



HUMAN APOLIPOPROTEIN E DETECTION KITS

PROTOCOL

Part # 64APOEPEG & 64APOEPEH

Test size#: 500 tests (64APOEPEG) and 10,000 tests (64APOEPEH) - assay volume: 20 μ L

Revision: 01 oct-21

Store at: -60°C or below (64APOEPEG); -60°C or below (64APOEPEH)

For research use only. Not for use in diagnostic procedures.

ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of human Apolipoprotein E (Apo-E) in cell-based formats and offers a fast alternative to ELISA.

The detection principle of this kit is based on HTRF® technology (Homogeneous Time-Resolved Fluorescence). As shown in Figure 1, Human Apolipoprotein E is detected in a sandwich assay by using anti Human Apolipoprotein E antibody labeled with Europium cryptate (donor), and anti Human Apolipoprotein E antibody labeled with d2 (acceptor).

When the dyes are in close proximity, the excitation of the donor with a light source (laser or flash lamp) triggers a Fluorescence Resonance Energy Transfer (FRET) towards the acceptor, which in turn fluoresces at a specific wavelength (665 nm). Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the Human Apolipoprotein E concentration.

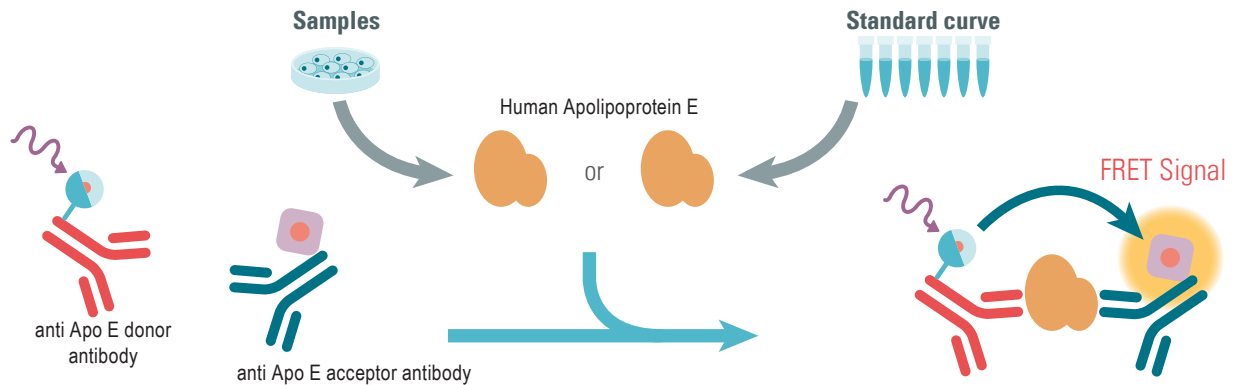
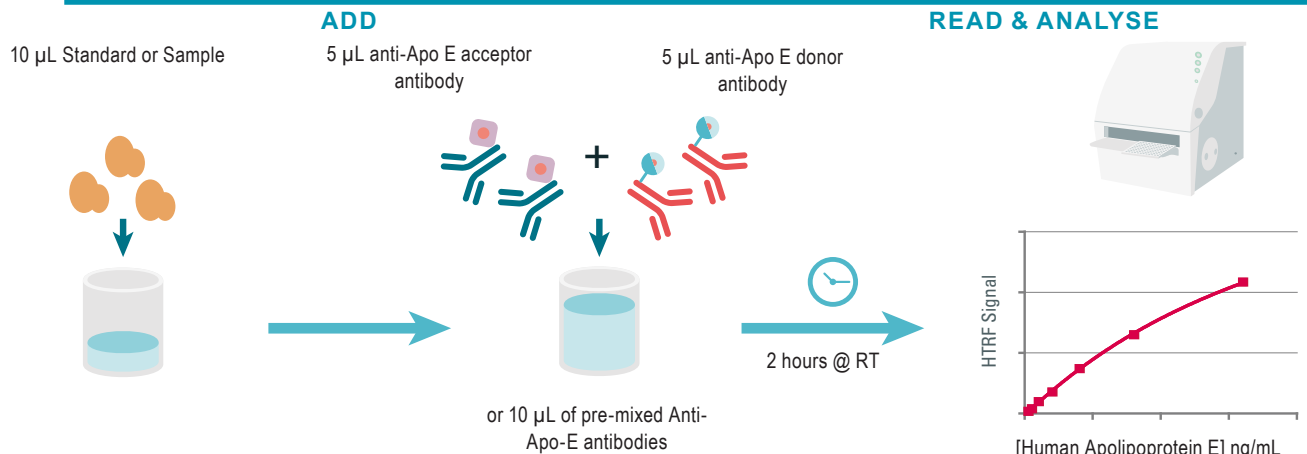


Figure 1: Principle of HTRF Human Apolipoprotein E sandwich assay.

PROTOCOL AT A GLANCE



Make sure to use the set-up for Eu Cryptate.

MATERIALS PROVIDED:

KIT COMPONENTS	500 TESTS * CAT # 64APOEPEG	10,000 TESTS * CAT # 64APOEPEH
Human Apolipoprotein E Standard Frozen	1 vial - 50 μ L 250 ng/mL	1 vial - 50 μ L 250 ng/mL
Human Apolipoprotein E Eu Cryptate Antibody	1 vial - 50 μ L Frozen - 50X	1 vial - 1 mL Frozen - 50X
Human Apolipoprotein E d2 Antibody	1 vial - 50 μ L Frozen - 50X	1 vial - 1 mL Frozen - 50X
Diluent #5 ** 5X	2 vials 2 mL	1 vial 10 mL
Detection buffer *** ready-to-use	1 vial 7 mL Detection Buffer #3	1 vial 105 mL Detection Buffer #3

* When used as advised, the two available kit sizes will provide sufficient reagents for 500 tests and 10,000 tests respectively in 20 μ L final volume..

Assay volumes can be adjusted proportionally to run the assay in 96 or 1536 well microplates.

** Medium like cell culture medium can be an alternative to the diluent.

*** The Detection buffer is used to prepare working solutions of acceptor and donor reagents.

PURCHASE SEPARATELY:

- HTRF[®]-Certified Reader. **Make sure the setup for Eu Cryptate is used.**

For a list of HTRF-compatible readers and set-up recommendations, please visit www.cisbio.com/compatible-readers

- Small volume (SV) detection microplates. For information about microplate recommendations, please visit our website at: cisbio.com/microplates-recommendations

STORAGE AND STABILITY

Store the kit at -60°C or below.

Under proper storage conditions, reagents are stable until the expiry date indicated on the label. Diluent and detection buffer are shipped frozen, but can be stored at $2-8^{\circ}\text{C}$ in your premises.



Reagents

If lyophilized, reconstituted reagents, antibodies, and standard stock solutions may be frozen and thawed only once. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below .

Volume of ApoE standard aliquots should not be under 10 μ L.




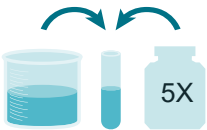
Thawed diluent and detection buffer can be stored at $2-8^{\circ}\text{C}$ in your premises.

REAGENT PREPARATION**BEFORE YOU BEGIN:**

- It is very important to prepare reagents in the specified buffers. The use of an incorrect diluent may affect reagent stability and assay results.
- Thaw the frozen reagents at room temperature, allow them to warm up to room temperature for at least 30 mins before use
- Before use, allow Diluent and Detection buffer to warm up at room temperature and homogenize them with a vortex.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.
- Human Apolipoprotein E standards (for standard curve) must be prepared in diluent or in the same medium as the samples.

TAKE CARE TO PREPARE STOCK AND WORKING SOLUTIONS ACCORDING TO THE DIRECTIONS FOR THE KIT SIZE YOU HAVE PURCHASED.

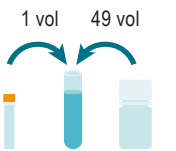
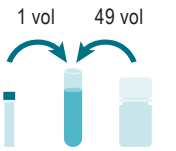

TO PREPARE REAGENT STOCK SOLUTIONS:

500 TESTS KIT - 64APOEPEG			10,000 TESTS KIT - 64APOEPEH
Anti-Human Apolipoprotein E Eu Cryptate antibody			
Thaw the Human Apolipoprotein E Eu Cryptate antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below..			Thaw the Human Apolipoprotein E Eu Cryptate antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below..
Anti-Human Apolipoprotein E d2 antibody			
Thaw the Human Apolipoprotein E d2 antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below..			Thaw the Human Apolipoprotein E d2 antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -60°C or below..
Human Apolipoprotein E Standard			
Thaw the ApoE standard stock solution (5,000 ng/mL) at RT. Mix gently. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solution into disposable plastic vials for storage at -60°C or below.			Thaw the ApoE standard stock solution (5,000 ng/mL) at RT. Mix gently. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solution into disposable plastic vials for storage at -60°C or below.
Diluent			
Dilute 5-fold the 5 X diluent #5 with distilled water: homogenize the 5 X diluent #5 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water (e.g., 1 mL of diluent + 4 mL of distilled water). Mix gently after dilution. This 1X diluent can be frozen and stored at -60°C or below.	4 vol 	1 vol	Dilute 5-fold the 5 X diluent #5 with distilled water: homogenize the 5 X diluent #5 with a vortex and add 1 volume of stock solution in 4 volumes of distilled water (e.g., 10 mL of diluent + 40 mL of distilled water). Mix gently after dilution. This 1X diluent can be frozen and stored at -60°C or below.
Detection buffer			
The Detection buffer is ready-to-use.			The Detection buffer is ready-to-use.

TO PREPARE ANTIBODY WORKING SOLUTIONS:

Each well requires 5 µL of Human Apolipoprotein E-Eu Cryptate Antibody and 5 µL of Human Apolipoprotein E-d2 Antibody.

Prepare the two antibody solutions in separate vials.

500 TESTS KIT - 64APOEPEG			10,000 TESTS KIT - 64APOEPEH
Human Apolipoprotein E Eu Cryptate antibody			
Dilute 50-fold the 50X stock solution (thawed reagent) of human Apo-E Eu Cryptate antibody stock solution with the Detection buffer #3 : add 1 volume of Eu Cryptate antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 0.05 mL of Eu Cryptate antibody stock solution + 2.45 mL of Detection Buffer #3).			Dilute 50-fold the 50X stock solution (thawed reagent) of human Apo-E Eu cryptate antibody stock solution with the Detection buffer #3 : add 1 volume of Eu Cryptate antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 1 mL of Eu Cryptate antibody stock solution + 49 mL of Detection Buffer #3).
Human Apolipoprotein E d2 antibody			
Dilute 50-fold the 50X stock solution (thawed reagent) of human Apo E-d2 antibody stock solution with the Detection buffer #3 : add 1 volume of d2-antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 0.05 mL of d2-antibody stock solution + 2.45 mL of Detection Buffer #3).			Dilute 50-fold the 50X stock solution (thawed reagent) of human Apo E d2 antibody stock solution with the Detection buffer #3 : add 1 volume of d2 antibody stock solution in 49 volumes of Detection buffer #3 (e.g., 1 mL of d2 antibody stock solution + 49 mL of Detection Buffer #3).
Antibody mix			
It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2 antibody solution to 1 volume of Cryptate antibody solution (e.g. 1 mL of d2 antibody + 1 mL of Cryptate antibody).			It is possible to pre-mix the two ready-to-use antibody solutions just prior to dispensing the reagents by adding 1 volume of d2 antibody solution to 1 volume of Cryptate antibody solution (e.g. 1 mL of d2 antibody + 1 mL of Cryptate antibody).

TO PREPARE STANDARD WORKING SOLUTIONS:

- Each well requires 10 μL of standard.
- Dilute the standard stock solution serially with diluent #5 (1X) or in the medium used for the preparation of the samples.
- If culture medium is used to dilute the standard, we recommend to supplement it with serum (2 to 10%) or BSA (0.2 to 1%) in order to avoid ApoE sticking to assay plates.
- In order to check for a potential interference effect from your own assay buffer when using the assay for the first time, we highly recommend the parallel preparation of a standard curve in your own supplemented cell culture medium and in diluent #5 (1X).
- In order to counteract any standard sticking, we recommend changing tips between each dilution.

A recommended standard dilution procedure is listed and illustrated below:

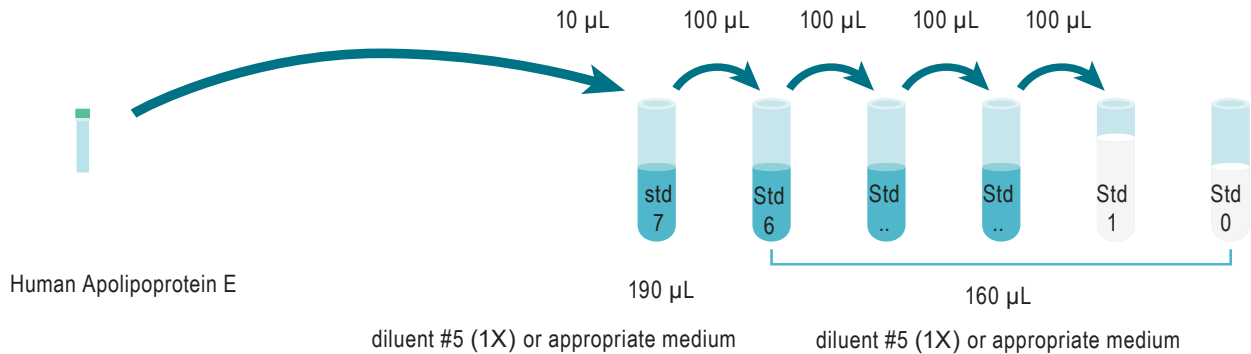
Dilute the standard stock solution 20-fold with diluent; this yields the Standard Max solution (250 ng/mL)

Dilute the standard stock solution 20-fold with diluent #5 (1X) to prepare high standard (std 7): e.g. take 10 μL of standard stock solution and add it to 190 μL of diluent #5 (1X). Mix gently.

Use the high standard (std 7) to prepare the standard curve using 1/2.6 serial dilutions as follows:

- Dispense 160 μL of diluent #5 (1X) in each vial from Std 6 to Std 0.
- Add 100 μL of standard to 160 μL of diluent #5 (1X), mix gently and repeat the 1/2.6 serial dilution to make standard solutions: std6, std5, std4, std3, std2, std1.

This will create 7 standards for the analyte. Std 0 (Negative control) is diluent #5 (1X) or appropriate culture medium alone.








STANDARD	SERIAL DILUTIONS	APOE WORKING SOLUTIONS (NG/ML)
Standard Stock solution	Thawed stock solution	5 000
Standard 7	10 μL Standard Stock Solution + 190 μL Diluent #5 (1X)	250
Standard 6	100 μL standard 7 + 160 μL Diluent #5 (1X)	96,15
Standard 5	100 μL standard 6 + 160 μL Diluent #5 (1X)	36,98
Standard 4	100 μL standard 5 + 160 μL Diluent #5 (1X)	14,22
Standard 3	100 μL standard 4 + 160 μL Diluent #5 (1X)	5,47
Standard 2	100 μL standard 3 + 160 μL Diluent #5 (1X)	2,10
Standard 1	100 μL standard 2 + 160 μL Diluent #5 (1X)	0,81
Standard 0	160 μL Diluent #5 (1X)	-

TO PREPARE SAMPLES:

- Each well requires 10 μ L of sample.
- Just after their collection, put the samples at 4°C and test them immediately. For later use, samples should be dispensed into disposable plastic vials and stored at -60°C or below. Avoid multiple freeze/thaw cycles.
- Samples with a concentration above the highest standard (std 7) must be diluted diluent #5 (1X) in your appropriate sample medium, prepared, as recommended above.
- To obtain additional information or support, please contact the HTRF technical support team at cisbio.com/contact-us

ASSAY PROTOCOL

	Standard (Std 0 - std 7)	Samples
Step 1 	Dispense 10 μ L of each Human Apolipoprotein E standard (Std 0 - std 7) into each standard well	Dispense 10 μ L of each sample into each sample well
Step 2 	Add 5 μ L of Human Apolipoprotein E d2 antibody working solution to all wells	
Step 3 	Add 5 μ L of Human Apolipoprotein E Eu Cryptate antibody working solution to all wells	
Step 4 	Seal the plate and incubate 2 hours @ RT Following incubation, the signal remains stable over a period of 48 hours.	
Step 5 	Remove the plate sealer and read on an HTRF® compatible reader	

DATA REDUCTION

1. Calculate the ratio of the acceptor and donor emission signals for each individual well.

$$\text{Ratio} = \frac{\text{Signal 665 nm}}{\text{Signal 620 nm}} \times 10^4$$

2. Calculate the % CVs. The mean and standard deviation can then be worked out from ratio replicates.

$$\text{CV (\%)} = \frac{\text{Standard deviation}}{\text{Mean Ratio}} \times 100$$

3. Calculate the % delta F which reflects the signal to background of the assay. The negative control (Standard 0) plays the role of an internal assay control. Delta F is used for the comparison of day to day runs of the same assay.

$$\text{delta F (\%)} = \frac{\text{Ratio Standard or sample} - \text{Ratio Negative Control}}{\text{Ratio Negative Control}} \times 100$$

For more information about data reduction, please visit <http://www.cisbio.com/htrf-ratio-and-data-reduction>

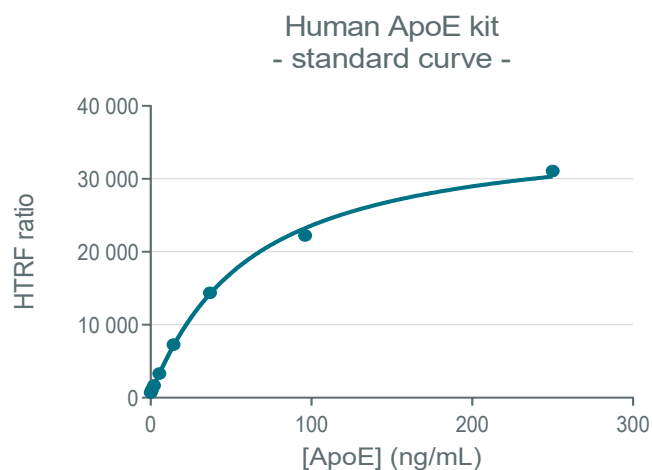
RESULTS

This data must not be substituted for the data obtained in the laboratory and should be considered only as an example.

Results may vary from one HTRF® compatible reader to another.

Standard curve fitting with the 4 Parameter Logistic (4PL) model (with 1/Y² weighting):

	Ratio (1)	CV (2)
Standard 0 - Negative control	732	0%
Standard 1 - 0.81 ng/mL	1089	4%
Standard 2 - 2.10 ng/mL	1696	8%
Standard 3 - 5.47 ng/mL	3304	5%
Standard 4 - 14.22 ng/mL	7298	10%
Standard 5 - 36.98 ng/mL	14383	7%
Standard 6 - 96.15 ng/mL	22239	3%
Standard 7 - 250 ng/mL	31093	2%



ANALYTICAL CHARACTERISTICS

ASSAY PERFORMANCES

Assay range (LOQ* to Std max)	0.32 - 250 ng/mL
Limit Of Detection (LOD)* = Mean Std 0 + 2 SD	0.17 ng/ml
Incubation time	2h at RT

*The LOD and LOQ were calculated from data obtained in diluent with the PHERAstar FS reader (flash lamp excitation) after overnight incubation. These values may vary from one HTRF compatible reader to another.

This product contains material of biologic origin. Use for research purposes only. Do not use in humans or for diagnostic purposes. The purchaser assumes all risk and responsibility concerning reception, handling and storage. The use of the cell line will be done with appropriate safety and handling precautions to minimize health and environmental impact. Remaining disclaimer.

Copyright 2020 Cisbio Bioassays. All rights reserved. HTRF and the HTRF logo are trademarks or registered trademarks of Cisbio Bioassays.

FOR MORE INFORMATION

Europe and other countries +33(0)466-796-705 U.S. and Canada 1-888-963-4567 China +86-21-5018-9880

Japan +81-(0)43-306-8712 Visit www.cisbio.com to find a list of our regional distributors