

AMPIGENE[®] Taq DNA Polymerase

Optimized standard PCR for enhanced speed, yield, and specificity

AMPIGENE[®] Taq DNA Polymerase uses the latest developments in polymerase technology and buffer chemistry to optimize PCR. The AMPIGENE[®] Taq DNA Polymerase and 5x AMPIGENE[®] reaction buffer are separate for flexibility and end-user customization. It can be used for routine PCR applications including genotyping, screening, and library construction.

Ordering Information

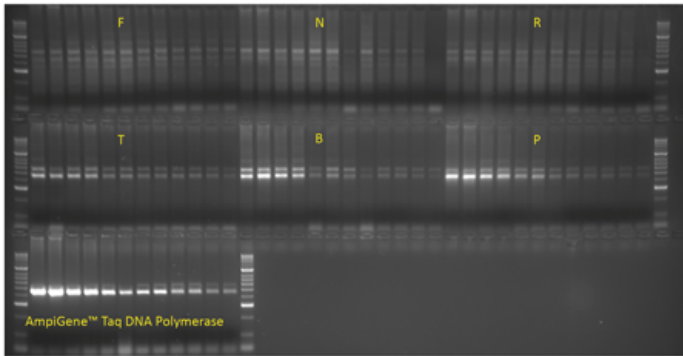
[Order Online »](#)

ENZ-PRT100-0500	500U
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

- For everyday PCR applications
- Consistent DNA polymerase
- Flexible format with separate buffer



Higher yield and sensitivity. 450bp fragment of the human myc gene was amplified with AmpGene™ Taq and compared with Taq Polymerases from other suppliers. A serial dilution of human genomic DNA was incubated under standard PCR conditions. AmpGene™ Taq delivers higher yield and sensitivity as compared with all six competing products.

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

Contents AMPIGENE[®] Taq DNA polymerase (5u/μl)
5x AMPIGENE[®] reaction buffer (15mM MgCl₂, 5mM dNTPs, enhancers and stabilizers)



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)