

# Saphir Bst2.0 Turbo GreenMaster, S pack

Master mix for isothermal DNA amplification  
with EvaGreen

Saphir Bst2.0 Turbo GreenMaster is a complete 2x conc. master mix for isothermal amplification of DNA. The mix is based on a genetically enhanced Bst polymerase with an additionally fused DNA-binding domain. The mix is the ideal choice for ultra-fast and robust 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 109 which is comparable to approx. 30 cycles in a PCR assay. The polymerase is 2-3x faster compared to Saphir Bst2.0 Polymerase (#PCR-389/#PCR-388/#PCR-387) and allows detection of a target gene within 5-10 minutes.

## Ordering Information

[Order Online »](#)

JBS-PCR-393S	2x1.25ml
--------------	----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance liquid

Concentration 2x conc.

Emission Maximum 530 nm (EvaGreen bound to DNA)

Excitation Maximum 500 nm (EvaGreen bound to DNA)

**Technical Info / Product Notes** For the Original Manufacturer's data sheet please [click here](#).



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)