

AMPIGENE® Bst Mix

Recombinant *Bst* DNA Polymerase mix

AMPIGENE® *Bst* Polymerase (recombinant fragment of *Geobacillus stearothermophilus* DNA polymerase) is ideal for nucleic acid amplification methods taking place at a single incubation temperature. such as whole genome amplification, multiple displacement amplification, loop mediated isothermal amplification (LAMP), and isothermal amplification. A number of applications use these properties in point-of-care diagnostics, field-based detection, and fast diagnostic methods.

Ordering Information

[Order Online »](#)

ENZ-NUC140-0100	100Reactions
-----------------	--------------

Manuals, SDS & CofA

[View Online »](#)

- Rapid and consistent results
- 30 minute protocol
- Operates over broad temperature range (optimum 65°C)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application PCR

Contents AMPIGENE® *Bst* Mix, 20x Fluorescent Dye

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)