

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.

Manuals, SDS & CofA

[View Online »](#)

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

