

Applied Biosystems® 3500/3500xL Genetic Analyzer

Publication Number 4401689 Revision C

This guide contains the information needed to prepare your site for installation of the Applied Biosystems® 3500/3500xL Genetic Analyzer (Cat. no. 4405186, 4405187).

Site preparation workflow 3500 series

A Life Technologies representative will contact you to schedule the installation. When the installation is scheduled:

1. Receive and inspect the system (page 11).
2. Move the crated instrument to the installation site (page 11).
3. Complete the site preparation checklist (page 2).
4. Ensure:
 - The site preparation checklist is complete.
 - The purchase order is complete.

Installation timeline and training

After the instrument (Cat. no. 4405186, 4405187) is uncrated, installation and testing of the Applied Biosystems® 3500/3500xL Genetic Analyzer takes about 1.5 business days.

During and/or after installation, the Life Technologies service representative reviews data and provides some basic operator training. For additional training and reference information, see the user documents provided with the instrument.

Site preparation checklist

IMPORTANT! Complete, date, and initial all items in the following checklist before the scheduled installation date. If the site preparation checklist is not complete when the Life Technologies service representative arrives, the scheduled installation may be postponed.

✓	Date	Initials	Site preparation requirement	See page
<input type="checkbox"/>			Customer responsibilities have been reviewed and personnel have been assigned.	3
			The installation site is identified and meets requirements:	
			<input type="checkbox"/> Space and clearance	6
			<input type="checkbox"/> Environmental	6
			<input type="checkbox"/> Electrical	8
			<input type="checkbox"/> Network	9
			<input type="checkbox"/> Safety	9
<input type="checkbox"/>			Antivirus software is available for installation on the computer provided with the instrument.	10
<input type="checkbox"/>			All materials needed for installation and operation are available.	10
<input type="checkbox"/>			The instrument was received and inspected:	
			<input type="checkbox"/> All items on the shipping list are the same items ordered at the time of purchase.	11
			<input type="checkbox"/> Any damage to shipping containers was reported to the shipping company that delivered the instrument.	
			<input type="checkbox"/> Any damage or mishandling was recorded on the shipping documents.	
<input type="checkbox"/>			The installation kit was unpacked and stored as specified.	11
<input type="checkbox"/>			The installation site is cleared and ready for instrument installation.	11
<input type="checkbox"/>			The crated instrument and other shipping containers are moved to the installation site.	

Customer responsibilities

Personnel	Responsibilities
Site preparation/ installation coordinator	<ul style="list-style-type: none"> • Reviews the site preparation guide for safety information and instrument requirements. • Coordinates personnel and tasks. • Chooses the site. • Reviews checklists with applicable personnel, then with the service representative to verify that the site is properly prepared. • Receives and inspects the instrument. • Stores the reagents box according to the specifications indicated in the product inserts. • Schedules the installation and informs personnel of the installation day. • Ensures that the site is clear of unnecessary material on the installation day. • Is available to assist the service representative throughout installation.
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 the customer's facility.
Laboratory personnel/ primary users	<ul style="list-style-type: none"> • Review safety information. • Ensures that all customer-provided materials for installation are present at the site. • Ensures that 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"> • Ensures that the 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 - Safety and installation materials • If possible, moves the crated instrument 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 instrument.

Personnel	Responsibilities
Network or IT specialist (if the instrument 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 instrument to the network. • If applicable, provides and installs a network or dedicated printer. <p> CAUTION! Do not attempt to connect the instrument components to the network before the service representative arrives.</p>

Site requirements

Dimensions and weights

To prepare for installation, provide space for receipt and configuration of the following components. This section provides dimensions and weights for the crates and packages you will receive, and it describes the dimensions of the instrument after it has been installed and configured.

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

Crates and packages

Ensure the building clearances allow for the passage of the instrument crates and packages.

Crate or package	Height	Length (depth)	Width	Weight
Instrument	114.8 cm (45.2 in.)	75.5 cm (29.7 in.)	79 cm (31.1 in.)	95 kg (210 lb)
Computer	71.1 cm (28 in.)	72 cm (28.5 in.)	43.2 cm (17 in.)	31.7 kg (70 lb)
Monitor	21.6 cm (8.5 in.)	44.4 cm (17.5 in.)	38.1 cm (15 in.)	6.35 kg (14 lb)

Components

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

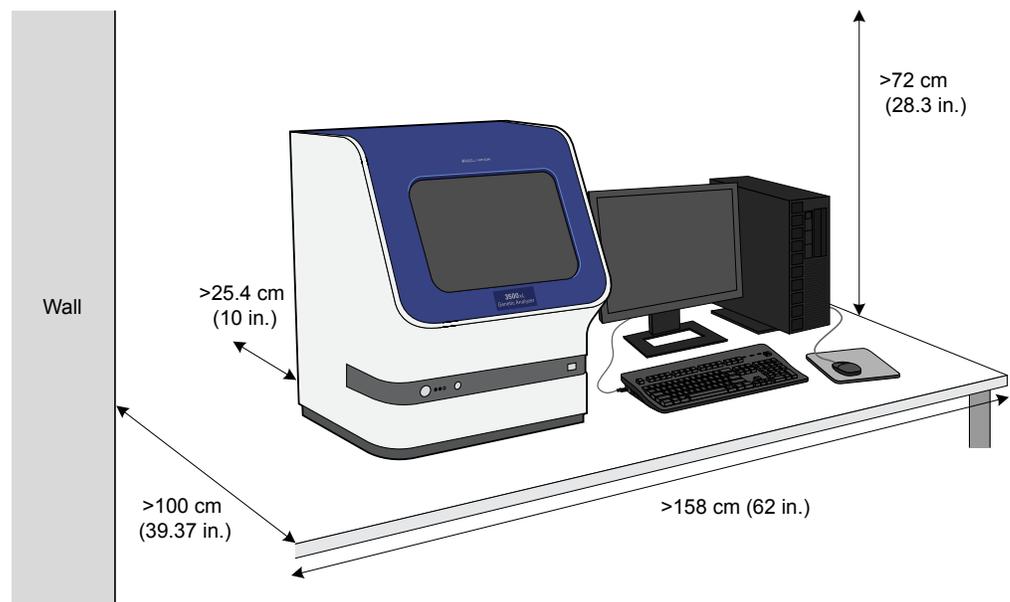
Component	Height	Length (depth)	Width	Weight
Instrument (Closed Door) (Open Door)	72 cm (28.3 in.)	61 cm (24 in.)	61 cm (24 in.) 122 cm (48 in.)	~82 kg (180 lbs)
Computer	34.91 cm (13.75 in.)	39.65 cm (15.61 in.)	10.93 cm (4.3 in.)	~9.4 kg (20.6 lbs)
Monitor	45.72 cm (18 in.)	18.29 cm (7.2 in.)	54.86 cm (21.6 in.)	~4.8 kg (10.4 lbs)
Keyboard	5 cm (2 in.)	15.25 cm (6 in.)	44.7 cm (17.5 in.)	.09 kg (0.2 lbs)

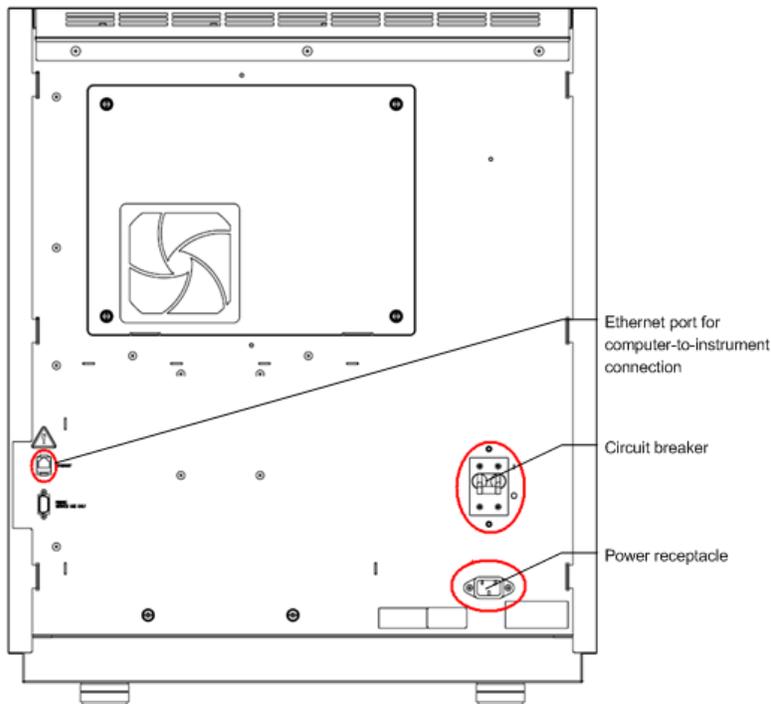


WARNING! PHYSICAL INJURY HAZARD.

Do not attempt to lift or move the instrument without professional assistance. The crated instrument is heavy. Any incorrect lifting or moving of the crated instrument can cause serious injury.

Configured system dimensions





Instrument clearances

During instrument setup and maintenance, it is necessary to access the back and sides of the instrument. If the back of the instrument faces a wall, it will be necessary to have enough space to rotate the instrument on the bench for access.

IMPORTANT! For safety, the power outlet used for powering the instrument must be accessible at all times.

Bench	Minimum clearance
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	>158 cm (62 in.) for the instrument, computer, and computer monitor.

Environmental requirements

Table 1 Environmental requirements

Condition	Requirement
Installation site	Indoor use only
Altitude	Safety tested up to 2000 m

Condition	Requirement
Electrical ratings	Power cord with ground pin required <ul style="list-style-type: none"> • Instrument - AC 100-240 V \pm10%, 50/60 Hz, 3.1 A, power rated 320 VA • Maximum current - 15 A • Maximum power dissipation - 417 VA, 371 W (approximately, not including computer and monitor)
	<ul style="list-style-type: none"> • Computer - AC 100-240 V \pm10%, 50/60 Hz , 2.1 A, power rated 125 VA
	<ul style="list-style-type: none"> • Monitor - AC 100-240 V \pm10%, 50/60 Hz , 1.5 A, power rated 65 VA
Mains AC line voltage tolerances	Up to \pm 10 percent of nominal voltage
Transient category	Installation categories II
Pollution degree	2
Operating conditions	15 to 30°C (59 to 86°F) (Room temperature should not fluctuate \pm 2°C during an instrument run) 20 to 80% relative humidity, noncondensing
Transport and storage conditions	– 30 to +60°C (– 22 to +140°F) Minimum 20% relative humidity, maximum 85% (non-condensing)

Electrical requirements

 **CAUTION!** Do not unpack or plug in any components until the Field Service Engineers (FSEs) have configured the system for the proper operating voltage.

 **WARNING!** For safety, the power outlet used for powering the instrument must be accessible at all times. See "Components" on page xx 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.

- Electric receptacle required: 2-prong with ground pin
- Maximum power dissipation: 417 VA, 371 W (approximately, not including computer and monitor)
- Mains AC line voltage tolerances must be up to ± 10 percent of nominal voltage

Device	Rated voltage	Circuit required	Rated frequency	Rated current	Rated power
Instrument	100–240 $\pm 10\%$ VAC ^[1]	10 A	50/60 Hz	3.1 Z A	320 VA
Computer	100–240 $\pm 10\%$ VAC	10 A	50/60 Hz	2.1 A	125 VA
Monitor				1.5 A	65 VA

^[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.

Electrical protective devices

We recommend several protective devices to protect the system in environments with large voltage and power fluctuations.

Device	Description
Power line regulator	<p>We recommend the use of a 1.5-kVA power line regulator 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> <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>

Device	Description
Uninterruptible power supply (UPS)	<p>We recommend the use of a 1.5-kVA uninterruptible power supply (UPS), especially 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.</p> <p> WARNING! PHYSICAL INJURY HAZARD. Do not attempt to lift the UPS unit without assistance of at least two people. Improper lifting can cause painful and permanent back injury. Refer to the UPS manufacturer user guide for more information.</p> <p>IMPORTANT! UPSs provide power for a limited time. They are meant to delay the effects of a power outage, not to serve as replacement power sources. 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.</p>
Surge protector	<p>We recommend the use of a 10-kVA surge protector (line conditioner) 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> <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>

Network requirements

The instrument is factory-configured for IPv4 TCP/IP communication and includes a fast Ethernet adapter (10/100 Mbps) with a RJ45-type connector and one 3 m (9.8 f) crossover Ethernet cable that connects the computer and the instrument.

If the instrument will be connected to a LAN, an active, tested network jack must be in place before the scheduled installation date. Due to differences in network connections, the Life Technologies service representative cannot configure the system to access a specific network.

You must supply a standard Category 5 Ethernet cable of the required length to connect the computer to your LAN.

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)

Antivirus software requirements

The computer provided with the instrument does not include antivirus software because customer preferences and network requirements vary.

We recommend Norton Antivirus, which has been tested and approved for use with the Applied Biosystems® 3500/3500xL Genetic Analyzer with 3500 Series Data Collection Software 3.

Materials for installation and operation

Installation materials

Have the following materials on hand before installation and operation of the instrument.

- Safety glasses, lab coats, and chemical-resistant, disposable gloves (powder-free)
- Glassware washing solution
- Lint-free tissues
- Mobile bench to allow access to the instrument for maintenance and service
- Easily accessible specified power outlet
- External network connection

- Mini vortexer, centrifuge, and sample tubes
- Freezer (–20°C)
- Optional electrical protective devices (universal power supply unit, surge protector, and/or power line regulator)
- Refrigerator or cold-room (4°C)
- Methanol or isopropanol, HPLC-grade or better
- Water
- Three sizes of micropipettors and tips:
 - 1- to 10- μ L
 - 10- to 100- μ L
 - 100- to 1000- μ L

Operation materials

Additional supplies and consumables are necessary for routine operation of the instrument. Contact a sales representative to order these additional supplies. Use only supplies as specified by Life Technologies.

Receive and inspect 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 and report any damage to the service representative. Record any damage or mishandling on the shipping documents.
3. Immediately unpack the reagents or installation kit box (boxed separately from the instrument components). Store the components as specified in “Installation kit contents” on page 13.

IMPORTANT! Do not unpack Applied Biosystems® 3500/3500xL Genetic Analyzer shipping containers, except for the reagents or installation kit box, to protect yourself from liability if any damage occurred during shipping.

Move the crated instrument to the installation site

1. Clear the installation site of all unnecessary materials.
2. If possible, move the crated instrument and other shipping containers to the installation site. Do not uncrate.



CAUTION! Do not tip the crated instrument on end. Tipping may damage the instrument hardware and electronics.

Note: After installation, retain the crate and instrument packaging in case you need to relocate the instrument.

Related documentation

Document	Publication number	Description
<i>3500/3500xL Genetic Analyzer with Data Collection Software 3 Quick Reference</i>	100026299	Provides abbreviated instructions for using the instrument and software.
<i>3500/3500xL Genetic Analyzer with 3500 Series Data Collection Software 3 User Guide</i>	100025036	Provides information needed to prepare your site for instrument installation.
<i>Polymer Delivery Pump Cleaning Kit Instructions</i>	4414004	Describes how to clean the pump.

Customer and technical support

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

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
- 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 SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

Installation kit contents

Sequencing chemistry installation kit

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4397722)	3500xL Kit (Cat. no. 4405774)
4393708	POP-7™ 3.5mL (384)	2	2 to 8	✓	✓
4404312	BigDye® Terminator (BDT) v3.1 Sequencing Standards, 3500/3500xL	1	2 to 8	✓	✓
4337035	Protocol, BigDye® Terminator v3.1 Cycle Sequencing Ready Reaction (RR-24) This is a protocol used by Field Application Specialists for customer training.	1	—	✓	✓
4404685	Capillary array 8-cap 50cm	1	15 to 30	✓	—
4404689	Capillary array 24-cap 50cm	1	15 to 30	—	✓
4401457	Hi-Di™ 5mL	2	-15 to -25	✓	✓
4393927	Anode buffer container (1x) (4pk)	1	2 to 8	✓	✓
4408256	Cathode buffer container (4pk)	1	2 to 8	✓	✓
4306737	96 TCII well plate w barcode (20pk)	1	—	✓	✓
4410228	96 well retainer and base std (4pk)	1	—	✓	✓
4412614	96 well septa (20pk)	1	—	✓	✓
4410715	Septa buffer/water/waste 3500 (20pk)	1	—	✓	✓
AM9930	Nuclease-free water 500ml (Buf Kit)	1	—	✓	✓
4393718	Conditioning reagent	3	2 to 8	✓	✓
4412619	Pouch cap for polymer pouch (4pk)	1	—	✓	✓

Fragment analysis chemistry installation kit

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4397785)	3500xL Kit (Cat. no. 4405776)
4393708	POP-7™ 3.5mL (384)	2	2 to 8	✓	✓
4345833	Multi-Capillary DS-33 (Dye Set G5) Matrix Std Kit ^[1] This is a sample kit used by Field Application Specialists for customer training.	1	2 to 8	✓	✓
4376911	Fragment Analysis Installation Standard (5-Dye) DS-33	1	2 to 8	✓	✓
4404685	Capillary array 8-cap 50cm	1	15 to 30	✓	—
4404689	Capillary array 24-cap 50cm	1	15 to 30	—	✓
4401457	Hi-Di™ 5mL	2	-15 to -25	✓	✓
4393927	Anode buffer container (1x) (4pk)	1	2 to 8	✓	✓
4408256	Cathode buffer container (4pk)	1	2 to 8	✓	✓
4306737	96 TCII well plate w barcode (20pk)	1	—	✓	✓
4410228	96 well retainer and base std (4pk)	1	—	✓	✓
4412614	96 well septa (20pk)	1	—	✓	✓
4410715	Septa buffer/water/waste 3500 (20pk)	1	—	✓	✓
AM9930	Nuclease-free water 500ml (Buf Kit)	1	—	✓	✓
4393718	Conditioning reagent	3	2 to 8	✓	✓
4412619	Pouch cap for polymer pouch (4pk)	1	—	✓	✓

^[1] Use DS-33 only to qualify 3500 and 3500xL instruments with GeneMapper® Software.

**Combined
sequencing and
fragment analysis
chemistry
installation kit**

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4405778)	3500xL Kit (Cat. no. 4406933)
4393708	POP-7™ 3.5mL (384)	2	2 to 8	✓	✓
4404312	BigDye® Terminator (BDT) v3.1 Sequencing Standards, 3500/3500xL	1	2 to 8	✓	✓
4337035	Protocol, BigDye® Terminator v3.1 Cycle Sequencing Ready Reaction (RR-24) This is a protocol used by Field Application Specialists for customer training.	1	—	✓	✓
4345833	Multi-Capillary DS-33 (Dye Set G5) Matrix Std Kit ^[1] This is a sample kit used by Field Application Specialists for customer training.	1	2 to 8	✓	✓
4376911	DS-33 GeneScan™ 600 LIZ® Installation	1	-15 to -25	✓	✓
4404685	Capillary array 8-cap 50cm	1	15 to 30	✓	—
4404689	Capillary array 24-cap 50cm	1	15 to 30	—	✓
4401457	Hi-Di™ 5mL	2	-15 to -25	✓	✓
4393927	Anode buffer container (1x) (4pk)	1	2 to 8	✓	✓
4408256	Cathode buffer container (4pk)	1	2 to 8	✓	✓
4306737	96 TCII well plate w barcode (20pk)	1	—	✓	✓
4410228	96 well retainer and base std (4pk)	1	—	✓	✓
4412614	96 well septa (20pk)	1	—	✓	✓
4410715	Septa buffer/water/waste 3500 (20pk)	1	—	✓	✓
AM9930	Nuclease-free water 500ml (Buf Kit)	1	—	✓	✓
4393718	Conditioning reagent	3	2 to 8	✓	✓
4412619	Pouch cap for polymer pouch (4pk)	1	—	✓	✓

^[1] Use DS-33 only to qualify 3500 and 3500xL instruments with GeneMapper® Software.

HID chemistry installation kit

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4397722)	3500xL Kit (Cat. no. 4405774)
4393715	POP-4 [®] 3.5mL (384)	2	2 to 8	✓	✓
4345833	Multi-Capillary DS-33 (Dye Set G5) Matrix Std Kit ^[1] This is a sample kit used by Field Application Specialists for customer training.	1	2 to 8	✓	✓
4396597	HID 5-Dye Installation Standard	1	2 to 8	✓	✓
4408399	GeneScan™ 600 LIZ [®] Size Standard v2.0	1	2 to 8	✓	✓
4404683	Capillary array 8-cap 36cm	1	15 to 30	✓	—
4404687	Capillary array 24-cap 36cm	1	15 to 30	—	✓
4401457	Hi-Di™ 5mL	2	-15 to -25	✓	✓
4393927	Anode buffer container (1x) (4pk)	1	2 to 8	✓	✓
4408256	Cathode buffer container (4pk)	1	2 to 8	✓	✓
4306737	96 TCII well plate w barcode (20pk)	1	—	✓	✓
4410228	96 well retainer and base std (4pk)	1	—	✓	✓
4412614	96 well septa (20pk)	1	—	✓	✓
4410715	Septa buffer/water/waste 3500 (20pk)	1	—	✓	✓
AM9930	Nuclease-free water 500ml (Buf Kit)	1	—	✓	✓
4393718	Conditioning reagent	3	2 to 8	✓	✓
4412619	Pouch cap for polymer pouch (4pk)	1	—	✓	✓

^[1] Use DS-33 only to qualify 3500 and 3500xL instruments with GeneMapper[®] Software.

HID and sequencing chemistry installation kit

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4397722)	3500xL Kit (Cat. no. 4405774)
4393708	POP-7™ 3.5mL (384)	2	2 to 8	✓	✓
4393715	POP-4® 3.5mL (384)	2	2 to 8	✓	✓
4404312	BigDye® Terminator (BDT) v3.1 Sequencing Standards, 3500/3500xL	1	2 to 8	✓	✓
4337035	Protocol, BigDye® Terminator v3.1 Cycle Sequencing Ready Reaction (RR-24) This is a protocol used by Field Application Specialists for customer training.	1	—	✓	✓
4345833	Multi-Capillary DS-33 (Dye Set G5) Matrix Std Kit ^[1] This is a sample kit used by Field Application Specialists for customer training.	1	2 to 8	✓	✓
4396597	HID 5-Dye Installation Standard	1	2 to 8	✓	✓
4408399	GeneScan™ 600 LIZ® Size Standard v2.0	1	2 to 8	✓	✓
4404685	Capillary array 8-cap 50cm	1	15 to 30	✓	—
4404689	Capillary array 24-cap 50cm	1	15 to 30	—	✓
4404683	Capillary array 8-cap 36cm	1	15 to 30	✓	—
4404687	Capillary array 24-cap 36cm	1	15 to 30	—	✓
4401457	Hi-Di™ 5mL	2	-15 to -25	✓	✓
4393927	Anode buffer container (1x) (4pk)	1	2 to 8	✓	✓
4408256	Cathode buffer container (4pk)	1	2 to 8	✓	✓
4306737	96 TCII well plate w barcode (20pk)	1	—	✓	✓
4410228	96 well retainer and base std (4pk)	1	—	✓	✓
4412614	96 well septa (20pk)	1	—	✓	✓
4410715	Septa buffer/water/waste 3500 (20pk)	1	—	✓	✓
AM9930	Nuclease-free water 500ml (Buf Kit)	1	—	✓	✓
4393718	Conditioning reagent	3	2 to 8	✓	✓
4412619	Pouch cap for polymer pouch (4pk)	1	—	✓	✓

^[1] Use DS-33 only to qualify 3500 and 3500xL instruments with GeneMapper® Software.

QST chemistry installation kit

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4407051)	3500xL Kit (Cat. no. 4413692)
4393717	POP-6™ 3.5mL (384)	2	2 to 8	✓	✓
4336824	BigDye® Terminator (BDT) v1.1 Matrix Standards Kit, 31xx and 3500	1	2 to 8	✓	✓
4404314	BigDye® Terminator v1.1 Sequencing Install Standard	1	-15 to -25	✓	✓
4404685	Capillary array 8-cap 50cm	1	15 to 30	✓	—
4404689	Capillary array 24-cap 50cm	1	15 to 30	—	✓
4401457	Hi-Di™ 5mL	2	-15 to -25	✓	✓
4393927	Anode buffer container (1x) (4pk)	1	2 to 8	✓	✓
4408256	Cathode buffer container (4pk)	1	2 to 8	✓	✓
4306737	96 TCII well plate w barcode (20pk)	1	—	✓	✓
4410228	96 well retainer and base std (4pk)	1	—	✓	✓
4412614	96 well septa (20pk)	1	—	✓	✓
4410715	Septa buffer/water/waste 3500 (20pk)	1	—	✓	✓
AM9930	Nuclease-free water 500ml (Buf Kit)	1	—	✓	✓
4393718	Conditioning reagent	3	2 to 8	✓	✓
4412619	Pouch cap for polymer pouch (4pk)	1	—	✓	✓

QST and fragment chemistry installation kit

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4413691)	3500xL Kit (Cat. no. 4413692)
4393708	POP-7™ 3.5mL (384)	2	2 to 8	✓	✓
4393717	POP-6™ 3.5mL (384)	2	2 to 8	✓	✓
4345833	Multi-Capillary DS-33 (Dye Set G5) Matrix Std Kit ^[1] This is a sample kit used by Field Application Specialists for customer training.	1	2 to 8	✓	✓
4376911	DS-33 GeneScan™ 600 LIZ® Installation	1	-15 to -25	✓	✓

Part Number	Name	Qty	Storage Temp. (°C)	3500 Kit (Cat. no. 4413691)	3500xL Kit (Cat. no. 4413692)
4404685	Capillary array 8-cap 50cm	1	15 to 30	✓	—
4404689	Capillary array 24-cap 50cm	1	15 to 30	—	✓
4401457	Hi-Di™ 5mL	2	-15 to -25	✓	✓
4393927	Anode buffer container (1x) (4pk)	1	2 to 8	✓	✓
4408256	Cathode buffer container (4pk)	1	2 to 8	✓	✓
4306737	96 TCII well plate w barcode (20pk)	1	—	✓	✓
4410228	96 well retainer and base std (4pk)	1	—	✓	✓
4412614	96 well septa (20pk)	1	—	✓	✓
4410715	Septa buffer/water/waste 3500 (20pk)	1	—	✓	✓
AM9930	Nuclease-free water 500ml (Buf Kit)	1	—	✓	✓
4393718	Conditioning reagent	3	2 to 8	✓	✓
4412619	Pouch cap for polymer pouch (4pk)	1	—	✓	✓

^[1] Use DS-33 only to qualify 3500 and 3500xL instruments with GeneMapper® Software.

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

DISCLAIMER

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, 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

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2015 Thermo Fisher Scientific Inc. All rights reserved.

For support visit lifetechnologies.com/support or email techsupport@lifetech.com
lifetechnologies.com

18 February 2015

