

inside the PerkinElmer LAMBDA 950/850/650 Series the highest performing UV/Vis/NIR

DEUTERIUM AND TUNGSTEN HALOGEN LIGHT SOURCES

Prealigned and prefocused for quick replacement and maximum uptime.

DOUBLE HOLOGRAPHIC GRATING MONOCHROMATORS For ultra-low stray light performance.

COMMON BEAM MASK Allows precise adjustment of beam height to match samples of different dimensions.

> COMMON BEAM DEPOLARIZER Corrects for inherent instrument polarization to allow accurate measurements of birefringent samples (optional)

Switches between sample and reference beam. Four-segment design provides individual blank readings for sample and reference, increasing



and UV/Vis available



LARGEST SAMPLE COMPARTMENT IN THE INDUSTRY Allows easy access to a wide variety of sampling accessories and sample types.

HIGH-SENSITIVITY PHOTOMULTIPLIER AND PELTIER-CONTROLLED PbS DETECTORS Provides full range UV/Vis/NIR coverage from 175 to 3,300 nm (LAMBDA 950 only).

Houses a range of snap-in sampling modules including transmission optics (shown) 60 mm and 150 mm integrating spheres and the Universal Reflectance Accessory (URA) for high-precision absolute reflectance measurements.

## Another PerkinElmer exclusive

All three of the LAMBDA™ Series instruments feature unmatched polarization measurement capabilities to match your analytical needs.

- Common beam depolarizer corrects for instrument bias.
- Automated polarizer/depolarizer drive in the large sample compartment provides further depolarization or allows study of oriented samples with polarized light.
- Compatible with transmission and reflectance measurements.

