

# AMPIGENE® Bst Polymerase with Dye

Recombinant *Bst* DNA Polymerase with 5'-3'  
activity and dye added

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-NUC139-1600	1600U
ENZ-NUC139-8000	8000U

## Manuals, SDS & CofA

[View Online »](#)

- Rapid and consistent results
- Flexible format – available with or without dye (for real-time detection with any thermocycler)
- 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<sup>®</sup> *Bst* Polymerase 8U/μL, 10x AMPIGENE<sup>®</sup> Buffer A, 5x AMPIGENE<sup>®</sup> Buffer B, 20x Fluorescent Dye



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)