

Importing Biomek® 2000 Programs



Importing a Method

1. The Beckman Biomek® 2000 program is zipped into a self-extracting program. Unzip the program by double-clicking on the program icon. This will bring up a “browse” window to define where the files will be downloaded. We suggest making a new file folder for each Biomek® 2000 method. This prevents confusion later when another method is imported. Once unzipped, about 42 files will appear. These comprise the Biomek® 2000 method. Make a note of where the files are downloaded (e.g., C:/Program Files/....., etc.).
2. Open the Bioworks™ software “B2kEdit”. Check to make sure that you are in the desired Lab Book. The current Lab Book is shown in the title bar of the Bioworks Method Edit window (e.g., <Lab Book> Bioworks Method Edit). If you are not in the desired Lab Book, close “B2KEdit” and open “Lab Book Manager” to select the desired Lab Book. Once you have done this, close “Lab Book Manager” and open “B2KEdit” again.
3. In “B2KEdit”, select “Import” from the “Method” menu (at the top of the page). When importing a new method, you cannot have any other methods open).
4. At the bottom of the “Import” window, click “Browse”.
5. Find and select the folder containing the files you downloaded in Step 1. Click “OK”.
6. The method name should appear in the Method window. Click on the method name, highlighting the method, and then click “OK”. The file will be imported to the Bioworks™ lab book. This step may take a few moments. You may be prompted with messages such as “Labware holder exists. Overwrite?” You can select either “yes” or “no” because Promega Biomek™ 2000 methods do not change the standard labware definitions. After the method is imported, you will need to close the Method window.
8. To open the new protocol, click on “method”, “open”. The newly imported method should appear in the list of available protocols.

Note: Make sure that you have first unzipped the method file (Step 1). The Bioworks Software will not recognize the method if it is zipped.

© 2003 Promega Corporation. All Rights Reserved.

Biomek is a registered trademark of Beckman Instruments, Inc.



Promega Corporation

2800 Woods Hollow Road
Madison, WI 53711-5399 USA
Telephone 608-274-4330
Fax 608-277-2516
Internet www.promega.com