

# FLUORESCENCE THAT OUTSHINES EXPECTATIONS

## Introducing the FL 6500™ and the FL 8500™ Fluorescence Spectrometers

We know that high performance and reliability are top priorities for labs everywhere. That's why we've developed a complete fluorescence solution that enables confident and precise sample testing. Whether you're looking for a high-energy pulsed Xenon light source that will preserve the integrity of your sample (FL 6500), or a high-sensitivity source for testing diluted or small samples (FL 8500), our revolutionary instruments give you the speed and accuracy you need right at your fingertips.

### FL 6500 Pulse Fluorescence Spectrometer



- Built for bioresearch, including cell-biology, molecular biology, immunology, enzymology, and protein analysis
- Unique adjustable power settings are ideal for phosphorescence and avoiding photobleaching

### FL 8500 Continuous Wave Fluorescence Spectrometer



- Made for material characterization; industrial tracing and manufacturing R&D; agricultural and environmental analysis; and analysis of LEDs, solar cells, and organic electroluminescent materials
- Fast scanning of 60,000 nm/min minimizes scan time

## Software Features and Functionality

- Spectra Scan
- 3D Spectra Scan
- Quantification
- Scanning Anisotropy/Polarization
- Single Read
- Anisotropy/Polarization
- Time Drive
- Kinetics
- Intracellular Ion Concentrations
- Quantum Yield (Relative and Absolute)
- Quenching
- Wavelength Program
- Lifetime
- Absorbance
- Validation
- Sipper Control
- Service Utility
- Sample Table
- Integrating Sphere for Absolute Quantum Yield and Total Fluorescence
- Pre-Scan Mode
- Synchronous Scan Mode (Constant Energy and Wavelength)
- 3D Synchronous Scan
- Read FLWinLab Data Files
- Read UVWinLab Data Files

## Interchangeable Accessories

Our broad accessories are specifically designed for the FL 6500 and FL8500 and versatile enough to handle virtually any type of sample.

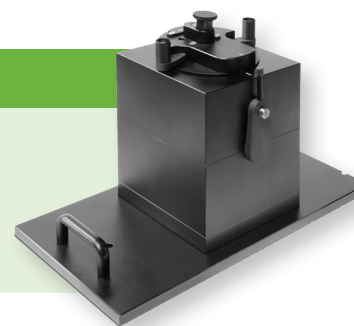
Accessories	Part Number
Single Cell Holder	N4201010
Single Cell Water-Jacket Holder	N4201011
Single Cell Water-Jacket Holder with Stirrer	N4201012
Micro Cell holder	N4201015
Micro Cell Water-Jacket Holder with Stirrer	N4201016
4 Position Multi-Cell holder	N4201025
4 Position Multi-Cell Water-Jacket Holder	N4201026
4 Position Multi-Cell Water-Jacket Holder with Stirrer	N4201027
Single Cell Peltier Holder	N4201029
4 Position Multi-Cell Peltier Holder	N4201030
Solid Sample Holder (for powders also require N4201032)	N4201013
Variable Angle Solid Sample Holder (for powders also require N4201032)	N4201014
Precision Cell for Powder Sample (used with N4201013 & N4201014)	N4201032

Accessories	Part Number
Vis Automated Polarizer	N4201022
UV-Vis Automated Polarizer	N4201019
Manual Polarizer Holder	N4201023
Rapid Mixing Accessory	N4104014
Rapid Mixing Accessory with Pneumatic Drive	N4104015
Rapid Mixing Accessory Cover (only needed with N4104015)	N4201020
Absorbance Module	N4201018
Fiber Optic Probe	N4201021
Microplate Reader	N4201024
Fast Filter (excluding filters)	N4201028
Auto Sipper (including Front Plate & FL Flow Cell)	N4201031
Auto Sampler S10	N2020004
Integrating Sphere	N4201017

## FEATURED ACCESSORY

### Integrating Sphere

Measure absolute quantum yield for liquid and solid samples. Achieve flexibility even with low absorbency and low quantum yield samples.



PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright © 2018, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.