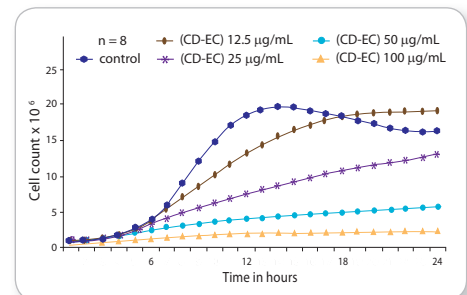




Laser-based microplate nephelometer for automated microbial growth monitoring

The NEPHELOstar is the world's only laser-based microplate nephelometer. It detects forward light scatter from particles in microplate-based samples using a sophisticated optical detection system and solid state laser.

- Monitoring of microbial growth kinetics
- High-throughput solubility screening
- Quantification of protein
- Monitoring of polymerization
- Stacker compatible
- Easy robotic integration



*Detecting the influence of antifungals on the growth of *Candida albicans* using nephelometry*

See more microplate readers on www.bmglabtech.com



PHERAstar FS



POLARstar



NOVOstar



FLUOstar



Stacker