

AMPIGENE® Taq Mix

Optimized standard PCR mix for enhanced speed, yield, and specificity

AMPIGENE® Taq Mix uses the latest developments in polymerase technology and buffer chemistry to optimize PCR. The AMPIGENE® Taq Mix is a robust mix for routine PCR applications including genotyping, screening, and library construction. An optimized buffer system allows efficient amplification under fast and standard cycling conditions.

Citations: 3

[View Online »](#)

Ordering Information

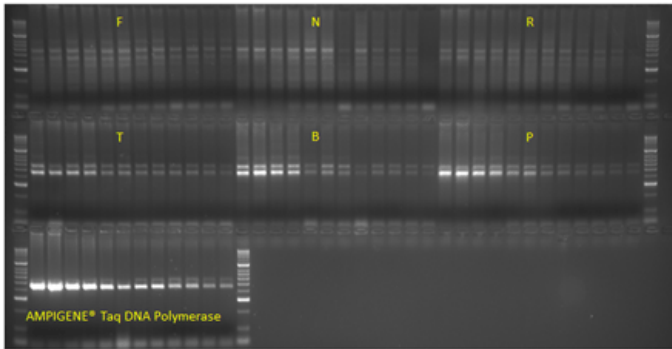
[Order Online »](#)

ENZ-NUC100-0040	40Reactions
ENZ-NUC100-0200	200Reactions
ENZ-NUC100-1000	1000Reactions

Manuals, SDS & CofA

[View Online »](#)

- For everyday PCR applications in an easy mix
- Consistent DNA polymerase
- Optimized buffer system for efficient amplification



Higher yield and sensitivity. 450bp fragment of the human myc gene was amplified with AMPIGENE® Taq and compared with Taq Polymerases from other suppliers. A serial dilution of human genomic DNA was incubated under standard PCR conditions. AMPIGENE® 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 2x AMPIGENE[®] Taq Mix (AMPIGENE[®] Taq Polymerase, 6mM MgCl₂, 2mM dNTPs, enhancer, 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)