

Stacker III

Mid-throughput
microplate handling system



Designed to optimize microplate handling
and increase sample throughput

- 20 and 50 microplate capacity magazines
- Accommodates all microplate formats
- Rapid plate loading, unloading and restacking
- Continuous load feature
- Barcode reader on left and front side of microplate
- User-definable advanced protocols

Stacker III

Software

BMG LABTECH's Control Software directly runs and coordinates both Stacker and plate reader. The script mode gives the user unlimited flexibility. It can be used to choose different test definitions for different plates in one batch run or to perform multiple measurements on one plate.



Technical specifications

Power requirements	85 to 240V AC, 50 to 60 Hz, 100 VA
Dimensions	Width 45 cm, depth 68 cm, height 19 cm Height with magazines 86 cm
Weight	15 kg
Barcode reader	Complies with Code 128, Code 39, Codabar, Code 11 UPC/EAN and 2/5 Interleaved and further barcodes

Compatible readers



PHERAstar® FSX



CLARIOstar®



Omega® Series



NEPHELOstar® Plus