

AMPIGENE[®] Hot Start Taq Mix (2X)

Convenient 2x Hot Start Taq DNA Mix for increased sensitivity for a broader range of samples

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

AMPIGENE[®] Hot Start *Taq* Mix has an error rate of approximately 1 error per 2.0×10^5 nucleotides incorporated. PCR products generated with AMPIGENE[®] Hot Start *Taq* Mix are A-tailed and may be cloned into TA cloning vectors.

Ordering Information

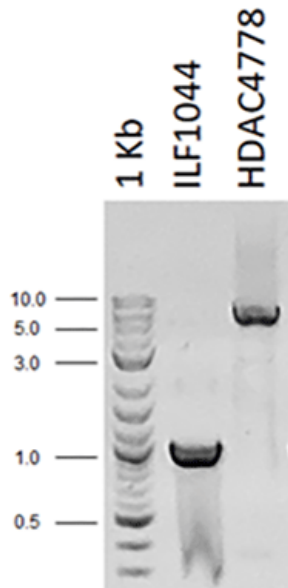
[Order Online »](#)

ENZ-NUC125-0200	200Reactions
-----------------	--------------

Manuals, SDS & CofA

[View Online »](#)

- For everyday PCR applications with increased sensitivity
- Advanced hot-start Taq DNA polymerase
- Convenient 2x Taq DNA polymerase mix



PCR amplification of human genomic DNA using AMPIGENE® Hot Start Taq Mix. PCR amplification was performed using two different primer sets. Amplification products were analyzed on 1.2% agarose gel. Human genomic DNA (25 ng) was used as template. Annealing temperature was 60°C. Left lane, 1 kb ladder.

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 2x AMPIGENE[®] Hot Start *Taq* Mix (AMPIGENE[®] Hot Start *Taq* DNA Polymerase, 3.5 mM MgCl₂, 0.4 mM dNTPs, enhancers and stabilizers)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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