



PR 3100 TSC

PR 3100 TSC Microplate Reader

The PR 3100 TSC is an intelligently designed compact, computer driven microplate reader which measures the absorbance of specimens in 96-well plates. This reader is configured with an on-board, fully functional computer that allows the user to manage the instrument operation and data reduction. Included in the configuration are five pre-installed filters (405, 450, 492, 550, 620 nm) which work directly with the assays that are validated on the reader. The broad colored touchscreen together with a wizard interface offers comprehensive data reduction with a streamlined operation for all assays.

Compact

- On-board computer
- Approximately 4000 plate results file storage

Streamlined Operation

- Comprehensive data reduction
- Validated assays on USB stick
- Rapid reading rates

Intelligent Design

- Broad colored touchscreen
- Wizard Interface
- 5 pre-installed filters
- Single and dual wavelength reading

