

Double your throughput...



...with a single PHERAstar microplate reader using simultaneous dual emission detection.

The Premier HTS Microplate Reader – The PHERAstar is worth a double take. Ease-of-use. Sensitivity. Speed. Flexibility. All qualities desired in an HTS microplate reader. This multidetection microplate reader performs all of the leading HTS assays in every plate format up to 1536 wells. A dedicated laser excitation source ensures high-end AlphaScreen® assay performance. Furthermore, its simultaneous dual emission detection technology substantially reduces plate read times and greatly

increases performance in fluorescence polarization, FRET, BRET and TR-FRET assays such as HTRF®. The PHERAstar's cutting-edge sensitivity is also based on pre-installed, application specific optic modules that are used in combination with five photomultiplier tubes, optimized for each detection mode. Because of continuous innovations, BMG LABTECH is a world leader in plate reading systems, offering a broad portfolio of solutions for your life science research.

Microplate readers for all your assay needs



FLUOstar

PHERAstar

NOVOstar

NEPHELOstar

Stacker

**BMG LABTECH**

www.bmglabtech.com