

FL 6500 Pulse Fluorescence Spectrometer and FL 8500 Continuous Wave Fluorescence Spectrometer



Fluorescence Spectroscopy

Preparation Checklist

- Order Overview
- Site Requirements
- Electrical Requirements
- Environmental Requirements
- PC Configuration
- Installation Overview

Order Overview

Upon receipt of the system, please review the order. Record any discrepancies between the PerkinElmer order and your Purchase Order, along with any agreements or commitments made by your PerkinElmer Sales Representative that are NOT listed on the order. Let your Customer Care Representative know about any discrepancies and/or commitments when you submit your Site Readiness confirmation.

Site Requirements

The FL 6500 and FL 8500 systems must be installed on a flat and level bench. The bench must be strong enough to support the weight specified below in addition to any ancillary equipment without warping, wobbling or swaying. Make sure that there are no overhanging shelves, and no water pipes or faucets that could leak onto the system.

Bench Space/Instrument Weights

Table 1. Bench Space/Instrument Weights.

Dimensions*				
Instrument	Width	Depth	Height	Weight
FL 6500	66 cm	64 cm	36 cm	47 kg
FL 8500	66 cm	64 cm	36 cm	47 kg

*The dimensions and weights listed above do not include the PC/Monitor.

The workbench where the FL 6500 and FL 8500 systems and PC are to be located should be at a suitable height to allow the user to work comfortably and must be strong and vibration free.

The bench space required should be slightly larger than the above dimensions to allow the user to work safely and comfortably when analyzing samples.

Specifically, there must be at least 13.5 in or 34.3 cm clear room above the instrument to allow the sample compartment lid to open.

Note: Floor vibrations or noise from heavy manufacturing equipment can affect the performance of the system. Avoid proximity to intense magnetic fields.

Electrical Requirements

Power Consumption

Max power consumption is 65 VA (approximately 40 W).

Maximum power consumption is experienced when the FL 6500 and FL 8500 systems are powering the (optional) wireless router.

In power save mode the mains power consumption is <10 VA (approximately 6 W) and the DC power consumption is 0.2 W.

Power Specifications

Any fluctuation in the line supply must not exceed $\pm 10\%$ of the nominal voltage.

The FL 6500 and FL 8500 Systems: 100 V to 230 V, 50 Hz or 60 Hz

Power Outlets:

Three (one for the FL 6500 and FL 8500, two for the PC/Monitor)

- If the electricity supply does not conform to the above specifications, please consult PerkinElmer prior to installation. If the supply of power is erratic please use an uninterruptible power supply (UPS). Incorrect shutdown, power fluctuations or brown-outs may damage the spectrometer or system.
- If possible, do not use photocopiers, discharge lamps, radio transmitters or other equipment with large or frequent transient loads on the same supply circuit as this may affect the performance of the system.
- Make sure that the power outlets at the electrical supply sockets to the system are not obstructed.

Environmental Requirements

Laboratory Environment

Operating Temperature Range: 0 °C to 50 °C*

Storage Temperature Range: -20 °C to 60 °C

Maximum Relative Humidity: 90% non-condensing

**If you intend to operate the instrument regularly at temperatures of 40 °C or higher, for optimum performance, it is recommended that you repeat the Wavelength Calibration at the appropriate operating temperature. Refer to the on-screen software help for more information.*

- The laboratory should be relatively free of dust, corrosive fumes and vibrations.
- The laboratory bench should be free from vibration and, if possible, isolated from other equipment that may induce vibration.
- Do not leave the instruments in direct sunlight.

- Make sure that there are no overhanging shelves, and no water pipes or faucets that could leak onto the instrument.
- Ensure there is sufficient clearance underneath the instrument to allow an adequate flow of cooling air.
- Major changes in temperature should be avoided – if the instrument is moved to a different environment which has a temperature difference, it should be left to stabilize before data collection can begin.

Safety Requirements

Electrical Safety

Connect the instrument power supply to a power outlet that includes a switch or other means of disconnection from the electricity supply.

Only plug the instrument power supply into a power outlet that is provided with a protective earth connection.

Do not operate the instrument with any covers or parts removed.

Servicing should be carried out only by a PerkinElmer service representative or similarly authorized and trained person.

Disconnect the instrument from all voltage sources before opening it for any adjustment, replacement, maintenance, or repair. If afterwards, the opened instrument must be operated for further adjustment, maintenance, or repair, this must only be done by a skilled person who is aware of the hazard involved.

Use only fuses with the required current rating and of the specified type for replacement. Do not use makeshift fuses or short-circuit the fuse holders.

Whenever it is likely that the instrument is no longer electrically safe for use, make the instrument inoperative and secure it against any unauthorized or unintentional operation.

The instrument is likely to be electrically unsafe when it:

- Shows visible damage.
- Fails to perform the intended measurement.
- Has been subjected to prolonged storage under unfavorable conditions.

PC Configuration

A pre-configured PC, which meets the system requirements, can be provided as part of the FL 6500 or FL8500 order, in which case, the system can be installed as normal.

If a customer-supplied PC is to be used, it must meet the following minimum specifications and hardware requirements:

Component	Requirement
Operating System	Windows® 10 (64-bit)
Processor	Intel Core i5-7500
Memory	8 GB (1x8G) 2400 MHz DDR4
HDD Size / HDD Spindle	1 TB 7.2k RPM SATA 6 Gbps
WWAN	Intel DB WLAN 8265 Software

Any extra time taken for the installation that is caused by issues encountered from trying to install the system on a PC which does not meet these requirements (or is otherwise unsuitable) will be billable at the current service rate.

PerkinElmer may not provide maintenance service on customer-supplied items.

Software Requirements

The FL 6500 and FL8500 instruments are operated with the Spectrum FL™ software.

NOTE: *It is important to note that you must be logged on at Administrator level before installing the software.*

The latest Microsoft Service Packs can be downloaded from: www.microsoft.com/msdownload/default.htm.

Installation Overview

The customer is recommended to remove the FL 6500 and FL 8500 instruments from the shipping box. The instrument should be placed on the bench along with associated components.

Customer to Hold Packing Material for CSE

When PerkinElmer is installing the instrument, the packing material must be retained until the installation is fully signed off (recommend to retain for the warranty period).

Physical Installation (Instrument Only)

The FL 6500 and FL 8500 instruments are small enough to be handled by a single person using the two handholds underneath on the ends of the instrument.

Physical Installation (Accessories)

As required.

Installation Test Standards

When PerkinElmer installs the instrument, the Service Engineer will test the instrument in order to ensure that its performance meets PerkinElmer's installation specifications.

Customer Orientation

When PerkinElmer installs the instrument, the service engineer will familiarize you with the basic instrument and software operation. If further training is required, please contact your local Sales representative.

Ensure that a staff member, who will have ultimate responsibility for the FL 6500 and FL 8500 System, is available during the installation.

Related Documents

The FL 6500/8500 instrument User's Guide informs you how to use the Fluorescence instrument. On-screen HTML Help Systems are supplied with all software applications.

Manuals for the PC and for the operating system of the PC are supplied with the PC (if purchased from PerkinElmer).