

Keep your cells happy ...



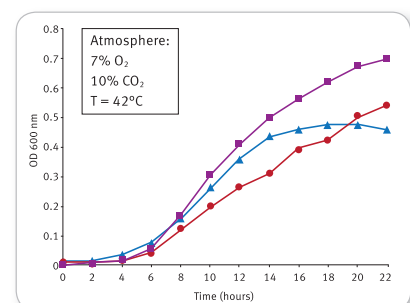
... with the **POLARstar Omega** microplate reader and the unique Atmospheric Control Unit (ACU).

With its independent control of O₂ and CO₂, the **POLARstar Omega** microplate reader with integrated ACU represents a true „walk away“ solution for long term cell based assays.

- Tandem Technology
- UV/Vis Absorbance Spectra (220-1000 nm)
- Fluorescence Intensity incl. FRET
- Fluorescence Polarization
- TRF, TR-FRET incl. HTRF® and LanthaScreen®
- Luminescence (flash and glow)
- AlphaScreen®/AlphaLISA®
- ACU – regulate O₂ (1-19%) and CO₂ (0-20%)

Visit us at **SLAS 2012** (booth #921).

Find further information on www.bmglabtech.com



Growth curve of three *Campylobacter* strains using the **POLARstar Omega** with the unique Atmospheric Control Unit.



HTRF is a registered trademark of Cisbio International.
LanthaScreen is a registered trademark of Invitrogen Corporation.
DLR is a trademark of Promega Corporation.
AlphaScreen and AlphaLISA are registered trademarks of PerkinElmer, Inc.
Transcreener is a registered trademark of BellBrook Labs.

BMG LABTECH

The Microplate Reader Company