

qPCR SybrMaster, S pack

Master mix for real-time qPCR with
SYBR®Green fluorescent DNA stain

qPCR SybrMaster is designed for quantitative real-time analysis of DNA samples. The mix contains all reagents required for qPCR (except template and primers) in a premixed 2x concentrated ready-to-use solution. It is recommended for routine PCR applications, high throughput PCR or genotyping and provides an improved specificity and sensitivity when amplifying low-copy-number targets or working with complex backgrounds. The mix is based on an optimized hot-start polymerase. Its activity is blocked by antibody at ambient temperature and switched on automatically at the onset of the initial denaturation. The thermal activation prevents the extension of nonspecifically annealed primers and primer-dimer formation at low temperatures during PCR setup. The fluorescent DNA stain SYBR®Green intercalates into the amplification product during the PCR process and allows the direct quantification of target DNA without the need to synthesize sequence-specific labeled probes (i.g. TaqMan® Probes). The mix can also be used in combination with ROX reference dye (#PCR-351) in PCR instruments that are compatible with the evaluation of the ROX signal. SYBR®Green Fluorescent DNA Stain SYBR®Green Fluorescent DNA Stain is a superior DNA intercalator dye specially developed for DNA analysis applications including real-time PCR (qPCR). Upon binding to DNA, the non-fluorescent dye becomes highly fluorescent while showing no detectable inhibition to the PCR process. The dye is extremely stable, providing convenience during routine handling. SYBR®Green is in contrast to EvaGreen® not recommended for high-resolution melting curve analysis (HRM). To perform the SYBR®Green-based assay simply select the optical setting for SYBR®Green on the detection instrument.

Ordering Information

[Order Online »](#)

JBS-PCR-372S	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

Emission Maximum 521 nm (bound to DNA)

Excitation Maximum 494 nm (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

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