

Importing Beckman Biomek® 2000 Programs

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, separate, file folder for each Biomek® 2000 method to unzip the files to. This prevents confusion later when another method is imported. Once unzipped, about 42 files will appear that comprise the Biomek® 2000 method. Make a note of where these files are downloaded. (For example: C:/Program Files/.....)
2. Create a new Bioworks™ notebook named PROMEGA.
 - a. Open BioWorks™ software and double-click on the Lab Book Manager icon.
 - b. Click on NEWBOOK to highlight.
 - c. Then click on the “Copy from list” button. This brings up a new window: “Copy Lab Book.”
 - d. Type “PROMEGA” in the new name box.
 - e. In the new path box highlight “NEWBOOK” between the forward slashes and replace with “PROMEGA”. The new path should look like “C:\BIOWORKS\PROMEGA\
 - f. Click “OK”
 - g. The new PROMEGA labbook should appear in the list of available labbooks.
 - h. Click on the PROMEGA labbook to highlight. Then click the close button. This will select the PROMEGA labbook as the labbook where you will import the new method.
3. Open “B2KEdit”.
4. Check to verify in the upper right hand corner, in the blue bar of the screen, that you are in the PROMEGA labbook. It should look like:
<PROMEGA> BioWorks Method Editor.
5. Click on “Method” at the top of the page.
6. Click on “Import...” (You can not have any methods open)
7. At the bottom of the “Import” window, click on “Browse”.
8. Find and select the folder that you unzipped the files into in step one above. Click “OK”.

9. The method name should appear in the “Method” window. Click on this method name, to highlight the method, and then click on “OK”. The file will be imported to the BioWorks™ lab book, and may take a few moments. You will probably be prompted with windows that say things like “Labware holder exists. Overwrite?” Promega Biomek™ 2000 methods do not change the standard labware definitions, so you can select either “yes” or “no”—it won’t matter. After the method is imported, you will need to close the “method” window.

10. To open the new protocol, click on “method”, “open”. The newly imported method should appear in the list of available protocols.

If you have any questions, please contact Promega Technical Services at techserv@promega.com or toll free in U.S. at 1-800-356-9526.

For Technical Service telephone numbers in other countries visit our website to view information for Worldwide contacts: <http://www.promega.com/worldwide/>