

USER GUIDE

applied
biosystems®
by *life* technologies™

SOLiD® EZ Bead™ System Software Installation User Guide

Software Installation v1.9

Publication Part Number 4472494 Rev. A
Revision Date October 2011



life
technologies™

For General Laboratory Use.

For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

Information in this document is subject to change without notice.

LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE. IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

TRADEMARKS

The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

© 2011 Life Technologies Corporation. All rights reserved.

Publication Part Number 4472494 Rev. A

Revision Date October 2011

Contents

Install Software Version 1.9	5
Purpose	5
Required materials	5
Obtain the software applications updates	5
Install the software applications updates	6
Reformat USB flash drive	6
SOLiD® EZ Bead™ Emulsifier install	6
SOLiD® EZ Bead™ Amplifier install	8
SOLiD® EZ Bead™ Enricher install	9
Documentation and Support	13
Related documentation	13
Obtaining support	14

Install Software Version 1.9

Purpose

The SOLiD® EZ Bead™ System Software Installation User Guide provides detailed information about downloading and installing software version 1.9 for the Emulsifier, Amplifier and Enricher systems.

Required materials

In order to update your SOLiD® EZ Bead™ System with the latest software version 1.9, you must have the following instrumentation and obtain their corresponding software application updates:

- SOLiD® EZ Bead™ Emulsifier
- SOLiD® EZ Bead™ Amplifier
- SOLiD® EZ Bead™ Enricher
- USB flash drive (must be reformatted)

Obtain the software applications updates

To obtain the software applications updates:

1. Send an email to EZBeadAssist@lifetech.com.
2. Respond to the instructions in the email reply.
3. Receive files.

You receive a single .ZIP file folder that contains the following files:

Name	Components
SOLiD® EZ Bead™ Emulsifier	3 files: <ul style="list-style-type: none">• EzBake.hex• firmware.bin• recipes.csv
SOLiD® EZ Bead™ Amplifier	1 file: <ul style="list-style-type: none">• Amplifier_v1_9.exe
SOLiD® EZ Bead™ Enricher	1 file: <ul style="list-style-type: none">• Enricher_v1_9.exe

IMPORTANT! Ensure that any USB sticks used for installation/updates are scanned for viruses using a computer that has the most up-to-date virus definitions.

4. Unzip the files.

Install the software applications updates

To install the software applications updates and to enable the instruments to prepare or sequence the nanobeads, follow the procedures indicated in each section below.

Reformat USB flash drive

Before performing the software installation you must reformat your USB flash drive.

1. Insert the USB flash drive into a USB port on your computer.
2. Select **My Computer** and in the list of **Devices with Removable Storage**, select and right-click on the USB flash drive.
3. Choose **Format**, then **Start**.
The USB flash drive is reformatted.

SOLiD® EZ Bead™ Emulsifier install

To install the software application updates for the SOLiD® EZ Bead™ Emulsifier follow these steps:

1. Save the following files to the reformatted USB stick:
 - EzBake.hex
 - firmware.bin
 - recipes.csv
2. Scan the USB stick for viruses, using a computer that has the most up-to-date virus definitions.
3. Update the Graphical User Interface (GUI) application by following these steps:
 - a. Power off the instrument.
 - b. Insert the USB stick into the USB port on the back of the instrument, power on the instrument, and wait for the GUI to display the following 3 buttons.

Note: This process may take a few minutes to complete.

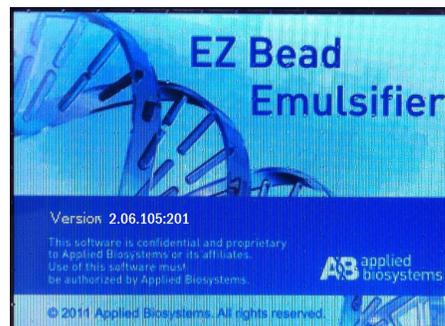
Note: If a USB error message appears, disconnect and reconnect the USB stick.
 - c. After data acquisition is complete, press the select button  and select **App**.
 - d. Press the play button  to initiate the application update.

- e. When prompted, confirm by pressing the multifunction button  and wait for the instrument to reboot and begin the update process.

Note: This process takes several minutes to complete. Do not remove the USB stick.

- f. When prompted on the GUI, remove the USB stick to allow the instrument to restart.
- g. Ensure that the GUI application is version 2.06.xxx.xxx.

Note: To verify that the correct version has been installed, power on the instrument and ensure that the splash screen displays the new version on the screen.



4. Update the IOC (Input/Output Controller) application by following these steps:

- a. Insert the USB stick into the USB port on the back of the instrument, power on the instrument, and wait for the GUI to display the following 3 buttons.



Note: This process may take a few minutes to complete.

Note: If a USB error message appears, disconnect and reconnect the USB stick, power off the instrument, reconnect the USB stick, and power on the instrument. If the USB error message appears again, reformat the USB stick, obtain the upgrade files, and proceed to update the IOC application.

- b. Using the GUI, press the select button  and select IOC.
- c. Press the play button  to initiate the application update.
- d. When prompted, confirm by pressing the multifunction button  and wait for the instrument to reboot and begin the update process.

Note: This process takes several minutes to complete. Do not remove the USB stick.

- e. When prompted on the GUI, remove the USB stick to allow the instrument to restart.
- f. Ensure that the IOC application is version 2.06.

5. Update the Rec (Recipes) application by following these steps:

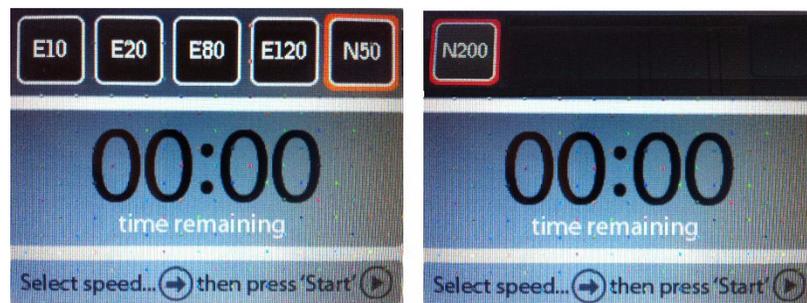
- a. Power off the instrument.
- b. Insert the USB stick into the USB port on the back of the instrument, power on the instrument, and wait for the GUI to display the following 3 buttons.



Note: This process may take a few minutes to complete.

Note: If a USB error message appears, disconnect and reconnect the USB stick, power off the instrument, reconnect the USB stick, and power on the instrument. If the USB error message appears again, reformat the USB stick, obtain the upgrade files, and proceed to update the IOC application.

- c. Using the GUI, press the select button  and select **REC**.
 - d. Press the play button  to initiate the application update.
 - e. When prompted, confirm by pressing the multifunction button  and wait for the instrument to reboot and begin the update process.
- Note:** This process takes several minutes to complete. Do not remove the USB stick.
- f. When prompted on the GUI, remove the USB stick to allow the instrument to restart.
 - g. Power off the instrument.
 - h. Power on the instrument.
 - i. Ensure that the GUI shows the correct scales. The correct scales are: E10, E20, E80, E120, N50, and N200.



Note: The N50 scale is not available at this time.

For any issues, please email EZBeadAssist@lifetech.com for assistance.

SOLiD® EZ Bead™ Amplifier install

To install the software application updates for the SOLiD® EZ Bead™ Amplifier follow these steps:

1. Save the “Amplifier_v1_9.exe” file to a USB stick.
2. Scan the USB stick for viruses, using a computer that has the most up-to-date virus definitions.
3. Power off the Amplifier.
4. Insert the USB stick into the Amplifier netbook computer, copy the “Amplifier_v1_9.exe” file from the USB stick, and paste it into the computer desktop.
5. Double-click **Amplifier_v1_9.exe** to begin installation.
6. Follow the prompts.
7. Select **I accept the License Agreement** and click **Next**.
8. Select **I accept the above 2 License Agreement(s)** and click **Next**.

9. Follow the prompts and click **Next** to complete the installation.
Note: The DOS window opens and closes when the upgrade is completed.

10. Close all the open windows and reboot the netbook.

11. Log in to the netbook computer.

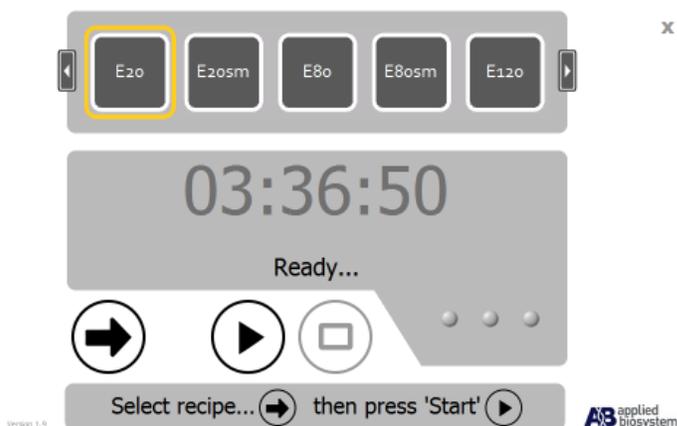
12. Power on the instrument.

13. From the desktop, double-click SOLiD® EZ Bead™ Amplifier icon  to launch the software.

14. Ensure that the splash screen displays the new version (1.9) on the lower right-hand corner of the screen.



15. Ensure that the GUI displays version 1.9 on the lower left-hand corner.



If you have any issues connecting to the instrument, please see the *Applied Biosystems SOLiD® EZ Bead™ Amplifier Getting Started Guide* (PN 4443494).

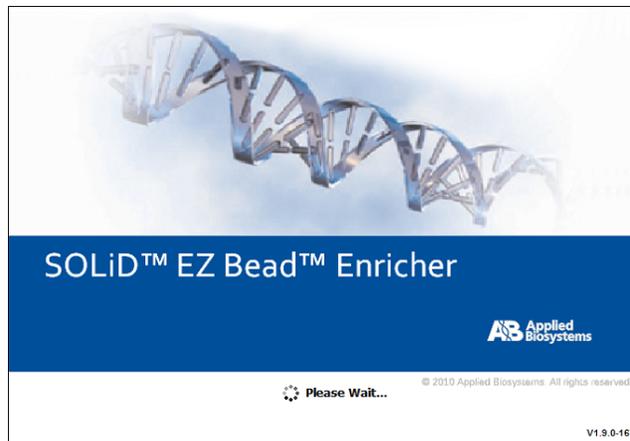
For any other issues, please email EZBeadAssist@lifetech.com for assistance.

SOLiD® EZ Bead™ Enricher install

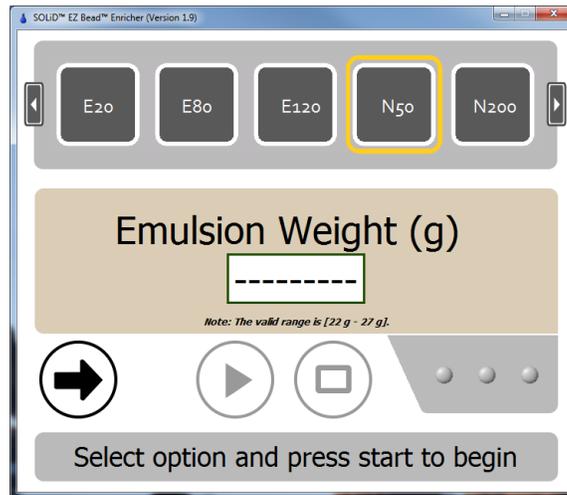
To install the software application updates for the SOLiD® EZ Bead™ Enricher follow these steps:

1. Save the “Enricher_v1_9.exe” file to a USB stick.

2. Scan the USB stick for viruses, using a computer that has the most up-to-date virus definitions.
3. Power off the Enricher.
4. Insert the USB stick into the Enricher laptop, copy the “Enricher.exe” file from the USB stick, and paste it into the laptop desktop.
5. Double-click **Enricher_v1_9.exe** to begin installation.
6. Follow the prompts.
7. Select **I accept the License Agreement** and click **Next**.
8. Select **I accept the above 2 License Agreement(s)** and click **Next**.
9. Follow the prompts and click **Next** to complete the installation.
Note: The DOS window opens and closes when the upgrade is completed.
10. Close all the open windows and reboot the laptop.
11. Log in to the laptop.
12. Power on the instrument.
13. From the desktop, double-click SOLiD® EZ Bead™ Enricher icon  to launch the software.
14. Ensure that the splash screen displays the new version (1.9) on the lower right-hand corner of the screen.



15. Ensure that the GUI displays version 1.9 in the top left corner.



If you have any issues connecting to the instrument, please see the *Applied Biosystems SOLiD® EZ Bead™ Enricher Getting Started Guide* (PN 4441486).

For any other issues, please email EZBeadAssist@lifetech.com for assistance.

Install Software Version 1.9
Install the software applications updates

Documentation and Support

Related documentation

To obtain the following related documents, contact your Life Technologies representative.

Document and part number	Description
<i>Applied Biosystems SOLiD[®] EZ Bead[™] Emulsifier Quick Reference Card (PN 4441487)</i>	Provides brief, step-by-step procedures on how to perform emulsion PCR. It is designed to help you quickly learn to use the SOLiD [®] EZ Bead [™] Emulsifier.
<i>Applied Biosystems SOLiD[®] EZ Bead[™] Emulsifier Getting Started Guide (PN 4441486)</i>	Describes how to perform emulsion PCR using the SOLiD [®] EZ Bead [™] Emulsifier.
<i>Applied Biosystems SOLiD[®] EZ Bead[™] Emulsifier Site Preparation Guide (PN 4452499)</i>	Provides information you need to fully prepare your site for the arrival and primary installation of the SOLiD [®] EZ Bead [™] Emulsifier.
<i>Applied Biosystems SOLiD[®] EZ Bead[™] Amplifier Quick Reference Card (PN 4443495)</i>	Provides brief, step-by-step procedures on how to perform emulsion PCR. It is designed to help you quickly learn to use the SOLiD [®] EZ Bead [™] Amplifier.
<i>Applied Biosystems SOLiD[®] EZ Bead[™] Amplifier Getting Started Guide (PN 4443494)</i>	Describes how to perform emulsion PCR using the SOLiD [®] EZ Bead [™] Amplifier.
<i>Applied Biosystems SOLiD[®] EZ Bead[™] Amplifier Site Preparation Guide (PN 4454035)</i>	Provides information you need to fully prepare your site for the arrival and primary installation of the SOLiD [®] EZ Bead [™] Amplifier.
<i>Applied Biosystems SOLiD[™] EZ Bead[™] Enricher Quick Reference Card (PN 4443497)</i>	Provides brief, step-by-step procedures on how to perform emulsion PCR. It is designed to help you quickly learn to use the SOLiD [®] EZ Bead [™] Enricher.
<i>Applied Biosystems SOLiD[™] EZ Bead[™] Enricher Getting Started Guide (PN 4443496)</i>	Describes how to perform emulsion PCR using the SOLiD [®] EZ Bead [™] Enricher.
<i>Applied Biosystems SOLiD[™] EZ Bead[™] Enricher Site Preparation Guide (PN 4454036)</i>	Provides information you need to fully prepare your site for the arrival and primary installation of the SOLiD [®] EZ Bead [™] Enricher.

Note:

For additional documentation, see [“Obtaining support” on page 14](#).

Obtaining support

For support for the *EZ Bead™ System Software v1.9* update, please email EZBeadAssist@lifetech.com for assistance.

For the latest services and support information for all locations, go to:

www.lifetechnologies.com

At the Life Technologies web site, you can:

- Access worldwide telephone and fax numbers to contact Life Technologies Technical Support and Sales facilities.
- Search through frequently asked questions (FAQs).
- Submit a question directly to Technical Support.
- Order Life Technologies user documents, SDSs, certificates of analysis, and other related documents.
- Download PDF documents.
- Obtain information about customer training.
- Download software updates and patches.

In addition, the Support page provides access to worldwide telephone and fax numbers to contact Life Technologies Technical Support and Sales facilities.

Headquarters

5791 Van Allen Way | Carlsbad, CA 92008 USA | Phone +1 760 603 7200 | Toll Free in USA 800 955 6288

For support visit www.appliedbiosystems.com/support

www.lifetechnologies.com

