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.

Ordering Information

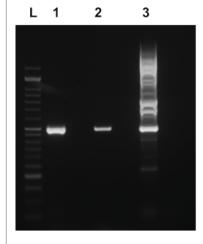
Order Online »

ENZ-NUC133-0100	100U
ENZ-NUC133-0500	500U

Manuals, SDS & CofA

View Online »

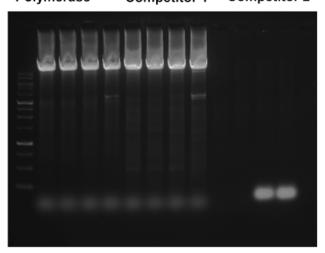
- 100x greater fidelity than Taq DNA polymerase
- Increased processivity (10-30s/kb)
- Amplification of challenging GC and AT-rich templates to over 17.5kb
- Hot start technology for maximized sensitivity and specificity



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

Lane 2 – Competitor 1 Lane 3 – Competitor 2

AMPIGENE® Hot Start High Fidelity Polymerase Competitor 1 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 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

AMPIGENE[®] Hot Start High Fidelity Polymerase (2U/μI), **Contents**

5x AMPIGENE® High Fidelity Rxn buffer, 10x AMPIGENE

® Enhancer

Last modified: May 29, 2024



info-

eu@enzolifesciences.com