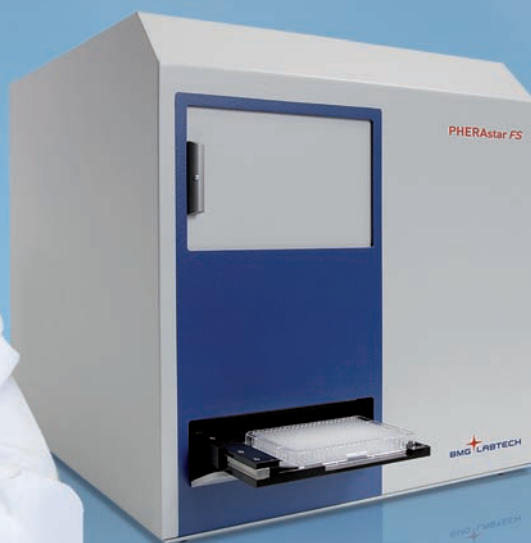


Simply the best.



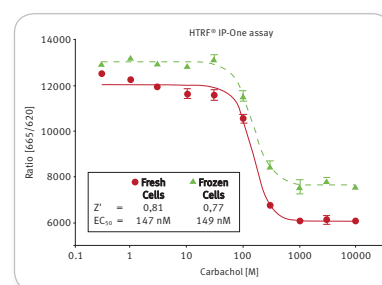
Reads plate formats up to
3456 well
BE ONE STEP AHEAD

PHERAstar FS - The Gold Standard Microplate Reader for High-Throughput Screening

High-end and unique features include:

- Laser for Alpha Technology
- Tandem Technology - Spectrometer / filter based detection
- Ultra fast UV/Vis spectrometer for full absorbance spectra
- Fluorescence Intensity, TRF, FP, Luminescence
- Dedicated UV-laser for TR-FRET measurements
- Advanced HTRF® performance
- Onboard reagent injection for precise kinetic measurements
- Advanced top and bottom reading
- High-resolution cell layer scanning
- Multi-user Control and Data Analysis Software MARS included

Find further information on www.bmglabtech.com



Dose response of carbachol from fresh and frozen cells using the IP-One assay

BMG LABTECH

The Microplate Reader Company

Australia · France · Germany · Japan · United Kingdom · United States

Alpha Technology includes AlphaScreen, AlphaLISA, and AlphaScreen SureFire. These assays are registered trademarks of PerkinElmer, Inc. HTRF is a registered trademark of Cisbio International.