

Love at first sight.



The new PHERAstar^{Plus} HTS microplate reader – the colleague you always dreamed of.

The PHERAstar Plus is a simultaneous dual emission multidetection HTS reader, offering Advanced HTRF® performance based on realtime ratio calculation. It is the only microplate reader featuring decay curve monitoring for assay optimization. A further hallmark is the fourfold increase in luminescence sensitivity.

The PHERAstar Plus provides FRET, TR-FRET, BRET, Fluorescence Polarization, and UV/Vis Absorbance measurements for plate formats up to 1536 wells. Proven assays include LanthaScreen™ and laser-based AlphaScreen®. Data analysis is achieved with the all new powerful MARS software.



HTRF is a registered trademark of Cisbio international.
Transreener is a patented technology of BellBrook Labs.
LanthaScreen is a registered trademark of Invitrogen Corp.
AlphaScreen is a registered trademark of PerkinElmer, Inc.



www.bmglabtech.com