



HUMAN FIBRONECTIN EDA DETECTION KITS

Part # 64HFNEDAPEG & 64HFNEDAPEH

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

Revision: #02 of Septmeber 2023

Store at: -16°C or below (64HFNEDAPEG); -16°C or below (64HFNEDAPEH)

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

ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of Human Fibronectin EDA in supernatant or cell lysate 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 Fibronectin EDA is detected in a sandwich assay by using anti Human Fibronectin EDA labeled with Europium cryptate (donor), and anti Human Fibronectin EDA 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 Fibronectin EDA concentration.

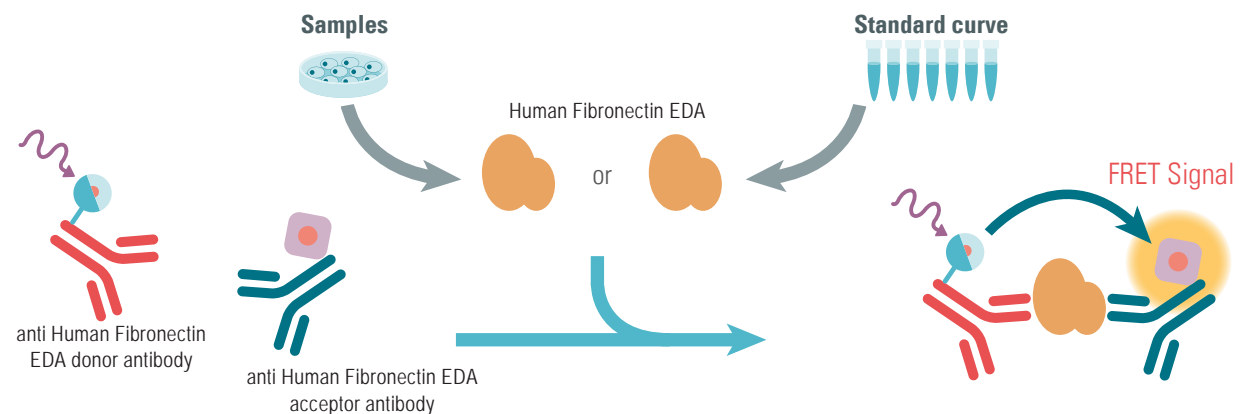
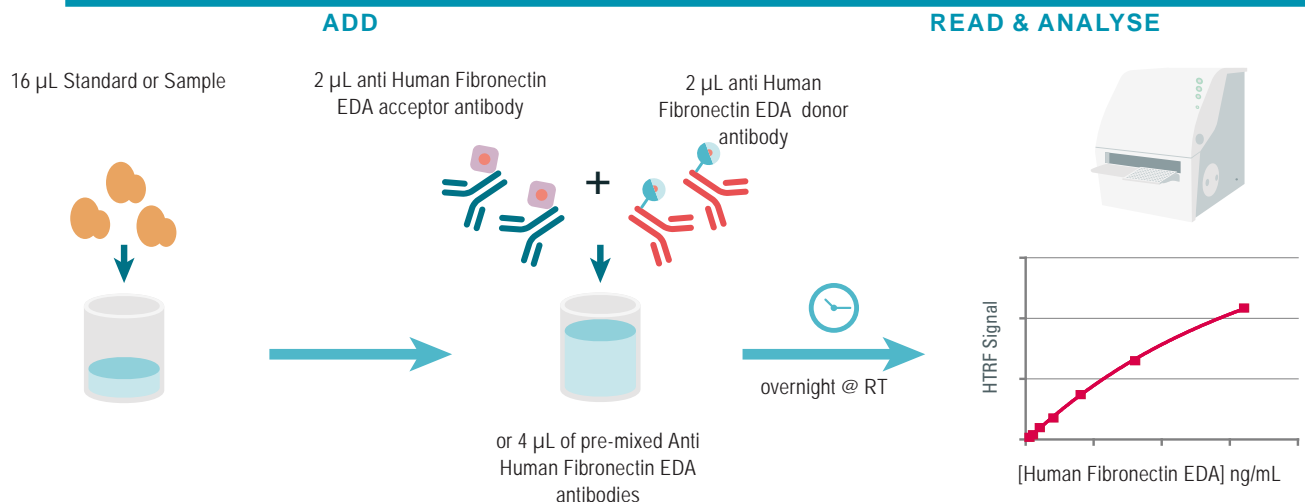


Figure 1: Principle of HTRF Human Fibronectin EDA sandwich assay.

MANUAL AT A GLANCE



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

MATERIALS PROVIDED:

KIT COMPONENTS	500 TESTS * CAT # 64HFNEDAPEG	10,000 TESTS * CAT # 64HFNEDAPEH
Human Fibronectin EDA Standard Lyophilized	1 vial 4,500 ng/mL	2 vials 4,500 ng/mL
Human Fibronectin EDA Eu Cryptate Antibody	1 vial - 50 µL Frozen - 20X	1 vial - 1 mL Frozen - 20X
Human Fibronectin EDA d2 Antibody	1 vial - 50 µL Frozen - 20X	1 vial - 1 mL Frozen - 20X
Lysis buffer #3 ** 4X	1 vial 130 mL	1 vial 130 mL
Detection buffer *** ready-to-use	2 vials 1.5 mL Detection Buffer #3	1 vial 50 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 µL 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.revvy.com

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

STORAGE AND STABILITY

Store the kit at -16°C or below.

Under proper storage conditions, reagents are stable until the expiry date indicated on the label. Lysis buffer and detection buffer are shipped frozen, but can be stored at 2-8°C in your premises.

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 -16°C or below. Volume of Human Fibronectin EDA standard aliquots should not be under 100 µL.







Thawed buffer and detection buffer can be stored at 2-8°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 lysis buffer may affect reagent stability and assay results.
- **Avoid freeze thawing cycles.**
- 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 Lysis buffer and Detection buffer to warm up at room temperature and homogenize them with a vortex.
- It is recommended to filter buffers.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.
- Human Fibronectin EDA standards (for standard curve) must be prepared in lysis buffer 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 - 64HFNEDAPEG			10,000 TESTS KIT - 64HFNEDAPEH
Anti-Human Fibronectin EDA Eu Cryptate antibody			
Thaw the Human Fibronectin EDA Eu Cryptate antibody . Mix gently. This 20X stock solution can be frozen and stored at -16°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -16°C or below.			Thaw the Human Fibronectin EDA Eu Cryptate antibody . Mix gently. This 20X stock solution can be frozen and stored at -16°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -16°C or below.
Anti-Human Fibronectin EDA d2 antibody			
Thaw the Human Fibronectin EDA d2 antibody . Mix gently. This 20X stock solution can be frozen and stored at -16°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -16°C or below.			Thaw the Human Fibronectin EDA d2 antibody . Mix gently. This 20X stock solution can be frozen and stored at -16°C or below. To avoid freeze/thaw cycles, it is recommended to dispense remaining stock solutions into disposable plastic vials for storage at -16°C or below.
Human Fibronectin EDA Standard			
Reconstitute the Human Fibronectin EDA Standard with distilled water in order to obtain a 4,500 ng/mL stock solution. See instructions on vial label for reconstitution volume. Mix gently after reconstitution. This stock solution can be frozen and stored at -16°C or below. After use, the reconstituted standard solution must be frozen and stored at -16°C or below			Reconstitute the Human Fibronectin EDA Standard with distilled water in order to obtain a 4,500 ng/mL stock solution. See instructions on vial label for reconstitution volume. Mix gently after reconstitution. This stock solution can be frozen and stored at -16°C or below. After use, the reconstituted standard solution must be frozen and stored at -16°C or below
Lysis buffer			
Dilute 4-fold the 4 X lysis buffer #3 with distilled water: homogenize the 4 X lysis buffer #3 with a vortex and add 1 volume of stock solution in 3 volumes of distilled water (e.g., 1 mL of lysis buffer + 3 mL of distilled water). Mix gently after dilution. This 1X buffer can be frozen and stored at -60°C or below.	3 vol		1 vol
Detection buffer			
The Detection buffer is ready-to-use.			The Detection buffer is ready-to-use.

TO PREPARE ANTIBODY WORKING SOLUTIONS:

Each well requires 2 µL of Human Fibronectin EDA-Eu Cryptate Antibody and 2 µL of Human Fibronectin EDA-d2 Antibody.

Prepare the two antibody solutions in separate vials.

500 TESTS KIT - 64HFNEDAPEG			10,000 TESTS KIT - 64HFNEDAPEH
Human Fibronectin EDA Eu Cryptate antibody			
Dilute 20-fold the 20X stock solution (thawed reagent) of Human Fibronectin EDA Eu Cryptate antibody with Detection buffer #3: add 1 volume of Eu Cryptate antibody stock solution in 19 volumes of detection buffer (e.g. 50 µL of Eu Cryptate antibody stock solution + 950 µL of detection buffer).	1 vol	19 vol	Dilute 20-fold the 20X stock solution (thawed reagent) of Human Fibronectin EDA Eu Cryptate antibody with Detection buffer #3: add 1 volume of Eu Cryptate antibody stock solution in 19 volumes of detection buffer (e.g. 1 mL of Eu Cryptate antibody stock solution + 19 mL of detection buffer).
Human Fibronectin EDA d2 antibody			
Dilute 20-fold the 20X stock solution (thawed reagent) of Human Fibronectin EDA d2 antibody with Detection buffer #3: add 1 volume of d2 antibody stock solution in 19 volumes of detection buffer (e.g. 50 µL of d2-antibody stock solution + 950 µL of detection buffer).	1 vol	19 vol	Dilute 20-fold the 20X stock solution (thawed reagent) of Human Fibronectin EDA d2 antibody with Detection buffer #3: add 1 volume of d2 antibody stock solution in 19 volumes of detection buffer (e.g. 1 mL of d2 antibody stock solution + 19 mL of detection buffer).
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. 20 mL of d2 antibody + 20 mL of Cryptate antibody).

TO PREPARE STANDARD WORKING SOLUTIONS:

- Each well requires 16 μL of standard.
- Dilute the standard stock solution serially with lysis buffer #3 (1X) . Please note: For all samples to test even if are cell supernatant, it is mandatory to use lysis buffer #3 only as diluent to allow better detection of Human Fibronectin EDA.
- For supernatant, we recommend to use medium supplemented with BSA (0.2 to 1%) instead of FCS, in order to avoid interferences or 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 lysis buffer #3 (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:

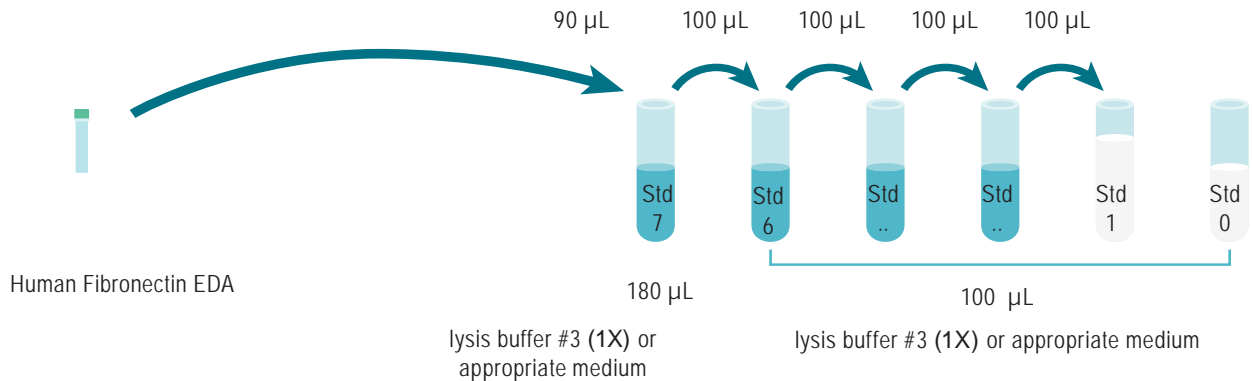
Dilute the standard stock solution 3-fold with diluent; this yields the Standard Max solution (1,500 ng/mL)

Dilute the standard stock solution 3-fold with lysis buffer #3 (1X) to prepare high standard (Std 7): e.g. take 90 μL of standard stock solution and add it to 180 μL of lysis buffer #3 (1X). Mix gently.

Use the high standard (Std 7) to prepare the standard curve using 1/2.0 serial dilutions as follows:

- Dispense 100 μL of lysis buffer #3 (1X) in each vial from Std 6 to Std 0.
- Add 100 μL of standard to 100 μL of lysis buffer #3 (1X), mix gently and repeat the 1/2.0 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 lysis buffer #3 (1X) or appropriate culture medium alone.








STANDARD	SERIAL DILUTIONS	FIBRONECTIN EDA WORKING SOLUTIONS (NG/ML)
Standard Stock solution	Reconstituted lyophilisate	4 500
Standard 7	90 μL stock solution + 180 μL lysis buffer #3 (1X)	1 500
Standard 6	100 μL standard 7 + 100 μL lysis buffer #3 (1X)	750
Standard 5	100 μL standard 6 + 100 μL lysis buffer #3 (1X)	375
Standard 4	100 μL standard 5 + 100 μL lysis buffer #3 (1X)	188
Standard 3	100 μL standard 4 + 100 μL lysis buffer #3 (1X)	94
Standard 2	100 μL standard 3 + 100 μL lysis buffer #3 (1X)	47
Standard 1	100 μL standard 2 + 100 μL lysis buffer #3 (1X)	23
Standard 0	100 μL Lysis buffer #3 (1X)	-

TO PREPARE SAMPLES:

- Each well requires 16 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.
- Cell supernatants must be prepared using a culture medium supplemented with serum (2 to 10%) or BSA (1%) to avoid Human Amyloid β 1-42 sticking to culture vessels. As Human Amyloid β 1-42 is prone to degradation, addition of antiprotease inhibitor cocktail can be beneficial.
- Samples with a concentration above the highest standard (Std 7) must be diluted lysis buffer #3 (1X) necessarily in lysis buffer#3 as diluent, prepared, as recommended above.
- In order to measure Human Fibronectin EDA in cell lysates, cells must be lyzed with Lysis Buffer #3 (1X) for 30 min at RT under gentle shaking. Then cell lysates and cell supernatant must be diluted ysis Buffer #3 (1X) as diluent to allow better detection of HUman Fibronectin EDA. Please note that the 4X stock solution of Lysis Buffer #3 must be ordered separately (Ref# 64KL3FDF, 130 mL) and 4-fold diluted with distilled water before use.
- To obtain additional information or support, please contact the HTRF technical support team at www.revvy.com

ASSAY MANUAL

	Standard (Std 0 - Std 7)	Samples
Step 1 	Dispense 16 μ L of each Human Fibronectin EDA standard (Std 0 - Std 7) into each standard well	Dispense 16 μ L of each sample into each sample well
Step 2 	Add 2 μ L of Human Fibronectin EDA d2 antibody working solution to all wells	
Step 3 	Add 2 μ L of Human Fibronectin EDA Eu Cryptate antibody working solution to all wells	
Step 4 	Seal the plate and incubate overnight @ RT The best signal is achieved after overnight incubation time but it can be read from 4h. 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.

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

For more information about data reduction, please visit www.revvity.com

RESULTS

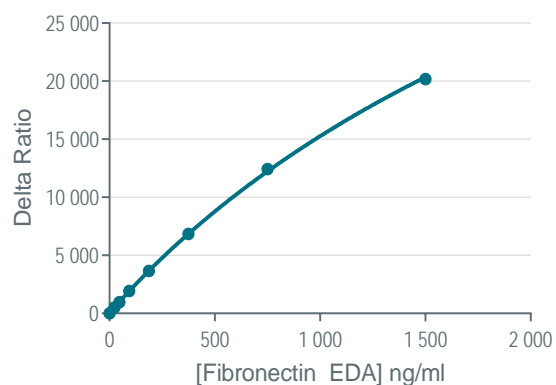
This data must not be substituted for the data obtained in the laboratory and should be considered only as an example.

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	CV	DELTA RATIO
Standard 0 - Negative control	856	4%	0
Standard 1 - 23 ng/ml	1371	3%	515
Standard 2 - 47 ng/ml	1825	3%	969
Standard 3 - 94 ng/ml	2791	3%	1 934
Standard 4 - 188 ng/ml	4505	0%	3 648
Standard 5 - 375 ng/ml	7697	3%	6 841
Standard 6 - 750 ng/ml	13278	1%	12 422
Standard 7 - 1500 ng/ml	21042	2%	20 185

Human Fibronectin EDA Standard Curve



ANALYTICAL CHARACTERISTICS

ASSAY PERFORMANCES

Assay range (LOQ* to Std max)	12 - 1,500 ng/mL
Limit Of Detection (LOD)* = Mean Std 0 + 2 SD	5.2 pg/ml
Incubation time	Overnight at RT (from 4h possible)

*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.



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
www.revvity.com

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