



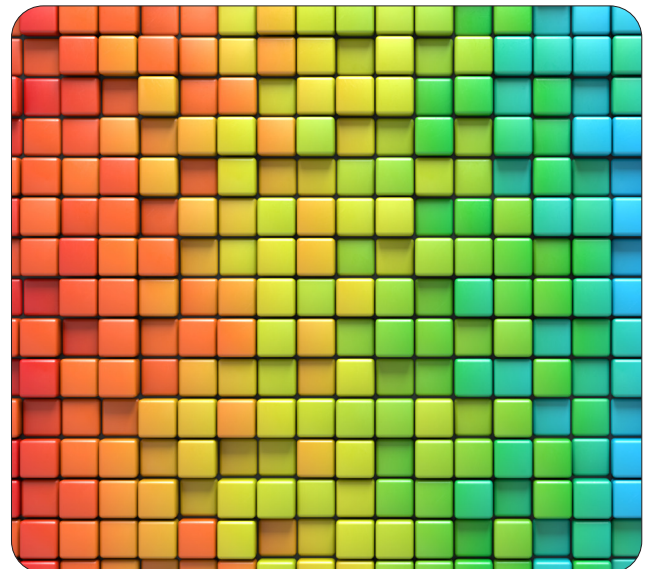
revvity

Greater sensitivity and flexibility with all the colors of the rainbow.

VICTOR Nivo rainbow filter exchanger:
Highly sensitive measurements and full spectrum flexibility for virtually any application

While being the industry's smallest multimode plate reader, the VICTOR® Nivo™ provides excellent performance, making it a great choice for multiuser labs that need a versatile workhorse instrument.

Every VICTOR Nivo comes rainbow-ready, giving you the flexibility of a monochromator, with the high sensitivity of a filter-based instrument.



At the heart of the VICTOR Nivo is a dynamic double-layered filter wheel system. With the capacity for up to 32 filters, all of the wavelengths you need for virtually any application are ready and available at any time.

With the rainbow filter exchanger, filters are:

- Exchanged between the inner and outer wheels
- Used for either Excitation or Emission lightpaths
- Locked in position so they don't get lost- ideal for multiuse labs

When fully loaded, the system is equipped to address virtually any application, with better sensitivity than a monochromator.



VICTOR Nivo multimode plate reader



Absorbance

Fluorescence intensity

Luminescence

TRF/TR-FRET

Fluorescence polarization

Alpha technology

Temperature control



Gas control



Dispenser



WiFi



GxP



Rainbow filter exchanger



To order, visit www.revvity.com or contact your sales representative.



revvity