

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.

Ordering Information

[Order Online »](#)

ENZ-KIT134-0001	96 wells
-----------------	----------

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

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

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

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

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