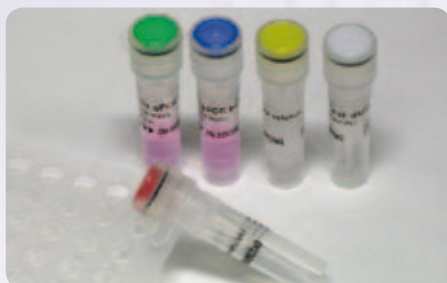


RT-PCR Kits and Mastermixes



All kits are based on SCRIPT Reverse Transcriptase, a genetically engineered M-MLV RTase with enhanced thermal stability. The enzyme shows increased specificity and improved efficiency for highly structured and long cDNA fragments. Corresponding premium quality components like ultrapure dNTPs and optimized reaction buffer ensure superior amplification results.

In particular, our **SCRIPT RT-PCR Kits and Mixes**

- guarantee highest cDNA yields
- are recommended for highly structured and/or long cDNA fragments
- allow fast and easy preparation set-ups to save time at the bench
- come at a price one can't ignore

Product	Cat.-No.	Size	Price
SCRIPT RT-PCR One-Step Master	PCR-508XS	20 reactions	90,- Euro
	PCR-508S	100 reactions	360,- Euro
	PCR-508L	500 reactions	1440,- Euro
SCRIPT RT-PCR One-Step Kit	PCR-507S	25 reactions	118,- Euro
	PCR-507L	200 reactions	670,- Euro
SCRIPT RT-PCR Two-Step Kit	PCR-506S	25 reactions	118,- Euro
	PCR-506L	200 reactions	670,- Euro
SCRIPT Reverse Transcriptase	PCR-505S	10,000 units	140,- Euro
	PCR-505L	50,000 units	560,- Euro
RNase Inhibitor	PCR-504S	2,000 units	45,- Euro
	PCR-504L	10,000 units	180,- Euro

Please contact us for inquiries and further information or request your customized quotation.

Find out more details at our website
www.jenabioscience.com



Jena Bioscience GmbH
 Loebstedter Str. 80
 07749 Jena
 Germany

Phone +49 (0)3641-62 85 000
 Fax +49 (0)3641-62 85 100
www.jenabioscience.com
info@jenabioscience.com

Jena Bioscience
www.jenabioscience.com