



# ESTRADIOL KITS

## PROTOCOL

### Part # 62ESTPEG & 62ESTPEH

**Test size#:** 500 tests (62ESTPEG) and 10,000 tests (62ESTPEH) - assay volume: 20  $\mu$ L

**Revision:** 05-July 2021

**Store at:** 2-8°C (62ESTPEG); -60°C or below (62ESTPEH)

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

### ASSAY PRINCIPLE

This kit is intended for the simple and rapid quantification of Native Estradiol produced by cells in buffered solution or in cell culture supernatants 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, Estradiol is detected in a competitive assay by using anti Estradiol antibody labeled with Europium cryptate (donor), and Estradiol labeled with XL665 (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 Estradiol present in the sample competes with the binding between the two HTRF detection solutions and thereby prevents FRET from occurring. The specific signal is inversely proportional to the Estradiol concentration.

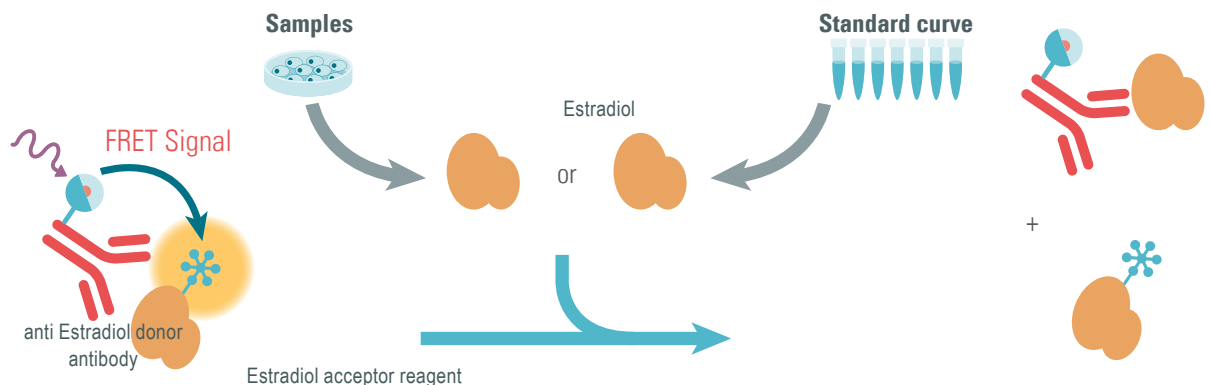
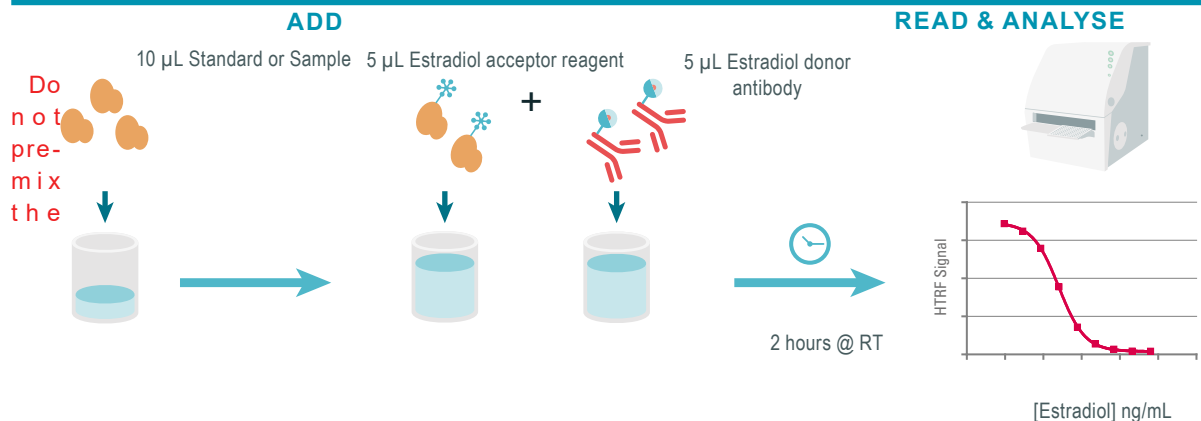


Figure 1: Principle of HTRF Estradiol competitive assay.

### PROTOCOL AT A GLANCE



XL665 and Cryptate solutions prior to dispensing.

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

**MATERIALS PROVIDED:**

| KIT COMPONENTS                               | 500 TESTS *<br>CAT # 62ESTPEG    | 10,000 TESTS *<br>CAT # 62ESTPEH |
|--|----------------------------------|----------------------------------|
| Estradiol Standard<br>Lyophilized            | 1 vial<br>Concentrated Estradiol | 1 vial<br>Concentrated Estradiol |
| anti Estradiol antibody Eu Cryptate antibody | 1 vial<br>Lyophilized            | 1 vial - 1 mL<br>Frozen - 50X    |
| Estradiol XL665 reagent                      | 1 vial<br>Lyophilized            | 1 vial - 1 mL<br>Frozen - 50X    |
| Diluent **<br>ready-to-use                   | 1 vial<br>20 mL                  | 1 vial<br>20 mL                  |
| Detection buffer ***<br>ready to use         | 1 vial<br>1 mL                   | 1 vial<br>5 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.

\*\* 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 [cisbio.com/compatible-readers](http://cisbio.com/compatible-readers)

- Small volume (SV) detection microplates - Use white plate only..

For more information about microplate recommendations, please visit our website at: [cisbio.com/microplates-recommendations](http://cisbio.com/microplates-recommendations)

**STORAGE AND STABILITY**

Store the kit 62ESTPEG at 2-8°C and the kit 62ESTPEH at -60°C or below.

Under proper storage conditions, reagents are stable until the expiry date indicated on the label.



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 (Can be stored 7 days at 4°C) .

**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.
- 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.
- Estradiol 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 - 62ESTPEG  |   |   | 10,000 TESTS KIT - 62ESTPEH   |
|---|---|---|---|
| anti Estradiol antibody Eu Cryptate antibody  |   |   |   |
| Reconstitute the anti Estradiol antibody Eu Cryptate antibody with 2.5 mL detection buffer. Mix gently. This ready to use 1X stock solution can be frozen and stored at -60°C or below. It can be stored unfrozen 7 days at 4°C.  |  |  | Thaw the anti Estradiol antibody Eu Cryptate antibody . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below. It can be stored unfrozen 7 days at 4°C.  |
| Estradiol XL665 reagent   |   |   |   |
| Reconstitute the Estradiol XL665 reagent with 2.5 mL detection buffer. Mix gently. This ready to use 1X stock solution can be frozen and stored at -60°C or below. It can be stored unfrozen 7 days at 4°C.   |  |  | Thaw the Estradiol XL665 reagent . Mix gently. This 50X stock solution can be frozen and stored at -60°C or below. It can be stored unfrozen 7 days at 4°C.   |
| Estradiol Standard  |   |   |   |
| Reconstitute the Estradiol Standard with distilled water in order to obtain a 100 ng/mL stock solution. See instructions on vial label for reconstitution volume. Mix gently after reconstitution. This stock solution is stable 2 days at 4°C. It can be frozen and stored at -60°C or below and thawed once only. |  |   | Reconstitute the Estradiol Standard with distilled water in order to obtain a 100 ng/mL stock solution. See instructions on vial label for reconstitution volume. Mix gently after reconstitution. This stock solution is stable 2 days at 4°C. It can be frozen and stored at -60°C or below and thawed once only. |
| Diluent   |   |   |   |
| The diluent is ready-to-use   |   |   | The diluent is ready-to-use   |
| 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 anti Estradiol antibody Eu Cryptate antibody and 5 µL Estradiol XL665 reagent.

Prepare the two solutions in separate vials.

| 500 TESTS KIT - 62ESTPEG   |   |   | 10,000 TESTS KIT - 62ESTPEH  |
|--|---|---|--|
| anti Estradiol antibody Eu Cryptate antibody                                       |   |   |  |
| After reconstitution, the Estradiol Eu Cryptate antibody is ready to use.          |  |  | Dilute 50-fold the stock solution of Estradiol Eu Cryptate antibody with detection buffer e.g. take 1 mL of Eu Cryptate antibody stock solution and add it to 49 mL of detection buffer. |
| Estradiol XL665 reagent  |   |   |  |
| After reconstitution, the Estradiol XL665 reagent is ready to use.                 |  |  | Dilute 50-fold the stock solution of Estradiol XL665 reagent with detection buffer e.g. take 1 mL of Eu Cryptate antibody stock solution and add it to 49 mL of detection buffer.        |
| Antibody mix   |   |   |  |
| <b>Do not pre-mix the XL665 and the Eu Cryptate solutions prior to dispensing.</b> |   |   |  |

## TO PREPARE STANDARD WORKING SOLUTIONS:

- Each well requires 10  $\mu\text{L}$  of standard.
- Dilute the standard stock solution serially with diluent
- 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 .
- In order to counteract any standard sticking, we recommend changing tips between each dilution.

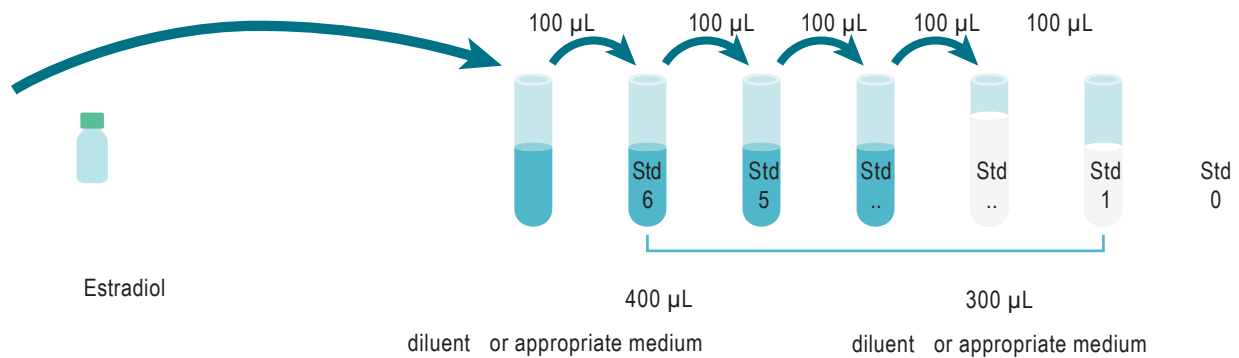
A recommended standard dilution procedure is listed and illustrated below:

Dilute the standard stock solution 5-fold with diluent to prepare high standard (Std 6): e.g. take 100  $\mu\text{L}$  of standard stock solution and add it to 400  $\mu\text{L}$  of diluent . Mix gently.

Use the high standard (Std 6) to prepare the standard curve using 1/4 serial dilutions as follows:

- Dispense 300  $\mu\text{L}$  of diluent in each vial from Std 5 to Std 0.
- Add 100  $\mu\text{L}$  of standard to 300  $\mu\text{L}$  of diluent , mix gently and repeat the 1/4 serial dilution to make standard solutions: std6, std5, std4, std3, std2, std1.

This will create 6 standards for the analyte. Std 0 (Positive control) is diluent or appropriate culture medium alone.








| STANDARD                | SERIAL DILUTIONS   | ESTRADIOL WORKING SOLUTION (ng/mL) |
|-------------------------|--|------------------------------------|
| Standard Stock solution | Reconstituted lyophilisate   | 100                                |
| Standard 6              | 100 $\mu\text{l}$ reconstituted standard + 400 $\mu\text{L}$ Diluent | 20                                 |
| Standard 5              | 100 $\mu\text{L}$ standard 6 + 300 $\mu\text{L}$ Diluent             | 5                                  |
| Standard 4              | 100 $\mu\text{L}$ standard 5 + 300 $\mu\text{L}$ Diluent             | 1.25                               |
| Standard 3              | 100 $\mu\text{L}$ standard 4 + 300 $\mu\text{L}$ Diluent             | 0.312                              |
| Standard 2              | 100 $\mu\text{L}$ standard 3 + 300 $\mu\text{L}$ Diluent             | 0.078                              |
| Standard 1              | 100 $\mu\text{L}$ standard 2 + 300 $\mu\text{L}$ Diluent             | 0.019                              |
| Standard 0              | 300 $\mu\text{L}$ Diluent  | 0                                  |

## TO PREPARE SAMPLES:

- Each well requires 10  $\mu$ 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 6) must be diluted diluent or in your appropriate sample medium.

## ASSAY PROTOCOL

|  |  | Negative control<br>or Cryptate control                              | Standard (Std 0 - Std 6)   | Samples  |
|--|--|--|--|--|
| <b>Step 1</b><br>   |  | Dispense 10 $\mu$ L of diluent into each negative control well       | Dispense 10 $\mu$ L of each Estradiol standard (Std 0 - Std 6) into each standard well | Dispense 10 $\mu$ L of each sample into each sample well |
| <b>Step 2</b><br>   |  | Add 5 $\mu$ L of Detection buffer to all negative control wells      | Add 5 $\mu$ L Estradiol acceptor reagent working solution to all wells                 |  |
| <b>Step 3</b><br>   |  | Add 5 $\mu$ L Estradiol donor antibody working solution to all wells |  |  |
| <b>Step 4</b><br>  |  | Seal the plate and incubate 2 hours @ RT                             |  |  |
| <b>Step 5</b><br> |  | 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 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>

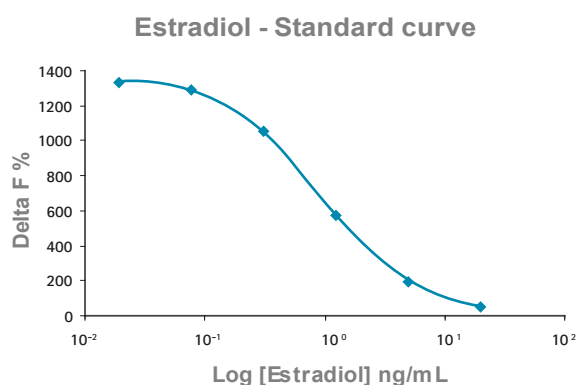
## RESULTS

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.

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

|                          | Ratio <sup>(1)</sup> | CV <sup>(2)</sup> | Delta F% <sup>(3)</sup> |
|--------------------------|----------------------|-------------------|-------------------------|
| Negative control         | 366                  | 1.5%              |                         |
| Std 0 – Positive control | 5,471                | 0.4%              | 1,395%                  |
| Std 1 - 0.019 ng/mL      | 5,250                | 0.8%              | 1,334%                  |
| Std 2 - 0.078 ng/mL      | 5,103                | 1.6%              | 1,294%                  |
| Std 3 - 0.312 ng/mL      | 4,226                | 1.2%              | 1,055%                  |
| Std 4 - 1.25 ng/mL       | 2,479                | 3%                | 577%                    |
| Std 5 - 5 ng/mL          | 1,073                | 0.9%              | 193%                    |
| Std 6 - 20 ng/mL         | 560                  | 4%                | 53%                     |



## ANALYTICAL CHARACTERISTICS

### DETECTION LIMIT AND CONVERSION TO NMOL/L

The minimum detectable dose of estradiol is 0.019 ng/ml (dose of mean zero - 2SD). Estradiol concentrations in ng/mL can easily be converted to nmol/l using the following formula : 1 ng/mL = 3.67 nmol/L.

### CROSS-REACTIVITY

|                         | Cross-reactivity (%) |                   | Cross-reactivity (%) |
|-------------------------|----------------------|-------------------|----------------------|
| Estradiol               | 100                  | Tamoxifen         | 0.007                |
| Estradiol 3 glucuronide | 58.1                 | 17 a Estradiol    | 0.006                |
| Estradiol 3 sulfate     | 12.3                 | Progesterone      | 0.002                |
| 17 a Ethynylestradiol   | 0.65                 | Estrone 3 sulfate | 0.002                |
| Estrone                 | 0.4                  | Corticosterone    | <0.001               |
| Estriol                 | 0.3                  | Cortisone         | <0.001               |
| estosterone             | 0.2                  | Cortisol          | <0.001               |
| Danazol                 | 0.02                 |                   |                      |

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](http://www.cisbio.com) to find a list of our regional distributors



cisbio