

Your research is too important ...

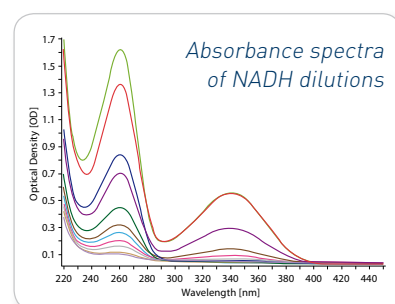


... to be left to others. Trust BMG LABTECH's
FLUOstar Omega microplate reader.

The FLUOstar Omega is based on BMG LABTECH's Tandem Technology: High-performance optical filters for exceptional detection limits, plus an ultra-fast UV/Vis absorbance spectrometer. The perfect solution for all your microplate based applications.

- Absorbance Spectra from 220 to 850 nm in 1s/well
- Fluorescence Intensity: < 0.2 fmol/well Fluorescein
- High-End Time-Resolved Fluorescence: < 3 amol/well Eu^{3+}
- Luminescence: 20 amol/well ATP
- HTRF®, LanthaScreen® certified*, DLReady™, AlphaScreen®

Visit our Applications Center at www.bmglabtech.com



Comprehensive MARS software
for powerful data analysis and
spectrum display.



HTRF is a registered trademark of Cisbio International.
LanthaScreen is a registered trademark of Invitrogen Corporation.
DLR is a trademark of Promega Corporation.
AlphaScreen is a registered trademark of PerkinElmer, Inc.



The Microplate Reader Company