AMPIGENE[®] Hot StartHigh Fidelity Red Mix

Convenient 2x Hot Start High Fidelity Red Mix with red dye for direct gel loading

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

Ordering Information

Order Online »

ENZ-NUC135-0050

50Reactions

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

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

2x AMPIGENE® Hot Start High Fidelity Red Mix **Contents**

Last modified: May 29, 2024



info-

eu@enzolifesciences.com

uk@enzolifesciences.com