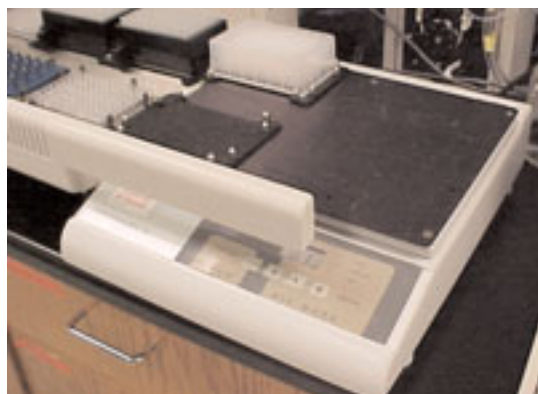


# DPC MicroMix<sup>®</sup> 5 Shaker Integration: Biomek<sup>®</sup> 2000

## DPC MicroMix<sup>®</sup> 5 Shaker Integration to the Biomek<sup>®</sup> 2000: Basic Installation Instructions

1. Unpack the shaker and integration kit.
2. Attach the black metal plate to the top of the shaker using the provided screws and spacers.
3. Attach the two black shaker labware holders on the left side of the shaker plate.
4. Attach the two shaker platform feet to the shaker platform.
5. Raise the deck of the Biomek<sup>®</sup> 2000 using the four screw-threaded perspex feet. Ensure that deck of the instrument is level.
6. Remove the right-side module of the Biomek<sup>®</sup> 2000.
7. Attach the shaker platform to the Biomek<sup>®</sup> 2000 base module.
8. Place the shaker onto the shaker platform.
9. Connect a free serial port on the back of the computer to the serial connection port on the back of the shaker.
10. Connect the power cord from the shaker to a power outlet.
11. Install the MicroMix<sup>®</sup> 5 software using all the default settings.



DPC MicroMix<sup>®</sup> 5 shaker integration to the Biomek<sup>®</sup> 2000 workstation.

© 2003 Promega Corporation. All Rights Reserved.

Biomek is a registered trademark of Beckman Coulter, Inc. MicroMix is a registered trademark of Clintec Nutrition, Co.



### Promega Corporation

2800 Woods Hollow Road  
Madison, WI 53711-5399 USA  
Telephone 608-274-4330  
Fax 608-277-2516  
Internet [www.promega.com](http://www.promega.com)