

RNA / DNA Purification and Cleanup Kits



- Excellent yields and recovery rates
- Highly pure RNA/DNA preparations for all down-stream applications
- Fast and easy preparation to save time at the bench
- Prices one can't ignore

Our RNA/DNA preparation section covers spin column based as well as solution based genomic DNA preparation kits for isolation of **RNA and DNA from blood, bacteria, plant cells, animal tissues and yeast.**

All kits are optimized for obtaining maximum yield of high quality DNA suitable for **automated sequencing, cloning, genotyping, real-time PCR, in vitro transcription and many more.**

Product	Cat.-No.	Size	Price
RNA Purification			
Total RNA Purification Kit	PP-210XS	10 preparations	24,- €
	PP-210S	50 preparations	90,- €
	PP-210L	250 preparations	360,- €
Genomic DNA Purification, spin column based			
Blood-Plant-Animal DNA Preparation Kit	PP-213S	50 preparations	70,- €
	PP-213L	250 preparations	280,- €
Bacteria DNA Preparation Kit	PP-214S	50 preparations	80,- €
	PP-214L	250 preparations	320,- €
Yeast DNA Preparation Kit	PP-215S	50 preparations	80,- €
	PP-215L	250 preparations	320,- €

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

For more information and for ordering, please visit the RNA / DNA Preparation and Cleanup section of our website at **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

RNA / DNA Purification and Cleanup Kits

Product	Cat.-No.	Size	Price
Genomic DNA Purification, solution based			
Blood DNA Preparation Kit	PP-205S	100 preparations	54,- €
	PP-205L	500 preparations	216,- €
Bacteria DNA Preparation Kit	PP-206S	100 preparations	61,- €
	PP-206L	500 preparations	245,- €
Plant DNA Preparation Kit	PP-207S	100 preparations	54,- €
	PP-207L	500 preparations	216,- €
Animal and Fungi DNA Preparation Kit	PP-208S	100 preparations	54,- €
	PP-208L	500 preparations	216,- €
Yeast DNA Preparation Kit	PP-209S	100 preparations	61,- €
	PP-209L	500 preparations	245,- €
Plasmid DNA Purification			
Fast-n-Easy Plasmid Mini-Prep Kit	PP-204XS	10 preparations	12,- €
	PP-204S	50 preparations	45,- €
	PP-204L	250 preparations	180,- €
DNA Cleanup			
PCR Purification Kit	PP-201XS	10 preparations	12,- €
	PP-201S	50 preparations	45,- €
	PP-201L	250 preparations	180,- €
Agarose Gel Extraction Kit	PP-202XS	10 preparations	12,- €
	PP-202S	50 preparations	45,- €
	PP-202L	250 preparations	180,- €

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

For more information and for ordering, please visit the RNA / DNA Preparation and Cleanup section of our website at 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