

# ApoE ELISA kit

**Highly sensitive ApoE ELISA kit enabling detection of ApoE in serum, plasma, cell culture supernates, and tissue homogenates in 3.5 hours.**

The ApoE ELISA kit is a colorimetric immunometric immunoassay kit with results in 3.5 hrs.

Apolipoprotein E (ApoE) transports lipoproteins, fat-soluble vitamins, and cholesterol into the lymph system and then into the blood. It is synthesized principally in the liver, but has also been found in other tissues such as the brain, kidneys, and spleen. In the nervous system, non-neuronal cell types, most notably astroglia and microglia, are the primary producers of ApoE, while neurons preferentially express the receptors for ApoE. ApoE was initially recognized for its importance in lipoprotein metabolism and cardiovascular disease. ApoE has been studied for its role in several biological processes including: type III hyperlipoproteinemia (HLP III), Alzheimer's disease (AD), immunoregulation, and cognition.

**Manuals, SDS & CofA**

[View Online »](#)

- Sensitive measurement of ApoE, detecting as little as 14.063 ng/ml
- Negligible reactivity to similar proteins.
- High throughput with results in 3.5 hours.
- Fully quantitative results that surpass semi-quantitative Western blot analysis

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Apolipoprotein E

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of ApoE in serum, plasma, cell culture supernates and tissue homogenates in human samples.

**Assay Time** 3.5 hours

**Compatibility** This product is compatible with the [Absorbance 96 Plate Reader](#).

**Contents** Microtiter Plate, Standard, Sample Buffer, Wash Buffer Concentrate, Antibody, Antibody Buffer, Conjugate, Conjugate Buffer Substrate, and Stop Solution

**Sensitivity** 14.063 ng/ml (23.438-1500 ng/ml)

**Species Reactivity** Human

**UniProt ID** P02649

**Wavelength** 450 nm

**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-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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