

AMPIGENE[®] HS Taq

DNA Polymerase

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

AMPIGENE[®] HS Taq DNA Polymerase uses the latest developments in polymerase technology and buffer chemistry to optimize and increase PCR sensitivity.

The AMPIGENE[®] HS Taq DNA Polymerase and 5x AMPIGENE[®] reaction buffer are separate for flexibility and end-user customization.

Hot-start technology is used to increase sensitivity for PCR applications including genotyping, multiplex PCR, screening, library construction, colony PCR and PCR direct from blood and urine.

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

Ordering Information

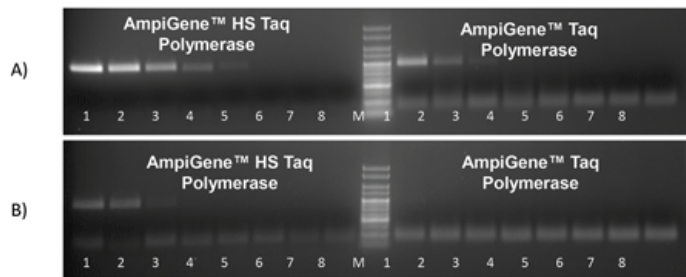
[Order Online »](#)

ENZ-PRT101-0500	500U
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- For everyday PCR applications with increased sensitivity
- Advanced hot-start DNA polymerase
- Flexible format with separate buffer
- Resistant to PCR inhibitors and suitable for unprocessed samples



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

A) β2MG (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 AMPIGENE[®] HS Taq DNA polymerase (5u/μl)
5x AMPIGENE[®] reaction buffer (15mM MgCl₂, 5mM dNTPs, 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)