

revvity

Introducing VICTOR Kira

multimode plate reader

Flexibility, performance, and endless possibilities.

This multimode reader offers detection for absorbance, fluorescence, and luminescence, and our monochromator-based technology allows for flexibility across multiple wavelengths.

With VICTOR Kira™, you'll enjoy a modern, easy to use software interface to support your complete workflow. The VICTOR Kira is a versatile, user-friendly solution that helps to empower your research.

Features:

- Versatile functionality across absorbance, fluorescence, and luminescence technologies for a wide variety of applications
- No need for exchanging filters when experimenting with different dyes
- Simple wavelength selection through intuitive software
- Characterize new fluorophores and analyze spectral shifts
- Quad monochromators help to greatly diminish stray light, allowing for more accurate and reliable data

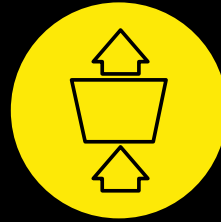


VICTOR Kira control

With the VICTOR Kira, you'll experience a sleek, intuitive software interface with built-in data analysis.

- **Versatile functionality** across absorbance, fluorescence, and luminescence technologies for a wide variety of applications
- **No need for exchanging filters** when experimenting with different dyes
- **Simple wavelength selection** through intuitive software
- Characterize new fluorophores and analyze spectral shifts
- **Quad monochromators** help to greatly diminish stray light, allowing for more accurate and reliable data
- **Versatile compatibility:** Operates on any device – laptop, tablet, etc. – and across all operating systems
- **Quick updates:** Stay current with the latest features through easy software updates
- **Ready out of the box:** Simple to set up and use from the start

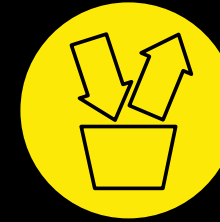
Absorbance



Absorbance applications

- ELISAs
- DNA/RNA quantification
- Protein quantification
- Bacterial growth assays

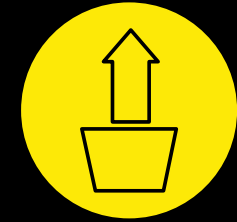
Fluorescence



Fluorescence applications

- Cell viability
- Cytotoxicity
- GPCRs
- Fluorescent protein quantification
- DNA/RNA quantification
- FRET

Luminescence



Luminescence applications

- Cell viability
- Reporter gene assays

Revvity's portfolio of plate readers is equipped with the latest detection technologies to meet the diverse assay requirements of labs today.

revvity