

Liquid Chromatography

Key Features and Benefits

- Install/change flow cell in one easy motion
- Flow cell design that sustains pressures up to 1500 psi
- Highest data acquisition rates of any PDA on the market
- Liquid core waveguide for optimal energy
- Best wavelength accuracy gives increased confidence in results
- Designed for minimal stray light to provide the widest linear dynamic range

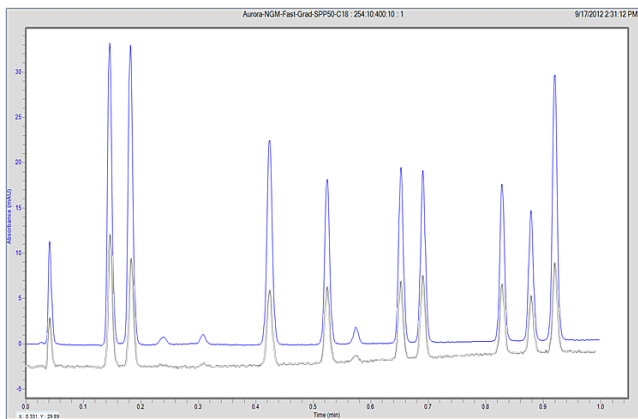


PDA Technology That Delivers New Levels of Performance and Productivity

With more stringent industry regulations and greater throughput demands, today's analytical laboratories are turning to PerkinElmer® for high performance liquid chromatography solutions. The newest addition to the LC family is the Flexar™ PDA Plus™ detector. This innovative technology brings the flexibility of modular design, the power of sensitivity and accuracy and the fast sampling speeds to the most optically advanced PDA detector on the market.

PerkinElmer's Flexar™ PDA Plus™ detector offers a unique design that includes two modular flow cells. The standard path length cell of 10 mm gives outstanding sensitivity, yet when very low levels of detection are needed, the 50 mm flow cell provides a five-fold increase in response levels.





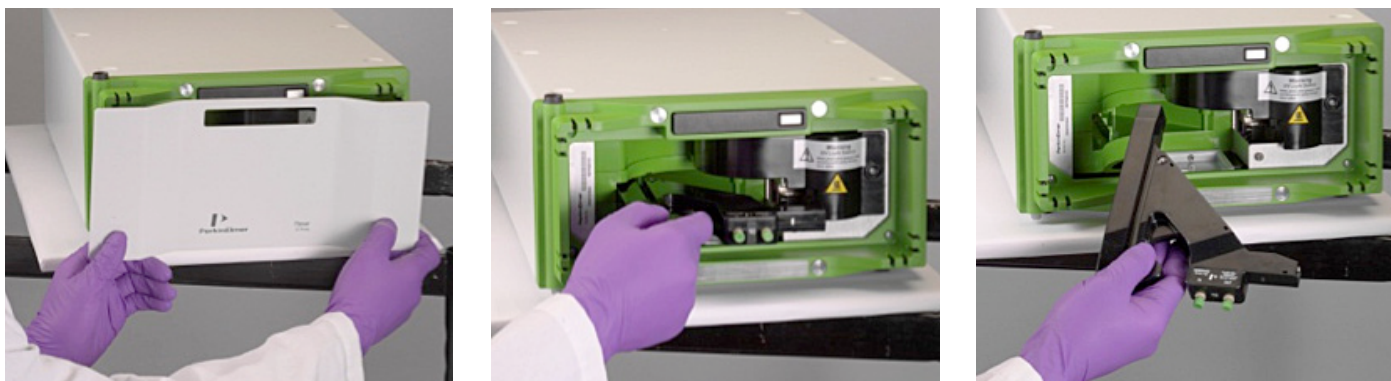
■ Flexar PDA Plus
■ Flexar EP

Overlay of Flexar EP w/67 mm 2.4 uL cell v. PDA Plus w/10 mm (1 uL) cell.

Easy to Install Flow cells for Even the Most Complex Analyses

The Flexar PDA Plus Detector is ideal for laboratories with high sample throughput requirements with multiple, highly sensitive methods. Its interchangeable modular flow cell design allows for one-motion installation and exchange, enabling greater flexibility and ultimately higher productivity in your lab.

Remove/replace flow cell in one simple movement. No tools or adjustments required.

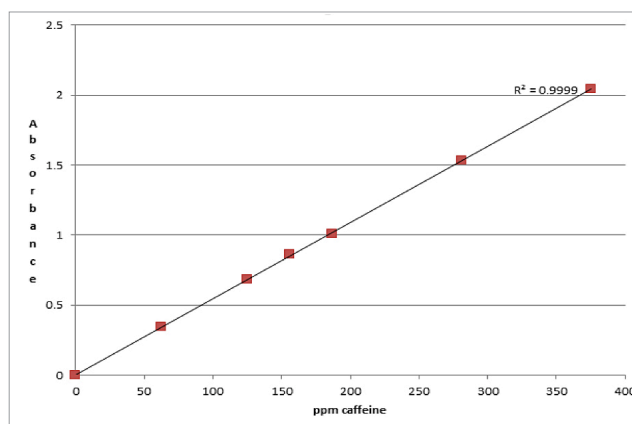


Unsurpassed Sensitivity and Flexibility

Additionally, the Flexar PDA Plus detector's unsurpassed linear dynamic range (up to 2.0 AU with less than 3% deviation) is the best of any PDA detector on the market and rivals or exceeds that of conventional UV detectors.

Low baseline noise coupled with the high linear dynamic range of the detector gives the user the ultimate in analyte concentration tolerance. This allows the user to detect from very low concentrations (those previously lost in the noise) to high concentrations with no need to dilute samples.

The combination of sampling rates up to 200 Hz, a flow cell capable of sustaining pressures up to 1500 psi with low optical dispersion, makes the Flexar PDA Plus the perfect system for any application.



Flexar PDA Plus calibration curve for caffeine indicates linearity over a wide dynamic range.

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



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright © 2013, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.