

# VICTOR Kira multimode plate reader



## Description

The VICTOR Kira™ multimode plate reader is designed with the intent to help streamline your lab work. 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.

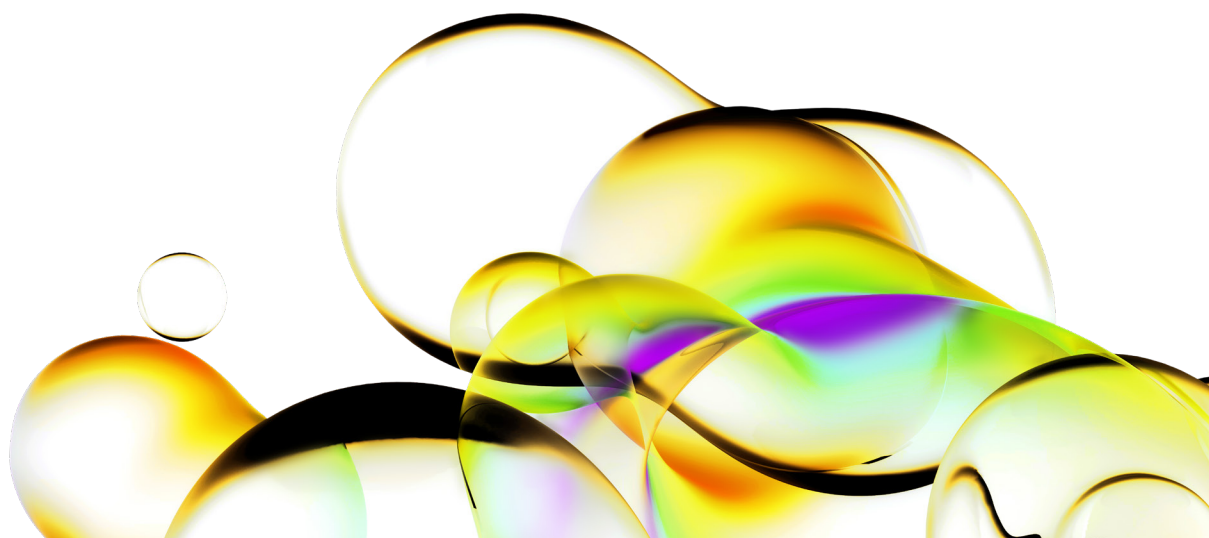
Complementing the VICTOR Kira, our comprehensive Revvity portfolio includes a wide array of assay kits and microplates. Additionally, our expert service and application support team is here to support you. The VICTOR Kira is a versatile, user-friendly solution that helps to empower your research.

## Detection technologies:

- Absorbance
- Fluorescence
- Luminescence

## Key features:

- **Quad monochromator detection:** Offers flexibility to run assays at multiple wavelengths
- **Three detection modes:** Absorbance, fluorescence, and luminescence for diverse applications
- **Modern software interface:** Intuitive and user-friendly supporting easy operation
- **Built-in calculations:** Provides analyzed data after measurement



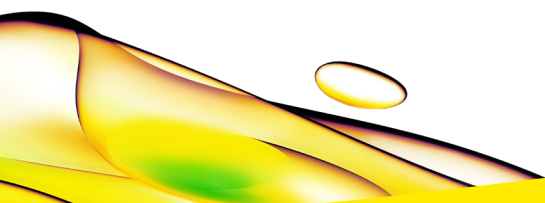
## Specifications

Physical Data		
Dimensions	Height	500 mm
	Width	330 mm
	Depth	550 mm
	Weight	45kg
Electrical requirements	Main voltage: 110-230V, 50/60 Hz	
Light source	Xenon flash lamp	
Plate formats	6-384 wp	
Temperature control	+3 °C above ambient up to 65 °C	
Shaking	Linear, orbital, double orbital	
Monochromator	Bandwidth	4 nm
	Wavelength range (increments)	230-850 nm (excitation), 1 nm increments; 350-900 nm (emission), 1 nm increments
Computer recommendations	Any operating system with common web browsers	

Typical Throughput (Time per Plate) (100ms exposure data)		
Detection	96-well	384-well
Absorbance	21s	60s
Fluorescence	22s	58s
Luminescence	22s	61s

Detection Limit Specifications		
<b>Absorbance</b>		
Dynamic range	0-4 OD	
Accuracy @ 2OD**	<1%	
Precision @ 2OD**	<0.2%	
Wavelength range	230-1000 nm	
Monochromator bandwidth	4 nm	
<b>Fluorescence</b>		
	<b>96-well</b>	<b>384-well</b>
Sensitivity (top reading)	0.6 fmol	0.2 fmol
Sensitivity (bottom reading)	<3 fmol	-
Wavelength range	Exc: 230-825 nm Ems: 245-850 nm	Exc: 230-825 nm Ems: 245-850 nm
Monochromator bandwidth (FWHM)	8 nm	8 nm
Dynamic range	>6 decades	>6 decades
<b>Luminescence</b>		
	<b>96-well</b>	<b>384-well</b>
Sensitivity (ATP) (top reading)	-	0.6 fmol/well (glow)
Wavelength range	400 nm-700 nm	

\*\* Measured at 490nm



revvity