



## 1.0 INTRODUCTION

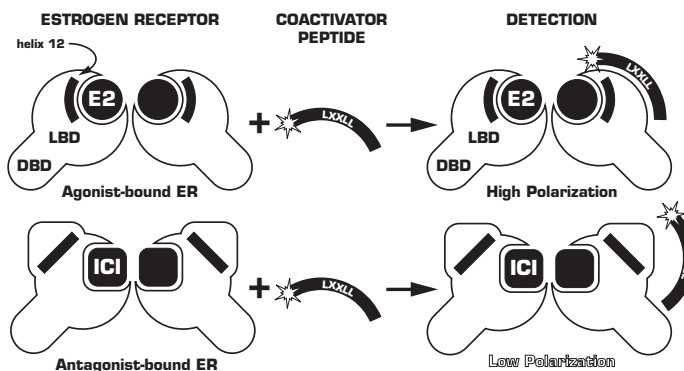
The Estrogen Receptor- $\alpha$  (ER $\alpha$ ) Coactivator Assay is designed as a screening assay for profiling known ER $\alpha$  ligands or test compounds. This kit contains the necessary reagents to perform assays:

1. To assess the ligand *dose dependency* for the recruitment of coactivator peptides to human ER $\alpha$  or
2. Alternatively, to perform *equilibrium binding* to measure the affinity of ER $\alpha$  for the coactivator peptide in the presence of saturating amounts of ligand.

The assays use a fluorescent coactivator peptide (D22), which is a rhodamine-labeled coactivator-like peptide containing an LXXLL motif and flanking sequences that resemble known coactivators (1).

When used to assess ligand *dose dependency*, ER $\alpha$  and D22 are added to increasing concentrations of ligand (or test compound), which results in either the formation or disruption of the ER $\alpha$ /D22 complex. Agonist-bound ER $\alpha$  can recruit D22, resulting in a larger fraction of bound D22 and therefore a higher polarization value (expressed as millipolarization units, mP) than the no-ligand control. In contrast, antagonist-bound ER $\alpha$  has a lower affinity for D22, yielding a smaller fraction of bound D22 and therefore a lower polarization value than the no-ligand control. The EC<sub>50</sub> value of the ligands, either to promote or to disrupt the ER $\alpha$ /D22 interaction, provides a means to classify test compounds as antagonists, agonists, or selective modulators.

An alternative protocol (for *equilibrium binding*) is provided to measure the affinity of the ER $\alpha$ /D22 interaction in the presence of saturating amounts of ligand (or test compound) with increasing concentrations of ER $\alpha$ . This equilibrium binding format offers another method to classify test compounds as antagonists, agonists, or selective modulators. This method provides a larger dynamic range for antagonist ligands.



**Figure 1.** Principle of the ER $\alpha$  Coactivator Assay. E2 = agonist; ICI = antagonist.

Please see PanVera's website for more information about:

- ER $\alpha$  Coactivator Assay Kit, Red ([www.panvera.com/catalog/P3071.html](http://www.panvera.com/catalog/P3071.html)).
- ER $\beta$  Coactivator Assay Kit, Red ([www.panvera.com/catalog/P2994.html](http://www.panvera.com/catalog/P2994.html)).
- Fluorescence polarization theory and techniques in PanVera's on-line Fluorescence Polarization Applications Guide (<http://www.panvera.com/tech/appguide/index.html>).

## 2.0 SAFETY PRECAUTIONS

Exercise normal precautions, such as using gloves, lab coats, eye protection and a fume hood, when working with any chemical reagents. Although the reagents in this kit are considered non-hazardous according to 29 CFR 1910.1200, the chemical, physical and toxicological properties of these products may not, as yet, have been thoroughly investigated.



### 3.0 DESCRIPTION

#### 3.1 Materials Supplied

Description	Composition	Amount	Part No.
10X D22 Coactivator peptide, Red (LPYEGSLLKLLRAPVEEV) (1)	10X in buffer (pH 7.5) containing peptide stabilizing agents	2 $\times$ 1 mL	P2993
Estrogen Receptor-alpha (ER $\alpha$ ), Human Recombinant	50 mM Tris•HCl (pH 8), 500 mM KCl, 2 mM DTT, 1 mM EDTA, 1 mM sodium orthovanadate and 10% glycerol	2 $\times$ 750 pmol	P2187
ER $\alpha$ Coactivator Assay Buffer	Buffer (pH 7.5) containing protein stabilizing agents	20 mL	P3070
1 M DTT	Aqueous	1 mL	P2325

#### 3.2 Materials Required But Not Supplied

- Fluorescence polarization (FP) instrument with suitable 535 nm excitation and 590 nm emission interference filters
- Black, round-bottom 384-well plates
- Red (FP) Standardization Kit (PanVera® Part No. P2888), recommended for verifying FP instrument performance. The Red Polarization Standard (RPS) may be used to determine whether the instrument is measuring polarization values accurately.
- Control Agonist: 17 $\beta$ -estradiol is recommended; 250  $\mu$ M stock in DMSO, ethanol (EtOH) or methanol (MeOH) suggested
- Control Antagonists: 4-OH-tamoxifen or ICI 182,780 is recommended; 250  $\mu$ M stock in DMSO, EtOH or MeOH suggested

### 4.0 STORAGE AND STABILITY

Description	Storage Temperature	Notes	Part No.
10X D22 Coactivator peptide, Red (1)	-20°C		P2993
Estrogen Receptor-alpha (ER $\alpha$ ), Human Recombinant	-80°C	Do not expose to more than four freeze/thaw cycles.	P2187
ER $\alpha$ Coactivator Assay Buffer	4°C	Store at 4°C after first use.	P3070
1 M DTT	-20°C		P2325

### 5.0 GENERAL CONSIDERATIONS FOR THE ASSAYS

- **Controls:** Use a control agonist, such as 17 $\beta$ -estradiol, and a control antagonist, such as 4-OH-tamoxifen or ICI 182,780, in each experiment for comparison to test compounds. In addition, include the controls specified under the specific assay experimental procedure for comparison to test compounds.
- **Solvents:** We recommend using a minimal amount of solvent in the assay. However, both assay formats will tolerate a final concentration of up to 5% DMSO, 4% EtOH or 4% MeOH.
- **Handle ER $\alpha$  gently:** For best results, thaw ER $\alpha$  on ice before use. **Never vortex ER $\alpha$ .**
- **Verify Instrument Calibration:** Verify that the instrument is calibrated properly and has suitable 535 nm excitation and 590 nm emission interference filters.

**6.0 DOSE DEPENDENCY ASSAY PROCEDURE****6.1 Introduction**

This kit provides sufficient reagents to perform 400 (40- $\mu$ L) reactions (dose dependency format) in a 384-well plate. Each well of a set contains a constant concentration of D22 (1X) and ER $\alpha$  (75 nM) with serially diluted ligand. Each set will assay a different ligand and should contain at least 16 wells (ligand dilutions). Alternatively, the assay may be performed in a 96-well plate (100- $\mu$ L reactions) by proportionally scaling up the volume of reagents used per well.

**6.2 Dose Dependency Protocol (384-well format)**

1. Place ER $\alpha$ , 10X D22, DTT and ER $\alpha$  Coactivator Assay Buffer on ice until ready to use.
2. Prepare the Complete ER $\alpha$  Coactivator Assay Buffer by adding sufficient 1 M DTT to the ER $\alpha$  Coactivator Assay Buffer to yield a final DTT concentration of 10 mM.
3. Dispense 20  $\mu$ L of Complete ER $\alpha$  Coactivator Assay Buffer into the wells of each ligand set in a 384-well plate, except for the first well. This first well will contain the highest ligand concentration. Use at least 16 wells (ligand dilutions) per set.
4. Prepare a 2X concentration of ligand in Complete ER $\alpha$  Coactivator Assay Buffer. Add 40  $\mu$ L of 2X ligand to the first well of the set.

**NOTE:** Although a recommended starting 2X concentration of 10  $\mu$ M is suggested for known ER $\alpha$  ligands, using higher concentrations of a ligand may be necessary if its affinity for ER $\alpha$  is low.

5. Perform two-fold serial dilutions of 20  $\mu$ L of ligand in Complete ER $\alpha$  Coactivator Assay Buffer from Well #1 through Well #16.
6. Prepare a 2X Master Mix containing 2X D22 and 150 nM ER $\alpha$ , diluted in Complete ER $\alpha$  Coactivator Assay Buffer, as shown below. The empty boxes underneath each example equation are for your calculations. The volumes given below are appropriate for 16 (40- $\mu$ L) reactions.

**First calculate the total volume of 2X Master Mix needed:**

$$\begin{array}{l} 20 \mu\text{L/well} \times 16 \text{ wells} \times 1.1 \text{ (for pipetting error)} = 352 \mu\text{L (total volume)} \\ \boxed{\phantom{000}} \mu\text{L/well} \times \boxed{\phantom{000}} \text{ wells} \times 1.1 \text{ (for pipetting error)} = \boxed{\phantom{000}} \mu\text{L (total volume)} \end{array}$$

**Calculate the volume of ER $\alpha$  stock needed:**

$$\begin{array}{l} [ 352 \mu\text{L} \times 150 \frac{\text{nM}}{1} \div 2000 \text{ nM ER}\alpha \text{ stock}^* = 26.4 \mu\text{L ER}\alpha \\ [ \boxed{\phantom{000}} \mu\text{L} \times 150 \frac{\text{nM}}{1} \div \boxed{\phantom{000}} \text{ nM ER}\alpha \text{ stock}^* = \boxed{\phantom{000}} \mu\text{L ER}\alpha \end{array}$$

\*The ER $\alpha$  concentration is lot specific. Use the concentration provided on the ER $\alpha$  tube or Certificate of Analysis to calculate the appropriate volume of concentrated ER $\alpha$ .

**Calculate the volume of D22 needed:**

$$\begin{array}{l} [ 352 \mu\text{L} \times 2\text{X Stock} ] \div 10\text{X stock} = 70.4 \mu\text{L D22} \\ [ \boxed{\phantom{000}} \mu\text{L} \times 2\text{X Stock} ] \div 10\text{X stock} = \boxed{\phantom{000}} \mu\text{L D22} \end{array}$$

**Calculate volume of Complete ER $\alpha$  Coactivator Assay Buffer needed:**

$$\begin{array}{l} [ 352 \mu\text{L} - 26.4 \mu\text{L ER}\alpha^* - 70.4 \mu\text{L D22} = 255.2 \mu\text{L Complete ER}\alpha \text{ Coactivator Assay Buffer} \\ [ \boxed{\phantom{000}} \mu\text{L} - \boxed{\phantom{000}} \mu\text{L ER}\alpha^* - \boxed{\phantom{000}} \mu\text{L D22} = \boxed{\phantom{000}} \mu\text{L Complete ER}\alpha \text{ Coactivator Assay Buffer} \end{array}$$

\*The ER $\alpha$  concentration is lot specific. Use the concentration provided on the ER $\alpha$  tube or Certificate of Analysis to calculate the appropriate volume of concentrated ER $\alpha$ .

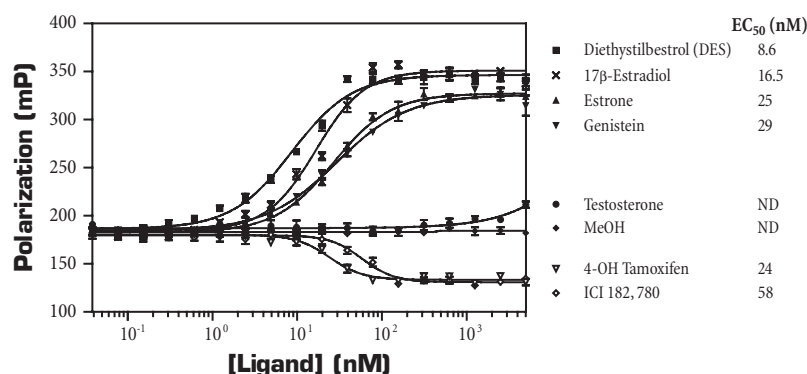
7. Gently mix the appropriate volumes of 10X D22, ER $\alpha$  and Complete ER $\alpha$  Coactivator Assay Buffer in a tube.
8. Dispense 20  $\mu$ L of 2X Master Mix into each well.



9. Prepare the following set of controls and standards (3–5 replicates per control or standard, 40  $\mu$ L per well), as follows:
  - 1X D22 in Complete ER $\alpha$  Coactivator Assay Buffer
  - 1X ER $\alpha$ /D22 Coactivator peptide (20  $\mu$ L of 2X Master Mix and 20  $\mu$ L of Complete ER $\alpha$  Coactivator Assay Buffer)
  - Complete ER $\alpha$  Coactivator Assay Buffer as a blank
  - 1:10 dilution of RPS in RPS buffer
  - RPS buffer as a blank for RPS
10. Mix components in 384-well plate by gentle agitation. Protect the plate from light and evaporation by covering with foil and then incubate for 1 to 5 hours at room temperature (20–25°C).
11. Measure polarization values.

## 7.0 RESULTS AND DISCUSSION – DOSE DEPENDENCY FORMAT

Below is an example of the dose dependency data generated using the ER $\alpha$  Coactivator Assay. In these experiments, 10  $\mu$ M (2X) of ligand was used as the starting concentration. The concentration of the ligand that results in a half-maximal increase or decrease in polarization equals the ligand EC<sub>50</sub> for the ER $\alpha$ /D22 interaction. The EC<sub>50</sub> values obtained in this example are listed after each ligand. The curve was fit using the sigmoidal dose response curve (variable slope) with Prism® software (GraphPad Software, Inc., San Diego, CA).



**Figure 2. Dose dependency of the ER $\alpha$ /D22 interaction.** Ligand was serially diluted in Complete ER $\alpha$  Coactivator Assay Buffer in a 384-well plate with final reaction conditions containing 75 nM ER $\alpha$  and 1X D22. The samples were incubated for 1 hour at room temperature. Values reported are ligand EC<sub>50</sub> values for the ER $\alpha$ /D22 interaction. ND = not determined; MeOH = unliganded ER $\alpha$  in solvent.

## 8.0 EQUILIBRIUM BINDING ASSAY PROCEDURE

### 8.1 Introduction

In this alternative protocol, an equilibrium binding curve can be generated by adding a constant concentration of D22 (1X) to increasing concentrations of ER $\alpha$  in the presence of a saturating concentration (typically 5  $\mu$ M) of ligand. Polarization values will be plotted against the concentration of ER $\alpha$ . The concentration of ER $\alpha$  that results in a half-maximal shift in polarization equals the EC<sub>50</sub> of the ER $\alpha$ /D22 interaction. The assay may be performed in a 96-well plate (100- $\mu$ L reactions) by proportionally scaling up the volume of reagents used per well.

### 8.2 Equilibrium Binding Protocol (384-well format)

1. Place ER $\alpha$ , 10X D22, DTT and ER $\alpha$  Coactivator Assay Buffer on ice until ready to use.
2. Prepare the Complete ER $\alpha$  Coactivator Assay Buffer by adding sufficient 1 M DTT to the ER $\alpha$  Coactivator Assay Buffer to yield a final DTT concentration of 10 mM.
3. Dispense 20  $\mu$ L of Complete ER $\alpha$  Coactivator Assay Buffer into each well set in a 384-well plate except for the first well of each ligand. This first well will contain the highest concentration of ER $\alpha$ . Use at least 16 wells (ER $\alpha$  dilutions) per set.
4. Prepare a 2X concentration of ER $\alpha$  (typically 1–4  $\mu$ M) in Complete ER $\alpha$  Coactivator Assay Buffer. Add 40  $\mu$ L of 2X ER $\alpha$  to the first well of the set.
5. Perform two-fold serial dilutions of 20  $\mu$ L of ER $\alpha$  from Well #1 through Well #16.



6. Prepare a 2X Master Mix containing a saturating concentration of ligand (2X, typically 10  $\mu$ M) and 2X D22, diluted in Complete ER $\alpha$  Coactivator Assay Buffer, as shown below. The empty boxes underneath each example equation are for your calculations. The volumes given in the table below are appropriate for 16 (40- $\mu$ L) reactions

**NOTE:** It may be necessary to use higher concentrations of a test compound if its affinity for ER $\alpha$  is low.

**First calculate the total volume of Master Mix needed:**

$$\begin{array}{l} 20 \text{ } \mu\text{L/well} \times 16 \text{ wells} \times 1.1 \text{ (for pipetting error)} = 352 \text{ } \mu\text{L (total volume)} \\ \boxed{\phantom{000}} \text{ } \mu\text{L/well} \times \boxed{\phantom{000}} \text{ wells} \times 1.1 \text{ (for pipetting error)} = \boxed{\phantom{000}} \text{ } \mu\text{L (total volume)} \end{array}$$

**Calculate the volume of D22 needed:**

$$\begin{array}{l} [ 352 \text{ } \mu\text{L} \times 2\text{X Stock} ] \div 10\text{X stock} = 70.4 \text{ } \mu\text{L D22} \\ [ \boxed{\phantom{000}} \text{ } \mu\text{L} \times 2\text{X Stock} ] \div 10\text{X stock} = \boxed{\phantom{000}} \text{ } \mu\text{L D22} \end{array}$$

**Calculate the volume of ligand needed:**

$$\begin{array}{l} [ 352 \text{ } \mu\text{L} \times 10 \text{ } \mu\text{M ligand*} ] \div 250 \text{ } \mu\text{M ligand*} = 14 \text{ } \mu\text{L ligand} \\ [ \boxed{\phantom{000}} \text{ } \mu\text{L} \times \boxed{\phantom{000}} \text{ } \mu\text{M ligand*} ] \div \boxed{\phantom{000}} \text{ } \mu\text{M ligand*} = \boxed{\phantom{000}} \text{ } \mu\text{L ligand} \end{array}$$

**NOTE:** Although a recommended starting 2X concentration of 10  $\mu$ M is suggested for known ER $\alpha$  ligands, using higher ligand concentrations may be necessary if the ligand has low affinity for ER $\alpha$ .

**Calculate volume of Complete ER $\alpha$  Coactivator Assay Buffer needed:**

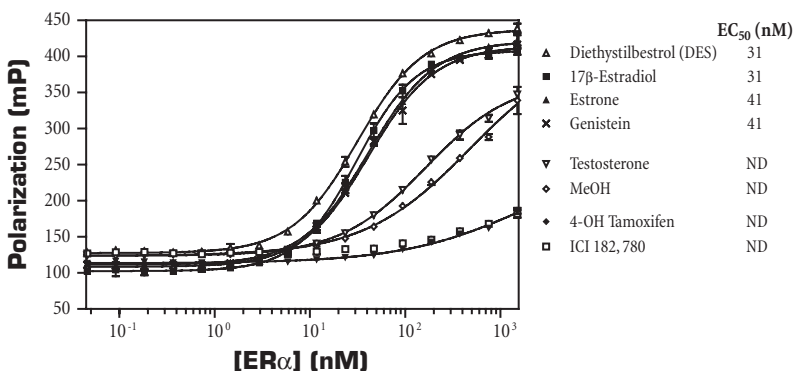
$$\begin{array}{l} 352 \text{ } \mu\text{L} - 70.4 \text{ } \mu\text{L D22} - 14 \text{ } \mu\text{L ligand} = 267.6 \text{ } \mu\text{L ER}\alpha \text{ Complete Coactivator Assay Buffer} \\ \boxed{\phantom{000}} \text{ } \mu\text{L} - \boxed{\phantom{000}} \text{ } \mu\text{L D22} - \boxed{\phantom{000}} \text{ } \mu\text{L ligand} = \boxed{\phantom{000}} \text{ } \mu\text{L ER}\alpha \text{ Complete Coactivator Assay Buffer} \end{array}$$

7. Gently mix the appropriate volumes of 10X D22, ligand and Complete ER $\alpha$  Coactivator Assay Buffer in a tube.
8. Dispense 20  $\mu$ L of 2X Master Mix into each well.
9. Prepare the following set of controls and standards (3–5 replicates per control or standard, 40  $\mu$ L per well), as follows:
- 1X D22 in Complete ER $\alpha$  Coactivator Assay Buffer
  - 1X ligand/D22 in Complete ER $\alpha$  Coactivator Assay Buffer (20  $\mu$ L of 2X Master Mix and 20  $\mu$ L of Complete ER $\alpha$  Coactivator Assay Buffer)
  - Complete ER $\alpha$  Coactivator Assay Buffer as a blank
  - 1:10 dilution of RPS in RPS buffer
  - RPS buffer as a blank for RPS
10. Mix components in 384-well plate by gentle agitation. Protect plate from light and evaporation by covering with foil and then incubate for 1 to 5 hours at room temperature (20–25°C).
11. Measure polarization values.



## 9.0 RESULTS AND DISCUSSION – EQUILIBRIUM BINDING FORMAT

Below is an example of the equilibrium binding data generated using the ER $\alpha$  Coactivator Assay with D22. In this experiment, the final concentration of each ligand was 5  $\mu$ M. The concentration of the ER $\alpha$  that results in a half-maximal shift in polarization equals the EC<sub>50</sub> of the ER $\alpha$ /D22 interaction. The EC<sub>50</sub> values for the ER $\alpha$ /D22 interaction are listed after each ligand. The EC<sub>50</sub> values can be converted mathematically to equilibrium binding constants ( $K_d$ ) by plotting bound ER $\alpha$  vs. log(free ER $\alpha$ ), also known as a Klotz plot. Agonists promote interaction of ER $\alpha$  with D22, but no interaction is detected in the presence of antagonists. The curve was fit using the sigmoidal dose response curve (variable slope) with Prism® software.



**Figure 3. Affinity of ER $\alpha$  for D22 in the presence of various ligands.** Purified ER $\alpha$  was serially diluted in Complete ER $\alpha$  Coactivator Assay Buffer in a 384-well plate with final reaction conditions containing 5  $\mu$ M ligand, 1X D22 and 0.5% MeOH. The samples were incubated for 1 hour at room temperature. Values reported are the EC<sub>50</sub> values for the interaction of ligand-occupied ER $\alpha$  and D22. ND = not determined; MeOH = unliganded ER $\alpha$  in solvent.

## 10.0 REFERENCES

1. Chang, C. *et al.* (1999) *Mol. Cell. Biol.* **19**:8226-39.

The performance of this product is guaranteed for six months from the date of purchase if stored and handled properly.

PanVera® is a registered trademark of PanVera® LLC.

Prism® is a registered trademark of GraphPad™ Software, Inc., San Diego, CA.

© 2002, PanVera® LLC. All rights reserved.

PURCHASER SHALL USE THE PRODUCTS PURCHASED HEREUNDER (the "Products") SOLELY FOR THE PURPOSE OF CONDUCTING INTERNAL RESEARCH AT ITS ORGANIZATION ("RESEARCH"). Purchaser will not sell, transfer, disclose or otherwise provide access to the Products to any third party. Purchaser agrees that Purchaser shall not have the right to authorize any third party to use or sell any Products or derivatives thereof. Unless a license has been executed between Purchaser and PanVera® LLC explicitly providing otherwise, Purchaser will not: (i) reformulate or create derivatives of the Products; or (ii) use the Products for providing services to a third party (e.g. screening or profiling); (iii) use the Products in a diagnostic process; (iv) use the Products in a quality control or quality assurance process for the manufacture of a product for sale; or (v) use the Products as a component of a kit.