors.

- For everyday PCR applications with increased sensitivity
- Advanced hot-start Taq DNA polymerase
- Flexible format with separate reaction buffer

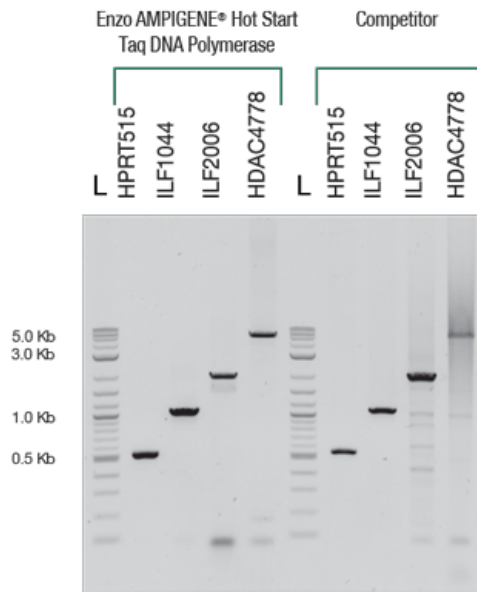
## Ordering Information

[Order Online »](#)

ENZ-NUC123-0200	200U
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)



**Performance of AMPIGENE® Hot Start Taq DNA Polymerase vs. Competitor.** AMPIGENE® Hot Start Taq DNA Polymerase shows strong signal and clear results. Human Genomic DNA (5ng) template was targeted with four separate primer sets in 20µl PCR volume. Annealing temperature was 60°C. 1 kb ladder (L).

## 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 that require higher sensitivity

Contents                              AMPIGENE<sup>®</sup> Hot Start *Taq* DNA Polymerase (5u/μL), 5x AMPIGENE<sup>®</sup> Reaction Buffer (8.75 mM MgCl<sub>2</sub>, 1 mM dNTPs, 250 mM KCl, 50 mM Tris, enhancers and stabilizers)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)