

INTRODUCING ENLIGHTENED UV/VIS



LAMBDA® 365+ UV/Vis System

The Perfect Balance of Versatility, Simplicity, and *Value*

QA/QC analysts and lab technicians in a wide variety of fields – pharma and biopharma, applied materials, optics, food, and academia among them – need to analyze samples quickly, accurately, and easily through their UV/Vis systems, with better turnaround times, confidence in results, and increased productivity. That's what drives our **LAMBDA® 365+ system**.

What's more, with the global workforce shifting to a hybrid work environment, the **LAMBDA 365+ system** can help labs securely coordinate and collaborate remotely across the globe, intuitively, with minimal training. All while offering real value for the long haul.

Performance at the Highest Level

Designed to the highest standards for speed and performance, the **LAMBDA 365+ UV/Vis** delivers variable spectral bandwidth capability from 0.5 nm to 20 nm, with a wavelength range of 190-1100 nm, and its innovative dual-beam optics provide high-quality data with low noise and stray light – so you benefit from accurate, repeatable results on a host of applications. The system enables:

- Measurement of higher absorbing liquids (up to 4A)
- Faster scan speeds to unravel chemical kinetics
- Higher wavelength reproducibility, photometric repeatability, and low stray light
- Built-in methods for fast, easy analysis



APPLICATIONS

Versatility for Any and All Applications

Pharmaceuticals, analytic chemists, geneticists, and manufacturing QA/QC analysts alike need quick, accurate analysis of their samples, with better turnaround times, confidence in results, and increased productivity. That's why versatility drives our **LAMBDA 365+ system**. It's UV/Vis that provides:

- A variety of sampling accessories with large sample compartments – and sampling accessories snap right in and self-calibrate, so you're productive right away
- Flexibility to choose between maximum compliance or productivity with an optional touchscreen
- Fast analysis and reproducible results with built-in workflows to analyze a wide variety of samples

The System That Drives Out Complexity

With a constantly changing laboratory landscape, high turnover, and business uncertainty, you need less complexity from your UV/Vis instruments and software. The **LAMBDA 365+** is designed to simplify your workflows and training needs so you can focus on what matters – your results. The system lets you:

- Simplify protocols and maximize productivity with out-of-the-box workflows
- Reduce setup time and effort with self-aligning plug-and-play accessories to minimize errors
- Maximize value with less maintenance and longer instrument life
- Collaborate effortlessly using touchscreen client-server software

APPLICATIONS



The Right Accessories for All Your Sampling Needs

The modular, flexible **LAMBDA 365+ system** accommodates a wide range of accessories that fit the type of samples your analysis requires. Most accessories snap right in – no tools necessary – and calibrate automatically, so everyone in your lab can be productive right from the start.

With the **LAMBDA 365+**, you get:

- Precision-engineered autosamplers to meet all your throughput needs
- Wider sample compartments to accommodate more sample types
- Plug-and-play accessories for maximum productivity

APPLICATIONS

Fixed-Angle Reflectance Holder

Designed for measuring specular reflectance spectra at a fixed angle



Advanced Transmission Holder

Used for measuring the transmittance of films, glass, plate, and so on, with adjustable fixing arms for samples of varying thicknesses



Autosipper Accessory

Sipper system with complete software control for automation of liquid sampling, including sample flow



Variable-Pathlength Cuvette Holder

An easy-to-install accessory for rectangular cells of different pathlengths



Water-Jacketed Single-Cell Holder with Stirrer

Software-recognized accessory with internal stirring functions for sample and reference positions



Water-Jacketed Microcell Holder with Stirrer

Software-recognized accessory with internal stirring functions for sample and reference positions, used with a water circulator



Auto-Peltier Controller

Comes with internal liquid-cooling system and N2 purge-gas capability, plus magnetic stirrer control



Cylindrical Cell Holder

Used for measuring the transmittance of cylindrically shaped samples



Faster Results with Simplified Workflows

Innovative, intuitive, easy-to-use **UV WinLab™ software** is designed for simplicity and efficiency in data collection and processing, while generating accurate results. It mirrors the way you work, guiding you through method development, analysis, reporting, and results in simple steps. And with advanced software capabilities such as macros and sample tables, it's easy to personalize all stages of your analysis.



APPLICATIONS

Autosamplers fully integrated with software

- Optimize automated sampling with the autosipper accessory and autosamplers
- Save parameters to your method for streamlined analysis workflow

Simplified data exports

- Data export options in the analysis method for easy exporting of sample data in common formats
- Export your data to Excel, an external server, or LIMS system

Built-in performance-validation methods

- Easily troubleshoot and resolve issues by making use of over 20 instrument performance validation (IPV) tests
- Monitor the performance of the instrument quickly and easily

Easy backup and data retrieval

- Manage your data effortlessly with the UV Winlab database tool to ensure your data is safe, secure, and 21 CFR Part 11 compliant
- Multilevel user permissions, password-protected access control, method-lock facility, and fully configurable e-signature points



S20 autosampler series



APPLICATIONS

FOCUS ON

Chemicals

With the **LAMBDA 365+ UV/Vis instrument**, you can monitor concentration changes of reactants to better understand rapid chemical processes and gain deeper insights into reaction kinetics, at exceptional scan speeds. The **LAMBDA 365+** also helps chemical manufacturers to identify, determine, characterize, and quantify their samples and help build better products that are stronger and better.

For example, sodium hypochlorite is included on the Environmental Protection Agency list of disinfectants recommended for use against COVID-19. The standard method for determination of sodium hypochlorite is an iodometric titration, which requires several chemical reagents and can be time consuming and costly when running many samples. The **LAMBDA 365+** provides a simple, fast solution for measurement of sodium hypochlorite in a QA/QC setting, enabling measurements in as little as 30 seconds using a single reagent.





APPLICATIONS

FOCUS ON

Coolants

Busy high-throughput analytical testing labs specializing in analysis of in-service lubricants and coolants need to quickly, easily, and accurately characterize the composition of additives to prevent equipment failure and costly repairs. Our **LAMBDA 365+ system** enables easy, fast, and automated analysis of nitrites, a common additive to coolant formulations. In-service lubricant and coolant testing labs also benefit from:

- Automation options, including an autosipper and autosamplers
- Touchscreen operation (with LAMBDA 365+ Touch/UVWinlab Touch)
- Built-in methods for fast workflow setup
- Large sample compartment for easy routing of flow cell tubing or manual sample changing
- Low maintenance – and greater confidence in instrument operation





APPLICATIONS

FOCUS ON

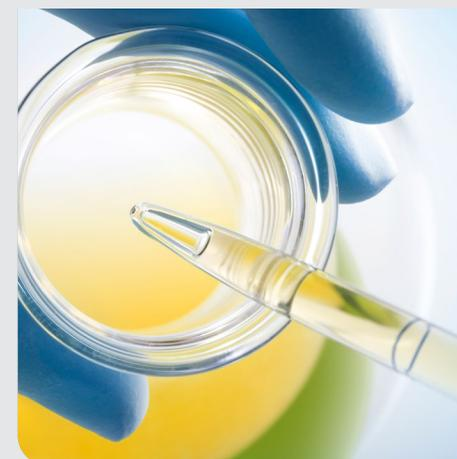
Food

Feed- and food-processing plants need quick, accurate, reliable ways to monitor ingredient quality, from incoming raw materials to finished products. Our **LAMBDA 365+ system** is the flexible, simple analytical solution for identification of key characteristics in foods. For example, it's ideal for analysis of:

- Brix/sugar in juices
- Ethanol, tannins, and sulfites in wines
- Olive oil acidity levels, for classification (olive pomace to extra virgin)
- Palm oil, to measure bleachability

The system is also perfect for more advanced methods that address food authenticity, provenance, and adulteration, including:

- Pomegranate juice adulteration (color changes and absence of antioxidants)
- Distilled spirits adulteration (water, ethanol, or methanol)
- Olive oil purity and degree of oxidation





APPLICATIONS

FOCUS ON

Pharmaceuticals

Pharmaceutical labs need the flexibility to perform a wide range of analyses, from analytical method development for QA/QC to basic research into new drug candidates to dissolution testing. The **LAMBDA 365+ system** shines in all these applications – and it's 21 CFR Part 11 compliant to meet your most stringent government and industry regulations.

Applications include:

- Quantification of DNA and proteins
- DNA melting experiments
- Enzyme kinetics
- USP method compliance
- Solvent analysis
- Biomolecule and biomedical implant characterization





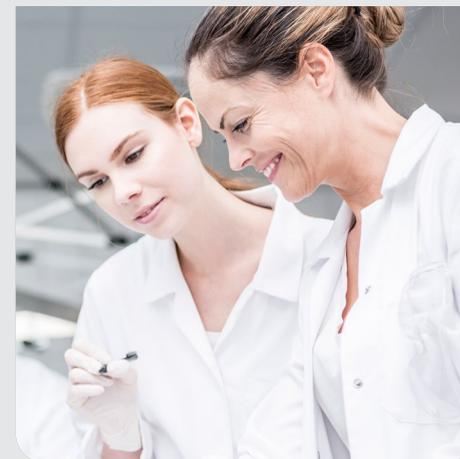
APPLICATIONS

FOCUS ON

Water

Drinking water analysis starts with confidence – and our established workflow solutions ensure you get accurate, timely results, day after day. And with faster turnarounds – and lab technician turnover – adding to your challenges, these simple, effective solutions instill a high level of confidence in your lab and help ensure your communities' water supplies are safe and unadulterated. Our **LAMBDA 365+ UV/Vis** spectrophotometer is the perfect solution for a whole range of water analysis challenges, including:

- Quantitative analysis of chemical oxygen demand (COD)
- Determination of nitrates, orthophosphate, and iron, using the ascorbic acid method for iron, the brucine method for nitrate-nitrogen, and more



Complete Services for Better Productivity and Efficiency



Today's lab leaders face daunting challenges, from tighter deadlines to increased budget scrutiny to teams with various degrees of comfort with lab equipment. Time that could be spent getting ahead is spent on noncore activities. To help you overcome barriers to success, OneSource® Laboratory Services has built a team of trained scientists and engineers who bring their real-life knowledge to you, helping increase your productivity with recommendations on how to best utilize your assets get you back to your core mission.

No matter what size of type of lab you're in, you need to know your equipment will work as expected every time you turn it on. From contracts and performance maintenance for our instruments (as well as other manufacturers' equipment) to full lab asset management delivered globally, we can help you make the most of your important lab assets. And for labs looking to introduce new equipment and techniques, we offer training at our facilities and at yours.

Compliance Services

Navigating the complexities of global regulatory entities and requirements can be difficult for even the best-run laboratories. OneSource Compliance Services can help get you back to focusing on your core activities.

With our consultative approach, we offer gap analysis or inspection readiness and provide comprehensive training, plus robust products and services to make the qualification process run smoothly. You simply review our standard recommended OQ protocols and customize to them to your specifications. Our highly trained service engineers then go to work, following the protocol to test instruments against approved specifications. Your final qualification report is delivered in a concise, easy-to-review layout in paper and secure electronic format. This is compliance made simple.

Education Services

Whether you're looking for a basic instrument refresher course, simple troubleshooting techniques, general application support, or method optimization, our field application scientists or service engineers can come directly to your lab. Through education, you'll gain knowledge and insights into the latest techniques, not only increasing your confidence, but also unlocking the full potential of your instrument.

APPLICATIONS



[For more information visit www.perkinelmer.com/lambda365plus](http://www.perkinelmer.com/lambda365plus)

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 ©2022, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

593359

PKI