

SCRIPT RT-PCR

Master (2x), L pack

One-Step RT-PCR Master Mix (2x conc.) for highly sensitive and specific amplification

SCRIPT RT-PCR Master is designed for performing highly sensitive and specific RT-PCR convenient in single tubes. The enzyme mix is based on a genetically engineered reverse transcriptase with enhanced thermal stability providing increased specificity, high cDNA yield and improved efficiency for highly structured and long cDNA fragments. The 2 x conc. mix contains all reagents required for RT-PCR (except template and primer) in one tube to ensure fast and easy preparation with a minimum of pipetting steps. Premium quality enzymes in combination with ultrapure dNTPs and an advanced buffer system ensure superior amplification results. RT-PCR is used to amplify double-stranded DNA from single-stranded RNA templates. In the RT step the reverse transcriptase synthesizes single-stranded DNA molecules (cDNA) complementary to the RNA template. In the first cycle of the PCR step hot start DNA polymerase synthesizes DNA molecules complementary to the cDNA, thus generating a double-stranded DNA template. During subsequent rounds of cycling the DNA polymerase exponentially amplifies this double-stranded DNA template. In one-step RT-PCR all components of RT and PCR are mixed in one tube so that the complete reaction can be performed without opening the tube. This offers tremendous convenience when applied in routine testing and minimizes the risk of contaminations. SCRIPT RT-PCR Master contains RNase inhibitor that is essential when working with low amounts of starting RNA.

Sensitivity: Targets can generally be detected from

Ordering Information

[Order Online »](#)

JBS-PCR-525L	10x1.25ml
--------------	-----------

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

Quantity 2500 reactions x 20 µl

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