





**Fig. 2:** Percentage reduction of alamarBlue® as a function of culture time, for substrates A and B.

## Conclusion

This application note shows that the reducing environment of cells can be accurately monitored using a BMG LABTECH microplate reader. The attraction of the alamarBlue® assay is that it incorporates a nontoxic reagent which allows continuous monitoring of cell proliferation on the same samples using either fluorescence or absorbance.

BMG LABTECH's readers are flexible multifunctional microplate readers that have six different measurement modes in one instrument: fluorescence intensity, time-resolved fluorescence, luminescence, fluorescence polarization, AlphaScreen® and absorbance. With two optional onboard injectors and a standard 45°C incubation chamber, BMG LABTECH microplate readers can easily become fully automated to perform any cell-based assays.



**PERAstar® FSX**

\*The PHERAstar FSX is the newest PHERAstar reader.



**CLARIOstar®**



**Omega Series**