

ICELL8® cx Single-Cell System Daily Maintenance Quick Start Guide

The following information is provided as a high-level introduction to the daily startup, clean-up, and system shutdown procedures of the ICELL8 cx Single-Cell System. For more detailed information, please refer to the [ICELL8 cx Single-Cell System User Manual](#).

Required Materials

- Deionized, filtered water
- 0.2% sodium hypochlorite solution
- Secondary containers (e.g., 1-L or 500-ml bottles)
- Kimwipes or similar dry wipes

Daily Startup

NOTE: If powering on the system, it takes ~10 min for the system to stabilize. Plan for this delay prior to proceeding to Step B.

If the system is already on, such as being left on overnight, skip to Step B.

A. Power on the system

1. If the instrument is powered down, turn it on.
2. If the system computer is shutdown, turn it on.
3. Start the CELLSTUDIO™ software on the computer.

B. Check and fill system containers

1. Ensure the water reservoir bottle and humidifier reservoir are each filled with at least 800 ml deionized, filtered water. Refill as needed. After the reservoirs have been refilled, screw the caps back on.
2. Check that the waste container is no more than half-full. If so, dispose of the liquid waste and reinstall the empty waste container.
3. Check the wash bottle. If the volume is <50% or if the solution is more than a week old, refill.
4. In the CELLSTUDIO interface in *Startup*, click the [Done] button for each of these four Daily Checklist categories when completed.

C. Run the [Daily Warmup]

1. In the Daily Checklist under *Startup*, click the “Daily warmup” [Start] button.
2. During warmup, monitor the syringes for trapped bubbles. If you see bubbles, repeat the daily warmup process up to two times.

NOTE: If large bubbles are still observed after three warmup cycles, refer to the User Manual.

D. Run the [Tip Clean] procedure


1. Click the *Maintenance* tab.
2. If the “Wash” well of the instrument trough is dry, click the [Wash Prime] button to fill it.
3. Click the [Tip Clean] button.

After the tip wash is complete, the system is ready for use.

Cleaning the ICELL8 cx system

1. After completing each dispense cycle, remove the 384-well source plate from the plate nest and properly dispose of it.
 2. Visually inspect the dispensing platform for any debris and remove carefully.
 3. Moisten a cloth with 70% isopropyl alcohol and wipe down the dispensing platform (chip nest, source plate nest, trough, and the areas on the base platform immediately surrounding them).
- NOTE: Do not wipe the plastic side panels.**
4. After the final dispense of the day, perform the [Tip Clean] procedure (Step D, above).

System Shutdown Procedure

1. Check that all nanowell chips and source plates have been removed from the instrument stage module.
2. Clean the ICELL8 cx instrument (see previous procedure).
3. Empty the contents of the water and humidifier reservoirs.
4. Press the [System Prime] button to empty the system tubing of water.
5. Dispose of the contents of the wash bottle.
6. In CELLSTUDIO software in the *Maintenance* tab, press [Wash Prime] until there is no new fluid coming up from the wash well.
7. Exit CELLSTUDIO software by clicking the [Close] button () at the top-right corner of the window.
8. Turn off the ICELL8 cx instrument by toggling the power switch on back of the system and power down the computer.
9. Properly dispose of contents of waste container, rinse, and allow it to dry.
10. Remove the back cover, connect a tube to the drain valve on the humidifier tank, and drain the humidifier.

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.

Takara Bio USA, Inc.

United States/Canada: +1.800.662.2566 • Asia Pacific: +1.650.919.7300 • Europe: +33.(0)1.3904.6880 • Japan: +81.(0)77.565.6999

© 2021 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.

07.21 US