

Attune™ NxT Flow Cytometer

Pub. No. 100024428 Rev. C.0

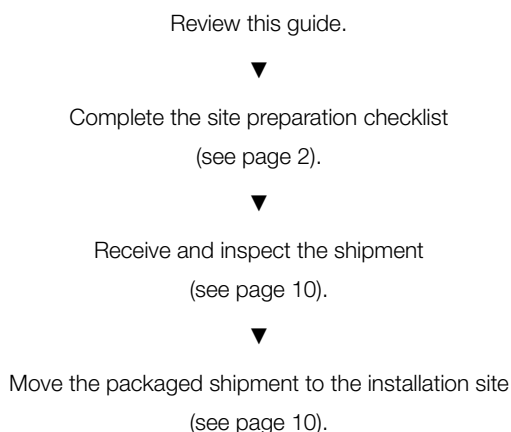
This guide contains the information needed to prepare your site for installation of the Attune™ NxT Acoustic Focusing Cytometer.

- Site preparation workflow 1
- Installation time and training 1
- Site preparation checklist 2
- Customer responsibilities 3
- Site requirements 4
- Materials for installation and operation 9
- Receive and inspect the shipment 10
- Move the packaged shipment to the installation site 10
- Register to access the Connect 11
- Related documentation 11

Site preparation workflow

IMPORTANT! Thermo Fisher Scientific does not install, service, or repair products in area designated BioSafety Level 3 (BSL-3) or BioSafety Level 4 (BSL-4).

A Thermo Fisher Scientific service representative will contact you to schedule the installation. When the installation date is scheduled, perform the following tasks.



Installation time and training

After the shipment is unpacked, the installation takes approximately four hours. When the system reaches proper operating status, the Thermo Fisher Scientific service representative returns to perform installation tests.


During and/or after installation, the service representative calibrates the instrument, performs run verification, reviews data, and provides some basic operator training. For additional training and reference information, see the user documents that are provided with the product.

Site preparation checklist

IMPORTANT! Complete, date, and initial all items in the following checklist before the scheduled installation date. After completion, return all pages of the completed checklist to: <mailto:concierge@thermofisher.com>. If the site preparation checklist is not complete when the service representative arrives, the scheduled installation may be postponed.

✓	Date	Initials	Site preparation requirement	See page
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	Customer responsibilities have been reviewed.	3
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	Personnel have been assigned tasks and responsibilities.	
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	The installation site is identified and meets the following requirements:	
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> Space and clearance	5
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> Environmental	6
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> Electrical	7
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> Network	8
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> Safety	8
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	If required, antivirus software is available for installation on the computer provided with the product.	9
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	The shipment was received and inspected as follows:	10
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> The items shown on the shipping list are the items that were ordered at the time of purchase.	
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> Damage to shipping containers was reported to the shipping company that delivered the shipment and to your service representative.	
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> Damage or mishandling was recorded on the shipping documents.	
	<input type="text"/>	<input type="text"/>	<input type="checkbox"/> If provided with the shipment, all reagents and plates are unpacked and stored as specified on package labels.	
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	The installation site is cleared and ready for the installation.	10
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	The packaged shipping containers are moved to the installation site.	
<input type="checkbox"/>	<input type="text"/>	<input type="text"/>	All materials for installation, qualification, and operation are available.	9

Customer responsibilities

Personnel	Responsibilities
Site preparation/installation coordinator	<ul style="list-style-type: none"> • Reviews the site preparation guide for safety information and system requirements. • Coordinates personnel and tasks. • Chooses the site. • Reviews checklists with applicable personnel, then with the Thermo Fisher Scientific service representative to verify that the site is properly prepared.^[1] • Receives and inspects the system. • Stores the Attune™ Starter Kit. • Schedules the installation and informs personnel of the installation date. • Ensures that the site is clear of unnecessary material on the installation day. • Is available to assist the service representative throughout installation.^[1]
Laboratory safety representative	<ul style="list-style-type: none"> • Reviews the site preparation guide for safety information. • Ensures that the required safety practices and equipment are in place. • Is in the vicinity and available to the service representative at all times while the service representative is at your facility. • Ensures that primary users (responsible for training other users) are available for training during the installation.
Laboratory personnel/primary users	<ul style="list-style-type: none"> • 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 be trained on the instrument.
Facilities personnel	<ul style="list-style-type: none"> • Ensure that installation requirements are met for: <ul style="list-style-type: none"> – Space at the installation site – Building clearances – Humidity and temperature – Waste collection – Electrical supply – Computer – Safety and installation materials • If possible, moves the packaged shipment to the site before the installation date. • Is available to assist service representative and laboratory personnel throughout installation. • If applicable, ensures that at least two people are available to help the service representative move and position the system.
Network or IT specialist (if the system will be connected to a network)	<ul style="list-style-type: none"> • Ensures that active, tested local area network (LAN) connections are in place before the scheduled installation date. • Ensures that network hardware is compatible with an RJ45-type connector. • If necessary, supplies additional cables. • Is available during installation to connect the system to the network. • If applicable, provides and installs a network or dedicated printer. <p data-bbox="329 1478 386 1528"></p> <p data-bbox="402 1478 1529 1528">CAUTION! Do not attempt to connect the system components to the network before the Thermo Fisher Scientific service representative arrives.</p>

^[1] Required for service representative installation of the instrument.

Site requirements

Dimensions and weights

To prepare for installation, provide space for receipt and configuration of the components listed in this section. This section provides dimensions and weights for the packages you will receive, and it describes the dimensions of the components after installation and configuration.

Components (packaged)

The Attune™ NxT Flow Cytometer and components are packed onto one pallet. Ensure that the building clearances allow for transport of the packaged components.

Table 1 Pallet dimensions

Height	Length (depth)	Width	Weight
69.8 cm (27.5 in.)	63.5 cm (25.0 in.)	76.2 cm (30.0 in.)	59.9 kg (132 lb)

CAUTION! PHYSICAL INJURY HAZARD. Do not attempt to lift or move the instrument without the assistance 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 people.

Typical system component layouts

IMPORTANT! We do not install, service, or repair Thermo Fisher Scientific instruments in areas designated BioSafety Level 3 (BSL-3) or BioSafety Level 4 (BSL-4).

The following images (not to scale) show typical system component layouts with various optional peripherals.

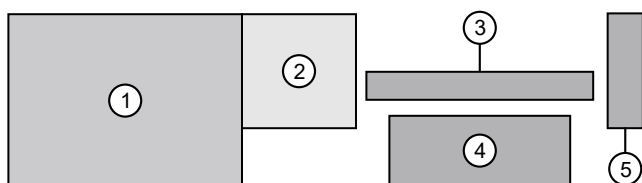


Figure 1 Attune™ NxT Flow Cytometer with Attune™ NxT Auto Sampler

- ① Attune™ NxT Flow Cytometer
- ② (Optional) Attune™ NxT Auto Sampler
- ③ Monitor
- ④ Keyboard
- ⑤ CPU tower

Note: When the system includes the optional Attune™ NxT Auto Sampler, the computer must be placed to the right of the cytometer and the auto sampler.

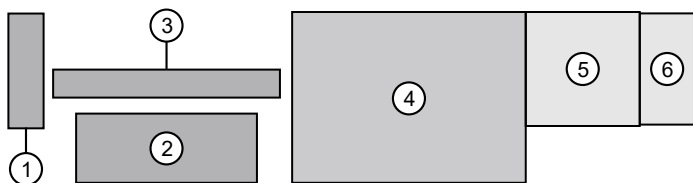


Figure 2 Attune™ NxT Flow Cytometer with CytKick™ / CytKick™ Max Autosampler

- ① CPU tower
- ② Monitor
- ③ Keyboard
- ④ Attune™ NxT Flow Cytometer
- ⑤ (Optional) CytKick™ / CytKick™ Max Autosampler
- ⑥ (Optional) CytKick™ / CytKick™ Max Autosampler Side Car bottle module (attached to the right side of the autosampler)

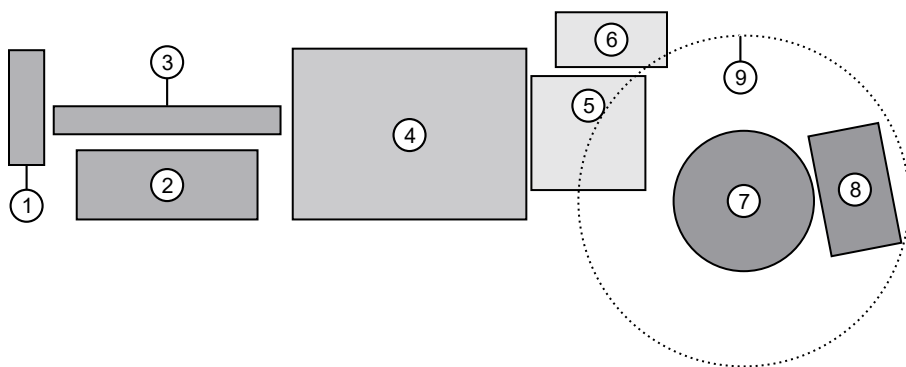


Figure 3 Attune™ NxT Flow Cytometer with CytKick™ / CytKick™ Max Autosampler and Orbitor™ RS2 Microplate Mover

- ① CPU tower
- ② Monitor
- ③ Keyboard
- ④ Attune™ NxT Flow Cytometer
- ⑤ (Optional) CytKick™ / CytKick™ Max Autosampler
- ⑥ (Optional) CytKick™ / CytKick™ Max Autosampler Side Car bottle module (must be placed in the back to accommodate the microplate mover)
- ⑦ Orbitor™ RS2 Microplate Mover
- ⑧ Microplate storage carousel for microplate mover
- ⑨ Range of movement for the bi-directional telescoping arm of the microplate mover

System component dimensions and weights

Ensure that the installation site bench space can accommodate the dimensions and support the weights of the components.

Component	Height	Length (depth)	Width	Weight
Instrument	40.0 cm (15.8 in.)	43.5 cm (17.1 in.)	58.1 cm (22.9 in.)	~27.3 kg (60 lbs) empty ~33.8 kg (74.5 lbs) with fluids
Computer	29.0 cm (11.4 in.)	29.2 cm (11.5 in.)	9.3 cm (3.7 in.)	~11.5 kg (25.4 lbs)
Monitor	41.9 cm (16.5 in.)	20.3 cm (8.0 in.)	57.2 cm (22.5 in.)	~4.8 kg (10.6 lbs)
Keyboard	5.1 cm (2.0 in.)	17.8 cm (7.0 in.)	45.7 cm (18.0 in.)	~0.9 kg (2.0 lbs)
(Optional) Attune™ NxT Auto Sampler	40.0 cm (15.8 in.)	29.0 cm (11.4 in.)	29.0 cm (11.4 in.)	15.9 kg (35.1 lbs)
(Optional) CytKick™ / CytKick™ Max Autosampler ^[1]	Autosampler: 40.6 cm (16.0 in.) Side Car: 29.0 cm (11.5 in.)	Autosampler: 28.5 cm (11.25 in.) Side Car: 28.0 cm (11.0 in.)	Autosampler: 29.0 cm (11.5 in.) Side Car: 15.0 cm (5.9 in.)	16.9 kg (37.2 lbs) empty 20.9 kg (46 lbs) with focus and waste bottles at full capacity
(Optional) Orbitor™ RS2 Microplate Mover ^[2]	72.0 cm (28.4 in.)	35.6 cm (14.0 in.) unit diameter 40.6 cm (16.0 in.) reach for telescoping arm 88.2 cm (34.7 in.) operational reach (bi-directional) ^[3]		25.0 kg (55.1 lbs)
(Optional) Microplate storage carousel assembly ^[4]	Dimensions and weight of the microplate storage carousel assembly depend on the combination of the stacks and hotels used. Refer to the Orbitor™ RS2 Microplate Mover product literature for the dimensions and weights of specific components.			

^[1] If using the optional Orbitor™ RS2 Microplate Mover, the Side Car bottle module must be placed in the back of the autosampler to allow automation access to the tray area.

^[2] The minimum bench space required to support the cytometer, autosampler, and the automated microplate mover with microplate storage carousel depend on the combination of the stacks and hotels used.

^[3] See Thermo Scientific™ Orbitor™ RS2 Microplate Mover product literature for more information.

^[4] Can consist of sequential access stacks, random access hotels, or a combination of both.

CAUTION! PHYSICAL INJURY HAZARD. Do not attempt to lift or move the instrument without the assistance 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 people.

Component clearances required for installation and maintenance

During instrument setup and maintenance, it is necessary to access the back of the Attune™ NxT Flow Cytometer and the optional Attune™ NxT Auto Sampler or the CytKick™ / CytKick™ Max Autosampler. If the back of the products face a wall, ensure that there is sufficient clearance on the bench to rotate the products for access.

IMPORTANT! For safety reasons, the power outlet for the product must be accessible at all times.


Bench	Minimum clearance ^[1]
Depth	<ul style="list-style-type: none"> >100 cm (39.37 in.) for a bench against a solid vertical surface >25.4 cm (10 in.) of clearance at the back of the instrument for air flow, service access, and cable routing. If the bench is at least 25.4 cm. (10 in.) from a wall, the bench can be 76.2 cm. (30 in.) deep. If the bench has wheels, it can facilitate access to the back of the instrument.
Width	<p>>158 cm (62 in.) for the instrument, computer, and computer monitor.</p> <p>IMPORTANT! If the system includes the optional Attune™ NxT Auto Sampler, then add 40 cm (15.8 in.) to the width of the bench, for a total bench width of ~ 198 cm (77.8 in.). If the system includes the optional CytKick™ / CytKick™ Max Autosampler, then add 55 cm (21.7 in.) to the width of the bench, for a total bench width of ~ 213 cm (83.9 in.)</p>
Height	74 cm (29 in.) allows the hinged lid to be placed in the vertical open position. This provides access to the optical compartment and allows changing of filters.


^[1] If you are using the optional Orbitor™ RS2 Microplate Mover, the minimum bench space required to support the cytometer, autosampler, and the automated microplate mover with microplate storage carousel depends on the combination of stacks and hotels used for the carousel assembly.

Environmental requirements

Condition	Requirement
Installation site	Indoor use only
Altitude	Safety tested up to 2000 m (6500 ft)
Humidity (instrument and computer)	Operation: 15%–80% (noncondensing)
Temperature (instrument and computer)	<p>15°C to 30°C (60°F to 85°F)</p> <p>Note: The room temperature must not fluctuate more than 2°C over a 2-hour period.</p> <p>Avoid placing the instrument and computer adjacent to heaters, cooling ducts, or in direct sunlight.</p>
Thermal Output	<p>Hot-air exhaust is vented from the Attune™ NxT Acoustic Focusing Cytometer through the hot-air waste port on the rear panel. The hot-air exhaust is designed to dissipate heat produced by the instrument.</p> <p>The maximum thermal output of the Attune™ NxT Flow Cytometer is ~ 500 Btu/h (150 W). Consult your facilities department to determine if the laboratory ventilation system can maintain room temperature with this level of thermal output. If it can maintain room temperature during instrument operation, the hot-air exhaust port can be vented directly to room air.</p>
Overvoltage category	Installation categories II
Vibration	Do not place the instrument adjacent to strong vibration sources, such as a centrifuge, pump, or compressor. Excessive vibration will affect instrument performance.
Pollution degree	<p>II</p> <p>Install the instrument in an environment that has nonconductive pollutants such as dust particles or wood chips. Typical environments with a Pollution Degree II rating are laboratories and sales and commercial areas.</p> <p>Ensure the room is away from any vents that could expel particulate material on the components.</p>
Liquid waste collection	Dispose of the buffer, reagents and any liquid waste as hazardous waste in compliance with local and national regulations.

Electrical requirements

 **CAUTION!** Do not unpack or plug in any components until they are configured for the proper operating voltage by the service representative.

 **WARNING!** For safety, the power outlet for the instrument must be accessible at all times. See “Component clearances required for installation and maintenance” on page 6 for information about the space needed between the wall and the instrument. In case of emergency, you must be able to immediately disconnect the main power supply to all the equipment. Allow adequate space between the wall and the equipment so that the power cords can be disconnected in case of emergency.

- Dedicated line and ground between the instrument and the main electrical service.
- Maximum power dissipation: ~150 W (not including computer, monitor, and the optional autosampler)
- Mains AC line voltage tolerances must be up to ± 10 percent of nominal voltage

Device	Rated voltage	Circuit required	Rated frequency	Rated power	Maximum power
Instrument	100–240 $\pm 10\%$ VAC ^[1]	15 A	50/60 Hz	960 W	150 W
Computer (desktop)	100–240 $\pm 10\%$ VAC	15 A	50/60 Hz	150 W	300 W
Monitor				60 W	100 W
Autosampler (optional)	100–240 $\pm 10\%$ VAC	15 A	50/60 Hz	25 W	30 W
Orbitor™ RS2 Microplate Mover	100–240 $\pm 10\%$ VAC	15 A	50/60 Hz	50 W	100 W

^[1] If the supplied power fluctuates beyond the rated voltage, a power line regulator may be required. High or low voltages can adversely affect the electronic components of the instrument.

The Attune™ NxT Flow Cytometer can be configured for operating voltages between 100 and 240 VAC at 50 or 60 Hz. The instrument is equipped with a universal power supply. The instrument requires a 15 A circuit for all indicated input voltages.

IMPORTANT! The Attune™ NxT Flow Cytometer and the optional autosampler are shipped with up to three power connectors. These connectors require standard 15 A wall receptacles with proper grounding. Do not use extension cords.

Electrical protective devices

We recommend several protective devices in environments with large voltage and power fluctuations.

Recommended devices
Power line regulator <ul style="list-style-type: none">• 1.5-kVA power line regulator• Use in areas where the supplied power fluctuates in excess of $\pm 10\%$ of the normal voltage.• Power fluctuations can adversely affect the function of the instrument and computer. <p>Note: A power line regulator monitors the input current and adjusts the power supplied to the instrument or computer. It does not protect against a power surge or failure.</p>
Surge protector <ul style="list-style-type: none">• 10-kVA surge protector (line conditioner)• Use in areas with frequent electrical storms or near devices that are electrically noisy, such as refrigerators, air conditioners, or centrifuges.• Short-duration, high-voltage power fluctuations can abruptly terminate the function of, and thereby damage the components of, the computer and the instrument. <p>Note: A dedicated line and ground between the instrument, computer, and the building's main electrical service can also prevent problems caused by power fluctuations.</p>

Recommended devices

Uninterruptible power supply (UPS)

- 1.5-kVA uninterruptible power supply (UPS)
- Use in areas prone to power failure.
- Power failures and other events that abruptly terminate the function of the instrument and computer can corrupt data and possibly damage the system.



CAUTION! PHYSICAL INJURY HAZARD. Do not attempt to lift or move the UPS unit without the assistance of at least two people. Improper lifting can cause painful and permanent back injury. See the UPS manufacturer user guide for more information.

IMPORTANT! A UPS provides power for a limited time. It is meant to delay the effects of a power outage, not to serve as a replacement power source. In the event of a power loss, power off the instrument and computer unless you expect to regain power within the battery life of the UPS.

Network requirements

The instrument is factory-configured for IPv4 TCP/IP communication and uses an Ethernet adapter (100/1,000 Mbps) with an RJ45-type connector for local area network (LAN) connection. Connect the instrument to the computer using the USB 3.0 connector.

For the instrument to be connected to a LAN, an active, tested LAN connection must be in place before the scheduled installation date.

The assigned IT or network specialist from your organization must be available during the installation to help connect the instrument to your network.

IMPORTANT! Do not use the instrument on a wireless network. A wireless network may interfere with data collection, resulting in data loss.

Safety requirements

Safety practices

A safety representative from your facility must ensure that:

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

Required safety equipment

Your laboratory has specific safety practices and policies designed to protect laboratory personnel from potential hazards that are present. Follow all applicable safety-related procedures at all times.

The following safety equipment and protection from hazards 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 service representative will work.
- Appropriate fire extinguisher:
 - You are responsible for providing an appropriate fire extinguisher for use on or near the equipment.
 - The types and sizes of fire extinguishers shall be suitable 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.
 - The installation of appropriate fire extinguishers shall be in addition to other fire-protection systems and not as a substitute or alternative to them.
- Eyewash
- Safety shower
- Eye and hand protection
- Adequate ventilation, including vent line/fume hood, if applicable
- Biohazard waste container, if applicable
- First-aid equipment
- Spill cleanup equipment
- Applicable Safety Data Sheets (SDSs)

Third-party software

Before installing third-party software on the computer running the product software, confirm that the third-party software will not:

- Restrict Ethernet communication.
- Interfere with instrument or computer operation.

Antivirus software requirements

The computer provided with the instrument does not include antivirus software because customer preferences and network requirements vary. You are responsible for installing antivirus software of your choice to protect the computer against viruses.

IMPORTANT! You must disable or deactivate antivirus software and antispyware during use of the instrument. Antivirus and antispyware monitoring can interfere with data collection, resulting in data loss.

Materials for installation and operation

Installation materials

Ensure that the following materials are available before installation of the product:

- Safety glasses, lab coats, and chemical-resistant, disposable gloves (powder-free)
- Easily accessible specified power outlet
- *(Optional)* Electrical protective devices (universal power supply unit, surge protector, and/or power line regulator)
- Lint-free tissues
- 10% bleach solution
- Methanol or isopropanol, HPLC-grade or better
- Water, Milli-Q™ grade
- Three sizes of micropipettors and tips:
 - 1- μ L to 10- μ L
 - 10- μ L to 100- μ L
 - 100- μ L to 1,000- μ L
- Mini vortexer, centrifuge equipped to accept 15-mL and 50-mL centrifuge tubes, and sample tubes

Operation materials

Additional supplies and consumables are necessary for routine operation. Contact a sales representative to order these additional supplies. Use only supplies as specified by Thermo Fisher Scientific.

Receive and inspect the shipment

1. Verify that the items shown on the shipping list are the same items that you ordered at the time of purchase.
2. Carefully inspect the shipping containers. Report any damage to the shipping company and to your service representative. Record any damage or mishandling on the shipping documents.
3. Immediately unpack the Attune™ NxT Starter Kit (boxed separately from the instrument components). Store the components as specified.


The Attune™ NxT Starter Kit contains the following items:

- Software Kit (includes a valid DESkey device, which is required for the operation of the instrument software)
- Reagents:

Reagent	Cat. No.	Storage conditions	Usage conditions	Stability
Attune™ Focusing Fluid, 6 x 1 L	4449791	15°C to 30°C	15°C to 30°C	The focusing fluid is stable on the instrument for 30 days after the bottle has been opened.
Attune™ Wash Solution	A24974	15°C to 30°C		The wash solution is stable on the instrument for 30 days after the bottle has been opened.
Attune™ Shutdown Solution	A24975	15°C to 30°C		The shutdown solution is stable on the instrument for 30 days after the bottle has been opened.
Attune™ Performance Tracking Beads	4449754	2°C to 8°C		The beads are stable for 1 year, when stored as directed.


IMPORTANT! Reagents can be stored at colder temperatures, but ensure that all reagents are at 15°C to 30°C before running the instrument.


IMPORTANT! Other than the Attune™ NxT Starter Kit, which contains reagents that require storage at specific conditions, do not unpack the shipping containers at this time. To protect yourself from liability for damage that occurred during shipping, inspect the shipping containers and report damage as described above.

 **WARNING! CHEMICAL HAZARD.** Some chemicals used with Thermo Fisher Scientific instruments are potentially hazardous and can cause injury, illness, or death. Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, work with, or dispose of any chemicals or hazardous materials

Move the packaged shipment to the installation site

1. Clear the installation site of all unnecessary materials.
2. Move the packaged shipment to the installation site.

 **CAUTION! PHYSICAL INJURY HAZARD.** Do not attempt to lift or move the instrument without the assistance 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 people.

 **CAUTION!** Do not tip the package on end. Tipping may damage the hardware and electronics.

Note: After installation, keep the packaging in case you need to relocate the components.

Register to access the Connect

The Connect cloud-based platform allows access to software applications and secure data storage. When signed to your Connect account, you can export and import of files to and from the Connect cloud-based platform to view performance data from a registered device.

Go to thermofisher.com/cloud, then follow the on-screen instructions to create a new account. .

Related documentation

Document	Publication number	Description
<i>Attune™ NxT Flow Cytometer User Guide</i>	100024235	This document describes how to operate the Attune™ NxT Flow Cytometer to acquire and analyze data.
<i>Attune™ Cytometric Software User Guide</i>	100024236	This document describes in detail the features of the Attune™ Cytometric Software that is used to control the Attune™ NxT System.
<i>Attune™ Flow Cytometry Maintenance and Troubleshooting Guide</i>	100024234	This document details the use, care, and maintenance for the Attune™ NxT Flow Cytometer.



Life Technologies Holdings Pte Ltd | Block 33 | Marsiling Industrial Estate Road 3 | #07-06, Singapore 739256
For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

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

Revision history: Pub. No. 100024428

Revision	Date	Description
C.0	06 July 2021	Add dimensions for the CytKick™ / CytKick™ Max Autosampler, update formatting.
B.0	14 December 2015	Basis for this revision.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.