

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

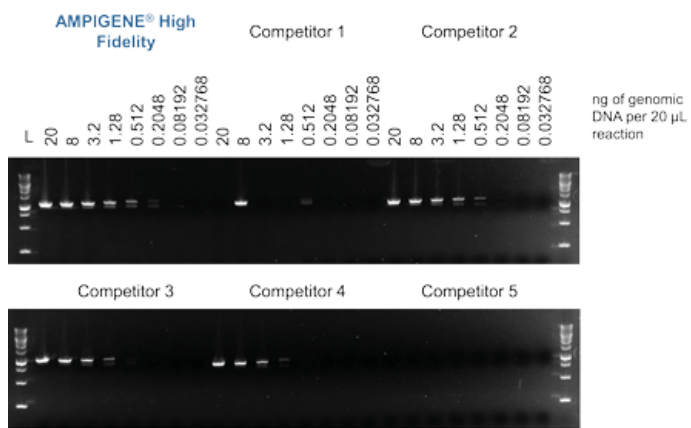
[Order Online »](#)

ENZ-NUC130-0500	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.

