

Absorbance 96 Plate Reader

Personal, Compact Plate Reader

The Absorbance 96 is a uniquely designed microplate reader for assays in 96-well format. 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 form, the Absorbance 96 delivers precise, accurate results at an affordable price. Together with its proprietary software, the Absorbance 96 is ideal for a variety of applications extending from ELISA to protein- and cell-based assays.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ENZ-INS-A96	1ea
-------------	-----

Manuals, SDS & CofA

[View Online »](#)

- Space Saving Design
- Fast Readout Speed
- Reliable Measurement Results
- Maintenance-free
- Solid State Technology



Handling & Storage

Shipping Ambient Temperature

Regulatory Status CE - CE Marked, RUO - Research Use Only

Product Details

Accuracy	405 nm: $\leq 1.5 \% + 0.010 \text{ OD (0–2 OD)}$; $\leq 3 \% + 0.010 \text{ OD (2–3 OD)}$ $\geq 450 \text{ nm: } \leq 1 \% + 0.010 \text{ OD (0–2 OD)}$ $\leq 1.5 \% + 0.010 \text{ OD (2–3 OD)}$
Connectivity/Communication	Through USB connection 5 V
Contents	Absorbance 96 Plate Reader Micro-USB Cable Hex Key Microfiber Cloth USB Stick with Software
Detection Method	Absorbance
Detection Mode	Endpoint, Kinetic
Dimensions (LxWxH)	15.4 x 9.6 x 5.5 cm
Light Source	4 x LEDs
Linearity	405 nm: $\leq 1.5 \% (0–2 \text{ OD})$; $\leq 3.0 \% (2–3 \text{ OD})$ $\geq 450 \text{ nm: } \leq 1 \% (0–2 \text{ OD})$; $\leq 1.5 \% (2–3 \text{ OD})$
Photometric Range	0.0–4.0 OD
Power	2.5 Watts
Reproducibility	$\leq 0.5 \% + 0.005 \text{ OD from } 0.0–2.0 \text{ OD}$ $\leq 1.0 \% + 0.010 \text{ OD from } 2.0–3.0 \text{ OD}$
Resolution	0.001 OD
Software	Compatible with Microsoft Windows and MacOS

Technical Info / Product Notes

Application Note:

[Benchmark Analysis of Novel Portable Absorbance Microplate Reader](#)

Throughput

96-well microplates

Voltage

5 V

Wavelength

405, 450, 492, 620 nm

Weight

0.9 kg

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)