

AMPIGENE® Hot Start Taq DNA Polymerase

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

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

AMPIGENE® Hot Start *Taq* DNA Polymerase has an error rate of approximately 1 error per 2.0×10^5 nucleotides incorporated. PCR products generated with AMPIGENE® 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

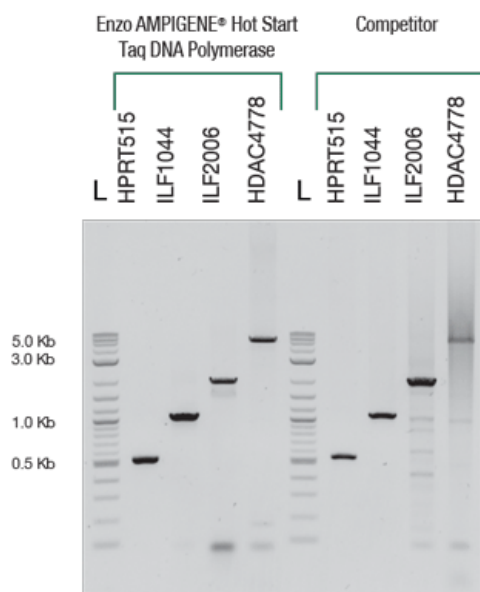
Ordering Information

[Order Online »](#)

| | |
|-----------------|------|
| ENZ-NUC123-0200 | 200U |
| ENZ-NUC123-0500 | 500U |

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[®] Hot Start *Taq* DNA Polymerase (5u/μL), 5x AMPIGENE[®] Reaction Buffer (8.75 mM MgCl₂, 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)