AMPIGENE® High Fidelity Polymerase

Optimized High Fidelity Polymerase for greater sequence accuracy

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

Ordering Information

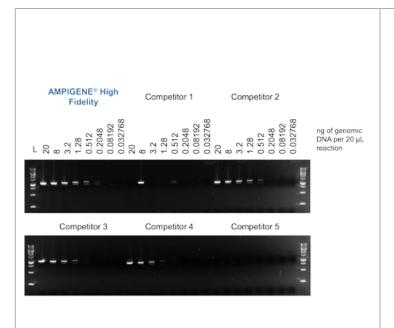
Order Online »

ENZ-NUC130-0100	100U
ENZ-NUC130-0500	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



AMPIGENE® High Fidelity Polymerase Provides Higher Yield and Sensitivity

AMPIGENE® High Fidelity Polymerase provides higher yield and sensitivity with low amount of target DNA compared to a competitor's product.

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

 $\mathsf{AMPIGENE}^{\circledR}$ High Fidelity Polymerase (2U/µI), 5x **Contents**

AMPIGENE® High Fidelity Rxn buffer, 10x AMPIGENE®

Enhancer

Last modified: May 29, 2024



info-

eu@enzolifesciences.com