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

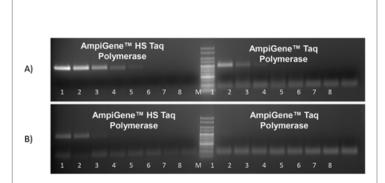
Order Online »

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) ß2MG (Beta-2 microglubulin) gene B) pgk gene.
AmpiGene™ 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

AMPIGENE® HS Taq DNA polymerase (5u/µl) **Contents**

5x AMPIGENE® reaction buffer (15mM MgCl₂, 5mM

dNTPs, enhancers and stabilizers)

Last modified: May 29, 2024



info-

eu@enzolifesciences.com