



## Saphir Bst2.0 Polymerase

Bst polymerase for isothermal DNA amplification  
Isothermal Amplification

Cat. No.	Amount
PCR-389S	2.000 Units
PCR-389L	10.000 Units

### For *in vitro* use only!

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Concentration:** 8 units/ $\mu$ l

### Description:

Saphir Bst2.0 Polymerase is a genetically improved Bst polymerase for rapid and specific amplification of DNA at constant temperature (60 to 65 °C). The enzyme shows high strand displacement activity and generates an amplification factor of up to  $10^9$  which is comparable to approx. 30 cycles in a PCR assay. This allows detection of a target gene within 10-30 minutes.

### Content:

#### Saphir Bst2.0 Polymerase

8 units/ $\mu$ l Bst DNA Polymerase in 10 mM Tris-HCl, 50 mM KCl, 50 % (v/v) Glycerol, pH 7.5 (25 °C) and stabilizers.

#### Saphir Bst2.0 Buffer

10 x conc. complete reaction buffer containing 200 mM Tris-HCl pH 8.8, KCl,  $(\text{NH}_4)_2\text{SO}_4$ , 20 mM  $\text{MgSO}_4$  and detergents.

#### $\text{MgSO}_4$ Stock Solution

25 mM  $\text{MgSO}_4$ .

### Detection

Although some methods have been developed to visualize DNA amplification by basic equipment or even the naked eye (increase of turbidity, color change of added dyes, hybridization to gold-bound ss-DNA) in general real-time detection of the DNA amplification by a fluorescent DNA-intercalator dye is recommended. Addition of EvaGreen Fluorescent DNA Stain (#PCR-379) to the assay allows a sensitive measurement of the increasing amount of DNA without influence on the reaction.

### Assay design

Isothermal amplification is an extremely sensitive detection method and care should be taken to avoid contamination of set-up areas and equipment with DNA of previous reactions. A common problem is amplification in no-template controls due to

1. carry-over contamination or
2. amplification of unspecifically annealed primers or primer dimer formations.

As sensitivity and non-template amplification of in-silico designed primers may vary, the evaluation of 2-4 real primer sets before choosing a final set is recommended.



## Saphir Bst2.0 Polymerase

Bst polymerase for isothermal DNA amplification  
Isothermal Amplification

### Assay set-up

Depending on the detection method and machine a reaction volume of 20–50 µl is recommended for most applications. Pipet with sterile filter tips and perform the set-up in an area separate from DNA preparation or analysis. No-template controls should be included in all amplifications.

First, prepare a 10x conc. primer pre-mix. Second, set-up the isothermal amplification assay:

component	stock conc.	final conc.	20 µl	50 µl
Saphir Bst2.0 Buffer	10x	1x	2 µl	5 µl
MgSO <sub>4</sub> Stock Solution	25 mM	6–8 mM	3.2–4.8 µl	8.0–12.0 µl
dNTP Mix	10 mM	1.4 mM	2.8 µl	7 µl
Primer Mix	10x	1x	2 µl	5 µl
Saphir Bst2.0 Polymerase	8 units/µl	0.32 units/µl	0.8 µl	2 µl
EvaGreen DNA Stain	100 µM	1.3 µM	0.26 µl	0.65 µl
Template DNA		<500 ng/assay	x µl	x µl
PCR-grade Water			fill up to 20 µl	fill up to 50 µl

- Use a specific detection instrument for isothermal amplification or a real-time PCR cycler to run the assays
- Set the instrument to a constant incubation temperature between 60 to 65 °C (depending on the primer annealing temperature)
- Measure the fluorescence intensity at an interval of 1 min for up to 30 min.

### Optimization of MgSO<sub>4</sub> concentration:

A final Mg<sup>2+</sup> concentration of 6.0 mM is optimal for most primer-template combinations. However, if an individual Mg<sup>2+</sup> optimization is essential add 25 mM MgSO<sub>4</sub> stock solution as shown in the table below.

final MgSO <sub>4</sub> conc.	20 µl final assay volume	50 µl final assay volume
6 mM	3.2 µl	8.0 µl
7 mM	4.0 µl	10.0 µl
8 mM	4.8 µl	12.0 µl

### Trouble shooting

If amplification in no-template controls occurs the following points should be reviewed.

#### Cross contamination from environments

- Clean equipment and areas with “DNA Away” solution
- Replace reagent stocks and pre-mixes with new components
- Stop reactions at an earlier point of time before non-template amplification occur

#### Carry-over contamination from previous reaction products

- Avoid opening reaction vessels after amplification
- Use separate preparation area and equipment if post-reaction processing is necessary

#### Non-template amplification from primers

- Increase incubation temperature stepwise by 1–2 °C
- Design a new set of primers for the target sequence

### Related Products:

EvaGreen DNA Stain, #PCR-379  
dNTP Mix / 10 mM, #NU-1006  
dNTP Mix / 25 mM, #NU-1023