

AMPIGENE[®] Hot Start High Fidelity Polymerase

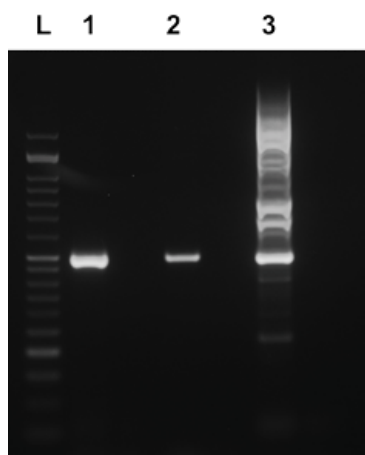
Optimized Hot Start High Fidelity Polymerase for greater sequence accuracy

AMPIGENE[®] Hot Start High Fidelity Polymerase was engineered to improve DNA binding resulting in amplification of difficult targets, greater yields, and shorter extension times.

Manuals, SDS & CofA

[View Online »](#)

- 100x greater fidelity than Taq DNA polymerase
- Increased processivity (10-30s/kb)
- Amplification of challenging GC and AT-rich templates to over 17.5kb
- Hot start technology for maximized sensitivity and specificity

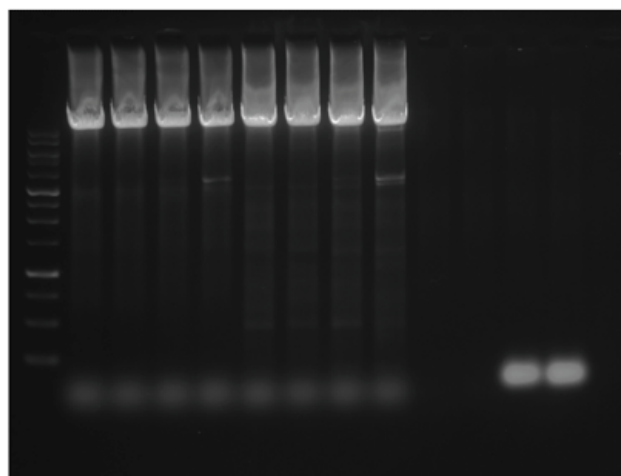


Lane L – Ladder
 Lane 1 – **AMPIGENE®**
Hot Start High Fidelity
Polymerase
 Lane 2 – Competitor 1
 Lane 3 – Competitor 2

AMPIGENE® Hot Start High Fidelity Polymerase Provides Enhanced Performance with GC-Rich Templates

AMPIGENE® Hot Start High Fidelity Polymerase exhibited superior performance with GC-rich template in comparison to the competitors' product. The template used was 500 pb long with 69% GC content.

AMPIGENE®
Hot Start High Fidelity
Polymerase **Competitor 1** **Competitor 2**



AMPIGENE® Hot Start High Fidelity Polymerase Provides Enhanced Performance with Difficult Templates

AMPIGENE® Hot Start High Fidelity Polymerase exhibited superior performance at different annealing temperatures (Ta1 = 68.5°C, Ta2 = 66.0°C, Ta3 = 63.0°C, Ta4 = 60.5°C) compared to the competitors' product. AMPIGENE® Hot Start High Fidelity Polymerase generated cleaner bands and no primer dimers. Human gDNA, 30 ng (8333 copies) and 13500 bp long was used as template.

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

Contents AMPIGENE[®] Hot Start High Fidelity Polymerase (2U/μl),
5x AMPIGENE[®] High Fidelity Rxn buffer, 10x AMPIGENE[®] Enhancer

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