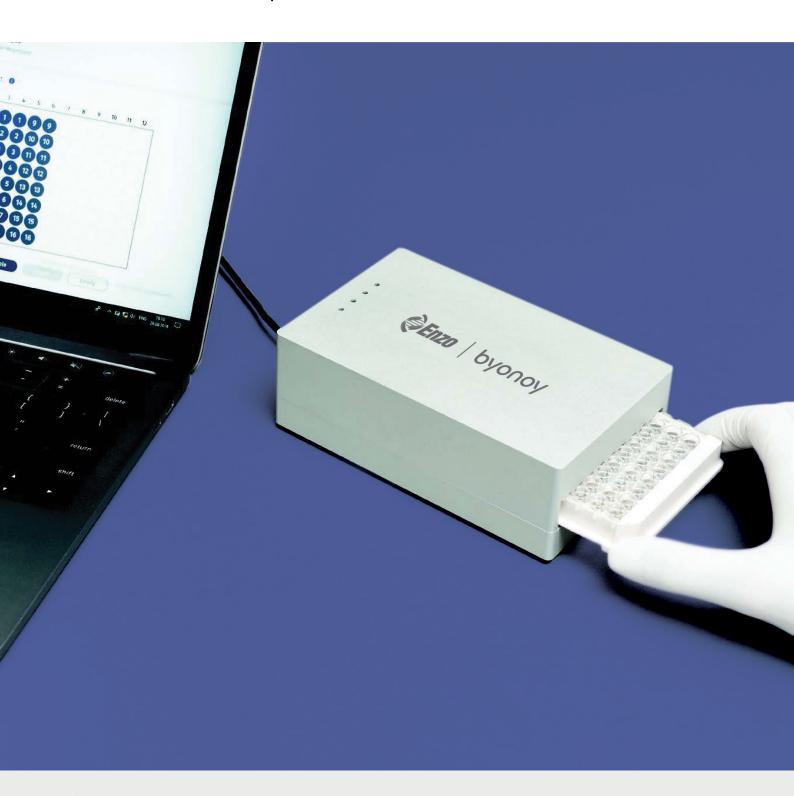
Absorbance 96

A Personal, Compact Plate Reader







Absorbance 96 – Innovative Compactness

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 reader is almost as small as the plate itself. Despite its size, this plate reader delivers precise and 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.

Utilize this plate reader with our catalog of nearly 300 ELISA kits, which includes sensitive, specific assays for relevant markers of cell viability, signaling pathways, steroid and peptide hormones, inflammation, and more. In addition to traditional antibody pair-based immunometric assays, our scientists specialize in the particularly challenging production of high-specificity competitive ELISAs for detection of small molecules. Enzo has over 20 years of assay development experience and state-of-the-art manufacturing facilities to ensure time-tested reproducibility.



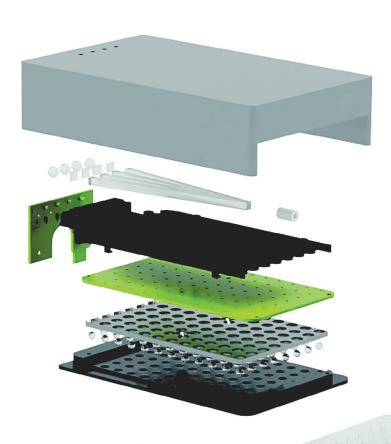
Usability

The goal was to develop a reader that would simplify the workflow in the laboratory. This philosophy of intuitive operation was translated into the Absorbance 96. The unique open design, in combination with the small size of the reader, leads to an entirely new user experience. A simple USB connection provides both the power supply and access to user-control software, which itself functions via plug-and-play technology. Easily transported within – and between – different workspaces and laboratories, the Absorbance 96 fits into every lab, saving bench space and providing unprecedented flexibility.

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



Advanced Technology



The Absorbance 96 is the first microplate reader on the market to contain 96 individual detection units. This means that all wells can be read simultaneously allowing for extremely fast photometric measurements. Without the need to scan across multiple wells, the Absorbance 96 has no moving parts. In combination with long-life LEDs, this solid-state design provides a maintenance-free user experience and ensures reliable, high quality results.

The four standard filters – 405, 450, 492, 620 nm – allow for ELISAs with a wide variety of substrates including PNPP, ABTS, OPD and TMB. Other filter combinations for protein (Bradford or BCA), cell-based (MTT, XTT), or cell density assays can also be provided.



- 96 detection units for fast reading
- Four-channel LED optics
- No scanning mechanics
- Maintenance-free
- 5-second read time capability

Software



The Absorbance 96 software is as flexible as the reader itself. It can be used on multiple devices and supports Windows as well as macOS. With plug-and-play technology, simply connect the Absorbance 96 and start your assay.



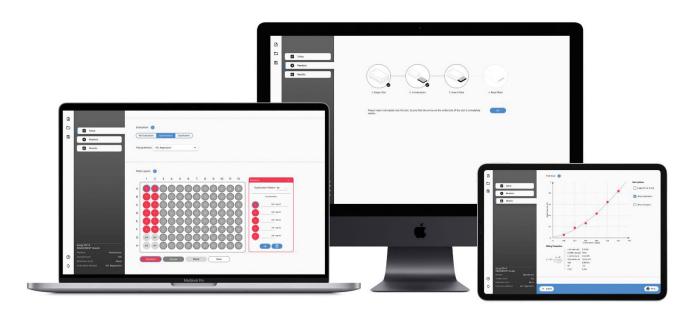


Windows

mac0S

MAIN FEATURES

- Endpoint and kinetic assays
- Quantitative and qualitative evaluation
- Custom protocols and reports
- Unlimited access across different devices



Pair with Quality ELISA Kits

We offer a large variety of immunoassay kits in both immunometric and competitive assay formats compatible with the Absorbance 96 Plate Reader. As scientists and manufacturers of kits, we understand the critical nature of your research. Each kit is put through rigorous testing to ensure high precision, accuracy, sensitivity, and specificity. You can be confident that you will obtain reproducible results, day-after-day and lot-after-lot.

- Consistent strict quality control guidelines ensure reproducible results, lot-after-lot
- Trusted reliability highlighted in peer-reviewed publications for three decades
- Flexible adaptable to automated equipment for high-throughput systems

EASY ANALYSIS

Absorbance 96 protocol files are available to download for selection of our ELISA kits.

Cyclic Nucleotides

The most sensitive colorimetric ELISA kits for detection of cAMP and cGMP, key second messenger signaling molecules.

Sensitive assays for steroid and peptide hormones

Steroid

Cortisol

DHEA

Estriol

FSH

LH

Melatonin

Progesterone

Testosterone

25(0H) Vitamin D

Aldosterone

Corticosterone

17β-Estradiol

that regulate stress, reproduction, metabolism,

cardiovascular physiology, and more.

Cyclic AMP Complete Cyclic GMP Complete Cyclic AMP Direct (Intracellular)

Cyclic GMP Direct (Intracellular)

Cyclic GMP (Extracellular) Cyclic GMP (Extracellular)

Hormones

Peptide

Gastrin

GLP-1

Insulin

Oxytocin

Proinsulin

Prolactin

Substance P

Vasopressin

Monoamine

Dopamine

Histamine

Serotonin

Angiotensin

Bradykinin

Endothelin-1

LVV-Hemorphin

Neuroscience

High-specificity, sensitive assays for biomarkers of neurodegeneration and neural signaling.

24(S) Hydroxycholesterol

Heat Shock & Oxidative Stress

The definitive collection of ELISA kits for heat shock and oxidative stress.

HSP60

HSP70 & HSP70 High Sensitivity

Grp94

Heme Oxygenase-1

Toxicoloay

Biomarker detection for assessing response to toxic compounds.

NGAL KIM-1

Wnt Pathway

Quantify key components of the Wnt pathway critical to cancer, development, bone remodeling, and neurodegeneration.

Dkk-1 **β-catenin**

24(S) Hydroxycholesterol

Dopamine

Serotonin

Survival Motor Neuron (SMN1)

APP∆C31

HSP27

Grp75

Grp78

DNA Damage (8-OHdG)

Bioprocess

Sensitive ELISA kits for biotherapeutic characterization and residual contaminant detection.

PEGylated Protein Protein A HEK²⁹³HCP CHO HCP

E. coli HCP

Programmed Cell Death

Quantify critical cell death pathway proteins. The perfect complement to our CELLESTIAL® cell-based assays.

p62 (Sequestosome 1)

Bax Bcl-2

Cytochrome C Survivin

XIAP

Immunology

Broad portfolio of sensitive, highspecificity ELISA kits for inflammation and immunological studies.

Adiponectin IL-33 CD40/CD40L Leptin **CRP** LTB_₄

Cysteinyl Methotrexate Nampt Leukotrienes **NGAL** FasL

Osteoprotegerin 12(S)-HETE 8-iso-PGF_{2a} 15(S)-HETE 6-keto-PGF_{1α} Histamine PGE₂ FPIA & CLIA 13(S)-HODE

PGE₂ IFN-γ $PGF_{2\alpha}$ IL-1B **PSA** IL-2 **RANKL** IL-4 Resistin IL-5

SARS-CoV-2 laG IL-6 TGF-β1 IL-8 TIMP-1 IL-10 TNF-α IL-12p70 Transferrin

IL-13 TXB₂ IL-17A **VEGF**

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: ≤ 1.5 % $(0.0-2.0 OD)$ /≥ $450 nm$: ≤ 1.0 % $(0.0-2.0 OD)$; ≤ 1.5 % $(2.0-3.0 OD)$)
Accuracy	405 nm: ≤ 1.5 % + 0.010 0D (0.0–2.0 0D)/≥ 450 nm: ≤ 1.0 % + 0.010 0D (0.0–2.0 0D) ≤ 1.5 % + 0.010 0D (2.0–3.0 0D)
Reproducibility	405 nm: ≤ 0.1 % + 0.005 0D (0.0–2.0 0D)/≥ 450 nm: ≤ 0.1 % + 0.005 0D (0.0–3.0 0D)
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

& Luxembourg Ph: +32 3 466 0420 info-eu@enzolifesciences.com info-be@enzolifesciences.com

Belgium, The Netherlands

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

Germany +49 7621 5500 526 Ph: info-de@enzolifesciences.com **UK & Ireland** Ph: 0845 601 1488 (UK customers)

Ph: +44 1392 825900 info-uk@enzolifesciences.com

For detailed product information visit us online: www.enzolifesciences.com All product names, logos, and brands are the property of their respective owners.

Put our experience to work for you!

Visit www.enzolifesciences.com to learn about our innovative tools for assay kits.