

Product Information & ELISA Manual

Non-human Primate Apolipoprotein B/ApoB
ELISA Kit
NBP3-11737

Enzyme-linked Immunosorbent Assay
for quantitative detection.

Contact

Web: www.bio-technne.com/brands/novus-biologicals/
Email: nb-customerservice@bio-technne.com
P: 888.506.6887 // P: 303.730.1950 // F: 303.730.1966

Novus kits are
guaranteed for 6 months
from date of receipt.

**For research use only.
Not for diagnostic or
therapeutic procedures.**

Table of Contents

Introduction	4
Shipping and storage	4
Contents.....	5
Preparation.....	6
Protocol	8
Performance.....	10

Introduction

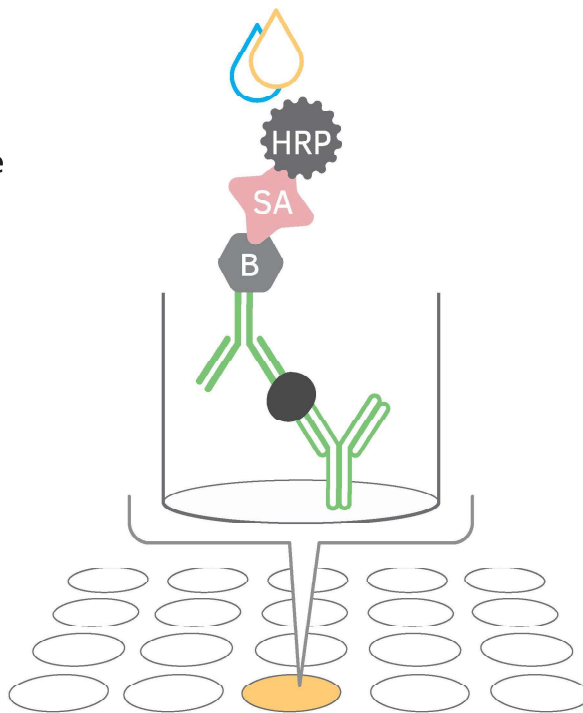
Novus Biologicals carefully validated ELISA Pro kits provide all the necessary reagents to conveniently quantify analytes in serum, plasma, and cell culture supernatants in a robust, sensitive, and specific manner. For Research Use Only (RUO). Not for use in diagnostic procedures.

ELISA assay principle

ELISA Pro kits are supplied with ELISA strip plates precoated with monoclonal antibody (mAb). Analyte in the sample is captured by the coated mAb and detected by the biotinylated detection mAb followed by Streptavidin-HRP (SA-HRP). Addition of TMB substrate will result in a colored substrate product. The reaction is stopped with sulfuric acid and the optical density can be quantified using an ELISA plate reader. The concentration of analyte is determined by comparison to a serial dilution of the ELISA standard analyzed in parallel.

Analysis of serum and plasma samples

The ELISA Pro kits include Apo ELISA buffer, a buffer that prevents false-positive signals. The buffer blocks heterophilic antibodies from cross-linking the assay antibodies. Heterophilic antibodies are commonly found in human serum/plasma and can also be present in other species. The buffer has been validated using serum/plasma samples from healthy human blood donors.



Shipping and storage

The kit is shipped at ambient temperature. All reagents should be stored at 4-8 °C upon receipt, except the standard, which should be stored at -20 °C. The expiry date indicates how long unopened products, stored according to instructions, are recommended for use. Do not combine components from different kit batches or components from other suppliers.

Contents

Component	1-plate kit		
Pre-coated ELISA strip plate: Anti-apoB mAbs LDL 20/17	1 x 96 wells		
Human apoB ELISA standard: Purified LDL in glycerol	1 vial		
Detection mAb LDL 11, biotin (1 mg/ml)	15 µl		
Streptavidin-HRP	15 µl		
Wash buffer concentrate (20x)	120 ml		
Apo ELISA buffer concentrate (5x)	60 ml		
Streptavidin-HRP diluent	15 ml		
Sample solvent	15 ml		
TMB substrate	15 ml		
Stop solution	15 ml		
Adhesive plate covers	3		

To ensure total recovery of the stated quantity, bottles and vials have been overfilled.

Materials required but not supplied

- Microplate reader capable of reading at 450 nm
- ELISA plate washer; automated or manual (e.g., multipipette or squirt bottle)
- Precision pipettes, tips, and graduated cylinders
- Tubes for standard and sample dilutions
- Distilled or deionized water

Safety information

The Stop solution, 0.18 M H₂SO₄ (< 1%), is irritating to eyes and skin and should be handled with care. The standard should also be handled carefully as the effects of exposure are unknown. Buffers and reagents in solution contain the preservative Kathon CG (0.002%), a potential allergen that may cause sensitization through skin contact. Human and animal samples should be treated as potentially hazardous biologic material. All material should be disposed of in accordance with local regulations. For further information please consult the Safety Data Sheet on our website.

Preparation

- Allow the plates and assay reagents to reach room temperature before starting the assay (except for the TMB substrate which should preferably be used cold).
- Plan the plate layout to include a standard curve, samples, and an assay background control, all in duplicate. The volume per well should not exceed 100 μ l. Include a plate blank (wells with only Substrate and Stop solution) to be used for subtraction before analysis.

Wash buffer

Add 50 ml Wash buffer concentrate to 950 ml distilled or deionized water (sufficient for all washing steps of 1 plate). If crystals have formed in the 20x concentrate, bring to room temperature and mix gently to dissolve.

Apo ELISA buffer

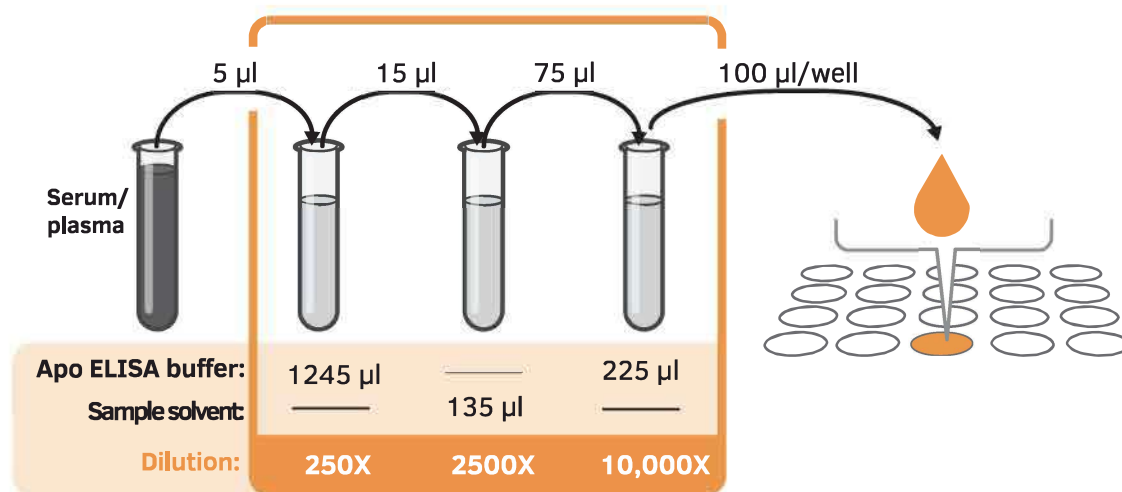
Prepare the required volume of Apo ELISA buffer by diluting Apo ELISA buffer concentrate 5-fold with distilled or deionized water. For each plate, add 30 ml Apo ELISA buffer concentrate to 120 ml water.

Samples

All samples should be diluted in Apo ELISA buffer in tubes or plates, adding the buffer before samples. Make sure to remove visible precipitates and avoid using strongly hemolyzed or hyperlipemic samples. To prevent interference by varying LDL particle sizes, all serum- and plasma-derived samples should be treated with Sample solvent, followed by vortex for 5 seconds. This treatment will not interfere with the analysis of other apolipoproteins, and it is not required when detecting apoB produced in cell lines. Avoid freeze-thawing of samples.

Dilution guidelines for human serum and plasma

For serum and plasma samples, we recommend the dilutions shown below. Note that an initial dilution in Apo ELISA buffer is required before the dilution in Sample solvent. Precise pipetting is important, change tips between dilution steps. Indicated volumes are sufficient for duplicates.

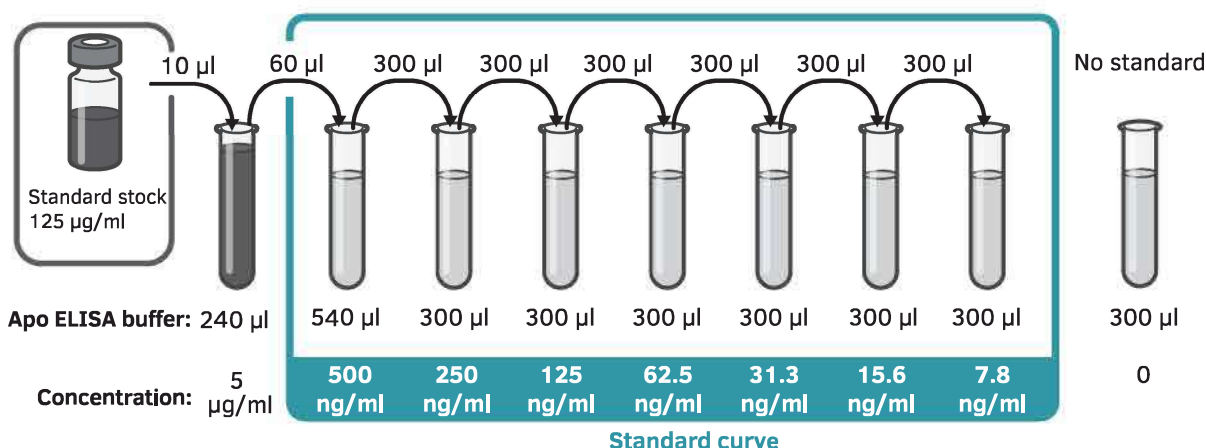


ELISA standard

The apoB standard is supplied as purified LDL stabilized by 50% glycerol. The concentration is 125 µg/ml. It is not necessary to aliquote the standard as the high content of glycerol keeps the standard in a liquid state. Store at -20 °C.

Preparation of standard curve

Dilute the standard stock solution to create a standard curve as shown. The indicated volumes are sufficient for duplicates. The last vial is used as an assay background control, i.e., the standard should be omitted. Prepare the standard curve within 30 minutes of use.



Detection antibody

Dilute the detection mAb in Apo ELISA buffer to a concentration of 1 µg/ml within 15 minutes of use. For each plate, add 12 µl detection mAb to 12 ml Apo ELISA buffer.

Streptavidin-HRP

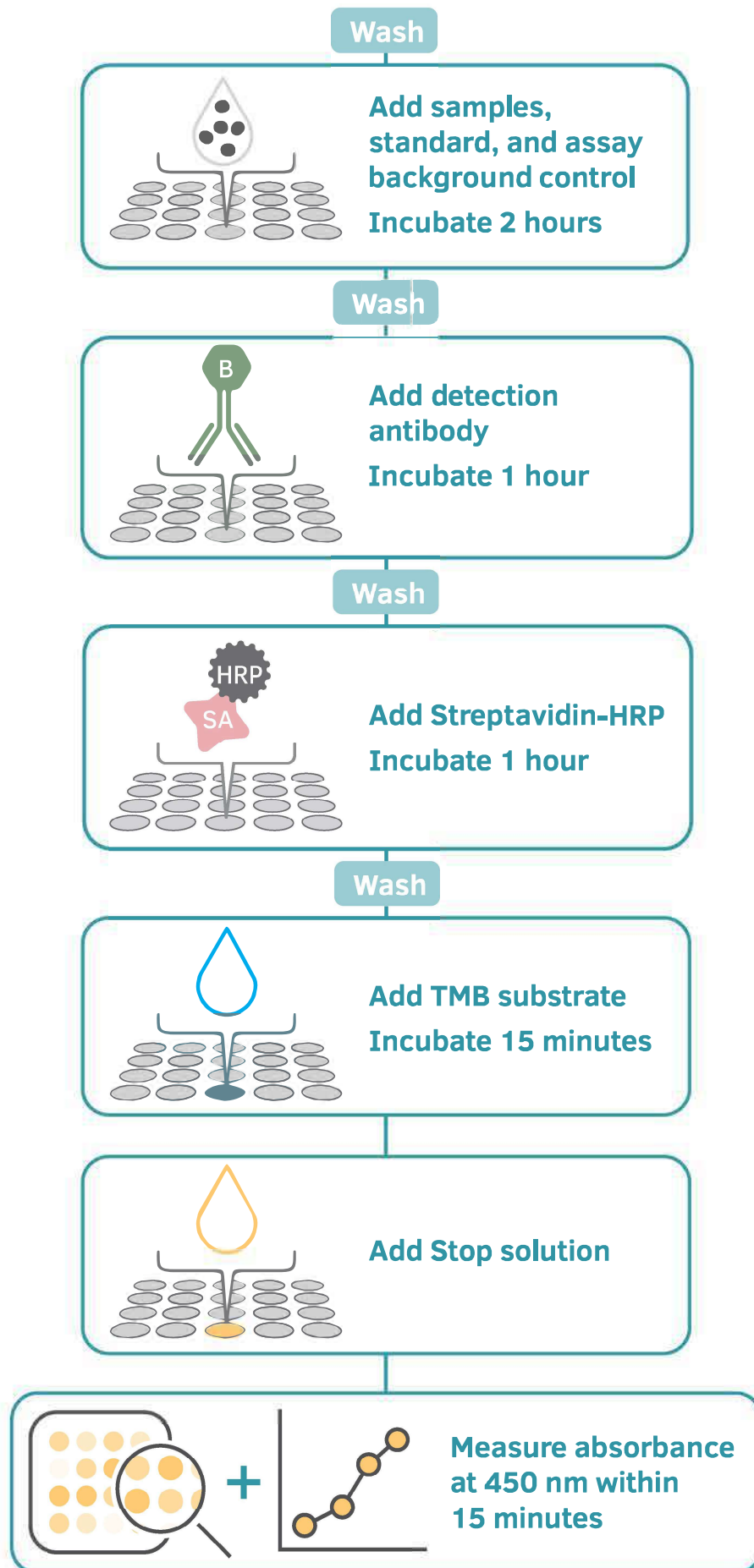
Dilute the Streptavidin-HRP 1:1000 in Streptavidin-HRP diluent within 15 minutes of use. For each plate, add 12 µl Streptavidin-HRP to 12 ml Streptavidin-HRP diluent.

Protocol

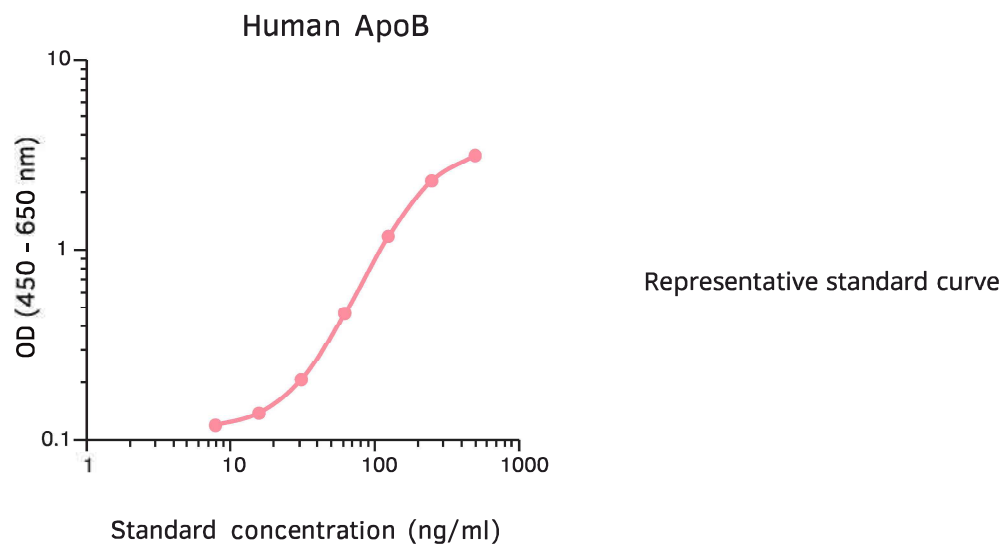
Prepare the reagents, standard curve, and samples as described in the Preparation section. Assemble the required number of strips in the plate frame and label the top of each strip. Store the remaining strips in the foil bag containing the desiccant at 4-8 °C.

1. Wash the plate 5 times with wash buffer, 300 µl per well. After the final wash, invert and tap the plate firmly against absorbent paper. Immediately proceed to the next step.
2. Add 100 µl per well of samples (diluted at least 2-fold), standard, and assay background control. Mix by tapping the plate. Cover the plate with an adhesive plate cover and incubate for 2 hours at room temperature.
3. Wash as in step 1.
4. Add 100 µl per well of detection mAb. Cover the plate and incubate for 1 hour at room temperature.
5. Wash as in step 1.
6. Add 100 µl per well of Streptavidin-HRP. Cover the plate and incubate for 1 hour at room temperature.
7. Wash as in step 1.
8. Add 100 µl of TMB substrate to each well. Incubate at room temperature, protected from direct light for 15 minutes.
9. Add 100 µl of Stop solution to each well to stop the color development.
10. Measure absorbance at 450 nm within 15 minutes. Preferably use a reader capable of subtracting a reference wavelength between 570 and 650 nm.

We recommend the use of an ELISA software utilizing a 4- or 5-parameter curve fit. Subtract the mean absorbance value of the blank from the samples, standard and assay background control prior to creating the standard curve and analyzing the results.



Performance



Standard range 7.8-500 ng/ml

Sensitivity 7 ng/ml

The lowest concentration that can be detected, but not necessarily quantified with precision and accuracy. This was determined by adding 2 standard deviations to the mean OD of background wells.

Calibration

No international standard exists for calibration of solubilized apoB. Please note that calibration is batch specific.

Precision

Sample	Intra-assay		Inter-assay	
	1	2	1	2
n	8	8	6	6
Mean (ng/ml)	268.3	146.8	261.0	136.6
SD	8.5	3.0	26.2	14.8
CV%	3.2	2.0	10.0	10.8

Intra-assay and inter-assay precision were determined at 2 different concentrations of analyte (8 replicates per concentration in 6 assays).

Linearity

Dilution of a human plasma gives a mid-curve recovery of 90-114% in repeated experiments (mean 99%).

Specificity

The kit is based on a matched pair of mAbs specific for human ApoB100. The antibodies cross-react with apoB from non-human primates.



The products are for research use only.

Novus Biologicals shall not be liable for the use or handling of the product or for consequential, special, indirect or incidental damages there from.
