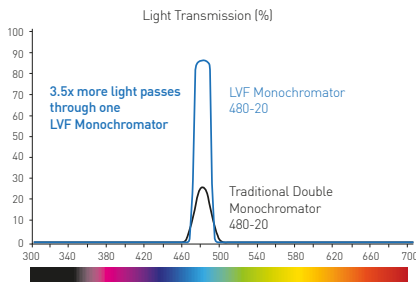


# 12 benefits of the CLARIOstar<sup>®</sup> Plus microplate reader

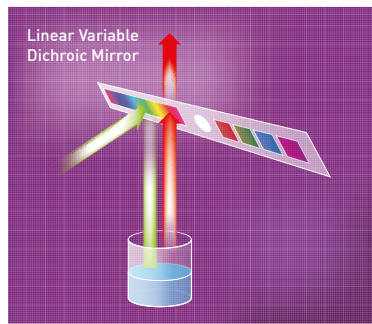
## LVF Monochromators™

The first and only monochromator using Linear Variable Filters, significantly increases light transmission and sensitivity over conventional monochromators based on diffraction gratings.



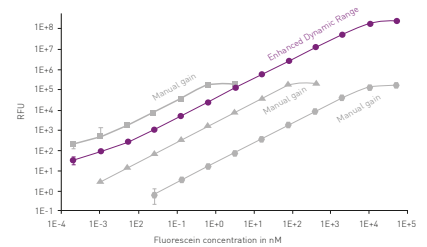
## Linear Variable Dichroic Mirror

A unique feature of the CLARIOstar Plus, it greatly reduces the background signal, improving sensitivity by more efficiently separating emission and excitation light.



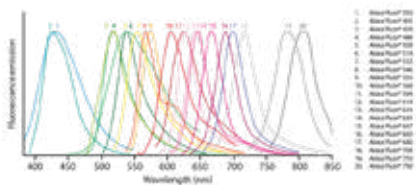
## Enhanced Dynamic Range (EDR)

EDR significantly simplifies measurement setup and provides an easier solution for assay development. It ensures reliable detection of samples at the largest range of concentrations and signal intensities (8 decades) in one measurement.



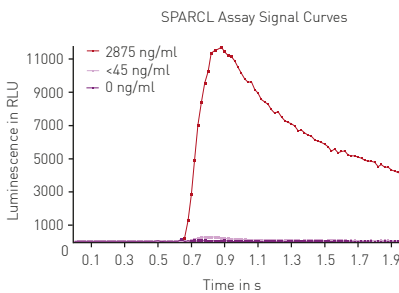
## Detect fluorophore emissions up to 900 nm

The extended wavelength PMT ensures high sensitivity detection across the entire spectrum. Superior performance with existing red fluorophores such as Alexa 647 and perfect for the ever expanding catalogue of red dyes.



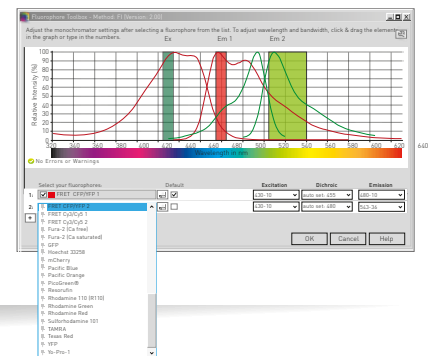
## Simultaneous injection and detection

An optimized injection system enables signal detection during injection. Even in the fastest of kinetic reactions, all data points will be reliably collected.



## Easy FRET and BRET detection

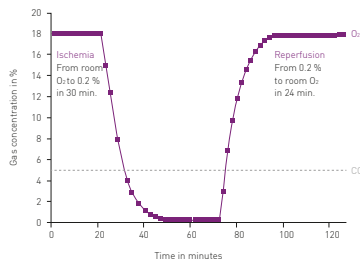
By displaying the spectra and optimizing settings by drag and drop, the Fluorophore-Toolbox significantly simplifies measurement definitions of dual emission assays.



# 12 benefits of the CLARIOstar<sup>®</sup> Plus microplate reader

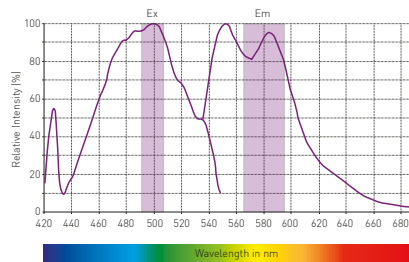
## Gas ramping

The ACU module offers the unique capability to run O<sub>2</sub> and CO<sub>2</sub> gas ramps. Samples can be maintained at constant CO<sub>2</sub> concentration during programmed de-oxygenation and rapid re-oxygenation to reflect ischemia/reperfusion for example.



## Customized fluorophore spectra

Choose one among more than 50 predefined settings. Alternatively, add fluorophores by scanning to the tool box and automatically determine optimal detection settings. This unique feature is made possible by the Fluorophore Toolbox.



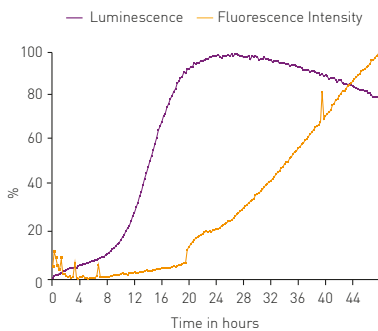
## Cross-talk reduction package

Hardware/software-based solutions help better discriminate between very low and high signals, even when they are proximal on the plate. Ideal for AlphaScreen<sup>®</sup> and luminescence based screening in 384 and 1536-well plates.



## Easy multiplex detection setup

Easily set up multiplex detection with only a few mouse clicks. Our Script Wizard enables a combination of up to 3 detection modes in one measurement run.



## Integrated barcode readers

Two optional barcode readers are positioned to read labels to the south and east of the microplate. Designed to read most common formats, they simplify data management in HTS or core labs.



## Ultra-fast spectrometer

Unlike conventional monochromators, our UV/vis absorbance spectrometer instantly captures an entire spectrum from 220 to 1000 nm in less than 1 second/well with no scanning needed.

