

## Absorbance 96 Plate Reader

As technology advances, scientific instruments have seen their use becoming more personal. No longer bulky, complicated machines, these devices are now small and portable. Who would imagine not having a personal mini-centrifuge on their benchtop?

Enzo is proud to introduce the next step of this revolution.

The Absorbance 96 Plate Reader is a personal 96-well format microplate reader that can be used for a variety of applications, including ELISA, protein quantification assays, cell-based assays, and endpoint and kinetic assays. By far the most compact reader on the market, the footprint of the plate reader is almost as small as the plate itself. Despite its scaled-down size, the Absorbance 96 delivers precise, accurate results at an affordable price. Together with proprietary software, the Absorbance 96 is ideal for a variety of applications extending from ELISA to protein- and cell-based assays.

### APPLICATIONS

- ELISAs – TMB, PNPP, ABTs, OPD
- Protein Assays – Bradford, BCA, Lowry
- Cell-based Assays – MTT, XTT, Cell Density

### Main Features

Space-saving design

Unique technical construction with 96 detection units

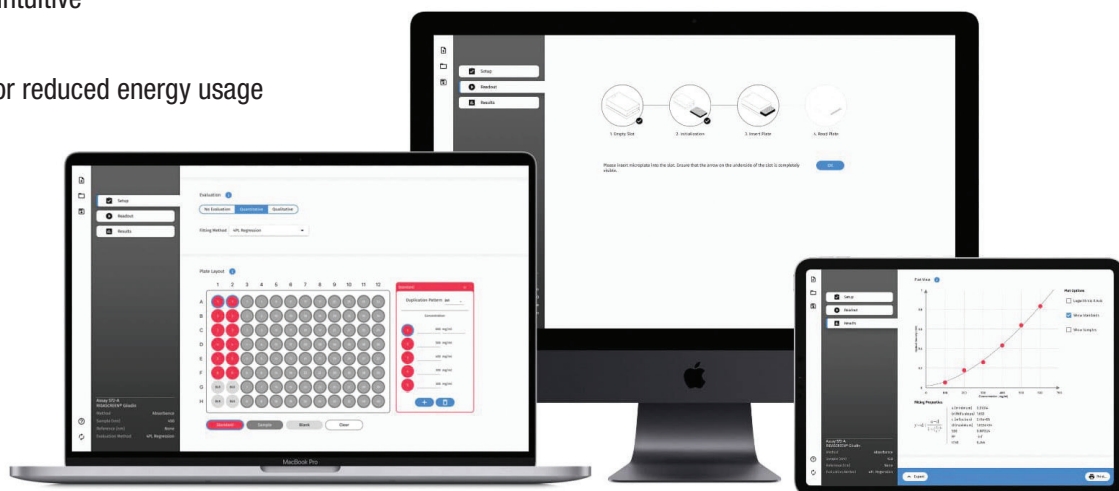
- » Enables rapid, reliable, and reproducible measurement
- » Maintenance-free

Proprietary software

- » User-friendly and intuitive

Affordable Price

Eco-friendly operation for reduced energy usage



# ABSORBANCE 96

## Specifications

### General

Part Number	ENZ-INS-A96
Detection method	Absorbance
Detection mode	Endpoint, kinetic
Microplate types	96-well microplates
Software	Custom software for external computer control

### Measurement

Light source	4 x LEDs
Detector	96 x Photodiodes
Wavelength selection	405, 450, 492, 620 nm
Photometric range	0.0-4.0 OD
Linearity	405 nm: $\leq 1.5\%$ (0-2 OD); $\leq 3.0\%$ (2-3 OD)/ $\geq 450$ nm: $\leq 1\%$ (0-2 OD); $\leq 1.5\%$ (2-3 OD)
Accuracy	405 nm: $\leq 1.5\% + 0.010$ OD (0-2 OD); $\leq 3\% + 0.010$ OD (2-3 OD)/ $\geq 450$ nm: $\leq 1\% + 0.010$ OD (0-2 OD) $\leq 1.5\% + 0.010$ OD (2-3 OD))
Reproducibility	$\leq 0.5\% + 0.005$ OD (0-2 OD); $\leq 1\% + 0.010$ OD (2-3 OD)
Resolution	0.001 OD
Read time	Down to 5 sec at single wavelength
Data output	USB 2.0 interface with PC

### Physical Characteristics

Dimensions	9.6 cm x 15.4 cm x 5.5 cm (W x L x H)
Weight	0.9 kg (2 lbs)
Power	Through USB connection 5V
Power consumption	2.5 Watts
Regulatory	CE marked

**Global Headquarters**  
ENZO LIFE SCIENCES, INC.  
Ph: 800.942.0430  
info-usa@enzolifesciences.com

**European Sales Office**  
ENZO LIFE SCIENCES (ELS) AG  
Ph: +41 61 926 8989  
info-eu@enzolifesciences.com

**Belgium, The Netherlands  
& Luxembourg**  
Ph: +32 3 466 0420  
info-be@enzolifesciences.com

**France**  
Ph: +33 472 440 655  
info-fr@enzolifesciences.com

**Germany**  
Ph: +49 7621 5500 526  
info-de@enzolifesciences.com

**UK & Ireland**  
Ph: 0845 601 1488 (UK customers)  
Ph: +44 1392 825900  
info-uk@enzolifesciences.com

For local distributors and detailed product information visit us online: [www.enzolifesciences.com](http://www.enzolifesciences.com)

All product names, logos, and brands are the property of their respective owners.