

One million times more with **quad4** monochromators™

Our exclusive quad4 monochromator™ technology, as featured in our Infinite™ M200 and Safire²™ monochromator microplate readers, guarantees improved sensitivity, accuracy, flexibility and speed. Unwanted stray light is the most critical factor affecting the performance and sensitivity of microplate readers, but the quad4 technology blocks stray light and the resulting background noise by a factor of 10^6 times more than typical dual monochromator-based plate readers.

The unique quad4 configuration consists of two monochromators for excitation and two monochromators for emission, connected in series. The technology eliminates the need for cut-off filters to reduce stray light, providing more accurate spectral scans and avoiding the artificial peak shifts seen when monochromators are used in combination with filters. Fluorescence scans can be performed up to six times faster with quad4, compared with instruments containing two monochromators combined with cut-off filters. The advanced technology ensures consistent performance across the entire wavelength spectrum and allows you to tune in any wavelength without having to optimize filters.



The Safire² microplate reader comes with quad4 monochromators

“Fluorescence scans can be performed up to six times faster with quad4”