

# History of ELISA

ELISA (enzyme-linked immunosorbent assay) is a method used to quantitatively detect an antigen within a sample. An antigen is a toxin or other foreign substance, for example a flu virus or environmental contaminant, that causes the vertebrate immune system to mount a defensive response. The range of potential antigens is vast, so ELISAs are used in many areas of research and drug discovery on a wide variety of sample types. Cell lysates, blood samples, food items, and more can be analyzed for specific substances of interest using ELISAs.

**1941**—Albert H. Coons and his colleagues are the first to label antibodies with a fluorescent dye, and use it to identify antigens in tissue sections. This method is known today as **immunofluorescence**.<sup>1</sup>

**1971**—Eva Engvall and Peter Perlman (independently) invent a method that revolutionized medicine called the **ELISA test**. The method uses antibodies to seek out the presence of hormones or viruses.<sup>3,4</sup>

**1977**—**Sandwich ELISA method**, in which the detection antibody is coated onto the plate surface before the protein of interest is added, is developed and tested on several substrates for proof-of-concept.<sup>6</sup>

**1985**—The ELISA test is the first screening test commonly employed for **HIV**. It was approved for use on March 2, 1985.<sup>9</sup>

**SpectraMax® iD3 Multi-Mode Microplate Reader**

Absorbance capable reader featuring a large, intuitive touchscreen and multimode detection abilities



**ABS and ABS Plus Absorbance ELISA Microplate Readers**

Absorbance readers providing consistent results with steady temperature regulation and tunable wavelength selection



**SoftMax® Pro Data Analysis Software**

Designed to provide the simplicity, flexibility and power required for advanced data analysis

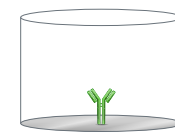


## References

1. Coons, A. H. The beginnings of immunofluorescence. *J. Immunol.* 87, 499–503 (1961).
2. Yalow, Rosalyn and Berson, Solomon. "Immunoassay of endogenous plasma insulin in man." *The Journal of Clinical Investigation.* 1960;39: 1157–75.
3. Perlmann, Peter et al. "Enzyme-linked immunosorbent assay (ELISA) quantitative assay of immunoglobulin G." *Immunochemistry.* 1971;8 (9): 871–4.
4. Schuur, A. "Immunoassay using antigen—enzyme conjugates." *FEBS Letters.* 1971;15 (3): 232–236.
5. Yorde, Donald et al. "Competitive Enzyme-Linked Immunoassay with Use of Soluble Enzyme/Antibody Immune Complexes for Labeling. I. Measurement of Human Chorionadotropin." *Clin. Chem.* 1976;22/8;1372–1377
6. Kato, K et al. "Use of rabbit antibody IgG bound onto plain and aminoalkylsilyl glass surface for the enzyme-linked sandwich immunoassay." *J Biochem.* 1977 Jul;82(1):261–6..
7. Lindström, P et al. "IgG autoantibody to human serum albumin studied by the ELISA-technique." *Scand J Immunol.* 1978;7(5):419–25.
8. Czerkinsky, C et al. "A solid-phase enzyme-linked immunospot (ELISPOT) assay for enumeration of specific antibody-secreting cells." *J Immunol Methods.* 1983;65 (1–2): 109–121.
9. Alexander, Thomas. "Human Immunodeficiency Virus Diagnostic Testing: 30 Years of Evolution." *Clinical and Vaccine Immunology.* 2016 Apr;23(4):249–253.

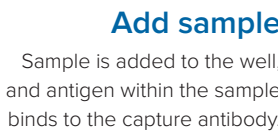
## Workflow of an ELISA protocol

The workflow of a typical sandwich ELISA protocol has multiple reagent addition, incubation and wash steps. Here we've highlighted each step and the instrumentation and tools needed to conduct the ELISA assay.

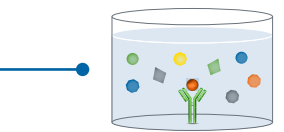


### 1 Capture antibody binds to wells

First, the capture antibody is bound to the bottom of the microplate well.

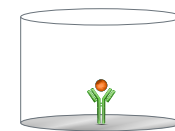


**Add sample**  
Sample is added to the well, and antigen within the sample binds to the capture antibody.



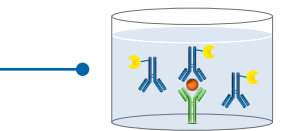
### 3 Wash microplate

Unbound material is washed away, leaving only the antigen of interest and minimizing the potential for high background signal.



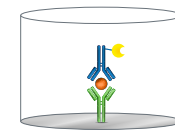
### 4 Add detection antibody

Enzyme-conjugated detection antibody binds to a second site on the antigen of interest, providing the means to detect the antigen.



### 5 Wash microplate

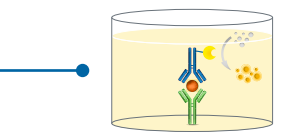
Unbound antibodies are washed away, leaving only those specific for the target of interest and again minimizing the potential for background signal.



### 6 Add substrate

Substrate is converted by the enzyme on the detection antibody, producing a color change, with intensity proportional to the amount of antigen present.

Depending on the enzyme and substrate used, the readout can also be fluorescent or luminescent.



### 7 Read plate

The microplate reader detects the colored reaction product and outputs optical density (OD) values that indicate how much light is absorbed by the contents of each well.



### 8 Calculate results

The amount of antigen in each sample is calculated, and different samples—for example, cells subjected to different treatment conditions—can be compared.

