

AMPIGENE® Hot Start High Fidelity Mix

Convenient 2x Hot Start High Fidelity Mix

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

Ordering Information

[Order Online »](#)

ENZ-NUC134-0050	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

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

Contents 2x AMPIGENE[®] Hot Start High Fidelity Mix

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)