

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.

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[®] *Bst* Polymerase 8U/μL, 10x AMPIGENE[®] Buffer A, 5x AMPIGENE[®] Buffer B, 20x Fluorescent Dye

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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