



## HTRF HUMAN ANDROGEN RECEPTOR DETECTION KIT

Part # 64ANDRPEG & 64ANDRPEH

Test Size#: 500 TESTS (64ANDRPEG), 10,000 TESTS (64ANDRPEH)

Revision: #04 of September 2023 Store at:  $\leq -60^{\circ}\text{C}$

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

### ASSAY PRINCIPLE

This assay is intended for the simple, rapid and direct detection of endogenous levels of human Androgen Receptor in cells. Androgen receptor is produced by cells and after lysis of the cell membrane, Total Androgen Receptor can be detected using the kit reagents. This total protein assay is used to monitor the steady state protein level in various cells.

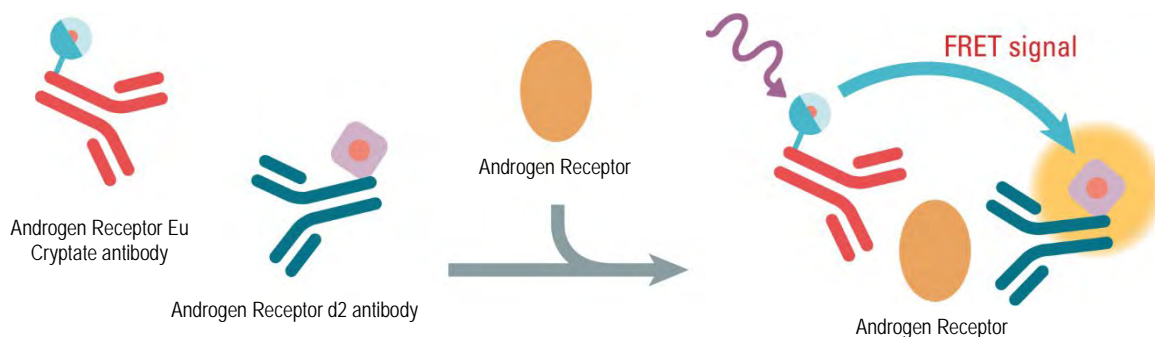


Figure 1: Principle of the HTRF sandwich assay

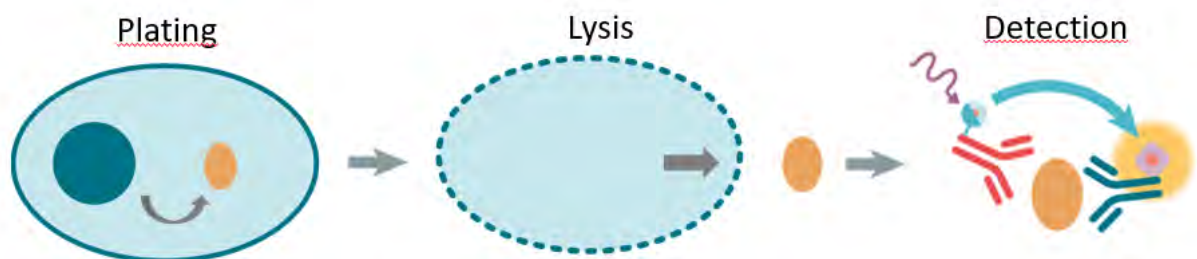
As shown here, Androgen Receptor is detected in a sandwich assay format using 2 different specific antibodies, one labeled with Eu<sup>3+</sup>-Cryptate (donor) and the second with d2 (acceptor).

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). The specific signal modulates positively in proportion to Androgen Receptor expression.

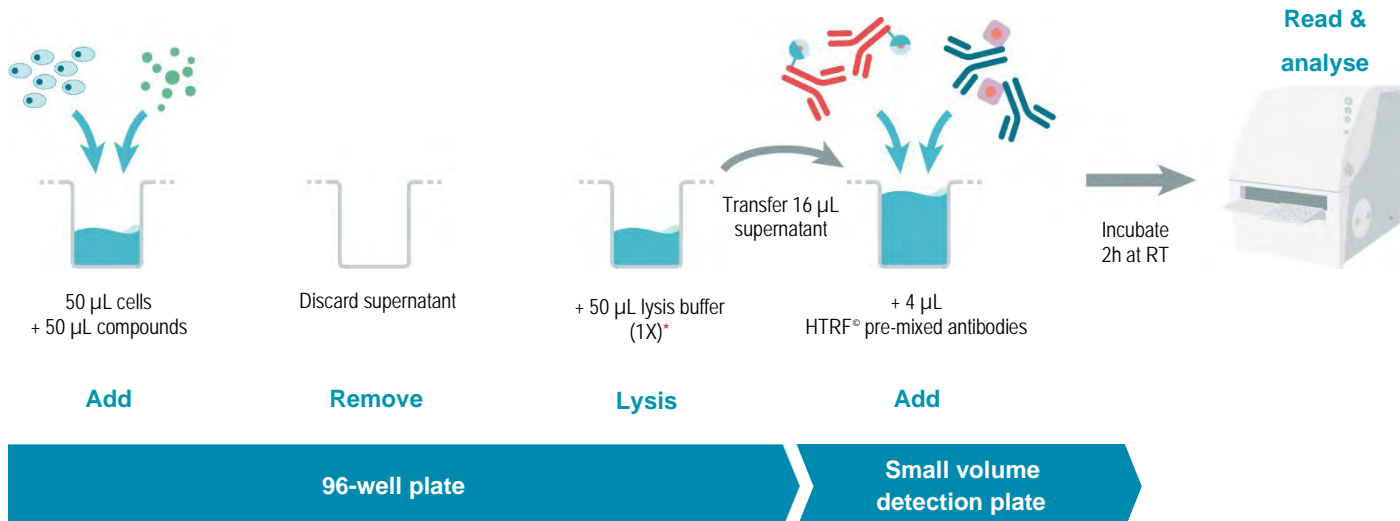
The assay can be run under a two-plate assay manual, where cells are plated and lysed in the same culture plate. Lysates are then transferred to the assay plate for the detection of Androgen Receptor by HTRF<sup>®</sup> reagents. This manual enables the cells viability and confluence to be monitored. It can also be further streamlined to a one-plate assay manual. Detection of Androgen Receptor with HTRF<sup>®</sup> reagents is performed in a single plate used for plating, stimulation and detection. No washing steps are required. This manual, HTS designed, allows miniaturization while maintaining HTRF<sup>®</sup> quality.

For tissue derived samples, please refer to the technical note: "Optimize your HTRF<sup>®</sup> cell signaling assays on tissues" on [www.revvity.com](http://www.revvity.com). Our technical support team can help you to set up this manual, or another. Please contact us at [www.revvity.com](http://www.revvity.com)

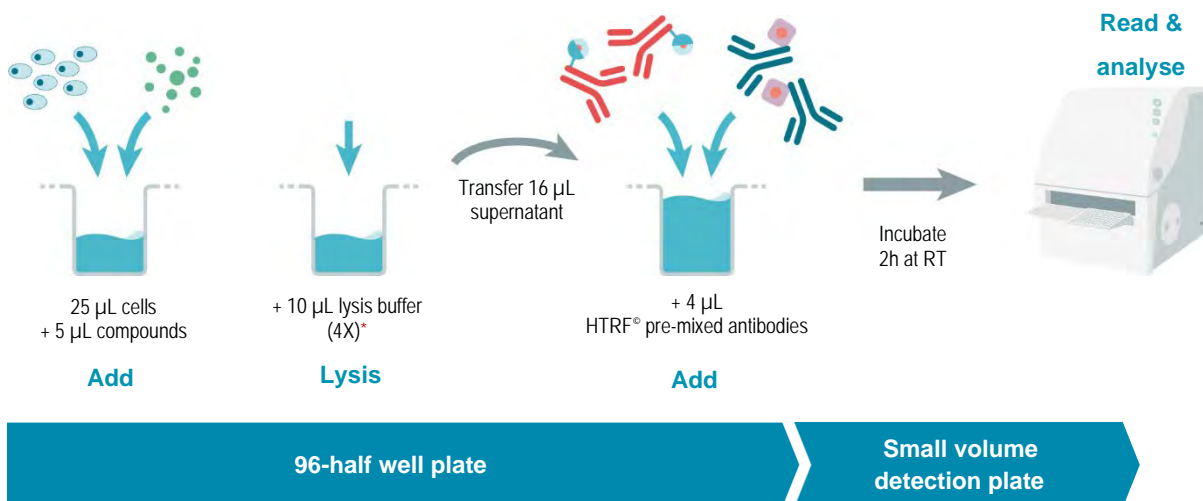
### MANUAL AT A GLANCE



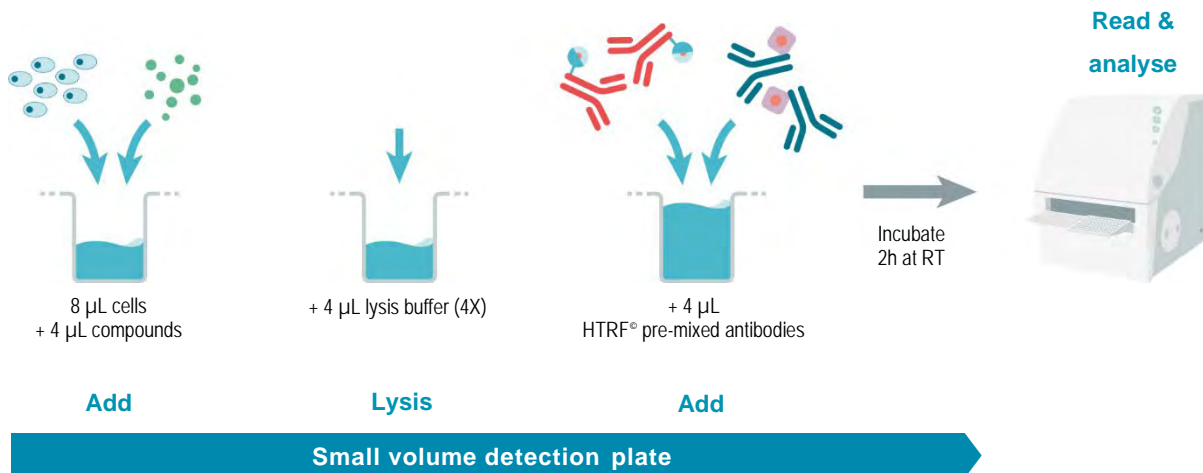
► **TWO-PLATE ASSAY MANUAL FOR ADHERENT CELLS:**



► **TWO-PLATE ASSAY MANUAL FOR SUSPENSION CELLS:**



► **ONE-PLATE ASSAY MANUAL:**















\*Depending on the cell lines used, the lysis volume should be optimized.  
Depending on the cell lines used, it may be necessary to dilute the cell lysate to ensure samples are within the assay linear range.

► **FOR HTRF CERTIFIED READER**

For more information about HTRF<sup>®</sup> compatible readers and for set-up recommendations, please visit our website at: [www.revvy.com](http://www.revvy.com)

**MATERIALS PROVIDED:**

KIT COMPONENTS	STORAGE	500 TESTS CAT# 64ANDRPEG		10,000 TESTS CAT# 64ANDRPEH	
		Green cap	1 vial - 150µL	Green cap	2 vials - 150µl
Androgen Receptor kit - Control lysate (ready-to-use)	≤-60°C	 Green cap	1 vial - 150µL	 Green cap	2 vials - 150µl
Androgen Receptor kit Eu cryptate antibody (20X)	≤-16°C	 Red cap	1 vial - 50µL	 Red cap	1 vial - 1mL
Androgen Receptor kit d2 antibody (20X)	≤-16°C	 Blue cap	1 vial - 50µL	 Blue cap	1 vial - 1mL
Blocking reagent* (stock solution 100X)	≤-16°C	 Purple cap	1 vial - 300µL	 Purple cap	3 vials - 2mL
Lysis buffer* #3 (4X)	≤-16°C	 Clear cap	4 vials - 2mL	 White cap	1 vial - 130mL
Phospho-total protein detection buffer** (ready-to-use)	≤-16°C	 Orange cap	2 vials - 2mL	 Red cap	1 vial - 50mL

\* Amounts of reagents provided are sufficient for generating 50 µL of cell lysate per well.

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

► **PURCHASE SEPARATELY**

96-well or 384-well small volume (SV) detection microplates - For more information about microplate recommendations, please visit our website at: [www.revvity.com](http://www.revvity.com)

## STORAGE AND STABILITY

### Storage upon reception:

Store the kit at -60°C or below until the expiration date indicated on the package.

### Storage and stability of thawed material:

When you are ready to use the kit, take the reagents out and prepare them following the manual provided in this document. Unused thawed reagents can be stored and conserved for future use. Refer to the table below for storage options and corresponding shelf life.

	Storage after Thawing/reconstitution
Lysis Buffer / Blocking Reagent / Detection buffer	2-8°C until the expiration date indicated on the package
Antibodies*	2-8°C for 48h or freeze at -16°C or below until the expiration date indicated on the package for long term storage
Protein/standard /Control Lysate*	freeze at -60°C or below until the expiration date indicated on the package for long term storage

\*For Antibodies, Protein, Standard & control lysate, Stock solutions may be thawed and frozen only once. Freeze in aliquots to avoid multiple freeze/thaw cycles (once aliquoted, single use of the reagent). Volume of antibodies aliquots should not be under 10µL. Volume of Protein, Standard & control lysate aliquots should not be under 20µL.

## REAGENT PREPARATION

Allow all reagents to thaw before use. We recommend centrifuging the vials gently after thawing, before pipetting the stock solutions. Prepare the working solutions from stock solutions by following the instructions below.

### TO PREPARE WORKING CONTROL LYSATE SOLUTION



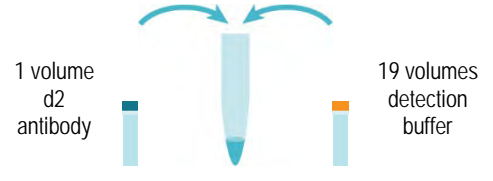
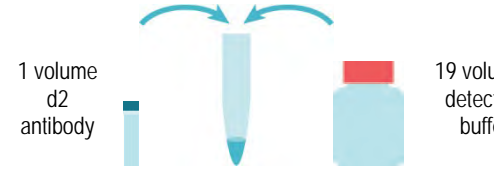

The control lysate is only provided as an internal assay control to check the quality of the results obtained. The window between the control lysate and the negative control should be greater than 2.

Thaw the control lysate. Mix gently, and then the control lysate is ready to use.

### TO PREPARE WORKING ANTIBODY SOLUTIONS:

HTRF® reagent concentrations have been set for optimal assay performances. Note that any dilution or improper use of the d2 or Europium Cryptate-antibodies will impair the assay's quality. Be careful, as the working solution preparation for the antibodies may differ between the 500 and 10,000 test data point kit.

Antibody working solutions are stable for 2 days at 2-8°C. Dilute the antibodies with detection buffer. In practice:

500 TESTS KIT CAT# 64ANDRPEG		10,000 TESTS KIT CAT# 64ANDRPEH	
<i>Androgen Receptor kit Eu cryptate antibody</i>			
 <p>1 volume Eu Cryptate antibody</p> <p>19 volumes detection buffer</p>	 <p>1 volume Eu Cryptate antibody</p> <p>19 volumes detection buffer</p>	<p>Dilute the frozen stock solution with detection buffer 20-fold, e.g. add 0.95 mL of detection buffer to 0.05 mL of Eu Cryptate-antibody stock solution.</p>	
<i>Androgen Receptor kit d2 antibody</i>			
 <p>1 volume d2 antibody</p> <p>19 volumes detection buffer</p>	 <p>1 volume d2 antibody</p> <p>19 volumes detection buffer</p>	<p>Dilute the frozen stock solution with detection buffer 20-fold, e.g. add 0.95 mL of detection buffer to 0.05 mL of d2-antibody stock solution.</p>	
<i>Antibody mix</i>			
<p>It is recommended 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 Eu Cryptate-antibody solution.</p>		<p>It is recommended 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 Eu Cryptate-antibody solution.</p>	

## TO PREPARE LYSIS BUFFER:

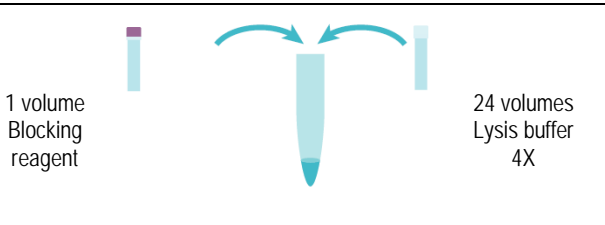
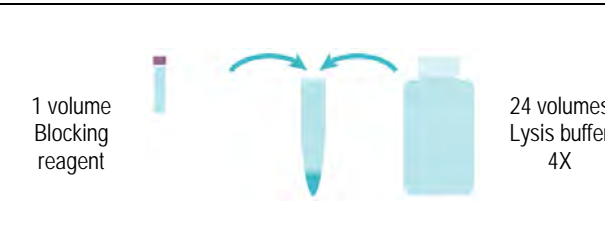
Make sure that the lysate has been generated by using the supplied kit lysis buffer.

Supplemented lysis buffer differs between the manuals. Make sure you use the appropriate supplemented lysis buffer depending on the chosen manual's specifications.

Prepare the required amount of lysis buffer before running the assay. Working solutions are stable for 2 days at 2-8°C.

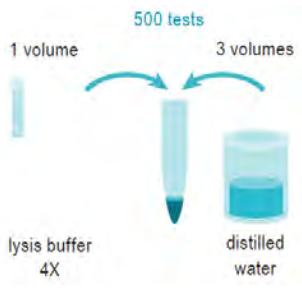
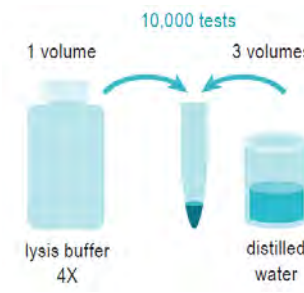
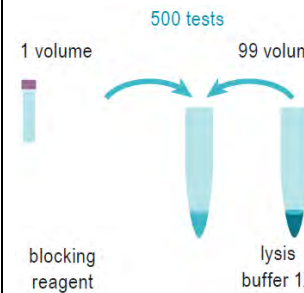
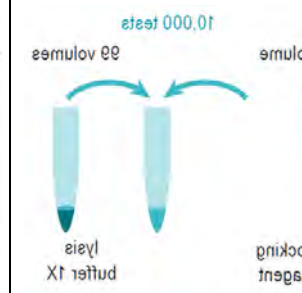
### ► *Supplemented Lysis buffer 4X for two-plate assay manual on suspension cells & one-plate assay manual*

Determine the amount of supplemented lysis buffer needed for the experiment. Each well requires 4 µL of supplemented lysis buffer for one-plate assay manual and 10 µL for two-plate assay manual on suspension cells. Dilute the blocking reagent stock solution 25-fold with lysis buffer 4X. In practice:

TWO-PLATE MANUAL ON SUSPENSION CELL & ONE-PLATE ASSAY MANUAL	
500 TESTS KIT 64ANDRPEG	10,000 TESTS KIT 64ANDRPEH
Preparation of Supplemented Lysis buffer 4X	
 <p>1 volume Blocking reagent</p> <p>24 volumes Lysis buffer 4X</p>	 <p>1 volume Blocking reagent</p> <p>24 volumes Lysis buffer 4X</p>
<p>Dilute the "blocking reagent stock solution" 25-fold with "lysis buffer 4X" e.g take 0.1 ml of "Blocking reagent stock solution" and add it to 2.4 ml of lysis buffer 4X. Mix gently.</p>	

### ► *Supplemented Lysis buffer 1X for two-plate assay manual on adherent cells*

Determine the amount of supplemented lysis buffer needed for the experiment. Each well generally requires 50µL of supplemented lysis buffer 1X. Prepare a lysis buffer solution 1X and then distribute on adherent cells. In practice:



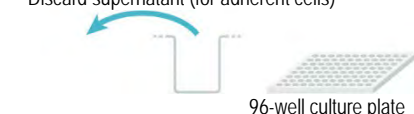
TWO-PLATE ASSAY MANUAL ON ADHERENT CELLS			
500 TESTS KIT 64ANDRPEG & 10,000 TESTS KIT 64ANDRPEH			
Preparation of lysis buffer 1X		Preparation of supplemented Lysis buffer 1X	
 <p>500 tests</p> <p>1 volume</p> <p>3 volumes</p> <p>lysis buffer 4X</p> <p>distilled water</p>	 <p>10,000 tests</p> <p>1 volume</p> <p>3 volumes</p> <p>lysis buffer 4X</p> <p>distilled water</p>	 <p>500 tests</p> <p>1 volume</p> <p>99 volumes</p> <p>blocking reagent</p> <p>lysis buffer 1X</p>	 <p>10,000 tests</p> <p>1 volume</p> <p>99 volumes</p> <p>blocking reagent</p> <p>lysis buffer 1X</p>
<p>Dilute the "lysis buffer 4X" 4-fold with distilled water to prepare lysis buffer 1X. e.g. take 1.25 mL of lysis buffer 4X and add it to 3.75 mL of distilled water. Mix gently.</p>		<p>Dilute the "blocking reagent" 100-fold with "Lysis buffer 1X". e.g. take 0.05 mL of "Blocking reagent stock solution" and add it to 4.95 mL of lysis buffer 1X. Mix gently.</p>	

## TWO PLATE ASSAY MANUAL

## GENERAL LAB WORK PRIOR USING REVVITY KIT: CELL PREPARATION

## FOR ADHERENT CELLS




## FOR SUSPENSION CELLS

1	Plate 50 $\mu\text{L}$ of cells in 96-well tissue-culture treated plate in appropriate growth medium and incubate overnight, at 37°C in CO <sub>2</sub> atmosphere.	Plate 25 $\mu\text{L}$ of cells in 96 half-well plate in your appropriate medium.	
	Cell seeding densities of 100K cells/well are generally sufficient for most cell lines, but the optimization of cell seeding densities is recommended. Depending on the receptor, a starving step with serum-free medium may be essential.		
2	Dispense 50 $\mu\text{L}$ of compound (2X) diluted in cell culture serum-free medium.	Dispense 5 $\mu\text{L}$ of compound (6X), diluted in your appropriate medium.	
	For most compounds, incubation time is 6h to 24h at 37°C. We recommend a time course study to determine the optimal stimulation time.		
3	Carefully remove cell supernatant either by aspirating the supernatant or by flicking the plate.	Do not remove your appropriate medium.	Discard supernatant (for adherent cells) 





## ANDROGEN RECEPTOR DETECTION USING REVVITY KIT

## FOR ADHERENT CELLS

## FOR SUSPENSION CELLS

4	Immediately add 50 $\mu\text{L}$ of lysis buffer (1X) and incubate for at least 30 minutes at room temperature under shaking.	Immediately add 10 $\mu\text{L}$ of supplemented lysis buffer (4X) and incubate for at least 30 minutes at room temperature under shaking.	
	Use the appropriate supplemented lysis buffer and incubate at room temperature under shaking. Lysis incubation time may be optimized. Lysis volume can be decreased down to 25 $\mu\text{L}$ .		
5	After homogenization by pipetting up and down, transfer 16 $\mu\text{L}$ of cell lysate from the 96-well cell-culture plate to a small volume (SV) white detection plate. <i>Depending on the cell lines used, it may be necessary to dilute the cell lysate to ensure samples are within the assay linear range.</i>		
6	Add 4 $\mu\text{L}$ of premixed antibody solutions (vol/vol) prepared in the detection buffer. Cover the plate with a plate sealer. Incubate for 2 hours at room temperature.		
	Set up your reader for Eu <sup>3+</sup> Cryptate and read the fluorescence emission at two different wavelengths (665nm and 620nm) on an HTRF <sup>®</sup> compatible reader.		

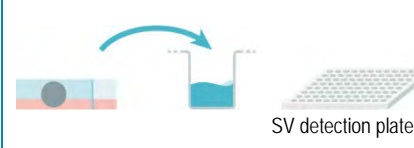
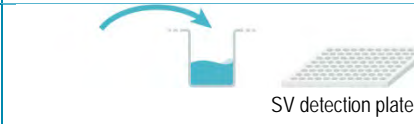
► Standard manual for two-plate assay manual in 20  $\mu\text{L}$  final volume (after lysis step)

	NON TREATED CELL LYSATE	TREATED CELL LYSATE	CONTROL LYSATE	NEGATIVE CONTROL
Step 1 	Dispense 16 $\mu\text{L}$ of non treated cell lysate	Dispense 16 $\mu\text{L}$ of treated cell lysate	Dispense 16 $\mu\text{L}$ of control lysate	Dispense 16 $\mu\text{L}$ of supplemented lysis buffer(1X)
Step 2 	Add 4 $\mu\text{L}$ of premixed Androgen Receptor kit d2 antibody + Androgen Receptor kit Eu cryptate antibody solution to all wells			
Step 3 	Cover the plate with a plate sealer. Incubate for 2h at room temperature.			
Step 4 	Remove the plate sealer and read on an HTRF compatible reader			



The Negative control is used to check the non-specific signal. The ratio between the control lysate signal / non-specific signal should be greater than 2.

## ONE PLATE ASSAY MANUAL







## GENERAL LAB WORK PRIOR USING REVVITY KIT: CELLS PREPARATION

<b>1</b>	Plate <b>8 <math>\mu</math>L of cells</b> in a small volume (SV) white detection plate in your appropriate medium. Cell seeding densities of 30K cells/well are generally sufficient for most cell lines, but optimization of cell seeding densities is recommended. A starving step with serum-free medium can be included.	
<b>2</b>	Dispense <b>4 <math>\mu</math>L of compounds (3X)</b> diluted in your appropriate medium. For most compounds, the incubation time is 4h at 37°C. We recommend a time course study to determine the optimal stimulation time.	

## ANDROGEN RECEPTOR DETECTION USING REVVITY KIT

<b>3</b>	Add <b>4 <math>\mu</math>L of supplemented lysis buffer (4X)</b> . Use the appropriate supplemented lysis buffer and incubate for at least 30 minutes at room temperature under shaking. Lysis incubation time may be optimized.	
<b>4</b>	Add <b>4 <math>\mu</math>L of premixed antibody solutions (vol/vol)</b> prepared in the detection buffer. Cover the plate with a plate sealer. Incubate for <b>2 hours</b> at room temperature. Set up your reader for <b>Eu<sup>3+</sup> Cryptate</b> and read the fluorescence emission at two different wavelengths (665nm and 620nm) on an HTRF <sup>®</sup> compatible reader.	

► Standard manual for one plate assay manual in 20  $\mu$ L final volume

		NON TREATED CELL LYSATE	TREATED CELL LYSATE	NEGATIVE CONTROL	CONTROL LYSATE
<b>GENERAL LAB WORK</b>	<b>Step 1</b> 	Dispense 8 $\mu$ L of cells			
	<b>Step 2</b> 	Add 4 $\mu$ L of your appropriate medium	Add 4 $\mu$ L of compound (3X)	Add 12 $\mu$ L of your appropriate medium	Dispense 16 $\mu$ L of control lysate
<b>TOTAL-ANDROGEN RECEPTOR DETECTION</b>	<b>Step 3</b> 	Add 4 $\mu$ L of lysis buffer (4X) - 30 min/RT.			
	<b>Step 4</b> 	Add 4 $\mu$ L of premixed Androgen Receptor kit d2 antibody + Androgen Receptor kit Eu cryptate antibody solution to all wells			
	<b>Step 5</b> 	Cover the plate with a plate sealer. Incubate for 2h at room temperature.			
	<b>Step 6</b> 	Remove the plate sealer and read on an HTRF compatible reader			

The Negative control is used to check the non-specific signal. The ratio between the control lysate signal / non-specific signal should be greater than 2.

## 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 % 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.revvy.com](http://www.revvy.com)

## RESULTS

These data below should only be considered as an example. Results may vary from one HTRF® compatible reader to another.

The results are drawn by plotting the HTRF® Ratio versus cell lysates from various cell lines in which Androgen Receptor is expressed.

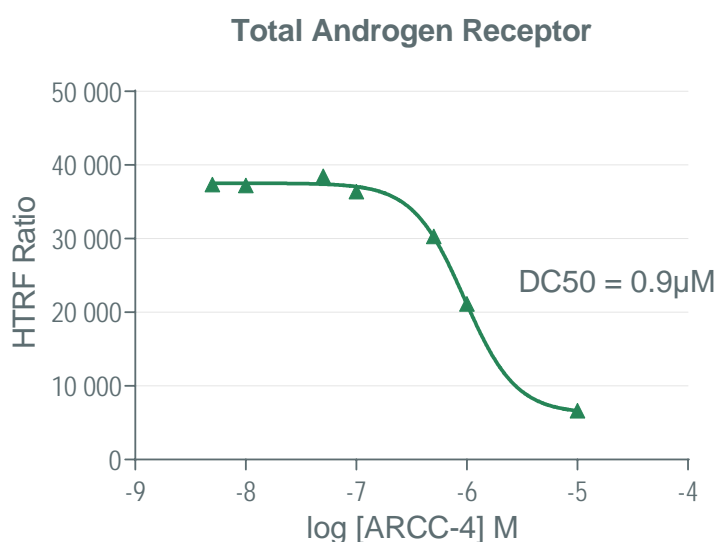
The curves are drawn up by plotting HTRF® Ratio versus the log [compound] concentrations.

The signal linearity is dependent both on the cell line and on the total protein detected. A cell density experiment is highly recommended to ensure working in optimal conditions.

Results on LNCaP cells (75,000 cells/well), using the two-plate assay manual for adherent cells.

Cells were treated for 24h with increasing concentrations of the PROTAC compound ARCC-4. Cells were then lysed with supplemented lysis buffer #3 (1X) for 30 minutes at room temperature. For each condition, a triplet was done by transferring 16µL of each cell lysate into a 384-well plate for the detection.

[ARCC-4] (nM)	Log [ARCC-4] (M)	Mean HTRF Ratio	CV%
5	-8.3	37371	2
10	-8	37226	2
50	-7.3	38478	2
100	-7	36414	1
500	-6.3	30300	3
1 000	-6	21140	1
10 000	-5	6630	2
Negative control		759	10





## GENERAL LAB WORK PRIOR USING REVVITY KIT: CELLS AND LYSATE PREPARATION FREQUENTLY ASKED QUESTIONS / TROUBLESHOOTING PARAMETERS

Using adherent cells, allow time for your cells to recover after plating.	Allow cells to regain full signaling capacity by plating them at least 6 hours before starting the pharmacological treatment.
Depending on the pathway, a serum starving step may be essential to reduce the basal level activity. This step should be optimized case-by-case.	Advice on cell culture conditions prior to using the Revvity kit: <ul style="list-style-type: none"> <li>- <u>For adherent cells</u> Before treating the cells with compounds, remove culture media from the plate and replace it with serum-free media before incubating from 2 hours up to overnight at 37°C.</li> <li>- <u>For suspension cells</u> The starvation step should be carried out in the flask. Harvest cells by centrifugation and re-suspend cells at a suitable cell density in serum-free media. Incubate from 2 hours up to overnight at 37°C.</li> </ul>
Generation of lysates	Ensure that the lysates used for the assay have been generated by using the HTRF <sup>®</sup> lysis buffer provided in the kit. Lysates generated with HTRF <sup>®</sup> buffers can be used in other technologies, like Western Blot. The lysis buffer is effective for creating cell extract under non denaturing conditions from both plated cells and cells pelleted from suspension cultures.
Using the two-plate assay manual, a low signal can often be improved by adjusting lysis volumes.	In most cases, a typical adherent cell line grown in 96-well plates is readily detected in a lysis volume of 50µL. However, the lysis volume can be adjusted from 25 µL to 200 µL.
Using an improper cell density can induce poor sensitivity and low signal.	Check that the cell density is correct. Too high or low cell numbers can affect assay performances.
Parameters such as cell density, stimulation time, and lysis incubation time should be optimized for each cell line used.	The assay can be used for many adherent and non-adherent cell types, including transfected cell lines and primary cells. However, the expression and phosphorylation of the readout of interest can vary from one cell line to another. Depending on the type of treatment, and the temperature, the stimulation time can vary widely. Because of this, we recommend a time course study to determine the optimal compound incubation time. Depending of the nature of your cells, lysis time may vary from 30' to 1h. Because of this, we also recommend determination of the optimal time.
Fluorescence reading	Using an inappropriate setup may seriously impair the results. For information about HTRF <sup>®</sup> compatible readers and for setup recommendations, please visit our website at: <a href="http://www.revvity.com">www.revvity.com</a>
Assaying for multiple targets from a single lysate	The two-plate assay manual indicates the use of 16µL of lysate per well, whereas the 96-well cell culture microplate would generate 50µL (or more) of lysate. Therefore, a typical cell lysate can be assayed for many targets, given that temporal and expression level constraints can vary from one target to another.
Batch production of cell lysates: example of T175 flask	General lab work - prior to using the HTRF Androgen Receptor Revvity kit: <b>Day1:</b> Dispense 10 million cells in T175cm2, add 25 mL of cell culture complete medium, and incubate 2 days at 37°C, with 5% CO <sub>2</sub> or without CO <sub>2</sub> . <b>Day3:</b> Cell stimulation Remove cell culture medium by aspiration, wash once (do not detach the cells), add 5 mL of agonist (1x) diluted in FCS free medium and incubate at 37°C, with 5% CO <sub>2</sub> or without CO <sub>2</sub> , for the optimized time <b>Day3:</b> Cell lysis Remove stimulation medium, wash once (do not detach the cells), add 3 ml of 1X HTRF <sup>®</sup> lysis buffer supplemented with the HTRF <sup>®</sup> blocking reagent for 30 min at room temperature under orbital shaking. Transfer the cell lysate to a 15 mL vial, centrifuge 10 min, 2400 rcf at RT, recover the cell lysate supernatant and store aliquots at -60°C or below. For long term conservation, aliquots should be stored in liquid nitrogen.

### 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.



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](http://www.revvity.com)

revvity

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA  
[www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.