

HTRF (h/m) Ataxin 2 Detection kit

Part # 64ATA2PEG & 64ATA2PEH

Test Size#: 500 tests (64ATA2PEG) and 10,000 tests (64ATA2PEH)

Assay volume: 20 μ L

Revision: #02 of September 2023

Store at: -60°C or below

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

ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of Ataxin 2 in 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, Ataxin 2 is detected in a sandwich assay by using anti Ataxin 2 antibody labeled with Europium cryptate (donor), and anti Ataxin 2 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 Ataxin 2 concentration.

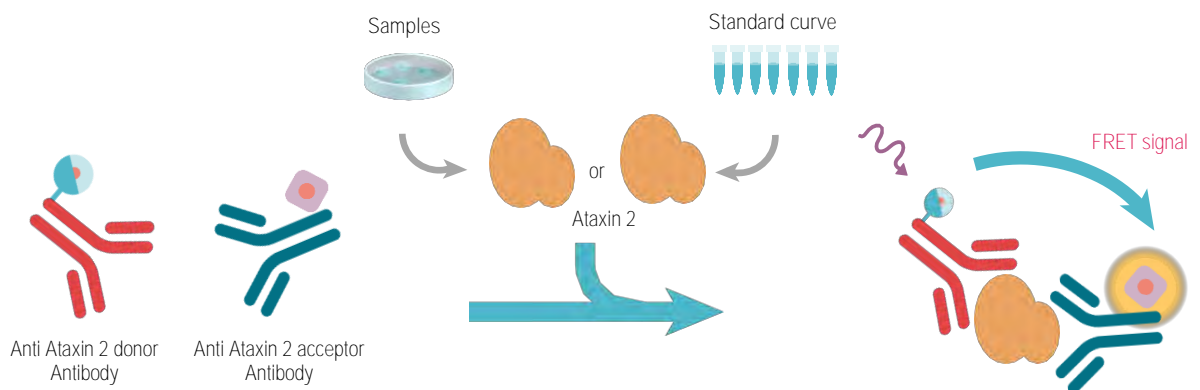
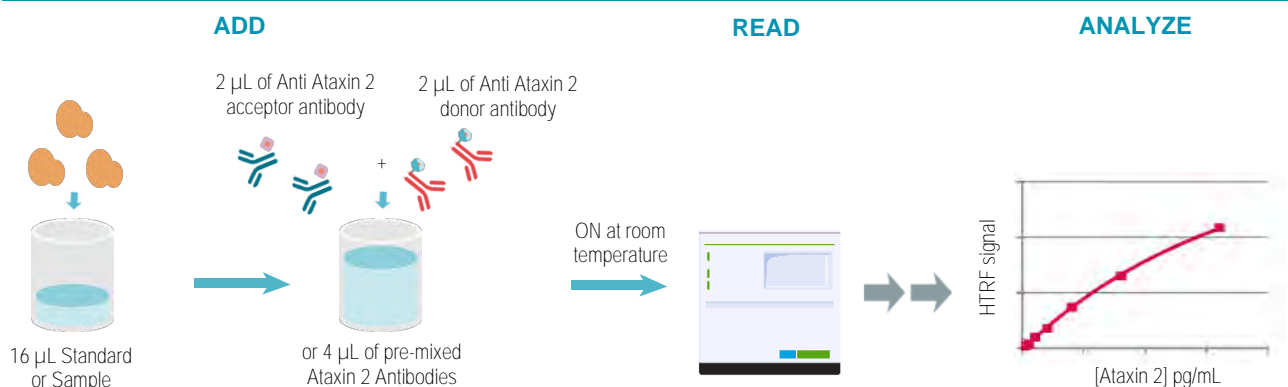


Figure 1. Principle of HTRF Ataxin 2 sandwich assay

MANUAL AT A GLANCE



Make sure to use the set-up for Eu Cryptate

MATERIALS

KIT COMPONENTS	500 TESTS*	10,000 TESTS*
Ataxin 2 Standard Frozen	1 vial – 150 µL 50 ng/mL	1 vial – 150 µL 50 ng/mL
Ataxin 2 Eu Cryptate Antibody Frozen 20X	1 vial 50 µL	1 vial 1 mL
Ataxin 2 d2 Antibody Frozen 20X	1 vial 50 µL	1 vial 1 mL
Lysis Buffer #1 4X	4 vials 2 mL	1 vial 130 mL
Detection Buffer** #3 Ready-to-use	2 vials 1.5 mL	1 vial 50 mL

* 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.

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

STORAGE AND STABILITY

KIT:

- Store the kit at -60°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.

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 -16°C or below.
- Volume of Human Ataxin 2 standard aliquots should not be under 10 µL.

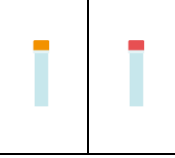







REAGENT PREPARATION

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.
- 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.
- Antibody solutions must be prepared in individual vials and can be mixed prior to dispensing.
- Ataxin 2 standards (for standard curve) must be prepared in lysis buffer.

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

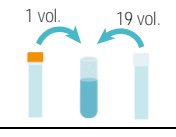
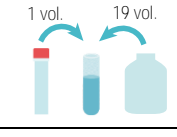
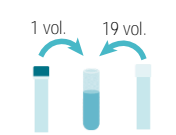



TO PREPARE REAGENT STOCK SOLUTIONS:

500 TESTS		10,000 TESTS	
Anti-Ataxin 2 Eu Cryptate antibody			
Thaw the Ataxin 2 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 Ataxin 2 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-Ataxin 2 d2 antibody			
Thaw the Ataxin 2 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 Ataxin 2 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.
Ataxin 2 Standard			
Thaw the Ataxin 2 standard stock solution (50 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 Ataxin 2 standard stock solution (50 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.
Lysis Buffer			
Dilute 5-fold the 4X Lysis Buffer #1 with distilled water: homogenize the 4X Lysis Buffer #1 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 lysis buffer can be frozen and stored at -16°C or below.			Dilute 5-fold the 4X Lysis Buffer #1 with distilled water: homogenize the 4X Lysis Buffer #1 with a vortex and add 1 volume of stock solution in 3 volumes of distilled water (e.g., 10 mL of lysis buffer + 30 mL of distilled water). Mix gently after dilution. This 1X lysis buffer can be frozen and stored at -16°C or below.
Detection buffer			
The Detection buffer is ready-to-use.			

TO PREPARE WORKING ANTIBODY SOLUTIONS:

Each well requires 2 µL of Ataxin 2-Eu Cryptate Antibody and 2 µL of Ataxin 2-d2 Antibody.

Prepare the two antibody solutions in separate vials

500 TESTS		10,000 TESTS	
Ataxin 2 Eu Cryptate antibody			
Dilute 20-fold the 20X stock solution (thawed reagent) of Ataxin 2 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).			Dilute 20-fold the 20X stock solution (thawed reagent) of Ataxin 2 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).
Ataxin 2 d2 antibody			
Dilute 20-fold the 20X stock solution (thawed reagent) of Ataxin 2 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).			Dilute 20-fold the 20X stock solution (thawed reagent) of Ataxin 2 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 WORKING STANDARDS SOLUTIONS:

- Each well requires 16 μL of standard.
- Dilute the standard stock solution serially with Lysis Buffer #1 (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:

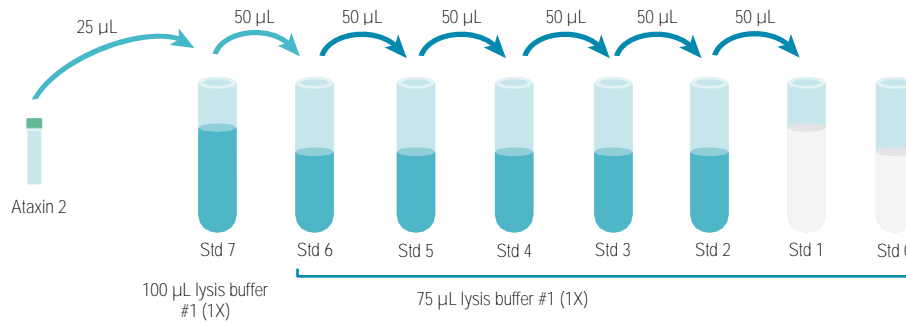
Dilute the standard stock solution 5-fold with lysis buffer; this yields the Standard Max solution (10,000 pg/mL)

Dilute the standard stock solution 5-fold with Lysis Buffer #1 (1X) to prepare high standard (Std 7): e.g. take 25 μL of standard stock solution and add it to 100 μL of Lysis Buffer #1 (1X). Mix gently.

Use the high standard (Std 7) to prepare the standard curve using 1/2.5 serial dilutions as follows:

- Dispense 75 μL of Lysis Buffer #1 (1X) in each vial from Std 6 to Std 0.
- Add 50 μL of standard to 75 μL of Lysis Buffer #1 (1X), mix gently and repeat the 1/2.5 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 #1 (1X) or appropriate culture medium alone



STANDARD	SERIAL DILUTIONS	WORKING SOLUTIONS
Standard Stock solution	Thaw stock solution	50 000 pg/mL
Standard 7	25 μL Standard stock Solution + 100 μL lysis buffer 1X	10 000 pg/mL
Standard 6	50 μL standard 7 + 75 μL lysis buffer 1X	4 000 pg/mL
Standard 5	50 μL standard 6 + 75 μL lysis buffer 1X	1 600 pg/mL
Standard 4	50 μL standard 5 + 75 μL lysis buffer 1X	640 pg/mL
Standard 3	50 μL standard 4 + 75 μL lysis buffer 1X	256 pg/mL
Standard 2	50 μL standard 3 + 75 μL lysis buffer 1X	102 pg/mL
Standard 1	50 μL standard 2 + 75 μL lysis buffer 1X	41 pg/mL
Standard 0	75 μL lysis buffer 1X	0

TO PREPARE SAMPLES

- Each well requires 16 μL 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 Lysis Buffer #1 (1X) in your appropriate sample medium, prepared, as recommended above.
- To obtain additional information or support, please contact the HTRF technical support team at www.revvy.com

DATA REDUCTION & INTERPRETATION

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 delta ratio of the acceptor and donor emission signals for each individual well. The Standard 0 (Negative control) plays the role of an internal assay control.

$$\text{delta Ratio} = \text{Ratio Standard or sample} - \text{Ratio Standard 0}$$

3. 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.revivity.com

RESULTS

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

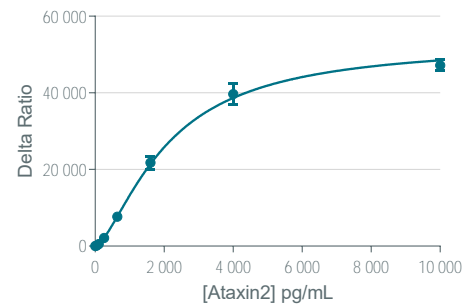
Results may vary from one HTRF® compatible reader to another.

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

(For more information about curve fitting please visit www.revivity.com)

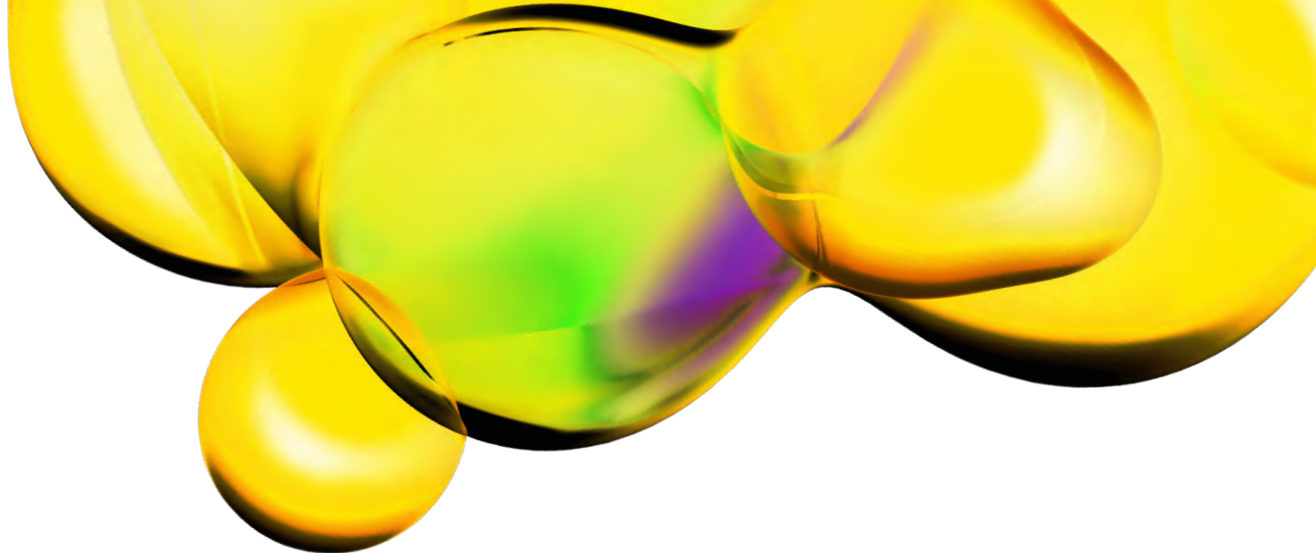
		Ratio (1)	delta R (2)	CV% (3)
Standard 0	Negative control	1272	0	1%
Standard 1	41 pg/mL	1419	147	8%
Standard 2	102 pg/mL	1802	530	14%
Standard 3	256 pg/mL	3420	2148	1%
Standard 4	640 pg/mL	8956	7684	3%
Standard 5	1,600 pg/mL	23025	21753	8%
Standard 6	4,000 pg/mL	40934	39662	7%
Standard 7	10,000 pg/mL	48432	47160	3%

Human/Mouse Ataxin2 Standard Curve in lysis buffer



REACH European regulations and compliance This product and/or some of its components include a Triton concentration of 0.1% or more and as such, it is concerned by the REACH European regulations. We recommend researchers using this product to act in compliance with REACH and in particular: to only use the product for in vitro research in appropriate and controlled premises by qualified researchers, ii) to ensure the collection and the treatment of subsequent waste, and iii) to make sure that the total amount of Triton handled does not exceed 1 ton per year

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
www.revvity.com

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