



EMnetik system

The EMnetik system can perform PCR and other enzymatic reaction cleanups and isolate plasmids from bacteria.

Users can:

- Perform 1-24 samples at a time
- Perform PCR cleanup in less than 16 minutes
- Expect at least 80% recovery with a DNA purity of (260/280) > 1.7 and (260/230) > 1.8
- Choose an elution volume between 20 µL and 50 µL



● **PHYSICAL DIMENSIONS**

17.15 cm W x 31.50 cm D x 17.02 cm H
6.75 in W x 12.40 in D x 6.70 in H

● **WEIGHT**

3.24 kg
7.15 lbs

● **OTHER**

KIT SIZE

EMnetik PCR Cleanup Kit: 500 preps for 50 µL reactions
EMnetik Plasmid Purification Kit: 96 preps

KIT SHELF LIFE

9 months from manufacturing

NOISE

< 52 dBA while in operation

● **ELECTRICAL REQUIREMENTS**

EXTERNAL POWER SUPPLY RATINGS

Input: AC voltage range from 100 to 240 VAC at 50/60 Hz, 10 A
Output: 24 VDC, 9.2 A

COMMUNICATIONS TO INTERNET (FOR SOFTWARE UPGRADES)

Ethernet



Not intended or validated for use in the diagnosis of disease or other conditions

© 2021 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com
21.08.2338.GEN