nucleic acid purification QUANTUM PREP® PLASMID MINIPREP KIT

FEATURES OF THE QUANTUM PREP PLASMID MINIPREP KIT:

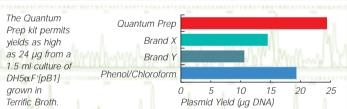
- YIELDS HIGHLY PURIFIED DNA, IDEAL FOR AUTOMATED FLUORESCENT SEQUENCING
- QC-SPECIFIED
 DNA SEQUENCING
 ACCURACY OF AT
 LEAST 98% OVER
 600 BASES
- FAST, FROM
 CELL CULTURE
 TO PURIFIED DNA
 IN LESS THAN
 15 MINUTES
- HIGH YIELD, WITH OVER
 20 µg PLASMID DNA FROM A
 1.5 ml OVERNIGHT CULTURE
- EASY SPIN
 COLUMN FORMAT
 ELIMINATES
 PHENOL
 EXTRACTIONS
 AND ETHANOL
 PRECIPITATION

A Patented Matrix for High Quality Sequencing Results

The advent of automated DNA sequencing and advances in genetic technology have brought a requirement for increased yields of ever more highly purified preparations of plasmid DNA. The Quantum Prep Plasmid Miniprep kit was developed to meet the stringent purity requirements of the DNA sequencer, while providing optimal yields of plasmid for gene transfer and cloning protocols. The key to the Quantum Prep matrix's consistently high purity yields lies in the unique Quantum Prep matrix: the highly-porous siliceous exoskeleton of the remains of prehistoric diatoms. Proprietary treatments produce a matrix with ideal DNA binding and purification characteristics.

Higher Yields

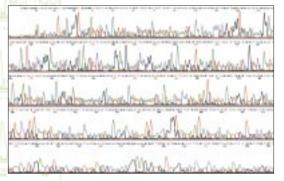
The unique 3-D structure of the diatom exoskeleton provides the greatest number of DNA binding sites available, so you are able to use enriched growth media such as Terrific Broth to obtain the highest yields on the market. Anion exchange and embedded silica resins don't come close to this kind of performance. The Quantum Prep plasmid miniprep kit is recommended for both high and low copy number plasmids.



u are to and ids.

Longer, More Accurate Sequencing Reads

To guarantee the highest quality DNA, each lot of Quantum Prep Plasmid Miniprep kits must pass stringent QC specifications (see table at right). This ensures that your DNA is suitable for the most demanding applications, including fluorescent sequencing, transfection and transformation protocols, *in vitro* transcription reactions, PCR* and subcloning experiments.



OC analysis of Quantum Prep plasmid miniprep kit-purified DNA. Fluorescent DNA sequence analysis using the AB™ PRISM™ Dye Terminator kit and analyzed on an ABI 377A DNA sequencer.



Lower Cost

The Quantum Prep Plasmid Miniprep kit has the lowest cost per microgram of purified DNA of any kit available. The patented matrix is tested to yield at least 20 µg of high copy plasmid from just 1.5 ml of culture, at a price you can afford. Each kit includes enough material to process 100 1.5 ml overnight cultures.

15-Minute Procedure

The Quantum Prep Plasmid Miniprep procedure takes less than 15 minutes from cell culture to purified plasmid. DNA is recovered in H₂O or TE for immediate use in sequencing, transfection, PCR, subcloning, and other protocols requiring highly purified DNA. Ready-to-use reagents do not require mixing or dilution; phenol extraction, CsCl₂, and ethanol precipitation are eliminated.

Specifications

PURITY >98% accuracy over

600 bases on ABI fluorescent sequencer

TIME 15 minutes, including

cell lysis

PLASMID YIELD $>20 \mu g$ (high copy

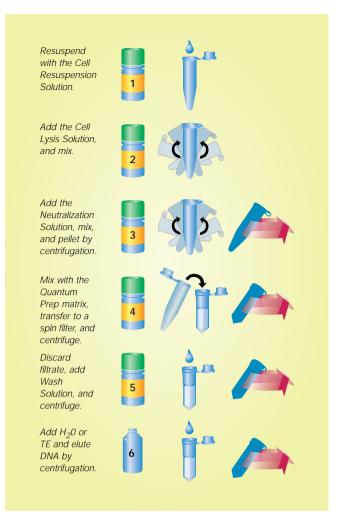
plasmids)

GROWTH MEDIA LB & TB overnight cultures **CULTURE VOLUME** 1-5 ml high-copy plasmids;

1-10 ml low-copy plasmids

METHOD Spin column/

microcentrifuge



^{*} The Polymerase Chain Reaction (PCR) process is covered by patents owned by Hoffman-LaRoche. Use of the PCR process requires a license.

ABI and PRISM are trademarks of The Perkin-Elmer Corp.

Ordering Information Catalog # Description

outuiog "	Bosonphon
732-6100	Quantum Prep Plasmid Miniprep Kit, 100 preps, includes 20 ml cell resuspension solution, 25 ml cell lysis solution,
	25 ml neutralization solution, 20 ml Quantum Prep matrix, 63 ml wash buffer, 100 spin filters, instructions
732-6120	Quantum Prep Plasmid Midiprep Kit, 20 preps
732-6130	Quantum Prep Plasmid Maxiprep Kit, 10 preps

REPLACEMENT ACCESSORIES

/32-0110	Quantum Prep Matrix, 20 mi
732-6115	Quantum Prep Neutralization Solution, 25 ml
732-6122	Quantum Prep Neutralization Solution, 125 ml
732-6027	Quantum Prep Mini Spin Filters, 100
732-6116	Quantum Prep Midi Spin Columns, 20
732-6137	Quantum Prep Midi Spin Baskets, 10



Bio-Rad Laboratories

Life Science Group Website www.bio-rad.com U.S. (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Canada (905) 712-2771 China 86-10-62051850 Denmark 45 39 17 99 47 Finland 358 (0)9 804 2200 France (01) 43 90 46 90 Germany 089 318 84-0 Hong Kong 852-2789-3300 India (91-11) 461-0103 Israel 03 951 4127 Italy 02-21609.1 Japan 03-5811-6270 Korea 82-2-3473-4460 The Netherlands 31 318-540666 New Zealand 64-9-4152280 Singapore 65-2729877 Spain (91) 661 70 85 Sweden 46 (0)8 627 50 00 Switzerland 01-809 55 55 United Kingdom 0800-181134

Bulletin 2277 US/EG Rev A 98-024 0498 Sig 120197