

AMPIGENE[®] HS Taq

Mix

Increased sensitivity standard PCR mix for a broader range of samples, with enhanced speed, yield, and specificity

AMPIGENE[®] HS Taq Mix uses the latest developments in polymerase technology and buffer chemistry to optimize and increase PCR sensitivity. The AMPIGENE[®] HS Taq Mix is a robust mix using hot-start technology to increase sensitivity for PCR applications including genotyping, multiplex PCR, screening, library construction, colony PCR and PCR direct from blood and urine.

An optimized buffer system allows efficient amplification under fast and standard cycling conditions.

The AMPIGENE[®] HS Taq Mix is resistant to PCR inhibitors and suitable for direct PCR from unprocessed samples: bacterial culture, bacterial colonies, blood, and urine.

Ordering Information

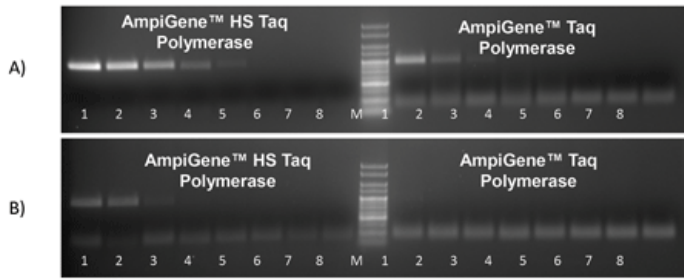
[Order Online »](#)

ENZ-NUC101-0040	40Reactions
ENZ-NUC101-0200	200Reactions

Manuals, SDS & CofA

[View Online »](#)

- For everyday PCR applications with increased sensitivity in an easy mix
- Advanced hot-start DNA polymerase
- Optimized buffer system for efficient amplification
- Resistant to PCR inhibitors and suitable for unprocessed samples



Human Genomic DNA 100ng/μl, then 1/3 serial dilution.

A) β 2MG (Beta-2 microglobulin) gene B) pgk gene.

AmpGene™ HS Taq DNA Polymerase does not produce primer-dimers and can amplify lower concentrations of DNA even from genes such as pgk that have high secondary structure levels.

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[®] HS Taq Mix (AMPIGENE[®] HS Taq Polymerase, 6mM MgCl₂, 2mM dNTPs, enhancer, 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