

iSort™ Automated Cell Sorter

Publication Number MAN0017638

Revision A.0

Site preparation workflow 1

Site preparation checklist 2

Pre-installation requirements..... 3

Operating environment and site requirements 4

Safety requirements 5

Materials required for installation..... 6

Prepare for installation 6

Related documentation and support 7

This guide contains the information to prepare your site for installation of the iSort™ Automated Cell Counter.

Site preparation workflow

IMPORTANT! Do not install the iSort™ Automated Cell Sorter yourself. A Thermo Fisher Scientific representative will contact you to schedule the installation.

When the installation is scheduled:

1. Receive and inspect the system (page 6).
2. Move the crated instrument to the installation site (page 6).
3. Complete the site preparation checklist before the installation date (page 2).

Timeline and training

Installation of the iSort™ Automated Cell Sorter is performed by a Thermo Fisher Scientific Field Service Engineer. Once successfully installed, the engineer performs the necessary installation tests. This entire process takes approximately two days.

During and/or after installation, the Thermo Fisher Scientific Service representative will walk through basic instrument maintenance functions. Additional training is provided separately by a Thermo Fisher Scientific Applications specialist.

Site preparation checklist

IMPORTANT! Complete all items in the checklist before the scheduled installation date. If the site preparation checklist is not complete when the Thermo Fisher Scientific service representative arrives, the scheduled installation may be delayed.

✓	Site preparation requirement	Page
	Customer responsibilities have been reviewed and personnel assigned.	3
	The installation site is identified and meets requirements: <input type="checkbox"/> Space and clearance <input type="checkbox"/> Environmental <input type="checkbox"/> Electrical <input type="checkbox"/> Network <input type="checkbox"/> Safety	4
	Materials that are needed for installation and operation are available.	6
	The system was received and inspected: <input type="checkbox"/> All items on the shipping list are the same items that were ordered at the time of purchase. <input type="checkbox"/> Any damage to shipping containers was recorded on the shipping containers	6
	The installation site is cleared and ready for instrument installation.	6
	The crated instrument and other shipping containers are moved to the installation site.	6

Pre-installation requirements

Personnel	Responsibilities
Instrument maintenance personnel	<ul style="list-style-type: none"> • Reviews the site preparation guide for safety information and system requirements. • Coordinates personnel and tasks. • Orders required materials. • Selects the site. • Reviews checklists with appropriate personnel, then with the Thermo Fisher Scientific service representative to ensure that the site is properly prepared. • Receives and inspects the system. • Schedules the installation and notifies personnel of the installation date. • Ensures that the site is clear of unnecessary material on the installation day. • Is available to help the service representative throughout installation. • Ensures that installation requirements are met for: <ul style="list-style-type: none"> – Space at the installation site – Building clearances – Temperature and humidity – Waste collection – Electrical supply – Computer – Required safety equipment (page 5) and installation materials (page 6) • If possible, move the crated system to the site before the installation date. • Available to help service representative and laboratory personnel throughout installation. • Review safety information. • Ensure that all customer-provided materials for installation are present at the site. • Primary users (responsible for training other users) are available during the installation, so that they can receive initial training on the instrument. • Reviews the user guide for safety information. • Ensures that the required safety practices and equipment are in place. • Is in the vicinity and available to the Thermo Fisher Scientific service representative when the service representative is at your facility.
Network or IT specialist	<ul style="list-style-type: none"> • Ensures that one active, tested wireless local area network (WLAN) is in place for Wi-Fi connection before the scheduled installation date. • Is available during installation to connect the system to the network. The Wi-Fi connections will also be used for remote troubleshooting if a problem arises. <p>Note: The Thermo Fisher Scientific Field Service Engineer will connect the Wireless Wi-Fi Dongle – Linksys™ AE3000 N900 Dual-Band Wireless-N USB Adapter during installation.</p>

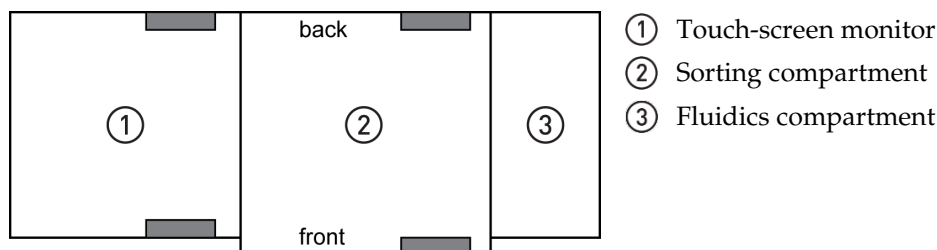
Operating environment and site requirements

- The iSort™ Automated Cell Sorter has a footprint (W × D × H) of approximately 67 cm × 41 cm × 49 cm (26.5 in × 16 in × 19 in) with the fluidics compartment door and the top cover closed.

With the top cover open, the height of the instrument is 74 cm (29 in).

With the fluidics compartment door open, the entire system requires a total bench width of ~135 cm (53.2 in).

- The setup for the iSort™ Automated Cell Sorter is shown below. The gray rectangles in the diagram represent the four handholds in the base that are used when lifting the instrument out of the box.



IMPORTANT! Do **not** lift the iSort™ Automated Cell Sorter from the fluidics compartment. Lifting from the fluidics compartment can damage the instrument. Lift the instrument using the designated handholds.

- The bench or table top must be able to accommodate 43 kg (95 lbs), the weight of the instrument plus the weight of the full 2-liter fluid bottles (sheath fluid and waste).
- Place the iSort™ Automated Cell Sorter on a level surface away from vibrations from other pieces of equipment. Tabletop centrifuges, vortex mixers, and other laboratory equipment can vibrate the instrument during a run and cause a decrease of instrument performance.
- The area should be free of excessive dust or moisture. Position the instrument so that the power button and main power supply can be reached easily.
- Allow at least 6.5 cm (2.5 in) free space at the back of the instrument to allow for proper ventilation and prevent overheating of electronic components.
- Ambient temperature range: 18°–25°C (64°–77°F)
- Storage temperature range: 5°–35°C (41°–95°F)
- Relative humidity range: 20–70%
- Electrical input: 100–240 VAC, 4.0 A
- Frequency: 50–60 Hz

Safety requirements

Safety practices

A safety representative from your facility must ensure that:

- Personnel establish and follow all appropriate safety practices and policies to protect laboratory personnel from potential hazards.
- All appropriate safety devices and equipment are always available.

Required safety equipment

The following safety protection and equipment must be available at the installation site:

- Protection from any sources of hazardous chemicals, radiation (for example, lasers, radioisotopes, radioactive wastes, and contaminated equipment), and potentially infectious biological material that may be present in the area where the Thermo Fisher Scientific service representative will work.
- Appropriate fire extinguisher appropriate for use on electrical and chemical fires as specified in current codes, regulations, and/or standards, and with approval of the Fire Marshall or other authority having jurisdiction.
- Eyewash
- Safety shower
- Eye and hand protection
- Sufficient ventilation, including vent line/fume hood, if applicable
- Biohazard waste container, if applicable
- First-aid equipment
- Spill cleanup equipment
- Applicable Safety Data Sheets (SDSs)



WARNING! BIOHAZARD. Cell sorting creates aerosols. Exposure to and inhalation of aerosols can potentially be hazardous depending on the cell types you are using. Consult your institution's safety officer for guidelines related to your individual samples.

The iSort™ Automated Cell Sorter may need to be placed in a biosafety cabinet and/or connected to an evacuator filter/pre-filter or liquid disinfectant trap, depending on your institutions safety requirements.

We recommend that you follow the International Society for the Advancement of Cytometry (ISAC) guidelines for cell sorter biosafety standards (www.ncbi.nlm.nih.gov/pmc/articles/PMC4117398/).

Materials required for installation

- Safety glasses, lab coats, and chemical-resistant, disposable gloves (powder-free)
- Lint-free tissues
- Isopropanol, HPLC-grade or better
- 10% bleach solution
- Deionized water
- PBS (1X), pH 7.4 (flow cytometry grade)
- PBS (10X), pH 7.4
- Tween™ 20 (50% solution)
- Attune™ Performance Tracking Beads
- Attune™ Wash Solution
- Decon™ Contrad™ 70 Liquid Detergent
- Texwipe™ TX761MD Microdenier Swabs
- 5-mL and 50-mL Falcon™ polypropylene tubes
- 1.5-mL centrifuge tubes
- 0–10 µL and 40–1000 µL pipettes (used in the instrument performance test)
- Analytical scale (with 3 to 4 decimal places) (to weigh polypropylene tubes)
- Ultrasonic cleaning bath (such as Fisherbrand™ CPX Series Digital Ultrasonic Cleaning Bath or similar) (used for nozzle tip cleaning)

Prepare for installation

- Receive and inspect the shipment** Carefully inspect the shipping containers and report any damage to the Thermo Fisher Scientific service representative. Record any damage or mishandling on the shipping documents.
- Move the crated instrument to the installation site**
1. Clear the installation site.
 2. If possible, move the crated instrument and other shipping containers to the installation site. **Do not uncrate.**



CAUTION! PHYSICAL INJURY HAZARD. The weight of the iSort™ Automated Cell Sorter is 41 kg (89 lbs). Do not attempt to lift or move the instrument without the help of others, the use of appropriate moving equipment, and proper lifting techniques. Improper lifting can cause painful and permanent back injury. Depending on the weight, moving or lifting an instrument may require two or more persons

IMPORTANT! Do not subject the iSort™ Automated Cell Sorter to sudden impact or excessive vibration. Handle the instrument with care to prevent damage.

Related documentation and support

Related documentation

The publication numbers in this section are for the latest product versions available at the time of publication. For documentation for newer product versions, go to thermofisher.com/support.

Document	Pub. No.
iSort™ Automated Cell Counter User Guide	MAN0016718
iSort™ Automated Cell Counter Quick Reference	MAN0017190

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation, including:
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)


Note: For the SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



IMPORTANT! Wiping the iSort™ Automated Cell Sorter computer (i.e., erasing the hard drive to remove all programs, files, and the operating system) voids the product warranty.

 **Manufacturer:** Cytonome | 9 Oak Park Dr. | Bedford, MA 01730

The information in this document is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: MAN0017638

Revision	Date	Description
A.0	19 October 2018	New site preparation guide for the iSort Automated Cell Sorter.

Important Licensing Information: This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

© 2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Kimwipes is a registered trademark of Kimberly-Clark Corporation. Linksys is a registered trademark of Belkin International, Inc.

thermofisher.com/support | thermofisher.com/askaquestion

thermofisher.com

ThermoFisher
S C I E N T I F I C