

Stacker



Mid-Throughput Microplate Handling System

- Increased throughput with microplate readers
- 50 plate capacity magazines
- Continuous load feature
- Barcode reader reads from several positions
- User defineable advanced protocols
- Compatible with PHERAstar *Plus* / *FS*, Omega series, OPTIMA series, NEPHELOstar, and RUBYstar microplate readers

