

I. Purpose

This document covers the procedure to install the SMART-Seq® Pro application workflow as a built-in option in the ICELL8® cx CELLSTUDIO™ software.

NOTE:

- It is strongly recommended that instead of running this protocol, you upgrade to the latest version of the ICELL8 cx CELLSTUDIO Software. As of **Version 2.5.27.0**, the SMART-Seq Pro application workflow comes preconfigured. If needed, please contact [technical support](#) for the instructions on how to upgrade.
- **If you are running Version 2.5.27.0 or later and are not seeing the SMART-Seq Pro tab**, refer to Section II.C, starting with [Step 6](#), for instructions on how to enable the application. If the SMART-Seq Pro application is not listed there, then start with [Section II.B](#), "Download Required Installation File".
- This procedure should only be used for versions of ICELL8 cx CELLSTUDIO software **2.5.6.0 OR LATER**. If you are running an earlier version of the CELLSTUDIO software, please upgrade by contacting technical support.

II. Protocol

NOTE: The ICELL8 cx instrument does not need to be turned off for this procedure.

A. Verify the ICELL8 cx CELLSTUDIO Software Version

1. Run the current ICELL8 cx CELLSTUDIO software (CELLSTUDIO software) on the system PC.
2. Navigate to **Help > About**

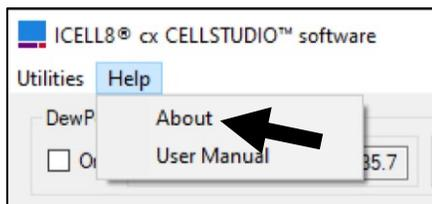


Figure 1. Finding Help > About in the ICELL8 cx CELLSTUDIO software interface.

3. Verify the version displayed in the *About* window is Version 2.5.6.0 or later.

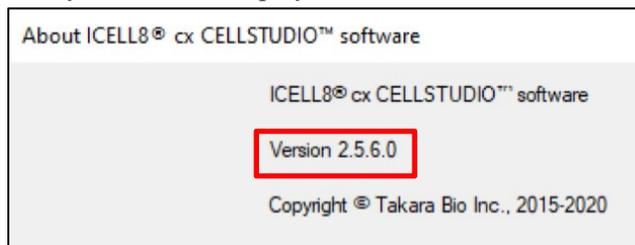


Figure 2. Partial view of the ICELL8 cx CELLSTUDIO software *About* window. The version information is highlighted by the red box.

4. If the version is earlier than 2.5.6.0, please stop process immediately and contact [technical support](#) for more instructions on the software upgrade process.

If the software version is 2.5.27.0 or later, proceed to Section II.C, [Step 6](#).

If the software version is between 2.5.6.0 and 2.5.27.0, close CELLSTUDIO software and proceed to the next section (Section II.B).

B. Download Required Installation File

If the ICELL8 cx computer is connected to the Internet:

1. Download the SMART-Seq_Pro_Application_XML_for_CELLSTUDIO_2.5.6.0.zip file from the Takara Bio website at the URL:

<https://www.takarabio.com/resourcedocument/x333121>

Save the zip file to a location on the instrument PC where you can easily find it.

If the ICELL8 cx computer is NOT connected to the Internet:

1. On a different computer with Internet connectivity, plug in a virus-free USB flash drive or external hard drive (ext-HD) that connects to the computer via a USB port.
2. Download the SMART-Seq_Pro_Application_XML_for_CELLSTUDIO_2.5.6.0.zip file from the Takara Bio website at the URL:

<https://www.takarabio.com/resourcedocument/x333121>

Save the zip file to a location on the external drive (flash drive or ext-HD) where you will be able to locate it easily later.

3. Transfer the USB flash drive or ext-HD to the ICELL8 cx instrument PC.
4. Copy the downloaded file from the external drive to a location on the instrument PC where you can easily find them.

C. Install the SMART-Seq Pro Application

1. From the location on the target computer where the zip file was saved (Section II.B), unzip the file. It should contain the file SMART-Seq Pro.xml.
2. Copy the SMART-Seq Pro.xml file, either by doing [CTRL] [C] or right-click on the file name in *File Explorer* and select **Copy**.

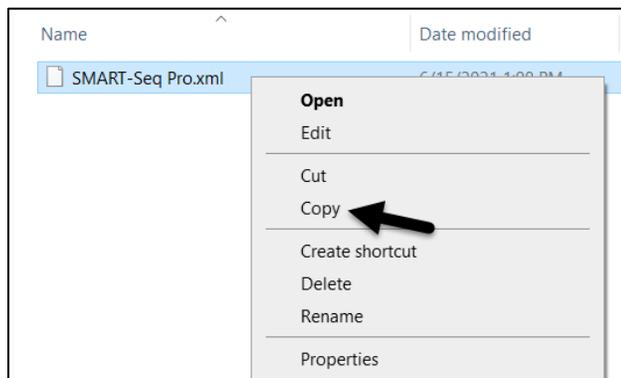


Figure 3. Right-mouse click menu to select Copy.

3. Navigate to the directory:

C:\ProgramData\takara\ICELL8 cx\Apps\

NOTE: By default, C:\ProgramData\ is a hidden directory that will not display in File Explorer. The easiest way to navigate to it is by copy and pasting or typing the file path into the input bar in File Explorer (see Figure 4).

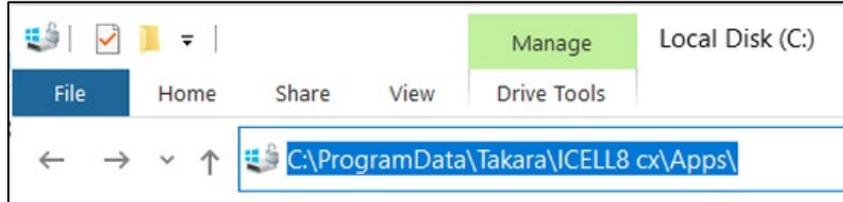


Figure 4. C:\ProgramData\takara\ICELL8 cx\Apps\ directory path in the File Explorer input field.

4. Paste the SMART-Seq Pro.xml file into the Apps\ directory, either by doing [CTRL] [V] or by right-mouse clicking in the File Explorer window and selecting **Paste**.

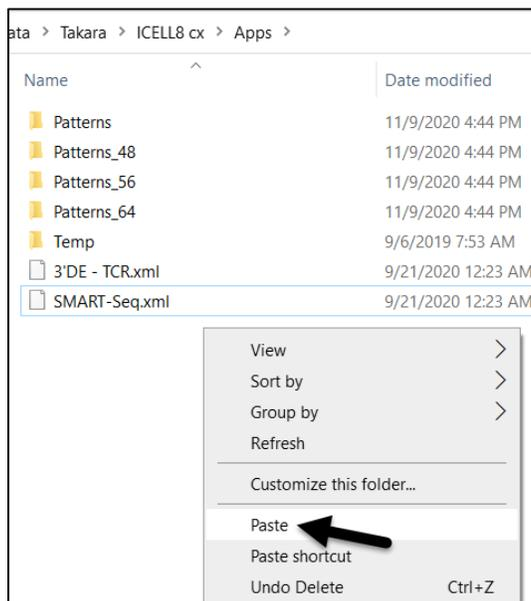


Figure 5. Right-mouse click menu to select Paste.

5. Restart CELLSTUDIO software.
6. Navigate to **Utilities > Application Manager**

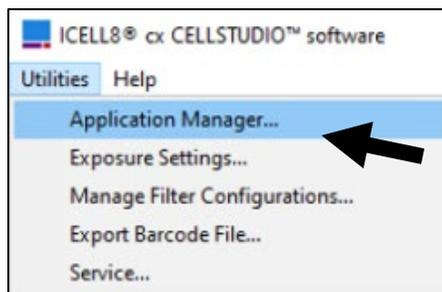


Figure 6. Finding Utilities > Application Manager in the CELLSTUDIO software interface.

SMART-Seq® Pro Application Installation Protocol-At-A-Glance

- In the *Application Manager* window, find SMART-Seq Pro in the list of applications and check the “Visible” box.

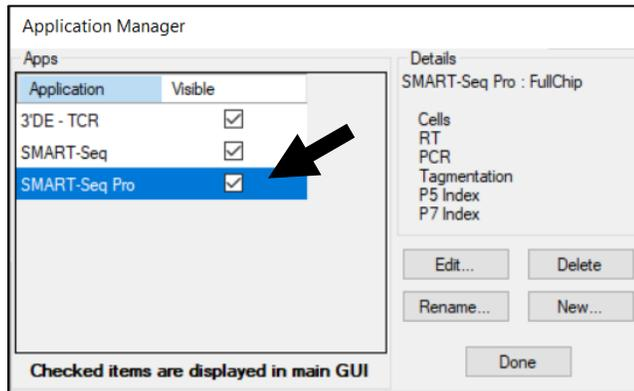


Figure 7. Check the “Visible” box for the SMART-Seq Pro application in the *Application Manager* apps list.

- Click the [Done] button.

A new tab, *SMART-Seq Pro*, should display in the main CELLSTUDIO software interface.

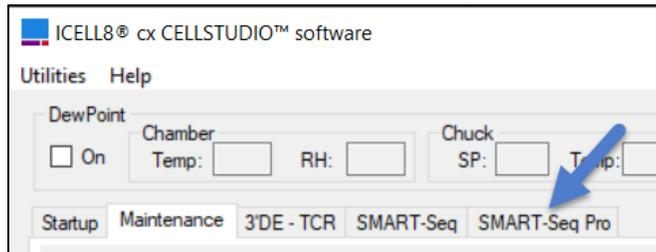


Figure 8. *SMART-Seq Pro* tab in the ICCELL8 cx CELLSTUDIO software after upgrade.

For more information about using the application in CELLSTUDIO software, refer to the [SMART-Seq Pro Application User Manual](#).

| Contact Us | |
|--|--|
| Customer Service/Ordering | Technical Support |
| tel: 800.662.2566 (toll-free) | tel: 800.662.2566 (toll-free) |
| fax: 800.424.1350 (toll-free) | fax: 800.424.1350 (toll-free) |
| web: takarabio.com/service | web: takarabio.com/support |
| e-mail: ordersUS@takarabio.com | e-mail: technical_support@takarabio.com |

Notice to Purchaser

Our products are to be used for **Research Use Only**. They may not be used for any other purpose, including, but not limited to, use in humans, therapeutic or diagnostic use, or commercial use of any kind. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without our prior written approval.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product’s web page at takarabio.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

© 2022 Takara Bio Inc. All Rights Reserved.

All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

This document has been reviewed and approved by the Quality Department.