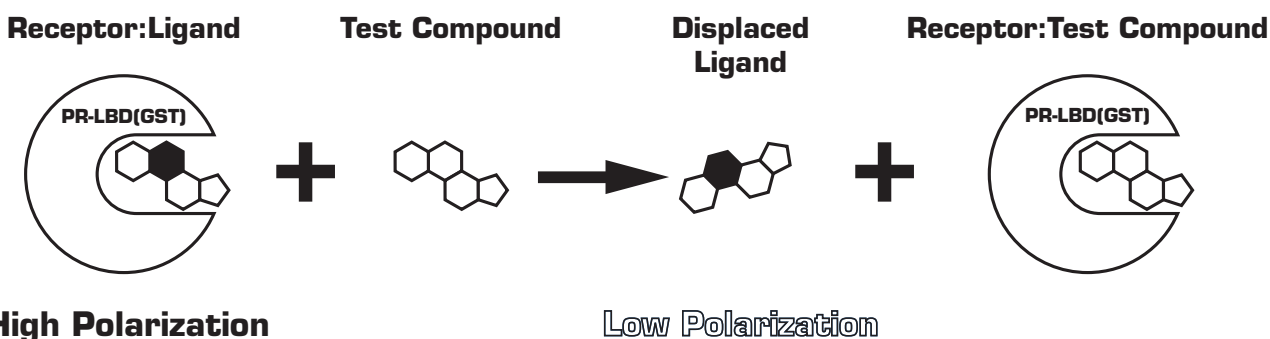


1.0 INTRODUCTION

This kit contains the necessary reagents to perform a competition assay for identification of progesterone receptor binding compounds. An N-terminal fusion of glutathione transferase to the ligand-binding domain of the human progesterone receptor (PR-LBD(GST)) and a proprietary, fluorescently-tagged progesterone ligand (Fluormone™ PL Green) are utilized. The presence of effective competitors prevents the formation of Fluormone™ PL Green/PR-LBD(GST) complex, which has high fluorescence polarization value. The relative binding affinities of test compounds can be determined by measuring the polarization value.



This kit contains enough reagents to perform 100 assays in a total volume of 100 µL in a 96-well plate.

If you would like more information, please see our on-line Fluorescence Polarization Applications Guide at:

<http://www.panvera.com/tech/fpguide/index.html>

If you would like more information about this specific product, please see our website at:

<http://www.panvera.com/catalog/P2895.html>

2.0 SAFETY PRECAUTIONS

Normal precautions exercised in handling laboratory reagents should be followed. All reagents in this kit are considered non-hazardous according to 29 CFR 1910.1200. The Fluormone™ PL Green may have steroid activity *in vivo* and therefore should be treated with caution. Methanol is toxic and may be absorbed through the skin. The chemical, physical, and toxicological properties of these products may not, as yet, have been thoroughly investigated. We recommend using gloves, lab coats, eye protection, and a fume hood when working with any chemical reagents.

3.0 DESCRIPTION

3.1 Materials Supplied

Description	Composition	Amount	Part No.
Fluormone™ PL Green	400 nM in 90% methanol	50 µL	P2897
PR-LBD(GST)	50 mM Tris (pH 8.0), 1M Urea, 500 mM KCl, 1 mM EDTA, 5 mM DTT and 50% glycerol	400 pmol	P2899
PR Screening Buffer, Green Assay	Buffer (pH 7.4) containing protein stabilizing agents and glycerol	20 mL	P2901
1 M DTT	in water	1 mL	P2325

3.2 Materials Required but Not Supplied

Multiwell fluorescence polarization instrument with suitable 485 nm excitation and 530 nm emission interference filters
Pipetting devices P20, P200, and P1000, suitable repeater pipettors, or multi-channel pipettors
Black multiwell plates for use in the fluorescence polarization instrument
Beacon® One-Step FP Standardization Kit (PanVera Part No. P2581): recommended for calibrating instrument

3.3 Storage and Stability

Description	Storage Temperature	Notes	Part No.
Fluormone™ PL Green	-20° C	PR-LBD(GST)/Fluormone™ PL Green complexes should be prepared in glass and stored on ice until dispensed.	P2897
PR-LBD(GST)	-80° C	Avoid repeated freeze-thaw cycles of PR-LBD(GST). PR-LBD(GST) must remain on ice once thawed.	P2899
PR Screening Buffer, Green Assay	Room Temperature	--	P2901
1 M DTT	-20° C	--	P2325

4.0 ASSAY CONSIDERATIONS

This kit contains enough reagents to perform 100, 100 µL assays in a 96-well plate. A competition curve is generated by adding Fluormone™ PL Green/PR-LBD(GST) complexes to a dilution series of the test compound and then plotting the polarization against the concentration of test compound. The test compound concentration resulting in a half-maximum polarization value shift equals its IC₅₀.

4.1 General Considerations

Controls: Wells that contain each of the following reagents should be included: 1X PR-LBD(GST), 1X Fluormone™ PL Green/PR-LBD(GST) complex, and PR Screening Buffer, Green Assay. Also, a control containing Fluormone™ PL Green/PR-LBD(GST) complex plus a competing compound, such as progesterone, may be included.

Fluormone™ PL Green/PR-LBD(GST) complex preparation: because of the hydrophobic nature of Fluormone™ PL Green, precautionary measures should be taken to minimize its adherence to plastic. We recommend preparing the Fluormone™ PL Green/PR-LBD(GST) complex in **glass**, or high-density polyethylene (HDPE) when glass is not available.

Handle PR-LBD(GST) gently: do not vortex. Mix by pipetting.

Solvent tolerances: Up to 0.5% ethanol, 1.25% methanol (in addition to the 0.45% methanol resulting from inclusion of 2 nM PL Green) or DMSO may be added to the assay without a reduction in dynamic range (•mP).

Note: The dynamic range will be decreased 10 to 15% when 1.25% ethanol, 5% methanol or 5% DMSO are included in the assay.

Standards: We recommend using 1 nM Low Polarization Solution and 1X High Polarization Solution from the Beacon® FP One-Step Standardization Kit (PanVera Part No. P2581) to determine if the instrument is measuring polarization accurately.

4.2 Competition Experiments

Design the fluorescence polarization competition experiments such that the starting polarization value will represent at least 50% of the maximum shift (*i.e.*, add enough PR-LBD(GST) to bind at least 50% of the Fluormone™ PL Green). The K_d of Fluormone™ PL Green with PR-LBD(GST) is approximately 10 nM. We recommend using 40 nM PR-LBD(GST) to achieve ~80% saturation with a final concentration of 2 nM Fluormone™ PL Green concentration. Fluormone™ PL Green concentrations greater than 2 nM may be required in instruments lacking wavelength-specific dichroic mirrors.

5.0 PROCEDURE

Remove PR-LBD(GST) from the -80°C freezer and Fluormone™ PL Green from the -20°C freezer and place on ice for at least 1 hour prior to use.

5.1 Prepare Test Reagents

1. Prepare serial dilutions of the test compounds in PR Screening Buffer, Green Assay directly in the multiwell plate. Prepare these dilutions in 50 μL volumes, so that the test compounds will be diluted two-fold when 50 μL PR-LBD(GST)/Fluormone™ PL Green complexes are added in the final reaction.

5.2 Prepare 4 nM Fluormone™ PL Green/PR-LBD(GST) Complex

1. Calculate the total volume of 2X Complex to be prepared: 50 μL of 2X Complex is needed per 100 μL assay.
2. Calculate the volume of each component required: PanVera recommends 2X Complexes containing 4 nM Fluormone™ PL Green and 80 nM PR-LBD(GST) in PR Screening Buffer, Green with 4 mM DTT. In the example shown below, 5 mL of 2X Complex is prepared using a 2000 nM PR-LBD(GST) stock. **Note that the concentration of PR-LBD(GST) is lot-dependent.** Five milliliters of 2X Complex will be sufficient for one 96-well plate. The empty rows of this table may be used as a work sheet for your experiments.

Total complex volume (μL)	[PR-LBD] (nM)	PR-LBD (μL) [80 nM @ 2X]	400 nM Fluormone™ PL Green (μL) [4 nM @ 2X]	1 M DTT (μL) [4 mM DTT @ 2X]	PR Screening Buffer, Green (μL)
5000	2000	200	50	20	4730

3. **Prepare the 2X Complex on ice:** Pipet the Green PR Screening Buffer into a **glass tube**. Add the required volume of DTT. Mix PR-LBD(GST) stock gently by pipetting; then add required volume to PR Screening Buffer, Green Assay/DTT mixture. Vortex 400 nM Fluormone™ PL Green. Pipette up and down to coat tip, then add needed volume into protein/buffer mixture. Mix well by pipetting.

5.3 Perform the Competition Assay

1. Add 50 μL of 2X Fluormone™ PL Green/PR-LBD(GST) Complex to the multiwell plate wells (already containing 50 μL of the test compound serial dilutions) and mix well by shaking on a plate shaker.
2. Make control wells. A “0% competition” negative control containing 1X Complex, without competitor is recommended. The polarization value for the negative control should be high. A “100% Competition” positive control containing 1X Complex with 10 μM Progesterone, may be used. The polarization value for the “100% Competition” positive control should be as low as the Fluormone™ PL Green without PR-LBD(GST); in other words, both should have polarization values equivalent to “free” Fluormone™ PL Green.
3. Incubate the plate in the dark at room temperature and read at 1-4 hours.

Note: After 4 hours at room temperature, polarization values will start to decrease. Plates may be stored at $+4^{\circ}\text{C}$ for up to 24 hours, then warmed to room temperature and read.

4. Measure polarization values using a plate reader equipped for measuring fluorescence polarization.

Note: All wells should be blanked against a control well containing 1X PR-LBD(GST) only. Alternatively, one can use PR Screening Buffer, Green Assay as the blank, since the inherent fluorescence of PR-LBD(GST) is very low.

6.0 RESULTS AND DISCUSSION

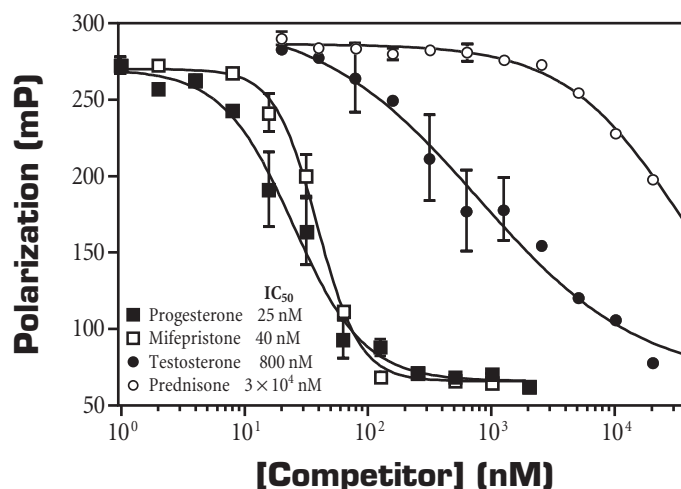
Below is an example of competition data generated on a 96-well plate. The concentration of test compound resulting in a half-maximum shift in polarization equals the IC_{50} of the compound.

This curve was plotted using the following equation:

$$Y = mP_{100\%} + (mP_{0\%} - mP_{100\%}) / (1 + 10^{((\text{Log}IC_{50} - X) \cdot \text{HillSlope})})$$

Where: Y = mP, X = Log [inhibitor], $mP_{100\%}$ = 100 % inhibition, and $mP_{0\%}$ = 0 % inhibition

Curve fitting was performed using Prism® software from GraphPad™ Software, Inc.



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

PanVera and CoreHTS are registered trademarks of Invitrogen Corporation.

Fluormone is a trademark of Invitrogen Corporation.

Prism is a registered trademark of GraphPad Software, Inc.

©2003, Invitrogen Corporation. All rights reserved. Reproduction with permission forbidden.

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 Invitrogen Corporation 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.