

AMPIGENE® Bst Polymerase

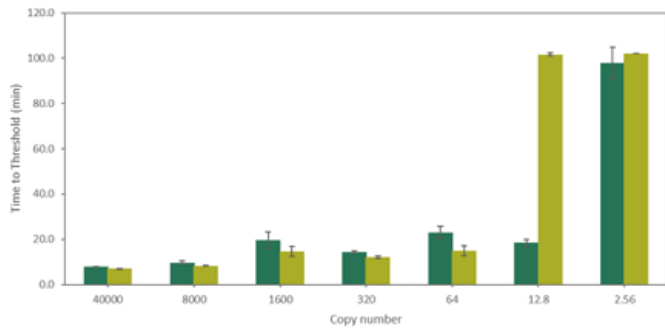
Recombinant *Bst* DNA Polymerase with 5'-3'
activity

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)



Isothermal amplification of SARS-CoV-2 N gene target using AMPIGENE® Bst Mix vs a competitor's product

AMPIGENE® Bst 1-Step Mix displayed earlier times to result and higher sensitivity compared to the competitor's product. Amplification of 13 copies of SARS-CoV-2 genome was obtained within 30 min.

