

# AMPIGENE® High Fidelity Red Mix

Convenient 2x High Fidelity DNA Polymerase Mix with red dye for direct gel loading

AMPIGENE® High Fidelity DNA Polymerase Red Mix 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
- Convenient 2x Mix format and red dye added

