

MANUAL

Technology: HTRF®

Biomarkers

HTRF Human Apolipoprotein A1 kit

Part number:	64APAPEG	64APAPEH
Test size	500 tests	10,000 tests

Storage: $\leq 60^{\circ}\text{C}$

Version: 05

Date: January 2024

ASSAY PRINCIPLE

This assay is intended for the quantitative determination of Human Apolipoprotein A1 (Apo-A1) using HTRF® technology.

As shown in the diagram to the right, Human Apolipoprotein A1 is detected in a sandwich assay format using two different specific antibodies, one labeled with Eu³⁺-Cryptate (donor) and the second with d2 (acceptor).

When these 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 (665nm).

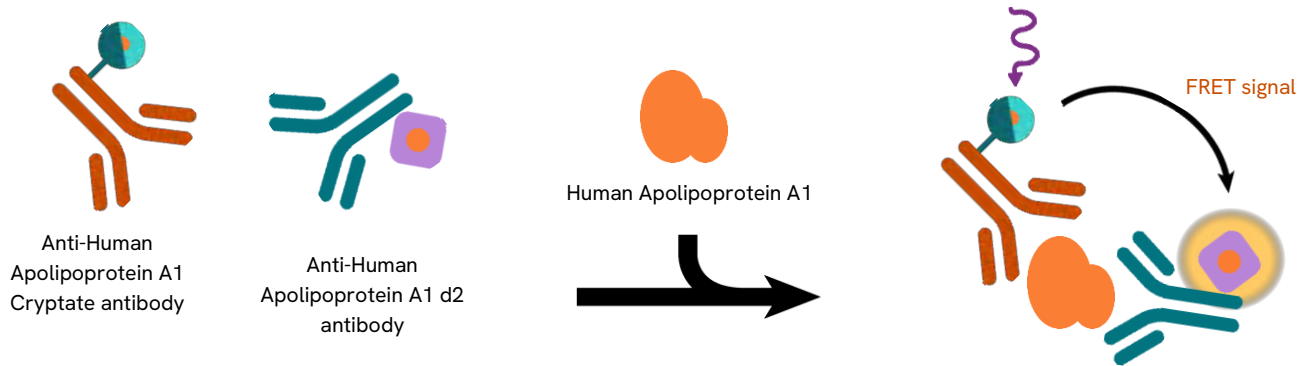
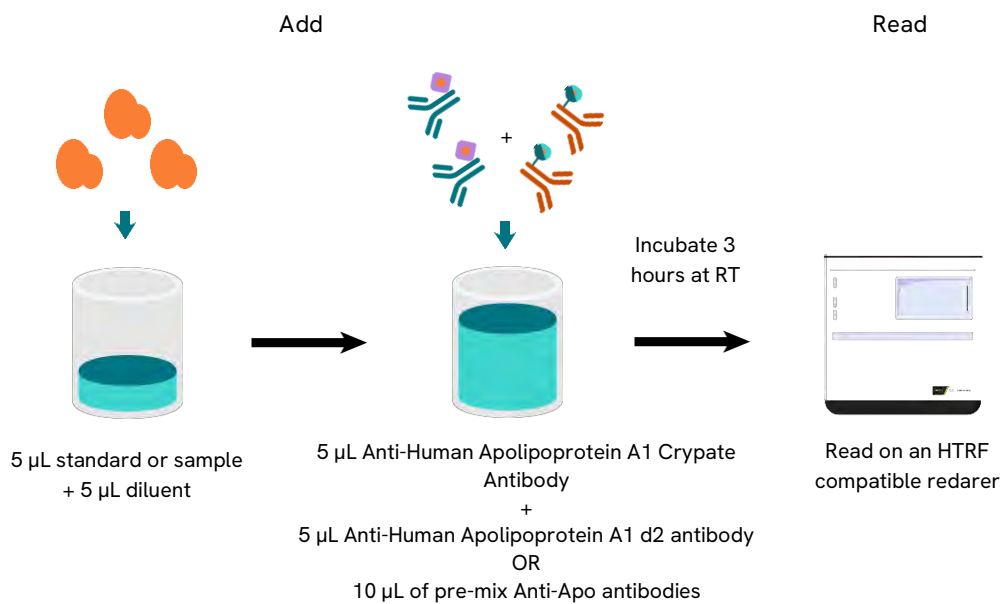


Figure 1: Principle of HTRF sandwich assay











The two antibodies bind to the antigen present in the sample, thereby generating FRET. Signal intensity is proportional to the number of antigen-antibody complexes formed and therefore to the Human Apolipoprotein A1 concentration.

PROTOCOL AT A GLANCE



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

MATERIAL PROVIDED

KIT COMPONENTS		Human Apo-A1 Standard		Anti-Human Apo-A1 d2 antibody		Anti-Human Apo-A1 Eu ³⁺ -Cryptate-antibody		Diluent		Detection buffer #3	
500 TESTS	Tubes		Green cap		Blue cap		Red cap		White cap		Transparent cap
	Stock solution	200 µL/vial 10 µg/mL		50 µL/vial		50 µL/vial		20 mL/vial		7 mL/vial 2 vials	
10,000 TESTS	Tubes		Green cap		Purple cap		Red cap		White cap		Red cap
	Stock solution	200 µL/vial 10 µg/mL		1 mL/vial		1 mL/vial		20 mL/vial		105 mL/vial 2 vials	
STORAGE		≤ 60°C		≤ 60°C		≤ 60°C		2-8°C to -20°C*		2-8°C to -20°C*	
CAUTION				Filter the Cryptate-antibody stock solution before use				Apolipoproteins are sticky components, we recommend to dilute them in the diluent provided in the kit, or in your own diluent containing over 0.5% BSA, or in cell culture medium supplemented with at least 2% FCS		Before use, we recommend to supplement the detection buffer with 0.02% triton X100.	
				Antibody stock solutions must be frozen in liquid nitrogen							
Use black plates only											

* Diluent and Detection buffer are shipped frozen, but can be stored at 2-8°C.

The Apo-A1 standard is prepared from human plasma. This was found to be negative for HBsAg, HIV-1, HIV-2, HIV-3, and HCV. As no test guarantees a product to be non-infectious, it is recommended to handle Apo-A1 solutions with the same precautions as potentially infectious specimens.

REAGENT PREPARATION

HTRF[®] reagent concentrations have been set for optimal assay performance. Any dilution or improper use of the d2 and Cryptate-antibodies will impair the quality of the assay.

For an accurate quantitative determination of sample, dilution must be carried out with the medium used for preparing the samples (i.e. diluent, culture medium or any other compatible medium).

Antibodies may be frozen and thawed once: to avoid freeze/thaw cycles it is recommended that you dispense remaining stock solutions of antibodies into disposable plastic vials. **Antibody stock solutions must be frozen in liquid nitrogen and stored at -60°C or below.**

Please note, working solution preparation may differ between the 500 and the 10,000 test size kits.

- Thaw all reagents at room temperature, allow them to warm up.
- Prepare the working solutions from stock solutions by following the instructions below.

One-plate assay protocol: cell-based

Dispense the reagents in the following order:



* Optimization of cell seeding densities is recommended.

** Time course study is recommended to determine the optimal stimulation time.

***The two working antibody solutions must be prepared in individual vials and can be mixed prior to dispense: add 1 volume of anti-Human Apolipoprotein A1-d2 antibody to 1 volume of anti-Human Apolipoprotein A1-Eu³⁺-cryptate antibody, and dispense 10 µL of the pre-mix anti-Human Apolipoprotein A1 antibodies.

- Cover the plate with a plate sealer.
- **Incubate at RT for 3 hours.**
- Remove the plate sealer and,
- Read the fluorescence emission at two different wavelengths (665nm and 620nm) on a compatible HTRF® reader.

For HTRF certified reader

For more information about HTRF® compatible readers and for set-up recommendations, please visit our website.

	Assays controls			
	Negative control	Cryptate control	Buffer control	Cells / Std
	<i>used to calculate the delta F%</i>	<i>used to check the Cryptate signal at 620 nm</i>	<i>used to check background fluorescence</i>	
Cells / Std	-	-	-	5 µL
Diluent	10 µL	10 µL	10 µL	5 µL
Anti-Human Apolipoprotein A1-d2 antibody	5 µL	-	-	5 µL
Anti-Human Apolipoprotein A1-Eu ³⁺ -Cryptate antibody	5 µL	5 µL	-	5 µL
Detection buffer#3	-	5 µL	10 µL	-

Preparation of antibody working solutions

Determine the amount of antibody needed for the experiment. Each well requires 5 µL of each antibody.

500 TESTS KIT	10,000 TESTS KIT
Anti-Human Apolipoprotein A1-d2 antibody	
Dilute 50-fold the stock solution of anti-human Apo-A1-d2 antibody with detection buffer: e.g. take 0.05 mL of d2-antibody stock solution and add it to 2.45 mL of detection buffer #3.	Dilute 50-fold the stock solution of anti-human Apo-A1-d2 antibody with detection buffer: e.g. take 1 mL of d2-antibody stock solution and add it to 49 mL of detection buffer #3.
Anti-Human Apolipoprotein A1-Eu³⁺-Cryptate-antibody	
Dilute 50-fold the stock solution of anti-human Apo-A1-cryptate antibody with detection buffer e.g. take 0.05 mL of cryptate-antibody stock solution and add it to 2.45 mL of detection buffer #3.	Dilute 50-fold the stock solution of anti-human Apo-A1-cryptate antibody with detection buffer e.g. take 1 mL of cryptate-antibody stock solution and add it to 49 mL of detection buffer #3.

Standard curve preparation

Determine how many samples and replicates will be tested.

Each well requires 5 µL of sample or standard.

Please note: If the sample to test is a cell supernatant, replace the diluent by culture medium. As apolipoproteins are sticky components, we recommend to dilute them using buffers containing over 0.5%BSA or cell culture medium supplemented with at least 2% FCS.

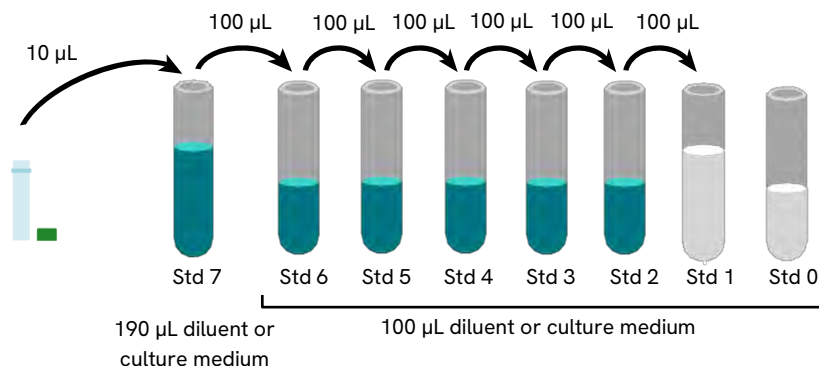
STANDARD	PREPARATION	WORKING CONCENTRATION (ng/mL)
Standard 7	200 µL standard stock solution	500
Standard 6	100 µL Std 7 + 100 µL diluent or culture medium	250
Standard 5	100 µL Std 6 + 100 µL diluent or culture medium	125
Standard 4	100 µL Std 5 + 100 µL diluent or culture medium	62.5
Standard 3	100 µL Std 4 + 100 µL diluent or culture medium	31.25
Standard 2	100 µL Std 3 + 100 µL diluent or culture medium	15.6
Standard 1	100 µL Std 2 + 100 µL diluent or culture medium	7.8
Standard 0	100 µL diluent or culture medium	0

A recommended standard dilution procedure is listed and illustrated below:

- 1) Pre-dilute the standard stock solution 20-fold with diluent. In practice: take 10 μ L of stock solution and add 190 μ L of diluent or culture medium in the standard stock solution. Mix gently. This yields the high standard (Std 7: 500 ng/mL) for the top of the curve.
- 2) Prepare the following serial dilutions:
 - Use the high standard (Std 7) to prepare the standard curve using 1/2 serial dilutions as follows:
 - Dispense 100 μ L of diluent in each vial from Std 6 to Std 1.
 - Add 100 μ L of standard to 100 μ L of diluent, mix gently and repeat the 1/2 serial dilution to make standard solutions: 250, 125, 62.5, 31.25, 15.6, 7.8 ng/mL.

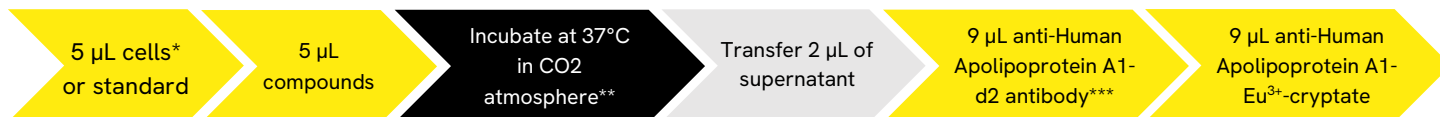
This will create 7 standards for the analyte. Std 0 (Negative control) is diluent or culture medium alone.

The standard dilution procedure is listed and illustrated below.



Two-plate assay protocol: supernatant

Dispense the reagents in the following order:



* Optimization of cell seeding densities is recommended.

** Time course study is recommended to determine the optimal stimulation time.

***The two working antibody solutions must be prepared in individual vials and can be mixed prior to dispense: add 1 volume of anti-Human Apolipoprotein A1-d2 antibody to 1 volume of anti-Human Apolipoprotein A1-Eu³⁺ cryptate antibody, and dispense 18 μ L of the pre-mix anti-Human Apolipoprotein A1 antibodies.

- Cover the plate with a plate sealer.
- **Incubate at RT for 3 hours.**
- Remove the plate sealer and,
- Read the fluorescence emission at two different wavelengths (665nm and 620nm) on a compatible HTRF® reader.

For HTRF certified reader

For more information about HTRF® compatible readers and for set-up recommendations, please visit our website.

	Assays controls			
	Negative control	Cryptate control	Buffer control	Cells / Std
	<i>used to calculate the delta F%</i>	<i>used to check the Cryptate signal at 620 nm</i>	<i>used to check background fluorescence</i>	
Supernatant / Std	-	-	-	2 µL
Diluent	2 µL	2 µL	2 µL	-
Anti-Human Apolipoprotein A1-d2 antibody	9 µL	-	-	9 µL
Anti-Human Apolipoprotein A1-Eu ³⁺ -Cryptate antibody	9 µL	9 µL	-	9 µL
Detection buffer#3	-	9 µL	18 µL	-

Preparation of antibody working solutions

Determine the amount of antibody needed for the experiment. Each well requires 5 µL of each antibody.

500 TESTS KIT	10,000 TESTS KIT
Anti-Human Apolipoprotein A1-d2 antibody	
Dilute 90-fold the stock solution of anti-human Apo-A1-d2 antibody with detection buffer: e.g. take 0.01 mL of d2-antibody stock solution and add it to 0.89 mL of detection buffer #3.	Dilute 90-fold the stock solution of anti-human Apo-A1-d2 antibody with detection buffer: e.g. take 0.01 mL of d2-antibody stock solution and add it to 0.89 mL of detection buffer #3.
Anti-Human Apolipoprotein A1-Eu³⁺-Cryptate-antibody	
Dilute 90-fold the stock solution of anti-human Apo-A1-cryptate antibody with detection buffer e.g. take 0.01 mL of cryptate-antibody stock solution and add it to 0.89 mL of detection buffer #3.	Dilute 90-fold the stock solution of anti-human Apo-A1-cryptate antibody with detection buffer e.g. take 0.01 mL of cryptate-antibody stock solution and add it to 0.89 mL of detection buffer #3.

Standard curve preparation

Determine how many samples and replicates will be tested.

Each well requires 5 μL of sample or standard.

Please note: If the sample to test is a cell supernatant, replace the diluent by culture medium. As apolipoproteins are sticky components, we recommend to dilute them using buffers containing over 0.5%BSA or cell culture medium supplemented with at least 2% FCS.

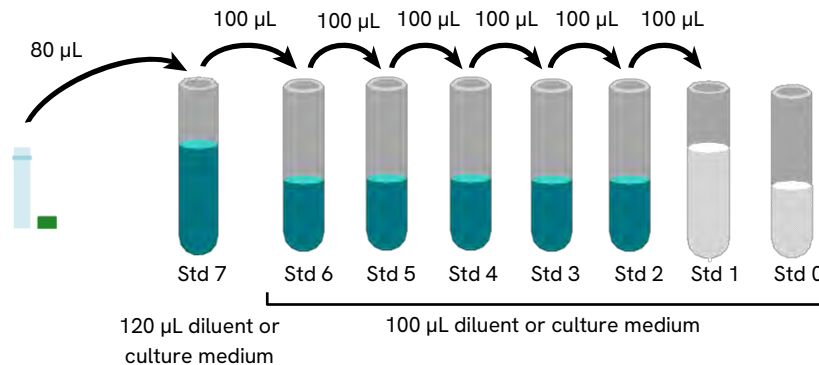
STANDARD	PREPARATION	WORKING CONCENTRATION (ng/mL)
Standard 7	200 μL standard stock solution	4 000
Standard 6	100 μL Std 7 + 100 μL diluent or culture medium	2 000
Standard 5	100 μL Std 6 + 100 μL diluent or culture medium	1 000
Standard 4	100 μL Std 5 + 100 μL diluent or culture medium	500
Standard 3	100 μL Std 4 + 100 μL diluent or culture medium	250
Standard 2	100 μL Std 3 + 100 μL diluent or culture medium	125
Standard 1	100 μL Std 2 + 100 μL diluent or culture medium	62.5
Standard 0	100 μL diluent or culture medium	0

A recommended standard dilution procedure is listed and illustrated below:

- 1) Pre-dilute the standard stock solution 2.5-fold with diluent. In practice: take 40 μL of stock solution and add 60 μL of diluent or culture medium in the standard stock solution. Mix gently. This yields the high standard (Std 7: 4,000 ng/mL) for the top of the curve.
- 2) Prepare the following serial dilutions:
 - Use the high standard (Std 7) to prepare the standard curve using 1/2 serial dilutions as follows:
 - Dispense 100 μL of diluent in each vial from Std 6 to Std 1.
 - Add 100 μL of standard to 100 μL of diluent, mix gently and repeat the 1/2 serial dilution to make standard solutions: 2,000 - 1,000 - 500 - 250 - 125 - 62.5 ng/mL.

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

The standard dilution procedure is listed and illustrated below.



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 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 our website.

RESULTS

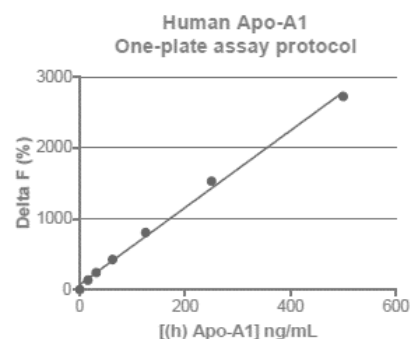
This data must not be substituted for that obtained in the laboratory and should be considered only as an example. Results may vary from one HTRF® compatible reader to another.

Standard curve were prepared in DMEM (GIBCO, low glucose), 2% FBS medium (penicillin, streptomycin).

The assay standard curve is created by plotting delta F% versus the analyte concentration:

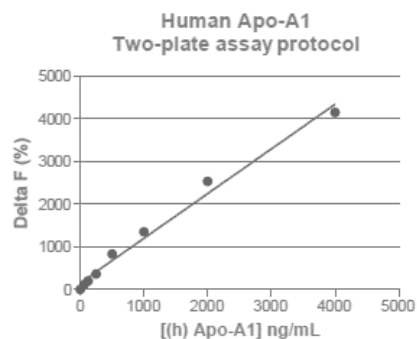
One-plate assay protocol

[Apo-A1 Std] (ng/mL)	Ratio (1)	CV % (2)	Delta F%
0	446	2%	0%
15.6	1 047	1%	135%
31.25	1 527	0%	242%
62.5	2 353	6%	427%
125	4 047	1%	806%
250	7 280	1%	1 531%
500	12 623	3%	2 727%



Two-plate assay protocol

[Apo-A1 Std] (ng/mL)	Ratio (1)	CV % (2)	Delta F%
0	479	3%	0%
62.5	997	3%	108%
125	1 422	4%	197%
250	2 218	3%	363%
500	4 468	7%	833%
1000	6 950	5%	1 351%
2 000	12 596	5%	2 530%
4 000	20 346	7%	4 148%



ANALYTICAL CHARACTERISTICS

Standardization

The HTRF® ApoA1 assay is calibrated against the ApoA1, A2 & B High Level Calibrator(Wako), based on IFCC Reference Preparation SP1-01.

Cross-reactivity

	Cross-reactivity %
Apo-A1	100%
Apo-B100	0.004%
Apo-A2	0.037%
Apo-C1	0.014%
Apo-C2	0.004%
Apo-C3	0.005%
Apo-E	0.004%

Detection limit

One-plate assay protocol: 0.3 ng/mL

Two-plate assay protocol (supernatant): 0.5 ng/mL

Extended range

One-plate assay protocol: 1,000 ng/mL

Two-plate assay protocol (supernatant): 8,000 ng/mL

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.



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

www.revvity.com

revvity

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000
www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.